



Unity™ Alert

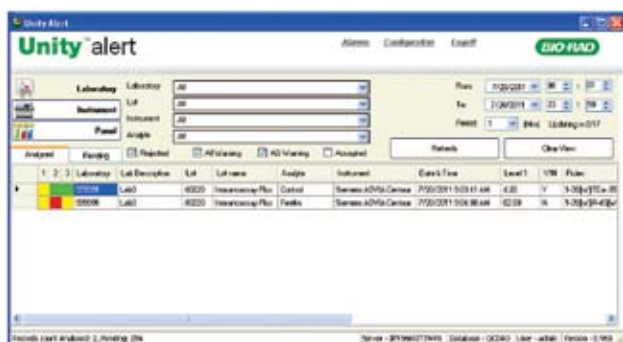
Automatic QC Status Notification Tool
to Help Monitor Instrument Performance



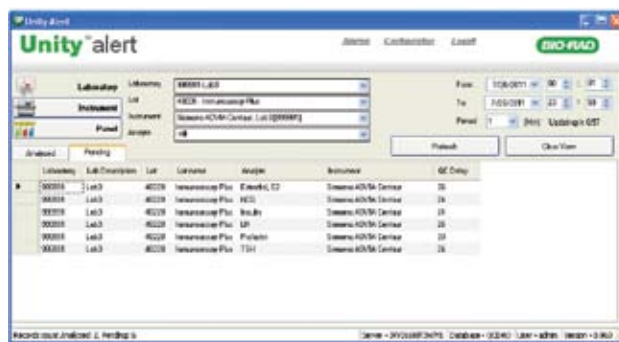
Unity™ Alert

Automatic QC Status Notification Tool

- 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.
- Setting up alert rules is easy and flexible, and can even be set up for specific instruments or tests.
- Lab supervisors and managers can receive emails remotely and stay informed regarding QC issues experienced in their laboratory even while they are out of the office.
- The color-coded displays on the Unity™ Alert client are helpful to draw attention to the most urgent QC issues.
- Unity™ Alert notifications are provided even when Unity Real Time® 2 is not in use.
- Unity™ Alert is available with Unity Real Time® 2.



Unity™ Alert provides a color-coded display for QC failures, helping draw attention to the most urgent issues.



Unity™ Alert can identify missing QC runs and indicate the delay since the test was last run, helping to ensure proper monitoring of instrument performance.



Unity™ Alert is part of the Unity™ Interlaboratory Program suite of data management products. Participation in the Unity™ Interlaboratory Program can help improve the quality of your laboratory test results and help you meet regulatory and accreditation requirements worldwide. You may also subscribe to one of our Unity™ QC Data Management Solutions designed to improve the effectiveness of your statistical process control. Ask your Bio-Rad Account Representative for more information.

Discover the power of Unity™ Solutions at www.QCNet.com.



**Bio-Rad
Laboratories**

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

**Clinical
Diagnostics Group**

Website www.bio-rad.com/diagnostics **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** 91-124-4029300 **Israel** 972-3-9636050 **Italy** +39-02-216091 **Japan** 81-3-6361-7070 **Korea** 82-2-3473-4460 **Mexico** 52(55)5200-0520 **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** 65-6415-3188 **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