



# 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 third party 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 care and cost effectiveness with quality control products, quality control data management solutions, industry-leading interlaboratory comparisons and personalized service from Bio-Rad.

*"The laboratory shall participate in interlaboratory comparisons. . ."*  
– ISO 15189:2007(E), Subclause 5.6.4

# 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 36,000 instruments in 67 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

## Receive On-Demand InstantQC™ Reports

### Get started at QCNet.com today!

InstantQC™ Reports permit 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.

Improved

## 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 can offer early awareness of shifts and trends to help avoid costly test repeats and unnecessary troubleshooting.

### Improving Outcomes on Proficiency Surveys

An interlaboratory program can also increase your confidence in proficiency surveys. If your laboratory compares well to other laboratories in an interlaboratory program, you are also likely to compare well to other laboratories participating in your proficiency testing program.

### Your Proficiency Program is Not Enough

Proficiency surveys merely 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 best way to be confident in test system performance every day, especially with an interlaboratory program that offers an option to receive instant interlaboratory feedback on-demand.

### Statistically Relevant Peer Data

Monthly Unity™ Interlaboratory Reports ensure that peer group sizes are maximized for regular, periodic review of your test systems. These reports are comprehensive and represent the largest peer groups in the industry.

*"For laboratory self-evaluation, peer-related bias and relative imprecision are useful parameters. Participation in an interlaboratory program provides an effective mechanism to complement external quality assessment (proficiency surveys) programs. Consequently, a laboratory should actively participate in interlaboratory QC programs when such programs are available."*

– CLSI (formerly NCCLS), C24-A3, Vol. 26, No 25.

# 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.

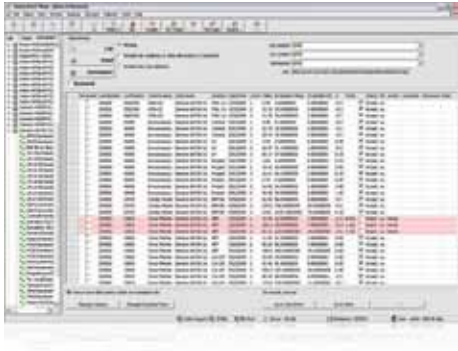
Want to know more?

Discover the power of Unity™ Solutions at [www.QCNet.com/Unity](http://www.QCNet.com/Unity)

	Basic User Options		Advanced User Options	
	 UnityWeb®	 Unity Real Time® LT	 Unity Real Time® online	 Unity Real Time®
<b>Type of Solution</b>				
Web Service (Internet-Based)	x		x	
Desktop Software		x		x
<b>Unity™ Interlaboratory Reports</b>				
Monthly Reports	x	x	x	x
InstantQC™ Reports	x	x	x	x
<b>Basic Intralaboratory Charts &amp; Reports</b>				
Westgard Rules	x	x	x	x
Various Charts and Reports	x	x	x	x
<b>Additional Features</b>				
Westgard Advisor™ (Subscription sold separately)	x		x	x
Bench Review	x		x	x
Bench & Supervisor Data Review			x	x
Analytical Goals			x	x
Dynamic Data Set Comparisons			x	x
<b>Database Platform</b>				
Bio-Rad Hosted	x		x	
Installed on Local Laboratory PC		x		x
<b>Connectivity (Sold separately)</b>				
WebConnect™	x		x	
UnityConnect™	x	x	x	x

*"The laboratory shall design internal quality control systems that verify the attainment of the intended quality of results. It is important that the control system provide staff members with clear and easily understood information on which to base technical and medical decisions. . ."*

– ISO 15189:2007(E), Subclause 5.6.1



## Unity Real Time®

### Expert QC Data Management Solution for Desktop or Online Users

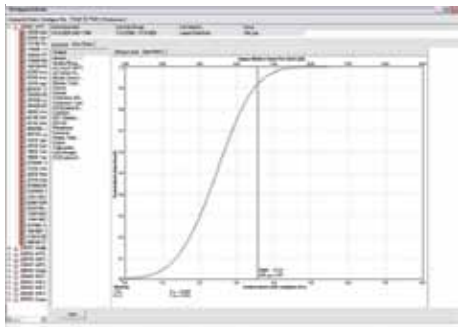
- Facilitate regulatory compliance under CLIA and ISO 15189
- Improve real-time bench and supervisor QC data review
- Implement best QC rules when used with Westgard Advisor™
- Run validation with comprehensive audit trails
- Advanced charts and reports for data analysis
- Reduce non-essential retests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



## UnityWeb® & Unity Real Time® LT

### Entry-level QC Data Management Solutions

- Basic QC rules, charts and reports
- Easily upgradable to Unity Real Time® for more advanced tools and features
- Upload QC data points from an LIS, middleware or instrument (optional)
- No software to install with UnityWeb®



## 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

## Automated Uploads to Unity™ Software

### 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.

- Eliminates manual keying of QC data
- Enables standard LIS QC reports to be used
- Provides automatic data import
- Provides real-time connection
- Transparent to the flow of data to the LIS
- Captures and analyzes QC data economically from laboratory instruments not connected to the LIS
- Easy-to-install optional interface hardware
- Choose from soft, serial, ethernet and wireless 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®.

- Eliminates manual keying of data
- Enables standard LIS QC reports to be used
- No software to install

# 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](http://www.QCNet.com).

**EQAS**

- View Reports
- Upload eForms
- Register
- View Product Information

**QCNet International**

**Local QCNet™**  
Visit your local QCNet™ portal, which will provide specific information related to your country\*.

Log In/Register | QCNet International | Contact Us

Home Unity™ Interlab LabTools™ EQAS QC Education QC Documents Support Lab Admin

Monday, November 08, 2010

**I'm a Quality System**

**QCNet Made Some Changes**

- Minimized Home Page**  
We have consolidated the content into less vertical space by creating the Documents & Resources area.
- Expanded Side Navigation**  
Get easier access to the most popular services through the expanded side navigation. One click and you're there!

**QCNet Made Some Changes What's New from Bio-Rad QC**

- Introducing the New & Improved Ligaschek™ Specialty Immunoassay Control**  
The enhancements include longer open vial stability for Anti-Tg & Anti-TPO and expanded testing methods for 25-OH Vitamin D.
- eInserts™ XML is Now Available From Bio-Rad**  
eInserts™ XML is an innovative solution to load assayed values directly into your Abbott ARCHITECTS Systems 7.0.
- Introducing the New Ligaschek™ Opiste Control from Bio-Rad**  
This third party control is designed to monitor urine toxicology testing procedures for a broad class of opiates.

**View My Unity Reports**

**Find Products & Information**

Search MSDS or search  
Search Bio-Rad Website

**Categories**

- Unity Interlaboratory Program**
  - My Unity Reports
  - InstantQC Reports
  - Unity Bulletins by Date Issued
  - Submission Calendar
- Data Management**
  - UnityWeb
  - UnityWeb 2
  - Unity Real Time online
  - Products & Store Info
- EQAS**
  - My EQAS Reports
  - User Documentation
  - eForms
  - Submission Calendar
  - Products & Store Info
- Training & Education**
  - QCNet University
  - Training Center
- LabTools**
  - Linearity
  - Urinalysis
  - Barcode Label Generator
  - eInserts XML
- Quality Controls**
  - Immunoassay
  - Serum Chemistry
  - Immunology & Protein

**Find More Online QC Tools on QCNet™**

- Linearity Web™
- eInserts™ XML
- Product Inserts (My eInserts™)
- Barcode Label Generator
- Urinalysis Tools
- QCNet™ University
- MSDS
- Code Lists for Unity™

\*Not all countries are available.



**Bio-Rad Laboratories**

For further information, please contact your local Bio-Rad office or learn more online at [www.QCNet.com/unity](http://www.QCNet.com/unity).

Clinical  
Diagnostics Group

Website [www.bio-rad.com/qualitycontrol](http://www.bio-rad.com/qualitycontrol) U.S. 1-800-2BIO-RAD Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium 32-9-385-5511 Brazil 5521-3237-9400 Canada 1-514-334-4372 China 86-21-64260808 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 1-800-180-1224 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 7-495-721-14-04 Singapore 65-6415-3170 South Africa 27-11-442-85-08 Spain 34-91-590-5200 Sweden 46-8-555-127-00 Switzerland 41-61-717-95-55 Thailand 662-651-8311 United Kingdom +44-(0)20-8328-2000