

Quality Control Products and Services



August 2025

This edition of the Quality Control Products and Services catalog features the full range of control families.

Recently added products to the **Immunoassay** family include:

NEW Liquichek and InteliQ Specialty Immunoassay Plus Controls

Specialty Immunoassay Plus is a four level, comprehensive, independent specialty immunoassay quality control designed to monitor assay precision of specialty analyte testing. Specialty Immunoassay Plus Control features high-demand, clinically relevant analytes such as Procalcitonin (PCT), IL-6, and Active Vitamin B₁₂. Available in multiple configurations to fit your laboratory's needs. Learn more:

- Liquichek Specialty Immunoassay Plus Control (page 62)
- InteliQ Specialty Immunoassay Plus Control (page 14)

NEW EQAS Specialty Immunoassay Program

The EQAS SPIA program is a liquid, independent, and confidential external assessment that consolidates 13 specialty immunoassay analytes, including procalcitonin, IL-6, active vitamin B₁₂, and fructosamine. Accredited to ISO/IEC 17043, this program offers labs a single shipment of 12 samples to test and report on monthly, with detailed performance reports provided. Learn more:

- EQAS Specialty Immunoassay Program (page 50)

Support for Bio-Rad controls and software products does not end after you receive your products

These support groups are available to help you with your questions and needs

By Phone within the U.S.

1-800-854-6737
Monday – Friday
5:00 AM to 4:00 PM
Pacific Time (PT)

By Fax

1-949-598-1531

Quality Controls Technical Support

When you purchase control products from Bio-Rad, you receive the backing of a full-service Technical Support Department. Bio-Rad Technical Service Representatives are trained to handle your inquiries in a prompt, knowledgeable and professional manner.

Unity™ and EQAS™ Program Support

The QC Program Department at Bio-Rad offers you expert technical support and information for all Unity and EQAS reports. Our QC Program staff answers inquiries about the Unity Interlaboratory Program, QCNet enrollments, as well as EQAS enrollments, submission of data and cycle dates.

Software and Connectivity Support

The Software and Connectivity Support Department is dedicated to helping your laboratory get up and running with the latest QC data management and connectivity solutions from Bio-Rad. With the help of our experts, we'll make sure your laboratory is able to capture, analyze and submit QC data successfully.

For Worldwide Customers

For service outside the United States, please refer to the inside back cover for the Bio-Rad location nearest you.

Online Resources

QCNet™

Learn more about the many QC tools and resources from Bio-Rad at www.QCNet.com

Unity™ Interlaboratory Reports

- Access your monthly Unity and InstantQC Reports in just a few quick clicks

Online Resources available on QCNet

- My eInserts
- QC Articles
- QC Webinars
- Training and Education
- Technical Documents
- Value Assignment Program
- Barcode Label Generator
- eInserts XML
- Linearity Web
- FAQs
- Link to YouTube channel

EQAS™ Resources on QCNet.com

Use the [EQAS page](#) for quick access to program essentials

Access program information, product inserts and user documentation through www.QCNet.com. This secure website allows registered participants access to training material, customer notifications and provides the ability to submit results with EQAS Online and view or print EQAS reports.

Access Bio-Rad product literature online

For more detailed product information, visit our [websites](#)

QCNet.com—Click on the Technical Documents page for access to Quality Control literature.

Bio-Rad.com—Click on the Literature Library link on the home page to search for a specific product or product family.



Powering Exceptional Labs

Bio-Rad harnesses independent QC to take your lab to the next level, so you can stand behind your results, minimize risk and costly errors, and improve your workflow to power exceptional performance for your laboratory.

We offer high-quality materials, insightful informatics systems and renowned customer support, including staff training, lot reservation and technical expertise. We go beyond just monitoring precision to help you improve performance in your laboratory. We demand and invest in quality—because patients deserve nothing less.

Table of Contents

Quality Control Products & Services

- QC Data Management Solutions 4
- IntelliQ Controls 10
- Infectious Disease: Molecular Quality Controls and Standards 15
- Blood Gas Controls 34
- Cardiac Assessment Controls 35
- Chemistry Controls & Serum Indices 37
- Coagulation Controls 44
- Diabetes/Hemoglobin Controls 45
- External Quality Assurance Services (EQAS) 47
- Hematology Controls 53
- Immunoassay Controls 59
- Immunology/Protein Controls 65
- Infectious Disease: Serology Controls 69
- Therapeutic Drug Monitoring Controls 77
- Toxicology: Drugs-of-Abuse Controls 79
- Toxicology: Specialty Controls 85
- Urinalysis Controls 87
- Accessories 88

Analyte Charts

- Control Analyte Index 89
- Blood Gas Controls Analyte Index 111
- Cardiac Assessment Controls Analyte Index 111
- Hematology Controls Analyte Index 112
- Infectious Disease: Molecular Quality Controls and Standards Analyte Index 113
- Infectious Disease: Serology Controls Analyte Index 117
- IntelliQ Analyte Index 119
- QC Products in Compact Vials Analyte Index 123
- External Quality Assurance Services (EQAS) Analyte Index 125

Index

- Alphabetical Index 130
- Catalog Number Index 133

Additional Information

- How to Place an Order 139
- Legal Notices 140
- Worldwide Contact Information Inside Back Cover

Unity™ QC Software Solutions

Your strategy for
smarter QC



With the world's largest QC peer group, Unity is the industry's leading software solution specifically designed to optimize laboratory QC practices and streamline workflows. By providing QC data in the form of real-time reports, your system will be better able to detect and correct problems before they impact patient results. Unity's robust desktop- or cloud-based interface and data management tools help improve your lab's analytical performance and attain a higher level of confidence in your test results. For a fully integrated QC system—from quality control to data analysis—choose Bio-Rad for independent QC.

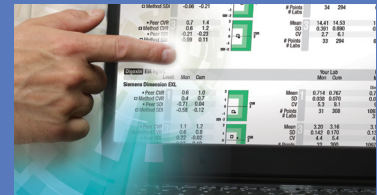
Powerful, real-time data management



Provides the largest community
of QC users



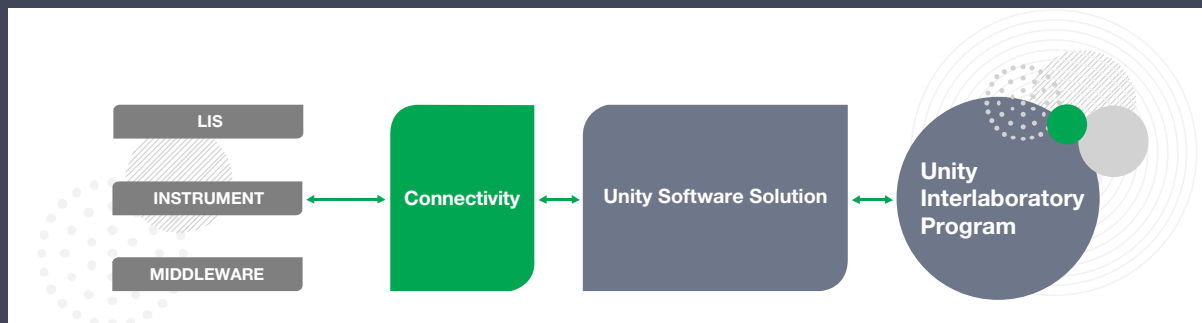
Streamlined workflow and
optimized performance



Increased confidence and
reliable patient results

Data Transfer and Connectivity Solutions

Bio-Rad offers several interoperability solutions to capture and process QC data directly from your laboratory devices or information systems. Bio-Rad connectivity solutions allow your lab to easily and automatically collate data, eliminate manual tasks, and efficiently troubleshoot problems. These solutions seamlessly connect a wide range of instruments and platforms to Bio-Rad's Unity software, helping you unlock the power of quality in your lab.



Join over 38,000 of your peers who already use the Unity Interlaboratory Program, the most robust program for clinical labs

Bio-Rad Quality Controls are part of the Unity Program with industry-leading features for benchmarking laboratory performance. Whether a lab produces five hundred or five million results a year, the lab director bears ultimate responsibility for the accuracy and timeliness of each test result.

A patient risk approach to QC means planning your QC based on the type or method of the test and its impact on patient outcomes when a failure is detected. As the world's leader in clinical laboratory QC, we've established the foundation to help you make the transition to patient-focused QC.

- Benefit from peer data generated from more than 59,000 instruments worldwide and 68 million monthly QC results
- Utilize real-time access to peer group comparisons to help with troubleshooting
- Consolidate your entire laboratory into one QC system

- Facilitate run-time decisions with optimized Westgard rules
- Facilitate compliance with simple tools to review rule violations and document corrective actions
- Define your own quality requirements and track key quality performance indicators
- Automate QC data from instruments, middleware and/or LIS systems with Unity connectivity

This sophisticated software includes best-in-class peer tools to support releasing results, benchmarking laboratory performance, and designing your QC system. Unity enables you to work smarter, making it possible to achieve optimum lab practice for your QC workflow. Reducing manual QC data management tasks with software automation frees your staff to focus on more critical tasks. Unity ultimately helps your team deliver objectives more efficiently and with greater ease.

Unity™ Interlaboratory Reports

There has never been a better time to be a member of the world's largest QC peer group program

Unity January 2025 • Lab 12345

Associated Regional Laboratory
123 Main Street
Anytown, NY 12345-6789
Attention: Lab Supervisor

Monthly Evaluation
Immunoassay Plus • Lot 12345 • Exp 01-Jan-2027

Please review your QC reports for January 2025.

! The tests listed below may require investigation or review !

Level	Lab	Peer	Method
1	2.15	2.80	3.00
	SD 0.14	0.170	0.253
	CV 6.5	6.0	8.4
	# Points 31	6797	20391
	# Labs 223		670
2	5.54	5.98	6.31
	SD 0.171	0.209	0.469
	CV 3.1	3.5	7.4
	# Points 32	6588	19764
	# Labs 220		659
3	7.94	7.87	8.13
	SD 0.215	0.273	0.575
	CV 2.7	3.5	7.1
	# Points 33	6797	20391
	# Labs 223		670

Data Exclusion: Lab Mean = 2.15
Acceptable values are 2.1941 to 3.5306
This data was not used as part of the Unity worldwide statistical database.

Warning: Acceptable values are above -2.0, below 2.0

This level is within established parameters

Unity January 2025 • Lab 12345

Associated Regional Laboratory
123 Main Street
Anytown, NY 12345-6789
Attention: Lab Supervisor

Laboratory Comparison Report
Immunoassay Plus • Lot 12345 • Exp 01-Jan-2027

Level	Mon	Cum	Year Lab Mon Cum	Peer Group Mon Cum	Method Group Mon Cum	
1	0.9	1.8	2.99	2.99	2.97	3.01
	SD 0.8	1.7	0.170	0.372	0.196	0.209
	CV 0.10	-0.12	5.7	12.5	6.6	7.0
	# Points 0.45	0.4	33	316	6794	53336
	# Labs 226				20382	160K
					677	925
2	0.9	0.9	8.75	8.66	8.73	8.88
	SD 0.6	0.8	0.314	0.366	0.369	0.402
	CV 0.05	-0.55	3.6	4.2	4.2	4.5
	# Points -0.06	-0.21	34	294	6412	50441
	# Labs 222				19236	151K
					666	909
3	0.7	1.4	14.41	14.53	14.54	14.68
	SD 0.6	1.2	0.391	0.890	0.588	0.661
	CV 0.21	-0.23	2.7	6.1	4.0	4.5
	# Points -0.09	0.11	33	294	6794	53336
	# Labs 226				20382	160K
					677	925

Unity Interlaboratory Reports provide visualizations to indicate your laboratory performance. With monthly evaluation reports, receive notifications of data exclusions and identify potential issues in your QC performance. Additional reports allow you to view your trending data over the past 12 months and vital statistics are shown in a comprehensive overview compared to peer and method consensus groups.

Unity Interlaboratory Program reports feature:

- Logical arrangement of data provides a quick and intuitive read
- Data consolidation and imagery simplifies interpretation of lab data

List of Available Reports

- Monthly Evaluation
- Laboratory Performance Overview
- Laboratory Comparison Report
- Laboratory Histogram
- Bias & Imprecision Histogram
- Worldwide Report
- Statistical Profile Report
- Affiliated Laboratory Comparison Report
- Affiliated Laboratory Comparison Report: Abbreviated Summary
- Affiliated Data Exception Report
- Urinalysis Reports
- Qualitative Worldwide Report
- Blood Typing/Serum Indices Report
- Instant QC Report

Unity™ Solutions

Bio-Rad Laboratories offers several options for participation in the Unity Interlaboratory Program. Refer to the product comparison grid below to help determine which solution is the most appropriate for your laboratory.

	Basic User Options			Advanced User Options	
	UnityWeb	Unity Real Time LT	Unity Next Peer QC	Unity Real Time online	Unity Real Time
Type of Solution					
Web Service (Web-based)	●		●	●	
Desktop Software		●			●
Unity Interlaboratory Reports					
Monthly Reports	●	●	●	●	●
Instant and real-time reports	accessible separately	accessible separately	●	accessible separately	accessible separately
Basic Intralaboratory Charts & Reports					
Westgard Rules	●	●	●	●	●
Various Charts and Reports	●	●	●	●	●
Additional Features					
Westgard Advisor (Subscription sold separately)	●			●	●
Bench Review with Data Review Report	●	●		●	●
Supervisor Data Review with Data Review Report				●	●
Analytical Goals				●	●
Dynamic Data Set Comparisons				●	●
Database Platform					
Bio-Rad Hosted	●		●	●	
Installed on Local Laboratory PC		●			●
Connectivity (Sold separately)					
WebConnect	●			●	
UnityConnect	●	●		●	●
Unity Next File Upload			●		

Want to know more? Discover the power of Unity Solutions at www.qcnet.com/qc-data-management



Customers agree:

“The documentation of Bio-Rad Controls using the Unity QC Program has been well accepted by CAP inspectors. Worry-free QC.”

— Laboratory Director,
Health Center

“To compare our data with a substantial peer group ensures that we are turning out reliable patient data.”

— Laboratory Supervisor,
School of Medicine

“The reports are clean and easy to read and understand. For each test, Unity Reports for our lab show a summary of our data, our instrument group, and our method group—reports are simple.”

— Clinical Chemist,
Hospital Laboratory

Unity Real Time™ LT

Entry-Level Desktop QC Data Management Solution

- Basic QC rules, charts and reports
- Configurable Levey-Jennings charts
- Upgrade easily to Unity Real Time for more advanced tools and features
- Upload QC data points from an LIS, middleware or instrument (optional)

Ordering Information

Cat #	Description
805-1	Unity Real Time LT Annual subscription
825i	Unity Remote Installation & Training Single use
858i	Additional Software Training Single use

UnityWeb™

Entry-Level Online QC Data Management Solution

- No software to install and maintain
- Use basic QC rules, charts and reports
- Upgrade easily to Unity Real Time online for more advanced tools and features
- Upload QC data points from an LIS, middleware or instrument (optional)

Ordering Information

Cat #	Description
870-1	UnityWeb Annual subscription
825i	Unity Remote Installation & Training Single use
858i	Additional Software Training Single use

Unity Next Peer QC

An entry-level online QC data management solution with instant access to peer comparison and the ability to perform basic QC processes

- Latest generation of Unity with user-friendly interface
- No software to install and maintain
- Use basic QC rules, Levey-Jennings charts and reports
- Instantly access peer reports for peer comparison and QC troubleshooting
- Upload QC data points from LIS, middleware, or instrument
- Submit QC data and download reports in one location

Ordering Information

Cat #	Description
11000517	Unity Next Peer QC (5 users) Annual subscription
11000506	Unity Next Onboarding Single use

WebConnect

A web-based connectivity solution that allows laboratories to easily upload QC data from LIS systems, middleware and/or instruments directly into Unity Real Time online or UnityWeb

- Eliminate manual keying of data
- Standard LIS QC reports can be used
- No software to install

Ordering Information

Cat #	Description
870-W1	WebConnect 2.0 Annual subscription
825i	Unity Remote Installation & Training Single use
858i	Additional Software Training Single use

Unity Next File Upload

A data upload solution allowing file uploads of QC data, specifically for Unity Next software

- Upload QC data from LIS, middleware, and/or instruments to Unity Next
- No software to install

Ordering Information

Cat #	Description
11001897	Unity Next File Upload Annual subscription

Unity Real Time™

Expert QC Data Management Solution for Desktop Users

- Facilitate compliance under CLIA and ISO 15189
- Improve real-time bench and supervisor QC data review
- Implement best QC rules when used with Westgard Advisor
- Run validation with comprehensive audit trails
- Advanced charts and reports for data analysis
- Reduce non-essential retests with Analytical Goal options
- Upload QC data points from an LIS, middleware or instrument (optional)
- RiLiBÄK module available to comply with German regulations
- GOST module available to comply with Russian regulations

Ordering Information

Cat #	Description
804-1	Unity Real Time Annual subscription
825i	Unity Remote Installation & Training. Single use
858i	Additional Software Training Single use

Unity Real Time™ online

Expert QC Data Management Solution for Online Users

- Facilitate compliance under CLIA and ISO 15189
- Improve real-time bench and supervisor QC data review
- Implement best QC rules when used with Westgard Advisor
- Run validation with comprehensive audit trails
- Advanced charts and reports for data analysis
- Reduce non-essential retests with Analytical Goal options
- Upload QC data points from an LIS, middleware or instrument (optional)

Ordering Information

Cat #	Description
804-W1	Unity Real Time online Annual subscription
825i	Unity Remote Installation & Training. Single use
858i	Additional Software Training Single use

UnityConnect

A connectivity solution that allows QC data from LIS systems, middleware and/or instruments to be quickly and easily imported into Unity software, enabling interlaboratory program comparison

- Eliminate manual keying of QC data
- Standard LIS QC reports can be used
- Automatic data import and real-time connection
- Capture and analyze QC data economically from laboratory instruments not connected to the LIS
- Easy-to-install optional interface hardware
- Choose between serial and ethernet connection

Ordering Information

Cat #	Description
11000009	UnityConnect Software Annual subscription
858i	Additional Software Training Single use

Hardware solutions available for serial and network devices. Contact your Bio-Rad sales representative for recommendation.

QC Lot Viewer

A cloud-based application that allows visibility into QC standing orders to help simplify order planning

- Standing QC order reports
- Shipment metric reports
- Multiple users
- Free and unlimited access

Ordering Information

Available to eligible customers. Contact sales for more information.

Available in the US only

Bio-Rad Mission: Control™

Take the guesswork out of your QC strategy

- Determine your laboratory’s Risk Management Index (RMI)
- Evaluate your test method performance based on your QC results
- Quantify the likelihood and severity of reporting incorrect patient results
- Assess how much you can lower your risk, based on instrument performance
- Create a QC Plan by confidently selecting rules and frequency that fit your risk tolerance
- Minimize your risk of reporting unreliable patient results

Ordering Information

Cat #	Description
11000002	Mission: Control. Annual Subscription
11000005	Mission: Control. Subscription and Service Bundle

Westgard Advisor

Automatic QC Rules Selection Engine

- Recommend and automatically apply best QC rules with patented technology
- Easy step-by-step automatic rule selection capabilities
- Reduce false rejections and desensitization to false error flags
- Save time and money by reducing unnecessary repeats and troubleshooting
- Improve laboratory test quality with optimally selected QC rules
- Available as an optional module with Unity QC data management solutions

Ordering Information

Cat #	Description
811-1	Westgard Advisor Annual subscription
825i	Unity Remote Installation & Training. Single use
858i	Additional Software Training Single use

Unity™ Alert

Continuously Monitor QC Status

- Notifications of missing QC runs can help to ensure proper monitoring of instrument performance
- Notifications of new violations can be useful for troubleshooting and monitoring problem analytes
- Alert rules are easy to set up and customize for specific instruments or tests
- Receive emails remotely to stay informed of QC issues in your laboratory
- Color-coded displays help draw attention to the most urgent QC issues
- Notifications are provided even when Unity Real Time is not in use
- Available as an optional module with Unity Real Time

Ordering Information

Cat #	Description
806-1	Unity Alert Annual subscription
825i	Unity Remote Installation & Training. Single use
858i	Additional Software Training Single use

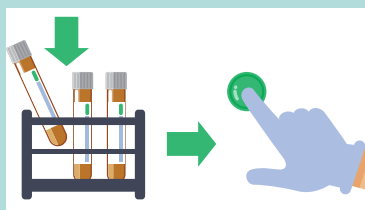
InteliQ™ Load-and-Go Quality Controls

Ready-to-use,
pre-barcoded tubes
help automate
your QC workflow

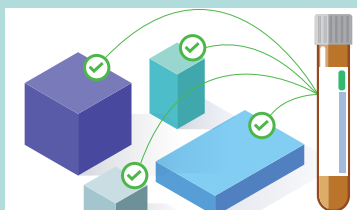


Transform your QC workflow with load-and-go efficiency and robust data management. Bio-Rad's InteliQ quality controls are the only barcoded, independent quality controls that are compatible with most major chemistry and immunoassay platforms. Because the pre-barcoded tubes are ready to load directly onto your instrument, InteliQ quality controls help automate your QC workflow—reducing turnaround time, limiting manual errors, and increasing your lab's overall efficiency. Combined with Unity data management, InteliQ provides a smarter way to manage your QC.

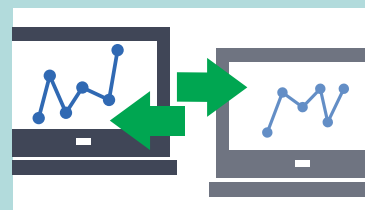
Transform your QC Workflow



Streamlined, automated workflow



Universal compatibility



Confidence in patient results

Benefits of InteliQ Quality Controls



Improve Turnaround Time

Pre-barcoded, ready-to-use quality controls reduce hands-on time and manual errors, providing faster turnaround from sample to result.



Enhance Staff Satisfaction

Automated QC set-up increases laboratory efficiency and promotes staff satisfaction.



Reduce Laboratory Waste

Eliminating pour-off steps reduces waste and saves money while open-vial storage options allow for repeated use of QC tubes.

InteliQ Assayed Multiquel Control



- Provides comprehensive range of 80 analytes to maximize testing efficiency for automated platforms
- Liquid, human serum based
- 3 year shelf life at -20°C to -70°C
- 7 day stability at 2-8°C for most analytes (refrigerated, off-board, all instruments)*
- 14 day stability at 2-8°C for most analytes (on-board Siemens Atellica Solution instrument)

Ordering Information

Cat #	Description
12008256	Level 1 12 x 3 mL
12008257	Level 2 12 x 3 mL
12008258	Level 3 12 x 3 mL
12008259	Trilevel MiniPak (1 of each level) 3 x 3 mL

Analytes

α-1-Antitrypsin	Bilirubin (Total)	Gentamicin	Potassium
α-1-Globulin	Calcium	Glucose	Prealbumin
α-2-Globulin	Calcium (Ionized)	Haptoglobin	Protein Serum (Total)
β-Globulin	Carbamazepine	Immunoglobulin A (IgA)	Salicylate
γ-Globulin	Carbon Dioxide (CO ₂)	Immunoglobulin G (IgG)	Sodium
γ-Glutamyltransferase (GGT)	Ceruloplasmin	Immunoglobulin M (IgM)	T3 (Total)
Acetaminophen	Chloride	Iron	T3 Uptake/T Uptake
Acid Phosphatase (Total)	Cholesterol (HDL)	Iron (TIBC)	T4 (Free)
Albumin	Cholesterol (LDL)	Iron (UIBC)	T4 (Total)
Albumin-to-Globulin Ratio	Cholesterol (Total)	Lactate (Lactic Acid)	Theophylline
Alkaline Phosphatase	Cholinesterase	Lactate Dehydrogenase (LDH)	Thyroid Stimulating Hormone (TSH)
ALT/SGPT	Complement C3	LAP Arylamidase	Tobramycin
Amikacin	Complement C4	Lipase	Transferrin
Amylase	Copper	Lithium	Triglycerides
Amylase (Pancreatic)	Cortisol	Magnesium	Urea
Apolipoprotein A-I	Creatine Kinase (CK)	Osmolality	Urea Nitrogen (BUN)
Apolipoprotein B	Creatinine	Phenobarbital	Uric Acid
AST/SGOT	Digoxin	Phenytoin	Valproic Acid
Bilirubin (Direct)	Ethanol	Phospholipids	Vitamin B ₁₂
Bilirubin (Neonatal)	Ferritin	Phosphorus	Zinc

Cardiac Advance

InteliQ Cardiac Advance Control



- Consolidated multi-analyte cardiac control optimized for high sensitivity Troponin and barcoded for load-and-go efficiency.
- Liquid, human serum based
- Optimized for high sensitivity troponin
- Level 1 Troponin values near limit of detection for most instruments.
- 28 month shelf life at -20°C to -70°C
- 10 day stability at 2-8°C for most analytes (refrigerated, off-board, all instruments)*
- 10 day stability at 2-8°C for most analytes (on-board Siemens Atellica Solution instrument)

Ordering Information

Cat #	Description
12016484	Level 1 6 x 3 mL
12016485	Level 2 6 x 3 mL
12016486	Level 3 6 x 3 mL
12016487	Level 4 6 x 3 mL

For Liquechek Cardiac Advance Control, see page 35

Analytes

BNP	Digitoxin	Myoglobin	Troponin T ¹
CK-MB Isoenzyme	Galectin-3	NT-proBNP	¹ Compatible with high sensitivity methods.
CRP ¹	Homocysteine	Troponin I ¹	

Compatible with Siemens and Roche platforms. Abbott users should use InteliQ Cardiac Markers Plus Control LT. Homocysteine not compatible with Siemens platforms. Siemens users should use InteliQ Homocysteine Control.

*Please refer to the package insert for platform-specific stability and performance claims. Refer to myInSerts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

InteliQ Cardiac Markers Plus Control LT



- Multi-analyte control designed to monitor precision of cardiac assessment assays
- Liquid, human serum based
- 3 year shelf life at -20°C to -70°C
- 10 day stability at 2-8°C with exceptions (refrigerated, off-board, all instruments)*
- 20 day stability at 2-8°C with exceptions (on-board Siemens Atellica Solution instrument)

Ordering Information

Cat #	Description	
12009955	Level 1	.6 x 3 mL
12009958	Level 1A	.6 x 3 mL
12009959	Level 1B	.6 x 3 mL
12009960	Level 1C	.6 x 3 mL
12009956	Level 2	.6 x 3 mL
12009957	Level 3	.6 x 3 mL
12009961	Level 1 MiniPak	.1 x 3 mL
12009964	Level 1A MiniPak	.1 x 3 mL
12009965	Level 1B MiniPak	.1 x 3 mL
12009966	Level 1C MiniPak	.1 x 3 mL
12009962	Level 2 MiniPak	.1 x 3 mL
12009963	Level 3 MiniPak	.1 x 3 mL

For Liquichek Cardiac Markers Plus Control LT, see page 35

Analytes

BNP
Creatine Kinase (CK)
CK-MB Isoenzyme
CRP¹
Digitoxin

Myoglobin
NT-proBNP
Troponin I¹
Troponin T¹

¹Compatible with high sensitivity methods.

InteliQ Diabetes Control



- Trilevel HbA1c control designed to monitor the precision of test procedures associated with diabetes monitoring for Hemoglobin A1C
- Liquid, human whole blood based
- 3 year shelf life at -20°C to -70°C
- 10 day stability at 2-8°C (refrigerated, off-board, all instruments)*
- 45 day stability at 2-8°C (on-board Siemens Atellica Solution instrument)

Ordering Information

Cat #	Description	
12008309	Level 1	.6 x 2 mL
12008310	Level 2	.6 x 2 mL
12008311	Level 3	.6 x 2 mL
12008312	Trilevel MiniPak (1 of each level)	.3 x 2 mL

Analyte

Hemoglobin A1C

InteliQ Ethanol/Ammonia Control



- Normal, elevated and toxic concentrations of Ethanol and Ammonia to monitor precision across the clinical range
- 2 year shelf life at 2-8°C
- Ammonia:
 - 4 day stability at 2-8°C (refrigerated, off-board, all instruments)*
 - 13 day stability at 2-8°C (on-board Siemens Atellica Solution instrument)
- Ethanol:
 - 3 day stability at 2-8°C (refrigerated, off-board, all instruments)*
 - 3 day stability at 2-8°C (on-board Siemens Atellica Solution instrument)

Ordering Information

Cat #	Description	
12008299	Level 1	.6 x 2 mL
12008300	Level 2	.6 x 2 mL
12008301	Level 3	.6 x 2 mL
12008302	Trilevel MiniPak (1 of each level)	.3 x 2 mL

Analytes

Ammonia
Ethanol

*Please refer to the package insert for platform-specific stability and performance claims. Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

InteliQ Homocysteine Control



- Monitors precision of Homocysteine assays
- Liquid, human serum based
- 30 day stability at 2-8°C (on-board Siemens Atellica Solution instrument)
- 5 day open-vial stability at 2-8°C (refrigerated, off-board, all instruments)*

Ordering Information

Cat #	Description
12004309	Level 16 x 1 mL
12004310	Level 26 x 1 mL
12004311	Level 36 x 1 mL
12004312	Trilevel MiniPak (1 of each level)3 x 1 mL

Analyte

Homocysteine

InteliQ Immunoassay Plus Control



- Offers a wide range of popular immunoassay analytes
- Provides an efficient solution for laboratories that focus on routine testing
- Liquid human serum based
- 28 month shelf life at -20°C to -70°C
- 10 day stability at 2-8°C for most analytes (refrigerated, off-board, all instruments)*
- 14 day stability at 2-8°C for most analytes (on-board Siemens Atellica Solution instrument)

Ordering Information

Cat #	Description
12009948	Level 112 x 4 mL
12009949	Level 212 x 4 mL
12009950	Level 312 x 4 mL
12009951	Trilevel MiniPak (1 of each level)3 x 4 mL

Analytes

Acetaminophen	Ferritin	Luteinizing Hormone (LH)	T3 (Total)
AFP	Folate	N-Acetyl Procainamide (NAPA)	T3 Uptake/T Uptake
Amikacin	FSH	Phenobarbital	T4 (Free)
Androstenedione	Gentamicin	Phenytoin	T4 (Total)
Caffeine	hCG	Primidone	Testosterone
Carbamazepine	hGH	Procainamide	Theophylline
CEA	Immunoglobulin A (IgA)	Progesterone	Thyroid Stimulating Hormone (TSH)
CK-MB Isoenzyme	Immunoglobulin E (IgE)	Prolactin	Thyroxine Binding Globulin (TBG)
Cortisol	Immunoglobulin G (IgG)	PSA (Free)	Tobramycin
Dehydroepiandrosterone-Sulfate (DHEA-Sulfate)	Immunoglobulin M (IgM)	PSA (Total)	Tricyclic Antidepressants (TCA)
Digoxin	Insulin	Quinidine	Valproic Acid
Estradiol	Iron	Salicylate	Vancomycin
Estril (Free)	Lidocaine	SHBG	Vitamin B ₁₂
	Lithium	T3 (Free)	

InteliQ Immunology Control



- Comprehensive menu of serum proteins and analytes, including Rheumatoid Factor
- Liquid human serum based
- 2 year shelf life at -20°C to -70°C
- 10 day stability at 2-8°C for most analytes (refrigerated, off-board, all instruments)*
- 30 day stability at 2-8°C for most analytes (on-board Siemens Atellica Solution instrument)

Ordering Information

Cat #	Description
12009941	Level 16 x 3 mL
12009942	Level 26 x 3 mL
12009943	Level 36 x 3 mL
12009944	Trilevel MiniPak (1 of each level)3 x 3 mL

Analytes

α-1-Acid Glycoprotein (AAG)	Antithrombin III	Ferritin	Lipoprotein (a)
α-1-Antitrypsin (AAT)	Apolipoprotein A-1	Haptoglobin	Prealbumin
α-2-Macroglobulin	Apolipoprotein B	Immunoglobulin A (IgA)	Properdin Factor B
β-2-Microglobulin	C1 Inhibitor	Immunoglobulin E (IgE)	Protein Serum (Total)
ADNase B	Ceruloplasmin	Immunoglobulin G (IgG)	Retinol Binding Protein (RBP)
Albumin	Complement C3	Immunoglobulin G Subclasses 1-4	Rheumatoid Factor
Antistreptolysin O (ASO)	Complement C4	Immunoglobulin M (IgM)	Transferrin
Anti-Cyclic Citrullinated Peptide (Anti-CCP)	C-Reactive Protein (CRP)	Kappa Light Chain	
	Cystatin C ¹	Lambda Light Chain	¹ Endogenous levels only.

*Please refer to the package insert for platform-specific stability and performance claims. Refer to myInSerts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

InteliQ Specialty Immunoassay Control



- Multi-analyte control for monitoring the precision of specialty immunoassay testing
- Liquid, human serum based
- 2 year shelf life at -20°C to -70°C
- 10 day stability at 2-8°C for most analytes (refrigerated, off-board, all instruments)*
- 30 day stability at 2-8°C for most analytes (on-board Siemens Atellica Solution instrument)

Ordering Information

Cat #	Description	
12008276	Level LTA	6 x 4 mL
12008277	Level 1	.6 x 4 mL
12008278	Level 2	.6 x 4 mL
12008279	Level 3	.6 x 4 mL
12008280	Four Level MiniPak (1 of each level)	.4 x 4 mL

Analytes

25-Hydroxy Vitamin D	Anti-Thyroperoxidase (Anti-TPO)	Erythropoietin (EPO)	Osteocalcin
Anti-Thyroglobulin (Anti-Tg)	C-Peptide	Insulin Growth Factor I (IGF-I)/Sorratomedin C	PTH (Intact)

NEW | InteliQ Specialty Immunoassay Plus Control



- Multi-analyte control for monitoring the precision of specialty immunoassay testing
- Liquid, human serum based
- 3 year shelf life at -20°C to -70°C
- 10 day stability at 2-8°C for most analytes (refrigerated, off-board, all instruments)*
- 30 day stability at 2-8°C for most analytes (on-board Siemens Atellica Solution instrument)

Ordering Information

Cat #	Description	
12015705	Level Low	6 x 3 mL
12015706	Level 1	.6 x 3 mL
12015707	Level 2	.6 x 3 mL
12015708	Level 3	.6 x 3 mL
12019558	Level Low MiniPak	.1 x 3 mL
12015709	Trilevel MiniPak	.3 x 3 mL

For Liquichek Specialty Immunoassay Plus Control, see page 62

Analytes

25-Hydroxy Vitamin D	C-Peptide	Interleukin-6 (IL-6)	Thyroid Stimulating Immunoglobulin (TSI)
Active Vitamin B ₁₂	Erythropoietin (EPO)	PTH (Intact)	Anti-TSH Receptor (TRAb)
Anti-Thyroglobulin (Anti-Tg)	Fructosamine	Osteocalcin	
Anti-Thyroperoxidase (Anti-TPO)	Insulin Growth Factor I (IGF-I)	Procalcitonin (PCT)	

InteliQ Tumor Marker Control



- Comprehensive, assayed, multi-analyte control
- Monitors precision of immunoassay-based oncology testing procedures
- Liquid, human serum based
- 28 month shelf life at -20°C to -70°C
- 10 day stability at 2-8°C for most analytes (refrigerated, off-board, all instruments)*
- 30 day stability at 2-8°C for most analytes (on-board Siemens Atellica Solution instrument)

Ordering Information

Cat #	Description	
12008289	Level 1	.6 x 2 mL
12008290	Level 2	.6 x 2 mL
12008291	Level 3	.6 x 2 mL
12008292	Trilevel MiniPak (1 of each level)	.3 x 2 mL

Analytes

α-Fetoprotein (AFP)	CA 27.29	Cyfra 21-1**	HER-2/neu	Prolactin	Thyroglobulin (Tg)
β-2-Microglobulin	CA 72-4**	Ferritin	Insulin Growth Factor I (IGF-I)	PSA (Free)	**International Use Only
CA 15-3	CA 125	hCG	PAP	PSA (Total)	
CA 19-9	Carcinoembryonic Antigen (CEA)	HE4	ProGRP**	SCC**	

InteliQ Urine Chemistry Control



- Liquid, bilevel human urine based control
- Ideal for monitoring urine chemistry test procedures
- Offers 19 analytes
- 2 year shelf life at 2-8°C
- 10 day stability at 2-8°C for most analytes (refrigerated, off-board, all instruments)*
- 30 day stability at 2-8°C for most analytes (on-board Siemens Atellica Solution instrument)

Ordering Information

Cat #	Description	
12009995	Level 1	12 x 3 mL
12009996	Level 2	12 x 3 mL
12009997	Bilevel Minipak (1 of each level)	.2 x 3 mL

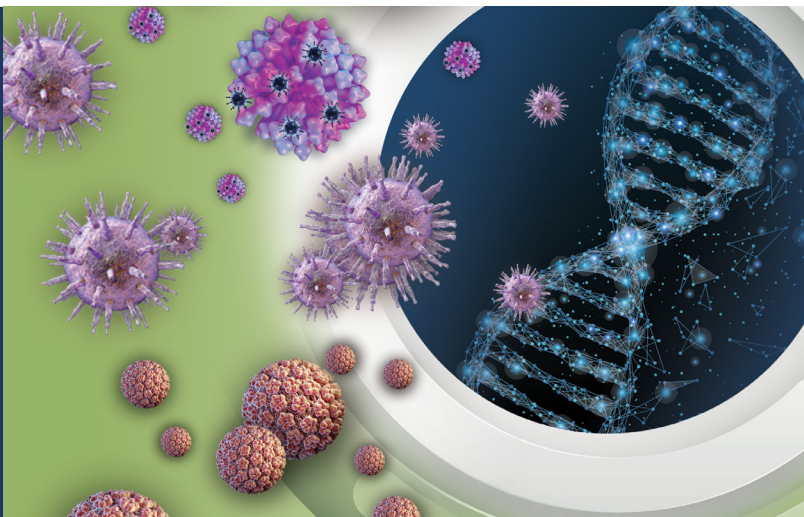
Analytes

Amylase	Creatinine	Osmolality	Pregnancy/hCG (Qualitative)	Urea
Calcium	Glucose	pH	Protein (Total)	Urea Nitrogen (BUN)
Chloride	Magnesium	Phosphorus	Sodium	Uric Acid
Cortisol	Microalbumin	Potassium	Specific Gravity	

*Please refer to the package insert for platform-specific stability and performance claims. Refer to myInserpts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Molecular Quality Controls and Panels

Setting a new standard in lab performance. Together.



With our expertise designing molecular quality controls and panels and the power of data management, we are helping you to advance the field of molecular diagnostics by driving quality and care.



Comprehensive portfolio of molecular diagnostic controls and panels

The Exact Diagnostics line of molecular controls and panels provide high-quality QC solutions for key diagnostic areas. These products feature unsurpassed quality, characterization, and consistency for which Bio-Rad is known. Our growing portfolio includes innovation in blood screening, syndromic panels and other molecular applications. Our key diagnostic areas include:

Transplantation

Virology

Sexually Transmitted Infections

Microbiology

Hospital Associated Infections

Respiratory

Tick-borne Diseases

Exact Diagnostics Quality Controls and Panels for Infectious Diseases



Patient-like Commutability
with simulation of patient specimens through extractable & inactivated whole organisms in high-quality clinically-relevant matrices



Standardization of Results
across laboratories through calibration to WHO International Standards (where applicable)



Independent Monitoring
of assay performance through long shelf-lives overlapping multiple assay reagent lots & assay-independent manufacturing



Consistency
through lot-to-lot reproducibility & concentration trueness using absolute quantification by droplet digital PCR (ddPCR)



Unity QC Data Management Ecosystem
for automated real-time result monitoring, QC rules customization & peer-grouping comparison

Ensuring Reliable Patient Results

Transplantation

Exact Diagnostics Adenovirus



- Verification Panel offers a wide dynamic range
- Manufacturing targets established using ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- True EDTA plasma matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
ADVL101*	Low Run Control	5 x 1.0 mL	IVD
ADVH102*	High Run Control	5 x 1.0 mL	IVD
Panel			
ADVP100	Verification Panel	5 x 1.0 mL	IVD

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analyte

Inactivated Adenovirus

Complementary product:

See page 26 for Negative Run Control (Cat# NEG000/NEG050)

See Page 26 for Bulk Urine Negative Run Control (Cat# 12019386)

Catalog #	Panel Member 1	Panel Member 2	Panel Member 3	Panel Member 4	Panel Member 5
ADVP100	200 cp/mL	2,000 cp/mL	20,000 cp/mL	200,000 cp/mL	2,000,000 cp/mL

Exact Diagnostics BKV



- Verification Panel offers a wide dynamic range
- Manufacturing targets established using ddPCR
- Calibrated against the 1st WHO International Standard for BK Virus
- High quality matrix to ensure lot-to-lot reproducibility
- True EDTA plasma matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
BKVL101*	Low Run Control	5 x 1.0 mL	IVD
BKVH102*	High Run Control	5 x 1.0 mL	IVD
Panels			
BKVP100	Verification Panel	6 x 1.0 mL	IVD
BKVP200	Verification Panel6 x 1.4 mL	IVD
BKVL201	1E3 IU/mL Precision9 x 1.4 mL	IVD
BKVH202	5E5 IU/mL Precision9 x 1.4 mL	IVD
BKVL0D	BKV LOD	10 x 1.0 mL	IVD
BKVUL0D	BKV Urine LOD	10 x 1.4 mL	IVD
Verification Bundle			
BKVB201	Verification Bundle . .	BKVP200 (4 kits), BKVL201 (2 kits), BKVH202 (2 kits), BKVL0D (2 kits), BKVUL0D (2 kits)	

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analyte

Inactivated BKV

Complementary product:

See page 26 for Negative Run Control (Cat# NEG000/NEG050)

See Page 26 for Bulk Urine Negative Run Control (Cat# 12019386)

Catalog #	Panel Member 1	Panel Member 2	Panel Member 3	Panel Member 4	Panel Member 5	Panel Member 6
BKVP100	200 IU/mL	2,000 IU/mL	20,000 IU/mL	200,000 IU/mL	2000,000 IU/mL	20,000,000 IU/mL
BKVP200	100 IU/mL	2,000 IU/mL	20,000 IU/mL	200,000 IU/mL	2000,000 IU/mL	20,000,000 IU/mL

Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Transplantation

Exact Diagnostics Parvovirus B19



- Verification Panel offers a wide dynamic range
- Manufacturing targets established using ddPCR
- Calibrated against the 3rd WHO International Standard for Parvovirus B19 (12/208)
- High quality matrix to ensure lot-to-lot reproducibility
- True EDTA plasma matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
B19L101*	Low Run Control	5 x 1.0 mL	IVD
B19H102*	High Run Control	5 x 1.0 mL	IVD
Panel			
B19P100	Verification Panel	5 x 1.0 mL	IVD

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analyte

Inactivated Parvovirus B19

Catalog #	Panel Member 1	Panel Member 2	Panel Member 3	Panel Member 4	Panel Member 5
B19P100	500 IU/mL	5,000 IU/mL	50,000 IU/mL	500,000 IU/mL	5,000,000 IU/mL

Exact Diagnostics CMV



- Verification Panel offers a wide dynamic range
- Manufacturing targets established using ddPCR
- Calibrated against the 1st WHO International Standard for Cytomegalovirus (09/162)
- High quality matrix to ensure lot-to-lot reproducibility
- True EDTA plasma matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
CMVL101*	Low Run Control	5 x 1.0 mL	IVD
CMVH102*	High Run Control	5 x 1.0 mL	IVD
Panels			
CMVP100	Verification Panel6 x 1.4 mL	IVD
CMVP200	Verification Panel6 x 1.4 mL	IVD
CMVL201	1E3 IU/mL Precision9 x 1.4 mL	IVD
CMVH202	5E5 IU/mL Precision9 x 1.4 mL	IVD
CMVLOD	CMV LOD	10 x 1.0 mL	IVD

Verification Bundle

CMVB200 Verification Bundle . . . CMVP200 (4 kits), CMVL201 (2 kits), CMVH202 (2 kits), CMVLOD (2 kits)

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analyte

Inactivated CMV

Complementary product:

See page 26 for Negative Run Control (Cat# NEG000/NEG050)

See Page 26 for Bulk Urine Negative Run Control (Cat# 12019386)

Catalog #	Panel Member 1	Panel Member 2	Panel Member 3	Panel Member 4	Panel Member 5	Panel Member 6
CMVP100	400 IU/mL	4,000 IU/mL	40,000 IU/mL	400,000 IU/mL	4,000,000 IU/mL	N/A
CMVP200	200 IU/mL	400 IU/mL	4,000 IU/mL	40,000 IU/mL	400,000 IU/mL	4,000,000 IU/mL

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Transplantation

Exact Diagnostics EBV



- Verification Panel offers a wide dynamic range
- Manufacturing targets established using ddPCR
- Calibrated against the 1st WHO International Standard for Epstein-Barr virus (09/260)
- High quality matrix to ensure lot-to-lot reproducibility
- True EDTA plasma matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
EBVL101*	Low Run Control	5 x 1.0 mL	IVD
EBVH102*	High Run Control	5 x 1.0 mL	IVD
Panels			
EBVP100	Verification Panel	5 x 1.0 mL	IVD
EBVP200	Verification Panel	7 x 1.4 mL	IVD
EBVL201	1E3 IU/mL Precision	9 x 1.4 mL	IVD
EBVH202	5E5 IU/mL Precision	9 x 1.4 mL	IVD
EBVL0D	EBV LOD	10 x 1.0 mL	IVD

Verification Bundle

EBVB200 Verification Bundle . . . EBVP200 (4 kits), EBVL201 (2 kits), EBVH202 (2 kits), EBVL0D (2 kits)

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analyte

Inactivated EBV

Complementary product:

See page 26 for Negative Run Control (Cat# NEG000/NEG050)

Catalog #	Panel Member 1	Panel Member 2	Panel Member 3	Panel Member 4	Panel Member 5	Panel Member 6	Panel Member 7
EBVP100	500 IU/mL	5,000 IU/mL	50,000 IU/mL	500,000 IU/mL	5,000,000 IU/mL		
EBVP200	100 IU/mL	500 IU/mL	5,000 IU/mL	50,000 IU/mL	500,000 IU/mL	5,000,000 IU/mL	50,000,000 IU/mL

Exact Diagnostics HHV-6



- Verification Panel offers a wide dynamic range
- Manufacturing targets established using ddPCR
- Calibrated against the 1st WHO International Standard for HHV-6B (15/266)
- High quality matrix to ensure lot-to-lot reproducibility
- True EDTA plasma matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
HV6L101*	Low Run Control	5 x 1.0 mL	IVD
HV6H102*	High Run Control	5 x 1.0 mL	IVD
Panels			
HV6AP100	Verification Panel	5 x 1.0 mL	IVD
HV6BP100	Verification Panel	5 x 1.0 mL	IVD

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analytes

Inactivated HHV-6A

Inactivated HHV-6B

Complementary product:

See page 26 for Negative Run Control (Cat# NEG000/NEG050)

Catalog #	Panel Member 1	Panel Member 2	Panel Member 3	Panel Member 4	Panel Member 5
HV6AP100	500 cp/mL	5,000 cp/mL	50,000 cp/mL	500,000 cp/mL	5,000,000 cp/mL
HV6BP100	500 cp/mL	5,000 cp/mL	50,000 cp/mL	500,000 cp/mL	5,000,000 cp/mL

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Transplantation

Exact Diagnostics JCV



- Calibrated against the 1st WHO International Standard for JC Virus (14/114)
- Manufacturing targets established using ddPCR
- Verification Panel offers a wide dynamic range
- High quality matrix to ensure lot-to-lot reproducibility
- True EDTA plasma matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
JCVL101*	Low Run Control	5 x 1.0 mL	IVD
JCVH102*	High Run Control	5 x 1.0 mL	IVD
Panels			
JCVP100	Verification Panel	6 x 1.0 mL	IVD

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analyte

Inactivated JCV

Complementary product:

See page 26 for Negative Run Control (Cat# NEG000/NEG050)

See Page 26 for Bulk Urine Negative Run Control (Cat# 12019386)

Catalog #	Panel Member 1	Panel Member 2	Panel Member 3	Panel Member 4	Panel Member 5	Panel Member 6
JCVP100	200 IU/mL	2,000 IU/mL	20,000 IU/mL	200,000 IU/mL	2,000,000 IU/mL	20,000,000 IU/mL

Sexually-Transmitted Infections

Exact Diagnostics CT/NG



- Manufacturing targets established using ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic Matrix
- Run Control has 30 day open vial stability when stored at 2-8°C
- Barcoded vials compatible with automated platforms

Ordering Information

Catalog #	Description	Format	Intended Use
Run Control			
CTNGPOS*	Positive Run Control	6 x 2.0 mL	IVD
Panels			
CTNGP200	Verification Panel	9 x 1.0 mL	IVD

Verification Bundle

CTNGB200 Verification Bundle CTNGP200 (5 kits)

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analytes

Chlamydia trachomatis (CT)

Neisseria gonorrhoeae (NG)

Complementary product:

See page 25 for STI Negative Run Control (Cat# NEGSTI)

See Page 26 for Bulk Urine Negative Run Control (Cat# 12019386)

Catalog #	Panel Members		
CTNGP200	CT (3X LOD)	CT (5X LOD)	CT (10X LOD)
	NG (3X LOD)	NG (5X LOD)	NG (10X LOD)
	CT/NG (3X LOD)	CT/NG (5X LOD)	CT/NG (10X LOD)

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Sexually-Transmitted Infections

Exact Diagnostics TV/MG



- Manufacturing targets established using ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic matrix
- 24 hour stability when stored at 2-8°C

Ordering Information

Catalog #	Description	Format	Intended Use
Panel			
TVMGP200	Verification Panel	9 x 1.0 mL	IVD
Verification Bundle			
TVMGB200	Verification Bundle	TVMGP200 (5 kits)	

Analytes

Trichomonas vaginalis
Mycoplasma genitalium

Complementary product:
 See Page 26 for Bulk Urine Negative Control (Cat# 12019386)

Catalog #	Panel Members		
TVMGP200	TV (3X LOD)	TV (5X LOD)	TV (10X LOD)
	MG (3X LOD)	MG (5X LOD)	MG (10X LOD)
	TV/MG (3X LOD)	TV/MG (5X LOD)	TV/MG (10X LOD)

Exact Diagnostics HPV



- Manufacturing targets established using ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- Formulated in PreservCyt
- Controls stable for 10 days when stored at room temperature
- Negative run control contains human cells to simulate adequate specimen sampling

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
HPV16C*	HPV Genotype 16 Control . . .	10 x 5.0 mL	IVD
HPV18C*	HPV Genotype 18 Control . . .	10 x 5.0 mL	IVD
HPV68C*	HPV Genotype 68 Control . . .	10 x 5.0 mL	IVD
HPVNEG*	HPV Negative Control	10 x 5.0 mL	IVD
Panels			
HPVP100	HPV Control Set		
	(1) HPV16C, (1) HPV18C, (1) HPV68C, (1) HPVNEG		
HPVP200	HPV Verification Panel	16 x 1.4 mL	IVD
Verification Bundle			
HPVB200	Verification Bundle	HPVP200 (5 kits)	

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analytes

Human Papillomavirus (HPV) Genotype 16
 Human Papillomavirus (HPV) Genotype 18
 Human Papillomavirus (HPV) Genotype 68
 Human Cell Line

Catalog #	Panel Members		
HPVP200	HPV 16 (3X LOD)	HPV 16 (5X LOD)	HPV 16 (10X LOD)
	HPV 18 (3X LOD)	HPV 18 (5X LOD)	HPV 18 (10X LOD)
	3 of HPV 16/18/68 (3X LOD)	HPV 16/18/68 (5X LOD)	HPV 16/18/68 (10X LOD)

Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Exact Diagnostics HSV-1



- Manufacturing targets established using ddPCR
- Verification Panel offers a wide dynamic range
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic CSF matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Complementary product:

See page 26 for Negative Run Control (Cat# NEG000/NEG050)

See Page 26 for Bulk Urine Negative Control (Cat# 12019386)

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
HSV1L101*	Low Run Control	5 x 1.0 mL	IVD
HSV1H102*	High Run Control	5 x 1.0 mL	IVD
Panels			
HSV1P100	Verification Panel	5 x 1.0 mL	IVD
HSVP200	Verification Panel	9 x 1.0 mL	IVD
Verification Bundle			
HSVB200	Verification Bundle	HSVP200 (5 kits)	

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analyte

Inactivated HSV-1

HSV-1

Catalog #	Panel Member 1	Panel Member 2	Panel Member 3	Panel Member 4	Panel Member 5
HSV1P100	200 cp/mL	2,000 cp/mL	20,000 cp/mL	200,000 cp/mL	2,000,000 cp/mL

HSV-1/HSV-2

Catalog #	Panel Members			
HSVP200	HSV-1 (3X LOD)	HSV-1 (5X LOD)	HSV-1 (10X LOD)	
	HSV-2 (3X LOD)	HSV-2 (5X LOD)	HSV-2 (10X LOD)	
	HSV-1/HSV-2 (3X LOD)	HSV-1/HSV-2 (5X LOD)	HSV-1/HSV-2 (10X LOD)	

Exact Diagnostics HSV-2



- Manufacturing targets established using ddPCR
- Verification Panel offers a wide dynamic range
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic CSF matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Complementary product:

See page 26 for Negative Run Control (Cat# NEG000/NEG050)

See Page 26 for Bulk Urine Negative Run Control (Cat# 12019386)

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
HSV2L101*	Low Run Control	5 x 1.0 mL	IVD
HSV2H102*	High Run Control	5 x 1.0 mL	IVD
Panels			
HSV2P100	Verification Panel	5 x 1.0 mL	IVD
HSVP200	Verification Panel	9 x 1.0 mL	IVD
Verification Bundle			
HSVB200	Verification Bundle	HSVP200 (5 kits)	

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analyte

Inactivated HSV-2

HSV-2

Catalog #	Panel Member 1	Panel Member 2	Panel Member 3	Panel Member 4	Panel Member 5
HSV2P100	200 cp/mL	2,000 cp/mL	20,000 cp/mL	200,000 cp/mL	2,000,000 cp/mL

HSV-1/HSV-2

Catalog #	Panel Members			
HSVP200	HSV-1 (3X LOD)	HSV-1 (5X LOD)	HSV-1 (10X LOD)	
	HSV-2 (3X LOD)	HSV-2 (5X LOD)	HSV-2 (10X LOD)	
	HSV-1/HSV-2 (3X LOD)	HSV-1/HSV-2 (5X LOD)	HSV-1/HSV-2 (10X LOD)	

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Virology

Exact Diagnostics HBV



- Calibrated against the 3rd WHO International Standard for Hepatitis B (10/264)
- Verification Panel offers a wide dynamic range
- Manufacturing targets established using ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- True EDTA plasma matrix
- Run controls stable for 5 days when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Analyte

Inactivated HBV

Complementary product:

See page 26 for Negative Run Control (Cat# NEG000/NEG050)

HBV

Catalog #	Panel Member 1	Panel Member 2	Panel Member 3	Panel Member 4	Panel Member 5	Panel Member 6	Panel Member 7
HBVP100	100 IU/mL	1,000 IU/mL	10,000 IU/mL	100,000 IU/mL	1,000,000 IU/mL	10,000,000 IU/mL	100,000,000 IU/mL
HBVP200	100 IU/mL	1,000 IU/mL	10,000 IU/mL	100,000 IU/mL	1,000,000 IU/mL	10,000,000 IU/mL	100,000,000 IU/mL

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
HBVL101*	HBV-1 Low Run Control	5 x 1.0 mL	IVD
HBVH102*	HBV-1 High Run Control	5 x 1.0 mL	IVD
Panels			
HBVP100	Verification Panel	7 x 1.0 mL	IVD
HBVP200	Verification Panel	7 x 1.4 mL	IVD
HBVL201	1E3 IU/mL Precision	9 x 1.4 mL	IVD
HBVH202	1E6 IU/mL Precision	9 x 1.4 mL	IVD
HBVLOD	HBV LOD	10 x 1.0 mL	IVD

Verification Bundle

HBVB200 Verification Bundle HBVP200 (5 kits)

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Exact Diagnostics HCV



- Calibrated against the 4th WHO International Standard for Hepatitis C (06/102)
- Verification Panel offers a wide dynamic range
- Manufacturing targets established using ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- True EDTA plasma matrix
- Run controls stable for 5 days when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Analyte

Recombinant HCV

Complementary product:

See page 26 for Negative Run Control (Cat# NEG000/NEG050)

HCV

Catalog #	Panel Member 1	Panel Member 2	Panel Member 3	Panel Member 4	Panel Member 5	Panel Member 6	Panel Member 7
HCVP100	100 IU/mL	300 IU/mL	3,000 IU/mL	30,000 IU/mL	300,000 IU/mL	3,000,000 IU/mL	30,000,000 IU/mL
HCVP200	100 IU/mL	300 IU/mL	3,000 IU/mL	30,000 IU/mL	300,000 IU/mL	3,000,000 IU/mL	30,000,000 IU/mL

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
HCVL101*	HCV-1 Low Run Control	5 x 1.0 mL	IVD
HCVH102*	HCV-1 High Run Control	5 x 1.0 mL	IVD
Panels			
HCVP100	Verification Panel	7 x 1.0 mL	IVD
HCVP200	HCV Verification Panel	7 x 1.4 mL	IVD
HCVL201	1E3 IU/mL Precision	9 x 1.4 mL	IVD
HCVH202	1E6 IU/mL Precision	9 x 1.4 mL	IVD
HCVLOD	HCV LOD	10 x 1.0 mL	IVD
HCVGTP†	Genotype Panel		
HCVGT 1a, HCVGT 1b, HCVGT 2, HCVGT 3, HCVGT 4	HCVGT 5		

Verification Bundle

HCVB200 Verification Bundle . . HCVP200 (4 kits), HCVL201 (2 kits), HCVH202 (2 kits), HCVLOD (2 kits)

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

†For research use only. Not for use in diagnostic procedures

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Virology

Exact Diagnostics HIV



- A family of unassayed external controls, verification panels and verification panel bundles used in the determination of values in the measurement of HIV-1 and HIV-2 RNA
- Verification Panel offers a wide dynamic range
- Calibrated against the 3rd WHO International Standard for HIV-1 (10/152)
- Calibrated against the 2nd WHO International Standard for HIV-2 (16/296)
- Calibrated with digital droplet PCR
- High quality matrix to ensure lot-to-lot reproducibility
- True EDTA Plasma
- Run controls stable for 5 days when stored at 2-8°C
- Controls can be frozen and reused a maximum of two (2) times

Analytes

Inactivated HIV-1
Inactivated HIV-2

Complementary product:

See page 26 for Negative Run Control (Cat# NEG000/NEG050)

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
HIVL101*	HIV-1 Low Run Control	5 x 1.0 mL	IVD
HIVH102*	HIV-1 High Run Control5 x 1.0 mL	IVD
Panels			
HIVP100	HIV-1 Verification Panel	7 x 1.0 mL	IVD
HIVP200	HIV-1 Verification Panel	7 x 1.4 mL	IVD
HIVL201	HIV-1 1E3 cp/mL Precision	9 x 1.4 mL	IVD
HIVH202	HIV-1 1E5 cp/mL Precision	9 x 1.4 mL	IVD
HIVLOD	HIV-1 LOD	10 x 1.0 mL	IVD
HIVQP200 [†]	HIV-1/HIV-2 Verification Panel	12 x 1.4 mL	RUO
Verification Bundles			
HIVB200	HIV-1 Verification Bundle...HIVP200 (4 kits), HIVL201 (2 kits), HIVH202 (2 kits), HIVLOD (2 kits)		
HIVQB200 [†]	HIV-1/HIV-2 Verification Bundle HIVQP200 (5 kits)		

*This product is CE Marked and is undergoing conversion to IVDR CE Marking
[†]For research use only. Not for use in diagnostic procedures

HIV-1

Catalog #	Panel Member 1	Panel Member 2	Panel Member 3	Panel Member 4	Panel Member 5	Panel Member 6	Panel Member 7
HIVP100	100 cp/mL	600 cp/mL	6,000 cp/mL	60,000 cp/mL	600,000 cp/mL	2,000,000 cp/mL	6,000,000 cp/mL
HIVP200	100 cp/mL	600 cp/mL	6,000 cp/mL	60,000 cp/mL	600,000 cp/mL	2,000,000 cp/mL	6,000,000 cp/mL

HIV-1/HIV-2

Catalog #	Panel Members		
HIVQP200	HIV-1 (3X LOD)	HIV-1 (5X LOD)	HIV-1 (10X LOD)
	HIV-2 (3X LOD)	HIV-2 (5X LOD)	HIV-2 (10X LOD)
	HIV-1/HIV-2 (2.5X LOD)	HIV-1/HIV-2 (5X LOD)	HIV-1/HIV-2 (10X LOD)

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Virology

Exact Diagnostics HBV, HCV and HIV Analytical Measurement Range



- HBV: Calibrated against the 3rd WHO International Standard for Hepatitis B (10/264)
- HCV: Calibrated against the 4th WHO International Standard for Hepatitis C (06/102)
- HIV: Calibrated against the 3rd WHO International Standard for HIV-1 (10/152)
- Manufacturing targets established using ddPCR
- AMR Panel offers a wide dynamic range
- High quality matrix to ensure lot-to-lot reproducibility
- True EDTA plasma matrix
- 24 hour stability when stored at 2-8°C

Complementary product:
See page 26 for Negative Run Control (Cat# NEG000/NEG050)

Ordering Information

Catalog #	Description	Format	Intended Use
Panels			
HBVAMR	HBV Analytical Measuring Range Panel . 3 x 1.0 mL	IVD	
HCVAMR	HCV Analytical Measuring Range Panel . 3 x 1.0 mL	IVD	
HIVAMR	HIV Analytical Measuring Range Panel . . 3 x 1.0 mL	IVD	
AMR Bundle			
AMRP100	Analytical Measuring Range BundleHBVAMR (1 kit), HCVAMR (1 kit), HIVAMR (1 kit)		

Analytes

Inactivated HBV
Recombinant HCV
Inactivated HIV-1

Catalog #	Panel Member 1	Panel Member 2	Panel Member 3
HBVAMR	100 IU/mL	100,000 IU/mL	10,000,000 IU/mL
HCVAMR	500 IU/mL	50,000 IU/mL	5,000,000 IU/mL
HIVAMR	600 IU/mL	60,000 IU/mL	6,000,000 IU/mL

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Negative Controls

Exact Diagnostics STI Negative



- Screened negative for 2 molecular targets
- High quality synthetic matrix to ensure lot-to-lot reproducibility
- Contains human cells to simulate adequate specimen sampling
- 30 day open vial stability when stored at 2-8°C
- Barcoded vials compatible with automated platforms

Ordering Information

Catalog #	Description	Format	Intended Use
-----------	-------------	--------	--------------

Run Control

NEGSTI*	STI Negative Run Control.	6 x 2.0 mL	IVD
---------	-----------------------------------	------------	-----

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analytes

Non-reactive for:

Chlamydia trachomatis (CT)
Neisseria gonorrhoeae (NG)

Exact Diagnostics Synthetic Negative



- Screened negative for 9 molecular targets
- High quality synthetic matrix to ensure lot-to-lot reproducibility
- 5 day stability when stored at 2-8°C

Ordering Information

Catalog #	Description	Format	Intended Use
-----------	-------------	--------	--------------

Run Control

NEGSYN*	Synthetic Negative Run Control . . .	5 x 1.0 mL	IVD
---------	--------------------------------------	------------	-----

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analytes

Non-reactive for:

<i>Anaplasma phagocytophilum</i>	Enterovirus (EV) Coxsackievirus A9
<i>Babesia microti</i>	Herpes Simplex Virus-1 (HSV-1)
<i>Bartonella quintana</i>	Herpes Simplex Virus-2 (HSV-2)
<i>Borrelia burgdorferi</i>	Varicella Zoster Virus (VZV)
<i>Ehrlichia chaffeensis</i>	

Exact Diagnostics HAI Negative



- Screened negative for 3 molecular targets
- High quality synthetic matrix to ensure lot-to-lot reproducibility
- 5 day stability when stored at 2-8°C

Ordering Information

Catalog #	Description	Format	Intended Use
-----------	-------------	--------	--------------

Run Control

NEGHA1*	HAI Negative Run Control.	5 x 1.0 mL	IVD
---------	-----------------------------------	------------	-----

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analytes

Non-reactive for:

Clostridium difficile (*C. difficile*)
Methicillin-resistant *Staphylococcus aureus* (MRSA)
Methicillin-susceptible *Staphylococcus aureus* (MSSA)

Negative Controls

Exact Diagnostics Negative



- Screened negative for 16 molecular targets
- High quality matrix to ensure lot-to-lot reproducibility
- True EDTA plasma matrix
- 5 day stability when stored at 2-8°C
- Control can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
NEG000*	Negative Run Control	5 x 1.0 mL	IVD
NEG050*	Bulk Negative Run Control	1 x 50 mL	IVD

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analytes

Non-reactive for:

Adenovirus DNA	HBsAg	HHV-6 DNA	JC Virus DNA
BKV DNA	HBV DNA	HIV-1/2 Antibodies	VZV DNA
CMV DNA	HCV Antibodies	HIV-1 RNA	Zika Virus RNA
EBV DNA	HCV RNA	HSV-1/2 DNA	

NEW | Exact Diagnostics Bulk CSF Negative



- Screened negative for 9 molecular targets
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic CSF matrix
- 5 day stability when stored at 2-8°C
- Control can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Control			
12019432*	Bulk CSF Negative Run Control . . .	1 x 50 mL	IVD

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analytes

Non-reactive for:

<i>Anaplasma phagocytophilum</i>	Coxsackievirus A9	Enterovirus (EV)
<i>Babesia microti</i>	Herpes Simplex Virus-1 (HSV-1)	
<i>Bartonella quintana</i>	Herpes Simplex Virus-2 (HSV-2)	
<i>Borrelia burgdorferi</i>	Varicella Zoster Virus (VZV)	
<i>Ehrlichia chaffeensis</i>		

NEW | Exact Diagnostics Bulk Urine Negative



- Screened negative for 10 molecular targets
- High quality matrix to ensure lot-to-lot reproducibility
- Pooled human urine
- 5 day stability when stored at 2-8°C
- Control can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Control			
12019386†	Bulk Urine Negative Run Control . . .	1 x 50 mL	RUO

†For research use only. Not for use in diagnostic procedures

Analytes

Non-reactive for:

Adenovirus (ADV)	John Cunningham Virus (JCV)
BK Virus (BKV)	<i>Mycoplasma genitalium</i> (MG)
<i>Candida auris</i> (<i>C. auris</i>)	<i>Neisseria gonorrhoeae</i> (NG)
<i>Chlamydia trachomatis</i> (CT)	<i>Trichomonas vaginalis</i> (TV)
Cytomegalovirus (CMV)	Zika Virus (ZKV)

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Meningitis and Encephalitis

Exact Diagnostics VZV



- Manufacturing targets established using ddPCR
- Verification Panel offers a wide dynamic range
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic CSF matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
VZVL101*	Low Run Control5 x 1.0 mL	IVD
VZVH102*	High Run Control5 x 1.0 mL	IVD
Panel			
VZVP100	Verification Panel	5 x 1.0 mL	IVD

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analyte

Inactivated Varicella Zoster Virus (VZV)

Complementary product:

See page 25 for Synthetic Negative Run Control (Cat# NEGSYN)
See Page 26 for Bulk CSF Negative Run Control (Cat# 12019432)

Catalog #	Panel Member 1	Panel Member 2	Panel Member 3	Panel Member 4	Panel Member 5
VZVP100	500 cp/mL	2,000 cp/mL	20,000 cp/mL	200,000 cp/mL	2,000,000 cp/mL

Exact Diagnostics Enterovirus



- Manufacturing targets established using ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Control			
EVRC*	Run Control	5 x 1.0 mL	IVD

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analyte

Inactivated Coxsackievirus A9

Complementary product:

See page 25 for Synthetic Negative Run Control (Cat# NEGSYN)
See Page 26 for Bulk CSF Negative Run Control (Cat# 12019432)

NEW | Exact Diagnostics GBS



- Internal value assignment using Droplet Digital PCR (ddPCR)
- High-quality clinically-relevant Lim broth matrix to ensure lot-to-lot reproducibility while monitoring pre-analytical workflow steps requiring the dipping of a swab in Lim broth
- Run controls have 5 days open vial stability when stored at 2-8°C or are single use (depending on the workflow)
- Negative run control contains human cells to simulate adequate specimen sampling

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
12020069*	Positive Run Control	5 x 0.25 mL	IVD
12020068*	Negative Run Control	5 x 0.25 mL	IVD
Panel			
12020070†	GBS Verification & Validation Panel .	5 x 0.25 mL	RUO
Verification Bundle			
17009668†	Verification Bundle	12020070 (4 kits)	

*This product is CE Marked and is undergoing conversion to IVDR CE Marking
†For research use only. Not for use in diagnostic procedures

Catalog #	Panel Members
12020070*	1,000 cp/mL

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Hospital-Associated Infections

Exact Diagnostics *C. difficile* Verification Bundle



- Manufacturing targets established using ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic matrix
- 24 hour stability when stored at 2-8°C

Ordering Information

Catalog #	Description	Format	Intended Use
Panel			
CDFP200	Verification Panel	3 x 1.0 mL	IVD
Verification Bundle			
CDFB200	Verification Bundle	CDFP200 (5 kits)	

Analyte

Inactivated *C. difficile*

Complementary product:

See page 25 for HAI Negative Run Control (Cat# NEGHA1)

Catalog #	Panel Members		
CDFP200	<i>C. difficile</i> (3X LOD)	<i>C. difficile</i> (5X LOD)	<i>C. difficile</i> (10X LOD)

Exact Diagnostics MRSA Verification Bundle



- Manufacturing targets established using ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic matrix
- 24 hour stability when stored at 2-8°C

Ordering Information

Catalog #	Description	Format	Intended Use
Panel			
MRSAP200	Verification Panel	6 x 1.0 mL	IVD
Verification Bundle			
MRSAB200	Verification Bundle	MRSAP200 (5 kits)	

Analytes

Inactivated Methicillin-resistant *Staphylococcus aureus* (MRSA)
 Inactivated Methicillin-susceptible *Staphylococcus aureus* (MSSA)

Complementary product:

See page 25 for HAI Negative Run Control (Cat# NEGHA1)

Catalog #	Panel Members		
MRSAP200	MRSA (3X LOD)	MRSA (5X LOD)	MRSA (10X LOD)
	MSSA (3X LOD)	MSSA (5X LOD)	MSSA (10X LOD)

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Hospital-Associated Infections

Amplichek II



- Liquid, multi-analyte quality control designed to monitor the performance of *in vitro* nucleic acid test systems for Healthcare Associated Infections (HAI)
- Use with multiple specimen types and sampling tools including swabs and transfer pipets
- Convenient single use configuration to minimize contamination
- Contains intact, inactivated bacteria to monitor the full NAT process from extraction to amplification and detection
- Negative, Levels 1 and 2 are suitable for Cepheid Xpert assays

Contact your local Bio-Rad office for additional assay compatibility and Level selection information.

Ordering Information

Catalog #	Description	Format	Intended Use
12000201	Negative	10 x 0.2 mL	IVD
12000202	Level 1	10 x 0.2 mL	IVD
12000203	Level 2	10 x 0.2 mL	IVD
12000204	Level 3	10 x 0.2 mL	IVD
12000575	Negative MiniPak	1 x 0.2 mL	IVD
12000576	Level 1 MiniPak	1 x 0.2 mL	IVD
12000577	Level 2 MiniPak	1 x 0.2 mL	IVD
12000578	Level 3 MiniPak	1 x 0.2 mL	IVD

Analytes

Clostridium difficile (*C diff.*)
 Methicillin-resistant *Staphylococcus aureus* (MRSA)
 Methicillin-sensitive *Staphylococcus aureus* (MSSA)
 Vancomycin-resistant *enterococci* (VRE or vanR)

Product Level Selection by Analyte

- Reactive
- Non-reactive

Analytes

Analyte	Level	Negative* (Non-reactive)	Level 1 (MSSA)	Level 2 (Low Positive)	Level 3 (High Positive)
<i>Clostridium difficile</i> (<i>C diff.</i>) genes: Toxin B, Binary Toxin, TcdC		●	●	●	●
Methicillin-Resistant <i>Staphylococcus aureus</i> (MRSA) genes: SPA, SCCmec, mecA		●	●	●	●
Methicillin-Sensitive <i>Staphylococcus aureus</i> (MSSA) gene: SPA		●	●		
Vancomycin-Resistant <i>enterococci</i> (VRE or vanR) genes: vanA, vanB**		●	●	●	●

Blood Screening

Exact Diagnostics HBV, HCV and HIV-1 Screen Controls



- Calibrated against the 3rd WHO International Standard for HBV (10/264)
- Calibrated against the 4th WHO International Standard for HCV (06/102)
- Calibrated against the 3rd WHO International Standard for HIV-1 (10/152)
- Citrate plasma matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
HBVSCREEN	HBV DNA Screen Control	10 x 1.4 mL	IVD
HCVSCREEN	HCV RNA Screen Control	10 x 1.4 mL	IVD
HIVSCREEN	HIV-1 RNA Screen Control	10 x 1.4 mL	IVD
VIRSCREEN	HIV-1, HCV, HBV Screen Control .	10 x 1.4 mL	IVD

Analytes

Inactivated HBV Inactivated HCV Inactivated HIV-1

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Respiratory

Exact Diagnostics SARS-CoV-2, Flu, RSV



- Whole organism SARS-CoV-2, Influenza A, Influenza B, Respiratory Syncytial Virus (A)
- Manufacturing targets established using ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic matrix
- Contains human genomic DNA (gDNA) to simulate adequate specimen sampling
- Controls can be frozen and reused a maximum of five (5) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
COVFLU*	Positive Run Control	5 x 1.0 mL	IVD
COVFLUNEG*	Negative Run Control.	5 x 1.0 mL	IVD
Panel			
COVFLUP200†	Verification Panel	15 x 1.0 mL	RUO
Verification Bundle			
COVFLUB200†	SARS-CoV-2, Flu, RSV Verification Bundle		
COVFLUP200	(5 kits)		

*This product is CE Marked and is undergoing conversion to IVDR CE Marking
 †For research use only. Not for use in diagnostic procedures

Targets

Whole Organism

- SARS-CoV-2
- Influenza A-H1N1
- Influenza B
- Respiratory Syncytial Virus (A)

Catalog #	Panel Members		
COVFLUP200	Flu A (3X LOD)	Flu A (5X LOD)	Flu A (10X LOD)
	Flu B (3X LOD)	Flu B (5X LOD)	Flu B (10X LOD)
	RSV A (3X LOD)	RSV A (5X LOD)	RSV A (10X LOD)
	SARS-CoV-2 (3X LOD)	SARS-CoV-2 (5X LOD)	SARS-CoV-2 (10X LOD)
	SARS-CoV-2, Flu, RSV (3X LOD)	SARS-CoV-2, Flu, RSV (5X LOD)	SARS-CoV-2, Flu, RSV (10X LOD)

Exact Diagnostics SARS-CoV-2



- Contains E, N, ORF1ab, RdRP and S genes
- Quantitation with ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic matrix
- Contains human genomic DNA (gDNA) to simulate adequate specimen sampling
- 24 hour stability when stored at 2-8°C

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
COV019CE*	Positive Run Control	5 x 0.3 mL	IVD
COV000CE*	Negative Run Control.	5 x 0.3 mL	IVD
Standards			
COV019†	Standard	5 x 0.3 mL	RUO
COV000†	Negative	5 x 0.3 mL	RUO

*This product is CE Marked and is undergoing conversion to IVDR CE Marking
 †For research use only. Not for use in diagnostic procedures

Targets

SARS-CoV-2 Genes

- | | |
|-----------------------|------------------------|
| Envelope (E) Gene | RdRP (RNA-dependent |
| Nucleocapsid (N) Gene | RNA polymerase) Gene |
| ORF1ab Gene | Spike Protein (S) Gene |

Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Respiratory

Exact Diagnostics SARS-CoV-2 S Gene Variant Controls



- Quantitation with ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic matrix
- Contains human genomic DNA (gDNA) to simulate adequate specimen sampling
- 7 day open-vial stability when stored at 2-8°C

Ordering Information

Catalog # Description

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

COVA	Alpha Variant Control	1 x 0.25 mL
COVB	Beta Variant Control.	1 x 0.25 mL
COVDK	Delta-Kappa Variant Control	1 x 0.25 mL
COVE	Epsilon Variant Control	1 x 0.25 mL
COVG	Gamma Variant Control	1 x 0.25 mL

Targets

Synthetic RNA Transcripts

Alpha Variant (B.1.1.7)	Delta-Kappa Variants
Beta Variant (B.1.351)	(B.1.617.1, B.1.617.2 and
Epsilon Variant (B.1.427, B.1.429)	AY sublineages, B.1.617.3)
Gamma Variant (P.1)	

Complementary product:

See page 30 for SARS-CoV-2, Negative (Cat# COV000)

Exact Diagnostics RP Positive & Negative



- 23 Analyte respiratory control
- Manufacturing targets established using ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic matrix

Ordering Information

Catalog #	Description	Format	Intended Use
Run Controls			
RPPOS*	RP Positive Control	6 x 0.3 mL	IVD
RPNEG*	RP Negative Control	6 x 0.3 mL	IVD

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analytes

Adenovirus	Coronavirus OC43	Middle Eastern Respiratory	Respiratory Syncytial Virus (A)
<i>Bordetella parapertussis</i>	Human metapneumovirus (RNA IVT)	Syndrome (RNA IVT)	Respiratory Syncytial Virus (B)
<i>Bordetella pertussis</i>	Influenza A H1N1	Parainfluenza 1	Severe Acute Respiratory Syndrome
<i>Chlamydia pneumoniae</i>	Influenza A H1N1-09 (RNA IVT)	Parainfluenza 2	Coronavirus 2
Coronavirus 229E	Influenza A H3N2 (RNA IVT)	Parainfluenza 3	
Coronavirus HKU1 (RNA IVT)	Influenza B	Parainfluenza 4	
Coronavirus NL63 (RNA IVT)	<i>Mycoplasma pneumoniae</i>	Rhinovirus	

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Vector-Borne

Exact Diagnostics Anaplasma



- Manufacturing targets established using ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Control			
ANARC*	Anaplasma Run Control	5 x 1.0 mL	IVD

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analyte

Anaplasma phagocytophilum

Complementary product:

See page 25 for Synthetic Negative Run Control (Cat# NEGSYN)

Exact Diagnostics Babesia



- Manufacturing targets established using ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Control			
BABRC*	Babesia Run Control	5 x 1.0 mL	IVD

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analyte

Babesia microti

Complementary product:

See page 25 for Synthetic Negative Run Control (Cat# NEGSYN)

Exact Diagnostics Bartonella



- Manufacturing targets established using ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Control			
BRTRC*	Bartonella Run Control	5 x 1.0 mL	IVD

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analyte

Bartonella quintana

Complementary product:

See page 25 for Synthetic Negative Run Control (Cat# NEGSYN)

See Page 26 for Bulk CSF Negative Run Control (Cat# 12019432)

Exact Diagnostics Borrelia



- Manufacturing targets established using ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Control			
BORRC*	Borrelia Run Control	5 x 1.0 mL	IVD

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analyte

Borrelia burgdorferi

Complementary product:

See page 25 for Synthetic Negative Run Control (Cat# NEGSYN)

See Page 26 for Bulk CSF Negative Run Control (Cat# 12019432)

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Vector Borne

Exact Diagnostics Ehrlichia



- Manufacturing targets established using ddPCR
- High quality matrix to ensure lot-to-lot reproducibility
- Synthetic matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Control			
ERLRC*	Ehrlichia Run Control	5 x 1.0 mL	IVD

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analyte

Ehrlichia chaffeensis

Complementary product:

See page 25 for Synthetic Negative Run Control (Cat# NEGSYN)

See Page 26 for Bulk CSF Negative Run Control (Cat# 12019432)

Exact Diagnostics ZKV



- Manufacturing targets established using ddPCR
- Verification Panel offers a wide dynamic range
- High quality matrix to ensure lot-to-lot reproducibility
- True EDTA plasma matrix
- 24 hour stability when stored at 2-8°C
- Controls can be frozen and reused a maximum of four (4) times

Ordering Information

Catalog #	Description	Format	Intended Use
Run Control			
ZKVL101*	Low Run Control.	5 x 1.0 mL	IVD
ZKVH102*	High Run Control	5 x 1.0 mL	IVD

Panels

ZKVP100	Verification Panel	5 x 1.0 mL	IVD
---------	------------------------------	------------	-----

*This product is CE Marked and is undergoing conversion to IVDR CE Marking

Analyte

Inactivated Zika Virus

Complementary product:

See page 26 for Negative Run Control (Cat# NEG000/NEG050)

See Page 26 for Bulk Urine Negative Run Control (Cat# 12019386)

Catalog #	Panel Member 1	Panel Member 2	Panel Member 3	Panel Member 4	Panel Member 5
ZKVP100	100 cp/mL	1,000 cp/mL	10,000 cp/mL	100,000 cp/mL	1,000,000 cp/mL

Liquichek Blood Gas Control



- Available in three distinct levels with values targeted at varying points within the clinical range
- Color coded matrix for easy identification
- Aqueous based matrix
- 4 year shelf life at 18-25°C

POCT methods listed

Ordering Information

Cat #	Description
571	Level 1 30 x 1.7 mL
572	Level 2 30 x 1.7 mL
573	Level 3 30 x 1.7 mL
570X	Trilevel MiniPak (3 of each level) 9 x 1.7 mL

Analytes

pCO₂
pH
pO₂

Liquichek Blood Gas Plus E Control



- Designed to monitor precision of laboratory measurements on Blood Gas and Ion Selective Electrode (ISE) instrumentation
- Clear aqueous matrix
- 4 year shelf life at 2-25°C

POCT methods listed

Ordering Information

Cat #	Description
576	Level 1 30 x 1.7 mL
577	Level 2 30 x 1.7 mL
578	Level 3 30 x 1.7 mL
575X	Trilevel MiniPak (3 of each level) 9 x 1.7 mL

Analytes

Calcium (Ionized) pO₂
Chloride Potassium
pCO₂ Sodium
pH

Liquichek Blood Gas Plus EGL Control



- Expanded test menu including pH, Blood Gas, Electrolytes and Metabolites
- Aqueous, clear matrix
- 3 year shelf life at 2-8°C
- Can be stored for up to 12 months at 20-25°C

POCT methods listed

Ordering Information

Cat #	Description
511	Level 1 30 x 1.7 mL
512	Level 2 30 x 1.7 mL
513	Level 3 30 x 1.7 mL
510X	Trilevel MiniPak (3 of each level) 9 x 1.7 mL

Analytes

Calcium (Ionized) pCO₂
Chloride pH
Glucose pO₂
Lactate (Lactic Acid) Potassium
Lithium Sodium
Magnesium

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Cardiac Advance

Liquichek Cardiac Advance Control



Available in Multiple Configurations

- A comprehensive, multi-analyte cardiac control optimized for high sensitivity Troponin testing
- Liquid, serum based

Standard Vials

- Level 1 Troponin values near limit of detection for most instruments.
- 28 month shelf life at -20°C to -70°C
- 20 day open-vial stability at 2-8°C, with exceptions

Compact Vials

- Level 1 Troponin values near limit of detection for most instruments.
- Barcoded vials for workflow optimization
- 28 month shelf life at -20°C to -50°C
- 5 day open-vial stability for all analytes when stored at 2-8°C



Compatible with Beckman Coulter, Siemens and Roche platforms. Abbott users should use Liquichek Cardiac Markers Plus Control LT.

Homocysteine not compatible with Siemens platforms. Siemens users should use Liquichek Homocysteine Control.

Ordering Information

Cat # Description

Standard Vials

12016447	Level 16 x 3 mL
12016448	Level 26 x 3 mL
12016449	Level 36 x 3 mL
12016450	Level 46 x 3 mL

Compact Vials

12016472	Level 1	6 x 2.5 mL
12016473	Level 2	6 x 2.5 mL
12016474	Level 3	6 x 2.5 mL
12016475	Level 4	6 x 2.5 mL

For InteliQ Cardiac Advance Control, see page 11

Analytes

BNP	Myoglobin
CK-MB Isoenzyme	NT-proBNP
CRP ¹	Troponin I ¹
Digitoxin	Troponin T* ¹
Galectin-3	
Homocysteine	*Not available in compact vial
	¹ Compatible with high sensitivity methods.

Liquichek Cardiac Markers Plus Control LT



Available in Multiple Configurations

- A comprehensive cardiac control with compatibility for high-sensitivity Troponin testing
- Liquid, human serum based

Standard Vials

- Choice of four Troponin levels (1, 1A, 1B, or 1C) for high-sensitivity Troponin methods
- 3 year shelf life at -20°C to -70°C
- 20 day open-vial stability at 2-8°C, with exceptions

Compact Vials

- Level 1C Troponin I values at or near the 99th percentile upper reference limit (URL) of Troponin I for TNIH method
- Barcoded vials for workflow optimization
- 5 day open-vial stability for all analytes when stored at 2-8°C
- 3 year shelf life when stored at -20°C to -50°C



POCT methods listed for standard vials

Ordering Information

Cat # Description

Standard Vials

146	Level 16 x 3 mL
27104	Level 1A6 x 3 mL
27105	Level 1B6 x 3 mL
27106	Level 1C6 x 3 mL
147	Level 26 x 3 mL
148	Level 36 x 3 mL
149	Level Low6 x 3 mL

146X	Level 1 MiniPak1 x 3 mL
27104X	Level 1A MiniPak1 x 3 mL
27105X	Level 1B MiniPak1 x 3 mL
27106X	Level 1C MiniPak1 x 3 mL
147X	Level 2 MiniPak1 x 3 mL
148X	Level 3 MiniPak1 x 3 mL
149X	Level Low MiniPak1 x 3 mL

Compact Vials

297	Level 1C	6 x 2.5 mL
298	Level 2	6 x 2.5 mL
299	Level 3	6 x 2.5 mL
298X	Trilevel MiniPak (1 of each level)	3 x 2.5 mL

For InteliQ Cardiac Markers Plus Control LT, see page 12

Analytes

BNP*	CRP ¹	Myoglobin	Troponin I ¹
Creatine Kinase (CK)	Digitoxin	NT-proBNP	Troponin T* ¹
CK-MB Isoenzyme	Homocysteine* ²		

*Available in standard vial configuration only.

¹Compatible with high sensitivity methods.

²No claim is made for expected value or stability.

Refer to myInsersts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Liquichek Cardiac Troponins Control



- Dedicated Troponins control
- Monitors the precision of high-sensitivity Troponin I and T assays near the 99th percentile upper reference limit and across the assay range
- Convenience of a liquid control
- 18 month shelf life at -20°C to -70°C
- Open-vial stability at 2-8°C
 - Troponin I: 14 days
 - Troponin T: 25 days

Ordering Information

Cat #	Description	
12001084	Level 1	6 x 3 mL
12006239	Level 2	6 x 3 mL
12006369	Level 3	6 x 3 mL
12001085	Level 1 MiniPak	1 x 3 mL
12006370	Level 2 MiniPak	1 x 3 mL
12006371	Level 3 MiniPak	1 x 3 mL

Analytes

Troponin I[†] Troponin T[†]

[†]Compatible with high sensitivity methods.

Liquichek Cardiac Markers Plus Control



- Comprehensive cardiac control monitoring the precision of non high-sensitivity Troponin assays
- Values provided for hs-CRP
- Liquid, human serum based
- 3 year shelf life at -20°C to -70°C
- 20 day open-vial stability at 2-8°C with exceptions

POCT methods listed

Ordering Information

Cat #	Description	
180	Trilevel (2 of each level)	6 x 3 mL
181	Level 1	6 x 3 mL
182	Level 2	6 x 3 mL
183	Level 3	6 x 3 mL
180X	Trilevel MiniPak (1 of each level)	3 x 3 mL

Analytes

BNP
Creatine Kinase (CK)
CK-MB Isoenzyme

CRP[†]
Digitoxin
Homocysteine*

Myoglobin
NT-proBNP

Troponin I
Troponin T

*No claim is made for expected values or stability.
[†]Compatible with high sensitivity methods.

Liquichek CK/LD Isoenzyme Control



- Monitors the precision of traditional and electrophoretic methods
- Liquid, human serum based
- 2 year shelf life at -20°C to -70°C
- 10 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description	
671	Level 1	6 x 1 mL
672	Level 2	6 x 1 mL
670X	Bilevel MiniPak (1 of each level)	2 x 1 mL

Analytes

Albumin
Creatine Kinase (CK)
CK Electrophoresis

CK-MB Isoenzyme
Globulin

LD Electrophoresis
LDH

Protein Electrophoresis
Protein Serum (Total)

Liquichek Homocysteine Control



- Designed to monitor Homocysteine precision across a wide range of target values
- Liquid, human serum based
- 3 year shelf life at -20°C to -70°C
- 30 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description	
287	Level 1	6 x 1 mL
288	Level 2	6 x 1 mL
289	Level 3	6 x 1 mL
288X	Trilevel MiniPak (1 of each level)	3 x 1 mL

Analyte

Homocysteine

For InteliQ Homocysteine Control, see page 13

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Liquid Assayed and Unassayed Multiquial



Available in Multiple Configurations

Provides a comprehensive range of chemistry analytes for monitoring the precision of general chemistry test procedures. Human serum based.

Liquid Assayed & Unassayed, Standard Vials

- 3 year shelf life at -20°C to -70°C
- 30 day unopened stability at 2-8°C for most analytes
- 14 day open-vial stability at 2-8°C for most analytes

Liquid Assayed, Compact Vials

- 5 day open-vial stability for all analytes when stored at 2-8°C
- 3 year shelf life when stored at -20°C to -50°C

POCT methods listed for standard vials

For InteliQ Assayed Multiquial Control, see page 11

Ordering Information

Cat #	Description	
Liquid Assayed Multiquial, Standard Vials		
694	Level 1	12 x 3 mL
695	Level 2	12 x 3 mL
696	Level 3	12 x 3 mL
695X	Trilevel MiniPak (1 of each level)	.3 x 3 mL
Liquid Assayed Multiquial, Compact Vials		
337	Level 1	12 x 2.5 mL
338	Level 2	12 x 2.5 mL
339	Level 3	12 x 2.5 mL
338X	Trilevel MiniPak (1 of each level)	3 x 2.5 mL
Liquid Unassayed Multiquial		
697	Level 1	12 x 10 mL
698	Level 2	12 x 10 mL
699	Level 3	12 x 10 mL
698X	Trilevel MiniPak (1 of each level)	3 x 10 mL

Analytes

	Standard Vials	Compact Vials		Standard Vials	Compact Vials		Standard Vials	Compact Vials
Acetaminophen	●	●	CO ₂	●	●	Osmolality	●	
Acid Phosphatase (Total)	●		Complement C3	●	●	PAP*	●	
Albumin	●	●	Complement C4	●	●	Phenobarbital	●	●
Albumin: Globulin Ratio	●		Copper	●		Phenytoin	●	●
Alkaline Phosphatase (ALP)	●	●	Cortisol	●		Phospholipids	●	
α-1-Antitrypsin	●	●	Creatine Kinase (CK)	●	●	Phosphorus	●	●
α-Hydroxybutyrate Dehydrogenase*	●		Creatinine	●	●	Potassium	●	●
ALT/SGPT	●	●	Digoxin	●	●	Prealbumin	●	●
Amikacin	●		Ethanol	●	●	Protein Serum (Total)	●	●
Amylase	●	●	Ferritin	●	●	Salicylate	●	●
Amylase (Pancreatic)	●		Gentamicin	●	●	Sodium	●	●
Apolipoprotein A-1	●	●	GGT	●	●	T3 (Total)	●	●
Apolipoprotein B	●	●	Globulin**	●		T3 Uptake/T-Uptake	●	●
AST/SGOT	●	●	Glucose	●	●	T4 (Free)	●	●
Bilirubin (Direct)	●	●	Haptoglobin	●	●	T4 (Total)	●	●
Bilirubin (Neonatal)	●		Immunoglobulin A (IgA)	●	●	Theophylline	●	●
Bilirubin (Total)	●	●	Immunoglobulin G (IgG)	●	●	Tobramycin	●	●
Calcium (Ionized)	●		Immunoglobulin M (IgM)	●	●	Transferrin	●	●
Calcium	●	●	Iron	●	●	Triglycerides	●	●
Carbamazepine	●	●	Iron (TIBC)	●	●	TSH	●	●
Ceruloplasmin	●	●	Iron (UIBC)	●		Urea	●	
Chloride	●	●	Lactate (Lactic Acid)	●	●	Urea Nitrogen (BUN)	●	●
Cholesterol (HDL)	●	●	Lactate Dehydrogenase (LDH)	●	●	Uric Acid	●	●
Cholesterol (LDL)	●	●	LAP Arylamidase	●		Valproic Acid	●	●
Cholesterol (Total)	●	●	Lipase	●	●	Vitamin B ₁₂	●	●
Cholinesterase	●	●	Lithium	●	●	Zinc	●	
CK-MB Isoenzyme*	●		Magnesium	●	●			

*No claim is made for expected value or stability.

**α-1-Globulin, α-2-Globulin, β-Globulin, γ-Globulin.

Lyphocek Assayed Chemistry Control and Lyphocek Unassayed Chemistry Control (Human)



- Includes a comprehensive list of claimed analytes for general chemistry testing
- Provides good stability for enzymes and CO₂
- No special diluents are required for reconstitution
- Human serum based
- Assayed: 40 month shelf life at 2-8°C
Unassayed: 37 month shelf life at 2-8°C
- 30 day reconstituted and frozen stability at -10°C to -20°C for most analytes
- 7 day open-vial reconstituted stability at 2-8°C for most analytes

POCT methods listed for
Lyphocek Assayed Chemistry Control

Ordering Information

Cat #	Description	
Assayed Chemistry Control		
C-310-5	Level 1	12 x 5 mL
C-315-5	Level 2	12 x 5 mL
313X	Bilevel MiniPak (1 of each level)	2 x 5 mL
Unassayed Chemistry Control (Human)		
731	Level 1	.25 x 5 mL
732	Level 2	.25 x 5 mL
730X	Bilevel MiniPak (1 of each level)	2 x 5 mL
C-320-10	Level 1	.50 x 10 mL
C-325-10	Level 2	.50 x 10 mL
323X	Bilevel MiniPak (1 of each level)	2 x 10 mL

Analytes

Acetaminophen	Ceruloplasmin	Immunoglobulin G (IgG)	T3 (Total)
Acid Phosphatase (Total)	Chloride	Immunoglobulin M (IgM)	T3 Uptake/T-Uptake
Albumin	Cholesterol (HDL)	Iron	T4 (Free)
Alkaline Phosphatase (ALP)	Cholesterol (LDL)	Iron (TIBC)	T4 (Total)
Alpha-1-Antitrypsin	Cholesterol (Total)	Iron (UIBC)	TBG
Alpha-Hydroxybutyrate Dehydrogenase (αHBDH)*	Cholinesterase	Lactate (Lactic Acid)	Theophylline
Alphafetoprotein (AFP)	CO ₂	Lactate Dehydrogenase (LDH)	Tobramycin
ALT/SGPT	Complement C3	LAP Arylamidase	Transferrin
Amylase	Complement C4*	Lipase	Triglycerides
Amylase (Alpha)	Copper	Lithium	TSH
Amylase (Pancreatic)	Cortisol	Magnesium	Urea
Apolipoprotein A-1	Creatine Kinase (CK)	Osmolality	Urea Nitrogen (BUN)
Apolipoprotein B	Creatinine	PAP	Uric Acid
AST/SGOT	Digoxin	Phenobarbital	Valproic Acid
Bilirubin (Direct)	Gentamicin	Phenytoin	Vancomycin
Bilirubin (Indirect)	GGT	Phosphorus	Vitamin B ₁₂
Bilirubin (Total)	GLDH	Potassium	Zinc
Calcium	Globulin	Protein Serum (Total)	
Calcium (Ionized)	Glucose	PSA (Total) ¹	
Carbamazepine	Haptoglobin	Salicylate	
CEA	hCG-β Subunit	Sodium	
	Immunoglobulin A (IgA)	T3 (Free)	

¹Available in Lyphocek Assayed Chemistry Control only.

*No claim is made for expected value or stability.

Liquicheck Unassayed Chemistry Control (Human)



- Unassayed control with larger 10 mL fill size for more cost effective serum chemistry testing
- Liquid, human serum based
- 2 year shelf life at -20°C to -70°C
- 15 day open-vial stability at 2-8°C for most analytes

Ordering Information

Cat #	Description
691	Level 1 25 x 10 mL
692	Level 2 25 x 10 mL
690X	Bilevel MiniPak (1 of each level) 2 x 10 mL

Analytes

Acetaminophen	Cholesterol (Total)	Iron (UIBC)*	Sodium
Acid Phosphatase (Total)	CO ₂	Lactate (Lactic Acid)	T3 (Free)*
Albumin	Copper	Lactate Dehydrogenase (LDH)	T3 (Total)
Aldolase*	Cortisol	LAP Arylamidase*	T3 Uptake/T-Uptake
Alkaline Phosphatase (ALP)	Creatine Kinase (CK)	Lipase	T4 (Free)*
Alpha-Hydroxybutyrate Dehydrogenase (αHBDH)*	Creatinine	Lithium	T4 (Total)
ALT/SGPT	Digoxin	Magnesium	Theophylline
Amylase	Ethanol	Osmolality	Tobramycin
AST/SGOT	Gentamicin	PAP*	Transferrin
Bilirubin (Direct)*	GGT	Phenobarbital	Triglycerides
Bilirubin (Indirect)*	GLDH*	Phenytoin	TSH
Bilirubin (Total)	Globulin	Phospholipids	Urea Nitrogen (BUN)
Calcium	Glucose	Phosphorus	Uric Acid
Calcium (Ionized)*	Immunoglobulin A (IgA)	Potassium	Valproic Acid
Carbamazepine	Immunoglobulin G (IgG)	Protein (Total)	Vancomycin
Chloride	Immunoglobulin M (IgM)	Protein Electrophoresis	Vitamin B ₁₂
Cholesterol (HDL)*	Iron	Pseudocholesterase	Zinc*
	Iron (TIBC)	Salicylate	

*No claims are made regarding performance or stability.

Liquicheck Ethanol/Ammonia Control



Available in Multiple Configurations

Offers normal, elevated and toxic concentrations of Ethanol and Ammonia.

Standard Vials

- 2 year shelf life at 2-8°C
- 20 day open-vial stability at 2-8°C

Compact Vials

- 20 day open-vial stability for both analytes when stored at 2-8°C
- 2 year shelf life when stored at 2-8°C

Ordering Information

Cat #	Description
Standard Vials	
544	Level 16 x 3 mL
545	Level 26 x 3 mL
546	Level 36 x 3 mL
545X	Trilevel MiniPak (1 of each level)3 x 3 mL
Compact Vials	
271	Level 16 x 2.5 mL
272	Level 26 x 2.5 mL
273	Level 36 x 2.5 mL
272X	Trilevel MiniPak (1 of each level)3 x 2.5 mL

For IntelliQ Ethanol/Ammonia Control, see page 12

Analytes

Ammonia
Ethanol

Liquichek Lipids Control



- Contains C-Reactive Protein (CRP) in levels optimized for hs-CRP assays
- Higher Triglycerides and Cholesterol values to monitor precision of the full assay range
- Provides values for direct HDL
- Liquid, human serum based
- 3 year shelf life at -20°C to -70°C
- 14 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description
641	Level 1 6 x 3 mL
642	Level 2 6 x 3 mL
640X	Bilevel MiniPak (1 of each level) 2 x 3 mL

Analytes

Apolipoprotein A-1	C-Reactive Protein (CRP)	Cholesterol (LDL)	Lipoprotein (a)
Apolipoprotein B	Cholesterol (HDL)	Cholesterol (Total)	Triglycerides

Liquichek Pediatric Control



Available in Multiple Configurations

- Suitable for monitoring high Bilirubin levels important for adult liver disease
- Liquid

Standard Vials

- Four Bilirubin parameters are available at two clinically significant levels for neonatal testing
- 3 year shelf life at -20°C to -70°C
- 3 month stability when stored unopened at 2-8°C
- 14 day open-vial stability for most analytes at 2-8°C

Compact Vials

- Monitors precision of analyte test procedures at higher concentrations important for neonatal testing
- 7 day open-vial stability for all analytes when stored at 2-8°C
- 3 year shelf life when stored at -20°C to -50°C

Ordering Information

Cat #	Description
Standard Vials	
354	Level 1 6 x 4 mL
355	Level 2 6 x 4 mL
353X	Bilevel MiniPak (1 of each level) 2 x 4 mL
Compact Vials	
176	Level 1 6 x 2.5 mL
177	Level 2 6 x 2.5 mL
178X	Bilevel MiniPak (1 of each level) 2 x 2.5 mL

Analytes

Bilirubin (Direct)*	Caffeine	Magnesium	Sodium
Bilirubin (Indirect)**	Calcium	Phenylalanine*	Theophylline
Bilirubin (Neonatal)**	Chloride	Potassium	
Bilirubin (Total)*	Glucose		*Available in Standard Vials only **7 day open-vial stability at 2-8°C

Refer to myInserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Liquid Assayed Multiquel Premium



- Combination of 95 analytes, including General Chemistry, Immunology and Therapeutic Drug Monitoring (TDM) analytes
- Introduces a new level of efficiency by offering analytes from multiple high-usage quality controls
- Liquid, human serum based
- 3 year shelf life at -20°C to -70°C
- 14 day frozen aliquot stability at 2-8°C for most analytes
- 30 day unopened stability at 2-8°C for most analytes

Ordering Information

Cat #	Description
284	Level 16 x 5 mL
285	Level 26 x 5 mL
286	Level 36 x 5 mL
285X	Trilevel MiniPak (1 of each level)3 x 5 mL

Analytes

General Chemistry

Acid Phosphatase (Total)
Albumin
Alkaline Phosphatase (ALP)
Alpha-Hydroxybutyrate
Dehydrogenase (α HBDH)¹
ALT/SGPT
Amylase
Amylase (Pancreatic)
AST/SGOT
Bile Acids
Bilirubin (Direct)
Bilirubin (Neonatal)
Bilirubin (Total)
Calcium
Calcium (Ionized)
Carbon Dioxide (CO₂)
Chloride
Cholesterol (HDL)
Cholesterol (LDL)
Cholesterol (Total)
Cholinesterase¹
Copper¹
Creatine Kinase (CK)
Creatinine
Ethanol
GGT

Glucose
Glutamate Dehydrogenase
(GLDH)**
Iron
Iron (TIBC)
Iron (UIBC)
Lactate (Lactic Acid)
Lactate Dehydrogenase (LDH)
LAP Arylamidase
Lipase
Magnesium
Osmolality
Phosphorus
Phospholipids
Potassium
Protein Serum (Total)
Sodium
Triglycerides
Urea
Urea Nitrogen (BUN)
Uric Acid
Zinc

Immunology

Alpha-1-Antitrypsin¹
Alpha-1-Acid Glycoprotein
(AAG)¹
Anti-CCP¹

ASO¹
Apolipoprotein A-1¹
Apolipoprotein B¹
 β -2-Microglobulin¹
Ceruloplasmin¹
C-Reactive Protein (CRP)
Complement C3¹
Complement C4¹
Cystatin C¹
Globulin¹
Haptoglobin¹
Immunoglobulin A (IgA)¹
Immunoglobulin G (IgG)¹
Immunoglobulin M (IgM)¹
Kappa Light Chain¹
Lambda Light Chain¹
Prealbumin¹
Transferrin¹

TDM

Acetaminophen
Amikacin
Caffeine
Carbamazepine
Digoxin
Disopyramide
Gentamicin
Lidocaine

Lithium
Methotrexate
NAPA
Phenobarbital
Phenytoin
Procainamide
Quinidine
Salicylate
Theophylline
Tobramycin
Valproic Acid
Vancomycin

Immunoassay

Cortisol
Ferritin¹
T3 Uptake / T Uptake
T3 (Total)
T4 (Free)
T4 (Total)
TSH
Vitamin B₁₂¹

*No claims are made regarding performance or stability.

**Not available for diagnostic use in the United States.

¹Endogenous levels only.

Liquichek Microalbumin Control



- Designed using human Albumin at levels found in early stage renal disease
- Values provided for a wide range of popular instrument platforms
- Liquid, human urine based
- 2 year shelf life at 2-8°C
- 90 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description
378	Level 112 x 10 mL
379	Level 212 x 10 mL
374X	Bilevel MiniPak (1 of each level)2 x 10 mL

Analytes

Creatinine
Microalbumin

Liquichek Urine Chemistry Control



Available in Multiple Configurations

- Monitors the precision of general urine chemistry testing procedures
- Includes Amylase and Creatinine
- Liquid, human urine based

Standard Vials

- 2 year shelf life at 2-8°C
- 30 day open-vial stability at 2-8°C

Compact Vials

- 30 day open-vial stability for all analytes when stored at 2-8°C
- 2 year shelf life when stored refrigerated at 2-8°C

POCT methods listed for standard vials

Ordering Information

Cat #	Description
Standard Vials	
397	Level 1 12 x 10 mL
398	Level 2 12 x 10 mL
395X	Bilevel MiniPak (1 of each level) 2 x 10 mL
Compact Vials	
195	Level 1 12 x 2.5 mL
196	Level 2 12 x 2.5 mL
198X	Bilevel MiniPak (1 of each level) 2 x 2.5 mL

For InteliQ Urine Chemistry Control, see page 14

Analytes

Amylase	Glucose	Phosphorus	Specific Gravity*
Calcium	Magnesium	Potassium	Urea*
Chloride	Microalbumin	Pregnancy/hCG (Qualitative)*	Urea Nitrogen (BUN)
Cortisol*	Osmolality*	Protein (Total)	Uric Acid
Creatinine	pH*	Sodium	

*Available in standard vial configuration only.

Lypochek Quantitative Urine Control



- Comprehensive analyte menu for quantitative assays that monitor the precision for general urine chemistry, hormones and esoteric testing including biogenic amines, metals and porphyrins
- Contains hCG
- Microalbumin available at levels found in early stages of proteinuria and associated nephropathy
- Value assigned for a comprehensive list of methods
- Human urine based
- 2 year shelf life at 2-8°C for most analytes
- 5 day reconstituted stability at 2-8°C for most analytes

Ordering Information

Cat #	Description
376	Level 1, Normal 12 x 10 mL
377	Level 2, Abnormal 12 x 10 mL
375X	Bilevel MiniPak (1 of each level) 2 x 10 mL

Analytes

3-Methoxytyramine*	Calcium	Hydroxyproline (Free)*	Potassium
5-Aminolevulinic Acid (ALA)	Chloride	Hydroxyproline (Total)*	Pregnancy (Qualitative)
5-HIAA	Copper*	Iron	Pregnanetriol*
11-β-Hydroxy-Androsterone*	Coproporphyrin	Lead	Protein (Total)
11-β-Hydroxy-Etiocholanolone*	Cortisol	Magnesium	Sodium
17-Hydroxycorticosteroids*	Creatine*	Mercury	Specific Gravity
11-Ketoandrosterone*	Creatinine	Metanephrine	Urea
11-Ketoetiocholanolone*	Cystine*	Metanephrine (Total)	Urea Nitrogen
17-Ketosteroids	Dehydroepiandrosterone (DHEA)*	Microalbumin	Uric Acid
17-Ketogenic Steroids	Dopamine	Norepinephrine	Uroporphyrins
Aldosterone	Epinephrine	Normetanephrine	Vanillylmandelic Acid (VMA)
AMP (Cyclic)*	Estriol (Total)*	Osmolality	Zinc*
Androsterone*	Etiocholanolone*	pH	
Arsenic	Glucose	Phosphorus	
Barbiturates*	Homovanillic Acid (HVA)	Porphyrins (Total)	

*No claims are made regarding performance or stability.

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Liquichek Serum Indices



- Intended for use as a pre-analytical step to determine if an instrument is detecting Hemolyzed, Icteric, or Lipemic (HIL) samples
- Prepared from human source material
- 2 year shelf life at -20°C to -70°C
- 14 day open-vial stability at 2-8°C
- 14 day closed-vial stability at 2-8°C
- 28 day frozen aliquot at -20°C to -70°C

Ordering Information

Cat #	Description
12012693	Hemolysis6 x 4 mL
12012694	Icterus6 x 4 mL
12012695	Lipemia6 x 4 mL
12012696	Non-Interfered6 x 4 mL
12012697	Hemolysis MiniPak.1 x 4 mL
12012698	Icterus MiniPak1 x 4 mL
12012699	Lipemia MiniPak1 x 4 mL
12012700	Non-Interfered MiniPak1 x 4 mL

Interferents

- Hemolysis
- Icterus
- Lipemia
- Non-Interfered

Lyphocek Coagulation Control



- Extended open-vial stability reduces the cost of daily QC
- Three year shelf life allows laboratories to save money by performing fewer crossover studies
- Human based
- 3 year shelf life at 2-8°C
- 48 hour open-vial stability at 2-25°C

Ordering Information

Cat #	Description
744	Level 1 12 x 1 mL
745	Level 2 12 x 1 mL
746	Level 3 12 x 1 mL
745X	Trilevel MiniPak (1 of each level) 3 x 1 mL

Analytes

Activated Partial Thromboplastin Time (APTT)	Fibrinogen
Antithrombin III (AT III)	Prothrombin Time (PT)
	Thrombin Time (TT)

Lyphocek Hemostasis Control



- Supplements basic parameters provided by Lyphocek Coagulation Control with additional factors
- Provides QC for more specialized diagnostic testing
- Human based
- 3 year shelf life at 2-8°C
- 8 hour open-vial stability at 2-25°C, except for Protein S at 2-8°C

Ordering Information

Cat #	Description
597	Level 1 12 x 1 mL
598	Level 2 12 x 1 mL
599	Level 3 12 x 1 mL
598X	Trilevel MiniPak (1 of each level) 3 x 1 mL

Analytes

Activated Partial Thromboplastin Time (APTT)	Protein C (Functional)
Antithrombin III (AT III)	Protein S (Functional)
Factors II, V, VII, VIII, IX, X, XI and XII	Prothrombin Time (PT)
Fibrinogen	Thrombin Time (TT)
Plasminogen	

Liquichek D-dimer Control



- Four level control monitoring assay precision across the critical range used in diagnosis of thrombotic events on multiple platforms
- Liquid, human-plasma based
- 40 month shelf life at -20°C to -70°C
- 15 day open-vial stability at 2-8°C

POCT methods listed

Ordering Information

Cat #	Description
27100	Level Low. 6 x 1 mL
27101	Level 1 6 x 1 mL
27102	Level 2 6 x 1 mL
27103	Level 3 6 x 1 mL
27102X	Four Level MiniPak (1 of each level) 4 x 1 mL

Analyte

D-dimer

Liquichek Diabetes Control



Available in Multiple Configurations

- Monitors the precision of test methods used for Hemoglobin A1C testing
- Liquid, human whole blood based

Standard Vials

- Higher third level HbA1c to monitor your assay range
- 3 year shelf life at -20°C to -70°C
- 45 day open-vial stability at 2-8°C

Compact Vials

- Compatible with applicable automated platforms such as Siemens Dimension Vista and Abbott Alinity
- 2D barcode for quick and easy QC set-up
- 3 year shelf life at -20°C to -50°C
- 30 day open-vial stability at 2-8°C

POCT methods listed for standard vials



Ordering Information

Cat #	Description
Standard Vials	
12011343	Level 16 x 1 mL
12011344	Level 26 x 1 mL
12011345	Level 36 x 1 mL
12011334	Trilevel MiniPak (1 of each level)3 x 1 mL
Compact Vials	
12011308	Level 16 x 2 mL
12011309	Level 26 x 2 mL
12011310	Level 36 x 2 mL
12011331	Trilevel MiniPak (1 of each level)3 x 2 mL

For IntelliQ Diabetes Control, see page 12

Analyte

Hemoglobin A1C

Lyphocheck Diabetes Control



- Lyophilized control available in two levels
- Human whole blood based control
- 3 year shelf life at 2-8°C
- 7 day reconstituted stability at 2-8°C

POCT methods listed

Ordering Information

Cat #	Description
740	Bilevel (3 of each level) 6 x 0.5 mL
740X	Bilevel MiniPak (1 of each level) 2 x 0.5 mL

Analytes

Hemoglobin (Total Glycated)*	Hemoglobin A1C
Hemoglobin (Total)	Hemoglobin F
Hemoglobin A1*	

*No claims are made regarding performance or stability.

Lyphocheck Hemoglobin A₂ Control



- Lyophilized
- Human whole blood based control
- 2 year shelf life at 2-8°C
- 21 day reconstituted stability at 2-8°C for most methods
- 14 day reconstituted stability at 2-8°C for Gel Electrophoresis

Assayed values available for common analyzers:

Bio-Rad

- D-10
- VARIANT™
- VARIANT™ II

Other

- SEBIA Hydragel
- SEBIA CAPILLARYS 2 FLEX PIERCING
- TOSOH G7

Ordering Information

Cat #	Description
553	Bilevel (2 of each level) 4 x 0.5 mL
553X	Bilevel MiniPak (1 of each level) 2 x 0.5 mL

Analytes

Hemoglobin A₂
Hemoglobin F
Hemoglobin S

Lymphochek Hemoglobin A1C Linearity Set



- Intended to verify linearity throughout the reportable range of Hemoglobin A1C assays
- Human whole blood based
- 3 year shelf life at 2-8°C
- 7 day reconstituted stability at 2-8°C

Linearity Web software available on QCNet.com

Ordering Information

Cat #	Description	Quantity
12000070	Set (1 of each level)	6 x 0.5 mL

Analyte

Hemoglobin A1C

Assayed values available for common analyzers:

Bio-Rad Analyzers

- D-10
- VARIANT™
- VARIANT™ II Turbo

Other Analyzers

- Beckman Coulter SYNCHRON Systems
- Siemens ADVIA
- Siemens Dimension Series
- Roche COBAS INTEGRA & Hitachi Systems
- Tosoh G7 & G8

Expected Values for Hemoglobin A1C (%NGSP)*

Level	Expected Value Range
Level 1	2.7 - 3.8%
Level 2	4.1 - 5.1%
Level 3	5.5 - 6.5%
Level 4	8.4 - 10%
Level 5	12 - 15%
Level 6	16 - 22%

*Ranges reflect typical recovery values for this product. Exact values may vary depending on test/assay methodology, instruments, reagent, system calibrators or laboratory technique.

Meter Trax Control



- Easy-to-use control for Glucose, Hemoglobin and Hematocrit testing on point-of-care instruments
- Glucose and Hemoglobin stable for 31 days at 2–25°C after opening
- Hematocrit stable for 31 days at 2-8°C after opening
- 6 month shelf life at 2-8°C

POCT methods listed

Assayed values available for common analyzers:

- | | |
|---|---|
| <ul style="list-style-type: none"> Abbott iSTAT BD Clay Adams Beckman StatSpin STI HemataSTAT HemoCue Hb | <ul style="list-style-type: none"> i-SENS CareSens LifeScan OneTouch Nova StatStrip Roche Accu-Chek |
|---|---|

Ordering Information

Cat #	Description	Quantity
970	Trilevel (2 of each level)	6 x 2 mL
971	Low Level	6 x 2 mL
972	Mid Level	6 x 2 mL
973	High Level	6 x 2 mL
970X	Trilevel MiniPak (1 of each level)	3 x 2 mL

Analytes

Glucose
Hematocrit (HCT)
Hemoglobin (Total)

Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Blood Gas Program



- Liquid, aqueous samples
- Comprehensive menu for external assessment of laboratory test performance of Blood Gas, Electrolytes, Glucose, Lactate and Magnesium
- 12-month cycle

Ordering Information

Catalog #
 BC31* 12 x 2.5 mL

Analytes

Calcium (Ionized)	pCO ₂
Chloride	pH
Glucose	pO ₂
Lactate (Lactic Acid)	Potassium
Magnesium	Sodium

Blood Typing Program



- Ideal for clinical lab, blood bank and transfusion center testing needs
- Barcoded and ready-to-use tubes for automated platforms
- 3 samples tested every 4 months
- Educational case studies with each shipment

Ordering Information

Catalog #	Description	Sample	Donor
12000837*	Shipment A	3 x 4 mL	+ 1 x 2 mL
12000838*	Shipment B	3 x 4 mL	+ 1 x 2 mL
12000839*	Shipment C	3 x 4 mL	+ 1 x 2 mL

Parameters

ABO Group	Compatibility (Cross-match)
Antibody Identification	DAT (Direct Antiglobulin Test)
Antibody Screen	Rh(D)
Antigen Typing	

Cardiac Markers Program



- Liquid, human plasma based
- Comprehensive menu of important cardiac assessment markers, plus D-dimer
- 12-month cycle

Ordering Information

Catalog #
 BC39* 12 x 1.5 mL

Analytes

BNP	Homocysteine	Myoglobin	Troponin I
CK-MB Isoenzyme (mass)	hs-CRP	NT-proBNP	Troponin T
D-dimer			

Coagulation Program



- Lyophilized, human plasma based
- 11 parameters for external assessment of laboratory coagulation test performance
- 12-month cycle

Ordering Information

Catalog #
 BC34* 12 x 1 mL

Analytes

Antithrombin III (ATIII)	Factor VIII	Protein C	Thrombin Time (TT)
APTT	Fibrinogen	Protein S	Von Willebrand Factor (vWF)
D-dimer	INR	PT	

*Aspects of the program are subcontracted from an approved supplier. Refer to myInSerts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Clinical Chemistry (Monthly) Program



- Lyophilized, human serum based
- 12-month cycle
- Submit results for the appropriate sample every month
- Reference Method Values provided for select analytes

Ordering Information

Catalog #
BC50 12 x 5 mL

Analytes

Acid Phosphatase (Total)	Calcium (Total)	Iron	Sodium
Alanine Aminotransferase (ALT/SGPT)	Carbon Dioxide (CO ₂) Chloride	Iron Binding Capacity, Total (TIBC)	Thyroid Stimulating Hormone (TSH)
Albumin	Cholesterol (HDL)	Iron Binding Capacity, Unsaturated (UIBC)	Thyroxine (T4), Free
Alkaline Phosphatase (ALP)	Cholesterol (LDL)	Lactate (Lactic Acid)	Thyroxine (T4), Total
Amylase (Pancreatic)	Cholesterol (Total)	Lactate Dehydrogenase (LDH)	Transferrin
Amylase (Total)	Cholinesterase	Lipase	Triglycerides
Aspartate Aminotransferase (AST/SGOT)	Copper	Lithium	Triiodothyronine (T3), Free
Bile Acids	Cortisol	Magnesium	Triiodothyronine (T3), Total
Bilirubin (Direct)	Creatine Kinase (CK)	Osmolality	Urea
Bilirubin (Indirect)	Creatinine	Phosphorus	Urea Nitrogen (BUN)
Bilirubin (Total)	Gamma Glutamyl Transferase (GGT)	Potassium	Uric Acid
Calcium (Ionized)	Glucose	Protein (Total)	Zinc

Ethanol/Ammonia Program



- Liquid, human serum based
- Clinically relevant levels of serum Ethanol and Ammonia
- 12-month cycle

Ordering Information

Catalog #
BC35* 12 x 3 mL

Analytes

Ammonia
Ethanol

Hematology Program



- Liquid product with human RBCs
- Choice of 11 basic hematology parameters
- For use on manual or automated analyzers
- Convenience of primary pierceable tubes
- 12-month cycle consisting of 4 separate shipments
- Submit results for the appropriate sample every month

Ordering Information

Catalog #	Description	
BC90A*	Shipment A 3 x 2 mL
BC90B*	Shipment B 3 x 2 mL
BC90C*	Shipment C 3 x 2 mL
BC90D*	Shipment D 3 x 2 mL

Parameters

Hematocrit (HCT)	Platelets (PLT)
Hemoglobin (HGB)	Red Blood Cells (RBC)
Mean Corpuscular Hemoglobin (MCH)	Red Blood Cell Distribution Width (RDW)
Mean Corpuscular Hemoglobin Concentration (MCHC)	Red Blood Cell Distribution Width-SD (RDW-SD)
Mean Corpuscular Volume (MCV)	White Blood Cells (WBC)
Mean Platelet Volume (MPV)	

*Aspects of the program are subcontracted from an approved supplier.

Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Hemoglobin Program



- Lyophilized, human whole blood based
- Includes HbA1C and HbA₂
- 12-month cycle
- Submit results for the appropriate sample every month
- **Reference Method Values** provided for Hemoglobin A1C

Ordering Information

Catalog #
BC80 12 x 0.5 mL

Analytes

Hemoglobin A1C Hemoglobin (Total Glycated)
Hemoglobin A₂

HIV/Hepatitis Program



- Liquid, human source material
- Suitable for serological tests used to detect key HIV and Hepatitis viral antibodies and antigens
- Qualitative and quantitative results reporting available
- Submit results and view reports each month through EQAS Online

Ordering Information

Catalog #
12000815 12 x 2 mL

Analytes

Anti-HAV	Anti-HCV	Anti-HTLV-I	HBeAg
Anti-HBc	Anti-HIV-1	HAV IgG	HBsAg
Anti-HBe	Anti-HIV-1/2	HAV IgM	HIV-1 Ag
Anti-HBs	Anti-HIV-2	HbC IgM	

Immunoassay (Monthly) Program



- Lyophilized, human serum based
- Comprehensive menu of the most commonly tested immunoassay analytes
- 12-month cycle
- Submit results for the appropriate sample every month

Ordering Information

Catalog #
BC75 12 x 5 mL

Analytes

11-Deoxycortisol	DHEA	NSE	Theophylline
17- α -OH-Progesterone	DHEA Sulfate	Phenobarbital	Thyroglobulin (Tg)
25-Hydroxy Vitamin D	Digoxin	Phenytoin	Thyroid Stimulating Hormone (TSH)
β -2 Microglobulin	Estradiol	Plasma Renin Activity	Thyroxine Binding Globulin (TBG)
ACE	Estradiol, Free (UE ₃)	Progesterone	Thyroxine, Free (FT ₄)
ACTH	Ferritin	Prolactin	Thyroxine (T4), Total
Aldosterone	Folate	PSA (Free)	Transferrin
Alphafetoprotein (AFP)	Follicle Stimulating Hormone (FSH)	PSA (Free/Total Ratio)	Triiodothyronine, Free (FT ₃)
Androstenedione	Fructosamine	PSA (Total)	Triiodothyronine (T3), Total
C-Peptide	Gastrin	Renin	Valproic Acid
CA 15-3	hCG	S100 Protein	Vitamin B ₁₂
CA 19-9	hGH	Sex Hormone Binding Globulin (SHBG)	<i>Pilot Analyte</i>
CA 27.29	Immunoglobulin E (IgE)	T3 Uptake	Calcitonin
CA 125	Insulin	T-Uptake	
Carbamazepine	Intact PTH	Testosterone	
Carcinoembryonic Antigen (CEA)	LH	Testosterone (Free)	
Cortisol			

Lipids Program



- Liquid, human serum based
- Comprehensive menu of lipid profile analytes
- 12-month cycle
- Submit results for the appropriate sample every month

Ordering Information

Catalog #
BC47* 12 x 3 mL

Analytes

Apolipoprotein A-1	Cholesterol (HDL)	Cholesterol (Total)	Triglycerides
Apolipoprotein B	Cholesterol (LDL)	Lipoprotein (a)	

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Serum Proteins Program



- Lyophilized, human serum based
- Comprehensive menu for external assessment of serum protein assays
- 12-month cycle
- Submit results for the appropriate sample every month

Ordering Information

Catalog #
BC23 12 x 1 mL

Analytes

α-1 Acid Glycoprotein	C-Reactive Protein (CRP)	Immunoglobulin E (IgE)	Protein (Total)
α-1 Antitrypsin	Ceruloplasmin	Immunoglobulin G (IgG)	Rheumatoid Factor (RF)
α-2 Macroglobulin	Complement C3	Immunoglobulin M (IgM)	Transferrin
β-2 Microglobulin	Complement C4	Kappa Light Chain	
Albumin	Haptoglobin	Lambda Light Chain	
Antistreptolysin O (ASO)	Immunoglobulin A (IgA)	Prealbumin	

NEW | Specialty Immunoassay Program



- Liquid, human serum based
- Comprehensive menu of complex and highly tested specialty immunoassay analytes
- 12-month cycle
- Submit results for the appropriate sample every month

Ordering Information

Cat # Description
12015696 12 x 3 mL

Analytes

25-Hydroxy Vitamin D	C-Peptide	Interleukin-6 (IL-6)*	TSH Receptor Antibody (TRAb)
Active Vitamin B ₁₂	Erythropoietin (EPO)	Osteocalcin	
Anti-Thyroglobulin (Anti-Tg)	Fructosamine	Procalcitonin (PCT)	
Anti-Thyroperoxidase (Anti-TPO)	Insulin Growth Factor I (IGF-I)	PTH (Intact)	*This analyte is not available for diagnostic use in the United States.

Syphilis/Chagas Program



- Liquid, human source material
- Suitable for serological tests detecting *Treponema pallidum*, Non-Treponemal and *Trypanosoma cruzi* antibodies
- 12-month cycle
- Qualitative and quantitative results reporting available
- Submit results and view reports each month through EQAS Online

Ordering Information

Catalog #
12000833 12 x 1.5 mL

Analytes

Non-Treponemal Antibodies	Anti- <i>Treponema pallidum</i> (Total)
<i>Treponema pallidum</i> IgG	<i>Trypanosoma cruzi</i> Antibodies

Therapeutic Drug Monitoring Program (TDM)



- Lyophilized, human serum based
- Analytes targeted at therapeutic, sub-therapeutic and toxic levels
- 12-month cycle
- Submit results for the appropriate sample every month
- **Weighed-in Values** provided for select analytes at end of cycle

Ordering Information

Catalog #
BC10 12 x 5 mL

Analytes

Acetaminophen	Digoxin	Phenobarbital	Tricyclic Antidepressants (TCA)
Amikacin	Ethosuximide	Phenytoin	Screen
Amitriptyline	Gentamicin	Primidone	Valproic Acid
Caffeine	Lithium	Salicylate	Vancomycin
Carbamazepine	Methotrexate	Theophylline	
Cyclosporine	Nortriptyline	Tobramycin	

*Aspects of the program are subcontracted from an approved supplier. Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

ToRCH/EBV/MuMZ Program



- Liquid, human source material
- Suitable for serological tests used to detect pediatric diseases and congenital infections
- Qualitative and quantitative results reporting available
- Submit results and view reports each month through EQAS Online

Ordering Information

Catalog #
12000825 12 x 2 mL

Analytes

Cytomegalovirus (CMV) IgG
Cytomegalovirus (CMV) IgM
Epstein-Barr Virus (EBNA) IgG
Epstein-Barr Virus (VCA) IgG
Epstein-Barr Virus (VCA) IgM
Herpes Simplex Virus (HSV-1) IgG
Herpes Simplex Virus (HSV-1/2) IgG
Herpes Simplex Virus (HSV-2) IgG

Measles Virus (Rubeola) IgG
Mumps Virus IgG
Rubella Virus IgG
Rubella Virus IgM
Toxoplasma gondii (Toxo) IgG
Toxoplasma gondii (Toxo) IgM
Varicella Zoster Virus (VZV) IgG

Urinalysis Program



- Liquid matrix
- Suitable for manual and automated dipstick analysis in both clinical and point-of-care testing (POCT)
- Submit qualitative results and view reports each month through EQAS Online

Ordering Information

Catalog #
12000810 12 x 12 mL

Analytes

Albumin
Albumin-to-Creatinine Ratio
Bilirubin
Blood/Hemoglobin
Creatinine
Glucose
Ketones
Leukocytes

Nitrite
pH
Pregnancy (hCG)
Protein (Total)
Protein-to-Creatinine Ratio
Specific Gravity
Urobilinogen

Urine Chemistry Program



- Lyophilized, human urine based
- Matrix appropriate samples include basic and specialty urine chemistry analytes
- 12-month cycle
- Submit results for the appropriate sample every month

Ordering Information

Catalog #
BC45 12 x 10 mL

Analytes

5-Hydroxyindoleacetic Acid (5-HIAA)
Calcium (Total)
Chloride
Cortisol (Free)
Creatinine
Dopamine

Epinephrine
Glucose
Homovanillic Acid (HVA)
Magnesium
Metanephrine
Microalbumin (Albumin)
Norepinephrine

Normetanephrine
Osmolality
Phosphorus
Potassium
Protein (Total)
Sodium

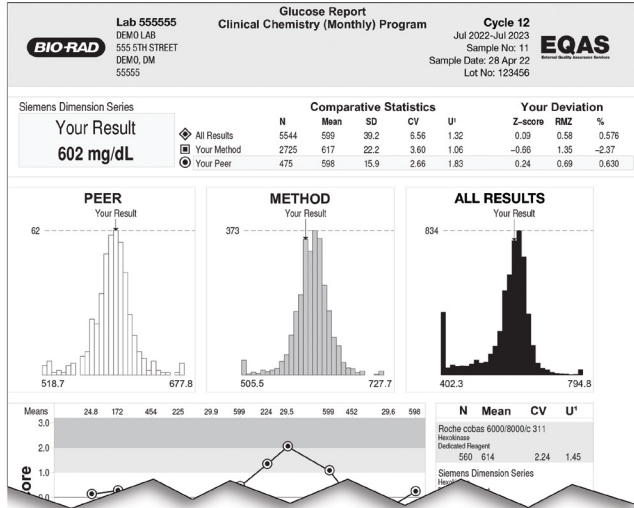
Urea Nitrogen (BUN)
Uric Acid
Vanillylmandelic Acid (VMA)
Pilot Analyte
Urea

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

EQAS™ Reports

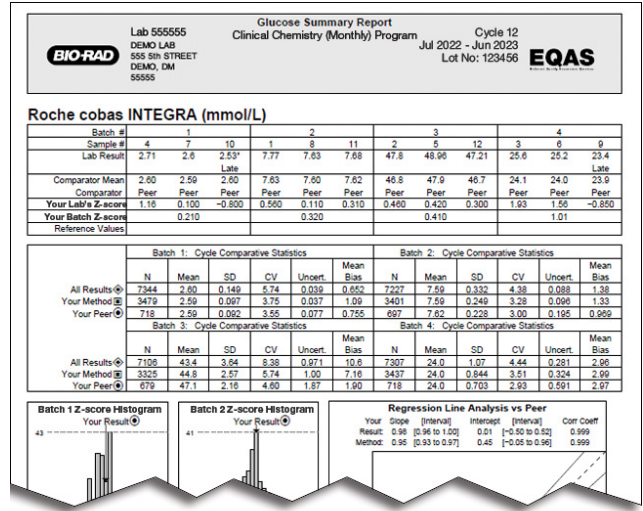
Participants receive twelve sample reports and one end-of-cycle report over each program cycle. These detailed reports are presented in an easy-to-read graphic format. Sample reports are available online within three days of the statistical analysis. Additional special reports (Quality Specification, Subgroup and Instrument/Method Summary) are available to registered participants.

Quantitative Reports



Sample Report

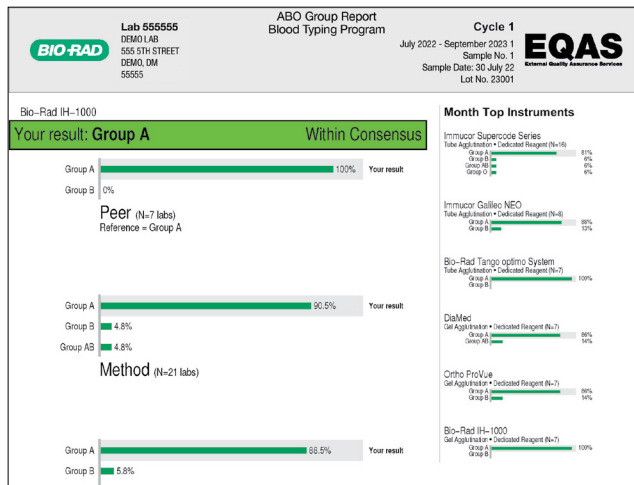
Detailed individual statistical analyses and peer group comparisons are provided.



End-of-Cycle Report

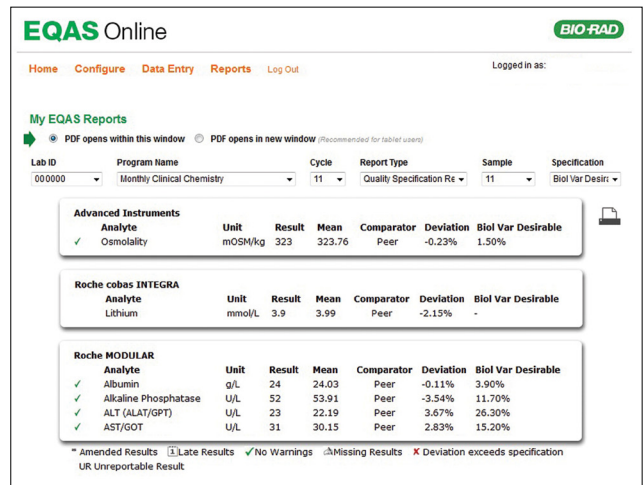
An overall assessment of laboratory performance over the entire cycle in terms of bias among participating laboratories is presented.

Qualitative Report



Sample Report

These reports provide clear and concise information about your qualitative test laboratory performance in the form of histograms and modified Levey-Jennings charts. They include detailed information about the distribution of the qualitative responses and the peer group consensus.



Quality Specification Report

An evaluation, in terms of percentage deviation, against set performance criteria from key regulatory and scientific organizations (ex. CLIA, BV, RCPA, IQMH, GOST and RiilBÄK), is available online for each sample analyte.

Refer to myInsersts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Liquichek Hematology Control (A)

For Abbott Systems



- Provides assayed values for the Abbott CELL-DYN 3200, Ruby and Sapphire analyzers
- Blood cells in plasma-like fluid
- 78 day shelf life at 2-8°C
- 8 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description
865	Trilevel (4 of each level) 12 x 3 mL
866	Low 4 x 3 mL
867	Normal. 4 x 3 mL
868	High 4 x 3 mL
865X	Trilevel MiniPak (1 of each level) 3 x 3 mL

Parameters

Basophils (BASO)
 % Basophils
 Eosinophils (EOS)
 % Eosinophils
 Granulocytes (GRAN)
 Hematocrit (HCT)
 Hemoglobin (Total)
 Lymphocytes (LYMPH)
 % Lymphocytes
 Mean Corpuscular Hemoglobin (MCH)
 Mean Corpuscular Hemoglobin Concentration (MCHC)
 Mean Corpuscular Volume (MCV)
 Mean Platelet Volume (MPV)
 MID
 Monocytes (MONO)
 % Monocytes

Neutrophils (NEUT)
 % Neutrophils
 Nucleated Red Blood Cells (NRBC)
 Nucleated Red Blood Cells/100 WBC (NRBC/WBC%)
 Platelecrit (PCT)*
 Platelet Volume Distribution Width (PDW)*
 Platelets (PLT)
 Platelets Impedance Method (PLTi)
 Red Blood Cell Distribution Width (RDW)
 Red Blood Cells (RBC)
 Red Blood Cells, Optical (RBCo)
 White Blood Cells (WBC)
 - White Optical Count (WOC)
 - White Impedance Count (WIC)
 - Nucleated Optical Count (NOC)

*Values reported by the analyzer are not for diagnostic use



Liquichek Reticulocyte Control (A)

For Abbott Systems

- Assayed whole blood reticulocyte control for use with Abbott CELL-DYN 3200 and Ruby analyzers
- Red cells suspended in plasma-like fluid
- 75 day shelf life at 2-8°C
- 16 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description
854	Level 1 4 x 3 mL
855	Level 2 4 x 3 mL
855X	Bilevel MiniPak (1 of each level) 2 x 3 mL



Liquichek Reticulocyte Control (A-I)

For Abbott Systems

- Assayed reticulocyte control for use with the CELL-DYN Sapphire hematology analyzer
- Red cells suspended in plasma-like fluid
- 78 day shelf life at 2-8°C
- 14 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description
916	Level 1 4 x 3 mL
917	Level 2 4 x 3 mL
915X	Bilevel MiniPak (1 of each level) 2 x 3 mL

Parameters for Liquichek™ Reticulocyte Control (A) and Liquichek™ Reticulocyte Control (A-I)

Immature Reticulocyte Fraction (IRF)*
 Red Blood Cells (RBC)*
 Reticulocyte (Retic)

*Available in Liquichek™ Reticulocyte Control (A-1) only.

Liquichek Hematology Control (C)

For Beckman Coulter Systems



- Assayed hematology control for monitoring values obtained on Beckman Coulter hematology analyzers with complete CBC and VCS 5-part differential technology
- Assayed values for the UniCel DxH 600/800, DxH900 and LH 700 analyzers
- Blood cells suspended in plasma-like fluid
- 105 day shelf life at 2-8°C
- 14 day open-vial stability at 2-8°C or 14 thermal cycles, whichever comes first

Ordering Information

Cat #	Description
904	Trilevel (4 of each level) 12 x 5 mL
905	Low 4 x 5 mL
906	Normal 4 x 5 mL
907	High 4 x 5 mL
904X	Trilevel MiniPak (1 of each level) 3 x 5 mL

Parameters

Basophils (BA)
 Eosinophils (EO)
 Hematocrit (HCT)
 Hemoglobin (HGB)
 Lymphocytes (LYMPH)
 Mean Corpuscular Hemoglobin (MCH)
 Mean Corpuscular Hemoglobin Concentration (MCHC)
 Mean Corpuscular Volume (MCV)
 Mean Platelet Volume (MPV)

Monocytes (MONO)
 Neutrophils (NEUT)
 Platelecrit (PCT)*
 Platelet Distribution Width (PDW)*
 Platelets (PLT)
 RDW-SD†
 Red Blood Cell Distribution Width (RDW)
 Red Blood Cells (RBC)
 White Blood Cells (WBC)

*Values reported by analyzer are not for diagnostic use.
 †For Beckman Coulter LH and Unicel DxH 600/800 only.

Liquichek Reticulocyte Control (C)

For Beckman Coulter Systems



- Assayed hematology control for monitoring values obtained on Beckman Coulter analyzers using VCS technology
- Red cells suspended in plasma-like fluid
- 14 day open-vial stability at 2-8°C
- 75 day shelf life at 2-8°C

Ordering Information

Cat #	Description
826	Level 1 4 x 4 mL
827	Level 2 4 x 4 mL
828	Level 3 4 x 4 mL
827X	Trilevel MiniPak (1 of each level) 3 x 4 mL

Parameters

Immature Reticulocyte Fraction (IRF)
 Mean Reticulocyte Volume (MRV)
 Red Blood Cells (RBC)
 Reticulocyte (Retic)

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Liquichek Hematology Control (S)

For Siemens Systems



- An assayed hematology control for monitoring values obtained on Siemens analyzers with 3-part and 5-part differential technology
- Blood cells suspended in plasma-like fluid
- 80 day shelf life at 2-8°C
- 14 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description
449	Trilevel (4 of each level) 12 x 3.5 mL
449X	Trilevel MiniPak (1 of each level) 3 x 3.5 mL

Parameters

Basophils (BASO)	Mean Platelet Volume (MPV)
Cellular Hemoglobin Concentration Mean (CHCM)	Monocytes (MONO)
Eosinophils (EOS)	NEUT X
Hematocrit (HCT)	NEUT Y
Hemoglobin (HGB)	Neutrophils (NEUT)
Hemoglobin Concentration Distribution Width (HDW)	Platelets (PLT)
Large Unstained Cells (LUC)	Red Blood Cell Distribution Width (RDW)
Lymphocytes (LYMPH)	Red Blood Cells (RBC)
Mean Corpuscular Hemoglobin (MCH)	White Blood Cells (WBC)
Mean Corpuscular Hemoglobin Concentration (MCHC)	White Blood Cells Counts Measured in Perox Channel (WBCP Perox)
Mean Corpuscular Volume (MCV)	

Liquichek Reticulocyte Control (S)

For Siemens Systems



- Assayed hematology control monitoring values obtained on Siemens analyzers with automated reticulocyte counting methods
- Red cells suspended in plasma-like fluid
- 80 day shelf life at 2-8°C
- 14 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description
921	Level 14 x 4 mL
922	Level 24 x 4 mL
923	Level 34 x 4 mL
922X	Trilevel MiniPak (1 of each level)3 x 4 mL

Parameters

Cellular Hemoglobin Concentration Mean of Gated Red Blood Cells (CHCMg)	Mean Hemoglobin Content of All Gated Red Cells (CHg)
Cellular Hemoglobin Concentration Mean of Reticulocytes (CHCMr)	Mean Hemoglobin Content of Reticulocytes (CHr)
Mean Corpuscular Volume of Gated Red Blood Cells (MCVg)	Reticulocyte (Retic)
Mean Corpuscular Volume of Reticulocytes (MCVr)	Total Red Blood Cell Counts in Retic Channel (Retic RBC)

Liquichek Hematology Control (X)

For Sysmex Systems



- An assayed hematology control for use with Sysmex hematology analyzers with 5-part differential technology
- Blood cells in plasma-like fluid
- 80 day shelf life at 2-8°C
- Stable for 15 days open-vial or 15 samplings, whichever comes first, when stored at 2-8°C

Ordering Information

Cat #	Description
487	Trilevel (4 of each level) 12 x 4.5 mL
487X	Trilevel MiniPak (1 of each level) 3 x 4.5 mL

Parameters

- | | |
|--|---|
| Basophils (BASO) | Monocytes (MONO) |
| Eosinophils (EOS) | Neutrophils (NEUT) |
| Hematocrit (HCT) | Nucleated Red Blood Cells (NRBC)* |
| Hemoglobin (HGB) | Platelets (PLT) |
| Lymphocytes (LYMPH) | Platelets, Optical Count (PLT-o)* |
| Mean Corpuscular Hemoglobin (MCH) | Red Blood Cell Distribution Width-CV (RDW-CV) |
| Mean Corpuscular Hemoglobin Concentration (MCHC) | Red Blood Cell Distribution Width-SD (RDW-SD) |
| Mean Corpuscular Volume (MCV) | Red Blood Cells (RBC) |
| Mean Platelet Volume (MPV) | White Blood Cells (WBC) |

*No claims made for expected value or stability.

Liquichek Reticulocyte Control (X)

For Sysmex Systems



- Assayed reticulocyte control for use with Sysmex hematology analyzers with automated reticulocyte counting methods
- Red cells suspended in plasma-like fluid
- 80 day shelf life at 2-8°C
- Stable for 15 days open-vial or 15 samplings, whichever comes first, when stored at 2-8°C

Ordering Information

Cat #	Description
926	Level 1 4 x 3 mL
927	Level 2 4 x 3 mL
928	Level 3 4 x 3 mL
927X	Trilevel MiniPak (1 of each level) 3 x 3 mL

Parameters

- Immature Reticulocyte Fraction (IRF)
- Red Blood Cells (RBC)
- Reticulocyte (Retic)

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Liquichek Hematology-16 Control



- Specifically designed for use on most hematology analyzers that measure up to 16 parameters and 3-part white blood cell differentials
- Can be used on either manual sampling or automatic cap-piercing instruments
- Assayed for most 2-16 parameter instruments
- Blood cell suspension
- 160 day shelf life at 2-8°C
- 21 day open-vial stability at 2-8°C

POCT methods listed

Ordering Information

Cat #	Description
760	Trilevel (2 of each level)6 x 3 mL
761	Low6 x 3 mL
762	Normal.6 x 3 mL
763	High.6 x 3 mL
760X	Trilevel MiniPak (1 of each level)3 x 3 mL

Parameters

- | | |
|--|---|
| Granulocytes (GRAN) | Mean Platelet Volume (MPV) |
| Hematocrit (HCT) | Mid-Sized Cells (MID) |
| Hemoglobin (HGB) | Monocytes (MONO) |
| Lymphocytes (LYMPH) | Platelets (PLT) |
| Mean Corpuscular Hemoglobin (MCH) | Red Blood Cell Distribution Width (RDW) |
| Mean Corpuscular Hemoglobin Concentration (MCHC) | Red Blood Cells (RBC) |
| Mean Corpuscular Volume (MCV) | White Blood Cells (WBC) |

Liquichek Hematology-16T Control



- A hematology reference control used in monitoring determinations of blood cell values on Sysmex and Abbott 3-part differential hematology analyzers
- Blood cells in plasma-like fluid
- 105 day shelf life at 2-8°C
- 14 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description
144	Trilevel (4 of each level). 12 x 2.5 mL
144X	Trilevel MiniPak (1 of each level) 3 x 2.5 mL

Parameters

- | | |
|--|---|
| Granulocytes (GRAN)* | MID* |
| Hematocrit (HCT) | MXD (W-MCC & W-MCR) |
| Hemoglobin (HGB) | NEUT (W-LCC & W-LCR) |
| LYMPH (W-SCC & W-SCR) | Platelets (PLT) |
| Lymphocytes (LYMPH)* | Red Blood Cell Distribution Width (RDW)* |
| Mean Corpuscular Hemoglobin (MCH) | Red Blood Cell Distribution Width-CV (RDW-CV) |
| Mean Corpuscular Hemoglobin Concentration (MCHC) | Red Blood Cell Distribution Width-SD (RDW-SD) |
| Mean Corpuscular Volume (MCV) | Red Blood Cells (RBC) |
| Mean Platelet Volume (MPV) | White Blood Cells (WBC) |

*No claims made for expected value or stability.

Liquichek Sedimentation Rate Control



- Bilevel, human whole blood control designed to monitor the precision of both manual and automated methods for measuring erythrocyte sedimentation rates
- Assayed for over 20 automated and manual methods, including Westergren and Wintrobe
- Stabilized human red blood cells in suspensions
- 540 day shelf life at 18-30°C
- 31 day open-vial stability at 18-30°C

Ordering Information

Cat #	Description
514	Level 1 4 x 9 mL
515	Level 2 4 x 9 mL
514X	Bilevel MiniPak (1 of each level) 2 x 9 mL

Parameter

Erythrocyte Sedimentation Rate (ESR)

Liquichek Sedimentation Rate Control may be used with the following instruments and test methods

- Diesse Mini-Cube
- Diesse Mini-Ves (Europe)
- Diesse Ves-matic 5
- Diesse Ves-Matic 10
- Diesse Ves-Matic 20/20 Plus New
- Diesse Ves-Matic 30/30 Plus
- Diesse Ves-Matic 60/60 Plus
- Diesse Ves-Matic Cube 30
- Diesse Ves-Matic Cube 30 Touch
- Diesse Ves-Matic Cube 80/200
- Diesse Ves-Matic Easy
- Diesse Ves-Matic Easy (Europe)
- Greiner Bio-One SRS100/II SRT10/II
- Kimble-Chase Disposable Westergren Tubes
- Vital Diagnostics Excyte M, 10, 20, 40
- Vital Diagnostics Excyte Mini
- Vital Diagnostics MICROsed-System
- Vital Diagnostics Monitor Series
- Polymedco Sedimat 60
- Polymedco Sediplast
- Sarstedt S-Sedivette
- Streck ESR-Auto Plus
- Westergren (Modified)

If your instrument model is not listed, please contact your nearest Bio-Rad office for current information.

Liquichek Immunoassay Plus Control



Available in Multiple Configurations

- Efficient solution for QC of routine Immunoassay and Therapeutic Drug Monitoring analytes
- Liquid, human serum based

Standard Vials

- Assigned values are available for major automated analyzers
- 2 year shelf life at -20°C to -70°C
- 14 day open-vial stability at 2-8°C for most analytes



Compact Vials

- 2 year shelf life when stored at -20°C to -50°C
- 4 day open-vial stability for all analytes when stored at 2-8°C

POCT methods listed

Ordering Information

Cat #	Description
Standard Vials	
360	Trilevel (4 of each level) 12 x 5 mL
361	Level 1 12 x 5 mL
362	Level 2 12 x 5 mL
363	Level 3 12 x 5 mL
360X	Trilevel MiniPak (1 of each level) 3 x 5 mL
Compact Vials	
267	Level 1 12 x 2.5 mL
268	Level 2 12 x 2.5 mL
269	Level 3 12 x 2.5 mL
268X	Trilevel MiniPak (1 of each level) 3 x 2.5 mL

For IntelliQ Immunoassay Plus Control, see page 13

Analytes

	Standard Vial	Compact Vial		Standard Vial	Compact Vial		Standard Vial	Compact Vial
11-Deoxycortisol*	●		Ethosuximide	●		Procainamide**	●	●
17- α -Hydroxyprogesterone**	●		Ferritin	●	●	Progesterone	●	
25-Hydroxy Vitamin D*	●		Flecainide**	●		Prolactin	●	●
Acetaminophen	●	●	Folate	●	●	Propranolol**	●	
Aldosterone	●		Fructosamine*	●		PSA	●	
Alphafetoprotein (AFP)	●	●	FSH	●	●	PSA (Free)	●	●
Amikacin	●	●	Gentamicin	●	●	PTH-MM*	●	
Amiodarone*	●		hCG	●		Quinidine	●	
Amitriptyline*	●		hCG- β Subunit	●	●	Salicylate	●	●
Androstenedione	●		hGH	●		SHBG (Sex Hormone Binding Globulin)	●	
Angiotensin I*	●		Ibuprofen**	●		Somatomedin-C*	●	
Anti-Thyroglobulin (Anti-Tg)*	●		Imipramine*	●		T3 (Free)	●	●
Anti-Thyroperoxidase (Anti-TPO)*	●		Immunglobulin A (IgA)	●	●	T3 (Total)	●	●
Caffeine	●	●	Immunglobulin E (IgE)	●	●	T3 Uptake/T-Uptake	●	●
Carbamazepine	●	●	Immunglobulin G (IgG)	●	●	T4 (Free)	●	●
Carbamazepine (Free)**	●		Immunglobulin M (IgM)	●	●	T4 (Total)	●	●
CEA	●	●	Insulin	●		TBG	●	
Chloramphenicol**	●		Iron	●	●	Testosterone	●	
CK-MB Isoenzyme	●	●	Iron (TIBC)*	●		Testosterone (Free)**	●	
Cortisol	●		LH	●	●	Theophylline	●	●
Cyclosporine*	●		Lidocaine	●	●	Thyroglobulin (Tg)*	●	
Dehydroepiandrosterone (DHEA)*	●		Lithium	●	●	Tobramycin	●	●
Desipramine**	●		N-Acetyl procainamide (NAPA)	●	●	Tricyclic Antidepressants (TCA)	●	●
DHEA Sulfate	●		Netilmicin*	●		TSH	●	●
Digoxin	●	●	Nortriptyline**	●		Valproic Acid (Free)**	●	
Disopyramide**	●		PAP	●		Valproic Acid**	●	●
Estradiol	●	●	Phenobarbital	●	●	Vancomycin	●	●
Estriol (Free)	●		Phenytoin	●	●	Vitamin B ₁₂	●	●
Estriol (Total)*	●		Phenytoin (Free)**	●				
Estrogen (Total)*	●		Primidone	●				

*No claims are made regarding performance or stability.
 **Values are not provided.

Lyphocek Immunoassay Plus Control



- Offers a vast array of popular routine immunoassay analytes
- Assigned values are available for all major automated analyzers
- Provides a highly efficient solution for laboratories that focus on routine tests
- Human serum based
- 3 year shelf life at 2-8°C
- 7 day reconstituted stability at 2-8°C for most analytes

POCT methods listed

Ordering Information

Cat #	Description
370	Trilevel (4 of each level) 12 x 5 mL
371	Level 1 12 x 5 mL
372	Level 2 12 x 5 mL
373	Level 3 12 x 5 mL
370X	Trilevel MiniPak (1 of each level) 3 x 5 mL

Analytes

25-Hydroxy Vitamin D	Desipramine**	Immunoglobulin M (IgM)	Salicylate
11-Deoxycortisol	DHEA Sulfate	Immunoreactive Trypsinogen (IRT)*	SHBG (Sex Hormone Binding Globulin)*
17- α -Hydroxyprogesterone	Digoxin	Insulin	Somatomedin-C
Acetaminophen	Disopyramide**	Iron	T3 (Free)
ACTH	Estradiol	Iron (TIBC)	T3 (Total)
Aldosterone	Estriol (Free)	LH	T3 Uptake/T-Uptake
Alphafetoprotein (AFP)	Estriol (Total)*	Lidocaine	T4 (Free)
Amikacin	Estrogen (Total)*	Lithium	T4 (Total)
Amiodarone*	Ethosuximide	N-Acetylprocainamide (NAPA)	TBG
Amitriptyline**	Ferritin	Netilmicin*	Testosterone
Androstenedione	Flecainide**	Nortriptyline**	Testosterone (Free)
Angiotensin I*	Folate	PAP	Theophylline
Anti-Thyroglobulin (Anti-Tg)*	Fructosamine**	Phenobarbital	Thyroglobulin (Tg)
Anti-Thyropoxidase (Anti-TPO)*	FSH	Phenytoin	Tobramycin
C-Peptide	Gastrin	Phenytoin (Free)**	Tricyclic Antidepressants (TCA)**
Caffeine	Gentamicin	Primidone	TSH
Calcitonin	Glucagon*	Procainamide	Valproic Acid
Carbamazepine	hCG	Progesterone	Valproic Acid (Free)**
Carbamazepine (Free)**	hCG- β Subunit	Prolactin	Vancomycin
CEA	hGH	Propranolol**	Vitamin B ₁₂
Chloramphenicol**	Imipramine**	PSA	
Cortisol	Immunoglobulin A (IgA)	PSA (Free)	
Cyclosporine*	Immunoglobulin E (IgE)	PTH (Intact)*	
Dehydroepiandrosterone (DHEA)*	Immunoglobulin G (IgG)	Quinidine	

*No claims are made regarding performance or stability.
 **Values are not provided.

Liquichek Tumor Marker Control



Available in Multiple Configurations

- Monitors the precision of tumor marker testing procedures
- Assayed, multi-analyte control
- Liquid, human based

Standard Vials

- Includes low PSA levels for high-sensitivity methods and low levels of cancer antigens and human-based Ferritin
- Human based
- 2 year shelf life when stored frozen at -20°C to -70°C
- 30 day open-vial stability at 2-8°C for most analytes



Compact Vials

- Monitors the precision of tumor marker testing procedures
- 2 year shelf life when stored at -20°C to -50°C
- 10 day open-vial stability for all analytes when stored at 2-8°C

Lyphocheck Tumor Marker Plus Control



- A comprehensive, assayed, multi-analyte control
- Ideal for monitoring both routine and esoteric tumor marker testing procedures
- Human serum based
- 3 year shelf life at 2-8°C
- 14 day reconstituted stability at 2-8°C for most analytes

For InteliQ Tumor Marker Control, see page 14

Analytes

Liquichek Tumor Marker and Lyphocheck Tumor Marker Plus Controls

	Liquid (Standard)	Liquid (Compact)	Lyophilized
β-2-Microglobulin	●	●	●
ACTH			●
Aldosterone			●
Alphafetoprotein (AFP)	●	●	●
CA 15-3	●	●	●
CA 19-9	●	●	●
CA 27.29	●		●
CA 50*1			●
CA 72-4*	●		●
CA 125	●	●	●
Calcitonin			●
Carcinoembryonic Antigen (CEA)	●	●	●
CASA*1			●
Cyfra 21-1*	●		●
Ferritin	●	●	●
hCG	●	●	●
hCG-β Subunit			●
HE4	●		
HER2/neu	●		
Insulin Growth Factor I (IGF-I)	●		
Neuron Specific Enolase (NSE)*			●
PAP	●		●
ProGRP*	●		
Prolactin	●	●	●
PSA (Free)	●	●	●
PSA (Total)	●	●	●
S-100*			●
SCC*	●		
Thyroglobulin (Tg)	●		●

*International use only. Not available for diagnostic use in the United States.
 1No claim is made for expected values or stability.

Ordering Information

Cat #	Description
Liquichek Tumor Marker Control, Standard Vials	
547	Level 16 x 2 mL
548	Level 26 x 2 mL
549	Level 36 x 2 mL
548X	Trilevel MiniPak (1 of each level)3 x 2 mL
Liquichek Tumor Marker Control, Compact Vials	
27114	Level 16 x 2 mL
27115	Level 26 x 2 mL
27116	Level 36 x 2 mL
27115X	Trilevel MiniPak (1 of each level)3 x 2 mL

Cat #	Description
Lyphocheck Tumor Marker Plus Control	
367	Level 16 x 2 mL
368	Level 26 x 2 mL
369	Level 36 x 2 mL
368X	Trilevel MiniPak (1 of each level)3 x 2 mL

Liquichek Specialty Immunoassay Control



- Offers complex, high usage specialty immunoassay analytes for laboratories conducting specialty testing in support of operating/emergency rooms, hospitals and physician offices
- Liquid, human serum based
- 3 year shelf life at -20°C to -70°C
- 30 day open-vial stability at 2-8°C for most analytes

Ordering Information

Cat #	Description	Volume
359	Level LTA	.6 x 5 mL
364	Level 1	.6 x 5 mL
365	Level 2	.6 x 5 mL
366	Level 3	.6 x 5 mL
359X	Four Level MiniPak (1 of each level)	.4 x 5 mL

For InteliQ Specialty Immunoassay Control, see page 14

Analytes

25-Hydroxy Vitamin D	Anti-Thyroperoxidase (Anti-TPO)	Erythropoietin (EPO)	PTH (Intact)
Anti-Thyroglobulin (Anti-Tg)	C-Peptide	Insulin Growth Factor I (IGF-I)	Osteocalcin

NEW | Liquichek Specialty Immunoassay Plus Control



- Enhanced formulation with high-usage analytes, including Procalcitonin, IL-6, and Active Vitamin B₁₂ for monitoring the precision of specialty immunoassay testing procedures
- Liquid, human serum based
- Multilevel control measure clinically relevant targets
- 3 year shelf life at -20°C to -70°C
- 30 day open-vial stability at 2-8°C for most analytes

Ordering Information

Cat #	Description	Volume
12015644	Level Low	.6 x 5 mL
12015645	Level 1	.6 x 5 mL
12015646	Level 2	.6 x 5 mL
12015647	Level 3	.6 x 5 mL
12019576	Level Low MiniPak	.1 x 5 mL
12015648	Trilevel MiniPak	.3 x 5 mL

For InteliQ Specialty Immunoassay Plus Control, see page 14

Analytes

25-Hydroxy Vitamin D	C-Peptide	Interleukin-6 (IL-6)	Thyroid Stimulating Immunoglobulin (TSI)
Active Vitamin B ₁₂	Erythropoietin (EPO)	PTH (Intact)	Anti-TSH Receptor (TRAb)
Anti-Thyroglobulin (Anti-Tg)	Fructosamine	Osteocalcin	
Anti-Thyroperoxidase (Anti-TPO)	Insulin Growth Factor I (IGF-I)	Procalcitonin (PCT)	

Lyphocheck Specialty Immunoassay Control



- Includes popular analytes with increasing clinical utilization
- Human serum based
- 3 year shelf life at 2-8°C
- 30 day reconstituted stability at 2-8°C for most analytes

Ordering Information

Cat #	Description	Volume
27124	Level 1	.6 x 2 mL
27125	Level 2	.6 x 2 mL
27126	Level 3	.6 x 2 mL
27125X	Trilevel MiniPak (1 of each level)	.3 x 2 mL

Analytes

25-Hydroxy Vitamin D	PTH (Intact)	Sex Hormone Binding Globulin (SHBG)
Erythropoietin (EPO)	Procalcitonin (PCT)	

Lyphocheck Anemia Control



- Monitors test precision at clinically significant low/deficient values for anemia-related analytes
- Human serum based
- 3 year shelf life at 2-8°C
- 14 day reconstituted stability at 2-8°C for most analytes

Ordering Information

Cat #	Description	Volume
500	Single Level	.6 x 3 mL
500X	MiniPak	.1 x 3 mL

Analytes

Ferritin	Iron	Iron (UIBC)	TSH	Prealbumin
Folate	Iron (TIBC)	Transferrin	Vitamin B ₁₂	

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Liquicheck AMH Control



- Liquid convenience to reduce handling errors
- Three levels targeted at low, medium and high clinically-relevant ranges
- Human base to more closely resemble a patient sample
- 26 month shelf life at -20°C to -70°C
- 14 day open-vial stability at 2-8°C

Ordering Information

Catalog #	Description
12006327	Trilevel (2 of each level)6 x 2 mL

Analyte

Anti-Mullerian Hormone

Lymphocheck Fertility Control



- Designed to monitor the precision of tests performed for fertility or pregnancy
- Elevated hCG and hCG-β Subunit levels allow laboratories to monitor their dilution procedure
- Human serum based
- 3 year shelf life at 2-8°C
- 7 day reconstituted stability at 2-8°C

POCT methods listed

Ordering Information

Cat #	Description
480	Trilevel (4 of each level).12 x 5 mL
481	Level 112 x 5 mL
482	Level 212 x 5 mL
483	Level 312 x 5 mL
480X	Trilevel MiniPak (1 of each level)3 x 5 mL

Analytes

Alphafetoprotein (AFP)	Estriol (Free)	FSH	LH
Digitoxin	Estriol (Total)*	hCG	Progesterone
Estradiol	Ferritin	hCG-β Subunit	Testosterone

*No claims are made regarding performance or stability.

Liquicheck Maternal Serum 1st Trimester Control



- Liquid, assayed quality control
- Designed to monitor the precision of immunoassay test methods used during maternal serum first trimester screening
- Includes clinically relevant levels of Free beta hCG and PAPP-A
- Human serum based
- 2 year shelf life at -20°C to -70°C
- 15 day open-vial stability at 2–8°C

Available outside the United States only

Ordering Information

Catalog #	Description
636	Trilevel (2 of each level)6 x 1 mL
636X	Trilevel MiniPak (1 of each level)3 x 1 mL

Analytes

hCG	PAPP-A
hCG-β Subunit (Free)	

Liquicheck Maternal Serum II Control



- Liquid, assayed quality control serum
- Analytes at clinically significant levels to monitor the precision of maternal serum analysis
- Human serum
- Includes the hormone Inhibin A
- Very elevated hCG
- Very low levels of Free Estriol and AFP
- 30 day open-vial stability at 2–8°C

Ordering Information

Catalog #	Description
402	Level 16 x 2.5 mL
403	Level 26 x 2.5 mL
404	Level 36 x 2.5 mL
403X	Trilevel MiniPak (1 of each level)3 x 2.5 mL

Analytes

AFP	Estriol (Free)
hCG	Inhibin A

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Liquichek Immunoassay Premium Control



- Combination of 62 analytes including routine Immunoassay, Specialty Immunoassay, Therapeutic Drug Monitoring (TDM) and Tumor Marker analytes
- Introduces a new level of efficiency by combining multiple high-usage quality controls
- Reduces QC set-up time, eliminates waste and provides cost and labor savings
- Liquid, human serum based
- 2 year shelf life at -20°C to -70°C
- 14 day open-vial stability at 2-8°C for most analytes
- 28 day unopened stability at 2-8°C for most analytes

Ordering Information

Cat #	Description
27110	Trilevel (2 of each level)6 x 5 mL
27111	Level 16 x 5 mL
27112	Level 26 x 5 mL
27113	Level 36 x 5 mL
27110X	MiniPak (1 of each level).3 x 5 mL

Analytes

Immunoassay	T4 (Free)	Carbamazepine	Valproic Acid
Acid Phosphatase (Prostatic)	T4 (Total)	Digoxin	Valproic Acid (Free)
AFP	TBG	Disopyramide	Vancomycin
Cortisol	Testosterone	Ethosuximide	Tumor Marker
DHEA-Sulfate	TSH	Gentamicin	β-2-Microglobulin
Estradiol	Vitamin B ₁₂	Lidocaine	CA 125
Estriol (Free)	Immunology	Lithium	CA 15-3
Ferritin	IgA	NAPA (N-Acetyl Procainamide)	CA 19-9
Folate	IgE	Phenobarbital	CEA (Carcinoembryonic Antigen)
FSH	IgG	Phenytoin	Cyfra 21-1*
hCG (Total Beta)	IgM	Phenytoin (Free)	PSA (Free)
hGH	Specialty Immunoassay	Primidone	PSA (Total)
Insulin	25-Hydroxy Vitamin D (Total)	Procainamide	
LH	SHBG	Quinidine	
Progesterone	TDM	Salicylate	
Prolactin	Acetaminophen	Theophylline	
T3 Uptake/T Uptake	Amikacin	Tobramycin	
T3 (Free)	Caffeine	Tricyclic Antidepressant (TCA)	
T3 (Total)		Screen	

*Not available for diagnostic use in the United States.

Lyphechek Hypertension Markers Control



- Provides relevant values in monitoring precision for analytes specific to hypertension disorders
- Human serum based
- 2 year shelf life at 2-8°C
- 21 day reconstituted stability at 2-8°C for most analytes

Ordering Information

Catalog #	Description
600	Trilevel (2 of each level)6 x 2 mL
600X	Trilevel MiniPak (1 of each level)3 x 2 mL

Analytes

ACTH	Cortisol
ADH*	Renin
Aldosterone	

*No claims are made regarding performance or stability.

Liquichek Immunology Control



Available in Multiple Configurations

- Comprehensive control to monitor the precision of serum proteins and related analyte test procedures
- Liquid, human serum based

Standard Vials

- Provides assayed values for most major integrated immunoassay platforms
- Contains more than 30 serum proteins and analytes, including Anti-Cyclic Citrullinated Peptide (Anti-CCP) and Rheumatoid Factor
- 2 year shelf life at -20°C to -70°C
- 30 day open-vial stability at 2-8°C for most analytes

Compact Vials

- 5 day open-vial stability for all analytes when stored at 2-8°C
- 2 year shelf life when stored at -20°C to -50°C

POCT methods listed for standard vials

Ordering Information

Cat #	Description
Standard Vials	
591	Level 1 6 x 1 mL
592	Level 2 6 x 1 mL
593	Level 3 6 x 1 mL
590X	Trilevel MiniPak (1 of each level) 3 x 1 mL
594	Level 1 6 x 3 mL
595	Level 2 6 x 3 mL
596	Level 3 6 x 3 mL
Compact Vials	
281	Level 1 6 x 2.5 mL
282	Level 2 6 x 2.5 mL
283	Level 3 6 x 2.5 mL
282X	Trilevel MiniPak (1 of each level) 3 x 2.5 mL

For InteliQ Immunology Control, see page 13

Analytes

α-1-Acid Glycoprotein	Apolipoprotein A-1	Immunoglobulin A (IgA)	Protein Serum (Total)
α-1-Antitrypsin	Apolipoprotein B	Immunoglobulin E (IgE)	Retinol Binding Protein (RBP)
α-2-Macroglobulin	C1 Inhibitor	Immunoglobulin G (IgG)	Rheumatoid Factor
β-2-Microglobulin	Ceruloplasmin	Immunoglobulin G Subclasses 1-4	Transferrin
ADNase B*	Complement C3	Immunoglobulin M (IgM)	
Albumin	Complement C4	Kappa Light Chain	
Anti-Cyclic Citrullinated Peptide (Anti-CCP)*	CRP	Lambda Light Chain	
Antistreptolysin O (ASO)	Cystatin C [†]	Lipoprotein (a)*	
Antithrombin III (AT III)*	Ferritin	Prealbumin	
	Haptoglobin	Properdin Factor B**/**	

*Available only in standard vials.
**Values are not provided.

Lyphochek Immunology Plus Control



- Contains 23 serum proteins and related analytes
- Lyophilized, human serum based
- 3 year shelf life at 2-8°C
- 15 day reconstituted stability at 2-8°C

POCT Methods Listed

Ordering Information

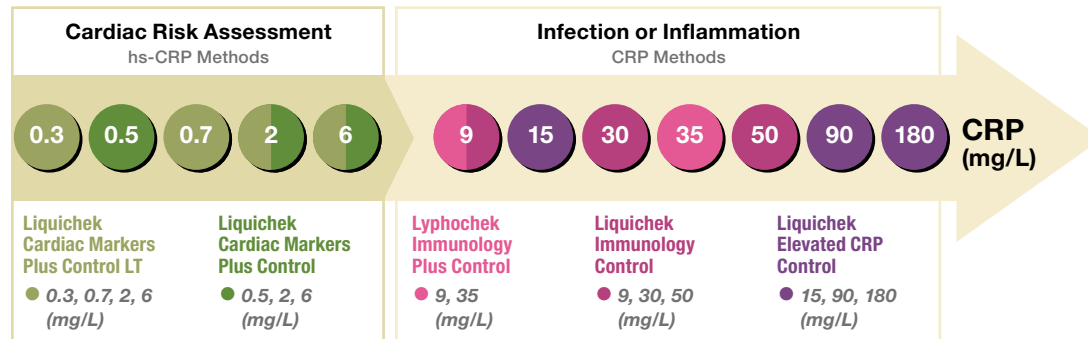
Cat #	Description
430	Bilevel (6 of each level) 12 x 1 mL
430X	Bilevel MiniPak (1 of each level) 2 x 1 mL

Analytes

α-1-Acid Glycoprotein	Ceruloplasmin	hs-CRP	Lambda Light Chain
α-1-Antitrypsin	Complement C1q*	Immunoglobulin A (IgA)	Prealbumin
α-2-Macroglobulin	Complement C3	Immunoglobulin E (IgE)	Retinol Binding Protein (RBP)
β-2-Microglobulin	Complement C4	Immunoglobulin G (IgG)	Rheumatoid Factor
Albumin	CRP	Immunoglobulin M (IgM)	Transferrin
Antistreptolysin O (ASO)	Haptoglobin	Kappa Light Chain	

*No claims are made regarding performance or stability.

Target CRP Levels in Bio-Rad Controls



Values reflect typical assayed values for these products. Recovered values may vary depending on methodology.

Liquichek Elevated CRP Control



- For monitoring the precision of C-Reactive Protein (CRP) assays as a general marker for infection and inflammation
- Target CRP levels of 15, 90 and 180 mg/L
- Liquid, human serum based
- 3 year shelf life at -20°C to -70°C
- 30 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description
254	Level 1 12 x 1 mL
255	Level 2 12 x 1 mL
256	Level 3 12 x 1 mL
255X	Trilevel MiniPak (1 of each level) 3 x 1 mL

Analyte

CRP

Liquichek Rheumatoid Factor Control



- For monitoring the precision of Rheumatoid Factor testing procedures including nephelometric and latex agglutination methods
- Target Rheumatoid Factor levels of 34, 79 and 150 U/mL
- Liquid, human serum based
- 2 year shelf life at -20°C to -70°C
- 30 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description
501	Level 1 6 x 2 mL
502	Level 2 6 x 2 mL
503	Level 3 6 x 2 mL
502X	Trilevel MiniPak (1 of each level) 3 x 2 mL

Analyte

Rheumatoid Factor

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Liquichek Spinal Fluid Control



Available in Multiple Configurations

- Multi-analyte control with relevant concentrations of analytes for cerebrospinal fluid tests
- Liquid, human based

Standard Vials

- 2 year shelf life at 2-8°C
- 30 day open-vial stability at 2-8°C

Compact Vials

- 2 year shelf life at 2-8°C
- 30 day open-vial stability for all analytes when stored at 2-8°C



Ordering Information

Cat #	Description	Volume
Standard Vials		
751	Level 1	.6 x 3 mL
752	Level 2	.6 x 3 mL
750X	Bilevel MiniPak (1 of each level)	.2 x 3 mL
Compact Vials		
303	Level 1	6 x 2.5 mL
304	Level 2	6 x 2.5 mL
302X	Bilevel MiniPak (1 of each level)	2 x 2.5 mL

Analytes

Albumin
Chloride
Glucose
Immunoglobulin A (IgA)*
Immunoglobulin G (IgG)
Immunoglobulin M (IgM)*
Lactate (Lactic Acid)

Lactate Dehydrogenase (LDH)
Protein (Total)
Protein Electrophoresis**
Sodium

*Available only in standard vials.

**Preatbumin, Albumin, α -1-Globulin, α -2-Globulin, β -Globulin, γ -Globulin.

Lyphocheck Allergen sIgE Control



- Multi-analyte control designed to monitor the precision of *in vitro* allergy test systems
- Provides assayed values for 15 commonly tested allergen-specific IgE antibodies
- Insert listings for fully-automated allergy diagnostic test systems
- Human sourced material
- 3 year shelf life at 2-8°C
- 28 day reconstituted stability at 2-8°C

Ordering Information

Cat #	Description	Volume
12000141	Negative	.6 x 2 mL
12000142	Panel A	.6 x 2 mL
12000143	Negative MiniPak	.1 x 2 mL
12000144	Panel A MiniPak	.1 x 2 mL

Analytes

D1 House Dust Mite (*Dermatophagoides pteronyssinus*)
D2 House Dust Mite (*Dermatophagoides farinae*)
E1 Cat Dander (*Felis domesticus*)
E3 Horse Dander (*Equus caballus*)
E5 Dog Dander (*Canis familiaris*)
F1 Egg White (*Gallus spp.*)
F2 Cow Milk (*Bos. spp.*)
F13 Peanut (*Arachis hypogaea*)

G2 Bermuda Grass (*Cynodon dactylon*)
G3 Cocksfoot/Orchard Grass (*Dactylis glomerata*)
G6 Timothy Grass (*Phleum pratense*)
M3 Mold (*Aspergillus fumigatus*)
M6 Mold (*Alternaria tenuis*)
T3 Birch (*Betula*)
W6 Mugwort (*Artemisia vulgaris*)

Liquichek Autoimmune Controls



- Individual unassayed pattern and detection controls for monitoring indirect immunofluorescent testing of antinuclear antibodies (ANA) and autoantibodies associated with systemic autoimmune diseases
- Liquid, human serum based
- 2 year shelf life at 2–8°C
- 6 month open-vial stability at 2–8°C

Package Description

- Negative and weak positive controls packaged separately
- 0.5 mL per vial

Ordering Information

Cat #	Description
107	ANA Control, Homogeneous Pattern. 3 x 0.5 mL
119	Anti-nDNA Control. 3 x 0.5 mL
127	Anti-Mitochondrial Control. 3 x 0.5 mL
129	Anti-Smooth Muscle Control 3 x 0.5 mL
130	Autoimmune Negative Control. 3 x 0.5 mL
131	ANA Controls Set, Positive 4 x 0.5 mL (Homogeneous, Speckled, Centromere and Nucleolar Patterns)

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Hepatitis and Retrovirus

VIROTROL I



- An unassayed quality control
- Human serum based
- 3 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C
- Available in multiple configurations and fill sizes

Ordering Information

Catalog #	Description
00100A	Single Level/Dropper Bottle 1 x 5 mL
00100B	Single Level/Dropper Bottle 1 x 5 mL
00100C	Single Level/Dropper Bottle 1 x 5 mL
00100E	Single Level/Dropper Bottle 1 x 5 mL
00100F	Single Level/Dropper Bottle 1 x 5 mL
00100H	Single Level/Dropper Bottle 1 x 5 mL
00101B	Single Level/Tube 10 x 4 mL
00101C	Single Level/Tube 10 x 4 mL
00101E	Single Level/Tube 10 x 4 mL
00101F	Single Level/Tube 10 x 4 mL

Analytes

Anti-CMV	Anti-HIV-1
Anti-HBc	Anti-HTLV-I
Anti-HCV	HBsAg

Assayed VIROTROL I-C | Assayed VIROTROL I-E | Assayed VIROTROL I-F



- Assayed quality controls
 - Human serum based
 - Values provided for popular automated methods
 - 3 year shelf life at 2-8°C
 - 60 day open-vial stability at 2-8°C
- Available outside the United States only.

Ordering Information

Cat #	Description
Assayed VIROTROL I-C	
00164	Single Level/Dropper Bottle 1 x 5 mL
Assayed VIROTROL I-E	
00166	Single Level/Dropper Bottle 1 x 5 mL
Assayed VIROTROL I-F	
00168	Single Level/Dropper Bottle 1 x 5 mL

Analytes

Anti-CMV*	Anti-HIV-1
Anti-HBc	Anti-HTLV-I*
Anti-HCV	HBsAg

*Values not provided.

VIROTROL II



- An unassayed quality control
- Human serum based
- 3 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C
- Available in multiple configurations and fill sizes

Ordering Information

Catalog #	Description
00104A	Single Level/Dropper Bottle 1 x 5 mL
00104B	Single Level/Dropper Bottle 1 x 5 mL
00115A	Single Level/Tube 10 x 4 mL
00115B	Single Level/Tube 10 x 4 mL

Analytes

Anti-HAV
Anti-HBs

Please note that not all products may be available in all countries. Refer to myelnserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Hepatitis and Retrovirus

Assayed VIROTROL II-B



- Assayed quality control
- Human serum based
- 3 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C

Available outside the United States only.

Ordering Information

Catalog #	Description
00152	Single Level/Dropper Bottle 1 x 5 mL

Analytes

Anti-HAV
Anti-HBs

VIROTROL III



- An unassayed quality control
- Human serum based
- 3 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C

Ordering Information

Catalog #	Description
00110A	Single Level/Dropper Bottle 1 x 5 mL
00110C	Single Level/Dropper Bottle 1 x 5 mL

Analytes

Anti-HAV
HAV-IgM
HBc-IgM

VIROTROL IV



- An unassayed quality control
- Human serum based
- 3 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C

Ordering Information

Catalog #	Description
00111	Single Level/Dropper Bottle 1 x 5 mL

Analyte

Anti-HBe

VIROTROL HIV-1 Ag



- An unassayed quality control
- Human serum based
- 2 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C

Ordering Information

Catalog #	Description
00108A	Single Level/Dropper Bottle 1 x 5 mL
00108B	Single Level/Dropper Bottle 1 x 5 mL

Analytes

HIV-1 Ag

Please note that not all products may be available in all countries. Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Hepatitis and Retrovirus

VIROTROL HIV-1 gO



- An unassayed control used to detect antibodies to HIV-1 group O
- Human serum based
- 2 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C

Ordering Information

Catalog #	Description
00113	Single Level/Tube5 x 4 mL
00113X	Single Level/Tube1 x 4 mL

VIROTROL HIV-2



- An unassayed quality control
- Human serum based
- 2 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C

Ordering Information

Catalog #	Description
00105A	Single Level/Dropper Bottle.1 x 5 mL
00105B	Single Level/Dropper Bottle.1 x 5 mL
00105C	Single Level/Dropper Bottle.1 x 5 mL

Analyte

Anti-HIV-2

VIROTROL PLUS-R



- An unassayed quality control
- Human serum based
- 2 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C

Ordering Information

Catalog #	Description
12000538	Single Level/Tube5 x 4 mL
12000539	Single Level/Tube1 x 4 mL

Analytes

Anti-CMV	Anti-HTLV I/II
Anti-HBc	Anti- <i>Treponema pallidum</i> (Syphilis)
Anti-HCV	HBsAg
Anti-HIV 1/2	

VIROTROL HAV-IgM



- An unassayed quality control
- Human serum based
- 3 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C

Ordering Information

Catalog #	Description
00142A	Single Level/Dropper Bottle.1 x 5 mL

VIROTROL HBeAg



- An unassayed quality control
- Human serum based
- 3 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C

Ordering Information

Catalog #	Description
00144	Single Level/Dropper Bottle.1 x 5 mL

Analyte

HBeAg

Please note that not all products may be available in all countries. Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Hepatitis and Retrovirus

VIROCLEAR



- An unassayed, non-reactive quality control
- Human serum based
- 3 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C
- Available in multiple configurations and fill sizes

Ordering Information

Cat #	Description
00106	Single Level/Dropper Bottle 1 x 5 mL
00112	Single Level/Tube 10 x 4 mL

Analytes

Non-reactive for:

Anti-CMV	Anti-HCV	HAV-IgM	Non-treponemal Antibodies (Syphilitic Reagin)
Anti-HAV	Anti-HIV-1	HBeAg	<i>Treponema pallidum</i> (Syphilis) IgG
Anti-HBc	Anti-HIV-2	HBsAg	<i>Treponema pallidum</i> (Syphilis) IgM
Anti-HBe	Anti-HTLV-I	HIV-1 Ag	<i>Treponema pallidum</i> (Syphilis) Total
Anti-HBs	Anti-HTLV-II		

Congenital / Pediatric Diseases

VIROTROL ToRCH



- An unassayed quality control
- Human serum based
- 2 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C

Ordering Information

Catalog #	Description
00109	Single Level/Dropper Bottle 1 x 5 mL

Analytes

Cytomegalovirus (CMV) IgG	Rubella Virus IgG
Herpes Simplex Virus Type 1 (HSV-1) IgG	<i>Toxoplasma gondii</i> IgG
Herpes Simplex Virus Type 2 (HSV-2) IgG	

VIROTROL ToRCH-M



- An unassayed quality control
- Human serum based
- 2 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description
00117A	Single Level/Dropper Bottle 1 x 5 mL
00117B	Single Level/Dropper Bottle 1 x 5 mL

Analytes

Cytomegalovirus (CMV) IgG	Herpes Simplex Virus Type 1 (HSV-1) IgG	Rubella Virus IgG	<i>Toxoplasma gondii</i> IgG
Cytomegalovirus (CMV) IgM	Herpes Simplex Virus Type 2 (HSV-2) IgG	Rubella Virus IgM	<i>Toxoplasma gondii</i> IgM

Please note that not all products may be available in all countries. Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Congenital / Pediatric Diseases

VIROCLEAR ToRCH



- An unassayed, non-reactive quality control
- Human serum based
- 2 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description
00118	Single Level/Dropper Bottle. 1 x 5 mL

Analytes

Non-reactive for:

Cytomegalovirus (CMV) IgG
 Cytomegalovirus (CMV) IgM
 Herpes Simplex Virus Type 1 (HSV-1) IgG

Herpes Simplex Virus Type 1 (HSV-1) IgM
 Herpes Simplex Virus Type 2 (HSV-2) IgG

Herpes Simplex Virus Type 2 (HSV-2) IgM
 Rubella Virus IgG
 Rubella Virus IgM

Toxoplasma gondii IgG
Toxoplasma gondii IgM

Liquichek ToRCH Plus Control



- A bilevel control used to monitor the precision of the most widely used serology test methods
- Liquid, human serum based
- 3 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C
- Positive and negative levels packaged separately

Ordering Information

Catalog #	Description
239	Positive 3 x 3 mL
239X	Positive MiniPak. 1 x 3 mL
227	Positive (Assayed)*. 3 x 3 mL
227X	Positive MiniPak (Assayed)*. 1 x 3 mL
228	Negative 3 x 3 mL
228X	Negative MiniPak. 1 x 3 mL

*Assayed control available outside the United States only.

Analytes

Cytomegalovirus (CMV) IgG
 Epstein-Barr Virus (EBV) IgG
Helicobacter pylori IgG
 Herpes Simplex Virus Type 1 (HSV-1) IgG
 Herpes Simplex Virus Type 2 (HSV-2) IgG

Rubella Virus IgG
 Syphilis RPR
Toxoplasma gondii IgG
Treponema pallidum (Syphilis) IgG
 Varicella Zoster Virus (VZV) IgG

VIROTROL MuMZ and VIROCLEAR MuMZ



- Unassayed quality controls
- Use VIROTROL MuMZ as a reactive control and VIROCLEAR MuMZ as a non-reactive control for the analytes listed at right
- Human serum based
- 2 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description
VIROTROL MuMZ	
00119	Single Level/Dropper Bottle. 1 x 5 mL
VIROCLEAR MuMZ	
00133	Single Level/Dropper Bottle. 1 x 5 mL

Analytes

Mumps Virus IgG
 Rubeola Virus (Measles) IgG
 Varicella Zoster Virus (VZV) IgG

Please note that not all products may be available in all countries. Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Respiratory

VIROTROL SARS-CoV-2



- An unassayed reactive control
- Human plasma based
- 2 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C

Ordering Information

200300A	Single Level/Tube	1 x 4 mL
200305A	Single Level/Tube5 x 4 mL

Analytes

SARS-CoV-2 Total IgG/IgM
SARS CoV-2 IgG

VIROCLEAR SARS-CoV-2



- An unassayed negative control
- Human plasma based
- 2 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C

Ordering Information

200500	Single Level/Tube	1 x 4 mL
200505	Single Level/Tube5 x 4 mL

Analytes

Non-reactive for:
SARS-CoV-2 Total IgG/IgM
SARS CoV-2 IgG

Please note that not all products may be available in all countries. Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Sexually Transmitted Diseases

VIROTROL Syphilis LR-A



- A low reactive (LR), unassayed quality control
- Human serum based
- 3 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description
00171	Single Level/Tube5 x 4 mL
00171X	Single Level/Tube1 x 4 mL

Analyte

Anti-*Treponema pallidum* (Syphilis)

VIROTROL RPR Panel



- An unassayed quality control
- VIROTROL RPR Panel consists of one non-reactive member, one minimally reactive member, and one moderately to highly reactive titratable member
- Human serum based
- 2 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C

Ordering Information

Catalog #	Description
01400	Three Levels/Vial 3 x 1.5 mL

Analyte

Non-treponemal Antibodies (Syphilitic Reagin)

Specialty Infectious Disease

VIROTROL Chagas



- An unassayed, reactive quality control
- Human serum based
- 2 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C

Ordering Information

Catalog #	Description
00125	Single Level/Tube5 x 4 mL
00125X	Single Level/Tube1 x 4 mL

Analyte

Trypanosoma cruzi (*T. cruzi*) Antibodies

VIROTROL EBV



- An unassayed quality control
- Human serum based
- 2 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C

Ordering Information

Catalog #	Description
00128A	Single Level/Dropper Bottle.1 x 5 mL

Analytes

EBNA IgG
Epstein-Barr Virus (VCA) IgG
Epstein-Barr Virus (VCA) IgM

Please note that not all products may be available in all countries. Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Specialty Infectious Disease

VIROCLEAR EBV



- An unassayed, non-reactive quality control
- Human serum based
- 2 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C

Ordering Information

Catalog #	Description
00131	Single Level/Dropper Bottle. 1 x 5 mL

Analytes

Non-reactive for:
 EBNA IgG
 EBV EA-D IgG
 Epstein-Barr Virus (VCA) IgG
 Epstein-Barr Virus (VCA) IgM

VIROTROL Lyme



- An unassayed quality control
- Human serum based
- 2 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C

Ordering Information

Catalog #	Description
00132	Single Level/Dropper Bottle. 1 x 5 mL

Analyte

Borrelia burgdorferi IgG

PyloriTrol



- An unassayed quality control
- Human serum based
- 2 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C

Ordering Information

Catalog #	Description
00121	Single Level/Dropper Bottle. 1 x 5 mL

Analyte

Helicobacter pylori IgG

Please note that not all products may be available in all countries. Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Liquichek Therapeutic Drug Monitoring Control (TDM)



- For labs that run a high volume of serum therapeutic drug tests or prefer dedicated TDM controls
- Includes markers for monitoring thyroid function and some esoteric drugs
- Liquid, human serum based
- 3 year shelf life at -20°C to -70°C
- 180 day unopened-vial stability at 2-8°C for most analytes
- 30 day open-vial stability at 2-8°C for most analytes

Ordering Information

Cat #	Description	Volume
724	Trilevel (4 of each level)	12 x 5 mL
725	Level 1	12 x 5 mL
726	Level 2	12 x 5 mL
727	Level 3	12 x 5 mL
724X	MiniPak (1 of each level)	.3 x 5 mL

Analytes

Acetaminophen	Digoxin	Nortriptyline*	T4 (Free)*
Amikacin	Disopyramide	Phenobarbital	T4 (Total)
Amitriptyline*	Estriol (Total)*	Phenytoin	Theophylline
Caffeine	Ethosuximide	Phenytoin (Free)	Thyroid Stimulating Hormone (TSH)
Carbamazepine	Flecainide*	Primidone	Tobramycin
Carbamazepine (Free)	Gentamicin	Procainamide	Tricyclic Antidepressants (TCA)
Chloramphenicol	Haloperidol*	Propranolol	Valproic Acid
Clonazepam*	Imipramine*	Quinidine	Valproic Acid (Free)
Cortisol	Lidocaine	Salicylate	Vancomycin
Cyclosporine*	Lithium	T3 (Free)	
Desipramine	Methotrexate	T3 (Total)	
Diazepam	N-Acetylprocainamide (NAPA)	T3 Uptake/T-Uptake	

*No claims are made regarding performance or stability.

Lyphocek Therapeutic Drug Monitoring Control (TDM)



- For labs that run a high volume of serum therapeutic drug tests or prefer dedicated TDM controls
- Contains over 40 analytes including markers for monitoring thyroid function
- Human serum based
- 3 year shelf life at 2-8°C
- 30 day reconstituted stability at 2-8°C for most analytes

Ordering Information

Cat #	Description	Volume
450	Trilevel (4 of each level)	12 x 5 mL
451	Level 1	12 x 5 mL
452	Level 2	12 x 5 mL
453	Level 3	12 x 5 mL
450X	MiniPak (1 of each level)	.3 x 5 mL

Analytes

Acetaminophen	Disopyramide	Nortriptyline	T4 (Total)
Amikacin	Estriol (Total)*	Phenobarbital	Theophylline
Amiodarone*	Ethosuximide	Phenytoin	Thyroid Stimulating Hormone (TSH)
Amitriptyline	Flecainide*	Phenytoin (Free)	Tobramycin
Caffeine	Gentamicin	Primidone	Tricyclic Antidepressants (TCA)
Carbamazepine	Haloperidol*	Procainamide	Valproic Acid
Carbamazepine (Free)	Imipramine	Propranolol	Valproic Acid (Free)
Chloramphenicol	Lidocaine	Quinidine	Vancomycin
Cortisol	Lithium	Salicylate	
Cyclosporine*	Methotrexate	T3 (Total)	
Desipramine	N-Acetylprocainamide (NAPA)	T3 Uptake/T-Uptake	
Digoxin	Netilmicin*	T4 (Free)	

*No claims are made regarding performance or stability.

Liquicheck Qualitative Urine Toxicology Control, Negative and Positive

Use with qualitative urine drug panels, including POCT methods such as the Bio-Rad TOX/See



- Human urine
- 3 year shelf life at 2-8°C
- 30 day open-vial stability at 2-25°C

POCT methods listed

Ordering Information

Cat #	Description
454	Negative Level6 x 3 mL
455	Positive Level6 x 3 mL
454X	Bilevel MiniPak (1 of each level)2 x 3 mL

Analyte Target Values

Analytes	Negative* (ng/mL)	Positive (ng/mL)
Amphetamines		
d-Amphetamine	<80	2,000
d-Methamphetamine	<80	2,000
3, 4-Methylenedioxymethamphetamine (MDMA)	<80	750
Barbiturates		
Secobarbital	<40	1,000
Benzodiazepines		
Nordiazepam	<25	3,000
Oxazepam	<25	1,000
Cannabinoids		
11-Nor-Δ-9-THC-9-COOH	<2	150
Cocaine		
Benzoylcegonine	<30	4,000
Ethanol (mg/dL)	<20	250
Lysergic Acid Diethylamide (LSD)	<0.1	2.0
Methadone	<30	750
Methaqualone	<25	750
Opiates		
Morphine (Free)	<50	3,000
Oxycodone	<25	150
Phencyclidine (PCP)	<2	200
Propoxyphene	<50	750
Tricyclic Antidepressants		
Nortriptyline	<25	2,000

*Negative level has been tested and found to be negative for the analytes listed in the package insert. Values above represent the testing method's Limit of Detection (LOD).

Liquicheck Urine Toxicology Negative Control

Use with Screening or Confirmatory methods



- Confirmed negative by GC-MS, LC-MS or other methods for 32 common drugs or drug metabolites
- Human urine
- 3 year shelf life at 2-8°C
- 60 day open-vial stability when stored refrigerated at 2-8°C

Ordering Information

Cat #	Description
460	Negative Control10 x 20 mL
460X	MiniPak1 x 20 mL

Liquichek Opiate Control



- Contains eight drugs or drug metabolites, at concentrations above and below assay cutoffs
- This control has been tested by GC/MS, LC/MS/MS and other methods to confirm the recovery of drug concentration
- Representative values for screening and confirmatory methods provided
- Human urine
- 3 year shelf life at -20°C to -70°C
- 30 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description
12005786	Level 1 6 x 5 mL
12005787	Level 2 6 x 5 mL
12002594	Level 1 MiniPak 1 x 5 mL
12002595	Level 2 MiniPak 1 x 5 mL

Gravimetric Reference Values

Analyte (ng/mL)	Level 1	Level 2	SAMHSA** Screen Cut-offs
6-Monoacetylmorphine (6-MAM, Heroin Metabolite)	7.5	12.5	10
Buprenorphine	5	15	
EDDP (Methadone Metabolite)	50	150	
Fentanyl	0.75	1.25	1
Meperidine*	100	300	
Methadone	150	450	
Oxycodone	50	150	
Tramadol	100	300	

*International use only.

**United States Substance Abuse and Mental Health Services Administration.

Liquichek Urine Toxicology Controls

SCREEN CONTROLS



Levels S1 & S2
Levels S1E & S2E

- Human urine
- 3 year shelf life at 2-8°C
- 30 day open-vial stability at 2-8°C

Ordering Information

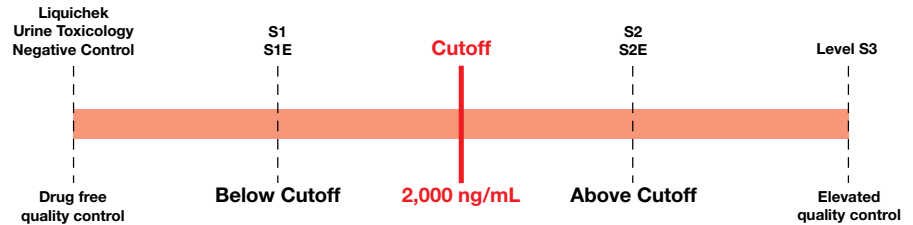
Cat #	Description
438	Level S1E 10 x 10 mL
439	Level S2E 10 x 10 mL
461	Level S1 10 x 10 mL
462	Level S2 10 x 10 mL

MiniPaks also available



Determine your cutoff

Controls on this page recommended for High Opiate cutoff (2,000 ng/mL)



Gravimetric Reference Values

Select the controls best suited for your Benzodiazepine assays

	Level S1E	Level S2E	Level S1	Level S2
<i>Example assays may include:</i>	<i>(e.g. Siemens EMIT II, EMIT DAT & Ortho VITROS)</i>		<i>(e.g. Roche OnLine)</i>	
	Lormetazepam		Nordiazepam	
Analytes (ng/mL)				
Amphetamines				
d-Amphetamine	–	–	750	1,250
d-Methamphetamine	750	1,250	–	–
MDMA (Ecstasy)	–	–	–	–
Barbiturates				
Secobarbital	150	375	150	375
Benzodiazepines				
Lormetazepam	150	375	–	–
Nordiazepam	–	–	75	375
Oxazepam	–	–	–	–
Cannabinoids				
11-Nor-Δ-9-THC-9-COOH	37.5	65	37.5	65
Cocaine				
Benzoylcegonine	225	375	225	375
Ethanol (mg/dL)	40	70	40	70
Lysergic Acid Diethylamide (LSD)	0.3	0.8	0.3	0.8
Methadone	225	375	225	375
Methaqualone	225	375	225	375
Opiates				
Morphine (Free)	1,500	2,500	1,500	2,500
Phencyclidine (PCP)	19	31	19	31
Propoxyphene	225	375	225	375
Tricyclic Antidepressants				
Nortriptyline	750	1,250	750	1,250

*United States Substance Abuse and Mental Health Services Administration. Refer to myInSerts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Liquichek Urine Toxicology Controls

SCREEN CONTROLS



Levels S1E & S2E Low Opiate

Levels S1 & S2 Low Opiate

- Human urine
- 3 year shelf life at 2-8°C
- 30 day open-vial stability at 2-8°C

Levels S10 & S20 Low Opiate

- Human urine
- 2 year shelf life at -20 to -70°C
- 30 day open-vial stability at 2-8°C
- 45 day closed-vial stability at 2-8°C

Ordering Information

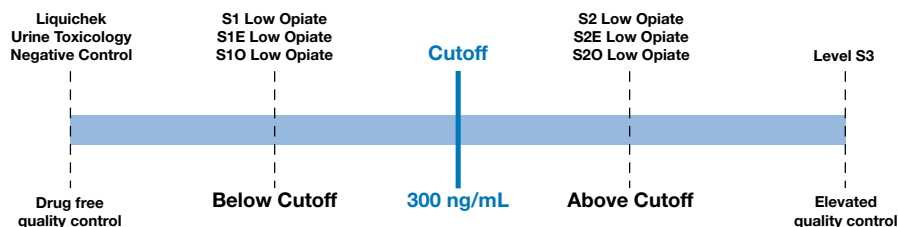
Cat #	Description
423	Level S1E Low Opiate 10 x 10 mL
424	Level S2E Low Opiate 10 x 10 mL
466	Level S1 Low Opiate 10 x 10 mL
467	Level S2 Low Opiate 10 x 10 mL
683	Level S10 Low Opiate 10 x 10 mL
706	Level S20 Low Opiate 10 x 10 mL

MiniPaks also available



Determine your cutoff

Controls on this page recommended for Low Opiate cutoff (300 ng/mL)



Gravimetric Reference Values

Select the controls best suited for your Benzodiazepine assays

	Level S1E Low Opiate	Level S2E Low Opiate	Level S1 Low Opiate	Level S2 Low Opiate	Level S10 Low Opiate	Level S20 Low Opiate
<i>Example assays may include:</i>	<i>(e.g.: Siemens EMIT II & EMIT DAT, Ortho VITROS)</i>		<i>(e.g.: Roche OnLine)</i>		<i>(e.g.: Abbott Multigent, Thermo DRI)</i>	
Analytes (ng/mL)	Lormetazepam		Nordiazepam		Oxazepam	
Amphetamines						
d-Amphetamine	–	–	750	1,250	–	–
d-Methamphetamine	750	1,250	–	–	750	1,250
MDMA (Ecstasy)	–	–	–	–	–	–
Barbiturates						
Secobarbital	150	375	150	375	150	375
Benzodiazepines						
Lormetazepam	150	375	–	–	–	–
Nordiazepam	–	–	75	375	–	–
Oxazepam	–	–	–	–	100	300
Cannabinoids						
11-Nor-Δ-9-THC-9-COOH	37.5	65	37.5	65	37.5	65
Cocaine						
Benzoyllecgonine	225	375	225	375	225	375
Ethanol (mg/dL)	40	70	40	70	40	70
Lysergic Acid Diethylamide (LSD)	0.3	0.8	0.3	0.8	0.3	0.8
Methadone	225	375	225	375	225	375
Methaqualone	225	375	225	375	225	375
Opiates						
Morphine (Free)	225	375	225	375	225	375
Phencyclidine (PCP)	19	31	19	31	19	31
Propoxyphene	225	375	225	375	225	375
Tricyclic Antidepressants						
Nortriptyline	750	1,250	750	1,250	–	–

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Liquichek Urine Toxicology Controls

SCREEN CONTROLS



Levels S1S & S2S

- Contains six drugs of abuse at concentrations 20-25% below and above SAMHSA's recommended cutoff levels
- Human urine
- 3 year shelf life at 2-8°C
- 30 day open-vial stability at 2-8°C

Ordering Information

Catalog #	Description
409	Level S1S. 10 x 10 mL
419	Level S2S. 10 x 10 mL

MiniPaks also available



Determine your cutoff

Levels S1S & S2S recommended for labs running SAMHSA* cutoffs



Level S3

- An elevated control for most urine Drugs-of-Abuse cutoffs
- For Low Opiate Cutoff
- Human urine
- 3 year shelf life at 2-8°C
- 30 day open-vial stability at 2-8°C

Ordering Information

Catalog #	Description
463	Level S3. 10 x 10 mL

MiniPaks also available



Determine your cutoff

Level S3 recommended for labs that need an Elevated/Strong Positive control

Gravimetric Reference Values

Analytes (ng/mL)	Level S1S	SAMHSA* Screen Cutoffs	Level S2S	Level S3
Amphetamines				
d-Amphetamine	–	–	–	2,000
d-Methamphetamine	375	500	625	–
MDMA (Ecstasy)	375	500	625	–
Barbiturates				
Secobarbital	–	–	–	1,000
Benzodiazepines				
Lormetazepam	–	–	–	–
Nordiazepam	–	–	–	1,000
Oxazepam	–	–	–	–
Cannabinoids				
11-Nor-Δ-9-THC-9-COOH	37.5	50	65	125
Cocaine				
Benzoyllecgonine	113	150	185	500
Ethanol (mg/dL)	–	–	–	250
Lysergic Acid Diethylamide (LSD)	–	–	–	2.0
Methadone	–	–	–	750
Methaqualone	–	–	–	750
Opiates				
Morphine (Free)	1,500	2,000	2,500	750
Phencyclidine (PCP)	19	25	31	200
Propoxyphene	–	–	–	750
Tricyclic Antidepressants				
Nortriptyline	–	–	–	2,000

*United States Substance Abuse and Mental Health Services Administration.

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Liquichek Urine Toxicology Controls

CONFIRM CONTROLS



- Human urine
- 3 year shelf life at 2-8°C
- 30 day open-vial stability at 2-8°C

Level C1

Contains drugs at the Limit of Quantitation (LOQ) or at very low concentrations

Levels C2 & C2 Low Opiate

Contains drugs at 20-25% below confirmatory drug cutoff levels

Levels C3 & C3 Low Opiate

Contains drugs at 20-25% above confirmatory drug cutoff levels

Level C4

Contains drugs at elevated concentrations

Ordering Information

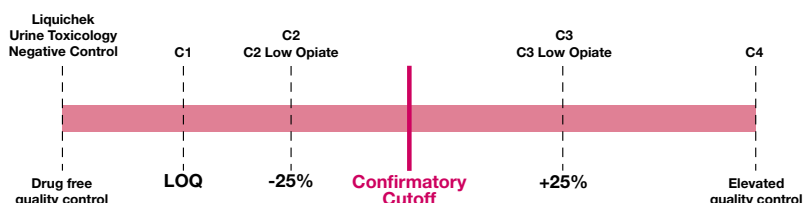
Catalog #	Description	Catalog #	Description
441	Level C1 10 x 20 mL	444	Level C4 10 x 20 mL
442	Level C2 10 x 20 mL	468	Level C2 Low Opiate 10 x 20 mL
443	Level C3 10 x 20 mL	469	Level C3 Low Opiate 10 x 20 mL

MiniPaks also available

Gravimetric Reference Values

Analytes (ng/mL)	Designed for SAMHSA* Cutoff				Level C2 Low Opiate	Level C3 Low Opiate
	Level C1	Level C2	Level C3	Level C4		
d-Amphetamine	100	188	312	750	375	625
d-Methamphetamine	100	188	312	750	375	625
3, 4-Methylenedioxymethamphetamine (MDMA)	100	188	312	750	-	-
3, 4-Methylenedioxyamphetamine (MDA)	100	188	312	750	-	-
3, 4-Methylenedioxy-N-ethylamphetamine (MDEA)	100	188	312	750	-	-
Amobarbital	80	150	250	1,200	150	250
Butalbital	80	150	250	1,200	150	250
Pentobarbital	80	150	250	1,200	150	250
Phenobarbital	80	150	250	1,200	150	250
Secobarbital	80	150	250	1,200	150	250
Alpha-Hydroxyalprazolam	80	225	375	1,000	225	375
Nordiazepam	80	225	375	1,000	225	375
Oxazepam**	**	**	**	**	**	**
11-Nor-Δ-9-THC-9-COOH	6	11.5	18.5	150	11.5	18.5
Benzoylcegonine	40	75	125	300	115	185
Ethanol (mg/dL)	20	40	70	250	40	70
Lysergic Acid Diethylamide (LSD)	0.2	0.35	0.65	4.0	0.35	0.65
Methadone	120	225	375	1,000	225	375
EDDP (Methadone Metabolite)**	**	**	**	**	**	**
Methaqualone	120	225	375	1,000	225	375
6-Monoacetylmorphine**	**	**	**	**	**	**
Codeine	120	1,500	2,500	4,000	225	375
Morphine-3-β-d-Glucuronide	120	1,500	2,500	4,000	225	375
Phencyclidine (PCP)	10	19	31	100	19	31
Norpropoxyphene	120	225	375	1,000	225	375
Propoxyphene	120	225	375	1,000	225	375

*United States Substance Abuse and Mental Health Services Administration.
 **Analyte present, no claims are made regarding performance or stability.



Refer to myelnserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Lyphocek Drug Free Serum



- Tested by HPLC and other methods
- Found to be drug free for more than 60 therapeutic drugs and illicit compounds
- Human serum
- 3 year shelf life at 2-8°C
- 10 day reconstituted stability at 2-8°C

Ordering Information

Catalog #	Description
456	Single Level 12 x 10 mL
456X	MiniPak 1 x 10 mL

Analytes

Acetaminophen	Desipramine	Meperidine	Phenytoin
Alprazolam	Desmethyldoxepin	Methadone	Primidone
Amikacin	Diazepam	Methamphetamine	Procainamide
Amitriptyline	Digitoxin	Methaqualone	Propoxyphene
Amphetamine	Digoxin	Methotrexate	Propranolol
Caffeine	Disopyramide*	Methprylon*	Pyrilamine
Cannabinoids	Doxepin	Morphine	Quinidine
Carbamazepine	Ethosuximide	N-Acetylprocainamide (NAPA)	Quinine
Chloramphenicol	Flecainide	Nicotine	Salicylate
Chlordiazepoxide	Gentamicin	Norclomipramine	Secobarbital
Cimetidine	Glutethimide	Nordiazepam	Theophylline
Clomipramine	Haloperidol	Nordoxepin	Tobramycin
Clonazepam	Hydromorphone	Nortriptyline	Trazodone
Cocaine	Ibuprofen	Oxazepam	Valproic Acid
Codeine	Imipramine	Phencyclidine (PCP)	Vancomycin
Cyclosporine	Lidocaine	Phenobarbital	
Desalkylflurazepam	Lithium	Phenothiazines	

*No claim is made for expected value or stability.

Lyphocek Benzo/TCA Control-Set B



- Used in conjunction with separation techniques, such as chromatography or mass spectrometry in identification and quantification of benzodiazepine and tricyclic antidepressant drugs
- Compatible with Roche COBAS INTEGRA Serum Benzodiazepines Assay for Alprazolam measurements
- Human serum based
- 3 year shelf life at 2-8°C
- 14 day reconstituted stability at 2-8°C

Ordering Information

Catalog #	Description
663	Level 16 x 5 mL
664	Level 26 x 5 mL
665X	Bilevel MiniPak (1 of each level)2 x 5 mL

Analytes

Alprazolam
Clomipramine
Norclomipramine
Trazodone

Clonazepam has been added to this product as a service to laboratories that wish to investigate this analyte. No claims are made regarding performance or stability.

Lyphocek Whole Blood Metals Control



- Contains major heavy metals of toxicological and environmental significance for whole blood analysis
- Assayed values for Lead include AA, ASV, ICP-MS and IDMS
- Human whole blood based
- 3 year shelf life at 2-8°C
- 14 day reconstituted stability at 2-8°C

Ordering Information

Catalog #	Description
527	Level 16 x 2 mL
528	Level 26 x 2 mL
529	Level 36 x 2 mL
528X	Trilevel MiniPak (1 of each level)3 x 2 mL

Analytes

Arsenic
Cadmium
Lead
Mercury
Thallium

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Liquichek Serum Volatiles Control



- Value assigned for volatile substances by multiple methods, including GC and clinical methods (Ethanol)
- Serum based
- 3 year shelf life at 2-8°C
- 10 day open-vial stability at 2-8°C

Ordering Information

Cat #	Description
383	Level 1 6 x 5 mL
384	Level 2 6 x 5 mL
385X	Bilevel MiniPak (1 of each level) 2 x 5 mL

Analytes

Acetone	Ethylene Glycol	Isopropanol (Isopropyl Alcohol)	Methanol (Methyl Alcohol)
Ethanol			

Lyphocheck Urine Metals Control



- Supports toxicological environmental and industrial testing by providing assayed values of critical trace elements, heavy metals and organic metabolites by AA, GC-MS, HPLC and ICP-MS
- Human urine
- 3 year shelf life at 2-8°C
- 5 day reconstituted stability at 2-8°C
- 30 day reconstituted stability at -10°C to -20°C

Ordering Information

Catalog #	Description
400	Level 1 10 x 25 mL
405	Level 2 10 x 25 mL
402X	Bilevel MiniPak (1 of each level) 2 x 25 mL

Analytes

5-Aminolevulinic Acid (ALA)	Cobalt	Mandelic Acid	Phenylglyoxylic Acid
Aluminum	Copper	Manganese	Selenium
Antimony	Fluoride	Mercury	Thallium
Arsenic	Hippuric Acid	Nickel	Trichloroacetic Acid
Cadmium	Lead	Pentachlorophenol	Zinc
Chromium	Magnesium	Phenol	

Lyphocheck Whole Blood Control



- Assayed values for Red Cell Folate for common clinical methods
- Assayed values for Lead include AA, ASV, ICP-MS
- Human whole blood based
- 3 year shelf life at 2-8°C
- 2 day reconstituted stability at 2-8°C (Folate)
- 14 day reconstituted stability at 2-8°C (Lead)
- 30 day reconstituted stability at -10° to -20°C

Ordering Information

Catalog #	Description
561	Level 1 6 x 2 mL
562	Level 2 6 x 2 mL
563	Level 3 6 x 2 mL
560X	Trilevel MiniPak (1 of each level) 3 x 2 mL

Analytes

Lead
Red Cell Folate
Serotonin*

*No claim is made for expected value or stability of this analyte.

qUAntify Advance Control



- Available in an easy-to-use plastic squeeze bottle
- Dipstick tests include Pregnancy and Microalbumin
- Human urine solution
- 18 month shelf life at 2-8°C
- 31 day open-vial stability at 2-25°C, including Ketones

Ordering Information

Cat #	Description
12007028	Bilevel (3 of each level) 6 x 12 mL
12007029	Bilevel MiniPak (1 of each level) 2 x 12 mL

Analytes

Albumin	Glucose	Microalbumin	Specific Gravity
Ascorbic Acid	hCG (Pregnancy)	Nitrite	Urobilinogen
Bilirubin	Hemoglobin (HGB)	pH	
Blood	Ketones	Protein (Total)	
Creatinine	Leukocytes	Protein-to-Creatinine Ratio	

qUAntify Plus Control



- Choose from a conical 12 mL tube or 120 mL flip top squeeze bottle
- Both options can be used for urine dipstick and microscopic testing
- Liquid
- 2 year shelf life at 2-8°C
- 31 day open-vial stability at 2-25°C for most analytes

POCT methods listed

Ordering Information

Catalog #	Description
995	Bilevel (5 of each level)/Tube 10 x 12 mL
995X	Bilevel MiniPak (1 of each level)/Tube. 2 x 12 mL
962	Bilevel (2 of each level)/Bottle 4 x 120 mL
962X	Bilevel MiniPak (1 of each level)/Bottle. 2 x 120 mL

Analytes

Albumin	Glucose	Microscopics (RBC, WBC, Crystals, Casts)	Specific Gravity
Albumin-to-Creatinine Ratio	hCG (Pregnancy)	Nitrite	Urobilinogen
Ascorbic Acid*	Hemoglobin (HGB)	pH	*Tested and found to be negative.
Bilirubin	Ketones	Protein (Total)	
Blood	Leukocytes	Protein-to-Creatinine Ratio	
Creatinine	Microalbumin		

Liquichek Urinalysis Control



- Ready-to-use urine control formulated to monitor the precision of dipstick and microscopic analyzers
- Available in 12 mL glass vials
- Liquid, human urine based
- 2.5 year shelf life at 2-8°C
- 30 day open-vial stability at 2-25°C

Ordering Information

Catalog #	Description
435	Bilevel (6 of each level) 12 x 12 mL
436	Level 1 12 x 12 mL
437	Level 2 12 x 12 mL
435X	Bilevel MiniPak (1 of each level) 2 x 12 mL

Analytes

Bilirubin	Glucose	Microscopics (RBC, WBC, Crystals, Casts)	Protein (Total)
Blood	hCG (Pregnancy)	Nitrite	Protein-to-Creatinine Ratio
Clarity	Ketones	Osmolality	Specific Gravity
Color	Leukocytes	pH	Urobilinogen
Creatinine	Microalbumin		

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Dispenser Tips for Screw Caps



- For controls fitted with a screw cap closure
- Disposable plastic tips enable fluid delivery directly from glass vial
- Reduces risk of product contamination that may be introduced via pipetting
- Easy and convenient to use

Ordering Information

Catalog #	Description
987000	Dispenser Tips for Screw Caps 100/Package

Validated for use with Liquichek Therapeutic Drug Monitoring Control (TDM) (catalog numbers 724-727 and 724X), Liquid Assayed Multiquel (catalog numbers 694-696 and 695X), Liquid Unassayed Multiquel (catalog numbers 697-699 and 698X) and Lyphocheck Immunology Plus Control (catalog numbers 430, 430X).

Dispenser Tips for Crimp Seals



- For controls fitted with a metal crimp seal and stopper
- Plastic tips allow the glass vial to be used as a dropper bottle
- Easy and convenient to use

Ordering Information

Catalog #	Description
588	Dispenser Tips for Crimp Seals 50/Package

Please use with caution; Cannabinoid Metabolites may bind to certain types of plastic and glass.

Blood Gas Ampule Breaker



- Designed to reduce the risk of injury when opening blood gas ampules
- Easy and convenient to use
- Reusable

Ordering Information

Catalog #	Description
587	Blood Gas Ampule Breaker 1 each

Quality Control Tray

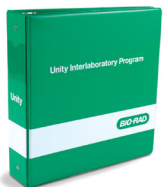


- 28-position durable acrylic tray to prepare or transport controls
- Accommodates up to 10 large and 18 standard sized control vials or dropper bottles
- Dimensions: 10" x 9.5" x 1.25" (25.4 cm x 24.1 cm x 3.2 cm)

Ordering Information

Catalog #	Description
12000846	Quality Control Tray 1 unit

Unity™ Binder



- Keep your Unity Interlaboratory Program reports organized and readily available
- 2 inch Capacity
- 3-Ring for Letter Size (8.5 x 11 inches)
- Monthly Tabs Provided

To request a binder, please call your local Bio-Rad office

Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

See page number

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ✱ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

	111	111	112	117	119	79	67	63	62	38	37	41	68	85	44	44	45	45	85	66	39	63	46	45	44	64	59	60	64	65	
	See Blood Gas Chart	See Cardiac Assessment Chart	See Hematology Chart	See Infectious Disease/Serology Chart	See IntelliQ Chart	Urine Toxicology (See Product Pages)	Allergen sIgE	AMH	Anemia	Assayed Chemistry	Assayed Multiqua (Liquid)	Assayed Multiqua Premium	Autoimmune	Benzo/TCA-Set B	Coagulation (Lyphocheck)	D-dimer	Diabetes (Liquicheck)	Diabetes (Lyphocheck)	Drug Free Serum	Elevated CRP	Ethano/Ammonia (Liquicheck)	Fertility	Hemoglobin A1C Linearity Set	Hemoglobin A ₂	Hemostasis	Hypertension Markers	Immunoassay Plus (Liquicheck)	Immunoassay Plus (Lyphocheck)	Immunoassay Premium	Immunology (Liquicheck)	
3-Methoxytyramine																															
5-Aminolevulinic Acid (ALA)																															
5-HIAA																															
6-Monoacetylmorphine						■																									
11-Ketoetiocholanolone																															
11-Nor-Δ-9-THC-9-COOH						●																									
11-Ketoandrosterone																															
11-Deoxycortisol																												▲	●		
11-β-Hydroxy-Androsterone																															
11-β-Hydroxy-Etiocholanolone																															
17-α-Hydroxyprogesterone																												●	●		
17-Hydroxycorticosteroids																															
17-Ketosteroids																															
25-Hydroxy Vitamin D						●																						▲	○	●	
α-1-Acid Glycoprotein						●						○																			●
α-1-Antitrypsin						○				○	○	○																			●
α-1-Globulin						○					○																				
α-2-Globulin						○					○																				
α-2-Macroglobulin																															●
αHBDH										□	□	□																			
β-2-Microglobulin						●						○																		●	●
β-Globulin						○																									
β-hCG						●																									
γ-Globulin						○					○																				
Acetaminophen						●				●	●	●								▶								●	●	●	
Acetone																															
Acid Phosphatase (Total)						●				●	●	●																			
ACTH																											●		●		
Activated Partial Thromboplastin Time															●											●					
Active Vitamin B ₁₂																															
ADH																											▲				
ADNase B																															●
ALA																															
Albumin		●				●				●	●	●																			●
Albumin (BCG)						●																									
Albumin (BCP)						●																									
Albumin-to-Creatinine Ratio																															
Albumin-to-Globulin Ratio						○																									
Alcohol (Ethanol)						●					●	●										●									
Alcohol (Isopropanol)																															
Alcohol (Methanol)																															
Aldolase																															
Aldosterone																											●	●	●		
Alkaline Phosphatase (ALP)						●				●	●	●																			
Alpha-Hydroxyalprazolam						●																									
Alphafetoprotein (AFP)						●				○													●					●	●	●	
Alprazolam														●							▶										
ALT/SGPT						●				●	●	●																			
Aluminum																															
Amikacin						●					●	●																●	●	●	
Amiodarone																												▲	▲		
Amitriptyline																												▲	●		

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Control Analyte Index

65	40	63	63	46	41	80	40	87	87	42	66	58	86	62	62	62	67	77	77	61	61	39	38	37	87	42	86	86	78	78	85						
Immunology Plus (Lyphocheck)	Lipids	Maternal Serum 1st Trimester	Maternal Serum II	Meter Trax	Microalbumin	Opiate	Pediatric	qAntify Advance	qAntify Plus	Quantitative Urine	Rheumatoid Factor	Sedimentation Rate	Serum Volatiles	Specialty Immunoassay (Liquichek)	Specialty Immunoassay Plus (Liquichek)	Specialty Immunoassay (Lyphocheck)	Spinal Fluid	TDM (Liquichek)	TDM (Lyphocheck)	Tumor Marker (Liquichek)	Tumor Marker Plus (Lyphocheck)	Unassayed Chemistry (Liquichek)	Unassayed Chemistry (Lyphocheck)	Unassayed Multiquel (Liquid)	Urinalysis (Liquichek)	Urine Chemistry (Liquichek)	Urine Metals	Whole Blood	Whole Blood Immunosuppressant (Liquid)	Whole Blood Immunosuppressant (Lyoph.)	Whole Blood Metals						
										●																							3-Methoxytyramine				
										●																		●					5-Aminolevulinic Acid (ALA)				
						●				●																							5-HIAA				
										▲																							6-Monoacetylmorphine				
										▲																								11-Ketoetiocolanolone			
										▲																								11-Nor-Δ-9-THC-9-COOH			
										▲																								11-Ketoandrosterone			
										▲																								11-Deoxycortisol			
										▲																								11-β-Hydroxy-Androsterone			
										▲																								11-β-Hydroxy-Etiocolanolone			
										▲																								17-α-Hydroxyprogesterone			
										▲																								17-Hydroxycorticosteroids			
										▲																								17-Ketosteroids			
														●	●	●																		25-Hydroxy Vitamin D			
●																																		α-1-Acid Glycoprotein			
●																								○	○									α-1-Antitrypsin			
																		●																	α-1-Globulin		
																		●																	α-2-Globulin		
●																																			α-2-Macroglobulin		
●																																			αHBDH		
																																				β-2-Microglobulin	
																																				β-Globulin	
																																				β-hCG	
																																				γ-Globulin	
																																				Acetaminophen	
														●																						Acetone	
																																				Acid Phosphatase (Total)	
																																				ACTH	
																																				Activated Partial Thromboplastin Time	
																																				Active Vitamin B ₁₂	
																																				ADH	
																																				ADNase B	
																																				ALA	
●																																				Albumin	
																																				Albumin (BCG)	
																																				Albumin (BCP)	
																																				Albumin-to-Creatinine Ratio	
																																				Albumin-to-Globulin Ratio	
																																				Alcohol (Ethanol)	
																																				Alcohol (Isopropanol)	
																																				Alcohol (Methanol)	
																																				Aldolase	
																																				Aldosterone	
																																					Alkaline Phosphatase (ALP)
																																				Alpha-Hydroxyalprazolam	
																																				Alphafetoprotein (AFP)	
																																				Alprazolam	
																																				ALT/SGPT	
																																				Aluminum	
																																				Amikacin	
																																				Amiodarone	
																																				Amitriptyline	

See page number

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ✦ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

Refer to myInserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

See page number

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ✱ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

	111	111	112	117	119	79	67	63	62	38	37	41	68	85	44	44	45	45	85	66	39	63	46	45	44	64	59	60	64	65	
	See Blood Gas Chart	See Cardiac Assessment Chart	See Hematology Chart	See Infectious Disease/Serology Chart	See IntellIQ Chart	Urine Toxicology (See Product Pages)	Allergen sigE	AMH	Anemia	Assayed Chemistry	Assayed Multiqua (Liquit)	Assayed Multiqua Premium	Autoimmune	Benzo/TCA-Set B	Coagulation (Lyphocheck)	D-dimer	Diabetes (Liquicheck)	Diabetes (Lyphocheck)	Drug Free Serum	Elevated CRP	Ethano/Ammonia (Liquicheck)	Fertility	Hemoglobin A1C Linearity Set	Hemoglobin A ₂	Hemostasis	Hypertension Markers	Immunoassay Plus (Liquicheck)	Immunoassay Plus (Lyphocheck)	Immunoassay Premium	Immunology (Liquicheck)	
Ammonia					●																●										
Amobarbital						●																									
AMP (Cyclic)																															
Amphetamines						●														▶											
Amylase					●					●	●	●																			
Amylase (Alpha)										●																					
Amylase (Pancreatic)					●					●	●	●																			
ANA, Homogeneous Pattern													●																		
Androstenedione																											●	●			
Androsterone																												▲	□		
Angiotensin I																															
Anti-CCP												○																			●
Anti-Cytomegalovirus (CMV)				●																											
Anti-HAV				●																											
Antidiuretic Hormone																											▲				
Anti-HBc				●																											
Anti-HBe				●																											
Anti-HBs				●																											
Anti-HCV				●																											
Anti-HIV-1				●																											
Anti-HIV-2				●																											
Anti-HTLV-I				●																											
Anti-HTLV-I/II				●																											
Anti-HTLV-II				●																											
Anti-Mitochondrial													●																		
Anti-Mullerian Hormone								●																							
Anti-nDNA													●																		
Anti-Smooth Muscle													●																		
Anti-Thyroglobulin (Anti-Tg)					●																						▲	▲			
Anti-Thyroperoxidase (Anti-TPO)					●																						▲	▲			
Anti-Treponema pallidum (Syphilis)				●																											
Anti-TSH Receptor (TRAb)																															
Antimony																															
Antistreptolysin O (ASO)					●							○																			●
Antithrombin III (AT III)															●										●						●
Apolipoprotein A-1					○					○	○	○																			●
Apolipoprotein B					○					○	○	○																			●
Arsenic																															
Ascorbic Acid																															
AST/SGOT					●					●	●	●																			
Autoimmune Negative													●																		
Barbiturates						●																									
Basophils (BASO)			●																												
Benzodiazepines						●																									
Benzoyllecgonine (Cocaine Metabolite)						●																									
Bile Acids												●																			
Bilirubin																															
Bilirubin (Direct)					●					●	●	●																			
Bilirubin (Indirect)										●																					
Bilirubin (Neonatal)					●						●	●																			
Bilirubin (Total)					●					●	●	●																			
Blood																															

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Control Analyte Index

65	40	63	63	46	41	80	40	87	87	42	66	58	86	62	62	62	67	77	77	61	61	39	38	37	87	42	86	86	78	78	85	See page number
Immunology Plus (Lyphocheck)	Lipids	Maternal Serum 1st Trimester	Maternal Serum II	Meter Trax	Microalbumin	Opiate	Pediatric	qUAntify Advance	qUAntify Plus	Quantitative Urine	Rheumatoid Factor	Sedimentation Rate	Serum Volatiles	Specialty Immunoassay (Liquichek)	Specialty Immunoassay Plus (Liquichek)	Specialty Immunoassay (Lyphocheck)	Spinal Fluid	TDM (Liquichek)	TDM (Lyphocheck)	Tumor Marker (Liquichek)	Tumor Marker Plus (Lyphocheck)	Unassayed Chemistry (Liquichek)	Unassayed Chemistry (Lyphocheck)	Unassayed Multiqua (Liquid)	Urinanalysis (Liquichek)	Urine Chemistry (Liquichek)	Urine Metals	Whole Blood	Whole Blood Immunosuppressant (Liquid)	Whole Blood Immunosuppressant (Lyoph.)	Whole Blood Metals	
																																Ammonia
																																Amobarbital
																																AMP (Cyclic)
																																Amphetamines
																																Amylase
																																Amylase (Alpha)
																																Amylase (Pancreatic)
																																ANA, Homogeneous Pattern
																																Androstenedione
																																Androsterone
																																Angiotensin I
																																Anti-CCP
																																Anti-Cytomegalovirus (CMV)
																																Anti-HAV
																																Anti-HBc
																																Anti-HBe
																																Anti-HBs
																																Anti-HCV
																																Anti-HIV-1
																																Anti-HIV-2
																																Anti-HTLV-I
																																Anti-HTLV-I/II
																																Anti-HTLV-II
																																Anti-Mitochondrial
																																Anti-Mullerian Hormone
																																Anti-nDNA
																																Anti-Smooth Muscle
																																Anti-Thyroglobulin (Anti-Tg)
																																Anti-Thyroperoxidase (Anti-TPO)
																																Anti-Treponema pallidum (Syphilis)
																																Anti-TSH Receptor (TRAb)
																																Antimony
																																Antistreptolysin O (ASO)
																																Anti-thrombin III (AT III)
																																Apolipoprotein A-1
																																Apolipoprotein B
																																Arsenic
																																Ascorbic Acid
																																AST/SGOT
																																Autoimmune Negative
																																Barbiturates
																																Basophils (BASO)
																																Benzodiazepines
																																Benzoylcegonine (Cocaine Metabolite)
																																Bile Acids
																																Bilirubin
																																Bilirubin (Direct)
																																Bilirubin (Indirect)
																																Bilirubin (Neonatal)
																																Bilirubin (Total)
																																Blood
																																BNP

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ✳ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

See page number

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ★ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

	111	111	112	117	119	79	67	63	62	38	37	41	68	85	44	44	45	45	85	66	39	63	46	45	44	64	59	60	64	65
	See Blood Gas Chart	See Cardiac Assessment Chart	See Hematology Chart	See Infectious Disease/Serology Chart	See IntelliQ Chart	Urine Toxicology (See Product Pages)	Allergen sigE	AMH	Anemia	Assayed Chemistry	Assayed Multiqua (Liquit)	Assayed Multiqua Premium	Autoimmune	Benzo/TCA-Set B	Coagulation (Lyphocheck)	D-dimer	Diabetes (Liquicheck)	Diabetes (Lyphocheck)	Drug Free Serum	Elevated CRP	Ethano/Ammonia (Liquicheck)	Fertility	Hemoglobin A1C Linearity Set	Hemoglobin A ₂	Hemostasis	Hypertension Markers	Immunoassay Plus (Liquicheck)	Immunoassay Plus (Lyphocheck)	Immunoassay Premium	Immunology (Liquicheck)
BNP		●			●																									
Borrelia burgdorferi IgG				●																										
Buprenorphine																														
Butalbital					●																									
C-Peptide					●																							●		
C1 Inhibitor																														●
CA 15-3					●																									●
CA 19-9					●																									●
CA 27.29					●																									●
CA 50*																														
CA 72-4*					●																									
CA 125					●																									●
Cadmium																														
Caffeine												●								▶								●	●	●
Calcitonin																														
Calcium					●					●	●	●																		
Calcium (GPC)					●																									
Calcium (Ionized)	●				●					○	●	●																		
Cannabinoid Metabolite						●																								
Cannabinoids						●														▶										
Carbamazepine					●					●	●	●								▶								●	●	●
Carbamazepine (Free)																												●	●	●
CASA*																														
Casts																														
CEA					●					○																		●	●	●
Cellular Hemoglobin Concentration Mean			●																											
Cellular Hemoglobin Concentration Mean of Gated Red Blood			●																											
Cellular Hemoglobin Concentration Mean of Reticulocytes			●																											
Ceruloplasmin					○					○	○	○																		●
Chlamydia trachomatis				●																										
Chlamydia trachomatis IgG				●																										
Chloramphenicol																				▶								●	●	
Chlordiazepoxide																				▶										
Chloride	●				●					●	●	●																		
Cholesterol (HDL)					●					●	●	●																		
Cholesterol (LDL)					●					●	●	●																		
Cholesterol (Total)					●					●	●	●																		
Cholinesterase					○					○	○	○																		
Chromium																														
Cimetidine																					▶									
CK (Total)					●																									
CK-MB Isoenzyme		●			●						▲																	●		
Clarity																														
Clomipramine															●					▶										
Clonazepam																				▶										
Clostridium difficile (C diff.)				●																										
CO ₂					●					●	●	●																		
Cobalt																														
Cocaine					●															▶										
Cocaine Metabolite					●																									

*International use only.
Refer to myInSerts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

65	40	63	63	46	41	80	40	87	87	42	66	58	86	62	62	62	67	77	77	61	61	39	38	37	87	42	86	86	78	78	85	See page number			
Immunology Plus (Lyphocheck)	Lipids	Maternal Serum 1st Trimester	Maternal Serum II	Meter Trax	Microalbumin	Opiate	Pediatric	qUAntify Advance	qUAntify Plus	Quantitative Urine	Rheumatoid Factor	Sedimentation Rate	Serum Volatiles	Specialty Immunoassay (Liquichek)	Specialty Immunoassay Plus (Liquichek)	Specialty Immunoassay (Lyphocheck)	Spinal Fluid	TDM (Liquichek)	TDM (Lyphocheck)	Tumor Marker (Liquichek)	Tumor Marker Plus (Lyphocheck)	Unassayed Chemistry (Liquichek)	Unassayed Chemistry (Lyphocheck)	Unassayed Multiqua (Liquid)	Urinalysis (Liquichek)	Urine Chemistry (Liquichek)	Urine Metals	Whole Blood	Whole Blood Immunosuppressant (Liquid)	Whole Blood Immunosuppressant (Lyphocheck)	Whole Blood Metals				
						●									●	●																Borrelia burgdorferi IgG			
																																Buprenorphine			
																																Butalbital			
																																C-Peptide			
																																C1 Inhibitor			
																						●	●									CA 15-3			
																					●	●										CA 19-9			
																					●	●										CA 27.29			
																					●	●										CA 50*			
																					●	●										CA 72-4*			
																					●	●										CA 125			
																												●				●	Cadmium		
							●											●	●														Caffeine		
																						●											Calcitonin		
											●												●	●	●		●						Calcium		
																																	Calcium (CPC)		
																							▲	○	●								Calcium (Ionized)		
																																	Cannabinoid Metabolite		
																																		Cannabinoids	
																							●	●	●									Carbamazepine	
																							●	●										Carbamazepine (Free)	
																																		CASA*	
																						▲												Casts	
																							●	●		○								CEA	
																																		Cellular Hemoglobin Concentration Mean	
																																		Cellular Hemoglobin Concentration Mean of Gated Red Blood	
																																		Cellular Hemoglobin Concentration Mean of Reticu-	
●																																		Ceruloplasmin	
																																			Chlamydia trachomatis
																																			Chlamydia trachomatis IgG
																																			Chloramphenicol
																																			Chlordiazepoxide
																																			Chloride
	●																																		Cholesterol (HDL)
	●																																		Cholesterol (LDL)
	●																																		Cholesterol (Total)
																																			Cholinesterase
																																			Chromium
																																			Cimetidine
																																			CK (Total)
																																			CK-MB Isoenzyme
																																			Clarity
																																			Clomipramine
																																			Clonazepam
																																			Clostridium difficile (C diff.)
																																			CO ₂
																																			Cobalt
																																			Cocaine
																																			Cocaine Metabolite
																																			Codeine

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- * Values not for diagnostic use
- ▲ No claim is made for expected value or stability

*International use only.

Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

See page number

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ✱ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

	111	111	112	117	119	79	67	63	62	38	37	41	68	85	44	44	45	45	85	66	39	63	46	45	44	64	59	60	64	65		
	See Blood Gas Chart	See Cardiac Assessment Chart	See Hematology Chart	See Infectious Disease/Serology Chart	See IntelliQ Chart	Urine Toxicology (See Product Pages)	Allergen sigE	AMH	Anemia	Assayed Chemistry	Assayed Multiqua (Liquit)	Assayed Multiqua Premium	Autoimmune	Benzo/TCA-Set B	Coagulation (Lyphocheck)	D-dimer	Diabetes (Liquicheck)	Diabetes (Lyphocheck)	Drug Free Serum	Elevated CRP	Ethano/Ammonia (Liquicheck)	Fertility	Hemoglobin A1C Linearity Set	Hemoglobin A ₂	Hemostasis	Hypertension Markers	Immunoassay Plus (Liquicheck)	Immunoassay Plus (Lyphocheck)	Immunoassay Premium	Immunology (Liquicheck)		
Codeine						●													▶													
Color																																
Complement C1q																																
Complement C3					○					○	○																				●	
Complement C4					○					□	○	○																			●	
Copper					○					○	○	○																				
Coproporphyrin																																
Cortisol						●				●	●	●															●	●	●	●		
Creatine Kinase (CK)		●			●					●	●	●																				
Creatinine					●	◆				●	●	●																				
CRP		●			●							●									●										●	
Crystals																																
Cyclosporine																				▶								▲	▲			
Cyfra 21-1*					●																											
Cystatin C					○							○																			○	
Cystine																																
Cytomegalovirus (CMV) IgG				●																												
Cytomegalovirus (CMV) IgM				●																												
D1 House Dust Mite (D. pteronyssinus)							●																									
D2 House Dust Mite (D. farinae)							●																									
d-Amphetamine						●																										
D-dimer																●																
d-Methamphetamine						●																										
Dehydroepiandrosterone (DHEA)																											▲	▲				
Desalkylflurazepam																				▶												
Desipramine																				▶								●	●			
Desmethyldoxepin																				▶												
DHEA Sulfate						●																						●	●	●		
Diazepam																				▶												
Digitoxin		●			●															▶		●										
Digoxin					●					●	●	●								▶								●	●	●		
Disopyramide												●								▶								●	●	●		
Dopamine																																
Doxepin																				▶												
E1 Cat Dander (Felis domesticus)							●																									
E3 Horse Dander (Equus caballus)							●																									
E5 Dog Dander (Canis familiaris)							●																									
EBNA IgG				●																												
EBV EA-D IgG				●																												
Ecstasy (MDMA)						●																										
EDDP (Methadone Metabolite)						■																										
Eosinophils (EOS)			●																													
Epinephrine																																
Epstein-Barr Virus (VCA) IgG				●																												
Epstein-Barr Virus (VCA) IgM				●																												
Erythrocyte Sedimentation Rate (ESR)			●																													
Erythropoietin (EPO)					●																											
Estradiol					●																		●					●	●	●		
Estriol (Free)																						●						●	●	●		
Estriol (Total)																						▲						▲	▲			

*International use only.
Refer to myInSerts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

65	40	63	63	46	41	80	40	87	87	42	66	58	86	62	62	62	67	77	77	61	61	39	38	37	87	42	86	86	78	78	85	See page number			
Immunology Plus (Lyphocheck)	Lipids	Maternal Serum 1st Trimester	Maternal Serum II	Meter Trax	Microalbumin	Opiate	Pediatric	qUAntify Advance	qUAntify Plus	Quantitative Urine	Rheumatoid Factor	Sedimentation Rate	Serum Volatiles	Specialty Immunoassay (Liquichek)	Specialty Immunoassay Plus (Liquichek)	Specialty Immunoassay (Lyphocheck)	Spinal Fluid	TDM (Liquichek)	TDM (Lyphocheck)	Tumor Marker (Liquichek)	Tumor Marker Plus (Lyphocheck)	Unassayed Chemistry (Liquichek)	Unassayed Chemistry (Lyphocheck)	Unassayed Multiqua (Liquid)	Urinalysis (Liquichek)	Urine Chemistry (Liquichek)	Urine Metals	Whole Blood	Whole Blood Immunosuppressant (Liquid)	Whole Blood Immunosuppressant (Lyoph.)	Whole Blood Metals				
																																Color			
▲																																	Complement C1q		
●																																	Complement C3		
●																																	Complement C4		
																																	Copper		
																																	Coproporphyrin		
																																	Cortisol		
																																	Creatine Kinase (CK)		
																																	Creatinine		
●	●				●			●	●	●																							CRP		
																																	Crystals		
																																	Cyclosporine		
																																		Cyfra 21-1*	
																																		Cystatin C	
																																		Cystine	
																																		Cytomegalovirus (CMV) IgG	
																																		Cytomegalovirus (CMV) IgM	
																																		D1 House Dust Mite (D. pteronyssinus)	
																																		D2 House Dust Mite (D. farinae)	
																																		d-Amphetamine	
																																		D-dimer	
																																		d-Methamphetamine	
																																		Dehydroepiandrosterone (DHEA)	
																																		Desalkylflurazepam	
																																		Desipramine	
																																		Desmethyldoxepin	
																																		DHEA Sulfate	
																																		Diazepam	
																																		Digitoxin	
																																		Digoxin	
																																		Disopyramide	
																																		Dopamine	
																																			Doxepin
																																		E1 Cat Dander (Felis domesticus)	
																																		E3 Horse Dander (Equus caballus)	
																																		E5 Dog Dander (Canis familiaris)	
																																		EBNA IgG	
																																		EBV EA-D IgG	
																																		Ecstasy (MDMA)	
																																		EDDP (Methadone Metabolite)	
																																		Eosinophils (EOS)	
																																		Epinephrine	
																																		Epstein-Barr Virus (VCA) IgG	
																																		Epstein-Barr Virus (VCA) IgM	
																																		Erythrocyte Sedimentation Rate (ESR)	
																																		Erythropoietin (EPO)	
																																		Estradiol	
																																		Estriol (Free)	
																																		Estriol (Total)	
																																		Estrogen (Total)	

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ✳ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

*International use only.

Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

See page number

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ✱ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

	111	111	112	117	119	79	67	63	62	38	37	41	68	85	44	44	45	45	85	66	39	63	46	45	44	64	59	60	64	65	
	See Blood Gas Chart	See Cardiac Assessment Chart	See Hematology Chart	See Infectious Disease/Serology Chart	See IntellIQ Chart	Urine Toxicology (See Product Pages)	Allergen sigE	AMH	Anemia	Assayed Chemistry	Assayed Multiqua (Liquit)	Assayed Multiqua Premium	Autoimmune	Benzo/TCA-Set B	Coagulation (Lyphocheck)	D-dimer	Diabetes (Liquicheck)	Diabetes (Lyphocheck)	Drug Free Serum	Elevated CRP	Ethano/Ammonia (Liquicheck)	Fertility	Hemoglobin A1C Linearity Set	Hemoglobin A ₂	Hemostasis	Hypertension Markers	Immunoassay Plus (Liquicheck)	Immunoassay Plus (Lyphocheck)	Immunoassay Premium	Immunology (Liquicheck)	
Estrogen (Total)																															
Ethanol					●	●						●	●									●									
Ethosuximide																			▶									●	●	●	
Ethylene Glycol																															
Etioccholanolone																															
Everolimus																															
F1 Egg White (Gallus spp.)							●																								
F2 Cow Milk (Bos. spp.)							●																								
F13 Peanut (Arachis hypogaea)							●																								
Factors II, V, VII, VIII, IX, X, XI and XII																									●						
Fentanyl																															
Ferritin					○				●		○	○											●					●	●	●	●
Fibrinogen															●										●						
Flecainide																				▶								●	●		
Fluoride																															
Folate					●				●																			●	●	●	
Folate (RBC)																															
Fructosamine																											□	○			
FSH					●																		●					●	●	●	
G2 Bermuda Grass (Cynodon dactylon)							●																								
G3 Cocksfoot/Orchard Grass (Dactylis)							●																								
G6 Timothy Grass (Phleum pratense)							●																								
Gastrin																													●		
Galectin-3		●			●																										
Gentamicin					●					●	●	●								▶								●	●	●	
GGT					●					●	●	●																			
GLDH*										○	●																				
Globulin		●								○	○	□																			
Glucagon																												▲			
Glucose	●									●	●	●																			
Glucose (Hexokinase)					●																										
Glutethimide						■															▶										
Granulocytes (GRAN)			●																												
Haloperidol																					▶										
Haptoglobin					○					○	○	○																		●	
HAV-IgM				●																											
HBc-IgM				●																											
HBeAg				●																											
HBsAg				●																											
hCG					●																		●					●	●		
hCG (Intact)					●																										
hCG (Pregnancy)																															
hCG (Total)					●																							●	●	●	
hCG-β Subunit										○													●					●	●		
hCG-β Subunit (Free)																												●	●		
HDL Cholesterol										●	●	●																			
HE4 (Human Epididymis Protein 4)					●																										
Helicobacter pylori IgG				●																											
Hematocrit (HCT)			●																												
Hemoglobin (HGB)			●																												

*International use only.
Refer to myInSerts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

65	40	63	63	46	41	80	40	87	87	42	66	58	86	62	62	62	67	77	77	61	61	39	38	37	87	42	86	86	78	78	85	See page number					
Immunology Plus (Lyphocheck)	Lipids	Maternal Serum 1st Trimester	Maternal Serum II	Meter Trax	Microalbumin	Opiate	Pediatric	qUAntify Advance	qUAntify Plus	Quantitative Urine	Rheumatoid Factor	Sedimentation Rate	Serum Volatiles	Specialty Immunoassay (Liquichek)	Specialty Immunoassay Plus (Liquichek)	Specialty Immunoassay (Lyphocheck)	Spinal Fluid	TDM (Liquichek)	TDM (Lyphocheck)	Tumor Marker (Liquichek)	Tumor Marker Plus (Lyphocheck)	Unassayed Chemistry (Liquichek)	Unassayed Chemistry (Lyphocheck)	Unassayed Multiqua (Liquid)	Urinalysis (Liquichek)	Urine Chemistry (Liquichek)	Urine Metals	Whole Blood	Whole Blood Immunosuppressant (Liquid)	Whole Blood Immunosuppressant (Lyphocheck)	Whole Blood Metals						
													●																				Ethanol				
													●					●	●														Ethosuximide				
																																	Ethylene Glycol				
										▲																							Etiocholanolone				
																																	Everolimus				
																																	F1 Egg White (Gallus spp.)				
																																	F2 Cow Milk (Bos. spp.)				
																																	F13 Peanut (Arachis hypogaea)				
																																	Factors II, V, VII, VIII, IX, X, XI and XII				
								●																									Fentanyl				
																					●	●			○								Ferritin				
																																		Fibrinogen			
																			▲	▲														Flecainide			
																																		Fluoride			
																																		Folate			
																																		Folate (RBC)			
																																		Fructosamine			
																																			FSH		
																																			G2 Bermuda Grass (Cynodon dactylon)		
																																			G3 Cocksfoot/Orchard Grass (Dactylis)		
																																			G6 Timothy Grass (Phleum pratense)		
																																			Gastrin		
																																			Galectin-3		
																																			Gentamicin		
																																				GGT	
																																				GLDH*	
																																				Globulin	
																																				Glucagon	
																																				Glucose	
																																				Glucose (Hexokinase)	
																																				Glutethimide	
																																				Granulocytes (GRAN)	
																																				Haloperidol	
																																					Haloproglobin
																																					HAV-IgM
																																					HBeAg
																																					HBeAg
																																					hCG
																																					hCG (Intact)
																																					hCG (Pregnancy)
																																					hCG (Total)
																																					hCG-β Subunit
																																					hCG-β Subunit (Free)
																																					HDL Cholesterol
																																					HE4 (Human Epididymis Protein 4)
																																					Helicobacter pylori IgG
																																					Hematocrit (HCT)
																																					Hemoglobin (HGB)
																																					Hemoglobin (Total)

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- * Values not for diagnostic use
- ▲ No claim is made for expected value or stability

*International use only.

Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

See page number

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ✱ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

	111	111	112	117	119	79	67	63	62	38	37	41	68	85	44	44	45	45	85	66	39	63	46	45	44	64	59	60	64	65		
	See Blood Gas Chart	See Cardiac Assessment Chart	See Hematology Chart	See Infectious Disease/Serology Chart	See IntellIQ Chart	Urine Toxicology (See Product Pages)	Allergen sigE	AMH	Anemia	Assayed Chemistry	Assayed Multiqua (Liquit)	Assayed Multiqua Premium	Autoimmune	Benzo/TCA-Set B	Coagulation (Lyphocheck)	D-dimer	Diabetes (Liquicheck)	Diabetes (Lyphocheck)	Drug Free Serum	Elevated CRP	Ethano/Ammonia (Liquicheck)	Fertility	Hemoglobin A1C Linearity Set	Hemoglobin A ₂	Hemostasis	Hypertension Markers	Immunoassay Plus (Liquicheck)	Immunoassay Plus (Lyphocheck)	Immunoassay Premium	Immunology (Liquicheck)		
Hemoglobin (Total)																		●														
Hemoglobin (Total Glycated)																		▲														
Hemoglobin A1																		▲														
Hemoglobin A1C																	●	●						●								
Hemoglobin A ₂																								●								
Hemoglobin Concentration Distribution			●																													
Hemoglobin F																		●						●								
Hemoglobin S																								●								
Hepatitis B Virus (HBV)				●																												
Hepatitis C Virus (HCV)				●																												
HER2/neu					●																											
Heroin Metabolite						●																										
Herpes Simplex Virus Type 1 (HSV-1) IgG				●																												
Herpes Simplex Virus Type 1 (HSV-1) IgM				●																												
Herpes Simplex Virus Type 2 (HSV-2) IgG				●																												
Herpes Simplex Virus Type 2 (HSV-2) IgM				●																												
hGH																											●	●	●			
Hippuric Acid																																
HIV-1 Ag				●																												
Homocysteine		●			●																											
Homogenous Pattern												●																				
hs-CRP		●																														
Human Immunodeficiency Virus Type 1				●																												
Human Papillomavirus (HPV)				●																												
HVA																																
Hydromorphone																				▶												
Hydroxyproline (Free)																																
Hydroxyproline (Total)																																
Ibuprofen																				▶								●				
Imipramine						●														▶								▲	●			
Immature Reticulocyte Fraction (IRF)			●																													
Immunoglobulin A (IgA)					○					○	○	○																○	○	○	●	
Immunoglobulin E (IgE)				●																								●	●	●	●	
Immunoglobulin G (IgG)					○					○	○	○																○	○	○	●	
Immunoglobulin G Subclasses 1-4																															●	
Immunoglobulin M (IgM)					○					○	○	○																○	○	○	●	
Immunoreactive Trypsinogen (IRT)																													▲			
Inhibin A																																
Insulin					●																							●	●	●		
Insulin Growth Factor I (IGF-I)/Somatomedin C					●																							▲	○			
Interleukin-6 (IL-6)																																
Iron					●				●	●	●	●																○	○			
Iron (TIBC)					○				●	○	○	○																▲	○			
Iron (UIBC)					○				●	○	○	○																				
IRT																													▲			
Isopropanol (Isopropyl Alcohol)																																
Kappa Light Chain					●							○																			●	
Ketones																																
Lactate (Lactic Acid)	●				●					●	●	●																				
Lactate Dehydrogenase (LDH)		●			●					●	●	●																				
Lambda Light Chain					●							○																			●	
LAP Arylamidase					●					○	○	□																				

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Control Analyte Index

65	40	63	63	46	41	80	40	87	87	42	66	58	86	62	62	62	67	77	77	61	61	39	38	37	87	42	86	86	78	78	85	See page number			
Immunology Plus (Lipchocek)	Lipids	Maternal Serum 1st Trimester	Maternal Serum II	Meter Trax	Microalbumin	Opiate	Pediatric	qUAntify Advance	qUAntify Plus	Quantitative Urine	Rheumatoid Factor	Sedimentation Rate	Serum Volatiles	Specialty Immunoassay (Lipchocek)	Specialty Immunoassay Plus (Lipchocek)	Specialty Immunoassay (Lipchocek)	Spinal Fluid	TDM (Lipchocek)	TDM (Lipchocek)	Tumor Marker (Lipchocek)	Tumor Marker Plus (Lipchocek)	Unassayed Chemistry (Lipchocek)	Unassayed Chemistry (Lipchocek)	Unassayed Multiqua (Liquid)	Urinanalysis (Lipchocek)	Urine Chemistry (Lipchocek)	Urine Metals	Whole Blood	Whole Blood Immunosuppressant (Liquid)	Whole Blood Immunosuppressant (Lyoph.)	Whole Blood Metals				
																																	Hemoglobin (Total Glycated)		
																																		Hemoglobin A1	
																																		Hemoglobin A1C	
																																		Hemoglobin A ₂	
																																		Hemoglobin Concentration Distribution	
																																		Hemoglobin F	
																																		Hemoglobin S	
																																		Hepatitis B Virus (HBV)	
																																		Hepatitis C Virus (HCV)	
																																		HER2/neu	
																																		Heroin Metabolite	
																																		Herpes Simplex Virus Type 1 (HSV-1) IgG	
																																		Herpes Simplex Virus Type 1 (HSV-1) IgM	
																																		Herpes Simplex Virus Type 2 (HSV-2) IgG	
																																		Herpes Simplex Virus Type 2 (HSV-2) IgM	
																																		hGH	
																																		Hippuric Acid	
																																		HIV-1 Ag	
																																		Homocysteine	
																																		Homogenous Pattern	
																																			hs-CRP
																																			Human Immunodeficiency Virus Type 1
																																			Human Papillomavirus (HPV)
																																			HVA
																																			Hydromorphone
																																			Hydroxyproline (Free)
																																			Hydroxyproline (Total)
																																			Ibuprofen
																																			Imipramine
																																			Immature Reticulocyte Fraction (IRF)
																																			Immunoglobulin A (IgA)
																																			Immunoglobulin E (IgE)
																																			Immunoglobulin G (IgG)
																																			Immunoglobulin G Subclasses 1-4
																																			Immunoglobulin M (IgM)
																																			Immunoreactive Trypsinogen (IRT)
																																			Inhibin A
																																			Insulin
																																			Insulin Growth Factor I (IGF-I)/Somatomedin C
																																			Interleukin-6 (IL-6)
																																			Iron
																																			Iron (TIBC)
																																			Iron (UIBC)
																																			IRT
																																			Isopropanol (Isopropyl Alcohol)
																																			Kappa Light Chain
																																			Ketones
																																			Lactate (Lactic Acid)
																																			Lactate Dehydrogenase (LDH)
																																			Lambda Light Chain
																																			LAP Arylamidase
																																			Large Unstained Cells (LUC)

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ✱ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

See page number

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ✱ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

	111	111	112	117	119	79	67	63	62	38	37	41	68	85	44	44	45	45	85	66	39	63	46	45	44	64	59	60	64	65	
	See Blood Gas Chart	See Cardiac Assessment Chart	See Hematology Chart	See Infectious Disease/Serology Chart	See IntellIQ Chart	Urine Toxicology (See Product Pages)	Allergen sigE	AMH	Anemia	Assayed Chemistry	Assayed Multiqua (Liquit)	Assayed Multiqua Premium	Autoimmune	Benzo/TCA-Set B	Coagulation (Lyphocheck)	D-dimer	Diabetes (Liquicheck)	Diabetes (Lyphocheck)	Drug Free Serum	Elevated CRP	Ethano/Ammonia (Liquicheck)	Fertility	Hemoglobin A1C Linearity Set	Hemoglobin A ₂	Hemostasis	Hypertension Markers	Immunoassay Plus (Liquicheck)	Immunoassay Plus (Lyphocheck)	Immunoassay Premium	Immunology (Liquicheck)	
Large Unstained Cells (LUC)			●																												
LDH (Total)																															
LDL Cholesterol										●	●	●																			
Lead																															
Leukocytes			●																												
Lidocaine												●							▶								●	●	●		
Lipase					●					●	●	●																			
Lipoprotein (a)					●																										●
Lithium	●				●					●	●	●							▶								●	●	●		
Lormetazepam						●																									
LSD						●																									
Luteinizing Hormone (LH)					●																	●						●	●	●	
LYMPH (W-SCC & W-SCR)			●																												
Lymphocytes (LYMPH)			●																												
Lysergic Acid Diethylamide (LSD)						●																									
M3 Mold (Aspergillus fumigatus)							●																								
M6 Mold (Alternaria tenuis)							●																								
Magnesium	●				●					●	●	●																			
Mandelic Acid																															
Manganese																															
Marijuana Metabolite						●																									
MDA						●																									
MDEA						●																									
MDMA						●																									
Mean Corpuscular Hemoglobin (MCH)			●																												
Mean Corpuscular Hemoglobin Concentration			●																												
Mean Corpuscular Volume (MCV)			●																												
Mean Corpuscular Volume of Gated Red Blood			●																												
Mean Corpuscular Volume of Reticulocytes			●																												
Mean Hemoglobin Content of All Gated Red			●																												
Mean Hemoglobin Content of Reticulocytes			●																												
Mean Platelet Volume (MPV)			●																												
Meperidine*																			▶												
Mercury																															
Metanephrine																															
Metanephrine (Total)																															
Methadone																				▶											
Methadone Metabolite (EDDP)						■																									
Methamphetamine						●														▶											
Methanol (Methyl Alcohol)																															
Methaqualone						●														▶											
Methicillin Resistant Staphylococcus aureus				●																											
Methicillin Sensitive Staphylococcus aureus				●																											
Methotrexate												●								▶											
Methyprylon																				▶											
Microalbumin					●																										
Microscopics (RBC, WBC, Crystals, Casts)																															
MID			●																												
Mitochondrial												●																			
Mixed Cell Percent (MXD)			●																												

*International use only.

Refer to myInSerts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Control Analyte Index

65	40	63	63	46	41	80	40	87	87	42	66	58	86	62	62	62	67	77	77	61	61	39	38	37	87	42	86	86	78	78	85				
Immunology Plus (Lyphocheck)	Lipids	Maternal Serum 1st Trimester	Maternal Serum II	Meter Trax	Microalbumin	Opiate	Pediatric	qUantify Advance	qUantify Plus	Quantitative Urine	Rheumatoid Factor	Sedimentation Rate	Serum Volatiles	Specialty Immunoassay (Liquichek)	Specialty Immunoassay Plus (Liquichek)	Specialty Immunoassay (Lyphocheck)	Spinal Fluid	TDM (Liquichek)	TDM (Lyphocheck)	Tumor Marker (Liquichek)	Tumor Marker Plus (Lyphocheck)	Unassayed Chemistry (Liquichek)	Unassayed Chemistry (Lyphocheck)	Unassayed Multiqua (Liquid)	Urinanalysis (Liquichek)	Urine Chemistry (Liquichek)	Urine Metals	Whole Blood	Whole Blood Immunosuppressant (Liquid)	Whole Blood Immunosuppressant (Lyoph.)	Whole Blood Metals				
	●																●																LDH (Total)		
										●													●	●									LDL Cholesterol		
								●	●																								Lead		
																			●	●													Leukocytes		
																																	Lidocaine		
																							●	●	●									Lipase	
	○																																Lipoprotein (a)		
																							●	●	●									Lithium	
																																		Lormetazepam	
																																		Luteinizing Hormone (LH)	
																																		LSD	
																																		LYMPH (W-SCC & W-SCR)	
																																		Lymphocytes (LYMPH)	
																																		Lysergic Acid Diethylamide (LSD)	
																																		M3 Mold (Aspergillus fumigatus)	
																																		M6 Mold (Alternaria tenuis)	
							●			●																								Magnesium	
																																		Mandelic Acid	
																																		Manganese	
																																			Marijuana Metabolite
																																			MDA
																																			MDEA
																																			MDMA
																																			Mean Corpuscular Hemoglobin (MCH)
																																			Mean Corpuscular Hemoglobin Concentra-
																																			Mean Corpuscular Volume (MCV)
																																			Mean Corpuscular Volume of Gated Red Blood
																																			Mean Corpuscular Volume of Reticulo-
																																			Mean Hemoglobin Content of All Gated Red
																																			Mean Hemoglobin Content of Reticulo-
																																			Mean Platelet Volume (MPV)
								●																											Meperidine*
										●																									Mercury
										●																									Metanephrine
										●																									Metanephrine (Total)
																																			Methadone
																																			Methadone Metabolite (EDDP)
																																			Methamphetamine
																																			Methanol (Methyl Alcohol)
																																			Methaqualone
																																			Methicillin Resistant Staphylococcus
																																			Methicillin Sensitive Staphylococcus
																																			Methotrexate
																																			Methypylon
					●			●	●	●																									Microalbumin
																																			Microscopics (RBC, WBC, Crystals,
																																			MID
																																			Mitochondrial
																																			Mixed Cell Percent (MXD)
																																			Monocytes (MONO)

- See page number
- Analyte present
 - Analyte present in endogenous levels
 - Analyte present and no claims are made regarding performance or stability
 - Analyte present in endogenous levels and no claims are made regarding performance or stability
 - ▶ Tested and found to be negative
 - ◆ Specimen validity criteria
 - ✳ Values not for diagnostic use
 - ▲ No claim is made for expected value or stability

*International use only.

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

See page number

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ✱ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

	111	111	112	117	119	79	67	63	62	38	37	41	68	85	44	44	45	45	85	66	39	63	46	45	44	64	59	60	64	65
	See Blood Gas Chart	See Cardiac Assessment Chart	See Hematology Chart	See Infectious Disease/Serology Chart	See IntelliQ Chart	Urine Toxicology (See Product Pages)	Allergen sigE	AMH	Anemia	Assayed Chemistry	Assayed Multiqua (Liquit)	Assayed Multiqua Premium	Autoimmune	Benzo/TCA-Set B	Coagulation (Lyphocheck)	D-dimer	Diabetes (Liquicheck)	Diabetes (Lyphocheck)	Drug Free Serum	Elevated CRP	Ethano/Ammonia (Liquicheck)	Fertility	Hemoglobin A1C Linearity Set	Hemoglobin A ₂	Hemostasis	Hypertension Markers	Immunoassay Plus (Liquicheck)	Immunoassay Plus (Lyphocheck)	Immunoassay Premium	Immunology (Liquicheck)
Monocytes (MONO)			●																											
Morphine						●													▶											
Morphine (Free)						●																								
Morphine-3-β-d-Glucuronide						●																								
Mumps Virus IgG			●																											
MXD (W-MCC & W-MCR)			●																											
Mycoplasma pneumoniae IgG				●																										
Mycoplasma pneumoniae IgM				●																										
Myoglobin		●			●																									
N-Acetyl Procainamide (NAPA)												●							▶								●	●	●	
nDNA													●																	
Neisseria gonorrhoeae (NG or GC)				●																										
Netilmicin																											▲	▲		
Neuron Specific Enolase (NSE)*																														
NEUT (W-LCC & W-LCR)			●																											
NEUT X			●																											
NEUT Y			●																											
Neutrophils (NEUT)			●																											
Nickel																														
Nicotine																				▶										
Nitrite																														
Non-treponemal Antibodies (Syphilitic)				●																										
Norclomipramine														●						▶										
Nordiazepam						●														▶										
Nordoxepin																				▶										
Norepinephrine																														
Normetanephrine																														
Norpropoxyphene						●																								
Nortriptyline						●														▶								●	●	
NT-proBNP		●			●																									
Nucleated Red Blood Cells (NRBC)			●																											
Nucleated Red Blood Cells/100 WBC (NRBC/			●																											
Opiates						●																								
Osmolality					●					○	●	●																		
Osteocalcin					●																									
Oxazepam						●														▶										
Oxycodone						●																								
PAP					●					●	▲																	●	●	●
PAPP-A																														
pCO ₂	●																													
PCT			✱																											
PDW			✱																											
Pentachlorophenol																														
Pentobarbital						●																								
pH	●					◆																								
Phencyclidine (PCP)						●																								
Phenobarbital					●	●				●	●	●								▶								●	●	●
Phenol																														
Phenothiazines																				▶										
Phenylalanine																														

*International use only.
Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Control Analyte Index

65	40	63	63	46	41	80	40	87	87	42	66	58	86	62	62	62	67	77	77	61	61	39	38	37	87	42	86	86	78	78	85	See page number						
Immunology Plus (Lyphocheck)	Lipids	Maternal Serum 1st Trimester	Maternal Serum II	Meter Trax	Microalbumin	Opiate	Pediatric	qUantify Advance	qUantify Plus	Quantitative Urine	Rheumatoid Factor	Sedimentation Rate	Serum Volatiles	Specialty Immunoassay (Liquichek)	Specialty Immunoassay Plus (Liquichek)	Specialty Immunoassay (Lyphocheck)	Spinal Fluid	TDM (Liquichek)	TDM (Lyphocheck)	Tumor Marker (Liquichek)	Tumor Marker Plus (Lyphocheck)	Unassayed Chemistry (Liquichek)	Unassayed Chemistry (Lyphocheck)	Unassayed Multiqua (Liquid)	Urinanalysis (Liquichek)	Urine Chemistry (Liquichek)	Urine Metals	Whole Blood	Whole Blood Immunosuppressant (Liquid)	Whole Blood Immunosuppressant (Lypho.)	Whole Blood Metals							
																																	Morphine					
																																		Morphine (Free)				
																																		Morphine-3-β-d-Glucuronide				
																																		Mumps Virus IgG				
																																		MXD				
																																		Mycoplasma pneumoniae IgG				
																																		Mycoplasma pneumoniae IgM				
																																		Myoglobin				
																																			N-Acetyl Procainamide (NAPA)			
																																			nDNA			
																																			Neisseria gonorrhoeae (NG or GC)			
																																			Netilmicin			
																																				Neuron Specific Enolase (NSE)*		
																																				NEUT (W-LCC & W-LCR)		
																																				NEUT X		
																																				NEUT Y		
																																				Neutrophils (NEUT)		
																																				Nickel		
																																					Nicotine	
																																					Nitrite	
																																					Non-treponemal Antibodies (Syphilitic)	
																																					Norclomipramine	
																																					Nordiazepam	
																																						Nordoxepin
																																						Norepinephrine
																																						Normetanephrine
																																						Norpropoxyphene
																																						Nortriptyline
																																						NT-proBNP
																																						Nucleated Red Blood Cells (NRBC)
																																						Nucleated Red Blood Cells/100 WBC
																																						Opiates
																																						Osmolality
																																						Osteocalcin
																																						Oxazepam
																																						Oxycodone
																																						PAP
																																						PAPP-A
																																						pCO ₂
																																						PCT
																																						PDW
																																						Pentachlorophenol
																																						Pentobarbital
																																						pH
																																						Phencyclidine (PCP)
																																						Phenobarbital
																																						Phenol
																																						Phenothiazines
																																						Phenylalanine
																																						Phenylglyoxylic Acid

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ✳ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

*International use only.

Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

See page number

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ✱ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

	111	111	112	117	119	79	67	63	62	38	37	41	68	85	44	44	45	45	85	66	39	63	46	45	44	64	59	60	64	65
	See Blood Gas Chart	See Cardiac Assessment Chart	See Hematology Chart	See Infectious Disease/Serology Chart	See IntelliQ Chart	Urine Toxicology (See Product Pages)	Allergen sIgE	AMH	Anemia	Assayed Chemistry	Assayed Multiqua (Liquit)	Assayed Multiqua Premium	Autoimmune	Benzo/TCA-Set B	Coagulation (Lyphocheck)	D-dimer	Diabetes (Liquicheck)	Diabetes (Lyphocheck)	Drug Free Serum	Elevated CRP	Ethano/Ammonia (Liquicheck)	Fertility	Hemoglobin A1C Linearity Set	Hemoglobin A ₂	Hemostasis	Hypertension Markers	Immunoassay Plus (Liquicheck)	Immunoassay Plus (Lyphocheck)	Immunoassay Premium	Immunology (Liquicheck)
Phenylglyoxylic Acid																														
Phenytoin					●					●	●	●							▶								●	●	●	
Phenytoin (Free)																											▲	●	●	
Phospholipids					●					●	●	●																		
Phosphorus					●					●	●	●																		
Plasminogen																									●					
Platecrit (PCT)			●																											
Platelet Distribution Width (PDW)			✱																											
Platelet Volume Distribution Width (PDW)			✱																											
Platelets (PLT)			●																											
Platelets Impedance Method (PLTi)			●																											
Platelets, Optical Count (PLTo)			●																											
pO ₂	●																													
Porphyrins (Total)																														
Potassium	●				●					●	●	●																		
Prealbumin					○				○	○	○																			●
Pregnancy/hCG (Qualitative)																														
Pregnanetriol																														
Primidone																			▶								●	●	●	
Procainamide												●							▶								●	●	●	
Procalcitonin																														
Progesterone					●																	●					●	●	●	
ProGRP*					●																									
Prolactin					●																						●	●	●	
Properdin Factor B																														●
Propoxyphene						●													▶											
Propranolol																			▶											
Prostatic Acid Phosphatase (PAP)										●	■																●	▲	●	
Protein (Total)																											●	●	●	
Protein (Total Urine)					●																									
Protein C (Functional)																								●						
Protein Electrophoresis																								●						
Protein S (Functional)																								●						
Protein Serum (Total)		●			●					●	●	●																		●
Protein-to-Creatinine Ratio																														
Prothrombin Time (PT)															●									●						
PSA (Free)					●																						●	●	●	
PSA (Total)					●					□																	●	●	●	
Pseudocholinesterase																														
PTH (Intact)					●																							▲		
PTH-MM																											▲			
Pyridamine																				▶										
Quinidine												●							▶								●	▲	●	
Quinine																			▶											
RDW-SD			●																											
Red Blood Cells (RBC)			●																											
Red Blood Cells, Optical (RBCo)			●																											
Red Cell Distribution Width (RDW)			●																											
Red Cell Distribution Width-CV (RDW-CV)			●																											
Red Cell Distribution Width-SD (RDW-SD)			●																											

*International use only.
Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Control Analyte Index

65	40	63	63	46	41	80	40	87	87	42	66	58	86	62	62	62	67	77	77	61	61	39	38	37	87	42	86	86	78	78	85	See page number		
Immunology Plus (Lyphocheck)	Lipids	Maternal Serum 1st Trimester	Maternal Serum II	Meter Trax	Microalbumin	Opiate	Pediatric	qUAntify Advance	qUAntify Plus	Quantitative Urine	Rheumatoid Factor	Sedimentation Rate	Serum Volatiles	Specialty Immunoassay (Liquichek)	Specialty Immunoassay Plus (Liquichek)	Specialty Immunoassay (Lyphocheck)	Spinal Fluid	TDM (Liquichek)	TDM (Lyphocheck)	Tumor Marker (Liquichek)	Tumor Marker Plus (Lyphocheck)	Unassayed Chemistry (Liquichek)	Unassayed Chemistry (Lyphocheck)	Unassayed Multiqua (Liquid)	Urinalysis (Liquichek)	Urine Chemistry (Liquichek)	Urine Metals	Whole Blood	Whole Blood Immunosuppressant (Liquid)	Whole Blood Immunosuppressant (Lyphocheck)	Whole Blood Metals			
																																	Phenytoin	
																																		Phenytoin (Free)
																																		Phospholipids
																																		Phosphorus
																																		Plasminogen
																																		Plateletcrit (PCT)
																																		Platelet Distribution Width (PDW)
																																		Platelet Volume Distribution Width (PDW)
																																		Platelets (PLT)
																																		Platelets Impedance Method (PLTi)
																																		Platelets, Optical Count (PLTo)
																																		pO ₂
																																		Porphyrins (Total)
																																		Potassium
																																		Prealbumin
																																		Pregnancy/hCG (Qualitative)
																																		Pregnanetriol
																																		Primidone
																																		Procainamide
																																		Procalcitonin
																																		Progesterone
																																		ProGRP*
																																		Prolactin
																																		Properdin Factor B
																																		Propoxyphene
																																		Propranolol
																																		Prostatic Acid Phosphatase (PAP)
																																		Protein (Total)
																																		Protein (Total Urine)
																																		Protein C (Functional)
																																		Protein Electrophoresis
																																		Protein S (Functional)
																																		Protein Serum (Total)
																																		Protein-to-Creatinine Ratio
																																		Prothrombin Time (PT)
																																		PSA (Free)
																																		PSA (Total)
																																		Pseudocholesterase
																																		PTH (Intact)
																																		PTH-MM
																																		Pyrimidine
																																		Quinidine
																																		Quinine
																																		RDW-SD
																																		Red Blood Cells (RBC)
																																		Red Blood Cells, Optical (RBCo)
																																		Red Cell Distribution Width (RDW)
																																		Red Cell Distribution Width-CV (RDW-CV)
																																		Red Cell Distribution Width-SD (RDW-SD)
																																		Red Cell Folate

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ✳ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

*International use only.

Refer to myInserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

See page number

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ▲ Values not for diagnostic use
- ✱ No claim is made for expected value or stability

	111	111	112	117	119	79	67	63	62	38	37	41	68	85	44	44	45	45	85	66	39	63	46	45	44	64	59	60	64	65	
	See Blood Gas Chart	See Cardiac Assessment Chart	See Hematology Chart	See Infectious Disease/Serology Chart	See IntellIQ Chart	Urine Toxicology (See Product Pages)	Allergen sigE	AMH	Anemia	Assayed Chemistry	Assayed Multiqua (Liquit)	Assayed Multiqua Premium	Autoimmune	Benzo/TCA-Set B	Coagulation (Lyphocheck)	D-dimer	Diabetes (Liquicheck)	Diabetes (Lyphocheck)	Drug Free Serum	Elevated CRP	Ethano/Ammonia (Liquicheck)	Fertility	Hemoglobin A1C Linearity Set	Hemoglobin A ₂	Hemostasis	Hypertension Markers	Immunoassay Plus (Liquicheck)	Immunoassay Plus (Lyphocheck)	Immunoassay Premium	Immunology (Liquicheck)	
Red Cell Folate																															
Renin																											●				
Reticulocyte (Retic)			●																												
Retinol Binding Protein (RBP)																														●	
Rheumatoid Factor				●																										●	
Rubella Virus IgG			●																												
Rubella Virus IgM			●																												
Rubeola Virus (Measles) IgG			●																												
S-100*																															
Salicylate				●						●	●	●							▶									●	●	●	
SARS-CoV-2 IgG			●																												
SARS-CoV-2 Total IgG/IgM			●																												
SCC*				●																											
Secobarbital					●														▶												
Sedimentation Rate (ESR)																															
Selenium																															
Serotonin																															
SHBG (Sex Hormone Binding Globulin)				●																							○	▲	●		
Sirolimus																															
Smooth Muscle												●																			
Sodium	●			●						●	●	●																			
Specific Gravity					◆																										
Spun Hematocrit (Spun HCT)		●																													
Syphilis RPR			●																												
Syphilitic Reagin			●																												
T3 Birch (Betula)					●																										
T3 (Free)				●						○																		●	●	●	
T3 (Total)				●						●	●	●															●	●	●		
T3 Uptake/T-Uptake				●						●	●	●															●	●	●		
T4 (Free)				●						○	●	●															●	●	●		
T4 (Total)				●						●	●	●															●	●	●		
T-Uptake				●																											
Tacrolimus (FK 506)																															
TBG										○																	●	○	●		
Testosterone				●																		●					●	●	●		
Testosterone (Free)																											●	●			
Thallium																															
THC Metabolite					●																										
Theophylline				●						●	●	●							▶								●	●	●		
Thrombin Time (TT)															●										●						
Thyroglobulin (Tg)																											▲	○			
Thyroid Stimulating Immunoglobulin (TSI)																															
Tobramycin				●						●	●	●							▶								●	●	●		
Total Red Blood Cell Counts in Retic Channel			●																												
Toxoplasma gondii IgG			●																												
Toxoplasma gondii IgM			●																												
Tramadol																															
Transferrin				●					●	○	○	○			●															●	
Trazodone															●				▶												
Treponema pallidum (Syphilis) IgG			●																												

*International use only.
Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

65	40	63	63	46	41	80	40	87	87	42	66	58	86	62	62	62	67	77	77	61	61	39	38	37	87	42	86	86	78	78	85		
Immunology Plus (Lyphocheck)	Lipids	Maternal Serum 1st Trimester	Maternal Serum II	Meter Trax	Microalbumin	Opiate	Pediatric	qUAntify Advance	qUAntify Plus	Quantitative Urine	Rheumatoid Factor	Sedimentation Rate	Serum Volatiles	Specialty Immunoassay (Liquichek)	Specialty Immunoassay Plus (Liquichek)	Specialty Immunoassay (Lyphocheck)	Spinal Fluid	TDM (Liquichek)	TDM (Lyphocheck)	Tumor Marker (Liquichek)	Tumor Marker Plus (Lyphocheck)	Unassayed Chemistry (Liquichek)	Unassayed Chemistry (Lyphocheck)	Unassayed Multiqua (Liquid)	Urinanalysis (Liquichek)	Urine Chemistry (Liquichek)	Urine Metals	Whole Blood	Whole Blood Immunosuppressant (Liquid)	Whole Blood Immunosuppressant (Lypho.)	Whole Blood Metals		
																																Renin	
																																Reticulocyte (Retic)	
●																																Retinol Binding Protein (RBP)	
●											●																					Rheumatoid Factor	
																																Rubella Virus IgG	
																																Rubella Virus IgM	
																																Rubeola Virus (Measles) IgG	
																						●										S-100*	
																			●	●			●	●	●							Salicylate	
																																SARS-CoV-2 IgG	
																																SARS-CoV-2 Total IgG/IgM	
																					●											SCC*	
																																Secobarbital	
												●																				Sedimentation Rate (ESR)	
																											●					Selenium	
																												▲				Serotonin	
																	●															SHBG (Sex Hormone Binding Globulin)	
																												●	●			Sirolimus	
																																Smooth Muscle	
								●			●												●	●	●		●					Sodium	
								●	●	●	●															●	●					Specific Gravity	
																																Spun Hematocrit (Spun HCT)	
																																Syphilis RPR	
																																Syphilitic Reagin	
																																T3 Birch (Betula)	
																			●				▲	○								T3 (Free)	
																			●	●			●	●	●							T3 (Total)	
																			●	●			●	●	●							T3 Uptake/T-Uptake	
																			▲	●			▲	○	●							T4 (Free)	
																			●	●			●	●	●							T4 (Total)	
																																	T-Uptake
																													●	●			Tacrolimus (FK 506)
																																	TBG
																																	Testosterone
																																	Testosterone (Free)
																											●						Thallium
																																	THC Metabolite
								●												●	●		●	●	●							Theophylline	
																																	Thrombin Time (TT)
																						●	●										Thyroglobulin (Tg)
																●																	Thyroid Stimulating Immunoglobulin (TSI)
																						●	●	●									Tobramycin
																																	Total Red Blood Cell Counts in Retic Channel
																																	Toxoplasma gondii IgG
																																	Toxoplasma gondii IgM
									●																								Tramadol
●																																	Transferrin
																																	Trazodone
																																	Treponema pallidum (Syphilis) IgG
																																	Treponema pallidum (Syphilis) IgM

- See page number
- Analyte present
 - Analyte present in endogenous levels
 - Analyte present and no claims are made regarding performance or stability
 - Analyte present in endogenous levels and no claims are made regarding performance or stability
 - ▶ Tested and found to be negative
 - ◆ Specimen validity criteria
 - ✳ Values not for diagnostic use
 - ▲ No claim is made for expected value or stability

*International use only.

Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

See page number

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ✖ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

	111	111	112	117	119	79	67	63	62	38	37	41	68	85	44	44	45	45	85	66	39	63	46	45	44	64	59	60	64	65	
	See Blood Gas Chart	See Cardiac Assessment Chart	See Hematology Chart	See Infectious Disease/Serology Chart	See IntellIQ Chart	Urine Toxicology (See Product Pages)	Allergen sIgE	AMH	Anemia	Assayed Chemistry	Assayed Multiqua (Liquit)	Assayed Multiqua Premium	Autoimmune	Benzo/TCA-Set B	Coagulation (Lyphocheck)	D-dimer	Diabetes (Liquicheck)	Diabetes (Lyphocheck)	Drug Free Serum	Elevated CRP	Ethano/Ammonia (Liquicheck)	Fertility	Hemoglobin A1C Linearity Set	Hemoglobin A ₂	Hemostasis	Hypertension Markers	Immunoassay Plus (Liquicheck)	Immunoassay Plus (Lyphocheck)	Immunoassay Premium	Immunology (Liquicheck)	
Treponema pallidum (Syphilis) IgM				●																											
Trichloroacetic Acid																															
Tricyclic Antidepressants						●																									
Tricyclic Antidepressants (TCA) Screen																											●	▲	●		
Triglycerides					●					●	●	●																			
Troponin I		●			●																										
hsTroponin I		●			●																										
Troponin T		●			●																										
hsTroponin T		●			●																										
Trypanosoma cruzi (T. cruzi)				●																											
TSH					●				●	●	●	●																●	●	●	
Urea					●					●	●	●																			
Urea Nitrogen (BUN)					●					●	●	●																			
Uric Acid					●					●	●	●																			
Urobilinogen																															
Uroporphyrin																															
Valproic Acid					●					●	●	●								▶								●	●	●	
Valproic Acid (Free)																												▲	▲	●	
Vancomycin					●					●		●								▶								●	●	●	
Vancomycin Resistant Enterococci (VRE)				●																											
Varicella Zoster Virus (VZV) IgG				●																											
Vitamin B ₁₂					●				●	○	○	○																●	●	●	
VMA																															
W6 Mugwort (Artemesia vulgaris)						●																									
West Nile Virus IgG				●																											
West Nile Virus IgM				●																											
White Blood Cells (WBC)		●																													
White Blood Cells Counts Measured in Perox Chan-		●																													
Zinc					●					○	●	●																			

*International use only.
Refer to myInSerts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

65	40	63	63	46	41	80	40	87	87	42	66	58	86	62	62	62	67	77	77	61	61	39	38	37	87	42	86	86	78	78	85	See page number	
Immunology Plus (Lyphocheck)	Lipids	Maternal Serum 1st Trimester	Maternal Serum II	Meter Trax	Microalbumin	Opiate	Pediatric	qUAntify Advance	qUAntify Plus	Quantitative Urine	Rheumatoid Factor	Sedimentation Rate	Serum Volatiles	Specialty Immunoassay (Liquichek)	Specialty Immunoassay Plus (Liquichek)	Specialty Immunoassay (Lyphocheck)	Spinal Fluid	TDM (Liquichek)	TDM (Lyphocheck)	Tumor Marker (Liquichek)	Tumor Marker Plus (Lyphocheck)	Unassayed Chemistry (Liquichek)	Unassayed Chemistry (Lyphocheck)	Unassayed Multiqua (Liquid)	Urinalysis (Liquichek)	Urine Chemistry (Liquichek)	Urine Metals	Whole Blood	Whole Blood Immunosuppressant (Liquid)	Whole Blood Immunosuppressant (Lyoph.)	Whole Blood Metals		
																																	Trichloroacetic Acid
																																	Tricyclic Antidepressants
																																	Tricyclic Antidepressants (TCA) Screen
	●																																Triglycerides
																																	Troponin I
																																	Troponin I Ultra
																																	Troponin T
																																	Trypanosoma cruzi (T. cruzi)
																																	TSH
																																	Urea
																																	Urea Nitrogen (BUN)
																																	Uric Acid
																																	Urobilinogen
																																	Uroporphyrin
																																	Valproic Acid
																																	Valproic Acid (Free)
																																	Vancomycin
																																	Vancomycin Resistant Enterococci (VRE)
																																	Varicella Zoster Virus (VZV) IgG
																																	Vitamin B ₁₂
																																	VMA
																																	W6 Mugwort (Artemisia vulgaris)
																																	West Nile Virus IgG
																																	West Nile Virus IgM
																																	White Blood Cells (WBC)
																																	White Blood Cells Counts Measured in Perox
																																	Zinc

- Analyte present
- Analyte present in endogenous levels
- Analyte present and no claims are made regarding performance or stability
- Analyte present in endogenous levels and no claims are made regarding performance or stability
- ▶ Tested and found to be negative
- ◆ Specimen validity criteria
- ✳ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

*International use only.

Refer to myInserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Blood Gas Controls

	34	34	34
● Analyte present	Blood Gas	Blood Gas Plus E	Blood Gas Plus EGL
Chloride		●	●
Calcium (Ionized)		●	●
Glucose			●
Lactate (Lactic Acid)			●
Lithium			●
Magnesium			●
pCO ₂	●	●	●
pH	●	●	●
pO ₂	●	●	●
Potassium		●	●
Sodium		●	●

Cardiac Assessment Controls

	35	35	36	36	36	36
● Analyte present ■ Analyte present and no claims are made regarding performance or stability	Cardiac Advance	Cardiac Markers Plus LT	Cardiac Markers Plus	Cardiac Troponins	CK/LD Isoenzyme	Homocysteine
Albumin					●	
B-type Natriuretic Peptide (BNP)	●	●	●			
Creatine Kinase (CK)		●	●		●	
CK Electrophoresis					●	
CK-MB Isoenzyme	●	●	●		●	
CRP		●	●			
hsCRP	●	●	●			
Digitoxin	●	●	●			
Galectin-3	●					
Globulin					●	
Homocysteine	●	■	■			●
LD Electrophoresis					●	
LD-1 Isoenzyme					●	
LDH					●	
Myoglobin	●	●	●			
NT-proBNP	●	●	●			
Protein Serum (Total)					●	
Protein Electrophoresis					●	
Troponin I			●			
hsTroponin I	●	●		●		
Troponin T			●			
hsTroponin T	●	●		●		

Refer to myInsersts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Hematology Controls

See page number

(A) for Abbot Systems
(C) for Beckman Coulter Systems
(S) for Siemens Systems
(X & 16T) for Sysmex Systems
(16) for 3 -part differentials

- Analyte present
- ✘ Values not for diagnostic use
- ▲ No claim is made for expected value or stability

	53	53	53	54	54	55	55	56	56	57	57	58
	Hematology (A)	Reticulocyte (A)	Reticulocyte (A-1)	Hematology (C)	Reticulocyte (C)	Hematology (S)	Reticulocyte (S)	Hematology (X)	Reticulocyte (X)	Hematology-16T	Hematology-16	Sedimentation Rate
Basophils (BASO or BA)	●			●		●		●				
Cellular Hemoglobin Concentration Mean (CHCM)						●						
Cellular Hemoglobin Concentration Mean of Gated Red Blood Cells (CHCMg)							●					
Cellular Hemoglobin Concentration Mean of Reticulocytes (CHCMr)							●					
Eosinophils (EOS or EO)	●			●		●		●				
Erythrocyte Sedimentation Rate (ESR)												●
Granulocytes (GRAN)	●									●	●	
Hematocrit (HCT)	●			●		●		●		●	●	
Hemoglobin (HGB)	●			●		●		●		●	●	
Hemoglobin Concentration Distribution Width (HDW)						●						
Immature Reticulocyte Fraction (IRF)			●		●				●			
Large Unstained Cells (LUC)						●						
LYMPH (W-SCC & W-SCR)										●		
Lymphocytes (LYMPH or LY)	●			●		●		●		●	●	
Mean Corpuscular Hemoglobin (MCH)	●			●		●		●		●	●	
Mean Corpuscular Hemoglobin Concentration (MCHC)	●			●		●		●		●	●	
Mean Corpuscular Volume (MCV)	●			●		●		●		●	●	
Mean Corpuscular Volume of Gated Red Blood Cells (MCVg)							●					
Mean Corpuscular Volume of Reticulocytes (MCVr)							●					
Mean Hemoglobin Content of All Gated Red Cells (CHg)							●					
Mean Hemoglobin Content of Reticulocytes (CHr)							●					
Mean Platelet Volume (MPV)	●			●		●		●		●	●	
Mean Reticulocyte Volume (MRV)					●							
MID	●									●	●	
Monocytes (MONO or MO)	●			●		●		●			●	
MXD (W-MCC & W-MCR)										●		
NEUT (W-LCC & W-LCR)										●		
Neutrophils (NEUT or NE)	●			●		●		●				
NEUTx						●						
NEUTy						●						
NRBC/100 WBC (NRBC percent of WBC Count)	●											
Nucleated Optical Count (NOC)	●											
Nucleated Red Blood Cells (NRBC)	●							▲				
Platelet Distribution Width (PDW)	✘			✘								
Plateletcrit (PCT)	✘			✘								
Platelets (PLT)	●			●		●		●		●	●	
Platelets Impedance Method (PLTi)	●											
Platelets, Optical Count (PLT-o)								▲				
Red Blood Cell Distribution Width (RDW)	●			●		●				●	●	
Red Blood Cell Distribution Width-CV (RDW-CV)								●		●		
Red Blood Cell Distribution Width-SD (RDW-SD)				●				●		●		
Red Blood Cells (RBC)	●	●	●	●	●	●	●	●	●	●	●	
Red Blood Cells, Optical (RBCo)	●											
Reticulocyte (Retic)		●	●		●		●		●			
Total Red Blood Cell Counts in Retic Channel (Retic RBC)							●					
White Blood Cells (WBC)	●			●		●		●		●	●	
White Blood Cells Counts Measured in Perox Channel (WBCP Perox)						●						
White Impedance Count (WIC)	●											
White Optical Count	●											

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

See page number

Infectious Disease: Molecular Quality Controls and Standards

- Analyte present
- Negative for analyte

	16	29	24	32	32	32	16	32	28	17	19	18	33	27	27	22	24	29	22	24	18	23
	Positive Controls																					
	Adenovirus	Amplichek II	AMR Bundle	Anaplasma	Babesia	Bartonella	BKV	Borrelia	C. difficile Verification Bundle	CMV	CT/NG	EBV	Ehrlichia	Enterovirus	GBS	HBV	HBV AMR Panel	HBV, HCV & HIV-1 Screen Controls	HCV	HCV AMR Panel	HHV-6	HIV
Adenovirus																						
Anaplasma phagocytophilum				●																		
Babesia microti					●																	
Bartonella quintana						●																
BKV							●															
Inactivated Bordetella parapertussis																						
Inactivated Bordetella pertussis																						
Borrelia burgdorferi								●														
Inactivated Chlamydia trachomatis (CT)											●											
Inactivated Chlamydomphila pneumoniae																						
Candida auris (C. auris)																						
Clostridium difficile (C. diff.)		●																				
CMV																						
Inactivated Coronavirus 229E																						
Coronavirus HKU1 (IVT)																						
Coronavirus NL63 (IVT)																						
Inactivated Coronavirus OC43																						
EBV																						
Ehrlichia chaffeensis													●									
Enterovirus (EV) Coxsackievirus A9																						
HBsAg																						
HBV																						
HCV Antibodies																						
HCV																						
HHV-6																						
HIV-1																						
HIV-1/2 Antibodies																						
HSV-1/2																						
Human Cell Line																						
Human metapneumovirus (IVT)																						
Human Papillomavirus (HPV) Genotype 16																						
Human Papillomavirus (HPV) Genotype 18																						
Human Papillomavirus (HPV) Genotype 68																						
Inactivated Adenovirus	●																					
Inactivated BKV							●															
Inactivated C. difficile								●														
Inactivated CMV									●													
Inactivated Coxsackievirus A9										●				●								
Inactivated EBV												●										
Inactivated HBV			●													●	●	●				
Inactivated HCV																		●				
Inactivated HHV-6A																					●	
Inactivated HHV-6B																					●	
Inactivated HIV-1			●															●				●
Inactivated HSV-1																						
Inactivated HSV-2																						
Inactivated JCV																						
Inactivated Methicillin-resistant Staphylococcus aureus (MRSA)																						
Inactivated Methicillin-susceptible Staphylococcus aureus (MSSA)																						
Inactivated Parvovirus B19																						
Inactivated Varicella Zoster Virus (VZV)																						
Inactivated Zika Virus																						
Influenza A H1N1																						
Influenza A H1N1-2009 (IVT)																						

Refer to myInsersts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

24	20	21	21	19	26	17	31	30	31	31	31	30	20	27	33	29	26	26	25	25	25	26	See page number	
Positive Controls																Negative						<h2>Infectious Disease: Molecular Quality Controls and Standards</h2> <ul style="list-style-type: none"> ● Analyte present ■ Negative for analyte 		
HIV-1 AMR Panel	HPV	HSV-1	HSV-2	JCV	MRSA Verification Bundle	Parvovirus B19	RP Positive & Negative	SARS-CoV-2	SARS-CoV-2 S Gene Alpha Variant	SARS-CoV-2 S Gene Beta Variant	SARS-CoV-2 S Gene Delta-Kappa Variant	SARS-CoV-2 S Gene Epsilon Variant	SARS-CoV-2 S Gene Gamma Variant	SARS-CoV-2, Flu, RSV Run Control	TV/MG	VZV	ZKV	Amplichek II	CSF Negative	Exact Diagnostics Negative	HAI Negative		Synthetic Negative	STI Negative
																				■			■	Adenovirus
																				■			■	Anaplasma phagocytophilum
																				■			■	Babesia microti
																				■			■	Bartonella quintana
																				■				BKV
							●																	Inactivated Bordetella parapertussis
							●																	Inactivated Bordetella pertussis
																			■			■		Borrelia burgdorferi
																							■	Inactivated Chlamydia trachomatis (CT)
							●																■	Inactivated Chlamydomphila pneumoniae
																								Candida auris (C. auris)
																		■			■			Clostridium difficile (C. diff.)
																				■				CMV
							●																	Inactivated Coronavirus 229E
							●																	Coronavirus HKU1 (IVT)
							●																	Coronavirus NL63 (IVT)
							●																	Inactivated Coronavirus OC43
																				■				EBV
																			■			■		Ehrlichia chaffeensis
																			■			■		Enterovirus (EV) Coxsackievirus A9
																				■				HBsAg
																				■				HBV
																				■				HCV Antibodies
																				■				HCV
																				■				HHV-6
																				■				HIV-1
																				■		■		HIV-1/2 Antibodies
																			■	■				HSV-1/2
																								Human Cell Line
							●																	Human metapneumovirus (IVT)
	●																							Human Papillomavirus (HPV) Genotype 16
	●																							Human Papillomavirus (HPV) Genotype 18
	●																							Human Papillomavirus (HPV) Genotype 68
							●																	Inactivated Adenovirus
																								Inactivated BKV
																								Inactivated C. difficile
																								Inactivated CMV
																								Inactivated Coxsackievirus A9
																								Inactivated EBV
																								Inactivated HBV
																								Inactivated HCV
																								Inactivated HHV-6A
																								Inactivated HHV-6B
	●																							Inactivated HIV-1
		●																						Inactivated HSV-1
			●																					Inactivated HSV-2
				●																				Inactivated JCV
					●																			Inactivated Methicillin-resistant Staphylococcus aureus (MRSA)
						●																		Inactivated Methicillin-susceptible Staphylococcus aureus (MSSA)
							●																	Inactivated Parvovirus B19
														●										Inactivated Varicella Zoster Virus (VZV)
															●									Inactivated Zika Virus
							●																	Influenza A H1N1
							●																	Influenza A H1N1-2009 (IVT)

Refer to myInserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

See page number

Infectious Disease: Molecular Quality Controls and Standards

- Analyte present
- Negative for analyte

	16	29	24	32	32	32	16	32	28	17	19	18	33	27	27	22	24	29	22	24	18	23	
	Positive Controls																						
	Adenovirus	Amplichek II	AMR Bundle	Anaplasma	Babesia	Bartonella	BKV	Borrelia	C. difficile Verification Bundle	CMV	CT/NG	EBV	Ehrlichia	Enterovirus	GBS	HBV	HBV AMR Panel	HBV, HCV & HIV-1 Screen Controls	HCV	HCV AMR Panel	HHV-6	HIV	
Influenza A H3N2 (IVT)																							
Influenza A-H1N1																							
Influenza B																							
JC Virus																							
Methicillin-resistant <i>Staphylococcus aureus</i> (MRSA)		●																					
Methicillin-sensitive <i>Staphylococcus aureus</i> (MSSA)		●																					
Middle Eastern Respiratory Syndrome (IVT)																							
<i>Mycoplasma genitalium</i>																							
<i>Mycoplasma pneumoniae</i>																							
<i>Neisseria gonorrhoeae</i> (NG)											●												
Parainfluenza 1																							
Parainfluenza 2																							
Parainfluenza 3																							
Parainfluenza 4																							
Recombinant HCV			●																	●	●		
Respiratory Syncytial Virus (A)																							
Respiratory Syncytial Virus (B)																							
Rhinovirus																							
SARS-CoV-2 (Whole Virus)																							
SARS-CoV-2 Genes (E, N, ORF1ab, RdRP, S)																							
SARS-CoV-2, Synthetic RNA Transcript: Alpha Variant (B.1.1.7)																							
SARS-CoV-2, Synthetic RNA Transcript: Beta Variant (B.1.351)																							
SARS-CoV-2, Synthetic RNA Transcript: Delta-Kappa Variants (B.1.617.1, B.1.617.2 and AY sublineages, B.1.617.3)																							
SARS-CoV-2, Synthetic RNA Transcript: Epsilon Variant (B.1.427, B.1.427.1)																							
SARS-CoV-2, Synthetic RNA Transcript: Gamma Variant (P.1)																							
<i>Trichomonas vaginalis</i>																							
Vancomycin-resistant <i>Enterococci</i> (VRE or vanR)		●																					
VZV																							
Zika Virus (ZKV)																							

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Positive Controls														Negative											
24	20	21	21	19	28	17	31	30	31	31	31	31	31	30	20	27	33	29	26	26	25	25	25	26	
HIV-1 AMR Panel	HPV	HSV-1	HSV-2	JCV	MRSA Verification Bundle	Parvovirus B19	RP Positive & Negative	SARS-CoV-2	SARS-CoV-2 S Gene Alpha Variant	SARS-CoV-2 S Gene Beta Variant	SARS-CoV-2 S Gene Delta-Kappa Variant	SARS-CoV-2 S Gene Epsilon Variant	SARS-CoV-2 S Gene Gamma Variant	SARS-CoV-2, Flu, RSV Run Control	TV/MG	VZV	ZKV	Amplichek II	CSF Negative	Exact Diagnostics Negative	HAI Negative	Synthetic Negative	STI Negative	Urine Negative	
							●																		Influenza A H3N2 (IVT)
							●								●										Influenza A-H1N1
							●								●										Influenza B
																					■				JC Virus (JCV)
																		■				■			Methicillin-resistant <i>Staphylococcus aureus</i> (MRSA)
																		■				■			Methicillin-sensitive <i>Staphylococcus aureus</i> (MSSA)
							●																		Middle Eastern Respiratory Syndrome (IVT)
							●								●										<i>Mycoplasma genitalium</i> (MG)
							●																		<i>Mycoplasma pneumoniae</i>
							●																■	■	<i>Neisseria gonorrhoeae</i> (NG)
							●																		Parainfluenza 1
							●																		Parainfluenza 2
							●																		Parainfluenza 3
							●																		Parainfluenza 4
							●																		Recombinant HCV
							●							●											Respiratory Syncytial Virus (A)
							●																		Respiratory Syncytial Virus (B)
							●																		Rhinovirus
							●							●											SARS-CoV-2 (Whole Virus)
								●																	SARS-CoV-2 Genes (E, N, ORF1ab, RdRP, S)
									●																SARS-CoV-2, Synthetic RNA Transcript: Alpha Variant (B.1.1.7)
										●															SARS-CoV-2, Synthetic RNA Transcript: Beta Variant (B.1.351)
											●														SARS-CoV-2, Synthetic RNA Transcript: Delta-Kappa Variants (B.1.617.1, B.1.617.2 and AY sublineages, B.1.617.3)
												●													SARS-CoV-2, Synthetic RNA Transcript: Epsilon Variant (B.1.427, B.1.617.1, B.1.617.2 and AY sublineages, B.1.617.3)
													●												SARS-CoV-2, Synthetic RNA Transcript: Gamma Variant (P.1)
														●											<i>Trichomonas vaginalis</i> (TV)
																		■							Vancomycin-resistant Enterococci (VRE or vanR)
																			■	■			■		VZV
																					■			■	Zika Virus (ZKV)

See page number

Infectious Disease: Molecular Quality Controls and Standards

- Analyte present
- Negative for analyte

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

See page number

Infectious Disease: Serology Controls

- Analyte present
- Negative for analyte

	69	69	69	70	73	76	75	75	71	71	70	71	71	69	69	70	70
	Positive Controls																
	Assayed VIROTROL I-C	Assayed VIROTROL I-E	Assayed VIROTROL I-F	Assayed VIROTROL II-B	Liq. ToRCH Plus	PyJoriTrol	VIROTROL Chagas	VIROTROL EBV	VIROTROL HAV-IgM	VIROTROL HBeAg	VIROTROL HIV-1 Ag	VIROTROL HIV-1 gO	VIROTROL HIV-2	VIROTROL I	VIROTROL II	VIROTROL III	VIROTROL IV
Anti-Cytomegalovirus (Anti-CMV)	●	●	●											●			
Anti-HAV				●											●	●	
Anti-HBc	●	●	●											●			
Anti-HBe																	●
Anti-HBs				●											●		
Anti-HCV	●	●	●											●			
Anti-HIV-1	●	●	●											●			
Anti HIV-1/2																	
Anti-HIV-1 gO												●					
Anti-HIV-2													●				
Anti-HTLV-I	●	●	●											●			
Anti-HTLV-I/II																	
Anti-HTLV-II																	
Anti-Treponema pallidum (Syphilis)																	
Anti-Trypanosoma cruzi (T. cruzi)							●										
Borrelia burgdorferi IgG																	
Cytomegalovirus (CMV) IgG					●												
Cytomegalovirus (CMV) IgM																	
EBNA IgG								●									
EBV EA-D IgG																	
Epstein-Barr Virus (VCA) IgG					●			●									
Epstein-Barr Virus (VCA) IgM								●									
HAV-IgM									●							●	
HBc-IgM																●	
HBeAg										●							
HBsAg	●	●	●											●			
Helicobacter pylori IgG					●	●											
Herpes Simplex Virus Type 1 (HSV-1) IgG					●												
Herpes Simplex Virus Type 1 (HSV-1) IgM																	
Herpes Simplex Virus Type 2 (HSV-2) IgG					●												
Herpes Simplex Virus Type 2 (HSV-2) IgM																	
HIV-1 Ag											●						
Measles Virus (Rubeola) IgG																	
Mumps Virus IgG																	
Non-treponemal Antibodies (Syphilitic Reagin)																	
Rubella Virus IgG					●												
Rubella Virus IgM																	
SARS-CoV-2 Total IgG/IgM and SARS-CoV-2 IgG																	
Syphilis RPR					●												
Toxoplasma gondii IgG					●												
Toxoplasma gondii IgM																	
Treponema pallidum (Syphilis) IgG					●												
Treponema pallidum (Syphilis) IgM																	
Varicella Zoster Virus (VZV) IgG					●												

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

76	73	71	75	74	75	72	72	73	72	76	73	74	73	See page number
Positive Controls								Negative Controls					Infectious Disease: Serology Controls	
VIROCONTROL Lyme	VIROCONTROL MuMmZ	VIROCONTROL PLUS-R	VIROCONTROL RPR Panel	VIROCONTROL SARS-CoV-2	VIROCONTROL Syphilis LR-A	VIROCONTROL ToRCH	VIROCONTROL ToRCH-M	Liq. ToRCH Plus	VIROCLEAR	VIROCLEAR EBV	VIROCLEAR MuMmZ	VIROCLEAR SARS-CoV-2		VIROCLEAR ToRCH
		●							■					Anti-Cytomegalovirus (Anti-CMV)
									■					Anti-HAV
		●							■					Anti-HBc
									■					Anti-HBe
		●							■					Anti-HBs
									■					Anti-HCV
		●							■					Anti-HIV-1
									■					Anti HIV-1/2
									■					Anti-HIV-1 gO
									■					Anti-HIV-2
		●							■					Anti-HTLV-I
									■					Anti-HTLV-I/II
									■					Anti-HTLV-II
									■					Anti-Treponema pallidum (Syphilis)
									■					Anti-Trypanosoma cruzi (T. cruzi)
●														Borrelia burgdorferi IgG
						●	●	■					■	Cytomegalovirus (CMV) IgG
							●						■	Cytomegalovirus (CMV) IgM
										■				EBNA IgG
										■				EBV EA-D IgG
								■		■				Epstein-Barr Virus (VCA) IgG
										■				Epstein-Barr Virus (VCA) IgM
										■				HAV-IgM
										■				HBc-IgM
										■				HBeAg
		●								■				HBsAg
									■					<i>Helicobacter pylori</i> IgG
						●	●	■					■	Herpes Simplex Virus Type 1 (HSV-1) IgG
													■	Herpes Simplex Virus Type 1 (HSV-1) IgM
						●	●	■					■	Herpes Simplex Virus Type 2 (HSV-2) IgG
													■	Herpes Simplex Virus Type 2 (HSV-2) IgM
										■				HIV-1 Ag
	●												■	Measles Virus (Rubeola) IgG
	●												■	Mumps Virus IgG
										■				Non-treponemal Antibodies (Syphilitic Reagin)
						●	●	■					■	Rubella Virus IgG
							●						■	Rubella Virus IgM
				●									■	SARS-CoV-2 Total IgG/IgM and SARS-CoV-2 IgG
								■						Syphilis RPR
						●	●	■					■	Toxoplasma gondii IgG
							●						■	Toxoplasma gondii IgM
		●	●		●			■	■					Treponema pallidum (Syphilis) IgG
		●	●		●				■					Treponema pallidum (Syphilis) IgM
●								■			■			Varicella Zoster Virus (VZV) IgG

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

InteliQ™

See page number

- Analyte available
- Analyte present in endogenous levels

Test Menu

Analytes

		11	11	12	12	12	13	13	13	14	14	14	14
		InteliQ Assayed Multiquial	InteliQ Cardiac Advance	InteliQ Cardiac Markers Plus LT	InteliQ Diabetes	InteliQ Ethanol/Ammonia	InteliQ Homocysteine	InteliQ Immunoassay Plus	InteliQ Immunology	InteliQ Specialty Immunoassay	InteliQ Specialty Immunoassay Plus	InteliQ Tumor Marker	InteliQ Urine Chemistry
Anemia	Ceruloplasmin	○							●				
	Erythropoietin (EPO)									●	●		
	Ferritin	○						●	●			●	
	Folate							●					
	Active Vitamin B ₁₂	○						●			●		
Bone Metabolism	25-Hydroxy Vitamin D									●	●		
	PTH (Intact)									●	●		
	Osteocalcin									●	●		
Cancer/Oncology	α-Fetoprotein (AFP)							●				●	
	CA 15-3											●	
	CA 19-9											●	
	CA 27.29											●	
	CA 72-4*											●	
	CA 125											●	
	Carcinoembryonic Antigen (CEA)							●				●	
	Cyfra 21-1*											●	
	HE4											●	
	HER-2/neu											●	
	PAP											●	
	ProGRP*											●	
	PSA (Free)								●			●	
	PSA (Total)								●			●	
SCC*											●		
Thyroglobulin											●		
Cardiac	BNP		●	●									
	Creatine Kinase (CK)			●									
	CK-MB Isoenzyme		●	●				●					
	hsCRP		●	●									
	Galectin-3		●										
	Homocysteine		●										
	Myoglobin		●	●									
	NT-proBNP		●	●									
	hsTroponin I		●	●									
hsTroponin T		●	●										
Diabetes	C-Peptide									●	●		
	Fructosamine										●		
	Hemoglobin A1C				●								
	Insulin							●					
Fertility/Pregnancy	Androstenedione							●					
	DHEA-S							●					
	Estradiol							●					
	Estriol (Free)							●					
	FSH							●					
	hCG (Total)							●				●	
	LH							●					
	Pregnancy/hCG (Qualitative)												●
	Progesterone							●					
	Prolactin							●				●	
	SHBG (Sex Hormone Binding Globulin)							○					
Testosterone							●						

* International use only.



See page number

- Analyte available
- Analyte present in endogenous levels

Test Menu	Analytes	11	11	12	12	12	13	13	13	14	14	14	14
		InteliQ Assayed Multiquel	InteliQ Cardiac Advance	InteliQ Cardiac Markers Plus LT	InteliQ Diabetes	InteliQ Ethanol/Ammonia	InteliQ Homocysteine	InteliQ Immunoassay Plus	InteliQ Immunology	InteliQ Specialty Immunoassay	InteliQ Specialty Immunoassay Plus	InteliQ Tumor Marker	InteliQ Urine Chemistry
Growth	Insulin Growth Factor I (IGF-I)/Somatomedin C									●	●	●	
	hGH							●					
Inflammation	Antistreptolysin O (ASO)								●				
	CRP			●					●				
	Immunoglobulin E (IgE)							●	●				
	Interleukin-6 (IL-6)										●		
	Procalcitonin (PCT)										●		
Metabolic	Cortisol	●						●					●
	Homocysteine		●				●						
Therapeutic Drug Monitoring	Amikacin	●						●					
	Caffeine							●					
	Carbamazepine	●						●					
	Digitoxin		●	●									
	Digoxin	●						●					
	Gentamicin	●						●					
	Lidocaine							●					
	Lithium	●						●					
	NAPA							●					
	Phenobarbital	●						●					
	Phenytoin	●						●					
	Primidone							●					
	Procainamide							●					
	Quinidine							●					
	Theophylline	●						●					
	Tobramycin	●						●					
	Valproic Acid	●						●					
Vancomycin							●						
Thyroid	Anti-Thyroglobulin (Anti-Tg)									●	●		
	Anti-Thyropoxidase (Anti-TPO)									●	●		
	Anti-TSH Receptor (TRAb)										●		
	T3 (Free)							●					
	T3 (Total)	●						●					
	T3 Uptake/T Uptake	●						●					
	T4 (Free)	●						●					
	T4 (Total)	●						●					
	Thyroglobulin											●	
	Thyroid Stimulating Hormone (TSH)	●						●					
	Thyroid Stimulating Immunoglobulin (TSI)										●		
Thyroxine Binding Globulin (TBG)							●						
Toxicology: DOA	Acetaminophen	●						●					
	Ethanol	●				●							
	Salicylate	●						●					
	Tricyclic Antidepressants (TCA)							●					
	Zinc	●											

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.



See page number

- Analyte available
- Analyte present in endogenous levels

Test Menu	Analytes	11	11	12	12	12	13	13	13	14	14	14	14
		InteliQ Assayed Multiquel	InteliQ Cardiac Advance	InteliQ Cardiac Markers Plus LT	InteliQ Diabetes	InteliQ Ethanol/Ammonia	InteliQ Homocysteine	InteliQ Immunoassay Plus	InteliQ Immunology	InteliQ Specialty Immunoassay	InteliQ Specialty Immunoassay Plus	InteliQ Tumor Marker	InteliQ Urine Chemistry
General Chemistry	α-1-Antitrypsin	○											
	α-1-Globulin	○											
	α-2-Globulin	○											
	β-Globulin	○											
	γ-Globulin	○											
	γ-Glutamyltransferase (GGT)	●											
	Acid Phosphatase (Total)	●											
	Albumin	●							●				
	Albumin-to-Globulin Ratio	○											
	Alkaline Phosphatase	●											
	ALT/SGPT	●											
	Ammonia						●						
	Amylase	●											●
	Amylase (Pancreatic)	●											
	AST/SGOT	●											
	Bilirubin (Direct)	●											
	Bilirubin (Neonatal)	●											
	Bilirubin (Total)	●											
	Calcium	●											●
	Calcium (Ionized)	●											
	Carbon Dioxide (CO ₂)	●											
	Chloride	●											●
	Cholesterol (HDL)	●											
	Cholesterol (LDL)	●											
	Cholesterol (Total)	●											
	Cholinesterase	○											
	Copper	○											
	Creatine Kinase (CK)	●		●									
	Creatinine	●											●
	Glucose	●											●
	Iron	●							○				
	Iron (TIBC)	○											
	Iron (UIBC)	○											
	Lactate (Lactic Acid)	●											
	Lactate Dehydrogenase (LDH)	●											
	LAP Arylamidase	○											
	Lipase	●											
	Magnesium	●											●
	Microalbumin												●
	Osmolality	●											●
Phospholipids	●											●	
Phosphorus	●											●	
Potassium	●											●	
Protein (Total)	●											●	
Sodium	●											●	
Triglycerides	●											●	
Urea	●											●	
Urea Nitrogen (BUN)	●											●	
Uric Acid	●											●	

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.



See page number

- Analyte available
- Analyte present in endogenous levels

Test Menu	Analytes	11	11	12	12	12	13	13	13	14	14	14	14
		InteliQ Assayed Multiquel	InteliQ Cardiac Advance	InteliQ Cardiac Markers Plus LT	InteliQ Diabetes	InteliQ Ethano//Ammonia	InteliQ Homocysteine	InteliQ Immunoassay Plus	InteliQ Immunology	InteliQ Specialty Immunoassay	InteliQ Specialty Immunoassay Plus	InteliQ Tumor Marker	InteliQ Urine Chemistry
Specific Proteins	α-1-Acid Glycoprotein (AAG)								●				
	α-1-Antitrypsin (AAT)	○							●				
	α-2-Macroglobulin								●				
	ADNase B								●				
	Anti-Cyclic Citrullinated Peptide (Anti-CCP)								●				
	Antithrombin III (AT III)								●				
	Apolipoprotein A-1	○							●				
	Apolipoprotein B	○							●				
	β-2-Microglobulin								●			●	
	C1 Inhibitor								●				
	Complement C3	○							●				
	Complement C4	○							●				
	Cystatin C								○				
	Haptoglobin	○							●				
	IgG Subclass 1								●				
	IgG Subclass 2								●				
	IgG Subclass 3								●				
	IgG Subclass 4								●				
	Immunoglobulin A (IgA)	○							○	●			
	Immunoglobulin G (IgG)	○							○	●			
	Immunoglobulin M (IgM)	○							○	●			
	Kappa Light Chain								●				
	Lambda Light Chain								●				
	Lipoprotein (a)								●				
	pH												●
	Prealbumin	○							●				
	Propertin Factor B								●				
	Protein Serum (Total)	●							●				
Retinol Binding Protein (RBP)								●					
Rheumatoid Factor								●					
Specific Gravity												●	
Transferrin	○							●					

Refer to myinserts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

See page number

QC Products in Compact Vials

- Analyte available
- Analyte present in endogenous levels

	37	35	35	45	39	59	65	40	67	61	42
	Assayed Multiquant (Liquid)	Cardiac Advance	Cardiac Markers Plus LT (Liquichek)	Diabetes (Liquichek)	Ethanol/Ammonia (Liquichek)	Immunosay Plus (Liquichek)	Immunology (Liquichek)	Pediatric (Liquichek)	Spinal Fluid (Liquichek)	Tumor Marker (Liquichek)	Urine Chemistry (Liquichek)
α-1-Acid Glycoprotein							●				
α-1-Antitrypsin	○						●				
α-2-Macroglobulin							●				
β-2-Microglobulin							●			●	
Acetaminophen	●					●					
Albumin	●						●		●		
Alkaline Phosphatase (ALP)	●										
Alphafetoprotein (AFP)						●				●	
ALT/SGPT	●										
Amikacin						●					
Ammonia					●						
Amylase	●										●
Antistreptolysin O (ASO)							●				
Apolipoprotein A-1	○						●				
Apolipoprotein B	○						●				
AST/SGOT	●										
Bilirubin (Direct)	●							●			
Bilirubin (Total)	●							●			
BNP		●									
C1 Inhibitor							●				
CA 15-3										●	
CA 19-9										●	
CA 125										●	
Caffeine						●		●			
Calcium	●							●			●
Carbamazepine	●					●					
CEA						●				●	
Ceruloplasmin	○						●				
Chloride	●							●	●		●
Cholesterol (HDL)	●										
Cholesterol (LDL)	●										
Cholesterol (Total)	●										
Cholinesterase	○										
CK (Total)			●								
CK-MB Isoenzyme		●	●			●					
CO ₂	●										
Complement C3	○						●				
Complement C4	○						●				
Creatine Kinase (CK)	●										
Creatinine	●										●
CRP		●	●				●				
Cystatin C							○				
Digitoxin		●	●								
Digoxin	●					●					
Estradiol						●					
Ethanol	●				●						
Ferritin	○					●	●			●	
Folate						●					
FSH						●					
Galectin-3		●									
Gentamicin	●					●					
GGT	●										
Glucose	●							●	●		●

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

See page number

QC Products in Compact Vials

- Analyte available
- Analyte present in endogenous levels

	37	35	35	45	39	59	65	40	67	61	42
	Assayed Multiquial (Liquid)	Cardiac Advance	Cardiac Markers Plus LT (Liquichek)	Diabetes (Liquichek)	Ethanol/Ammonia (Liquichek)	Immunoassay Plus (Liquichek)	Immunology (Liquichek)	Pediatric (Liquichek)	Spinal Fluid (Liquichek)	Tumor Marker (Liquichek)	Urine Chemistry (Liquichek)
Haptoglobin	○						●				
hCG										●	
hCG-β Subunit						●					
Hemoglobin A1C				●							
Homocysteine		●									
Immunoglobulin A (IgA)	○					○	●				
Immunoglobulin E (IgE)						●	●				
Immunoglobulin G (IgG)	○					○	●		●		
Immunoglobulin G Subclasses 1–4							●				
Immunoglobulin M (IgM)	○					○	●				
Iron	●					○					
Iron (TIBC)	○										
Kappa Light Chain							●				
Lactate (Lactic Acid)	●								●		
Lactate Dehydrogenase (LDH)	●								●		
Lambda Light Chain							●				
Lidocaine						●					
Lipase	●										
Lithium	●					●					
Luteinizing Hormone (LH)						●					
Magnesium	●							●			●
Microalbumin											●
Myoglobin		●	●								
N-Acetylprocainamide (NAPA)						●					
NT-proBNP		●	●								
Phenobarbital	●					●					
Phenytoin	●					●					
Phosphorus	●										●
Potassium	●							●			●
Prealbumin	○						●				
Procainamide						●					
Prolactin						●				●	
Protein (Total)	●						●		●		●
PSA						●				●	
PSA (Free)						●				●	
Retinol Binding Protein (RBP)							●				
Rheumatoid Factor							●				
Salicylate	●					●					
Sodium	●							●	●		●
T3 (Free)						●					
T3 (Total)	●					●					
T3 Uptake/T Uptake	●					●					
T4 (Free)	●					●					
T4 (Total)	●					●					
Theophylline	●					●		●			
Tobramycin	●					●					
Transferrin	○						●				
Tricyclic Antidepressants (TCA)						●					
Triglycerides	●										
Troponin I		●	●								
TSH	●					●					
Urea Nitrogen (BUN)	●										●
Uric Acid	●										●
Valproic Acid	●					●					
Vancomycin						●					
Vitamin B ₁₂	○					●					

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

EQAS™ Programs

See page number

- Analyte present
- ▶ Pilot analyte

	47	47	47	48	47	48	48	49	49	49	49	50	50	50	50	51	51	51	
	Blood Gas	Blood Typing	Cardiac Markers	Clin. Chemistry (Monthly)	Coagulation	Ethanol/Ammonia	Hematology	Hemoglobin	HIV/Hepatitis	Immunoassay (Monthly)	Lipids	Serum Proteins	Specialty Immunoassay	Syphilis/Chagas	TDM	ToRCH/EBV/MuM/Z	Urinalysis	Urine Chemistry	
5-Hydroxyindoleacetic Acid (HIAA)																			●
11-Deoxycortisol										●									
17-α-OH-Progesterone										●									
25-Hydroxy Vitamin D										●									
α-1 Acid Glycoprotein												●							
α-1 Antitrypsin												●							
α-2 Macroglobulin												●							
β-2 Microglobulin										●		●							
ABO Group		●																	
Acetaminophen															●				
Acid Phosphatase (Total)				●															
ACTH										●									
Active Vitamin B ₁₂													●						
Alanine Aminotransferase (ALT/SGPT)				●															
Albumin				●								●						●	
Albumin-to-Creatinine Ratio																		●	
Aldosterone										●									
Alkaline Phosphatase (ALP)				●															
Alphafetoprotein (AFP)										●									
Amikacin															●				
Amitriptyline															●				
Ammonia						●													
Amylase (Pancreatic)				●															
Amylase (Total)				●															
Androstenedione										●									
Angiotensin Converting Enzyme (ACE)										●									
Anti-HAV																			
Anti-HBc										●									
Anti-HBe										●									
Anti-HBs										●									
Anti-HCV										●									
Anti-HIV-1										●									
Anti-HIV-1/2										●									
Anti-HIV-2										●									
Anti-HTLV-1										●									
Antibody Identification		●																	
Antibody Screen		●																	
Antigen Typing		●																	
Antistreptolysin O (ASO)												●							
Antithrombin III (ATIII)					●														
Anti-Thyroglobulin (Anti-Tg)													●						
Anti-Tyropoxidase (Anti-TPO)													●						
Apolipoprotein A-1											●								
Apolipoprotein B											●								
APTT					●														
Aspartate Aminotransferase (AST/SGOT)				●															
Bile Acids				●															
Bilirubin																		●	
Bilirubin (Direct)				●															
Bilirubin (Indirect)				●															
Bilirubin (Total)				●															
Blood/Hemoglobin																		●	
BNP			●																
C-Peptide										●			●						
C-Reactive Protein (CRP)												●							
CA 15-3										●									
CA 19-9										●									
CA 27.29										●									
CA 125										●									

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Control Analyte Index

EQAS™ Programs

See page number

- Analyte present
- ▶ Pilot analyte

	47	47	47	48	47	48	48	49	49	49	49	50	50	50	50	51	51	51
	Blood Gas	Blood Typing	Cardiac Markers	Clin. Chemistry (Monthly)	Coagulation	Ethanol/Ammonia	Hematology	Hemoglobin	HIV/Hepatitis	Immunoassay (Monthly)	Lipids	Serum Proteins	Specialty Immunoassay	Syphilis/Chagas	TDM	ToRCH/EBV/MuMz	Urinalysis	Urine Chemistry
Caffeine															●			
Calcitonin										▶								
Calcium (Ionized)	●			●														
Calcium (Total)				●														●
Carbamazepine										●					●			
Carbon Dioxide (CO ₂)				●														
Carcinoembryonic Antigen (CEA)										●								
Ceruloplasmin												●						
Chloride	●			●														●
Cholesterol (HDL)				●							●							
Cholesterol (LDL)				●							●							
Cholesterol (Total)				●							●							
Cholinesterase				●														
CK-MB			●															
Compatibility (Cross-match)		●																
Complement C3												●						
Complement C4												●						
Copper				●														
Cortisol				●						●								
Cortisol (Free)																		●
Creatine Kinase (CK)				●														
Creatinine				●												●		●
Cyclosporine															●			
Cytomegalovirus (CMV) IgG																●		
Cytomegalovirus (CMV) IgM																●		
DAT (Direct Antiglobulin Test)		●																
D-dimer			●		●													
DHEA										●								
DHEA Sulfate										●								
Digoxin										●					●			
Dopamine																		●
Epinephrine																		●
Epstein-Barr Virus (EBNA) IgG																●		
Epstein-Barr Virus (VCA) IgG																●		
Epstein-Barr Virus (VCA) IgM																●		
Erythropoietin (EPO)												●						
Estradiol										●								
Estradiol, Free (UE ₃)										●								
Ethanol						●												
Ethosuximide															●			
Factor VIII					●													
Ferritin										●								
Fibrinogen					●													
Folate										●								
Follicle Stimulating Hormone (FSH)										●								
Fructosamine										●		●						
Gamma Glutamyltransferase (GGT)				●														
Gastrin										●								
Gentamicin															●			
Glucose	●			●													●	●
Haptoglobin												●						
HAV IgG									●									
HAV IgM									●									
HBc IgM									●									
HBeAg									●									
HBSAg									●									
hCG									●									
Hematocrit (HCT)							●											
Hemoglobin A1C								●										

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

EQAS™ Programs

See page number

- Analyte present
- ▶ Pilot analyte

	47	47	47	48	47	48	48	49	49	49	49	50	50	50	50	51	51	51
	Blood Gas	Blood Typing	Cardiac Markers	Clin. Chemistry (Monthly)	Coagulation	Ethanol/Ammonia	Hematology	Hemoglobin	HIV/Hepatitis	Immunoassay (Monthly)	Lipids	Serum Proteins	Specialty Immunoassay	Syphilis/Chagas	TDM	ToRCH/EBV/MuM/Z	Urinalysis	Urine Chemistry
Hemoglobin A ₂								●										
Hemoglobin							●											
Hemoglobin (Total Glycated)								●										
Herpes Simplex Virus (HSV-1) IgG																		●
Herpes Simplex Virus (HSV-1/2) IgG																		●
Herpes Simplex Virus (HSV-2) IgG																		●
hGH										●								
HIV-1 Ag									●									
Homocysteine			●															
Homovanillic Acid (HVA)																		●
hs-CRP			●															
IgA												●						
IgE										●		●						
IgG										●		●						
IgM												●						
INR					●													
Insulin										●								
Insulin Growth Factor (IGF-1)													●					
Intact PTH										●			●					
Interleukin-6 (IL-6)													●					
Iron				●														
Iron Binding Capacity, Total (TIBC)				●														
Iron Binding Capacity, Unsaturated (UIBC)				●														
Kappa Light Chain												●						
Ketones																		●
Lactate (Lactic Acid)	●			●														
Lactate Dehydrogenase (LDH)				●														
Lambda Light Chain												●						
Leukocytes																		●
LH										●								
Lipase				●														
Lipoprotein (a)											●							
Lithium				●											●			
Magnesium	●			●														●
Mean Corpuscular Hemoglobin (MCH)							●											
Mean Corpuscular Hemoglobin Concentration (MCHC)							●											
Mean Corpuscular Volume (MCV)							●											
Mean Platelet Volume (MPV)							●											
Measles Virus (Rubeola) IgG																		●
Metanephrine																		●
Methotrexate															●			
Microalbumin																		●
Mumps Virus IgG																		●
Myoglobin			●															
Nitrite																		●
Non-Treponemal antibodies														●				
Norepinephrine																		●
Normetanephrine																		●
Nortriptyline															●			
NSE										●								
NT-proBNP			●															
Osmolality				●														●
Osteocalcin													●					
pCO ₂	●																	
pH	●																	●
pO ₂	●																	
Phenobarbital										●					●			
Phenytoin										●					●			
Phosphorus				●														●

Refer to myInsersts.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

Control Analyte Index

EQAS™ Programs

See page number

- Analyte present
- ▶ Pilot analyte

	47	47	47	48	47	48	48	49	49	49	49	50	50	50	50	51	51	51
	Blood Gas	Blood Typing	Cardiac Markers	Clin. Chemistry (Monthly)	Coagulation	Ethanol/Ammonia	Hematology	Hemoglobin	HIV/Hepatitis	Immunoassay (Monthly)	Lipids	Serum Proteins	Specialty Immunoassay	Syphilis/Chagas	TDM	ToRCH/EBV/MuMZ	Urinalysis	Urine Chemistry
Plasma Renin Activity (PRA)										●								
Platelets (PLT)							●											
Potassium	●			●														●
Pregnancy (hCG)																	●	
Prealbumin												●						
Primidone															●			
Procalcitonin													●					
Progesterone										●								
Prolactin										●								
Protein C					●													
Protein S					●													
Protein (Total)				●								●						●
Protein-to-Creatinine Ratio																	●	
PSA (Free)										●								
PSA (Free/Total Ratio)										●								
PSA (Total)										●								
PT					●													
Red Blood Cells (RBC)							●											
Red Blood Cell Distribution Width (RDW)							●											
Red Blood Cell Distribution Width (RDW-SD)							●											
Renin										●								
Rh(D)		●																
Rheumatoid Factor (RF)												●						
Rubella Virus IgG																	●	
Rubella Virus IgM																	●	
S100 Protein										●								
Salicylate															●			
Sex Hormone Binding Globulin (SHBG)										●								
Sodium	●			●														●
Specific Gravity																	●	
T3 Uptake										●								
T-Uptake										●								
Testosterone										●								
Testosterone (Free)										●								
Theophylline										●					●			
Thrombin Time (TT)					●													
Thyroglobulin (Tg)										●								
Thyroid Stimulating Hormone (TSH)				●						●								
Thyroxine Binding Globulin (TBG)										●								
Thyroxine, Free (T ₄)				●						●								
Thyroxine (T ₄), Total				●						●								
Tobramycin															●			
Toxoplasma gondii (Toxo) IgG																	●	
Toxoplasma gondii (Toxo) IgM																	●	
Transferrin				●						●		●						
Treponema pallidum IgG													●					
Treponema pallidum Total Antibody													●					
Tricyclic Antidepressants (TCA) Screen															●			
Triglycerides				●							●							
Triiodothyronine, Free (T ₃)				●						●								
Triiodothyronine (T ₃), Total				●						●								
Troponin I			●															
Troponin T			●															
Trypanosoma cruzi Total Antibody													●					
TSH Receptor Antibody (TRAb)												●						
Urea				●														▶
Urea Nitrogen (BUN)				●														●
Uric Acid				●														●
Urobilinogen																	●	

Refer to myInserTs.com or the product insert (IFU) of currently available lots for specific analyte and stability claims.

EQAS™ Programs

See page number

- Analyte present
- ▶ Pilot analyte

	47	47	47	48	47	48	48	49	49	49	49	50	50	50	50	51	51	51
	Blood Gas	Blood Typing	Cardiac Markers	Clin. Chemistry (Monthly)	Coagulation	Ethanol/Ammonia	Hematology	Hemoglobin	HIV/Hepatitis	Immunoassay (Monthly)	Lipids	Serum Proteins	Specialty Immunoassay	Syphilis/Chagas	TDM	ToRCH/EBV/MUMZ	Urinalysis	Urine Chemistry
Valproic Acid										●					●			
Vancomycin															●			
Vanillylmandelic Acid (VMA)																		●
Varicella Zoster Virus (VZV) IgG																●		
Vitamin B ₁₂										●								
von Willebrand Factor (vWF)					●													
White Blood Cells (WBC)							●											
Zinc				●														

A		F	
Additional Software Training, Unity	7, 8, 9	FAQs (QCNet)	2
Affiliated Data Exception Report, Unity	5	Fertility Control (Lyphocheck)	63
Affiliated Laboratory Comparison Reports, Unity	5	H	
Allergen sIgE Control (Lyphocheck)	67	Hematology Control (A) (Liquichek)	53
AMH Control (Liquichek)	63	Hematology Control (C) (Liquichek)	54
Amplichek II	29	Hematology Control (S) (Liquichek)	55
ANA Control, Homogeneous Pattern (Liquichek)	68	Hematology Control (X) (Liquichek)	56
ANA Controls Set, Positive: Homogeneous, Speckled, Centromere and Nucleolar Patterns (Liquichek)	68	Hematology Program (EQAS)	48
Anemia Control (Lyphocheck)	62	Hematology-16 Control (Liquichek)	57
Anti-Mitochondrial Control (Liquichek)	68	Hematology-16T Control (Liquichek)	57
Anti-nDNA Control (Liquichek)	68	Hemoglobin A1C Linearity Set (Lyphocheck)	46
Anti-Smooth Muscle Control (Liquichek)	68	Hemoglobin A ₂ Control (Lyphocheck)	45
Assayed Chemistry Control (Lyphocheck)	38	Hemoglobin Program (EQAS)	49
Assayed Multiquel Control (InteliQ)	11	Hemostasis Control (Lyphocheck)	44
Assayed VIROTROL I-C	69	HIV/Hepatitis Program (EQAS)	49
Assayed VIROTROL I-E	69	Homocysteine Control (InteliQ)	13
Assayed VIROTROL I-F	69	Homocysteine Control (Liquichek)	36
Assayed VIROTROL II-B	70	Hypertension Markers Control (Lyphocheck)	64
Autoimmune Negative Control (Liquichek)	68	I	
B		Immunoassay (Monthly) Program (EQAS)	49
Barcode Label Generator	2	Immunoassay Plus Control (InteliQ)	13
Benzo/TCA Control-Set B (Lyphocheck)	85	Immunoassay Plus Control (Liquichek)	59
Bias and Imprecision Histogram Report, Unity	5	Immunoassay Plus Control (Lyphocheck)	60
Bio-Rad Mission: Control	9	Immunoassay Premium Control (Liquichek)	64
Blood Gas Ampule Breaker	88	Immunology Control (InteliQ)	13
Blood Gas Control (Liquichek)	34	Immunology Control (Liquichek)	65
Blood Gas Plus E Control (Liquichek)	34	Immunology Plus Control (Lyphocheck)	65
Blood Gas Plus EGL Control (Liquichek)	34	InstantQC Report, Unity	5
Blood Gas Program (EQAS)	47	L	
Blood Typing Program (EQAS)	47	Laboratory Comparison Report, Unity	5
Blood Typing/Serum Indices Report, Unity	5	Laboratory Histogram Report, Unity	5
C		Laboratory Performance Overview Report, Unity	5
Cardiac Advance Control (Liquichek)	35	Linearity Web	2
Cardiac Advance Control (InteliQ)	11	Lipids Control (Liquichek)	40
Cardiac Markers Plus Control (Liquichek)	36	Lipids Program (EQAS)	49
Cardiac Markers Plus Control LT (InteliQ)	12	Liquid Assayed Multiquel	37
Cardiac Markers Plus Control LT (Liquichek)	35	Liquid Assayed Multiquel Premium	41
Cardiac Markers Program (EQAS)	47	Liquid Unassayed Multiquel	37
Cardiac Troponins Control (Liquichek)	36	M	
CK/LD Isoenzyme Control (Liquichek)	36	Maternal Serum 1st Trimester Control (Liquichek)	63
Clinical Chemistry (Monthly) Program (EQAS)	48	Maternal Serum II Control (Liquichek)	63
Coagulation Control (Lyphocheck)	44	Meter Trax Control	46
Coagulation Program (EQAS)	47	Microalbumin Control (Liquichek)	41
D		Monthly Evaluation Report, Unity	5
D-dimer Control (Liquichek)	44	My elnserts	2
Diabetes Control (Liquichek)	45	O	
Diabetes Control (InteliQ)	12	Opiate Control (Liquichek)	80
Diabetes Control (Lyphocheck)	45	P	
Dispenser Tips for Crimp Seals	88	Pediatric Control (Liquichek)	40
Dispenser Tips for Screw Caps	88	PyloriTrol	76
Drug Free Serum (Lyphocheck)	85	E	
E		elnserts XML	2
Elevated CRP Control (Liquichek)	66	Elevated CRP Control (Liquichek)	66
End-of-Cycle Report, EQAS	52	End-of-Cycle Report, EQAS	52
EQAS Resources on QCNet	2	EQAS Resources on QCNet	2
Ethanol/Ammonia Control (InteliQ)	12	Ethanol/Ammonia Control (InteliQ)	12
Ethanol/Ammonia Control (Liquichek)	39	Ethanol/Ammonia Control (Liquichek)	39
Ethanol/Ammonia Program (EQAS)	48	Ethanol/Ammonia Program (EQAS)	48

	Q	
QC Articles	2	
QC Lot Viewer	8	
QCNet	2	
QCNet International	2	
QC Webinars	2	
Qualitative Urine Toxicology Control (Liquichek)	79	
Qualitative Worldwide Report, Unity	5	
Quality Control Tray	88	
Quality Specification Report, EQAS	52	
qUAntify Advance Control	87	
qUAntify Plus Control	87	
Quantitative Urine Control (Lyphocheck)	42	
	R	
Reticulocyte Control (A-I) (Liquichek)	53	
Reticulocyte Control (A) (Liquichek)	53	
Reticulocyte Control (C) (Liquichek)	54	
Reticulocyte Control (S) (Liquichek)	55	
Reticulocyte Control (X) (Liquichek)	56	
Rheumatoid Factor Control (Liquichek)	66	
	S	
Sample Report, EQAS	52	
Sedimentation Rate Control (Liquichek)	58	
Serum Indices (Liquichek)	43	
Serum Proteins Program (EQAS)	50	
Serum Volatiles Control (Liquichek)	86	
Specialty Immunoassay Control (InteliQ)	14	
Specialty Immunoassay Plus Control (InteliQ)	14	
Specialty Immunoassay Control (Liquichek)	62	
Specialty Immunoassay Plus Control (Liquichek)	62	
Specialty Immunoassay Control (Lyphocheck)	62	
Specialty Immunoassay Program (EQAS)	50	
Spinal Fluid Control (Liquichek)	67	
Statistical Profile Report, Unity	5	
Syphilis/Chagas Program (EQAS)	50	
	T	
Target CRP Levels in Bio-Rad Controls (Chart)	66	
Therapeutic Drug Monitoring Program (TDM) (EQAS)	50	
Technical Documents	2	
Therapeutic Drug Monitoring Control (TDM) (Liquichek)	77	
Therapeutic Drug Monitoring Control (TDM) (Lyphocheck)	77	
ToRCH/EBV/MuMZ Program (EQAS)	51	
ToRCH Plus Control (Liquichek)	73	
Training and Education	2	
Tumor Marker Control (InteliQ)	14	
Tumor Marker Control (Liquichek)	61	
Tumor Marker Plus Control (Lyphocheck)	61	
	U	
Unassayed Chemistry Control (Human) (Liquichek)	39	
Unassayed Chemistry Control (Human) (Lyphocheck)	38	
Unity Alert	9	
Unity Binder	88	
UnityConnect	6, 8	
Unity Interlaboratory Reports	6	
Unity Real Time	6, 8	
Unity Real Time LT	6, 7	
Unity Real Time online	6, 8	
UnityWeb	6, 7	
Urinalysis Control (Liquichek)	87	
Urinalysis Program (EQAS)	51	
Urinalysis Reports, Unity	5	
Urinalysis Tools	2	
Urine Chemistry Control (InteliQ)	14	
Urine Chemistry Control (Liquichek)	42	
Urine Chemistry Program (EQAS)	48	
Urine Metals Control (Lyphocheck)	86	
Urine Toxicology Control (Liquichek), Level C1	84	
Urine Toxicology Control (Liquichek), Level C2	84	
Urine Toxicology Control (Liquichek), Level C2 Low Opiate	84	
Urine Toxicology Control (Liquichek), Level C3	84	
Urine Toxicology Control (Liquichek), Level C3 Low Opiate	84	
Urine Toxicology Control (Liquichek), Level C4	84	
Urine Toxicology Control (Liquichek), Level S1	81	
Urine Toxicology Control (Liquichek), Level S1 Low Opiate	82	
Urine Toxicology Control (Liquichek), Level S1E	81	
Urine Toxicology Control (Liquichek), Level S1E Low Opiate	82	
Urine Toxicology Control (Liquichek), Level S1O Low Opiate	82	
Urine Toxicology Control (Liquichek), Level S1S	83	
Urine Toxicology Control (Liquichek), Level S2	81	
Urine Toxicology Control (Liquichek), Level S2 Low Opiate	82	
Urine Toxicology Control (Liquichek), Level S2E	81	
Urine Toxicology Control (Liquichek), Level S2E Low Opiate	82	
Urine Toxicology Control (Liquichek), Level S2O Low Opiate	82	
Urine Toxicology Control (Liquichek), Level S2S	83	
Urine Toxicology Control (Liquichek), Level S3	83	
Urine Toxicology Negative Control (Liquichek)	79	
	V	
Value Assignment Program	2	
VIROCLEAR	72	
VIROCLEAR EBV	76	
VIROCLEAR MuMZ	73	
VIROCLEAR SARS-CoV-2	74	
VIROCLEAR ToRCH	73	
VIROTROL Chagas	75	
VIROTROL EBV	75	
VIROTROL HAV-IgM	71	
VIROTROL HBeAg	71	
VIROTROL HIV-1 Ag	70	
VIROTROL HIV-1 gO	71	
VIROTROL HIV-2	71	
VIROTROL I	69	
VIROTROL II	69	
VIROTROL III	70	
VIROTROL IV	70	
VIROTROL Lyme	76	
VIROTROL MuMZ	73	
VIROTROL PLUS-R	71	
VIROTROL RPR Panel	75	
VIROTROL SARS-CoV-2	74	
VIROTROL Syphilis LR-A	75	
VIROTROL ToRCH	72	
VIROTROL ToRCH-M	72	

W

Westgard Advisor	6, 9
WebConnect	6, 7
Whole Blood Control (Lyphochek)	86
Whole Blood Immunosuppressant Control, (Liquichek)	78
Whole Blood Immunosuppressant Control (Lyphochek)	78
Whole Blood Metals Control (Lyphochek)	85
Worldwide Report, Unity	5

Y

YouTube Channel (youtube.com/bioradqc)	2
---	---

Exact Diagnostics

Adenovirus (Exact Diagnostics)	16
Analytical Measurement Range Bundle (Exact Diagnostics)	24
Anaplasma (Exact Diagnostics)	32
Babesia (Exact Diagnostics)	32
Bartonella (Exact Diagnostics)	32
BKV (Exact Diagnostics)	16
Borrelia (Exact Diagnostics)	32
Bulk CSF Negative (Exact Diagnostics)	26
Bulk Urine Negative (Exact Diagnostics)	26
<i>C. difficile</i> Verification Bundle (Exact Diagnostics)	28
CMV (Exact Diagnostics)	17
CT/NG (Exact Diagnostics)	19
EBV (Exact Diagnostics)	18
Ehrlichia (Exact Diagnostics)	33
Enterovirus (Exact Diagnostics)	27
GBS (Exact Diagnostics)	27
HAI Negative (Exact Diagnostics)	25
HBV (Exact Diagnostics)	22
HBV Analytical Measurement Range Panel (Exact Diagnostics)	24
HBV, HCV and HIV-1 Screen Controls (Exact Diagnostics)	29
HCV (Exact Diagnostics)	22
HCV Analytical Measurement Range Panel (Exact Diagnostics)	24
HHV-6 (Exact Diagnostics)	18
HIV (Exact Diagnostics)	23
HIV Analytical Measurement Range Panel (Exact Diagnostics)	24
HPV (Exact Diagnostics)	20
HSV-1 (Exact Diagnostics)	21
HSV-2 (Exact Diagnostics)	21
JCV (Exact Diagnostics)	19
MRSA Verification Bundle (Exact Diagnostics)	28
Negative (Exact Diagnostics)	26
Parvovirus B19 (Exact Diagnostics)	17
RP Positive & Negative (Exact Diagnostics)	31
SARS-CoV-2, Flu, RSV (Exact Diagnostics)	30
SARS-CoV-2 S Gene Variant (Exact Diagnostics)	31
SARS-CoV-2 (Exact Diagnostics)	30
STI Negative (Exact Diagnostics)	25
Synthetic Negative (Exact Diagnostics)	25
TV/MG (Exact Diagnostics)	20
VZV (Exact Diagnostics)	27
ZKV (Exact Diagnostics)	33

100

144 Hematology-16T Control (Liquichek), Trilevel. 57

144X Hematology-16T Control (Liquichek), Trilevel MiniPak. 57

146 Cardiac Markers Plus Control LT (Liquichek),
Level 1 (3 mL) 35

146X Cardiac Markers Plus Control LT (Liquichek),
Level 1 MiniPak (3 mL) 35

147 Cardiac Markers Plus Control LT (Liquichek),
Level 2 (3 mL) 35

147X Cardiac Markers Plus Control LT (Liquichek),
Level 2 MiniPak (3 mL) 35

148 Cardiac Markers Plus Control LT (Liquichek),
Level 3 (3 mL) 35

148X Cardiac Markers Plus Control LT (Liquichek),
Level 3 MiniPak (3 mL) 35

149 Cardiac Markers Plus Control LT (Liquichek),
Level Low (3 mL) 35

149X Cardiac Markers Plus Control LT (Liquichek),
Level Low MiniPak (3 mL) 35

176 Pediatric Control (Liquichek), Level 1 (2.5 mL) 40

177 Pediatric Control (Liquichek), Level 2 (2.5 mL) 40

178X Pediatric Control (Liquichek), Bilevel MiniPak (2.5 mL) 40

180 Cardiac Markers Plus Control (Liquichek), Trilevel. 36

180X Cardiac Markers Plus Control (Liquichek),
Trilevel MiniPak 36

181 Cardiac Markers Plus Control (Liquichek), Level 1 36

182 Cardiac Markers Plus Control (Liquichek), Level 2 36

183 Cardiac Markers Plus Control (Liquichek), Level 3 36

195 Urine Chemistry Control (Liquichek), Level 1 (2.5 mL) 42

196 Urine Chemistry Control (Liquichek), Level 2 (2.5 mL) 42

198X Urine Chemistry Control (Liquichek),
Bilevel MiniPak (2.5 mL) 42

200

227 ToRCH Plus Control (Liquichek), Positive 73

227X ToRCH Plus Control (Liquichek), Positive, MiniPak 73

228 ToRCH Plus Control (Liquichek), Negative 73

228X ToRCH Plus Control (Liquichek), Negative MiniPak 73

239 ToRCH Plus Control (Liquichek), Positive 73

239X ToRCH Plus Control (Liquichek), Positive MiniPak 73

254 Elevated CRP Control (Liquichek), Level 1 66

255 Elevated CRP Control (Liquichek), Level 2 66

255X Elevated CRP Control (Liquichek), Trilevel MiniPak 66

256 Elevated CRP Control (Liquichek), Level 3 66

267 Immunoassay Plus Control (Liquichek), Level 1 (2 mL) 59

268 Immunoassay Plus Control (Liquichek), Level 2 (2 mL) 59

268X Immunoassay Plus Control (Liquichek),
Trilevel MiniPak (2 mL) 59

269 Immunoassay Plus Control (Liquichek), Level 3 (2 mL) 59

271 Ethanol/Ammonia Control (Liquichek), Level 1 (2.5 mL) 39

272 Ethanol/Ammonia Control (Liquichek), Level 2 (2.5 mL) 39

272X Ethanol/Ammonia Control (Liquichek),
Trilevel MiniPak (2.5 mL) 39

273 Ethanol/Ammonia Control (Liquichek),
Level 3 (2.5 mL) 39

274 Whole Blood Immunosuppressant Control
(Lyphocek), Level 1 78

275 Whole Blood Immunosuppressant Control (Lyphocek),
Level 2 78

276 Whole Blood Immunosuppressant Control (Lyphocek),
Level 3 78

277 Whole Blood Immunosuppressant Control (Lyphocek),
Level 4 78

278 Whole Blood Immunosuppressant Control (Lyphocek),
Level 5 78

279X Whole Blood Immunosuppressant Control (Lyphocek),
Five Level MiniPak 78

281 Immunology Control (Liquichek), Level 1 (2.5 mL) 65

282 Immunology Control (Liquichek), Level 2 (2.5 mL) 65

282X Immunology Control (Liquichek),
Trilevel MiniPak (2.5 mL) 65

283 Immunology Control (Liquichek), Level 3 (2.5 mL) 65

284 Liquid Assayed Multiquel Premium, Level 1 41

285 Liquid Assayed Multiquel Premium, Level 2 41

285X Liquid Assayed Multiquel Premium, Trilevel MiniPak 41

286 Liquid Assayed Multiquel Premium, Level 3 41

287 Homocysteine Control (Liquichek), Level 1 36

288 Homocysteine Control (Liquichek), Level 2 36

288X Homocysteine Control (Liquichek), Trilevel MiniPak 36

289 Homocysteine Control (Liquichek), Level 3 36

297 Cardiac Markers Plus Control LT (Liquichek),
Level 1C (2.5 mL) 35

298 Cardiac Markers Plus Control LT (Liquichek),
Level 2 (2.5 mL) 35

298X Cardiac Markers Plus Control LT (Liquichek),
Trilevel MiniPak (2.5 mL) 35

299 Cardiac Markers Plus Control LT (Liquichek),
Level 3 (2.5 mL) 35

300

302X Spinal Fluid Control (Liquichek), Bilevel MiniPak (2.5 mL) 67

303 Spinal Fluid Control (Liquichek), Level 1 (2.5 mL) 67

304 Spinal Fluid Control (Liquichek), Level 2 (2.5 mL) 67

313X Assayed Chemistry Control (Lyphocek), Bilevel MiniPak. 38

323X Unassayed Chemistry Control (Human) (Lyphocek)
Bilevel MiniPak (10 mL) 38

337 Liquid Assayed Multiquel, Level 1 (2.5 mL) 37

338 Liquid Assayed Multiquel, Level 2 (2.5 mL) 37

338X Liquid Assayed Multiquel, Trilevel MiniPak (2.5 mL) 37

339 Liquid Assayed Multiquel, Level 3 (2.5 mL) 37

353X Pediatric Control (Liquichek), Bilevel MiniPak (4 mL) 40

354 Pediatric Control (Liquichek), Level 1 (4 mL) 40

355 Pediatric Control (Liquichek), Level 2 (4 mL) 40

359 Specialty Immunoassay Control (Liquichek), Level LTA 62

359X Specialty Immunoassay Control (Liquichek),
Four Level MiniPak 62

360 Immunoassay Plus Control (Liquichek), Trilevel (5 mL) 59

360X Immunoassay Plus Control (Liquichek),
Trilevel MiniPak (5 mL) 59

361 Immunoassay Plus Control (Liquichek), Level 1 (5 mL) 59

362 Immunoassay Plus Control (Liquichek), Level 2 (5 mL) 59

363 Immunoassay Plus Control (Liquichek), Level 3 (5 mL) 59

364 Specialty Immunoassay Control (Liquichek), Level 1 62

365 Specialty Immunoassay Control (Liquichek), Level 2 62

366 Specialty Immunoassay Control (Liquichek), Level 3 62

367 Tumor Marker Plus Control (Lyphocek), Level 1 61

368 Tumor Marker Plus Control (Lyphocek), Level 2 61

368X Tumor Marker Plus Control (Lyphocek), Trilevel MiniPak. 61

369 Tumor Marker Plus Control (Lyphocek), Level 3 61

370 Immunoassay Plus Control (Lyphocek), Trilevel. 60

370X Immunoassay Plus Control (Lyphocek), Trilevel MiniPak. 60

371 Immunoassay Plus Control (Lyphocek), Level 1 60

372 Immunoassay Plus Control (Lyphocek), Level 2 60

373 Immunoassay Plus Control (Lyphocek), Level 3 60

374X Microalbumin Control (Liquichek), Bilevel MiniPak 41

375X Quantitative Urine Control (Lyphocek), Bilevel MiniPak 42

376 Quantitative Urine Control (Lyphocek), Level 1, Normal 42

377 Quantitative Urine Control (Lyphocek),
Level 2, Abnormal 42

378 Microalbumin Control (Liquichek), Level 1 41

379	Microalbumin Control (Liquichek), Level 2	41	469	Urine Toxicology Control (Liquichek), Level C3 Low Opiate	84
383	Serum Volatiles Control (Liquichek), Level 1	86	480	Fertility Control (Lyphocek), Trilevel	63
384	Serum Volatiles Control (Liquichek), Level 2	86	480X	Fertility Control (Lyphocek), Trilevel MiniPak	63
385X	Serum Volatiles Control (Liquichek), Bilevel MiniPak	86	481	Fertility Control (Lyphocek), Level 1	63
395X	Urine Chemistry Control (Liquichek), Bilevel MiniPak (10 mL)	42	482	Fertility Control (Lyphocek), Level 2	63
397	Urine Chemistry Control (Liquichek), Level 1 (10 mL)	42	483	Fertility Control (Lyphocek), Level 3	63
398	Urine Chemistry Control (Liquichek), Level 2 (10 mL)	42	487	Hematology Control (X) (Liquichek), Trilevel	56
	400		487X	Hematology Control (X) (Liquichek), Trilevel MiniPak	56
400	Urine Metals Control (Lyphocek), Level 1	86		500	
402	Maternal Serum II (Liquichek), Level 1	63	500	Anemia Control (Lyphocek), Single Level	62
402X	Urine Metals Control (Lyphocek), Bilevel MiniPak	86	500X	Anemia Control (Lyphocek), MiniPak	62
403	Maternal Serum II Control (Liquichek), Level 2	63	501	Rheumatoid Factor Control (Liquichek), Level 1	66
403X	Maternal Serum II Control (Liquichek), Trilevel MiniPak	63	502	Rheumatoid Factor Control (Liquichek), Level 2	66
404	Maternal Serum II Control (Liquichek), Level 3	63	502X	Rheumatoid Factor Control (Liquichek), Trilevel MiniPak	66
405	Urine Metals Control (Lyphocek), Level 2	86	503	Rheumatoid Factor Control (Liquichek), Level 3	66
409	Urine Toxicology Control (Liquichek), Level S1S	83	510X	Blood Gas Plus EGL Control (Liquichek), Trilevel MiniPak	34
419	Urine Toxicology Control (Liquichek), Level S2S	83	511	Blood Gas Plus EGL Control (Liquichek), Level 1	34
423	Urine Toxicology Control (Liquichek), Level S1E Low Opiate	82	512	Blood Gas Plus EGL Control (Liquichek), Level 2	34
424	Urine Toxicology Control (Liquichek), Level S2E Low Opiate	82	513	Blood Gas Plus EGL Control (Liquichek), Level 3	34
430	Immunology Plus Control (Lyphocek), Bilevel	65	514	Sedimentation Rate Control (Liquichek), Level 1	58
430X	Immunology Plus Control (Lyphocek), Bilevel MiniPak	65	514X	Sedimentation Rate Control (Liquichek), Bilevel MiniPak	58
435	Urinalysis Control (Liquichek), Bilevel	87	515	Sedimentation Rate Control (Liquichek), Level 2	58
435X	Urinalysis Control (Liquichek), Bilevel MiniPak	87	527	Whole Blood Metals Control (Lyphocek), Level 1	85
436	Urinalysis Control (Liquichek), Level 1	87	528	Whole Blood Metals Control (Lyphocek), Level 2	85
437	Urinalysis Control (Liquichek), Level 2	87	528X	Whole Blood Metals Control (Lyphocek), Trilevel MiniPak	85
438	Urine Toxicology Control (Liquichek), Level S1E	81	529	Whole Blood Metals Control (Lyphocek), Level 3	85
439	Urine Toxicology Control (Liquichek), Level S2E	81	544	Ethanol/Ammonia Control (Liquichek), Level 1 (3 mL)	39
441	Urine Toxicology Control (Liquichek), Level C1	84	545	Ethanol/Ammonia Control (Liquichek), Level 2 (3 mL)	39
442	Urine Toxicology Control (Liquichek), Level C2	84	545X	Ethanol/Ammonia Control (Liquichek), Trilevel MiniPak (3 mL)	39
443	Urine Toxicology Control (Liquichek), Level C3	84	546	Ethanol/Ammonia Control (Liquichek), Level 3 (3 mL)	39
444	Urine Toxicology Control (Liquichek), Level C4	84	547	Tumor Marker Control (Liquichek), Level 1 (2 mL)	61
449	Hematology Control (S) (Liquichek), Trilevel	55	548	Tumor Marker Control (Liquichek), Level 2 (2 mL)	61
449X	Hematology Control (S) (Liquichek), Trilevel MiniPak	55	548X	Tumor Marker Control (Liquichek), Trilevel MiniPak (2 mL)	61
450	Therapeutic Drug Monitoring Control (TDM) (Lyphocek), Trilevel	77	549	Tumor Marker Control (Liquichek), Level 3 (2 mL)	61
450X	Therapeutic Drug Monitoring Control (TDM) (Lyphocek), Trilevel MiniPak	77	553	Hemoglobin A ₂ Control (Lyphocek), Bilevel	45
451	Therapeutic Drug Monitoring Control (TDM) (Lyphocek), Level 1	77	553X	Hemoglobin A ₂ Control (Lyphocek), Bilevel MiniPak	45
452	Therapeutic Drug Monitoring Control (TDM) (Lyphocek), Level 2	77	560X	Whole Blood Control (Lyphocek), Trilevel MiniPak	86
453	Therapeutic Drug Monitoring Control (TDM) (Lyphocek), Level 3	77	561	Whole Blood Control (Lyphocek), Level 1	86
454	Qualitative Urine Toxicology Control (Liquichek), Negative Level	79	562	Whole Blood Control (Lyphocek), Level 2	86
454X	Qualitative Urine Toxicology Control (Liquichek), Bilevel MiniPak	79	563	Whole Blood Control (Lyphocek), Level 3	86
455	Qualitative Urine Toxicology Control (Liquichek), Positive Level	79	570X	Blood Gas Control (Liquichek), Trilevel MiniPak	34
456	Drug Free Serum (Lyphocek)	85	571	Blood Gas Control (Liquichek), Level 1	34
456X	Drug Free Serum (Lyphocek), MiniPak	85	572	Blood Gas Control (Liquichek), Level 2	34
460	Urine Toxicology Negative Control (Liquichek)	79	573	Blood Gas Control (Liquichek), Level 3	34
460X	Urine Toxicology Negative Control (Liquichek), MiniPak	79	575X	Blood Gas Plus E Control (Liquichek), Trilevel MiniPak	34
461	Urine Toxicology Control (Liquichek), Level S1	81	576	Blood Gas Plus E Control (Liquichek), Level 1	34
462	Urine Toxicology Control (Liquichek), Level S2	81	577	Blood Gas Plus E Control (Liquichek), Level 2	34
463	Urine Toxicology Control (Liquichek), Level S3	83	578	Blood Gas Plus E Control (Liquichek), Level 3	34
466	Urine Toxicology Control (Liquichek), Level S1 Low Opiate	82	587	Blood Gas Ampule Breaker	88
467	Urine Toxicology Control (Liquichek), Level S2 Low Opiate	82	588	Dispenser Tips for Crimp Seals	88
468	Urine Toxicology Control (Liquichek), Level C2 Low Opiate	84	590X	Immunology Control (Liquichek), Trilevel MiniPak (1 mL)	65
			591	Immunology Control (Liquichek), Level 1 (1 mL)	65
			592	Immunology Control (Liquichek), Level 2 (1 mL)	65
			593	Immunology Control (Liquichek), Level 3 (1 mL)	65
			594	Immunology Control (Liquichek), Level 1 (3 mL)	65
			595	Immunology Control (Liquichek), Level 2 (3 mL)	65
			596	Immunology Control (Liquichek), Level 3 (3 mL)	65
			597	Hemostasis Control (Lyphocek), Level 1	44
			598	Hemostasis Control (Lyphocek), Level 2	44
			598X	Hemostasis Control (Lyphocek), Trilevel MiniPak	44
			599	Hemostasis Control (Lyphocek), Level 3	44

600

600	Hypertension Markers Control (Lyphocheck), Trilevel	64
600X	Hypertension Markers Control (Lyphocheck), Trilevel MiniPak	64
636	Maternal Serum 1st Trimester Control (Liquichek), Trilevel . .	63
636X	Maternal Serum 1st Trimester Control (Liquichek), Trilevel MiniPak	63
640X	Lipids Control (Liquichek), Bilevel MiniPak	40
641	Lipids Control (Liquichek), Level 1	40
642	Lipids Control (Liquichek), Level 2	40
663	Benzo/TCA Control Set B (Lyphocheck), Level 1	85
664	Benzo/TCA Control Set B (Lyphocheck), Level 2	85
665X	Benzo/TCA Control Set B (Lyphocheck), Bilevel MiniPak	85
670X	CK/LD Isoenzyme Control (Liquichek), Bilevel MiniPak	36
671	CK/LD Isoenzyme Control (Liquichek), Level 1	36
672	CK/LD Isoenzyme Control (Liquichek), Level 2	36
683	Urine Toxicology Control (Liquichek), Level S10 Low Opiate	82
690X	Unassayed Chemistry Control (Human) (Liquichek), Bilevel MiniPak	39
691	Unassayed Chemistry Control (Human) (Liquichek), Level 1	39
692	Unassayed Chemistry Control (Human) (Liquichek), Level 2	39
694	Liquid Assayed Multiqual, Level 1 (3 mL)	37
695	Liquid Assayed Multiqual, Level 2 (3 mL)	37
695X	Liquid Assayed Multiqual, Trilevel MiniPak (3 mL)	37
696	Liquid Assayed Multiqual, Level 3 (3 mL)	37
697	Liquid Unassayed Multiqual, Level 1	37
698	Liquid Unassayed Multiqual, Level 2	37
698X	Liquid Unassayed Multiqual, Trilevel MiniPak	37
699	Liquid Unassayed Multiqual, Level 3	37
700		
706	Urine Toxicology Control (Liquichek), Level S20 Low Opiate	82
724	Therapeutic Drug Monitoring Control (TDM) (Liquichek), Trilevel	77
724X	Therapeutic Drug Monitoring Control (TDM) (Liquichek), Trilevel MiniPak	77
725	Therapeutic Drug Monitoring Control (TDM) (Liquichek), Level 1	77
726	Therapeutic Drug Monitoring Control (TDM) (Liquichek), Level 2	77
727	Therapeutic Drug Monitoring Control (TDM) (Liquichek), Level 3	77
730X	Unassayed Chemistry Control (Human) (Lyphocheck), Bilevel MiniPak (5 mL)	38
731	Unassayed Chemistry Control (Human) (Lyphocheck), Level 1 (5 mL)	38
732	Unassayed Chemistry Control (Human) (Lyphocheck), Level 2 (5 mL)	38
740	Diabetes Control (Lyphocheck), Bilevel	45
740X	Diabetes Control (Lyphocheck), Bilevel MiniPak	45
744	Coagulation Control (Lyphocheck), Level 1	44
745	Coagulation Control (Lyphocheck), Level 2	44
745X	Coagulation Control (Lyphocheck), Trilevel MiniPak	44
746	Coagulation Control (Lyphocheck), Level 3	44
750X	Spinal Fluid Control (Liquichek), Bilevel MiniPak (3 mL)	67
751	Spinal Fluid Control (Liquichek), Level 1 (3 mL)	67
752	Spinal Fluid Control (Liquichek), Level 2 (3 mL)	67
760	Hematology-16 Control (Liquichek), Trilevel	57
760X	Hematology-16 Control (Liquichek), Trilevel MiniPak	57
761	Hematology-16 Control (Liquichek), Low	57
762	Hematology-16 Control (Liquichek), Normal	57
763	Hematology-16 Control (Liquichek), High	57

800

804-1	Unity Real Time (Annual Subscription)	8
804-W1	Unity Real Time online (Annual Subscription)	8
805-1	Unity Real Time LT	7
806-1	Unity Alert (Annual Subscription)	9
811-1	Westgard Advisor (Annual Subscription)	9
825i	Unity Remote Installation & Training	7, 8, 9
826	Reticulocyte Control (C) (Liquichek), Level 1	54
827	Reticulocyte Control (C) (Liquichek), Level 2	54
827X	Reticulocyte Control (C) (Liquichek), Trilevel MiniPak	54
828	Reticulocyte Control (C) (Liquichek), Level 3	54
854	Reticulocyte Control (A) (Liquichek), Level 1	53
855	Reticulocyte Control (A) (Liquichek), Level 2	53
855X	Reticulocyte Control (A) (Liquichek), Bilevel MiniPak	53
858i	Additional Software Training	7, 8, 9
865	Hematology Control (A) (Liquichek), Trilevel	53
865X	Hematology Control (A) (Liquichek), Trilevel MiniPak	53
866	Hematology Control (A) (Liquichek), Low	53
867	Hematology Control (A) (Liquichek), Normal	53
868	Hematology Control (A) (Liquichek), High	53
870-1	UnityWeb (Annual Subscription)	7
870-W1	WebConnect (Annual Subscription)	7

900

904	Hematology Control (C) (Liquichek), Trilevel	54
904X	Hematology Control (C) (Liquichek), Trilevel MiniPak	54
905	Hematology Control (C) (Liquichek), Low	54
906	Hematology Control (C) (Liquichek), Normal	54
907	Hematology Control (C) (Liquichek), High	54
915X	Reticulocyte Control (A-I) (Liquichek), Bilevel MiniPak	53
916	Reticulocyte Control (A-I) (Liquichek), Level 1	53
917	Reticulocyte Control (A-I) (Liquichek), Level 2	53
921	Reticulocyte Control (S) (Liquichek), Level 1	55
922	Reticulocyte Control (S) (Liquichek), Level 2	55
922X	Reticulocyte Control (S) (Liquichek), Trilevel MiniPak	55
923	Reticulocyte Control (S) (Liquichek), Level 3	55
926	Reticulocyte Control (X) (Liquichek), Level 1	56
927	Reticulocyte Control (X) (Liquichek), Level 2	56
927X	Reticulocyte Control (X) (Liquichek), Trilevel MiniPak	56
928	Reticulocyte Control (X) (Liquichek), Level 3	56
962	qUAntify Plus Control, Bilevel, Bottle	87
962X	qUAntify Plus Control, Bilevel MiniPak, Bottle	87
970	Meter Trax Control, Trilevel	46
970X	Meter Trax Control, Trilevel MiniPak	46
971	Meter Trax Control, Low Level	46
972	Meter Trax Control, Mid Level	46
973	Meter Trax Control, High Level	46
995	qUAntify Plus Control, Bilevel	87
995X	qUAntify Plus Control, Bilevel MiniPak	87
987000	Dispenser Tips for Screw Caps	88

	00100	
00100A	VIROTROL I (5 mL)	69
00100B	VIROTROL I (5 mL)	69
00100C	VIROTROL I (5 mL)	69
00100E	VIROTROL I (5 mL)	69
00100F	VIROTROL I (5 mL)	69
00100H	VIROTROL I (5 mL)	69
00101B	VIROTROL I (4 mL)	69
00101C	VIROTROL I (4 mL)	69
00101E	VIROTROL I (4 mL)	69
00101F	VIROTROL I (4 mL)	69
00104A	VIROTROL II (5 mL)	69
00104B	VIROTROL II (5 mL)	69
00105A	VIROTROL HIV-2	71
00105B	VIROTROL HIV-2	71
00105C	VIROTROL HIV-2	71
00106	VIROCLEAR (5 mL)	72
00108A	VIROTROL HIV-1 Ag	70
00108B	VIROTROL HIV-1 Ag	70
00109	VIROTROL ToRCH	70
00110A	VIROTROL III	70
00110C	VIROTROL III	70
00111	VIROTROL IV	70
00112	VIROCLEAR (4 mL)	72
00113	VIROTROL HIV-1 gO (4 tubes)	71
00113X	VIROTROL HIV-1 gO (1 tube)	71
00115A	VIROTROL II (4 mL)	66
00115B	VIROTROL II (4 mL)	69
00117A	VIROTROL ToRCH-M	72
00117B	VIROTROL ToRCH-M	72
00118	VIROCLEAR ToRCH	73
00119	VIROTROL MuMZ	73
00121	PyloriTrol	76
00125	VIROTROL Chagas	75
00125X	VIROTROL Chagas	75
00128A	VIROTROL EBV	75
00131	VIROCLEAR EBV	76
00132	VIROTROL Lyme	76
00133	VIROCLEAR MuMZ	70
00142A	VIROTROL HAV-IgM	71
00144	VIROTROL HBeAg	71
00152	Assayed VIROTROL II-B	70
00164	Assayed VIROTROL I-C	69
00166	Assayed VIROTROL I-E	69
00168	Assayed VIROTROL I-F	69
00171	VIROTROL Syphilis LR-A	75
00171X	VIROTROL Syphilis LR-A	75
01400	VIROTROL RPR Panel	75
	27000	
27100	D-dimer Control (Liquichek), Level Low	44
27101	D-dimer Control (Liquichek), Level 1	44
27102	D-dimer Control (Liquichek), Level 2	44
27102X	D-dimer Control (Liquichek), Four Level MiniPak	44
27103	D-dimer Control (Liquichek), Level 3	44
27104	Cardiac Markers Plus Control LT (Liquichek), Level 1A (3 mL)	35
27104X	Cardiac Markers Plus Control LT (Liquichek), Level 1A MiniPak (3 mL)	35
27105	Cardiac Markers Plus Control LT (Liquichek), Level 1B (3 mL)	35
27105X	Cardiac Markers Plus Control LT (Liquichek), Level 1B MiniPak (3 mL)	35
27106	Cardiac Markers Plus Control LT (Liquichek), Level 1C (3 mL)	35
27106X	Cardiac Markers Plus Control LT (Liquichek), Level 1C MiniPak (3 mL)	35
27110	Immunoassay Premium Control (Liquichek), Trilevel	64
27110X	Immunoassay Premium Control (Liquichek), Level 1	64
27111	Immunoassay Premium Control (Liquichek), Level 2	64
27112	Immunoassay Premium Control (Liquichek), Level 3	64
27113	Immunoassay Premium Control (Liquichek), Trilevel MiniPak	64
27114	Tumor Marker Control (Liquichek), Level 1 (2 mL)	61
27115	Tumor Marker Control (Liquichek), Level 2 (2 mL)	61
27115X	Tumor Marker Control (Liquichek), Trilevel MiniPak (2 mL)	61
27116	Tumor Marker Control (Liquichek), Level 3 (2 mL)	61
27124	Specialty Immunoassay Control (Lyphocheck), Level 1	62
27125	Specialty Immunoassay Control (Lyphocheck), Level 2	62
27125X	Specialty Immunoassay Control (Lyphocheck), Trilevel MiniPak	62
27126	Specialty Immunoassay Control (Lyphocheck), Level 3	62
	200000	
200300A	VIROTROL SARS-CoV-2	74
200305A	VIROTROL SARS-CoV-2	74
200500	VIROCLEAR SARS-CoV-2	74
200505	VIROCLEAR SARS-CoV-2	74
	1100000	
1100002	Bio-Rad Mission: Control, Annual Subscription	9
1100005	Bio-Rad Mission: Control, Subscription and Service Bundle	9
	11000000	
11000009	UnityConnect (Annual Subscription)	8
11000506	Unity Next Onboarding	7
11000517	Unity Next Peer QC (Annual Subscription)	7
11001897	Unity Next File Upload (Annual Subscription)	7
	12000000	
12000070	Hemoglobin A1C Linearity Set (Lyphocheck)	46
12000141	Allergen sIgE Control (Lyphocheck), Negative	67
12000142	Allergen sIgE Control (Lyphocheck), Panel A	67
12000143	Allergen sIgE Control (Lyphocheck), Negative MiniPak	67
12000144	Allergen sIgE Control (Lyphocheck), Panel A MiniPak	67
12000201	Amplichek II, Negative	29
12000202	Amplichek II, Level 1	29
12000203	Amplichek II, Level 2	29
12000204	Amplichek II, Level 3	29
12000404	Whole Blood Immunosuppressant Control (Liquichek), Level 1	78
12000405	Whole Blood Immunosuppressant Control (Liquichek), Level 2	78
12000406	Whole Blood Immunosuppressant Control (Liquichek), Level 3	78
12000407	Whole Blood Immunosuppressant Control (Liquichek), Level 4	78
12000408	Whole Blood Immunosuppressant Control (Liquichek), Four Level MiniPak	78
12000538	VIROTROL PLUS-R (5 x 4 mL)	71
12000539	VIROTROL PLUS-R (1 x 4 mL)	71
12000575	Amplichek II, Negative MiniPak	29
12000576	Amplichek II, Level 1 MiniPak	29
12000577	Amplichek II, Level 2 MiniPak	29
12000578	Amplichek II, Level 3 MiniPak	29
12000810	Urinalysis Program (EQAS)	51
12000815	HIV/Hepatitis Program (EQAS)	49
12000825	ToRCH/EBV/MuMZ Program (EQAS)	51
12000833	Syphilis/Chagas Program (EQAS)	50
12000837	Blood Typing Program (EQAS), Shipment A	47
12000838	Blood Typing Program (EQAS), Shipment B	47
12000839	Blood Typing Program (EQAS), Shipment C	47
12000846	Quality Control Tray	88
12001084	Cardiac Troponins Control (Liquichek), Level 1	36
12001085	Cardiac Troponins Control (Liquichek), Level 1 MiniPak	36

12002594	Opiate Control (Liquichek), Level 1 MiniPak	80
12002595	Opiate Control (Liquichek), Level 2 MiniPak	80
12004309	Homocysteine Control (InteliQ), Level 1	13
12004310	Homocysteine Control (InteliQ), Level 2	13
12004311	Homocysteine Control (InteliQ), Level 3	13
12004312	Homocysteine Control (InteliQ), Trilevel MiniPak	13
12005786	Opiate Control (Liquichek), Level 1	80
12005787	Opiate Control (Liquichek), Level 2	80
12006239	Cardiac Troponins Control (Liquichek), Level 2	36
12006327	AMH Control (Liquichek), Trilevel	63
12006369	Cardiac Troponins Control (Liquichek), Level 3	36
12006370	Cardiac Troponins Control (Liquichek), Level 2 MiniPak	36
12006371	Cardiac Troponins Control (Liquichek), Level 3 MiniPak	36
12007028	qUAntify Advance, Bilevel	87
12007029	qUAntify Advance, Bilevel MiniPak	87
12008256	Assayed Multiqual Control (InteliQ), Level 1	11
12008257	Assayed Multiqual Control (InteliQ), Level 2	11
12008258	Assayed Multiqual Control (InteliQ), Level 3	11
12008259	Assayed Multiqual Control (InteliQ), Trilevel MiniPak	11
12008276	Specialty Immunoassay Control (InteliQ), Level LTA	14
12008277	Specialty Immunoassay Control (InteliQ), Level 1	14
12008278	Specialty Immunoassay Control (InteliQ), Level 2	14
12008279	Specialty Immunoassay Control (InteliQ), Level 3	14
12008280	Specialty Immunoassay Control (InteliQ), Four Level MiniPak	14
12008289	Tumor Marker Control (InteliQ), Level 1	14
12008290	Tumor Marker Control (InteliQ), Level 2	14
12008291	Tumor Marker Control (InteliQ), Level 3	14
12008292	Tumor Marker Control (InteliQ), Trilevel MiniPak	14
12008299	Ethanol/Ammonial Control (InteliQ), Level 1	12
12008300	Ethanol/Ammonial Control (InteliQ), Level 2	12
12008301	Ethanol/Ammonial Control (InteliQ), Level 3	12
12008302	Ethanol/Ammonial Control (InteliQ), Trilevel MiniPak	12
12008309	Diabetes Control (InteliQ), Level 1	12
12008310	Diabetes Control (InteliQ), Level 2	12
12008311	Diabetes Control (InteliQ), Level 3	12
12008312	Diabetes Control (InteliQ), Trilevel MiniPak	12
12009941	Immunology Control (InteliQ), Level 1	13
12009942	Immunology Control (InteliQ), Level 2	13
12009943	Immunology Control (InteliQ), Level 3	13
12009944	Immunology Control (InteliQ), Trilevel MiniPak	13
12009948	Immunoassay Plus Control (InteliQ), Level 1	13
12009949	Immunoassay Plus Control (InteliQ), Level 2	13
12009950	Immunoassay Plus Control (InteliQ), Level 3	13
12009941	Immunoassay Plus Control (InteliQ), Trilevel MiniPak	13
12009955	Cardiac Markers Plus Control LT, (InteliQ), Level 1	12
12009956	Cardiac Markers Plus Control LT, (InteliQ), Level 2	12
12009957	Cardiac Markers Plus Control LT, (InteliQ), Level 3	12
12009958	Cardiac Markers Plus Control LT, (InteliQ), Level 1A	12
12009959	Cardiac Markers Plus Control LT, (InteliQ), Level 1B	12
12009960	Cardiac Markers Plus Control LT, (InteliQ), Level 1C	12
12009961	Cardiac Markers Plus Control LT, (InteliQ), Level 1 MiniPak	12
12009962	Cardiac Markers Plus Control LT, (InteliQ), Level 2 MiniPak	12
12009963	Cardiac Markers Plus Control LT, (InteliQ), Level 3 MiniPak	12
12009964	Cardiac Markers Plus Control LT, (InteliQ), Level 1A MiniPak	12
12009965	Cardiac Markers Plus Control LT, (InteliQ), Level 1B MiniPak	12
12009966	Cardiac Markers Plus Control LT, (InteliQ), Level 1C MiniPak	12
12009995	Urine Chemistry Control (InteliQ), Level 1	14
12009996	Urine Chemistry Control (InteliQ), Level 2	14
12009997	Urine Chemistry Control (InteliQ), Bilevel MiniPak	14
12011308	Diabetes Control (Liquichek), Level 1 (2 mL)	45
12011309	Diabetes Control (Liquichek), Level 2 (2 mL)	45
12011310	Diabetes Control (Liquichek), Level 3 (2 mL)	45
12011331	Diabetes Control (Liquichek), Trilevel MiniPak (2 mL)	45
12011334	Diabetes Control (Liquichek), Trilevel MiniPak (1 mL)	45
12011343	Diabetes Control (Liquichek), Level 1 (1 mL)	45
12011344	Diabetes Control (Liquichek), Level 2 (1 mL)	45
12011345	Diabetes Control (Liquichek), Level 3 (1 mL)	45
12012693	Serum Indices (Liquichek), Hemolysis	43
12012694	Serum Indices (Liquichek), Icterus	43
12012695	Serum Indices (Liquichek), Lipemia	43
12012696	Serum Indices (Liquichek), Non-interfered	43
12012697	Serum Indices (Liquichek), Hemolysis MiniPak	43
12012698	Serum Indices (Liquichek), Icterus MiniPak	43
12012699	Serum Indices (Liquichek), Lipemia MiniPak	43
12012700	Serum Indices (Liquichek), Non-interfered MiniPak	43
12015644	Specialty Immunoassay Plus Control (Liquichek), Level Low	62
12015645	Specialty Immunoassay Plus Control (Liquichek), Level 1	62
12015646	Specialty Immunoassay Plus Control (Liquichek), Level 2	62
12015647	Specialty Immunoassay Plus Control (Liquichek), Level 3	62
12015648	Specialty Immunoassay Plus Control (Liquichek), Level 1-3 MiniPak	62
12015696	Specialty Immunoassay Program (EQAS)	50
12015705	Specialty Immunoassay Plus Control (InteliQ), Level Low	14
12015706	Specialty Immunoassay Plus Control (InteliQ), Level 1	14
12015707	Specialty Immunoassay Plus Control (InteliQ), Level 2	14
12015708	Specialty Immunoassay Plus Control (InteliQ), Level 3	14
12015709	Specialty Immunoassay Plus Control (InteliQ), Level 1-3 MiniPak	14
12019558	Specialty Immunoassay Plus Control (InteliQ), Level Low MiniPak	14
12019576	Specialty Immunoassay Plus Control (Liquichek), Level Low MiniPak	62
B		
BC10	TDM Program (EQAS)	50
BC23	Serum Proteins Program (EQAS)	50
BC31	Blood Gas Program (EQAS)	47
BC34	Coagulation Program (EQAS)	47
BC35	Ethanol/Ammonia Program (EQAS)	48
BC39	Cardiac Markers Program (EQAS)	47
BC45	Urine Chemistry Program (EQAS)	51
BC47	Lipids Program (EQAS)	49
BC50	Clinical Chemistry (Monthly) Program (EQAS)	48
BC75	Immunoassay (Monthly) Program (EQAS)	49
BC80	Hemoglobin Program (EQAS)	44
BC90A	Hematology Program (EQAS), Shipment A	48
BC90B	Hematology Program (EQAS), Shipment B	48
BC90C	Hematology Program (EQAS), Shipment C	48
BC90D	Hematology Program (EQAS), Shipment D	48
C		
C-310-5	Assayed Chemistry Control (Lyphochek), Level 1	38
C-315-5	Assayed Chemistry Control (Lyphochek), Level 2	38
C-320-10	Unassayed Chemistry Control (Human) (Lyphochek), Level 1 (10 mL)	38
C-325-10	Unassayed Chemistry Control (Human) (Lyphochek), Level 2 (10 mL)	38

Exact Diagnostics

ADVH102	Adenovirus High Run Control (Exact Diagnostics)	16	HCV200	HCV Verification Panel (Exact Diagnostics)	22
ADVL101	Adenovirus Low Run Control (Exact Diagnostics)	16	HCVSCREEN	HCV RNA Screen Control (Exact Diagnostics)	29
ADVP100	Adenovirus Verification Panel (Exact Diagnostics)	16	HIVAMR	HIV-1 Analytical Measurement Range Panel (Exact Diagnostics)	24
AMRP100	Analytical Measurement Range Bundle (Exact Diagnostics)	24	HIVB200	HIV-1 Verification Bundle (Exact Diagnostics)	23
ANARC	Anaplasma Run Control (Exact Diagnostics)	32	HIVH102	HIV-1 High Run Control (Exact Diagnostics)	23
B19H102	Parvovirus B19 High Run Control (Exact Diagnostics)	17	HIVL101	HIV-1 Low Run Control (Exact Diagnostics)	23
B19L101	Parvovirus B19 Low Run Control (Exact Diagnostics)	17	HIVP100	HIV-1 Verification Panel (Exact Diagnostics)	23
B19P100	Parvovirus B19 Verification Panel (Exact Diagnostics)	17	HIVP200	HIV-1 Verification Panel (Exact Diagnostics)	23
BABRC	Babesia Run Control (Exact Diagnostics)	32	HIVQP200	Verification Panel (Exact Diagnostics)	23
BKVB201	BKV Verification Bundle (Exact Diagnostics)	16	HIVQB200	Verification Bundle (Exact Diagnostics)	23
BKVH102	BKV High Run Control (Exact Diagnostics)	16	HIVSCREEN	HIV-1 RNA Screen Control (Exact Diagnostics)	29
BKVL101	BKV Low Run Control (Exact Diagnostics)	16	HPV16C	HPV Genotype 16 (Exact Diagnostics)	20
BKVP100	BKV Verification Panel (Exact Diagnostics)	16	HPV18C	HPV Genotype 18 (Exact Diagnostics)	20
BKVP200	BKV Verification Panel (Exact Diagnostics)	16	HPV68C	HPV Genotype 68 (Exact Diagnostics)	20
BORRC	Borrelia Run Control (Exact Diagnostics)	32	HPVB200	HPV Verification Bundle (Exact Diagnostics)	20
BRTRC	Bartonella Run Control (Exact Diagnostics)	32	HPVNEG	HPV Negative Control (Exact Diagnostics)	20
CDFP200	<i>C. difficile</i> Verification Panel (Exact Diagnostics)	28	HPVP100	HPV Control Set (Exact Diagnostics)	20
CDFB200	<i>C. difficile</i> Verification Bundle (Exact Diagnostics)	28	HPVP200	HPV Verification Panel (Exact Diagnostics)	20
CMVB200	CMV Verification Bundle (Exact Diagnostics)	17	HSV1H102	HSV-1 High Run Control (Exact Diagnostics)	21
CMVH102	CMV High Run Control (Exact Diagnostics)	17	HSV1L101	HSV-1 Low Run Control (Exact Diagnostics)	21
CMVL101	CMV Low Run Control (Exact Diagnostics)	17	HSV1P100	HSV-1 Verification Panel (Exact Diagnostics)	21
CMVP100	CMV Verification Panel (Exact Diagnostics)	17	HSV2H102	HSV-2 High Run Control (Exact Diagnostics)	21
CMVP200	CMV Verification Panel (Exact Diagnostics)	17	HSV2L101	HSV-2 Low Run Control (Exact Diagnostics)	21
COV000	SARS-CoV-2 (Negative) (Exact Diagnostics)	30	HSV2P100	HSV-2 Verification Bundle (Exact Diagnostics)	21
COV00CE	SARS-CoV-2 (Negative Run Control) (Exact Diagnostics)	30	HSVB200	HSV-1 Verification Bundle (Exact Diagnostics)	21
COV019	SARS-CoV-2 (Standard) (Exact Diagnostics)	30	HV6AP100	HHV-6A Verification Panel (Exact Diagnostics)	18
COV019CE	SARS-CoV-2 (Positive Run Control) (Exact Diagnostics)	30	HV6BP100	HHV-6B Verification Panel (Exact Diagnostics)	18
COVA	SARS-CoV-2 S Gene Alpha Variant Control	31	HV6H102	HHV-6 High Run Control (Exact Diagnostics)	18
COVB	SARS-CoV-2 S Gene Beta Variant Control	31	HV6L101	HHV-6 Low Run Control (Exact Diagnostics)	18
COVDK	SARS-CoV-2 S Gene Delta-Kappa Variant Control	31	JCVH102	JCV Verification Panel (Exact Diagnostics)	19
COVE	SARS-CoV-2 S Gene Epsilon Variant Control	31	JCVL101	JCV Low Run Control (Exact Diagnostics)	19
COVG	SARS-CoV-2 S Gene Gamma Variant Control	31	JCVP100	JCV Verification Panel (Exact Diagnostics)	19
COVFLU	SARS-CoV-2, Flu, RSV Run Control (Positive) (Exact Diagnostics)	30	MRSAP200	MRSA Verification Panel (Exact Diagnostics)	28
COVFLUNEG	SARS-CoV-2, Flu, RSV Run Control (Negative) (Exact Diagnostics)	30	MRSAB200	MRSA Verification Bundle (Exact Diagnostics)	28
CTNGPOS	CT/NG Positive Run Control (Exact Diagnostics)	19	NEG000	Negative Run Control (Exact Diagnostics)	26
CTNGP200	CT/NG Verification Panel (Exact Diagnostics)	19	NEG050	Negative Control (Bulk) (Exact Diagnostics)	26
CTNGB200	CT/NG Verification Bundle (Exact Diagnostics)	19	NEGHA1	HAI Negative Run Control (Exact Diagnostics)	25
EBVB200	EBV Verification Bundle (Exact Diagnostics)	18	NEGSTI	STI Negative Run Control (Exact Diagnostics)	25
EBVH102	EBV High Run Control (Exact Diagnostics)	18	NEGSYN	Synthetic Negative Run Control (Exact Diagnostics)	25
EBVL101	EBV Low Run Control (Exact Diagnostics)	18	RPNEG	RP Negative Run Control (Exact Diagnostics)	31
EBVP100	EBV Verification Panel (Exact Diagnostics)	18	RPPOS	RP Positive Run Control (Exact Diagnostics)	31
EBVP200	EBV Verification Panel (Exact Diagnostics)	18	TVMGP200	TV/MG Verification Panel (Exact Diagnostics)	20
ERLRC	Ehrlichia Run Control (Exact Diagnostics)	33	TVMGB200	TV/MG Verification Bundle (Exact Diagnostics)	20
EVRC	Enterovirus Run Control (Exact Diagnostics)	27	VIRSCREEN	HBV, HCV and HIV-1 Screen Control (Exact Diagnostics)	29
HBVAMR	HBV Analytical Measurement Range Panel (Exact Diagnostics)	24	VZVH102	VZV High Run Control (Exact Diagnostics)	27
HBVB200	HBV Verification Bundle (Exact Diagnostics)	22	VZVL101	VZV Low Run Control (Exact Diagnostics)	27
HBVH102	HBV High Run Control (Exact Diagnostics)	22	VZVP100	VZV Verification Panel (Exact Diagnostics)	27
HBVL101	HBV Low Run Control (Exact Diagnostics)	22	ZKVH102	ZKV High Run Control (Exact Diagnostics)	33
HBVP100	HBV Verification Panel (Exact Diagnostics)	22	ZKVL101	ZKV Low Run Control (Exact Diagnostics)	33
HBVP200	HBV Verification Panel (Exact Diagnostics)	22	ZKVP100	ZKV Verification Panel (Exact Diagnostics)	33
HBVSCREEN	HBV DNA Screen Control (Exact Diagnostics)	29	12019432	Bulk CSF Negative Control (Exact Diagnostics)	26
HCVAMR	HCV Analytical Measurement Range Panel (Exact Diagnostics)	24	12019386	Bulk Urine Negative Control (Exact Diagnostics)	26
HCVB200	HCV Verification Bundle (Exact Diagnostics)	22	12020068	GBS Negative Run Control (Exact Diagnostics)	27
HCVGTP	HCV Genotype Panel (Exact Diagnostics)	22	12020069	GBS Positive Run Control (Exact Diagnostics)	27
HCVH102	HCV High Run Control (Exact Diagnostics)	22	12020070	GBS Verification & Validation Panel (Exact Diagnostics)	27
HCVL101	HCV Low Run Control (Exact Diagnostics)	22	17009668	GBS Verification Bundle (Exact Diagnostics)	27
HCV200	HCV Verification Panel (Exact Diagnostics)	22			

To Order Within the United States

By Phone

1-800-224-6723
Monday – Friday
7:00 AM to 2:30 PM
Pacific Standard Time (PST)

By Fax

1-510-741-6373
or 1-800-883-1887

By Mail

Bio-Rad Laboratories, Inc.
Clinical Diagnostics Group
Customer Service
4000 Alfred Nobel Drive
Hercules, CA 94547

By Email

ask_customer_care@bio-rad.com

Be Sure Your Order Includes

Purchase order number, “ship to” address and “bill to” address, catalog number, product name, quantity required (unit and size), desired lot number and the required delivery date. Desired lot number requests can be shipped up to 30 days prior to the lot expiration. To avoid duplicate shipments, confirming orders must be clearly marked “Confirming Order/Do Not Ship.”

Standing Orders

As a convenience to our customers, we can set up standing orders for a specific lot of product to be shipped on scheduled dates. Shipment dates with respective quantities must be specified at the time of order. Restocking charges will apply to canceled standing orders.

Terms of Sale

Our terms are net due 30 days from the date of invoice. Prices are FOB shipping point. For a complete listing of terms and conditions regarding quantity, pricing, applicable taxes, delivery, discrepancies, returned goods and cancellation policies, please contact our Customer Service Department.

**For Worldwide Customers**

For service outside the United States, please refer to the inside back cover for the Bio-Rad location nearest you.

Notice of Rights

Copyright © 2022 Bio-Rad Laboratories, Inc. All rights reserved. Printed in the United States of America. No part of this book may be reproduced, stored or transmitted in any form by any means, electronic, mechanical, optical, photocopying, recording or otherwise, without prior permission of the publisher.

Timeliness, Errors and Omissions

The information provided in this catalog carries no express or implied warranties as to accuracy, timeliness, or appropriateness for a particular purpose. Bio-Rad Laboratories disclaims responsibility for content errors and omissions, for any allegedly infringing material, and for any damages whatsoever associated with user reliance on information provided in this catalog.

Applicable Laws

This was published by Bio-Rad Laboratories from its offices in Irvine, California. Bio-Rad makes no representation that the information and products in this catalog are appropriate or available for use in your country, and access to such information and products from territories where such information and products are illegal is prohibited. If you choose to use this catalog, you do so on your own initiative and at your own risk, and are responsible for compliance with applicable local laws. Contact your local Bio-Rad office if you have questions regarding the availability of products listed in this catalog. These terms will be governed by and construed in accordance with the laws of the State of California, without giving effect to any principles of conflicts of laws.

Trademarks

Bio-Rad Laboratories

The following are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.

Amplichek
Bio-Rad
EQAS
IntelliQ

Liquichek
Lyphocheck
Mission: Control
QCNet

qUAntify
Unity
Unity Real Time
UnityWeb

VARIANT
VIROCLEAR
VIROTROL

All trademarks used herein are the property of their respective owner.

Worldwide Contact Information

www.bio-rad.com/qualitycontrol



UNITED STATES OF AMERICA

Bio-Rad Laboratories, Inc.
Clinical Diagnostics
4000 Alfred Nobel Drive
Hercules, CA 94547
Phone: (510) 724-7000
Toll Free (US): 1 (800) 2 BIORAD (1-800-224-6723)
Fax: (510) 741-5800/1 (800) 879-2289 (US)
www.bio-rad.com/qualitycontrol



AUSTRALIA

Bio-Rad Laboratories Pty. Ltd.
South Granville, New South Wales
Phone: +61 (2) 9914 2800
Toll Free: 1800 224 354 (within Australia only)
Fax: +61 (2) 9914 2888
Email: Sales.Australia@bio-rad.com



AUSTRIA

Bio-Rad Laboratories Ges.m.b.H.
Wien (Vienna)
Phone: +43 (0) 1 877 89 01 9
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +43 (0) 1 876 56 29
Email: Contact_CentralEurope@bio-rad.com



BELGIUM

Bio-Rad Laboratories N.V.
Temse
Phone: +32 (0) 3 710 53 00
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +32 (0) 3 710 53 01
Email: Benelux_Contact@bio-rad.com



BRAZIL

Bio-Rad Laboratórios Brasil Ltda
Santo Amaro, São Paulo
Phone: +55 11 3065 7550
Email: atendimento@bio-rad.com



CANADA

Bio-Rad Laboratories, Ltd.
Montréal, Québec
Phone: (514) 334 4372
Freephone: 1 (800) 361 1808
Fax: (514) 334 0872
Email: cdg_canada_salesmarketing@bio-rad.com



CHINA

Bio-Rad Laboratories (Shanghai) Co. Ltd.
Shanghai
Phone: +86 21 6169 8500
Fax: +86 21 6169 8599
Email: Sales.China@bio-rad.com



CZECH REPUBLIC

Bio-Rad spol, s.r.o.
Prague
Phone: +36 1 459 6192
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +420 241 431 6602
Email: Logistika_CZ@bio-rad.com



DENMARK

Bio-Rad Laboratories
Copenhagen
Phone: +46 844 98051 / +45 44 52 10 00
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +45 44 52 10 01
Email: denmark@bio-rad.com



FINLAND

Bio-Rad Finland Oy
Helsinki
Phone: 00 800 00 246 723 / 09 804 22 00
Fax: +358 (0)9 7597 5010
Email: finland@bio-rad.com



FRANCE

Bio-Rad
Marnes-la-Coquette
Phone: +33 (0) 1 47 95 93 00
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +33 (0) 1 47 95 61 45
Email: Contact_France@bio-rad.com



GERMANY

Bio-Rad Laboratories GmbH
Feldkirchen
Phone: +49 89 3188 4393
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +49 (0)89 31884 100
Contact_CentralEurope@bio-rad.com



GREECE

Bio-Rad Laboratories M E.P.E.
Athens
Phone: +30 210 7774396
Fax: +30 210 7774376



HONG KONG

Bio-Rad Pacific Ltd.
Quarry Bay
Phone: +85 2 2789 3300
Fax: +85 2 2789 1257
Email: Sales.HongKong@bio-rad.com



HUNGARY

Bio-Rad Hungary Ltd.
Budapest
Phone: +36 1 459 6190
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +36 1 459 6101
Email: Info_Hungary@bio-rad.com



INDIA

Bio-Rad Laboratories (India) Pvt. Ltd.
Gurugram, Haryana
Phone: 91 124 4029300
Fax: +91 124 2398115
Email: sales.india@bio-rad.com



ISRAEL

Bio-Rad Laboratories Ltd.
Rishon Le Zion
Phone: +972 3 963 6050
Fax: +972 3 951 4129
Email: isreal_cdg@bio-rad.com



ITALY

Bio-Rad Laboratories S.r.l.
Milan
Phone: +39 02 216091
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +39 02 21609553
Email: contact_italy@bio-rad.com



JAPAN

Bio-Rad Laboratories K.K.
Tokyo
Phone: +81 3 6361 7070
Fax: +81 3 5463 8481
Email: diag-jp@bio-rad.com



KOREA

Bio-Rad Korea Ltd.
Seoul
Phone: +82 080 007 73730
Fax: +82 (2) 3472 7003
Email: Sales.Korea@bio-rad.com



MEXICO

Bio-Rad, S.A.
Mexico City
Phone: +52 (55) 5488-7670
Fax: +52 (55) 1107-7246
Email: cdg_mexico@bio-rad.com



THE NETHERLANDS

Bio-Rad Laboratories B.V.
Veenendaal
Phone: +31 (0) 318 540 666
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +31 (0)318 542216
Email: Benelux_Contact@bio-rad.com



NEW ZEALAND

Bio-Rad New Zealand
Auckland
Phone: +64 (9) 415 2280
Toll Free: 0508 805 500 (within New Zealand only)
Fax: +64 (9) 415 2284
Email: Sales.NZ@bio-rad.com



NORWAY

Bio-Rad Laboratories
Oslo
Phone: +46 844 98052 / +47 23 38 41 30
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +47 23 38 41 39
Email: norway@bio-rad.com



POLAND

Bio-Rad Polska Sp. z o.o.
Warsaw
Phone: +48 22 331 99 81
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +48 22 331 99 88
Email: Logistyka_PL@bio-rad.com



PORTUGAL

Bio-Rad Laboratories, Lda.
Amadora
Phone: +351 214 727 700
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +351 214 727 7777
Email: customerservice_portugal@bio-rad.com



RUSSIA

Bio-Rad Laboratorii
Moscow
Phone: +7 495 721 1404
Fax: +7 495 721 1412
Email: info_russia@bio-rad.com



SINGAPORE

Bio-Rad Laboratories (Singapore) Pte. Ltd.
Phone: +65 6415 3170
Fax: +65 6415 3189
Email: sales.singapore@bio-rad.com



SOUTH AFRICA

Bio-Rad Laboratories (Pty) Ltd.
Johannesburg
Phone: +27 11 442 8508 / +27 10 055 8040
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +27 11 442 8525
Email: CDG_TechSupport_EEMEA@bio-rad.com



SPAIN

Bio-Rad Laboratories, S.A.
Madrid
Phone: + 34 91 49 06 580
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +34 91 590 52 11
Email: customerservice_spain@bio-rad.com



SWEDEN

Bio-Rad Laboratories A.B.
Solna
Phone: +46 844 98053
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +46 8 55 51 27 80
Email: sweden@bio-rad.com



SWITZERLAND

Bio-Rad Laboratories AG
Cressier
Phone: +41 (0) 61 717 9555
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +41-(0) 26 674 52 19
Email: Contact_CentralEurope@bio-rad.com



TAIWAN

Bio-Rad Laboratories Taiwan Ltd.
Taipei
Phone: +886 (2) 2578 7189
Fax: +886 (2) 2578 6890



THAILAND

Bio-Rad Laboratories Ltd.
Bangkok
Phone: 662 651 8311
Fax: 662 651 8312
Email: Sales.Thailand@bio-rad.com



UNITED KINGDOM

Bio-Rad Laboratories Ltd.
Watford, Hertfordshire
Phone: +44 01 923 47 1301
Freephone: 00 800 00 BIORAD (00 800 00 24 67 23)
Fax: +44 01 923 47 1340
Email: UK_contact@bio-rad.com

EQAS

An independent, external assessment of laboratory performance in comparison to your peers.

QCNet.com/eqas

Independent QC

Ongoing, proactive, unbiased daily QC that helps identify errors as they occur or begin to trend.

QCNet.com/quality-controls

Unity™

QC Data Management tools that help you create a strategy to reduce risk and streamline QC workflow.

QCNet.com/qc-data-management

Want to know more? Visit us online at www.bio-rad.com/qualitycontrol



Connect with us

@BioRadQC | Twitter

LinkedIn/company/Bio-Rad | LinkedIn

BioRadQC | YouTube

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Clinical
Diagnostics Group

Website www.bio-rad.com/diagnostics **U.S.** 1 800 224 6723 **Australia** +61 (2) 9914 2800 **Austria** +43 (0) 1 877 89 01 9 **Belgium** +32 (0) 3 710 53 00
Brazil +55 11 3065 7550 **Canada** +1 514 334 4372 **China** +86 21 6169 8500 **Czech Republic** +420 241 431 660 **Denmark** +45 44 52 10 00
Finland +358 9 804 22 00 **France** +33 (0) 1 47 95 60 00 **Germany** +49 (0) 89 31884 393 **Greece** +30 210 7774396 **Hong Kong** +85 2 2789 3300
Hungary +36 1 459 6190 **India** +91 124 4029300 **Israel** +972 03 963 6025 **Italy** +39 024 94 86 600 **Japan** +81 3 6361 7070 **Mexico** +52 (55) 5488 7670
The Netherlands +31 (0) 318 540 666 **New Zealand** +64 (9)415 2280 **Norway** +47 23 38 41 30 **Poland** +48 22 331 99 99 **Portugal** +351 21 47 27 700
Republic of Korea +82 080 007 7373 **Russia** +7 (495) 721-14-04 **Singapore** +65 6415 3170 **South Africa** +27 11 442 8508 **Spain** +34 91 490 6580
Sweden +46 844 98053 **Switzerland** +41 (0) 61 717 9555 **Taiwan** +886 (2) 2578-7189 **Thailand** (662) 651 8311 **United Kingdom** +44 (0) 1923 471301

