



Liquichek™ ToRCH Plus Controls



Liquichek™ ToRCH Plus Control

A bilevel, liquid control used to monitor the performance of immunoassay methods to detect IgG antibodies with the most widely used serology instruments and methods. Both positive and negative controls allow you to monitor the performance of your test system at the critical decision making levels.



- Liquid, human serum based
- 3 year shelf life at 2-8°C
- 60 day open-vial stability at 2-8°C
- Positive and negative controls packaged separately
- Assayed and unassayed versions available

Liquichek™ ToRCH Plus IgM Control

A bilevel, unassayed liquid control used to monitor the performance of immunoassay methods to detect IgM antibodies to ToRCH panel analytes.



- Liquid, human serum based
- Positive and negative controls packaged separately
- 3 year shelf life when stored at -20°C to -70°C
- 60 day open-vial stability at 2-8°C

Analytes

Toxoplasma gondii IgG
Rubella Virus IgG
Cytomegalovirus (CMV) IgG
Herpes Simplex Virus
Type 1 (HSV-1) IgG
Herpes Simplex Virus
Type 2 (HSV-2) IgG

Epstein-Barr Virus (VCA) IgG
Varicella Zoster Virus (VZV) IgG
Helicobacter pylori IgG
Treponema pallidum (Syphilis) IgG
Syphilis RPR

Refer to the package insert of currently available lots for performance and stability claims

Analytes

Toxoplasma gondii IgM
Rubella Virus IgM
Cytomegalovirus (CMV) IgM
Herpes Simplex Virus Type 1 (HSV-1) IgM
Herpes Simplex Virus Type 2 (HSV-2) IgM

Refer to the package insert of currently available lots for performance and stability claims

Ordering Information

Cat #	Liquichek™ ToRCH Plus Control	Quantity
239	Positive	3 x 3 mL
239X	Positive MiniPak.	1 x 3 mL
227	Positive, Assayed (available outside the United States only)	3 x 3 mL
227X	Positive, Assayed MiniPak (available outside the United States only)	1 x 3 mL
228	Negative	3 x 3 mL
228X	Negative MiniPak.	1 x 3 mL

Ordering Information

Cat #	Liquichek™ ToRCH Plus IgM Control	Quantity
229	Positive	3 x 2 mL
229PX	Positive MiniPak.	1 x 2 mL
230	Negative	3 x 2 mL
230NX	Negative MiniPak.	1 x 2 mL



These controls are 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