



EQAS Urine Chemistry Program

External Quality Assessment Suited for Your Laboratory

- Compare your laboratory's results to an international database of participants
- Evaluate and report sample results each month
- All 12 samples for each cycle are provided in a single shipment
- Human urine based samples
- Select from a large menu of basic and specialty urine chemistry analytes
- Laboratory specific reports in an easy-to-read format with online viewing
- Electronic data entry is available

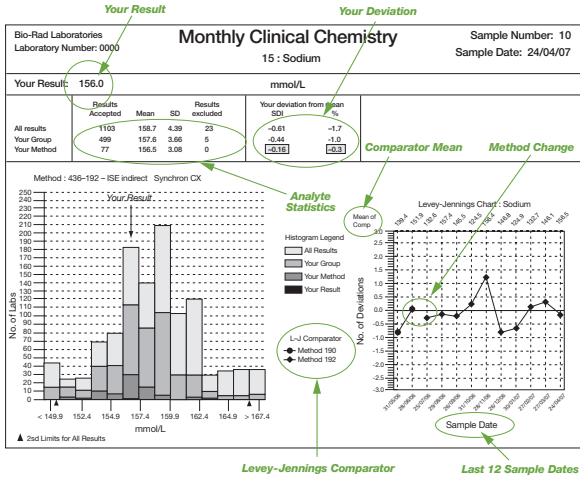
BIO-RAD

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.

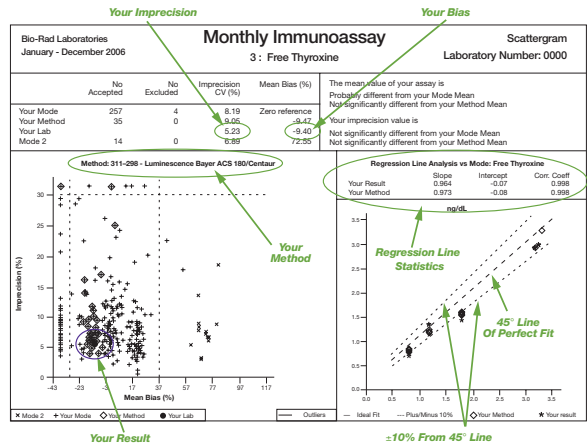
Monthly Report (Example)

Detailed individual statistical analyses and peer group comparisons are provided.



End-of-Cycle Report (Example)

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

Analytes for EQAS Urine Chemistry Program

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

Refer to the package insert of currently available lots for specific analyte and stability claims

Ordering Information (Additional Programs Available)

Cat #	EQAS Programs	Quantity
BC40	Urine Chemistry Program (12 analytes)	12 x 10 mL
BC45	Urine Chemistry Program (24 analytes)	12 x 10 mL
CS40	Urine Chemistry Program (12 analytes)	Report Only
CS45	Urine Chemistry Program (24 analytes)	Report Only
QC40	Urine Chemistry Program (specimen pack only)	12 x 10 mL



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 81-3-6361-6902 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