



VIROTROL Infectious Disease Controls

For use with the Ortho Clinical Diagnostics VITROS® HIV Combo Assay

Bio-Rad controls contain suitable levels of Anti-HIV-1, Anti-HIV-2, Anti-HIV-1 group O, and HIV-1 Ag, which are recommended for use with the Ortho VITROS® HIV Combo Assay. Improve your lab's efficiency and confidence in reporting patient test results with a comprehensive line of human matrix, independent quality controls that offer:

- An unbiased assessment
- Liquid, ready to use convenience
- Multiple packaging configurations designed to suit your lab's specific needs
- Worldwide peer group comparison available through Unity Software



VIROTROL I

Anti-HIV-1 Positive Control

An unassayed quality control for assays detecting antibodies to Human Immunodeficiency Virus Type 1.

- Human serum based
- 3 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C
- Available in multiple configurations and fill sizes

VIROTROL HIV-1 Ag

HIV-1 Antigen Positive Control

An unassayed quality control for assays detecting Human Immunodeficiency Virus Type 1 antigens.

- Human serum based
- 2 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C

VIROTROL HIV-1 gO

Anti-HIV-1 gO Positive Control

An unassayed control for the Ortho VITROS® HIV Combo Assay used to detect antibodies to HIV-1 group O.

- Human serum based
- 2 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C

VIROTROL HIV-2

Anti-HIV-2 Positive Control

An unassayed quality control for assays detecting antibodies to Human Immunodeficiency Virus Type 2.

- Human serum based
- 2 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C

VIROCLEAR

Negative Control

An unassayed, non-reactive quality control. Non-reactive for commonly tested Retrovirus analytes.

- Human serum based
- 3 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C
- Available in multiple configurations and fill sizes

Ordering Information

Cat # Description

VIROTROL I

00100E Single Level/Dropper Bottle 1 x 5 mL
 00101E Single Level/Tube 10 x 4 mL

VIROTROL HIV-1 Ag

00108A Single Level/Dropper Bottle 1 x 5 mL

VIROTROL HIV-1 gO

00113 Single Level/Tube 5 x 4 mL
 00113X Single Level/Tube 1 x 4 mL

VIROTROL HIV-2

00105C Single Level/Dropper Bottle 1 x 5 mL

VIROCLEAR

00106 Single Level/Dropper Bottle 1 x 5 mL
 00112 Single Level/Tube 10 x 4 mL

Refer to myinserts.com or the package insert of currently available lots for specific analyte and stability claims.

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

Product availability in other countries subject to local regulatory requirements.



EQAS

An independent, external assessment of performance in comparison to your peers.



Independent QC

Ongoing, proactive, unbiased daily QC that helps identify errors as they occur or begin to trend.



Unity

QC Data Management tools that help you create a strategy to reduce risk and streamline QC workflow.



**Bio-Rad
Laboratories, Inc.**

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com/idcontrols

Clinical
Diagnostics Group

Website www.bio-rad.com/qualitycontrol **USA** 1 800 224 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 8901 **Belgium** 32 03 710 53 00
Brazil 55 31 3689 6600 **Canada** 1 514 334 4372 **China** 86 21 61698500 **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** 1800 180 1224
Israel 972 3 9636050 **Italy** 39 02 216091 **Japan** 81 3 6361 7070 **Korea** 82 2 3473 4460 **Mexico** 52 55 5488 7670 **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 1404 **Singapore** 65 6415 3170
South Africa 27 11 442 85 08 **Spain** 34 91 590 5200 **Sweden** 46 8 555 127 00 **Switzerland** 41 0 26 674 55 05 06 **Taiwan** 886 2 2578 7189
Thailand 662 651 8311 **United Kingdom** 44 0 20 8328 2000

The Bio-Rad logo or Bio-Rad Laboratories products and services are not affiliated with Ortho Clinical Diagnostics. Information is subject to change without notice. EQAS, Unity and VIROTROL are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. VITROS is a registered trademark of Ortho Clinical Diagnostics.

© 2017 Bio-Rad Laboratories, Inc. Printed in USA QSD17-272 Q-1224 Rev. 09/2017

