



EQAS

External Quality Assurance Services





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.

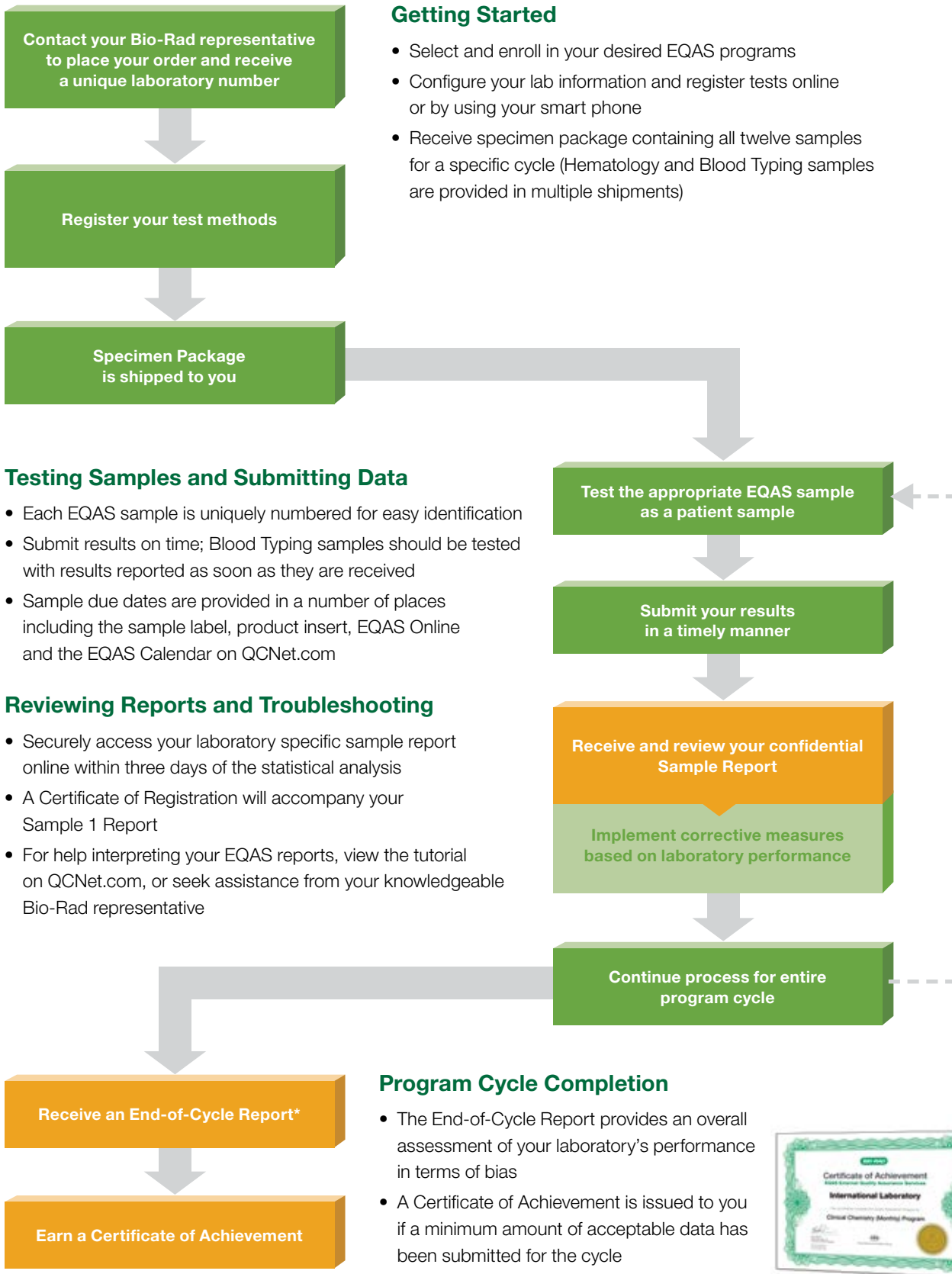
External quality assessment programs are accepted around the world as invaluable tools used by laboratories to periodically assess the performance of their test systems. Results are objectively compared to other laboratories using the same methodologies, instruments and reagents. When used in conjunction 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 35 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



*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.



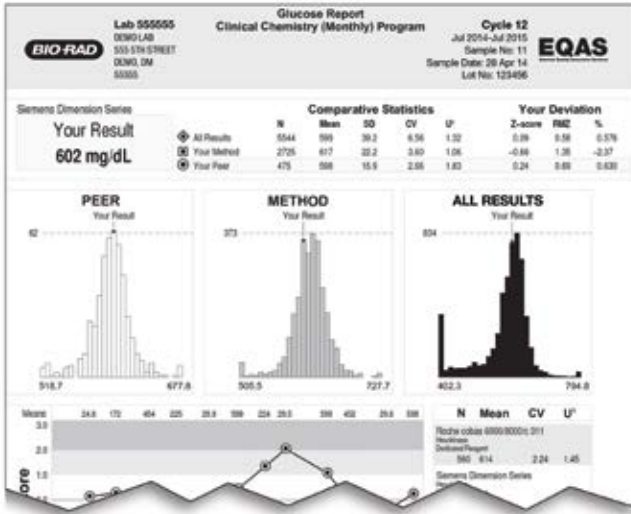
Submitting Results via EQAS Online/Mobile

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

EQAS Reports

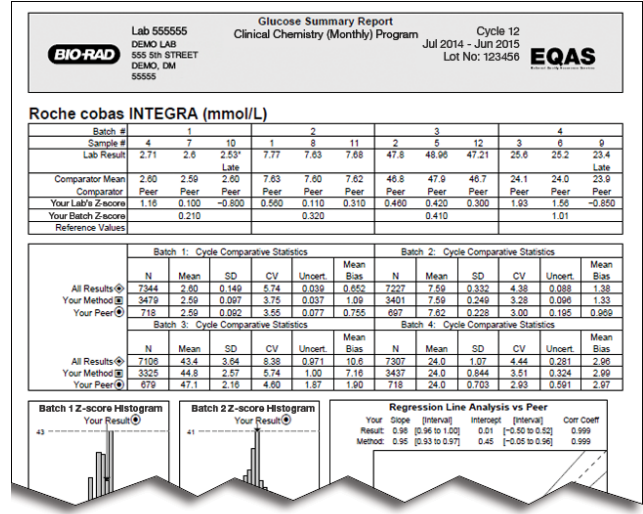
Participants receive twelve sample reports and one end-of-cycle report over each program cycle. These detailed reports are presented in an easy-to-read graphic format. Sample reports are available online within three days of the statistical analysis. Additional special reports (Quality Specification,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

Quantitative Reports



Sample Report

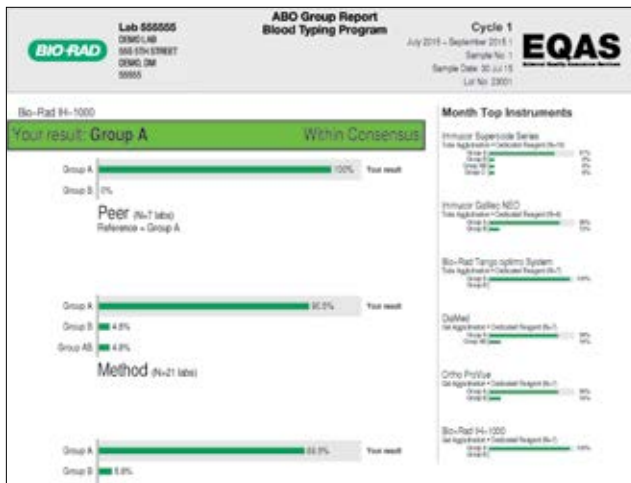
Detailed individual statistical analyses and peer group comparisons are provided.



End-of-Cycle Report

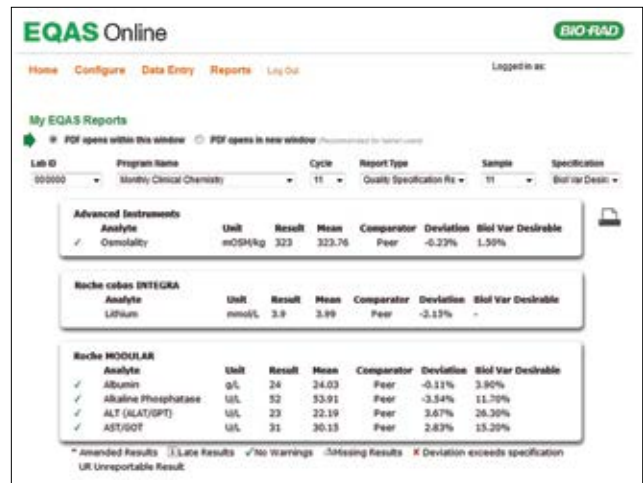
An overall assessment of laboratory performance over the entire cycle in terms of bias among participating laboratories is presented.

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.



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
- 12 x 2.5 mL

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
- 3 x 4 mL Sample and 1 x 2 mL Donor (each shipment)

Parameters

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

Cardiac Markers Program

- Liquid, human plasma based
- Comprehensive menu of important cardiac assessment markers, plus D-dimer
- 12-month cycle
- 12 x 1.5 mL

Analytes

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



Refer to myinserts.com or the package insert for current analyte 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
- 12 x 5 mL

Analytes

Acid Phosphatase (Total)	Carbon Dioxide (CO ₂)	Iron	Protein (Total)
Alanine Aminotransferase (ALT/SGPT)	Chloride	Iron Binding Capacity, Total (TIBC)	Sodium
Albumin	Cholesterol (HDL)	Iron Binding Capacity, Unsaturated (UIBC)	Thyroid Stimulating Hormone (TSH)
Alkaline Phosphatase (ALP)	Cholesterol (LDL)	Lactate (Lactic Acid)	Thyroxine (FT ₄), Free
Amylase (Pancreatic)	Cholesterol (Total)	Lactate Dehydrogenase (LDH)	Thyroxine (T4), Total
Amylase (Total)	Cholinesterase	Lipase	Triglycerides
Aspartate Aminotransferase (AST/SGOT)	Copper	Lithium	Triiodothyronine (FT ₃), Free
Bilirubin (Direct)	Cortisol	Magnesium	Triiodothyronine (T3), Total
Bilirubin (Total)	Creatine Kinase (CK)	Osmolality	Urea
Calcium (Ionized)	Creatinine	Phosphorus	Urea Nitrogen
Calcium (Total)	Gamma Glutamyl Transferase (GGT)	Potassium	Uric Acid
	Glucose		Zinc

Coagulation Program

- Lyophilized, human plasma based
- 9 parameters for external assessment of laboratory coagulation test performance
- 12-month cycle
- 12 x 1 mL

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
- 12 x 3 mL

Analytes

Ammonia
Ethanol



Refer to myinserts.com or the package insert for current analyte 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
- 3 x 2 mL (each shipment)

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 HbA1C and HbA₂
- 12-month cycle
- Submit results for the appropriate sample every month
- Reference Method Values provided for Hemoglobin A1C
- 12 x 0.5 mL

Analytes

Hemoglobin A1C
 Hemoglobin A₂
 Hemoglobin (Total Glycated)

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



Refer to myinserts.com or the package insert for current analyte 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
- 12 x 5 mL

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 (T ₄), Total
Aldosterone	Ferritin	Prolactin	Transferrin
Alphafetoprotein (AFP)	Folate	PSA (Free)	Triiodothyronine, Free (FT ₃)
Androstenedione	Follicle Stimulating Hormone (FSH)	PSA (Free/Total Ratio)	Triiodothyronine (T ₃), Total
C-Peptide	Gastrin	PSA (Total)	Valproic Acid
CA 15-3	hCG	Renin	Vitamin B ₁₂
CA 19-9	hGH	S100 Protein	<i>Pilot Analyte</i>
CA 27.29	Immunoglobulin E (IgE)	Sex Hormone Binding Globulin (SHBG)	Thyroglobulin (Tg)
CA 125	Insulin	T3 Uptake	
Carbamazepine	Intact PTH	T-Uptake	

Lipids Program

- Liquid, human serum based
- Comprehensive menu of lipid profile analytes
- 12-month cycle
- Submit results for the appropriate sample every month
- 12 x 3 mL

Analytes

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



Refer to myinserts.com or the package insert for current analyte 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
- 12 x 1 mL

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
- 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

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 at end of cycle
- 12 x 5 mL

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 for current analyte claims.

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

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
- 12 x 12 mL

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
- 12 x 10 mL

Analytes

5-Hydroxyindoleacetic Acid
(5-HIAA)
Aldosterone
Calcium (Total)
Chloride
Cortisol (Free)

Creatinine
Dopamine
Epinephrine
Glucose
Homovanillic Acid (HVA)
Hydroxyproline (Total)

Magnesium
Metanephrine
Microalbumin
Norepinephrine
Normetanephrine
Osmolality

Phosphorus
Potassium
Protein (Total)
Sodium
Urea Nitrogen
Uric Acid
Vanillylmandelic Acid (VMA)



Refer to myinserts.com or the package insert for current analyte claims.

EQAS Analyte Chart

See page number

- Analyte present
- ▶ Pilot analyte

	4	5	6	7	8	9											
	Blood Gas	Blood Typing	Cardiac Markers	Clinical Chemistry (Monthly)	Coagulation	Ethanol/Ammonia	Hematology	Hemoglobin	HIV/Hepatitis	Immunoassay (Monthly)	Lipids	Serum Proteins	Syphilis	Therapeutic Drug Monitoring	ToRCH/EBV/MUMZ	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)										●							
Antistreptolysin O (ASO)												●					
Antithrombin III (ATIII)					●												
Anti-HAV									●								
HAV IgG									●								
HAV IgM									●								
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		▶															
Apolipoprotein A-1												●					
Apolipoprotein B												●					
APTT					●												
Aspartate Aminotransferase (AST/SGOT)				●													
Bilirubin																	●
Bilirubin (Direct)				●													
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)	●			●													

EQAS Analyte Chart

- Analyte present
- ▶ Pilot analyte

	4	5	6	7	8	9											
	Blood Gas	Blood Typing	Cardiac Markers	Clinical Chemistry (Monthly)	Coagulation	Ethanol/Ammonia	Hematology	Hemoglobin	HIV/Hepatitis	Immunoassay (Monthly)	Lipids	Serum Proteins	Syphilis	Therapeutic Drug Monitoring	ToRCH/EBV/MUMZ	Urinalysis	Urine Chemistry
Calcium (Total)				●													●
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														●			
DAT (Direct Antiglobulin Test)		●															
Cytomegalovirus (CMV) IgG																●	
Cytomegalovirus (CMV) IgM																●	
D-dimer			●		●												
DHEA										●							
DHEA Sulfate										●							
Digoxin										●				●			
Dopamine																	●
Epinephrine																	●
Epstein-Barr Virus (EBNA) IgG																●	
Epstein-Barr Virus (VCA) IgG																●	
Epstein-Barr Virus (VCA) IgM																●	
Estradiol										●							
Estradiol, Free (UE ₃)										●							
Ethanol						●											
Ethosuximide														●			
Ferritin										●							
Fibrinogen					●												
Folate										●							
Follicle Stimulating Hormone (FSH)										●							
Gamma Glutamyltransferase (GGT)				●													
Gastrin										●							
Gentamicin														●			
Glucose	●			●												●	●
Haptoglobin												●					
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 Analyte Chart

See page number

- Analyte present
- Pilot analyte

	4	5	6	7	8	9											
	Blood Gas	Blood Typing	Cardiac Markers	Clinical Chemistry (Monthly)	Coagulation	Ethanol/Ammonia	Hematology	Hemoglobin	HIV/Hepatitis	Immunoassay (Monthly)	Lipids	Serum Proteins	Syphilis	Therapeutic Drug Monitoring	ToRCH/EBV/MUMZ	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														●			

EQAS Analyte Chart

See page number

- Analyte present
- ▶ Pilot analyte

	4	5	6	7	8	9											
	Blood Gas	Blood Typing	Cardiac Markers	Clinical Chemistry (Monthly)	Coagulation	Ethanol/Ammonia	Hematology	Hemoglobin	HIV/Hepatitis	Immunoassay (Monthly)	Lipids	Serum Proteins	Syphilis	Therapeutic Drug Monitoring	ToRCH/EBV/MUMZ	Urinalysis	Urine Chemistry
Progesterone										●							
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										●							
Theophylline										●				●			
Thrombin Time (TT)					●												
Thyroid Stimulating Hormone (TSH)				●						●							
Thyroglobulin										▶							
Thyroxine Binding Globulin (TBG)										●							
Thyroxine, Free (FT ₄)				●						●							
Thyroxine (T ₄), Total				●						●							
Tobramycin														●			
Transferrin										●		●					
<i>Toxoplasma gondii</i> (Toxo) IgG															●		
<i>Toxoplasma gondii</i> (Toxo) IgM															●		
<i>Treponema pallidum</i> IgG													●				
<i>Treponema pallidum</i> IgM													●				
<i>Treponema pallidum</i> Total Antibody													●				
Tricyclic Antidepressants (TCA) Screen														●			
Triglycerides				●							●						
Triiodothyronine, Free (FT 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 #	Configuration
Blood Gas Program*	
BC31	12 x 2.5 mL
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
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 myeinserts.com or the package insert for current analyte claims.

Cat #	Configuration
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



**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