

EQAS Hematology Program

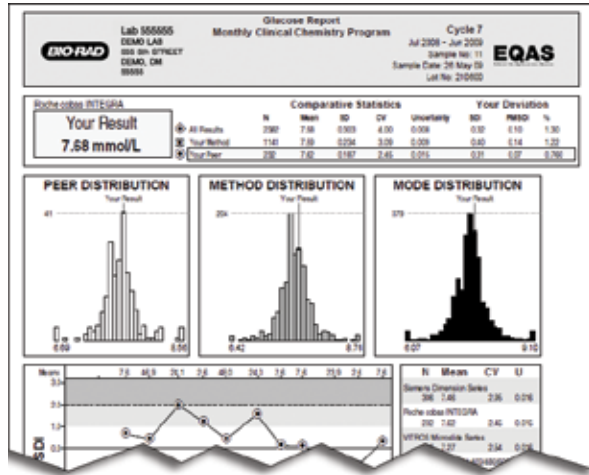
External Quality Assessment Suited for Your Laboratory

- Compare your laboratory's results to an international database of participants
- Convenience of primary piercable tubes
- 12 samples per cycle provided in 4 separate shipments
- Evaluate and report sample results on a monthly basis
- Select from 11 basic parameters (non-differential)
- For use on manual or automated analyzers
- Electronic data entry available
- Laboratory specific reports in an easy-to-read format with online viewing

BIO-RAD

EQAS Reports

Participants receive 12 sample reports over each program cycle. All reports are available within three days of the statistical analysis.



Sample Report (Example)

Detailed individual statistical analyses and peer group comparisons are provided.

Additional Reports (available on request): Sub-Group Reports

Analytes for EQAS Hematology Program

- | | | |
|--|--|-------------------------|
| Hematocrit (HCT) | Mean Corpuscular Volume (MCV) | Red Blood Cells (RBC) |
| Hemoglobin (Total) | Mean Platelet Volume (MPV) | White Blood Cells (WBC) |
| Mean Corpuscular Hemoglobin (MCH) | Platelets (PLT) | |
| Mean Corpuscular Hemoglobin Concentration (MCHC) | Red Blood Cell Distribution Width (RDW) | |
| | Red Blood Cell Distribution Width (RDW-SD) | |

Refer to the package insert for current parameter claims.

Ordering Information

Cat #	EQAS Programs	Quantity
BC90A	Hematology Program, Shipment A	3 x 2 mL
BC90B	Hematology Program, Shipment B	3 x 2 mL
BC90C	Hematology Program, Shipment C	3 x 2 mL
BC90D	Hematology Program, Shipment D	3 x 2 mL
CS90	Hematology Program	Report Only
QC90A	Hematology Program, Shipment A (specimen pack only)	3 x 2 mL
QC90B	Hematology Program, Shipment B (specimen pack only)	3 x 2 mL
QC90C	Hematology Program, Shipment C (specimen pack only)	3 x 2 mL
QC90D	Hematology Program, Shipment D (specimen pack only)	3 x 2 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