



Bio-Rad Quality Control Products in Dimension Vista® Vials





Bio-Rad Controls in Dimension Vista® Vials

Our comprehensive line of liquid, multi-analyte quality controls are filled in Vista vials, ready for use on Dimension Vista® instruments. These assayed controls, with verified QC stability claims for on-board instrument storage, allow for routine quality control of Vista® assays.

Controls in Vista vials allow Dimension Vista® customers to:

- Rely on Bio-Rad controls and QC performance
- Take advantage of shared control lot numbers with existing product configurations
- Access extensive Unity™ peer group reports and comparisons
- Scan 2D barcodes to automatically set-up QC with Vista® assigned values to prevent transcription errors
- Alleviate manual pour-off tasks to minimize handling errors and reduce the cost of extra consumables
- Monitor analytes that are specific to the Dimension Vista® test menu



Siemens recommends the family of Bio-Rad Controls in Dimension Vista® vials for your Quality Control needs.

Bio-Rad Controls make using your Dimension Vista® Even More Efficient



2D barcodes are scanned from the product insert to automatically set-up QC with Dimension Vista® assigned values to prevent transcription errors.



QC vials are inserted into the vial carrier and loaded onto the instrument for automated scanning to identify the specific lot and catalog numbers.



QC vials are loaded and stored refrigerated on the Dimension Vista®, providing automated on board QC. Based on the number of analytes tested, an appropriate aliquot is taken from the vial and QC is automatically performed.



On board QC is automatically monitored and the Dimension Vista® alerts the technologist when new QC vials are required for analysis.

Unity
Worldwide Report
Multiquel 1, 2, 3 Unassayed

Abbott ARCHITECT (e, I, G)
Mean: 2.60
SD: 0.032
CV: 1.3
Points: 2346
Labs: 4859

Beckman Coulter CX Series
Mean: 2.44
SD: 0.051
CV: 2.1
Points: 88
Labs: 4

Beckman Coulter LX20, LX1725, Dx
Mean: 2.44
SD: 0.045
CV: 1.8
Points: 436
Labs: 10

Beckman Coulter UniCel DxC Series
Mean: 2.43

Unity
Laboratory Comparison Report
Immunoassay Plus • Lot 12345 • Exp 01-Jan-2016

Level	Your Lab		Peer Group		Method Group	
	Mean	CV	Mean	CV	Mean	CV
0.9	1.8	2.99	2.97	3.01	2.90	2.90
0.8	1.7	0.170	0.170	0.209	0.201	0.219
0.10	-0.12	5.7	12.5	6.6	7.0	6.9
0.45	0.4	6794	53336	226	308	20382
						1606
						677
						925

January 2014
Conventional Units

January 2014 • Lab 12345
Associated Regional Laboratory
123 Main Street
Anytown, NY 12345-6789
Attention: Lab Supervisor

Level	Your Lab		Peer Group		Method Group	
	Mean	CV	Mean	CV	Mean	CV
0.9	0.9	8.75	8.86	8.73	8.88	8.77
0.8	0.8	0.014	0.366	0.369	0.622	0.414
0.05	-0.55	3.6	4.2	4.2	4.5	4.7
-0.06	-0.21	34	294	642	50841	19236
						151K
						666
						969

The Unity™ Interlaboratory Program provides easily interpreted reports and worldwide peer group comparisons essential for more reliable patient test results.

Liquid Assayed Multiquel® Control

A trilevel control for monitoring general chemistry test procedures.

- Offers more than 60 analytes
- 5 day open-vial (punctured) stability for all analytes when stored at 2–8°C
- 3 year shelf life when stored at -20°C to -50°C

Analytes

General Chemistry

Albumin	GGT
Alkaline Phosphatase (ALP)	Glucose
ALT/SGPT	Iron
Amylase	Iron (TIBC)
AST/SGOT	Lactate (Lactic Acid)
Bilirubin (Direct)	Lactate Dehydrogenase (LDH)
Bilirubin (Total)	Lipase
Calcium (Total)	Magnesium
Chloride	Phosphorus
Cholesterol (HDL)	Potassium
Cholesterol (LDL)	Protein (Total)
Cholesterol (Total)	Pseudocholesterase
CO ₂	Sodium
Creatine Kinase	Triglycerides
Creatinine	Urea Nitrogen
Ethanol	Uric Acid

Immunology

Alpha-1-Antitrypsin
Apolipoprotein A-1
Apolipoprotein B
Ceruloplasmin
Complement C3
Complement C4
Haptoglobin
Immunoglobulin A (IgA)
Immunoglobulin G (IgG)
Immunoglobulin M (IgM)
Prealbumin
Transferrin

Therapeutic Drug Monitoring

Acetaminophen
Carbamazepine
Digoxin

Gentamicin
Lithium
Phenobarbital
Phenytoin
Salicylate
Theophylline
Tobramycin
Valproic Acid

Immunoassay

Ferritin
T3 (Total)
T3 Uptake/T-Uptake
T4 (Free)
T4 (Total)
TSH
Vitamin B ₁₂

Liquichek™ Cardiac Markers Plus Control LT

A comprehensive cardiac assessment control designed to monitor precision of Troponin I at lower levels.

- Offers 7 analytes
- 10 day open-vial (punctured) stability for all analytes when stored at 2–8°C
- 3 year shelf life when stored at -20°C to -50°C.

Analytes

CK (Total)	Myoglobin
CK-MB Isoenzyme	NT-proBNP
CRP	Troponin I
Digitoxin	

Liquichek™ Diabetes Control

A liquid, human whole blood based control to monitor the precision of Hemoglobin A1C. Available in three critical levels.

- 14 day open-vial (punctured) stability for all analytes when stored at 2–8°C
- 3 year shelf life when stored at -10°C to -50°C

Analytes

Hemoglobin A1C



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

Liquichek™ Ethanol/Ammonia Control

Designed to monitor the precision of Ethanol and Ammonia test procedures across the clinical range.

- 20 day open-vial (punctured) stability for all analytes when stored at 2–8°C
- 2 year shelf life when stored at 2–8°C

Analytes

Ammonia
Ethanol

Liquichek™ Immunoassay Plus Control

An efficient solution for QC of routine Immunoassay and Therapeutic Drug Monitoring analytes.

- Offers more than 40 analytes
- 4 day open-vial (punctured) stability for all analytes when stored at 2–8°C
- 2 year shelf life when stored at -20°C to -50°C

Analytes

Therapeutic Drug Monitoring

Acetaminophen
Amikacin
Caffeine
Carbamazepine
Digoxin
Gentamicin
Lidocaine
Lithium
N-Acetylprocainamide (NAPA)
Phenobarbital
Phenytoin
Procainamide

Salicylate
Theophylline
Tobramycin
Tricyclic Antidepressants (TCA)
Screen
Valproic Acid
Vancomycin

Immunology

Immunoglobulin A (IgA)
Immunoglobulin E (IgE)
Immunoglobulin G (IgG)
Immunoglobulin M (IgM)

Immunoassay

Alphafetoprotein (AFP)
CEA
Estradiol
Ferritin
Folate
FSH
HCG-β Subunit
LH
Prolactin
PSA
PSA (Free)
T3 (Free)

T3 (Total)
T3 Uptake/T-Uptake
T4 (Free)
T4 (Total)
TSH
Vitamin B₁₂

Cardiac

CK-MB Isoenzyme

General Chemistry

Iron

Liquichek™ Immunology Control

A liquid, trilevel control to monitor the precision of serum proteins and related analytes.

- Offers 31 analytes
- 5 day open-vial (punctured) stability for all analytes when stored at 2–8°C
- 2 year shelf life when stored at -20°C to -50°C

Analytes

α-1-Acid Glycoprotein
α-1-Antitrypsin
α-2-Macroglobulin
β-2-Microglobulin
Albumin
Antistreptolysin O (ASO)
Apolipoprotein A-1

Apolipoprotein B
C1 Inhibitor
Ceruloplasmin
Complement C3
Complement C4
CRP
Cystatin C

Ferritin
Haptoglobin
Immunoglobulin A (IgA)
Immunoglobulin E (IgE)
Immunoglobulin G (IgG)
Immunoglobulin G Subclasses 1-4
Immunoglobulin M (IgM)

Kappa Light Chain
Lambda Light Chain
Prealbumin
Protein (Total)
Retinol Binding Protein
Rheumatoid Factor
Transferrin



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

Liquichek™ Pediatric Control

A liquid, bilevel control to monitor the precision of analytes important to neonatal testing. Four Bilirubin parameters are available at two clinically significant levels.

- Offers 10 analytes
- 14 day open-vial (punctured) stability for all analytes when stored at 2–8°C
- 3 year shelf life when stored at -20°C to -50°C

Analytes

Bilirubin (Direct)	Glucose
Bilirubin (Total)	Magnesium
Caffeine	Potassium
Calcium	Sodium
Chloride	Theophylline

Liquichek™ Spinal Fluid Control

A liquid, bilevel control to monitor the precision of assays for proteins and other critical analytes in cerebrospinal fluid.

- Offers 8 analytes
- 30 day open-vial (punctured) stability for all analytes when stored at 2–8°C
- 2 year shelf life at 2–8°C

Analytes

Albumin	Lactate (Lactic Acid)
Chloride	Lactate Dehydrogenase (LDH)
Glucose	Protein (Total)
Immunoglobulin G (IgG)	Sodium

Liquichek™ Tumor Marker Control

A liquid, trilevel human based control designed to monitor the precision of tumor marker testing procedures.

- Offers 11 analytes, including CA 125, CA 15-3 and CA-19-9
- 10 day open-vial (punctured) stability for all analytes when stored at 2–8°C
- 2 year shelf life when stored at -20°C to -50°C.

Analytes

Alphafetoprotein (AFP)	Ferritin
β-2-Microglobulin	hCG
CA 15-3	Prolactin
CA 19-9	PSA (Free)
CA 125	PSA (Total)
Carcinoembryonic Antigen (CEA)	

Liquichek™ Urine Chemistry Control

A liquid, bilevel human urine based control for monitoring urine chemistry test procedures.

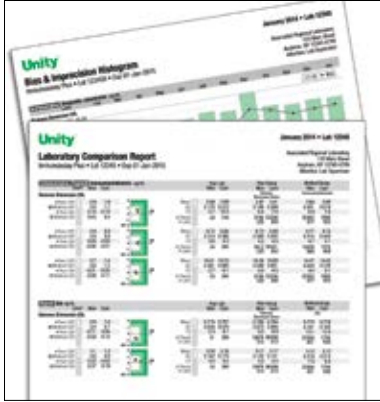
- Offers 13 analytes
- 30 day open-vial (punctured) stability for all analytes when stored at 2–8°C
- 2 year shelf life when stored refrigerated at 2–8°C

Analytes

Amylase	Phosphorus
Calcium	Potassium
Chloride	Protein (Total)
Creatinine	Sodium
Glucose	Urea Nitrogen
Magnesium	Uric Acid
Microalbumin	



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



Unity™ QC Data Management Solutions

Bio-Rad Quality Controls are part of the Unity™ Program with its industry-leading features for benchmarking laboratory performance.

- Benefit from peer data generated from more than 40,000 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

You can consolidate your entire laboratory into one QC system through our Unity™ Program. This sophisticated software includes best-in-class tools to support releasing results, benchmarking laboratory performance, and designing your QC system.

Unity™ Software Solutions

Contact your local sales representative for more information about all of our QC Data Management Solutions.



Unity Real Time®



Expert QC Data Management Solution for Desktop or Online Users

- Facilitate regulatory 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 Advisor™ module available to comply with German 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)
- Access UnityWeb® via mobile device



Basic



Advanced



Desktop based



Web based



Mobile

Want to know more? Discover the power of Unity™ Solutions at www.bio-rad.com/qc-datamangement

Ordering Information

Cat #	Product Name	Quantity	Cat #	Product Name	Quantity
Liquid Assayed MultiquaI® Control			Liquichek™ Immunology Control		
337	Level 1	12 x 2.5 mL	281	Level 1	6 x 2.5 mL
338	Level 2	12 x 2.5 mL	282	Level 2	6 x 2.5 mL
339	Level 3	12 x 2.5 mL	283	Level 3	6 x 2.5 mL
338X	Trilevel MiniPak (1 of each level)	3 x 2.5 mL	282X	Trilevel MiniPak (1 of each level)	3 x 2.5 mL
Liquichek™ Cardiac Markers Plus Control LT			Liquichek™ Pediatric Control		
297	Level 1C	6 x 2.5 mL	176	Level 1	6 x 2.5 mL
298	Level 2	6 x 2.5 mL	177	Level 2	6 x 2.5 mL
299	Level 3	6 x 2.5 mL	178X	Bilevel MiniPak (1 of each level)	2 x 2.5 mL
298X	Trilevel MiniPak (1 of each level)	3 x 2.5 mL	Liquichek™ Spinal Fluid Control		
Liquichek™ Diabetes Control			303	Level 1	6 x 2.5 mL
291	Level 16 x 2 mL	304	Level 2	6 x 2.5 mL
292	Level 26 x 2 mL	302X	Bilevel MiniPak (1 of each level)	2 x 2.5 mL
293	Level 36 x 2 mL	Liquichek™ Tumor Marker Control		
292X	Trilevel MiniPak (1 of each level)3 x 2 mL	27114	Level 16 x 2 mL
Liquichek™ Ethanol/Ammonia Control			27115	Level 26 x 2 mL
271	Level 1	6 x 2.5 mL	27116	Level 36 x 2 mL
272	Level 2	6 x 2.5 mL	27115X	Trilevel MiniPak (1 of each level)3 x 2 mL
273	Level 3	6 x 2.5 mL	Liquichek™ Urine Chemistry Control		
272X	Trilevel MiniPak (1 of each level)	3 x 2.5 mL	195	Level 1	12 x 2.5 mL
Liquichek™ Immunoassay Plus Control			196	Level 2	12 x 2.5 mL
267	Level 1	12 x 2.5 mL	198X	Bilevel MiniPak (1 of each level)	2 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			

Bio-Rad is Dedicated to Advancing QC Solutions Worldwide

Quality Controls

Broad portfolio of independent controls provides an unbiased assessment of methods and system performance.

QC Data Management Solutions

Manage and interpret your QC results using the industry-leading Unity™ Interlaboratory Program and benefit from data generated by users worldwide.

Customer Support

Receive the backing of full-service technical and customer support trained to handle your inquiries in a prompt, knowledgeable manner.

Education

We offer the educational tools and resources to help improve your ability to produce reliable test results.

Efficiently manage your QC product inserts with My eInserts™. Visit www.myeinserts.com for more information.



**Bio-Rad
Laboratories**

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

Clinical
Diagnostics Group

Website www.bio-rad.com/qualitycontrol Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium +32 (3)710-53-00 Brazil +55 (31)3689-6600 Canada 1-514-334-4372 China 86-21-61698500 Czech Republic 420-241-430-532 Denmark +45-4452-1000 Finland 358-9-804-22-00 France 33-1-47-95-60-00 Germany +49 (0)89-318-840 Greece 30-210-7774396 Hong Kong 852-2789-3300 Hungary +36-1-459-6100 India 1800-180-1224 Israel 972-3-9636050 Italy +39-02-216091 Japan 81-3-6361-7070 Korea 82-2-3473-4460 Mexico +52 (55)5488-7670 The Netherlands +31-318-540666 New Zealand 64-9-415-2280 Norway +47-23-38-41-30 Poland 48-22-3319999 Portugal 351-21-472-7700 Russia +7-495-721-1404 Singapore 65-6415-3170 South Africa 27-11-442-85-08 Spain 34-91-590-5200 Sweden +46-8-555-127-00 Switzerland +41 (0)26-674-55-05/06 Taiwan 886-2-2578-7189 Thailand 662-651-8311 United Kingdom +44 (0)20-8328-2000