

Infectious Disease Controls for Serology Testing

A comprehensive guide for monitoring testing for Hepatitis, Retrovirus, Respiratory, Sexually Transmitted and Congenital Diseases



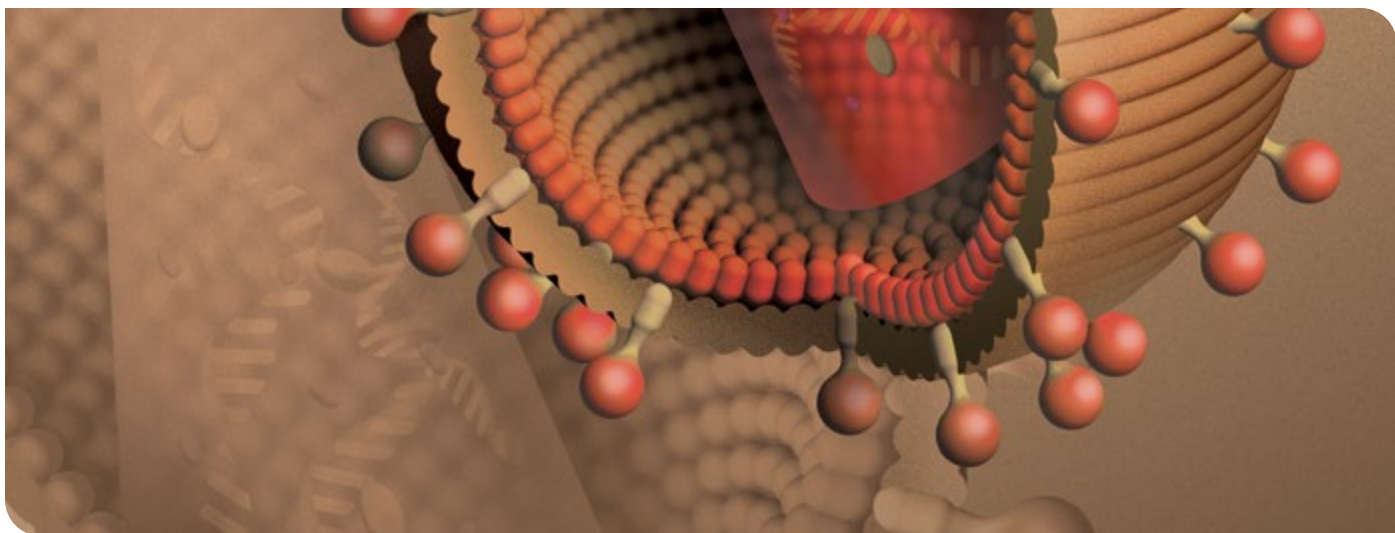


Table of Contents

For Your Consideration

The Impact of Inaccurate Results to Patients and Your Laboratory	2
Why Choose Infectious Disease Quality Controls from Bio-Rad?	3
Assuring Quality in the Laboratory	4
Effective Document Control with Unity QC Data Management Systems	4
Important Differences Between Independent and In-kit Controls	5

Reference

List of Popular Instruments, Assays and Bio-Rad Controls	6
--	---

Quality Control Products and Services

Hepatitis and Retrovirus Controls	7
Congenital/Pediatric Disease Controls	10
Respiratory Disease Controls	11
Sexually Transmitted Disease Controls	12
Specialty Infectious Disease Controls	13
QC Data Management Solutions	14



The Impact of Inaccurate Results to Patients and Your Laboratory

Each year, approximately 16 to 22 million people in the United States are tested for HIV. The Centers for Disease Control and Prevention (CDC) reports an estimated 39,000 new HIV infections annually. Around the world, the impact of all infectious diseases is even greater. The role of the laboratory in providing accurate test results is a key component in controlling the spread of infection.

“Results of control materials must meet the laboratory’s and, as applicable, the manufacturer’s test system criteria for acceptability before reporting patient test results.”

CLIA 493.1256 Standard: Control procedures (f)

The implications of reporting a false positive result are obvious, but the consequences of reporting a false negative can be just as devastating in terms of necessary treatment and personal life style changes for the patient. Using Bio-Rad infectious disease controls as an independent assessment of your testing systems helps to provide additional confidence in your test results.

Source: Centers for Disease Control and Prevention



Why Choose Infectious Disease Quality Controls from Bio-Rad?

We Speak From More Than 30 Years of Experience in Clinical Diagnostics

Bio-Rad Laboratories offers the most comprehensive line of infectious disease quality controls. Choose from over 25 infectious disease control products that cover a wide-range of testing, including hepatitis and retrovirus, sexually transmitted diseases, congenital diseases, panels and other specialty testing. Our infectious disease controls are used on a wide-range of test systems and methods, which include automated immunoassay systems, EIA processors and manual EIA methods.

Where Can Our Controls Be Used?

- Blood Banks
- Public Health Laboratories
- Virology Laboratories
- Reference Laboratories
- Microbiology Departments
- Serology Departments
- Immunology Departments
- Special Chemistry

How Do Bio-Rad Infectious Disease Quality Control Products and Services Stand Out From the Competition?

- Unbiased independent assessment of test systems
- Worldwide peer group comparisons available
- Liquid stable, ready-to-use, multi-analyte
- Basic to advanced QC data management solutions
- Assayed virology products available outside the US only

Assuring Quality in the Laboratory

Regulations often provide plenty of general information on quality control, but may not always offer specific information regarding quality control especially for infectious disease testing.

Without this specific direction, laboratories may unknowingly perform the minimum amount of quality control, or use only the manufacturer provided (in-kit) controls.

A responsible laboratory understands the critical nature of infectious disease testing. Establishing an internal QC system that incorporates independent controls and comprehensive data management will help you assure the quality and reliability of your infectious disease testing.

“The laboratory shall design quality control procedures that verify the attainment of the intended quality of results.”

ISO 15189:2012 5.6.2.1

Effective Document Control With Unity QC Data Management Systems

Laboratory inspectors often need to see documentation of QC data as evidence of proper quality control procedures. Yet documentation and monitoring of quality controls can become a huge burden on a busy laboratory. Bio-Rad provides QC data management solutions that allow you to optimize your laboratory performance. Choose from basic to advanced statistical rules – available as a web-based or desktop software solution.

Unity Interlaboratory Program

Participation in an interlaboratory program is extremely important for laboratories testing for hepatitis, retrovirus, STDs and other infectious diseases. Join the Unity Interlaboratory Program – the world’s largest community of quality control users.

Compare to peer group data to increase confidence in your patient test results.

▼ Example Unity Report

Unity January 2020

Worldwide Report Conventional Units

VIROTROL III • Lot 23456 • Exp 01-Jan-2022

Anti-HAV, IgM Chemiluminescence Sample/Cutoff (Index, ISR)

Level	Mon	Cum
Mean	24.8	2.42
SD	0.181	0.230
CV	7.3	9.51
# Points	792	5490
# Labs	24	35

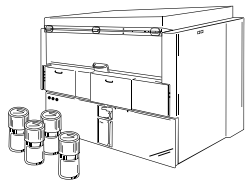
Siemens Diagnostics Centaur/XP Series
Dedicated Reagent

Anti-HBc, IgM Chemiluminescence Sample/Cutoff (Index, ISR)

Level	Mon	Cum
Mean	1.96	1.98
SD	0.134	0.214
CV	6.8	10.8
# Points	642	4885
# Labs	18	30

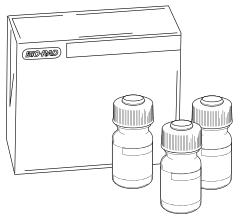
Siemens Diagnostics Centaur/XP Series
Dedicated Reagent

Important Differences Between Independent and In-kit Controls



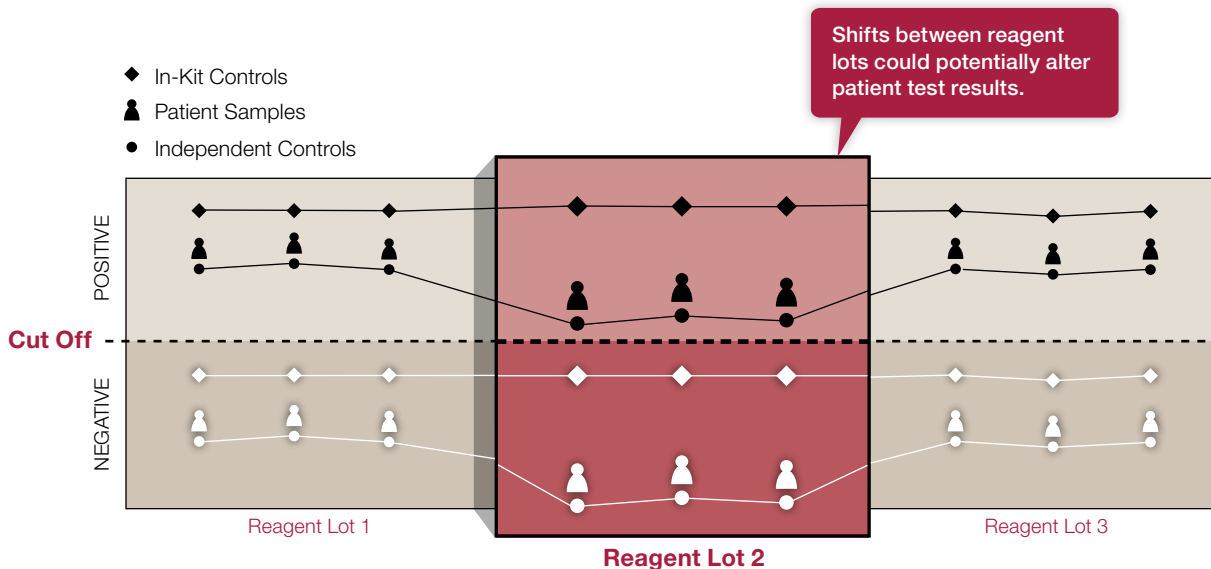
In-kit Controls

- May be same material as calibrators
- Specific to each kit/reagent lot
- May not identify shifts between reagent lots
- Can be used to calculate assay cut-off
- Levels typically well above or below assay cut-off
- Frequent lot changes don't allow for long-term monitoring
- May not be similar to patient sample(s)



Independent Controls

- Different from kit calibrators
- Independent of kit/reagent lot
- Identify shifts between reagent lots
- Designed to monitor the precision of the test system
- Levels are typically targeted to be reactive within a relevant range
- Same control lot can be used across multiple reagent lots allowing long-term monitoring of assay performance
- Control may be treated the same as a patient sample



Controls should always be run with new reagent lots. However, it's important to be aware that in-kit controls are typically used as part of the process to determine assay cut-off, and are specific to each new lot of reagents. Independent controls are independent of reagent lots, and therefore, may detect shifts or changes that may not be identified by the in-kit controls.

In the illustration above, the in-kit controls matched to Reagent Lot 2, did not identify a change in sensitivity that resulted in a shift in patient sample results and the independent controls.

Without independent controls, a situation like this could potentially result in reporting incorrect results.

List of Popular Instruments, Assays, Methods and Bio-Rad Controls

Please contact your Bio-Rad representative for additional details.

Manufacturer/Instrument/Assay/Method	▼ Positive Controls														▼ Negative Controls									
	VIROTROL I	VIROTROL II	VIROTROL III	VIROTROL IV	VIROTROL HIV-1 Ag	VIROTROL HIV-1 gO	VIROTROL HIV-2	VIROTROL HAV-IgM	VIROTROL HBeAg	VIROTROL HBc-IgM	VIROTROL PLUS-R	VIROTROL ToRCH	VIROTROL ToRCH-M	Liquichek ToRCH Plus (Positive)	VIROTROL Chagas	VIROTROL MuMZ	VIROTROL EBV	PyloriTrol	VIROTROL Lyme	VIROCLEAR	VIROCLEAR ToRCH	Liquichek ToRCH Plus (Negative)	VIROCLEAR MuMZ	VIROCLEAR EBV
Abbott Diagnostics																								
Alinity	●	●	●	●	●		●		●	●		●	●	●			●			●	●			●
Architect Series	●	●	●	●	●		●	●	●	●		●	●	●	●					●	●	●		
ARCHITECT i System																								
PRISM	●				●		●								●					●				
Beckman Coulter																								
Access Series/LXi/UniCel DxI	●	●	●				●				●	●	●							●	●	●		
BioMérieux																								
VIDAS	●	●	●	●	●		●	●	●	●		●	●	●		●		●		●	●	●	●	●
Bio-Rad Laboratories																								
BioPlex 2200	●				●		●					●	●			●	●			●	●		●	●
Evolis/EIA (GS, Genscreen, Monolisa, Platelia)	●	●	●	●	●		●	●	●	●		●		●		●	●	●	●	●	●	●	●	●
Geenius HIV 1/2	●						●													●				
Platelia SARS Screen																								
Cellex																								
Cellex Rapid Test																								
DiaSorin																								
LIAISON/ LIAISONXL	●	●	●	●	●		●	●	●	●		●	●	●		●	●		●	●	●		●	●
ETI-MAX Systems	●	●	●	●				●	●	●		●		●		●	●	●		●	●		●	●
Diesse																								
Chorus Line												●	●								●	●		
Ortho Clinical Diagnostics																								
VITROS	●	●	●	●		●	●	●	●	●		●	●	●						●	●	●		
VITROS ECi/ECiQ/3600/5600/XT/7600																								
Roche Diagnostics																								
cobas 4000/6000/8000 Series		●	●	●	●		●	●	●	●	●	●	●	●						●	●	●		
Cobas e411/e601/e602/e801																								
Elecsys Systems		●						●	●	●	●	●	●	●						●	●	●		
Siemens																								
ADVIA Centaur Series	●	●	●				●	●		●		●	●	●						●	●			
Atellica	●	●	●		●		●		●				●	●						●	●			
IMMULITE Series	●	●	●				●					●	●	●						●	●	●		

The controls listed here include analyte(s) that may be appropriate for use on the test systems noted.

Content is provided as a guide and for informational purposes only. Results and specific performance characteristics may vary with different manufacturers' test kits and assay procedures, different lots of the same test kit, and different lots of the control material. Please note that not all products may be available in all countries. Refer to the product insert for current product claims.

Hepatitis & 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

Analytes

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

Analytes

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

Analytes

Anti-HAV
Anti-HBs

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.

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

Analytes

Anti-HAV
HAV-IgM
HBc-IgM

Hepatitis & Retrovirus

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

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

Analyte

HIV-1 Ag

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

Analytes

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

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

Analyte

Anti-HIV-2

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

Analyte

Anti-HIV-1 group O

Refer to myInserTs.com or the package insert of currently available lots for specific analyte and stability claims.

Hepatitis & 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

Analytes

Non-reactive for:

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

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

Analyte

HAV-IgM

VIROTROL HBc-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

Analyte

HBc-IgM

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

Analyte

HBeAg

Congenital/Pediatric Disease

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

Analytes

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

VIROTROL ToRCH-M



- An unassayed quality control
- Human serum based
- 2 year shelf life at 2–8°C (Cat #00117A) or 18 month shelf life at 2–8°C (Cat #00117B)
- 60 day open-vial stability at 2–8°C

Analytes

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

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

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

Analytes

Non-reactive for:

Toxoplasma gondii IgG
Toxoplasma gondii IgM
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

Liquichek ToRCH Plus Control



- An unassayed, 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

Analytes

Helicobacter pylori IgG
Toxoplasma gondii IgG
Treponema pallidum (Syphilis) IgG
Cytomegalovirus (CMV) IgG
Epstein-Barr Virus (EBV) IgG
Herpes Simplex Virus Type 1 (HSV-1) IgG
Herpes Simplex Virus Type 2 (HSV-2) IgG
Rubella Virus IgG
Syphilis RPR
Varicella Zoster Virus (VZV) IgG

Refer to myInserTs.com or the package insert of currently available lots for specific analyte and stability claims.

Congenital/Pediatric Disease

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

Analytes

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

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

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

Analytes

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

PneumoTrol



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

Analytes

Chlamydia trachomatis IgG
Mycoplasma pneumoniae IgG
Mycoplasma pneumoniae IgM

Sexually Transmitted Disease

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

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 titerable member
- Human serum based
- 2 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C

Analyte

Non-treponemal Antibodies (Syphilitic Reagin)

Specialty Infectious Disease

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

Analytes

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

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

Analytes

Non-reactive for:

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

Refer to myInserTs.com or the package insert of currently available lots for specific analyte and stability claims.

Specialty Infectious Disease

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

Analyte

Helicobacter pylori IgG

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

Analyte

Trypanosoma cruzi (*T. cruzi*)

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

Analyte

Borrelia burgdorferi IgG

VIROTROL WNV



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

Analytes

West Nile Virus IgG
West Nile Virus IgM

QC Data Management Solutions

Join over 29,000 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.¹

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 50,000 instruments worldwide
- 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 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.

Choose From Four 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 Real Time online	Unity Real Time
Type of Solution				
Web Service (Web-based)	X		X	
Desktop Software		X		X
Unity Interlaboratory Reports				
Monthly Reports	X	X	X	X
InstantQC Reports	X	X	X	X
Basic Intralaboratory Charts & Reports				
Westgard Rules	X	X	X	X
Various Charts and Reports	X	X	X	X
Additional Features				
Westgard Advisor (Subscription sold separately)	X		X	X
Bench Review with Data Review Report	X	X	X	X
Supervisor Data Review with Data Review Report			X	X
Analytical Goals			X	X
Dynamic Data Set Comparisons			X	X
Database Platform				
Bio-Rad Hosted	X		X	
Installed on Local Laboratory PC		X		X
Connectivity (Sold separately)				
WebConnect	X		X	
UnityConnect	X	X	X	X

Want to know more? Discover the power of Unity Solutions at www.qcnet.com/datamanagement/

¹Malone B. A New Approach to Quality Control. AACC. 2011;37:11. Available at: <http://www.aacc.org/publications/cln/2011/november/Pages/ANewApproachtoQualityControl.aspx#>. Accessed June 14, 2013.



Basic



Advanced



Desktop based



Web based



Mobile

Premier Unity Software Solutions

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

UnityWeb

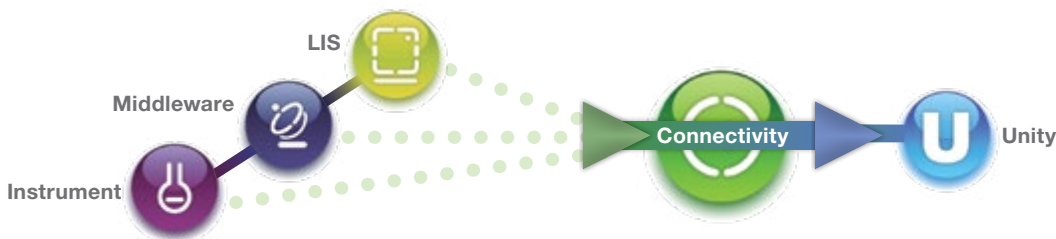


Entry-Level Online QC Data Management Solution

- 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)

Connectivity Solutions for Automated Uploads to Unity Software

Simple automated uploading of your data from your instruments, middleware and LIS to Unity software solutions



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
- Transparent to the flow of data to the LIS
- Capture and analyze QC data economically from laboratory instruments not connected to the LIS
- Easy-to-install optional interface hardware
- Choose from soft, serial and ethernet connections

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

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



Regulatory Requirements

"The laboratory shall design quality control procedures that verify the attainment of the intended quality of results."

– ISO 15189:2012(E), Subclause 5.6.2.1

Unity Interlaboratory Reports

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

Bio-Rad drives innovation in the development of Quality Control for the clinical laboratory and we're raising the bar again. It's time to take a fresh look at your Unity peer reports. Redesigned formats make reports easier than ever to read. The addition of color, graphics and statistical information allows you to interpret your data with ease, enabling prompt action to potential problems. Condensed report formats require fewer pages and view equally well printed in color or black and white.

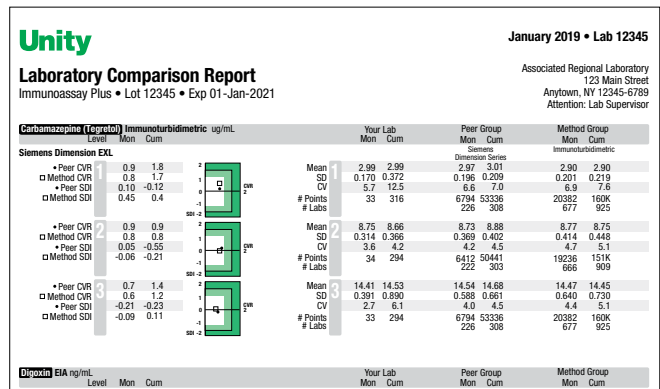
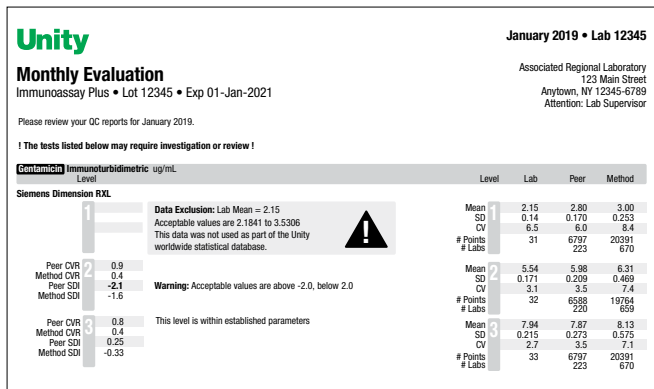
Unity Interlaboratory Program Reports Feature:

- 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
- InstantQC Report

Contact your Bio-Rad Sales Representative or visit QCNet.com for more information.



Receive on-demand InstantQC Reports via QCNet today!

InstantQC Reports allow you to compare your QC results to other laboratories at any time, with real-time peer information. On-demand reports are particularly useful for troubleshooting problems with test system performance.



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

Ordering Information

Cat #	Description
Hepatitis & Retrovirus	
VIROTROL I	
00100A / 00100B / 00100C / 00100E / 00100F / 00100G / 00100H	Single Level/Dropper Bottle1 x 5 mL
00101B / 00101C / 00101E / 00101F	Single Level/Tube10 x 4 mL
00102B / 00102C	Single Level/Dropper Bottle2 x 20 mL
Assayed VIROTROL I-C*	
00164	Single Level/Dropper Bottle1 x 5 mL
Assayed VIROTROL I-E*	
00166	Single Level/Dropper Bottle1 x 5 mL
Assayed VIROTROL I-F*	
00168	Single Level/Dropper Bottle1 x 5 mL
VIROTROL II	
00104A / 00104B	Single Level/Dropper Bottle1 x 5 mL
00115A / 00115B	Single Level/Tube10 x 4 mL
Assayed VIROTROL II-B	
00152	Single Level/Dropper Bottle1 x 5 mL
VIROTROL III	
00110A / 00110C	Single Level/Dropper Bottle1 x 5 mL
VIROTROL IV	
00111	Single Level/Dropper Bottle1 x 5 mL
VIROTROL HIV-1 Ag	
00108A / 00108B	Single Level/Dropper Bottle1 x 5 mL
VIROTROL PLUS-R	
12000538	Single Level/Tube5 x 4 mL
12000539	Single Level/Tube1 x 4 mL
VIROTROL HIV-2	
00105A / 00105B / 00105C	Single Level/Dropper Bottle1 x 5 mL
VIROTROL HIV-1 gO	
00113	Single Level/Tube5 x 4 mL
00113X	Single Level/Tube1 x 4 mL
VIROCLEAR	
00106	Single Level/Dropper Bottle1 x 5 mL
00107	Single Level/Dropper Bottle2 x 20 mL
00112	Single Level/Tube10 x 4 mL
VIROTROL HAV-IgM	
00142A	Single Level/Dropper Bottle1 x 5 mL
VIROTROL HBc-IgM	
00143	Single Level/Dropper Bottle1 x 5 mL
VIROTROL HBsAg	
00144	Single Level/Dropper Bottle1 x 5 mL
Congenital Disease	
VIROTROL ToRCH	
00109	Single Level/Dropper Bottle1 x 5 mL
VIROTROL ToRCH-M	
00117A	Single Level/Dropper Bottle1 x 5 mL
00117B	Single Level/Dropper Bottle1 x 5 mL

Cat #	Description
Congenital Disease (continued)	
VIROCLEAR ToRCH	
00118	Single Level/Dropper Bottle1 x 5 mL
Liquichek ToRCH Plus Control	
239	Positive3 x 3 mL
239X	Positive MiniPak1 x 3 mL
227	Positive, Assayed*3 x 3 mL
227X	Positive Assayed MiniPak*1 x 3 mL
228	Negative3 x 3 mL
228X	Negative MiniPak1 x 3 mL
VIROTROL MuMZ	
00119	Single Level/Dropper Bottle1 x 5 mL
VIROCLEAR MuMZ	
00133	Single Level/Dropper Bottle1 x 5 mL
Respiratory	
VIROTROL SARS-CoV-2	
200300A	Single Level/Tube1 x 4 mL
200305A	Single Level/Tube5 x 4 mL
200300B	Single Level/Tube1 x 4 mL
200305B	Single Level/Tube5 x 4 mL
200300C	Single Level/Tube1 x 4 mL
200305C	Single Level/Tube5 x 4 mL
VIROCLEAR SARS-CoV-2	
200500	Single Level/Tube1 x 4 mL
200505	Single Level/Tube5 x 4 mL
PneumoTrol	
00129	Single Level/Dropper Bottle1 x 5 mL
Sexually Transmitted Disease	
VIROTROL Syphilis LR-A	
00171	Single Level/Tube5 x 4 mL
00171X	Single Level/Tube1 x 4 mL
VIROTROL RPR Panel	
01400	Three Levels/Dropper Bottle3 x 1.5 mL
Specialty Infectious Disease	
VIROTROL EBV	
00128A	Single Level/Dropper Bottle1 x 5 mL
VIROCLEAR EBV	
00131	Single Level/Dropper Bottle1 x 5 mL
PyloriTrol	
00121	Single Level/Dropper Bottle1 x 5 mL
VIROTROL Chagas	
00125	Single Level/Tube5 x 4 mL
00125X	Single Level MiniPak1 x 4 mL
VIROTROL Lyme	
00132	Single Level/Dropper Bottle1 x 5 mL
VIROTROL WNV	
00116	Single Level/Dropper Bottle1 x 5 mL

(Continued on Back Cover)

*Available outside the United States only.
Please note that not all products may be available in all countries.

Ordering Information

(Continued from Inside Back Cover)

Cat # Description

QC Data Management Solutions

Unity Real Time

804	Installation Package	Single use
804-1	Unity Real Time	Annual subscription
825i	Unity Remote Installation & Training	Single use
858i	Additional Software Training	Single use

Unity Web

870-1	UnityWeb	Annual subscription
825i	Unity Remote Installation & Training	Single use

UnityConnect

820-1	UnityConnect Software	Annual subscription
858i	Additional Software Training	Single use

WebConnect

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

EQAS

An independent, external assessment of 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/independentqc

Unity

QC Data Management tools that help you create a strategy to reduce risk and streamline QC workflow.

QCNet.com/datamanagement



**Bio-Rad
Laboratories, Inc.**

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com/idcontrols

Clinical
Diagnostics Group

Website www.bio-rad.com/diagnostics **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 **Korea** +82 080 007 7373 **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 **Russia** +7 495 721 1404 **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

BIOPLEX, BIO-RAD, LIQUICHEK, UNITY REAL TIME, UNITYWEB, VIROCLEAR and VIROTROL are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.
All trademarks used herein are the property of their respective owner.

© 2021 Bio-Rad Laboratories, Inc. Printed in USA QSD21-0091 Q-1565 Rev. 04/2021

Printed on recycled paper with soy-based inks