



# VIROCLEAR® ToRCH



## VIROCLEAR® ToRCH

For use as an unassayed, non-reactive quality control with assays detecting IgG and IgM antibodies to *Toxoplasma gondii*, IgG and IgM antibodies to Rubella Virus, IgG and IgM antibodies to Cytomegalovirus, and IgG and IgM 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 positive controls VIROTROL® ToRCH and VIROTROL® ToRCH-M

### Analytes

**Non-reactive for:**

*Toxoplasma gondii* IgG  
*Toxoplasma gondii* IgM  
 Rubella IgG  
 Rubella IgM  
 Cytomegalovirus IgG

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

Refer to the package insert for performance and stability claims

### Ordering Information

Cat #	Product Name	Quantity
<b>VIROCLEAR® ToRCH</b>		
00118	Single Level/Dropper Bottle . . . . .	1 x 5 mL
<b>VIROTROL® ToRCH</b>		
00109	Single Level/Dropper Bottle . . . . .	1 x 5 mL
<b>VIROTROL® ToRCH-M</b>		
00117	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](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/qualitycontrol](http://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