



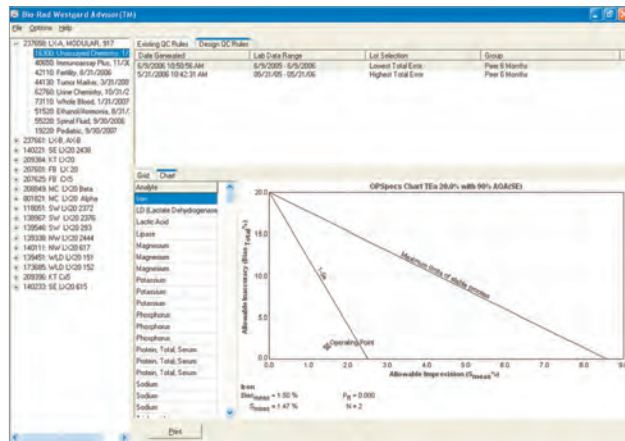
# Westgard Advisor™



# Westgard Advisor™

## Automatic QC Rules Selection Engine

- Recommend and automatically apply best QC rules with patented technology
- Easy step-by-step automatic rule selection capabilities
- Reduce false rejections and desensitization to false error flags
- Save time and money by reducing unnecessary repeats and troubleshooting
- Improve laboratory test quality with optimally selected QC rules
- Available as an optional module with Unity™ QC Data Management Solutions




Utilize OPSpecs charts to document test quality for each of your tests.

Level	TEa Selected	Bias %	CV	Sigma	Existing Revert Rule	Suggested Rule
CLIA	20.0	1.36	2.50	7.45		1-5s
3SD	14.4	0.879	4.66	2.95		1-3s/2-4s/1-10s
3V Mean	95.5	0.950	4.69	11.6		1-5s
CLIA	25.0	0.750	2.16	11.2		1-5s
CLIA	25.0	0.646	1.87	13.0		1-5s
CLIA	25.0	1.05	2.94	7.95		1-5s
CLIA	12.8	0.107	1.96	6.45		1-4s
CLIA	12.8	0.740	1.41	8.42		1-5s
CLIA	12.8	1.00	1.54	7.60		1-5s
3SD	11.5	0.016	3.11	3.77		1-3s/2-4s/1-10s
3SD	22.8	3.01	5.06	3.98		1-3s/2-4s/1-10s
3SD	19.9	2.28	2.12	4.52		1-2.5s
CLIA	10.0	0.036	1.24	7.40		1-5s
CLIA	10.0	0.180	2.38	4.13		1-2.5s
CLIA	10.0	0.438	1.33	7.21		1-4s
CLIA	10.0	0.213	0.960	2.68		1-3s/2-4s/1-10s
CLIA	2.99	1.33	1.11	1.49		1-3s/2-4s/1-10s

Automatically set up Westgard rules to evaluate each data point using a simple 4-step wizard.

## Ordering Information

Contact your local Bio-Rad Account Representative for more information.



Westgard Advisor™ is part of the Unity™ Interlaboratory Program. Participation in the Unity™ Interlaboratory Program can help improve the quality of your laboratory test results and help you meet regulatory and accreditation requirements worldwide. You may also subscribe to one of our Unity™ QC Data Management Solutions designed to improve the effectiveness of your statistical process control. Ask your Bio-Rad Account Representative for more information.

**Discover the power of Unity™ Solutions at [www.QCNet.com](http://www.QCNet.com).**



**Bio-Rad Laboratories**

For further information, please contact the Bio-Rad office nearest you or visit our website at [www.bio-rad.com/qualitycontrol](http://www.bio-rad.com/qualitycontrol)

Clinical Diagnostics Group

Website [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics) 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 86-21-64268080 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 91-124-4029300 Israel 972-3-9636050 Italy +39-02-216091 Japan 81-3-6361-7070 Korea 82-2-3473-4460 Mexico 52(55)5200-0520 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