

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 assess 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.



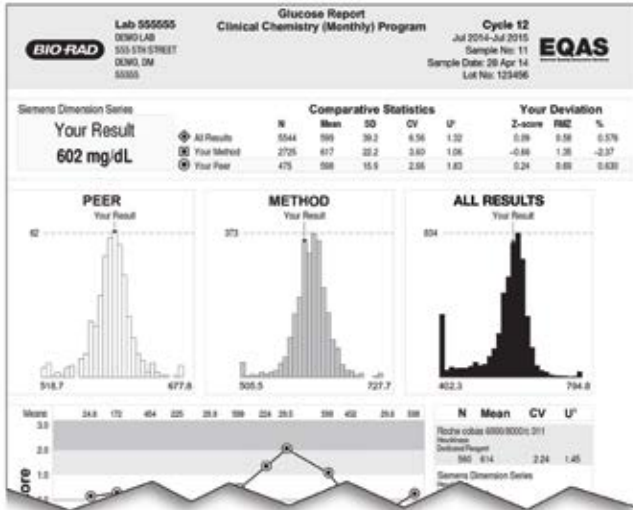
Submitting Results via EQAS Online/Mobile

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.

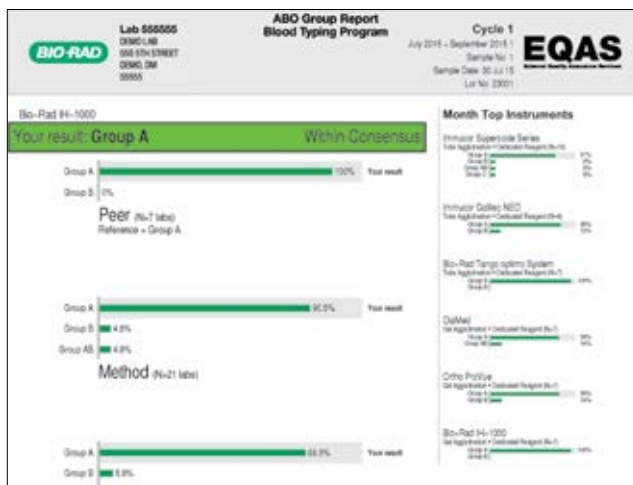
Quantitative Reports



Sample Report

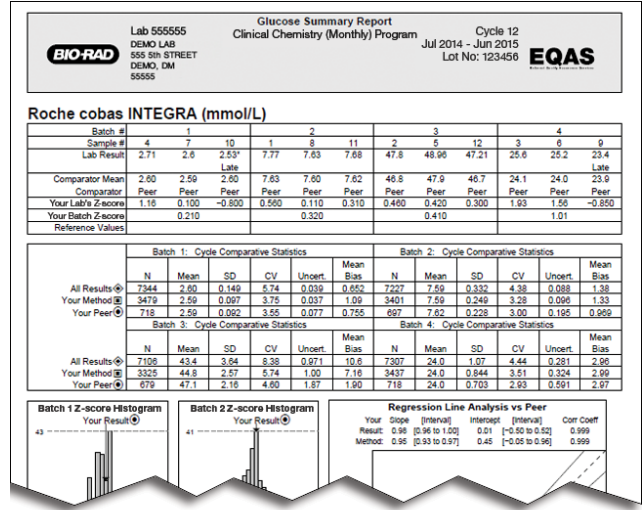
Detailed individual statistical analyses and peer group comparisons are provided.

Qualitative Report



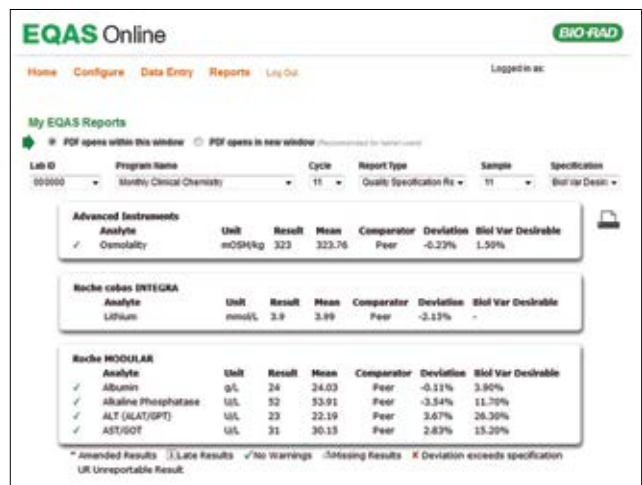
Sample Report

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 RiliBÄK), 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	<i>Pilot Analyte</i>
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	

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

Coagulation Program



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

Analytes

Antithrombin III (ATIII)	Protein C
APTT	Protein S
D-dimer	PT
Fibrinogen	Thrombin Time (TT)
INR	

Ethanol/Ammonia Program



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

Analytes

Ammonia
Ethanol

*Aspects of the program are subcontracted from an approved supplier.
Refer to myelinserts.com or the package insert of currently available lots for specific analyte and stability claims.

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	

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

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	<i>Pilot Analyte</i>
Apolipoprotein B	Cholesterol (non-HDL)
Cholesterol (HDL)	
Cholesterol (LDL)	
Cholesterol (Total)	
Lipoprotein (a)	
Triglycerides	

*Aspects of the program are subcontracted from an approved supplier.

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

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

Syphilis Program



- Liquid, human source material
- Suitable for serological tests detecting treponemal and non-treponemal 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 IgM
Treponema pallidum 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
 Cytomegalovirus (CMV) IgM
 Epstein-Barr Virus (EBNA) IgG
 Epstein-Barr Virus (VCA) IgG
 Epstein-Barr Virus (VCA) IgM
 Herpes Simplex Virus (HSV-1) IgG
 Herpes Simplex Virus (HSV-1/2) IgG
 Herpes Simplex Virus (HSV-2) IgG

Measles Virus (Rubeola) IgG
 Mumps Virus IgG
 Rubella Virus IgG
 Rubella Virus IgM
Toxoplasma gondii (Toxo) IgG
Toxoplasma gondii (Toxo) IgM
 Varicella Zoster Virus (VZV) IgG

Urinalysis Program



- Liquid
- 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
 Albumin-to-Creatinine Ratio
 Bilirubin
 Blood/Hemoglobin
 Creatinine
 Glucose
 Ketones
 Leukocytes

Nitrite
 pH
 Pregnancy (hCG)
 Protein (Total)
 Protein-to-Creatinine Ratio
 Specific Gravity
 Urobilinogen

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)
 Aldosterone
 Calcium (Total)
 Chloride
 Cortisol (Free)
 Creatinine

Dopamine
 Epinephrine
 Glucose
 Homovanillic Acid (HVA)
 Hydroxyproline (Total)
 Magnesium
 Metanephrine

Microalbumin (Albumin)
 Norepinephrine
 Normetanephrine
 Osmolality
 Phosphorus
 Potassium
 Protein (Total)

Sodium
 Urea Nitrogen
 Uric Acid
 Vanillylmandelic Acid (VMA)
Pilot Analyte
 Urea

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

EQAS Programs

- Analyte present
- ▶ Pilot analyte

	4	5	6	7	8	9											
	Blood Gas	Blood Typing	Cardiac Markers	Clin. Chemistry (Monthly)	Coagulation	Ethanol/Ammonia	Hematology	Hemoglobin	HIV/Hepatitis	Immunoassay (Monthly)	Lipids	Serum Proteins	Syphilis	TDM	ToRCH/EBV/MeuNZ	Urinalysis	Urine Chemistry
5-Hydroxyindoleacetic Acid (HIAA)																	●
11-Deoxycortisol										●							
17-α-OH-Progesterone										●							
25-OH Vitamin D										●							
α-1 Acid Glycoprotein												●					
α-1 Antitrypsin												●					
α-2 Macroglobulin												●					
β-2 Microglobulin										●		●					
ABO Group		●															
Acetaminophen														●			
Acid Phosphatase (Total)				●													
ACTH										●							
Alanine Aminotransferase (ALT/SGPT)				●													
Albumin				●								●					●
Albumin-to-Creatinine Ratio																●	
Aldosterone										●							●
Alkaline Phosphatase (ALP)				●													
Alphafetoprotein (AFP)										●							
Amikacin																●	
Amitriptyline																●	
Ammonia						●											
Amylase (Pancreatic)				●													
Amylase (Total)				●													
Androstenedione										●							
Angiotensin Converting Enzyme (ACE)										●							
Anti-HAV									●								
Anti-HBc									●								
Anti-HBe									●								
Anti-HBs									●								
Anti-HCV									●								
Anti-HIV-1									●								
Anti-HIV-1/2									●								
Anti-HIV-2									●								
Anti-HTLV-I									●								
Antibody Identification		●															
Antibody Screen		▶															
Antigen Typing		▶															
Antistreptolysin O (ASO)												●					
Antithrombin III (ATIII)					●												
Apolipoprotein A-1											●						
Apolipoprotein B											●						
APTT					●												
Aspartate Aminotransferase (AST/SGOT)				●													
Bile Acid				▶													
Bilirubin																	●
Bilirubin (Direct)				●													
Bilirubin (Indirect)				▶													
Bilirubin (Total)				●													
Blood/Hemoglobin																	●
BNP			●														
C-Peptide										●							
C-Reactive Protein (CRP)												●					
CA 15-3										●							
CA 19-9										●							
CA 27.29										●							
CA 125										●							
Caffeine																●	
Calcium (Ionized)	●			●													
Calcium (Total)				●													●

EQAS Programs

- Analyte present
- ▶ Pilot analyte

	4	5	6	7	8	9											
	Blood Gas	Blood Typing	Cardiac Markers	Clin. Chemistry (Monthly)	Coagulation	Ethanol/Ammonia	Hematology	Hemoglobin	HIV/Hepatitis	Immunoassay (Monthly)	Lipids	Serum Proteins	Syphilis	TDM	ToRCH/EBV/nuMZ	Urinalysis	Urine Chemistry
Carbamazepine																	
Carbon Dioxide (CO ₂)				●													
Carcinoembryonic Antigen (CEA)										●							
Ceruloplasmin												●					
Chloride	●			●													●
Cholesterol (HDL)				●							●						
Cholesterol (LDL)				●							●						
Cholesterol (non-HDL)											▶						
Cholesterol (Total)				●							●						
Cholinesterase				●													
CK-MB			●														
Compatibility (Cross-match)		●															
Complement C3												●					
Complement C4												●					
Copper				●													
Cortisol				●						●							
Cortisol (Free)																	●
Creatine Kinase (CK)				●													
Creatinine				●												●	●
Cyclosporine															●		
Cytomegalovirus (CMV) IgG																●	
Cytomegalovirus (CMV) IgM																●	
DAT (Direct Antiglobulin Test)		●															
D-dimer			●		●												
DHEA										●							
DHEA Sulfate										●							
Digoxin										●							
Dopamine																	●
Epinephrine																	●
Epstein-Barr Virus (EBNA) IgG																●	
Epstein-Barr Virus (VCA) IgG																●	
Epstein-Barr Virus (VCA) IgM																●	
Estradiol										●							
Estriol, Free (UE ₃)										●							
Ethanol						●											
Ethosuximide															●		
Ferritin										●							
Fibrinogen					●												
Folate										●							
Follicle Stimulating Hormone (FSH)										●							
Gamma Glutamyltransferase (GGT)				●													
Gastrin										●							
Gentamicin															●		
Glucose	●			●												●	●
Haptoglobin												●					
HAV IgG									●								
HAV IgM									●								
HBc IgM									●								
HBeAg									●								
HBSAg									●								
hCG										●							
Hematocrit (HCT)							●										
Hemoglobin A1C							●										
Hemoglobin A ₂							●										
Hemoglobin							●										
Hemoglobin (Total Glycated)							●										
Herpes Simplex Virus (HSV-1) IgG																●	
Herpes Simplex Virus (HSV-1/2) IgG																●	
Herpes Simplex Virus (HSV-2) IgG																●	

EQAS Programs

- Analyte present
- Pilot analyte

	4		5		6		7	8		9							
	Blood Gas	Blood Typing	Cardiac Markers	Clin. Chemistry (Monthly)	Coagulation	Ethanol/Ammonia	Hematology	Hemoglobin	HIV/Hepatitis	Immunoassay (Monthly)	Lipids	Serum Proteins	Syphilis	TDM	ToRCH/EBV/nuMZ	Urinalysis	Urine Chemistry
hGH																	
HIV-1 Ag									●								
Homocysteine			●														
Homovanillic Acid (HVA)																	●
hs-CRP			●														
Hydroxyproline (Total)																	●
IgA												●					
IgE												●					
IgG												●					
IgM												●					
INR					●												
Insulin												●					
Intact PTH												●					
Iron					●												
Iron Binding Capacity, Total (TIBC)					●												
Iron Binding Capacity, Unsaturated (UIBC)					●												
Kappa Light Chain												●					
Ketones																	●
Lactate (Lactic Acid)	●			●													
Lactate Dehydrogenase (LDH)				●													
Lambda Light Chain												●					
Leukocytes																	●
LH												●					
Lipase				●													
Lipoprotein (a)												●					
Lithium				●										●			
Magnesium	●			●													●
Mean Corpuscular Hemoglobin (MCH)							●										
Mean Corpuscular Hemoglobin Concentration (MCHC)							●										
Mean Corpuscular Volume (MCV)							●										
Mean Platelet Volume (MPV)							●										
Measles Virus (Rubeola) IgG																●	
Metanephrine																	●
Methotrexate														●			
Microalbumin																	●
Mumps Virus IgG																●	
Myoglobin			●														
Nitrite																	●
Non-Treponemal antibodies													●				
Norepinephrine																	●
Normetanephrine																	●
Nortriptyline														●			
NSE												●					
NT-proBNP			●														
Osmolality				●													●
pCO ₂	●																
pH	●																●
pO ₂	●																
Phenobarbital												●		●			
Phenytoin												●		●			
Phosphorus				●													●
Plasma Renin Activity (PRA)												●					
Platelets (PLT)							●										
Potassium	●			●													●
Pregnancy (hCG)																	●
Prealbumin												●					
Primidone														●			
Progesterone										●							

EQAS Programs

- Analyte present
- ▶ Pilot analyte

	4	5	6	7	8	9											
	Blood Gas	Blood Typing	Cardiac Markers	Clin. Chemistry (Monthly)	Coagulation	Ethanol/Ammonia	Hematology	Hemoglobin	HIV/Hepatitis	Immunoassay (Monthly)	Lipids	Serum Proteins	Syphilis	TDM	ToRCH/EBV/nuMZ	Urinalysis	Urine Chemistry
Prolactin										●							
Protein C					●												
Protein S					●												
Protein (Total)				●								●					●
Protein-to-Creatinine Ratio																●	
PSA (Free)										●							
PSA (Free/Total Ratio)										●							
PSA (Total)										●							
PT					●												
Red Blood Cells (RBC)							●										
Red Blood Cell Distribution Width (RDW)							●										
Red Blood Cell Distribution Width (RDW-SD)							●										
Renin										●							
Rh(D)		●															
Rheumatoid Factor (RF)												●					
Rubella Virus IgG																●	
Rubella Virus IgM																●	
S100 Protein										●							
Salicylate													●				
Sex Hormone Binding Globulin (SHBG)										●							
Sodium	●			●													●
Specific Gravity																●	
T3 Uptake										●							
T-Uptake										●							
Testosterone										●							
Testosterone (Free)										▶							
Theophylline										●			●				
Thrombin Time (TT)					●												
Thyroglobulin (Tg)										▶							
Thyroid Stimulating Hormone (TSH)				●						●							
Thyroxine Binding Globulin (TBG)										●							
Thyroxine, Free (T ₄)				●						●							
Thyroxine (T ₄), Total				●						●							
Tobramycin																●	
<i>Toxoplasma gondii</i> (Toxo) IgG																●	
<i>Toxoplasma gondii</i> (Toxo) IgM																●	
Transferrin				▶						●		●					
<i>Treponema pallidum</i> IgG												●					
<i>Treponema pallidum</i> IgM												●					
<i>Treponema pallidum</i> Total Antibody												●					
Tricyclic Antidepressants (TCA) Screen														●			
Triglycerides				●							●						
Triiodothyronine, Free (T ₃)				●						●							
Triiodothyronine (T ₃), Total				●						●							
Troponin I			●														
Troponin T			●														
Urea				●													▶
Urea Nitrogen				●													●
Uric Acid				●													●
Urobilinogen																●	
Valproic Acid										●				●			
Vancomycin														●			
Vanillylmandelic Acid (VMA)																	●
Varicella Zoster Virus (VZV) IgG																●	
Vitamin B ₁₂										●							
White Blood Cells (WBC)							●										
Zinc				●													

Ordering Information

Cat #	Description	Sample	Donor
Blood Gas Program*			
BC31	12 x 2.5 mL	
Blood Typing Program*			
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 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.



Independent QC

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



Unity

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

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

Web site www.bio-rad.com/qualitycontrol USA 1 800 224 6723 Australia 61 2 9914 2800 Austria 43 1 877 8901 Belgium 32 03 710 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