



EQAS eForms 2

**Automate Your EQAS Configuration, Results Entry
and Data Transmission With Our Easy-to-Use Software**

- No more handwritten result forms
- Eliminate late data caused by postal delays
- Eliminate transcription errors due to poor fax quality
- No more Method Questionnaires
- Codes automatically selected when you choose your instrument
- Reduce the use of paper to help preserve our planet



EQAS eForms 2

eForms 2 is available free of charge to all Bio-Rad EQAS participants. Please contact your local Bio-Rad representative for applicable programs.

Customize Your Lab Profile, Test Setup and Data Entry Options

- Eliminate the need to manually complete and mail in Change Notification Forms
- Quick and easy instrument setup tool
- Update your codes with one simple click
- Enter and save your lab's results at any time
- Sort your test display order for easy data entry
- Automatically generate a transmission file to email your data or upload it through QCNet.com
- Data transmission dates recorded for traceability

Analyte	Method	Instrument	Reagent	Unit
Ferritin	Chemiluminescence	Siemens ADVIA Centaur XP	Dedicated Reagent	µg/L
Folate	Chemiluminescence	Siemens ADVIA Centaur XP	Dedicated Reagent	nmol/L
HCG	Chemiluminescence	Siemens ADVIA Centaur XP	Dedicated Reagent	mIU/mL
T3, Free	Chemiluminescence	Siemens ADVIA Centaur XP	Dedicated Reagent	pmol/L
T3, Total	Chemiluminescence	Siemens ADVIA Centaur XP	Dedicated Reagent	nmol/L
T3/T Uptake	Chemiluminescence	Siemens ADVIA Centaur XP	Dedicated Reagent	% Uptake
T4, Free	Chemiluminescence	Siemens ADVIA Centaur XP	Dedicated Reagent	pmol/L
T4, Total	Chemiluminescence	Siemens ADVIA Centaur XP	Dedicated Reagent	nmol/L
TSH	Chemiluminescence	Siemens ADVIA Centaur XP	Dedicated Reagent	mIU/L



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 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 +36-1-459-6100 India 1-800-180-1224 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 7-495-721-14-04 Singapore 65-6415-3170 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