



EQAS

External Quality Assessment Suited for the Small Laboratory

- 12 samples per program cycle
- Evaluate and report sample results each month
- All samples provided in a single shipment
- Human serum based
- Select from a large menu of available analytes
- Laboratory specific reports in an easy-to-read format

Clinical Chemistry (Monthly) Program

Select Up to 10 Analytes

Analytes

Alanine Aminotransferase (ALT/SGPT)	Iron
Albumin	Iron Binding Capacity, Total (TIBC)
Alkaline Phosphatase (ALP)	Lactate Dehydrogenase (LDH)
Amylase, Pancreatic	Lipase
Amylase, Total	Lithium
Aspartate Aminotransferase (AST/SGOT)	Magnesium
Bicarbonate (CO ₂)	Osmolality
Bilirubin, Direct	Phosphate
Bilirubin, Total	Potassium
Calcium, Ionized	Protein, Total
Calcium, Total	Sodium
Chloride	Thyroid Stimulating Hormone (TSH)
Cholesterol, Total	Thyroxine (T ₄), Total
Creatine Kinase (CK)	Triglycerides
Creatinine	Triiodothyronine (T ₃), Total
γ-Glutamyl Transferase (GGT)	Urea
Glucose	Uric Acid
HDL Cholesterol	

Refer to the package insert for current analyte claims

Immunoassay (Monthly) Program

Select Up to 8 Analytes

Analytes

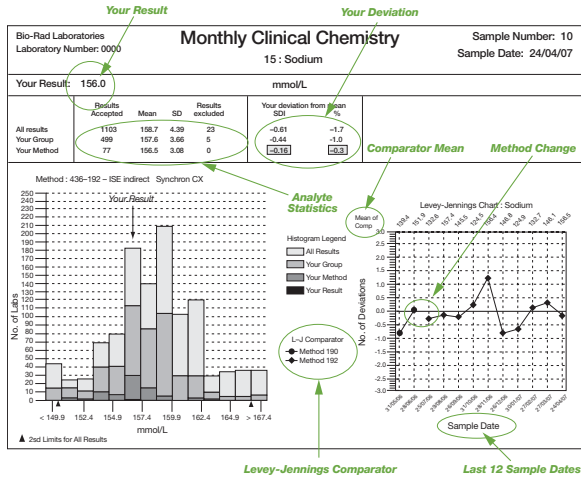
17-α-OH-Progesterone	Immunoglobulin E (IgE)
Alphafetoprotein (AFP)	Luteinizing Hormone (LH)
β-2 Microglobulin	Parathyroid Hormone (PTH), Intact
CA 15-3	Phenobarbital
CA 19-9	Phenytoin
CA 125	Progesterone
Carbamazepine	Prolactin
Carcinoembryonic Antigen (CEA)	Prostate Specific Antigen (PSA), Free
Chorionic Gonadotropin (hCG)	Prostate Specific Antigen (PSA), Total
Cortisol	Testosterone
DHEA-Sulfate	Theophylline
Digoxin	Thyroid Stimulating Hormone (TSH)
Estradiol	Thyroxine (T ₄), Total
Ferritin	Triiodothyronine (T ₃), Total
Folate	Valproic Acid
Follicle Stimulating Hormone (FSH)	Vitamin B ₁₂
Free Thyroxine (FT ₄)	
Free Triiodothyronine (FT ₃)	
Growth Hormone (hGH)	

EQAS Reports

Participants receive 12 monthly reports and one end-of-cycle report over each program cycle. All reports are available within three days of the statistical analysis.

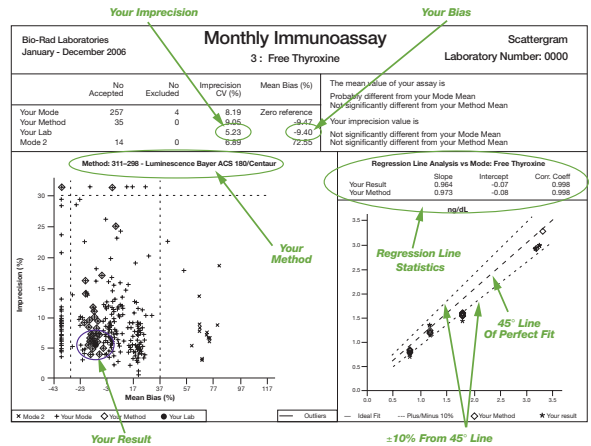
Monthly Report

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 upon request)

- Sub-Group Reports
- Instrument/Method Summary Report

Ordering Information

Cat #	EQAS Programs	Quantity
BC5L	Clinical Chemistry (Monthly) Program (10 analytes)	12 x 5 mL
BC7L	Immunoassay (Monthly) Program (8 analytes)	12 x 5 mL

Additional Programs Available



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