

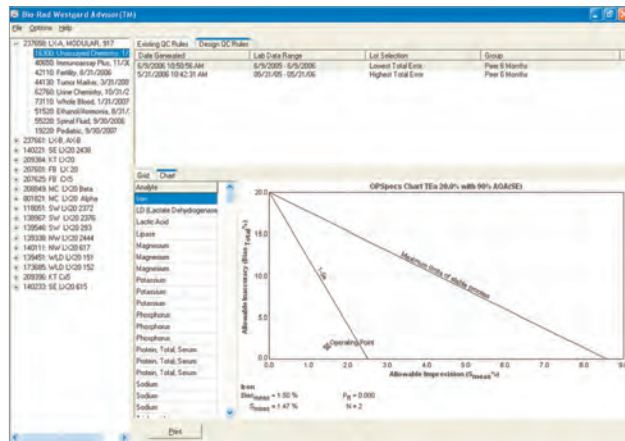
Westgard Advisor™



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	TE & Select	TE _A	Bias, %	CV	Sigma	Existing Revert Rule	Suggested Rule
1	CLIA	20.0	1.36	2.50	7.45		1-5s
1	3SD	14.4	0.679	4.66	2.96		1-3s/2-4s/1-10s
1	3V Mean/SD	95.5	0.950	4.69	11.6		1-5s
1	CLIA	25.0	0.750	2.16	11.2		1-5s
1	CLIA	25.0	0.646	1.87	13.0		1-5s
2	CLIA	25.0	-1.95	2.94	7.95		1-5s
1	CLIA	12.8	0.107	1.96	6.45		1-4s
1	CLIA	12.8	-0.740	1.41	8.42		1-5s
1	CLIA	12.8	-1.00	1.54	7.60		1-5s
1	3SD	11.5	0.016	3.11	3.77		1-3s/2-4s/1-10s
1	3SD	22.8	-3.01	5.06	3.98		1-3s/2-4s/1-10s
1	3SD	19.9	2.28	2.12	4.52		1-2.5s
2	CLIA	10.0	-0.036	1.24	7.40		1-5s
1	CLIA	18.0	0.180	2.38	4.13		1-2.5s
2	CLIA	16.0	-0.408	1.33	7.21		1-4s
1	CLIA	3.00	0.213	0.960	2.88		1-3s/2-4s/1-10s
1	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.



Bio-Rad Laboratories

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

Clinical Diagnostics Group

Website 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** 86-21-64268088 **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