



EQAS[®] Programs for Blood Banks

Donating blood saves lives – but so does reporting reliable patient results

Recognized as the worldwide leader in providing laboratory quality controls, Bio-Rad is committed to providing a range of External Quality Assessment (EQA) programs to assist blood banks and transfusion centers increase their confidence when reporting test results. When used in conjunction with daily quality controls, these external programs give laboratories added confidence in reporting their test results, ensuring the safety of patients and blood supplies.

Bio-Rad EQAS[®] Programs:

- Are developed and maintained under a compliant and comprehensive quality management system
- Provide participants with the confidence and assurance of high quality programs supported by human based samples of measured stability and homogeneity
- Offer comprehensive test menus that cover a wide range of parameters suitable for manual and automated methods
- Objectively compare results to other laboratories using the same instruments and reagents

Enroll today to utilize convenient and easy to interpret comparison reports, international peer groups, and unparalleled expert customer service and support to take your laboratory's test system performance to the next level.

BIO-RAD

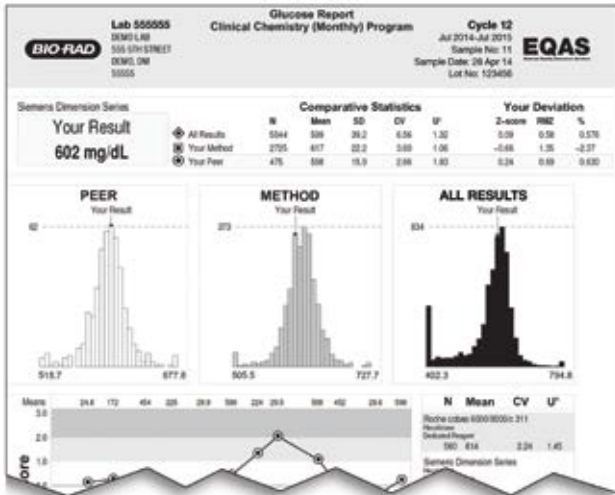
Online EQAS® Resources

Easily access product information, product inserts and user documentation through www.QCNet.com. This secure website allows registered participants access to training material, including a Troubleshooting Checklist, reference method values and customer notifications. It also provides entry to EQAS® Online, where you can manage your contact details and test configurations, submit your results and view or print all your EQAS® reports.

EQAS® Reports

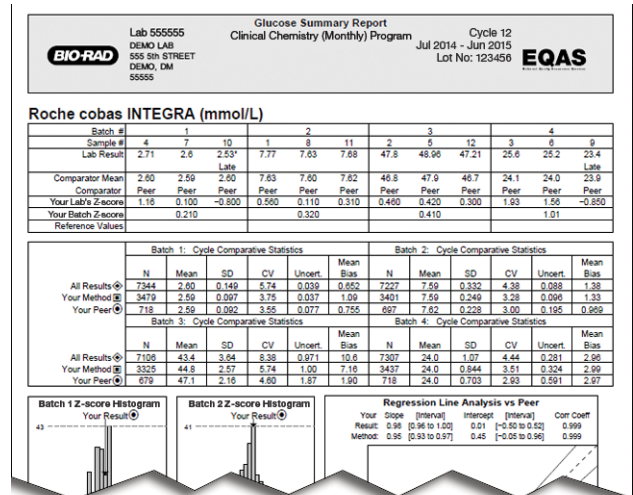
Participants receive regular sample reports once enrolled in a program. These detailed reports are presented in an easy-to-read graphic format. Sample reports are available online within three days of the statistical analysis.

Quantitative Reports



Sample Report

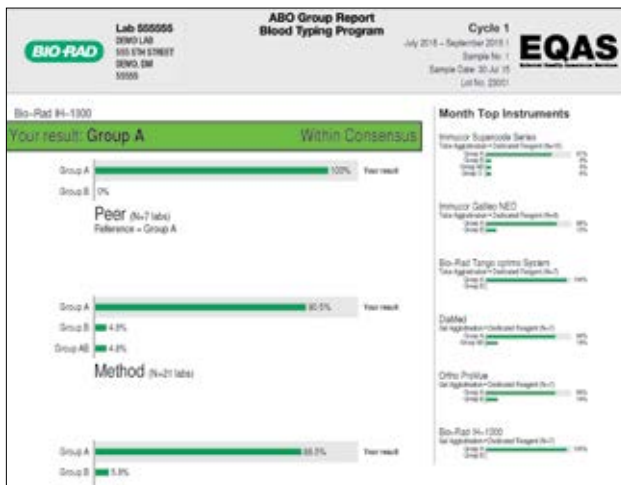
Detailed individual statistical analyses and peer group comparisons are provided.



End-of-Cycle Report

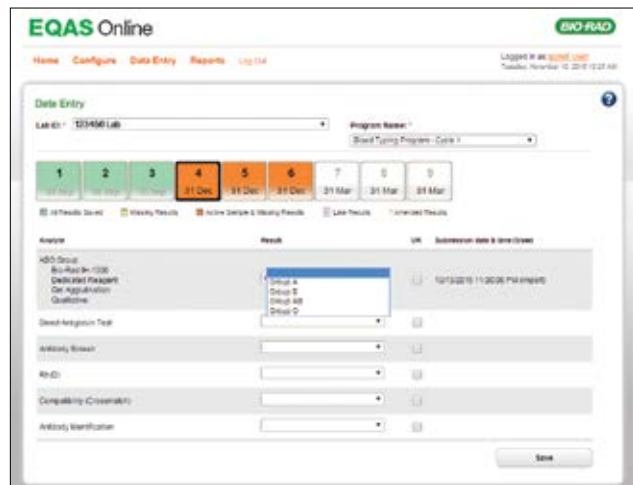
An overall assessment of laboratory performance over the entire cycle in terms of bias among participating laboratories is presented.

Qualitative Report



Sample Report

These reports provide clear and concise information about your qualitative test laboratory performance in the form of histograms and modified Levey-Jennings charts. They include detailed information about the distribution of the qualitative responses and the peer group consensus.



EQAS® Online

Conveniently manage test configurations, submit results and view reports for all programs you participate in. Set up notifications and get email reminders of upcoming due dates and when reports are available for viewing and printing. Accessible through www.QCNet.com.

EQAS® Programs

Blood Typing Program

- Ideal for clinical lab, blood bank and transfusion center testing needs
- Barcoded and ready-to-use tubes for automated platforms
- 3 samples tested every 4 months
- Submit results and view reports through EQAS Online
- 3 x 4 mL Sample and 1 x 2 mL Donor (each shipment)

Parameters

ABO Group	Compatibility (Cross-match)
Antibody Identification	DAT (Direct Antiglobulin Test)
Antibody Screen	Rh(D)

HIV/Hepatitis Program

- Liquid, human source material
- Suitable for serological tests used to detect key HIV and Hepatitis viral antibodies and antigens
- A qualitative and quantitative program is available
- Submit results and view reports each month through EQAS Online
- 12 x 2 mL

Analytes

Anti-HAV	Anti-HBe	Anti-HIV-2
Anti-HAV IgG	Anti-HBs	Anti-HTLV-1
Anti-HAV IgM	Anti-HCV	HBsAg
Anti-HBc	Anti-HIV-1	HBsAg
Anti-HBc IgM	Anti-HIV-1/2	HIV-1 Ag

Syphilis Program

- Liquid, human source material
- Suitable for serological tests detecting treponemal and non-treponemal antibodies
- 12-month cycle
- A qualitative and quantitative program is available
- Submit results and view reports each month through EQAS Online
- 12 x 1.5 mL

Analytes

Non-Treponemal antibodies
Treponema pallidum IgG
Treponema pallidum IgM
Treponema pallidum Antibodies

ToRCH/EBV/MuMZ Program

- Liquid, human source material
- Suitable for serological tests used to detect pediatric diseases and congenital infections
- A qualitative and quantitative program is available
- Submit results and view reports each month through EQAS Online
- 12 x 2 mL

Analytes

Anti-Cytomegalovirus (CMV) IgG
Anti-Cytomegalovirus (CMV) IgM
Anti-Epstein-Barr Virus (EBNA) IgG
Anti-Epstein-Barr Virus (VCA) IgG
Anti-Epstein-Barr Virus (VCA) IgM

Anti-Herpes Simplex Virus (HSV-1) IgG
Anti-Herpes Simplex Virus (HSV-1/2) IgG
Anti-Herpes Simplex Virus (HSV-2) IgG
Anti-Measles Virus (Rubeola) IgG
Anti Mumps Virus IgG


Anti-Rubella Virus IgG
Anti-Rubella Virus IgM
Anti-*Toxoplasma gondii* (Toxo) IgG
Anti-*Toxoplasma gondii* (Toxo) IgM
Anti-Varicella Zoster Virus (VZV) IgG



Ordering Information

Cat #	Configuration
Blood Typing Program*	
	<i>Sample Donor</i>
12000837	Shipment A3 x 4 mL + 1 x 2 mL
12000838	Shipment B3 x 4 mL + 1 x 2 mL
12000839	Shipment C3 x 4 mL + 1 x 2 mL
HIV/Hepatitis Program	
12000815 12 x 2 mL
Syphilis Program	
12000833 12 x 1.5 mL
ToRCH/EBV/MuMZ Program	
12000825 12 x 2 mL

*Aspects of the program are subcontracted from an approved supplier. Refer to myinserts.com or the package insert for current analyte claims.



For more information, contact your local Bio-Rad representative or send an email to eqas_support@bio-rad.com



**Bio-Rad
Laboratories**

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

*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