



VIROTROL® HIV-1 Ag Control



VIROTROL® HIV-1 Ag

For use as an unassayed reactive quality control with assays detecting Human Immunodeficiency Virus Type 1 antigens.

- Human serum based
- 60 day open-vial stability at 2–8°C
- 18 month shelf life at 2–8°C
- Convenient liquid formulation eliminates potential reconstitution variability
- Recommended for Bio-Rad GENSCREEN® ULTRA HIV Ag-Ab Assay which uses HIV p24 antigen.



Use with negative control VIROCLEAR®

Analyte

HIV-1 Ag

Refer to the package insert for performance and stability claims

Ordering Information

Cat #	Product Name	Quantity
VIROTROL® HIV-1 Ag		
00108	Single Level/Dropper Bottle (Available outside the United States only)	1 x 5 mL
VIROCLEAR®		
00106	Single Level/Dropper Bottle	1 x 5 mL
00107	Single Level/Dropper Bottle	2 x 20 mL
00112	Single Level/Tube	10 x 4 mL

Please note that not all products may be available in all countries.



This control is featured in 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.



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/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 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