



VIROTROL[®] HAV IgM



VIROTROL® HAV IgM

For use as an unassayed quality control with assays detecting Immunoglobulin M (IgM) antibodies to Hepatitis A Virus.

- Human serum based
- Convenient liquid formulation eliminates potential reconstitution variability
- 3 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C



Use with negative control VIROCLEAR®

Analyte

HAV IgM

Refer to the package insert for performance and stability claims

Ordering Information

| Cat # | Product Name | Quantity |
|--------------------------|---------------------------------------|-----------|
| VIROTROL® HAV IgM | | |
| 00142 | Single Level/Dropper Bottle | 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.



**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** 86-21-64260808 **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