

Infectious Disease Panels

Ordering Information

Catalog No. Description

BioPlex® 2200 EBV IgG Panel

665-1250	BioPlex®2200 EBV IgG Reagent Pack.....	1 pack
663-1200	BioPlex®2200 EBV IgG Calibrator Set.....	1 set
663-1230	BioPlex®2200 EBV IgG Control Set.....	1 set

BioPlex® 2200 EBV IgM Panel

665-1350	BioPlex®2200 EBV IgM Reagent Pack.....	1 pack
663-1300	BioPlex®2200 EBV IgM Calibrator Set.....	1 set
663-1330	BioPlex®2200 EBV IgM Control Set.....	1 set

BioPlex® 2200 Syphilis IgG Panel

665-1450	BioPlex®2200 Syphilis IgG Reagent Pack.....	1 pack
663-1400	BioPlex®2200 Syphilis IgG Calibrator Set.....	1 set
663-1430	BioPlex®2200 Syphilis IgG Control Set.....	1 set

BioPlex® 2200 Syphilis IgM Panel‡

665-1550	BioPlex®2200 Syphilis IgM Reagent Pack.....	1 pack
663-1500	BioPlex®2200 Syphilis IgM Calibrator Set.....	1 set
663-1530	BioPlex®2200 Syphilis IgM Control Set.....	1 set

BioPlex® 2200 ToRC IgG Panel

665-1650	BioPlex®2200 ToRC IgG Reagent Pack.....	1 pack
663-1600	BioPlex®2200 ToRC IgG Calibrator Set.....	1 set
663-1630	BioPlex®2200 ToRC IgG Control Set.....	1 set

BioPlex® 2200 ToRC IgM Panel‡

665-1750	BioPlex®2200 ToRC IgM Reagent Pack.....	1 pack
663-1700	BioPlex®2200 ToRC IgM Calibrator Set.....	1 set
663-1730	BioPlex®2200 ToRC IgM Control Set.....	1 set

BioPlex® 2200 Rubella & CMV IgM Panel*

665-3850	BioPlex®2200 Rubella & CMV IgM Reagent Pack.....	1 pack
663-3800	BioPlex®2200 Rubella & CMV IgM Calibrator Set.....	1 set
663-3830	BioPlex®2200 Rubella & CMV IgM Control Set.....	1 set

BioPlex® 2200 HSV-1 & HSV-2 IgG Panel

665-3350	BioPlex®2200 HSV-1 & HSV-2 IgG Reagent Pack.....	1 pack
663-3300	BioPlex®2200 HSV-1 & HSV-2 IgG Calibrator Set.....	1 set
663-3330	BioPlex®2200 HSV-1 & HSV-2 IgG Control Set.....	1 set

BioPlex® 2200 MMRV IgG Panel

665-2450	BioPlex®2200 MMRV IgG IgG Reagent Pack.....	1 pack
663-2400	BioPlex®2200 MMRV IgG IgG Calibrator Set.....	1 set
663-2430	BioPlex®2200 MMRV IgG IgG Control Set.....	1 set

Additional BioPlex® 2200 Panels in Development

BioPlex®2200 HIV	BioPlex®2200 Hepatitis
------------------	------------------------

‡ Available Outside US Only
* Available in US Only



BioPlex® 2200 Infectious Disease Panels

Advancing productivity in infectious disease testing with a broad menu of assays and multiplex panels

EBV IgG and IgM
Syphilis IgG and IgM†
ToRC IgG and IgM†
Rubella & CMV IgM*

HSV-1 & HSV-2 IgG
MMRV IgG
HIV†
Hepatitis†

† In development
‡ Available outside US only
* Available in US only




Bio-Rad
Laboratories

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

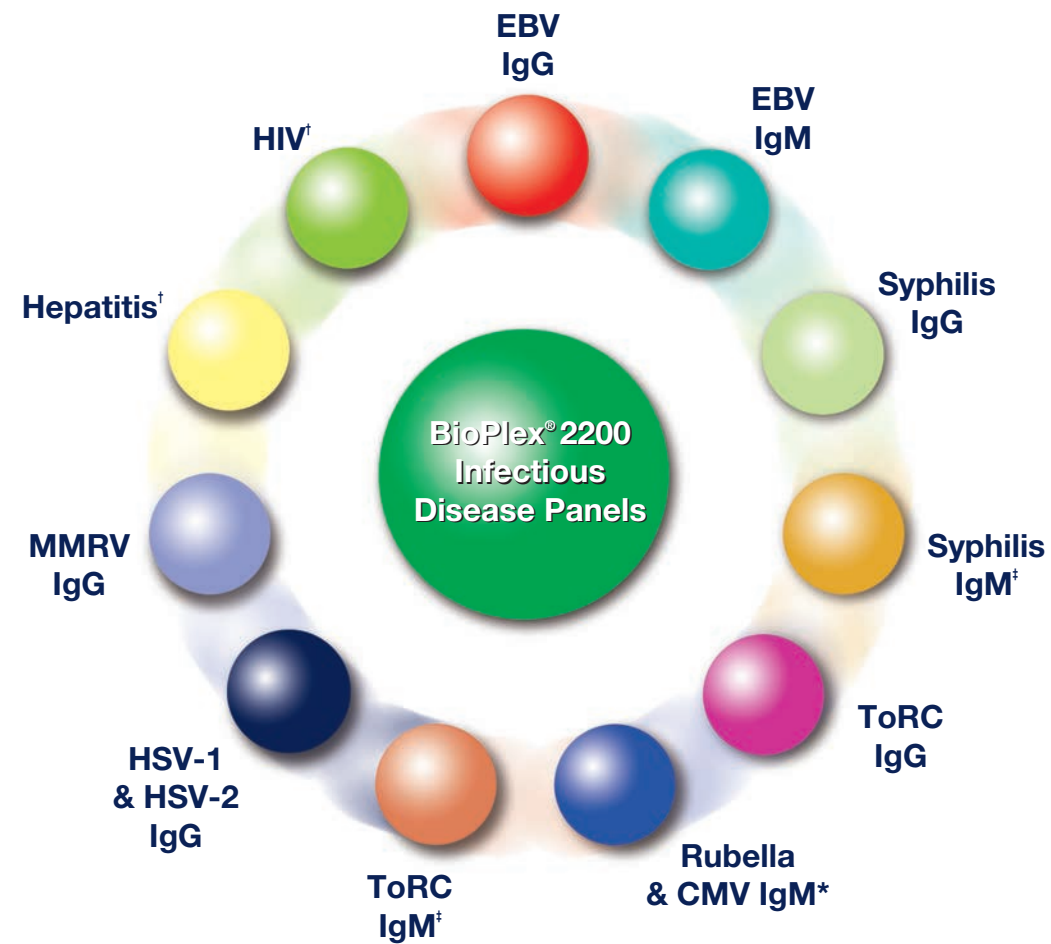
Clinical
Diagnostics Group

Website www.bio-rad.com/diagnostics **U.S.** 1-800-2BIO-RAD (1-800-224-6723) **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** 1-800-180-1224 **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** 7-495-721-14-04 **Singapore** 65-6415-3170 **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

 Printed on recycled paper with soy-based inks

© 2012 Bio-Rad Laboratories, Inc. Printed in USA 03/12 DG11-1258 T-167





BioPlex® 2200 Infectious Disease Panels

Leading the way with an expanding multiplexed assay menu

Comprehensive assay menu offering multiplexed solutions for simultaneous detection of important infectious disease markers.

Offering the benefits of multiplex testing:

- High throughput multiplex testing improves workflow efficiencies and eliminates hands-on steps
- Consolidates traditional single-analyte tests into one panel
- Time-to-first result in approximately 45 minutes
- Ready-to-use multiplexed calibrators and controls, with curve stability up to 30 days

† In development
‡ Available outside US only
* Available in US only

Epstein-Barr Virus

BioPlex® 2200 EBV IgG and IgM Panels



The BioPlex® 2200 EBV IgG and IgM kits provide laboratories with more relevant information for improved diagnosis of Infectious Mononucleosis and for management and follow-up of EBV infected patients.

BioPlex® 2200 EBV IgG

Analytes

- EBV VCA IgG
- EBV NA-1 IgG
- EBV EA-D IgG

BioPlex® 2200 EBV IgM

Analytes

- EBV VCA IgM
- Heterophile IgM

Herpes Simplex Virus

BioPlex® 2200 HSV-1 & HSV-2 Panel



The BioPlex® 2200 HSV-1 and HSV-2 kit simultaneously detects and differentiates the two most clinically relevant IgG antibodies in serum and plasma.

BioPlex® 2200 HSV-1 & HSV-2 IgG

Analytes

- HSV-1 IgG
- HSV-2 IgG

Syphilis

BioPlex® 2200 Syphilis IgG and IgM Panels



The BioPlex® 2200 Syphilis IgG and IgM kits detect the five most clinically relevant antibodies against individual recombinant proteins to *Treponema pallidum* in a multiplex format.

BioPlex® 2200 Syphilis IgG

Analytes

- Syphilis IgG Composite[‡]
- T. pallidum* 15kDa**
- T. pallidum* 17kDa**
- T. pallidum* 47kDa**

BioPlex® 2200 Syphilis IgM[†]

Analytes

- Syphilis IgM Composite[‡]
- T. pallidum* 17kDa**
- T. pallidum* 47kDa**

‡ Available outside US only
‡ Syphilis IgG and IgM are a rollup of the individual results.
** Individual results available outside US

Measles, Mumps, Rubella and Varicella-zoster virus

BioPlex® 2200 MMRV IgG Panel



The BioPlex® 2200 MMRV IgG kit provides a fast, accurate and efficient method to determine serological status.

Toxoplasmosis, Rubella and Cytomegalovirus Infections

BioPlex® 2200 ToRC IgG and IgM[†] Panels



The BioPlex® 2200 ToRC IgG kit offers a simple, centralized, cost-effective method for the simultaneous identification of antibodies associated with the most commonly tested infectious diseases.

BioPlex® 2200 ToRC IgG

Analytes

- Toxoplasma gondii* (quantitative)
- Rubella (quantitative)
- Cytomegalovirus (qualitative)

BioPlex® 2200 ToRC IgM[†]

Analytes

- Toxoplasma gondii* (qualitative)
- Rubella (qualitative)
- Cytomegalovirus (qualitative)

BioPlex® 2200 Rubella & CMV IgM^{*}

Analytes

- Rubella (qualitative)
- Cytomegalovirus (qualitative)

‡ Available outside US only
* Available in US only

BioPlex® 2200 MMRV IgG

Analytes

- Measles IgG
- Mumps IgG
- Rubella IgG (German Measles)
- Varicella-zoster virus IgG