



EQAS

Participate in an internationally recognized quality assessment program





Bio-Rad EQAS programs are fully accredited to help meet the regulatory needs of today's clinical laboratories.

External quality assessment programs are accepted around the world as invaluable tools used by laboratories to periodically monitor the performance of their test systems. Results are objectively compared to other laboratories using the same methodologies, instruments and reagents. When used in conjunction with daily quality controls, these external programs can give laboratories added confidence in reporting their patient test results.

As the worldwide leader in the development of quality control systems, participants in EQAS can benefit from the experience, knowledge and reliability that Bio-Rad has offered for more than 40 years.

Bio-Rad EQAS programs provide an independent and confidential external assessment of individual laboratory performance.

- Large international database with participants from over 110 countries
- Fully accredited to ISO/IEC 17043:2010 "Conformity assessment—General requirements for proficiency testing"
- Comprehensive, easy-to-read reports
- Technical support from experienced professionals
- Convenient electronic reporting options: EQAS Online and EQAS Mobile
- Each cycle includes 12 high quality samples together in one convenient shipment (except Hematology and Blood Typing Programs)
- Comprehensive analyte menus covering a wide range of diagnostic tests suitable for clinical labs and blood banks
- Analyte levels that challenge both normal and abnormal ranges
- Reference Method Values provided for select programs

EQAS Program Overview

Contact your Bio-Rad representative to place your order and receive a unique laboratory number



Register your test methods



Specimen Package is shipped to you

Getting Started

- Select and enroll in your desired EQAS programs
- Configure your lab information and register tests online or by using your smart phone
- Receive specimen package containing all twelve samples for a specific cycle (Hematology and Blood Typing samples are provided in multiple shipments)

Testing Samples and Submitting Data

- Each EQAS sample is uniquely numbered for easy identification
- Submit results on time; Blood Typing samples should be tested with results reported as soon as they are received
- Sample due dates are provided in a number of places including the sample label, product insert, EQAS Online and the EQAS Calendar on QCNet.com

Test the appropriate EQAS sample as a patient sample



Submit your results in a timely manner



Receive and review your confidential Sample Report

Implement corrective measures based on laboratory performance



Continue process for entire program cycle

Reviewing Reports and Troubleshooting

- Securely access your laboratory specific sample report online within three days of the statistical analysis
- A Certificate of Registration will accompany your Sample 1 Report
- For help interpreting your EQAS reports, view the tutorial on QCNet.com, or seek assistance from your knowledgeable Bio-Rad representative

Receive an End-of-Cycle Report*



Earn a Certificate of Performance

Earn a Certificate of Achievement

Program Cycle Completion

- The End-of-Cycle Report provides an overall assessment of your laboratory's performance in terms of bias
- A Certificate of Performance lists all analytes that passed the performance criteria for the cycle
- A Certificate of Achievement is issued to you if a minimum amount of acceptable data has been submitted for the cycle



*Provided for quantitative programs only.



Online EQAS Resources

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.

Become a Registered QCNet User

Go to www.QCNet.com today and click on the Log In/Register link. Once you've completed the new user form, you will have access to all available EQAS tools.



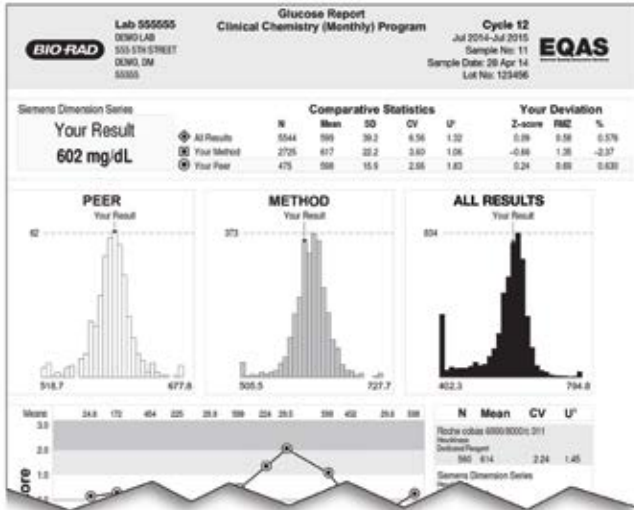
Submitting Results via EQAS Online/Mobile

Conveniently manage your laboratory's test configuration and submit results from any device with an internet connection.

EQAS Reports

Participants receive twelve sample reports and one end-of-cycle report over each program cycle. These detailed reports are presented in an easy-to-read graphic format. Sample reports are available online within three days of the statistical analysis. Additional special reports (Quality Specification, Subgroup and Instrument/Method Summary) are available to registered participants.

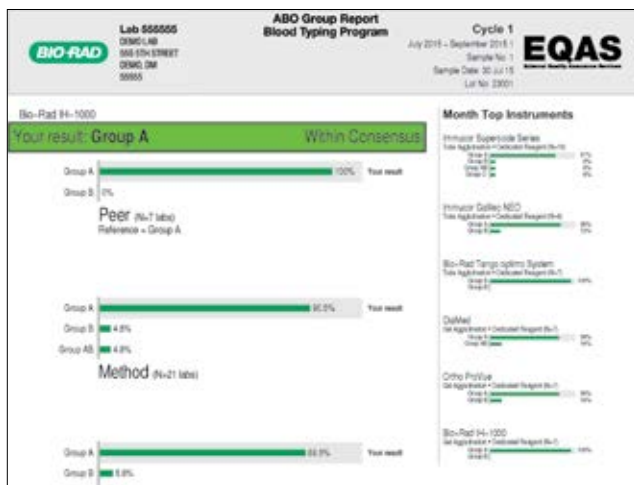
Quantitative Reports



Sample Report

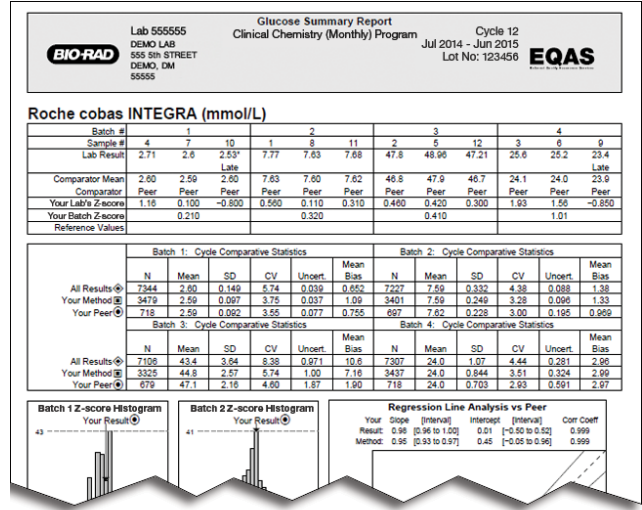
Detailed individual statistical analyses and peer group comparisons are provided.

Qualitative Report



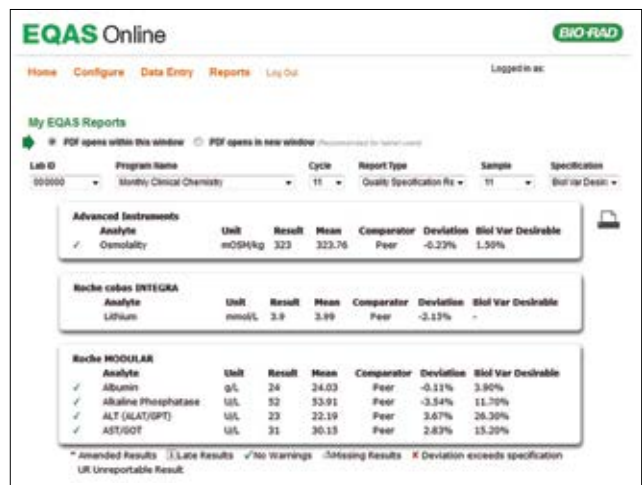
Sample Report

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.



End-of-Cycle Report

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



Quality Specification Report

An evaluation, in terms of percentage deviation, against set performance criteria from key regulatory and scientific organizations (ex. CLIA, BV, RCPA, IQMH, GOST and RiikbÄK), is available online for each sample analyte.



EQAS Programs

Blood Gas Program



- Liquid, aqueous samples
- Comprehensive menu for external assessment of laboratory test performance of Blood Gas, Electrolytes, Glucose, Lactate and Magnesium
- 12-month cycle

Analytes

Calcium (Ionized)	pCO ₂
Chloride	pH
Glucose	pO ₂
Lactate (Lactic Acid)	Potassium
Magnesium	Sodium

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
- Includes 3 case studies per year

Parameters

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

Cardiac Markers Program



- Liquid, human plasma based
- Comprehensive menu of important cardiac assessment markers, plus D-dimer
- 12-month cycle

Analytes

BNP	Myoglobin
CK-MB	NT-proBNP
D-dimer	Troponin I
Homocysteine	Troponin T
hs-CRP	

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

Clinical Chemistry (Monthly) Program



- Lyophilized, human serum based
- 12-month cycle
- Submit results for the appropriate sample every month
- **Reference Method Values** provided for select analytes

Analytes

Acid Phosphatase (Total)	Chloride	Iron Binding Capacity, Unsaturated (UIBC)	Thyroxine (T4), Free
Alanine Aminotransferase (ALT/SGPT)	Cholesterol (HDL)	Lactate (Lactic Acid)	Thyroxine (T4), Total
Albumin	Cholesterol (LDL)	Lactate Dehydrogenase (LDH)	Transferrin
Alkaline Phosphatase (ALP)	Cholesterol (Total)	Lipase	Triglycerides
Amylase (Pancreatic)	Cholinesterase	Lithium	Triiodothyronine (T3), Free
Amylase (Total)	Copper	Magnesium	Triiodothyronine (T3), Total
Aspartate Aminotransferase (AST/SGOT)	Cortisol	Osmolality	Urea
Bilirubin (Direct)	Creatine Kinase (CK)	Phosphorus	Urea Nitrogen
Bilirubin (Indirect)	Creatinine	Potassium	Uric Acid
Bilirubin (Total)	Gamma Glutamyl Transferase (GGT)	Protein (Total)	Zinc
Calcium (Ionized)	Glucose	Sodium	Pilot Analytes
Calcium (Total)	Iron	Thyroid Stimulating Hormone (TSH)	Bile Acid
Carbon Dioxide (CO ₂)	Iron Binding Capacity, Total (TIBC)		

Coagulation Program



- Lyophilized, human plasma based
- 11 parameters for external assessment of laboratory coagulation test performance
- 12-month cycle

Analytes

Antithrombin III (ATIII)	Protein C
APTT	Protein S
D-dimer	PT
Factor VIII	Thrombin Time (TT)
Fibrinogen	Von Willebrand Factor (vWF)
INR	

Ethanol/Ammonia Program



- Liquid, human serum based
- Clinically relevant levels of serum Ethanol and Ammonia
- 12-month cycle

Analytes

Ammonia
Ethanol

Hematology Program*



- Liquid product with human RBCs
- Choice of 11 basic hematology parameters
- For use on manual or automated analyzers
- Convenience of primary pierceable tubes
- 12-month cycle consisting of 4 separate shipments
- Submit results for the appropriate sample every month

Parameters

Hematocrit (HCT)	Platelets (PLT)
Hemoglobin	Red Blood Cells (RBC)
Mean Corpuscular Hemoglobin (MCH)	Red Blood Cell Distribution Width (RDW)
Mean Corpuscular Hemoglobin Concentration (MCHC)	Red Blood Cell Distribution Width-SD (RDW-SD)
Mean Corpuscular Volume (MCV)	White Blood Cells (WBC)
Mean Platelet Volume (MPV)	

Hemoglobin Program



- Lyophilized, human whole blood based
- Includes HbA_{1C} and HbA₂
- 12-month cycle
- Submit results for the appropriate sample every month
- **Reference Method Values** provided for Hemoglobin A_{1C}

Analytes

Hemoglobin A _{1C}	Hemoglobin (Total Glycated)
Hemoglobin A ₂	

HIV/Hepatitis Program



- Liquid, human source material
- Suitable for serological tests used to detect key HIV and Hepatitis viral antibodies and antigens
- Qualitative and quantitative results reporting available
- Submit results and view reports each month through EQAS Online

Analytes

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

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

Immunoassay (Monthly) Program



- Lyophilized, human serum based
- Comprehensive menu of the most commonly tested immunoassay analytes
- 12-month cycle
- Submit results for the appropriate sample every month

Analytes

11-Deoxycortisol	DHEA	Phenobarbital	Thyroid Stimulating Hormone (TSH)
17- α -OH-Progesterone	DHEA Sulfate	Phenytoin	Thyroxine Binding Globulin (TBG)
25-OH Vitamin D	Digoxin	Plasma Renin Activity	Thyroxine, Free (FT ₄)
β -2 Microglobulin	Estradiol	Progesterone	Thyroxine (T4), Total
ACE	Estriol, Free (UE ₃)	Prolactin	Transferrin
ACTH	Ferritin	PSA (Free)	Triiodothyronine, Free (FT ₃)
Aldosterone	Folate	PSA (Free/Total Ratio)	Triiodothyronine (T3), Total
Alphafetoprotein (AFP)	Follicle Stimulating Hormone (FSH)	PSA (Total)	Valproic Acid
Androstenedione	Gastrin	Renin	Vitamin B ₁₂
C-Peptide	hCG	S100 Protein	Pilot Analytes
CA 15-3	hGH	Sex Hormone Binding Globulin (SHBG)	Calcitonin
CA 19-9	Immunoglobulin E (IgE)	T3 Uptake	Fructosamine
CA 27.29	Insulin	T-Uptake	Testosterone (Free)
CA 125	Intact PTH	Testosterone	
Carbamazepine	LH	Theophylline	
Carcinoembryonic Antigen (CEA)	NSE	Thyroglobulin (Tg)	
Cortisol			

Lipids Program



- Liquid, human serum based
- Comprehensive menu of lipid profile analytes
- 12-month cycle
- Submit results for the appropriate sample every month

Analytes

Apolipoprotein A-1	Pilot Analyte
Apolipoprotein B	Cholesterol (non-HDL)
Cholesterol (HDL)	
Cholesterol (LDL)	
Cholesterol (Total)	
Lipoprotein (a)	
Triglycerides	

Serum Proteins Program



- Lyophilized, human serum based
- Comprehensive menu for external assessment of serum protein assays
- 12-month cycle
- Submit results for the appropriate sample every month

Analytes

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

C-Reactive Protein (CRP)
 Ceruloplasmin
 Complement C3
 Complement C4
 Haptoglobin
 Immunoglobulin A (IgA)

Immunoglobulin E (IgE)
 Immunoglobulin G (IgG)
 Immunoglobulin M (IgM)
 Kappa Light Chain
 Lambda Light Chain
 Prealbumin

Protein (Total)
 Rheumatoid Factor (RF)
 Transferrin
Pilot Analyte
 Anti-DNase B

Syphilis/Chagas Program



- Liquid, human source material
- Suitable for serological tests detecting *Treponema pallidum*, Non-Treponemal and *Trypanosoma cruzi* antibodies
- 12-month cycle
- Qualitative and quantitative results reporting available
- Submit results and view reports each month through EQAS Online

Analytes

Non-Treponemal Antibodies
Treponema pallidum IgG

Treponema pallidum Antibodies
Trypanosoma cruzi Antibodies

Therapeutic Drug Monitoring Program



- Lyophilized, human serum based
- Analytes targeted at therapeutic, sub-therapeutic and toxic levels
- 12-month cycle
- Submit results for the appropriate sample every month
- **Weighed-in Values** provided for select analytes

Analytes

Acetaminophen
 Amikacin
 Amitriptyline
 Caffeine
 Carbamazepine
 Cyclosporine
 Digoxin
 Ethosuximide
 Gentamicin
 Lithium
 Methotrexate

Nortriptyline
 Phenobarbital
 Phenytoin
 Primidone
 Salicylate
 Theophylline
 Tobramycin
 Tricyclic Antidepressants (TCA) Screen
 Valproic Acid
 Vancomycin

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

ToRCH/EBV/MuMZ Program



- Liquid, human source material
- Suitable for serological tests used to detect pediatric diseases and congenital infections
- Qualitative and quantitative results reporting available
- Submit results and view reports each month through EQAS Online

Analytes

Cytomegalovirus (CMV) IgG	Measles Virus (Rubeola) IgG
Cytomegalovirus (CMV) IgM	Mumps Virus IgG
Epstein-Barr Virus (EBNA) IgG	Rubella Virus IgG
Epstein-Barr Virus (VCA) IgG	Rubella Virus IgM
Epstein-Barr Virus (VCA) IgM	<i>Toxoplasma gondii</i> (Toxo) IgG
Herpes Simplex Virus (HSV-1) IgG	<i>Toxoplasma gondii</i> (Toxo) IgM
Herpes Simplex Virus (HSV-1/2) IgG	Varicella Zoster Virus (VZV) IgG
Herpes Simplex Virus (HSV-2) IgG	

Urinalysis Program



- Liquid matrix
- Suitable for manual and automated dipstick analysis in both clinical and point-of-care testing (POCT)
- Submit qualitative results and view reports each month through EQAS Online

Analytes

Albumin	Nitrite
Albumin-to-Creatinine Ratio	pH
Bilirubin	Pregnancy (hCG)
Blood/Hemoglobin	Protein (Total)
Creatinine	Protein-to-Creatinine Ratio
Glucose	Specific Gravity
Ketones	Urobilinogen
Leukocytes	

Urine Chemistry Program



- Lyophilized, human urine based
- Matrix appropriate samples include basic and specialty urine chemistry analytes
- 12-month cycle
- Submit results for the appropriate sample every month

Analytes

5-Hydroxyindoleacetic Acid (5-HIAA)	Epinephrine	Normetanephrine	Urea Nitrogen
Calcium (Total)	Glucose	Osmolality	Uric Acid
Chloride	Homovanillic Acid (HVA)	Phosphorus	Vanillylmandelic Acid (VMA)
Cortisol (Free)	Magnesium	Potassium	Pilot Analyte
Creatinine	Metanephrine	Protein (Total)	Urea
Dopamine	Microalbumin (Albumin)	Sodium	
	Norepinephrine		

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

Ordering Information

Cat #	Description		
Blood Gas Program*			
BC31	12 x 2.5 mL	
Blood Typing Program*			
		<i>Sample</i>	<i>Donor</i>
12000837	Shipment A	3 x 4 mL + 1 x 2 mL
12000838	Shipment B	3 x 4 mL + 1 x 2 mL
12000839	Shipment C	3 x 4 mL + 1 x 2 mL
Cardiac Markers Program*			
BC39	12 x 1.5 mL	
Clinical Chemistry (Monthly) Program			
BC50	12 x 5 mL	
Coagulation Program*			
BC34	12 x 1 mL	
Ethanol/Ammonia Program*			
BC35	12 x 3 mL	
Hematology Program*			
BC90A	Shipment A	3 x 2 mL
BC90B	Shipment B	3 x 2 mL
BC90C	Shipment C	3 x 2 mL
BC90D	Shipment D	3 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.

Cat #	Description	
Hemoglobin Program		
BC80	12 x 0.5 mL
HIV/Hepatitis Program		
12000815	12 x 2 mL
Immunoassay (Monthly) Program		
BC75	12 x 5 mL
Lipids Program*		
BC47	12 x 3 mL
Serum Proteins Program		
BC23	12 x 1 mL
Syphilis/Chagas Program		
12000833	12 x 1.5 mL
Therapeutic Drug Monitoring Program		
BC10	12 x 5 mL
ToRCH/EBV/MuMZ Program		
12000825	12 x 2 mL
Urinalysis Program		
12000810	12 x 12 mL
Urine Chemistry Program		
BC45	12 x 10 mL

EQAS

An independent, external assessment of performance in comparison to your peers.

QCNet.com/eqas

Independent QC

Ongoing, proactive, unbiased daily QC that helps identify errors as they occur or begin to trend.

QCNet.com/independentqc

Unity

QC Data Management tools that help you create a strategy to reduce risk and streamline QC workflow.

QCNet.com/datamanagement

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

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/diagnostics 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 6169 8500 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