



QC Data Management Solutions

Introduce your laboratory to a whole new world of Unity data management solutions





Commitment to Quality Control

Using our comprehensive product line of independent controls is an important step toward helping you improve the reliability of your laboratory results. The ability to successfully manage and interpret your quality control results is also essential to producing solid, dependable laboratory outcomes.

Bio-Rad Laboratories offers a vision and a commitment to quality control that is unsurpassed in the clinical laboratory industry. Today, your laboratory can be connected to the Unity Interlaboratory Program, the world's largest community of quality control users. Improve patient test results and cost effectiveness with quality control products, quality control data management solutions, industry-leading interlaboratory comparisons and personalized service from Bio-Rad.



Regulatory Requirements

"The laboratory shall participate in an interlaboratory comparison programme(s) . . ."

– ISO 15189:2012(E), Subclause 5.6.3.1



Unity Interlaboratory Program

Participation in an interlaboratory program is extremely important for laboratories to help ensure the reliability and precision of test systems. No program in the world is more powerful and effective at providing information and helping to improve laboratory analytical performance than the Unity Interlaboratory Program.

- Benefit from peer QC data generated from more than 59,000 instruments in 100+ countries
- Recognize analytical process improvement opportunities
- Increase confidence in proficiency survey outcomes
- Identify unrecognized trends or shifts that may occur between proficiency surveys
- Receive on-demand interlaboratory comparison reports
- Receive monthly reports with regular data submissions
- Submit data in real time using one of our Unity software solutions
- Detect and identify potentially important analytical errors
- Meet accreditation and regulatory requirements

Unity January 2018

InstantQC Laboratory Comparison Report
 Unassayed Chemistry • Lot 12345 • Exp 01-Dec-2020

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

Parameter	ISE Indirect method			Level	Year	Lab	Peer Group		Method Group		
	Level	Min	Com				Min	Com	Min	Com	
Roche cobas 8000											
Peer CR	0.2	0.4			Mean	4.11	4.08	4.04	4.05	4.04	4.04
Method CR	0.1	0.5			SD	0.021	0.040	0.05	0.037	0.024	0.041
Peer SD	0.85	0.34			CV	0.5	1.0	2.2	2.7	1.8	2.0
Method SD	0.39	0.52			# Points	7	24	107	100	102	100
					# Labs			108	106	102	100
Peer CR	0.4	0.7			Mean	6.39	6.33	6.34	6.35	6.31	6.31
Method CR	0.1	0.8			SD	0.021	0.027	0.025	0.032	0.014	0.019
Peer SD	0.50	0.21			CV	0.3	1.1	1.3	1.6	1.8	1.9
Method SD	0.40	0.10			# Points	7	24	112	114	110	106
					# Labs			103	103	815	878
Roche cobas 8000					Mean	83.00	83.09	82.48	82.32	85.09	85.66
Peer CR	0.3	0.5			SD	0.400	0.014	1.19	1.36	2.14	2.26
Method CR	0.2	0.2									

Receive On-demand InstantQC Reports

Get started at QCNet.com today!

InstantQC Reports allow you to compare your control results to other laboratories at any time. These on-demand reports are particularly useful for troubleshooting problems with test system performance.



The Advantages of Participating in the Unity Interlaboratory Program

Detecting Errors in Your Test System

Without a high quality interlaboratory program, your laboratory may not be aware of gradual or sudden changes in your test system that may be caused by events such as reagent or calibrator reformulations, standardization changes, or instrument software changes. An interlaboratory program may offer early awareness of shifts and trends to help avoid costly test repeats and unnecessary troubleshooting.

Your Proficiency Program is Not Enough

Proficiency surveys provide a snapshot in time and should not be used as a substitute for daily quality control with an interlaboratory program.

- Acceptable performance on the day of the proficiency testing does not guarantee testing reliability every day
- Participation in an interlaboratory program is the recommended way to improve confidence in test system performance every day, especially with an interlaboratory program that offers an option to receive instant interlaboratory feedback on-demand

Some advantages for participating in an interlaboratory QC program are:

- Verifying that a laboratory is producing QC results that are consistent with other laboratories using the same measurement procedure, and thus demonstrating that the laboratory is using the measurement procedure correctly

Detecting and identifying bias in a measurement procedure

- Bias can be caused by events such as reagent or calibrator lot changes or reformulations, changes in calibration traceability to reference systems, or instrument software changes. Comparison of an individual laboratory's QC result to a peer group mean value can identify a trend or shift, or ascertain if other laboratories are experiencing the same changes (see Subchapter 3.3.1)."

Supplementing PT/EQA programs

- PT/EQA programs verify performance at a point in time. Acceptable performance on the day of PT/EQA testing does not guarantee testing reliability every day because errors in a measurement procedure can occur at any time. In addition, the interlaboratory QC data can be used to investigate a failed or questionable PT/EQA result.

– CLSI, C24, 4th ed. Statistical Quality Control for Quantitative Measurement Procedures: Principles and Definitions. 7.2 Using Interlaboratory Quality Control to Assess a Quality Control Program.



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.

Internet-based solutions such as UnityWeb and Unity Real Time online eliminate the need to install and update software locally in your laboratory, and reduce the amount of support necessary from on-site IT staff.

Desktop software solutions, like Unity Real Time, are appropriate if your internet connection is not sufficient, or if you prefer maintaining local software.

	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



Basic



Advanced



Desktop based



Web based



Mobile

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

Basic User Options

Test Information	Date & Time	Level 1	Level 2	Level 3
U88	03/02/2012 12:00:00 A	0.24	0.24	0.24
U89	03/02/2012 12:00:00 A	0.24	0.24	0.24
U90	03/02/2012 12:00:00 A	0.24	0.24	0.24
U91	03/02/2012 12:00:00 A	0.24	0.24	0.24
U92	03/02/2012 12:00:00 A	0.24	0.24	0.24
U93	03/02/2012 12:00:00 A	0.24	0.24	0.24
U94	03/02/2012 12:00:00 A	0.24	0.24	0.24
U95	03/02/2012 12:00:00 A	0.24	0.24	0.24
U96	03/02/2012 12:00:00 A	0.24	0.24	0.24
U97	03/02/2012 12:00:00 A	0.24	0.24	0.24
U98	03/02/2012 12:00:00 A	0.24	0.24	0.24
U99	03/02/2012 12:00:00 A	0.24	0.24	0.24
U100	03/02/2012 12:00:00 A	0.24	0.24	0.24

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)

Test Information	Date & Time	Level 1	Level 2	Level 3
U88	03/02/2012 12:00:00 A	0.24	0.24	0.24
U89	03/02/2012 12:00:00 A	0.24	0.24	0.24
U90	03/02/2012 12:00:00 A	0.24	0.24	0.24
U91	03/02/2012 12:00:00 A	0.24	0.24	0.24
U92	03/02/2012 12:00:00 A	0.24	0.24	0.24
U93	03/02/2012 12:00:00 A	0.24	0.24	0.24
U94	03/02/2012 12:00:00 A	0.24	0.24	0.24
U95	03/02/2012 12:00:00 A	0.24	0.24	0.24
U96	03/02/2012 12:00:00 A	0.24	0.24	0.24
U97	03/02/2012 12:00:00 A	0.24	0.24	0.24
U98	03/02/2012 12:00:00 A	0.24	0.24	0.24
U99	03/02/2012 12:00:00 A	0.24	0.24	0.24
U100	03/02/2012 12:00:00 A	0.24	0.24	0.24

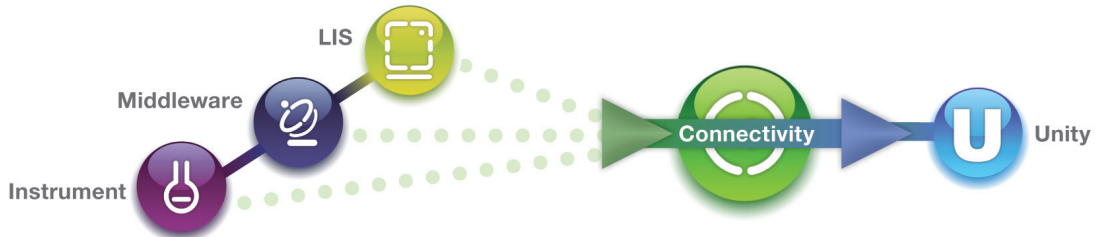
UnityWeb



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



Basic

WebConnect



Automated Uploads to Unity Software

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

Advanced

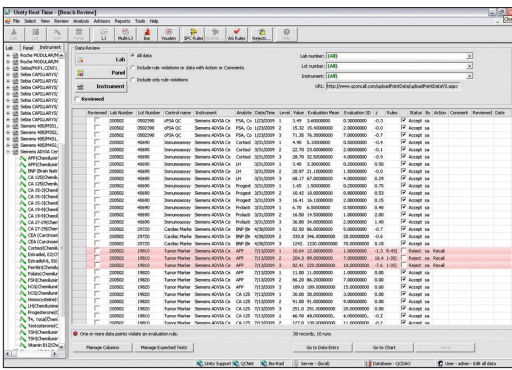
UnityConnect



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

- 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

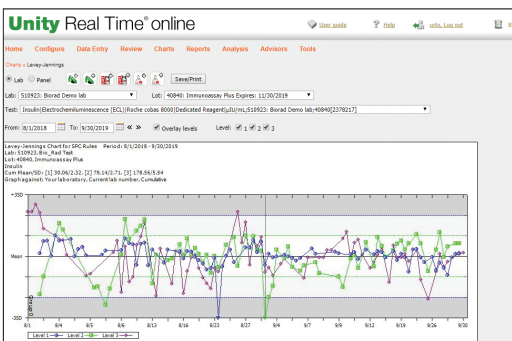
Advanced User Options



Unity Real Time

Expert QC Data Management Solution for Desktop Users

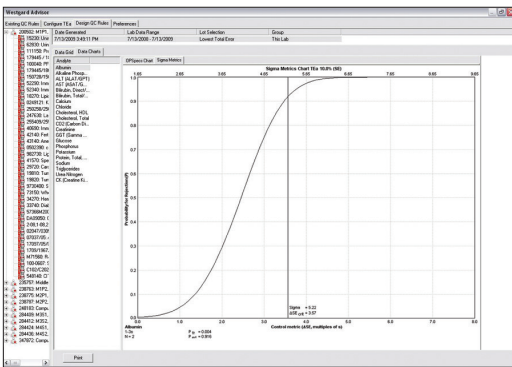
- Facilitate regulatory compliance under CLIA and ISO 15189
- Improve real-time bench and supervisor QC data review
- Implement best QC rules when used with Westgard Advisor
- Run validation with comprehensive audit trails
- Advanced charts and reports for data analysis
- Reduce non-essential retests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



Unity Real Time online

Expert QC Data Management Solution for Online Users

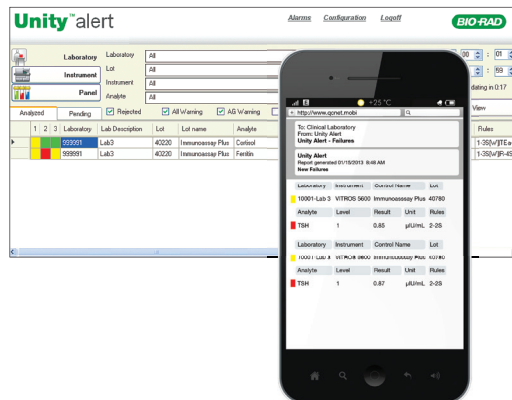
- Facilitate regulatory compliance under CLIA and ISO 15189
- Improve real-time bench and supervisor QC data review
- Implement best QC rules when used with Westgard Advisor
- Run validation with comprehensive audit trails
- Advanced charts and reports for data analysis
- Reduce non-essential retests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



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



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





Bio-Rad Mission: Control



Take the guesswork out of your QC strategy

Employing a robust quality control system is an essential step for any laboratory hoping to provide reliable results for appropriate patient care. Unfortunately for clinical laboratories, selecting the right QC rules and frequency can be challenging, and questions often arise.

- How often should controls be run for specific assays— is there an optimal frequency?
- How can QC be managed to minimize the risk of releasing incorrect patient results?

Bio-Rad Mission: Control is the first objective risk management solution for quality control in laboratories. The intuitive interface allows users to focus on patient safety by assessing the risk of reporting incorrect patient results in their laboratory.

Bio-Rad Mission: Control allows you to:

- 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

A patient risk approach to Quality Control is a progressive step toward improving quality assurance and patient care. Bio-Rad can help you understand what this means for your laboratory and how to take steps to implement an effective approach for Quality Control within your analytical test phase.

For more information, visit www.qcnet.com/missioncontrol



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

QCNet is the Leading QC Portal on the Internet

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

Quick Links to the most-used features

Unity Reports

- Get your Unity Reports in just a few clicks.

My elnserts

- Access product inserts

EQAS

- View Reports
- Access EQAS Online
- Training/Tutorials
- View Product Information

Access country-specific information

At the bottom of the QCNet home page, click on the country name shown to access the pop-up menu, then select the desired country*.

*Not all countries are available.

Explore the drop-down menus to find more online QC tools on QCNet

Quality Controls	QC Data Management	Resources	Support
COVID-91	UnityWeb	Training & Education	My Account - Log in
InteliQ Load and Go	Unity Real Time online	Technical Documents	Lab Admin
Cardiac Marker	Unity Software Solutions	Frequently Asked Questions	Contact us
Infectious Disease	Peer Reporting Program	Barcode Label Generator	
Immunoassay	Mission: Control	Value Assignment Program	
Chemistry	EQAS		
Diabetes & Hemoglobin	Linearity Web		
Immunology	Data Submission Calendar		
Blood Gas			
Toxicology			
Tumor Markers			
Therapeutic Drug Monitoring			
Hematology			
Coagulation			
Urinalysis			

Ordering Information

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

Cat #	Description
UnityWeb	
870-1	UnityWeb Annual subscription

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

Cat #	Description
UnityConnect	
820-1	UnityConnect Software. Annual subscription
825i	Unity Remote Installation & Training Single use

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

Cat #	Description
Unity Real Time online	
804-W1	Unity Real Time online Annual subscription

Cat #	Description
Westgard Advisor	
811-1	Westgard Advisor Annual subscription

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

Cat #	Description
Bio-Rad Mission: Control	
1100002	Mission: Control Annual Subscription
1100005	Mission: Control Subscription and Service Bundle

EQAS	Independent QC	Unity
An independent, external assessment of laboratory performance in comparison to your peers.	Ongoing, proactive, unbiased daily QC that helps identify errors as they occur or begin to trend.	QC Data Management tools that help you create a strategy to reduce risk and streamline QC workflow.
QCNet.com/eqas	QCNet.com/independentqc	QCNet.com/datamanagement



**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 **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) 721-14-04 **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