



EQAS

External Quality Assurance Services

BIO-RAD



Participate in an internationally recognized quality assessment program

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 control, these external programs can give laboratories added confidence in their patient test results.

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

- Compare your results to a large international database with participants from over 90 countries
- Evaluate your results using comprehensive, easy-to-read reports
- Benefit from the support of experienced professionals
- eForms electronic data entry (optional)
- All 12 samples for each cycle are shipped at the same time
- Extensive analyte menus spanning normal and abnormal levels

Our EQAS programs are accredited, providing participant confidence in quality assurance.



Immunoassay (Monthly) Program

- 12 x 5 mL human serum based samples
- 1 sample is tested every month/12-month cycle

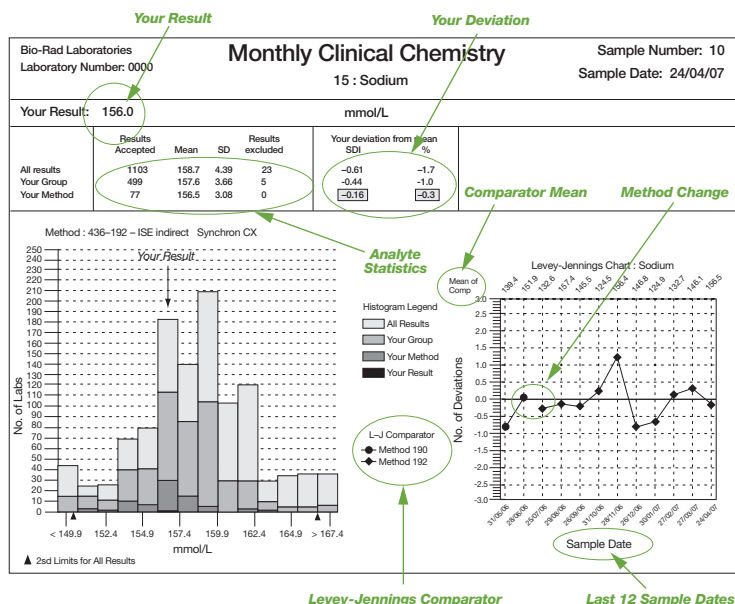
Analytes

17- α -OH-Progesterone
 β -2 Microglobulin
 Alphafetoprotein (AFP)
 CA 15-3
 CA 19-9
 CA 125
 Carbamazepine
 Carcinoembryonic Antigen (CEA)
 Chorionic Gonadotropin (hCG)
 Cortisol
 DHEA-Sulfate
 Digoxin

Estradiol
 Ferritin
 Folate
 Follicle Stimulating Hormone (FSH)
 Free Thyroxine (FT₄)
 Free Triiodothyronine (FT₃)
 Growth Hormone (hGH)
 Immunoglobulin E (IgE)
 Luteinizing Hormone (LH)
 Parathyroid Hormone (PTH), Intact
 Phenobarbital
 Phenytoin

Progesterone
 Prolactin
 Prostate Specific Antigen (PSA), Free
 Prostate Specific Antigen (PSA), Total
 Testosterone
 Theophylline
 Thyroid Stimulating Hormone (TSH)
 Thyroxine (T₄), Total
 Triiodothyronine (T₃), Total
 Valproic Acid
 Vitamin B₁₂

Refer to the package insert for current analyte claims



EQAS Reports

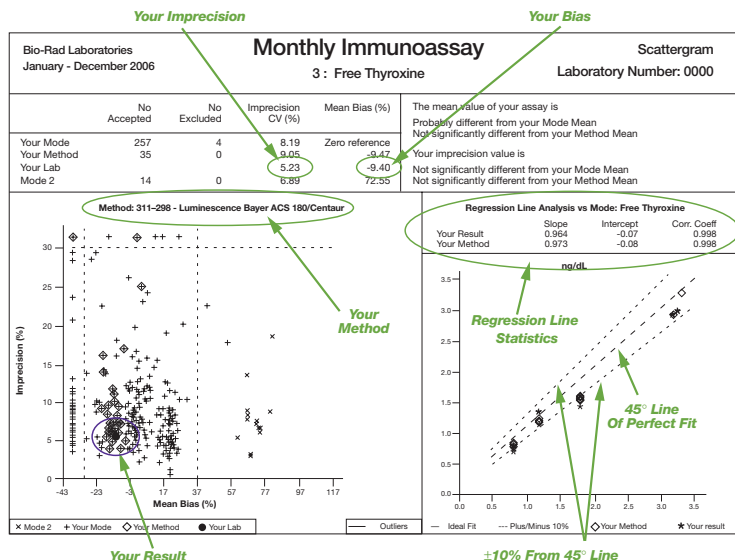
Participants receive twelve bi-weekly or monthly reports and one end-of-cycle report over each program cycle. These detailed reports are presented in an easy-to-read graphic format. All reports are available online within three days of the statistical analysis.

Bi-Weekly/Monthly 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 and imprecision among participating laboratories is presented.



Additional reports (available on request):

- Sub-Group Reports
- Instrument/Method Summary Report



Clinical Chemistry Program

- 12 x 5 mL human serum based samples
- 1 sample is tested every 2 weeks / 6-month cycle

Analytes

α -Hydroxybutyrate Dehydrogenase (HBDH)
 γ -Glutamyl Transferase (GGT)
 Acid Phosphatase, Total
 Alanine Aminotransferase (ALT/SGPT)
 Albumin
 Alkaline Phosphatase (ALP)
 Aluminum
 Amylase, Total
 Aspartate Aminotransferase (AST/SGOT)
 Bicarbonate (CO₂)
 Bilirubin, Total
 Calcium, Total

Chloride
 Cholesterol, Total
 Copper
 Cortisol
 Creatine Kinase (CK)
 Creatinine
 Glucose
 HDL Cholesterol
 Iron
 Iron Binding Capacity, Total (TIBC)
 Lactate Dehydrogenase (LDH)
 Lithium

Magnesium
 Osmolality
 Phosphate
 Potassium
 Prostatic Acid Phosphatase (PAP)
 Protein, Total
 Sodium
 Triglycerides
 Urea
 Uric Acid
 Zinc

Refer to the package insert for current analyte claims



Clinical Chemistry (Monthly) Program

- 12 x 5 mL human serum based samples
- 1 sample is tested every month / 12-month cycle

Analytes

γ -Glutamyl Transferase (GGT)
 Alanine Aminotransferase (ALT/SGPT)
 Albumin
 Alkaline Phosphatase (ALP)
 Amylase, Pancreatic
 Amylase, Total
 Aspartate Aminotransferase (AST/SGOT)
 Bicarbonate (CO₂)
 Bilirubin, Direct
 Bilirubin, Total
 Calcium, Ionized
 Calcium, Total

Chloride
 Cholesterol, Total
 Creatine Kinase (CK)
 Creatinine
 Glucose
 HDL Cholesterol
 Iron
 Iron Binding Capacity, Total (TIBC)
 Lactate Dehydrogenase (LDH)
 Lipase
 Lithium
 Magnesium

Osmolality
 Phosphate
 Potassium
 Protein, Total
 Sodium
 Thyroid Stimulating Hormone (TSH)
 Thyroxine (T₄), Total
 Triglycerides
 Triiodothyronine (T₃), Total
 Urea
 Uric Acid

Refer to the package insert for current analyte claims



Therapeutic Drug Monitoring Program

- 12 x 5 mL human serum based samples
- 1 sample is tested every 2 weeks / 6-month cycle

Analytes

Acetaminophen
 Amikacin
 Amitriptyline
 Caffeine
 Carbamazepine
 Cyclosporine
 Digoxin
 Ethosuximide

Gentamicin
 Lithium
 Methotrexate
 Nortriptyline
 Phenobarbital
 Phenytoin
 Primidone
 Quinidine

Salicylate
 Theophylline
 Tobramycin
 Tricyclic Screen
 Valproic Acid
 Vancomycin

Refer to the package insert for current analyte claims



Hemoglobin Program

- 12 x 0.5 mL human whole blood based samples
- 1 sample is tested every month / 12-month cycle
- End-of-cycle comparison to reference method values

Analytes

Hemoglobin A1c
Hemoglobin A2
Hemoglobin, Total Glycated (GHb)

Refer to the package insert for current analyte claims



Immunoassay Programs

- 12 x 5 mL (Programs 1 and 2) and 1 mL (Programs 3 and 4) human serum based samples
- 1 sample is tested every 2 weeks / 6-month cycle
- Each program may be purchased separately

Analytes

Program 1

Thyroid Function

Free Thyroxine (FT₄)
Free Triiodothyronine (FT₃)
T3 Uptake/T-Uptake
Thyroid Stimulating Hormone (TSH)
Thyroxine (T₄), Total
Thyroxine Binding Globulin (TBG)
Triiodothyronine (T₃), Total

Down Syndrome/Pregnancy

Alphafetoprotein (AFP)
Chorionic Gonadotropin (hCG)
Estril (E₃), Total
Unconjugated Estril (UE₃)

Iron Metabolism

Ferritin
Folate
Transferrin
Vitamin B₁₂

Program 2

Reproductive Endocrinology

Chorionic Gonadotropin (hCG)
DHEA-Sulfate
Estradiol
Follicle Stimulating Hormone (FSH)
Luteinizing Hormone (LH)
Progesterone
Prolactin
Sex Hormone Binding Globulin (SHBG)
Testosterone

General Endocrinology

17- α -OH-Progesterone
Aldosterone
Cortisol
Gastrin
Growth Hormone (hGH)
Insulin
Parathyroid Hormone (PTH), Intact

Others

Angiotensin Converting Enzyme (ACE)
Digoxin
Renin (Angiotensin 1)
Renin, Direct

Program 3

Specific Proteins

α -1 Acid Glycoprotein
 α -1 Antitrypsin
 α -2 Macroglobulin
 β -2 Microglobulin
Albumin
C-Reactive Protein (CRP)
Ceruloplasmin
Complement C3

Complement C4

Haptoglobin
Immunoglobulin A (IgA)
Immunoglobulin E (IgE)
Immunoglobulin G (IgG)
Immunoglobulin M (IgM)
Prealbumin
Protein, Total

Pathological Proteins

Antistreptolysin O (ASO)
Rheumatoid Factor (RF)

Program 4

Oncology

Alphafetoprotein (AFP)
CA 15-3
CA 19-9
CA 125

Carcinoembryonic Antigen (CEA)
Chorionic Gonadotropin (hCG)
Prostate Specific Antigen (PSA), Free
Prostate Specific Antigen (PSA), Free/Total Ratio

Prostate Specific Antigen (PSA), Total
Prostatic Acid Phosphatase (PAP)

Refer to the package insert for current analyte claims





QCNet™

EQAS participants have easy access to product information, product inserts and user documentation through QCNet™ (www.QCNet.com). This secure website allows registered participants access to upload their eForms data files and view or print their EQAS reports.

Sign up on QCNet™

In the interest of confidentiality, only registered users have access to their EQAS reports.

1. Type in www.QCNet.com.
The QCNet™ home page appears.
2. Click on Register.
3. Complete the New User Signup information.
4. Click on Submit.
5. A successful registration screen appears.
Click OK to proceed to login.

Register your EQAS Lab Number

1. Enter your QCNet™ Login ID and Password and click on Login.
2. Click on the EQAS tab or link.
3. Click on the EQAS Registration.
4. Enter your EQAS lab number and select your country from the drop-down menu.
5. Click on Register.

Please allow up to 2 business days to process your registration. Once your laboratory number has been verified by Bio-Rad, your EQAS reports for the current cycle will be posted.

Ordering Information

Cat #	Product Name	Quantity	Cat #	Product Name	Quantity
Clinical Chemistry Program			Immunoassay Programs		
BC36	Clinical Chemistry Program (17 analytes)	12 x 5 mL	BC21	Immunoassay Program 1	12 x 5 mL
BC37	Clinical Chemistry Program (35 analytes)	12 x 5 mL	BC22	Immunoassay Program 2	12 x 5 mL
CS36	Clinical Chemistry Program (17 analytes)	Report Only	BC23	Immunoassay Program 3	12 x 1 mL
CS37	Clinical Chemistry Program (35 analytes)	Report Only	BC24	Immunoassay Program 4	12 x 1 mL
QC37	Clinical Chemistry Program	12 x 5 mL (specimen pack only)	CS21	Immunoassay Program 1	Report Only
Clinical Chemistry (Monthly) Program			CS22	Immunoassay Program 2	Report Only
BC5L	Clinical Chemistry (Monthly) Program (10 analytes)	12 x 5 mL	CS23	Immunoassay Program 3	Report Only
BC50	Clinical Chemistry (Monthly) Program (35 analytes)	12 x 5 mL	CS24	Immunoassay Program 4	Report Only
CS50	Clinical Chemistry (Monthly) Program	Report Only	QC21	Immunoassay Program 1	12 x 5 mL (specimen pack only)
QC50	Clinical Chemistry (Monthly) Program	12 x 5 mL (specimen pack only)	QC22	Immunoassay Program 2	12 x 5 mL (specimen pack only)
Therapeutic Drug Monitoring Program			QC23	Immunoassay Program 3	12 x 1 mL (specimen pack only)
BC05	Therapeutic Drug Monitoring Program	12 x 5 mL (12 analytes)	QC24	Immunoassay Program 4	12 x 1 mL (specimen pack only)
BC10	Therapeutic Drug Monitoring Program	12 x 5 mL (22 analytes)	Immunoassay (Monthly) Program		
CS05	Therapeutic Drug Monitoring Program	Report Only	BC7L	Immunoassay (Monthly) Program	12 x 5 mL (8 analytes)
QC05	Therapeutic Drug Monitoring Program	12 x 5 mL (specimen pack only)	BC70	Immunoassay (Monthly) Program	12 x 5 mL (17 analytes)
Hemoglobin Program			BC75	Immunoassay (Monthly) Program	12 x 5 mL (35 analytes)
BC80	Hemoglobin Program	12 x 0.5 mL	CS70	Immunoassay (Monthly) Program	Report Only (17 analytes)
CS80	Hemoglobin Program	Report Only	CS75	Immunoassay (Monthly) Program	Report Only (35 analytes)
QC80	Hemoglobin Program	12 x 0.5 mL (specimen pack only)	QC75	Immunoassay (Monthly) Program	12 x 5 mL (specimen pack only)



**Bio-Rad
Laboratories**

**For further information, please contact your local Bio-Rad
office or learn more online at www.bio-rad.com/qualitycontrol**

**Clinical
Diagnostics Group**

Website www.bio-rad.com/qualitycontrol **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** 36-1-455-8800 **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)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** 7-495-721-14-04 **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