



VIROTROL® ToRCH



VIROTROL® ToRCH

For use as an unassayed, reactive quality control with assays detecting IgG antibodies to *Toxoplasma gondii*, IgG antibodies to Rubella Virus, IgG antibodies to Cytomegalovirus and IgG antibodies to Herpes Simplex Virus Types 1 and 2.

- Human serum based
- Multi-analyte control allows for product consolidation
- Convenient liquid formulation eliminates potential reconstitution variability
- 2 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C



Use with negative control VIROCLEAR® ToRCH

Analytes

Toxoplasma gondii IgG
 Rubella IgG
 Cytomegalovirus IgG

Herpes Simplex Virus Type 1 (HSV-1) IgG
 Herpes Simplex Virus Type 2 (HSV-2) IgG

Refer to the package insert for performance and stability claims

Ordering Information

Cat #	Product Name	Quantity
VIROTROL® ToRCH		
00109	Single Level/Dropper Bottle	1 x 5 mL
VIROCLEAR® ToRCH		
00118	Single Level/Dropper Bottle	1 x 5 mL



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/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