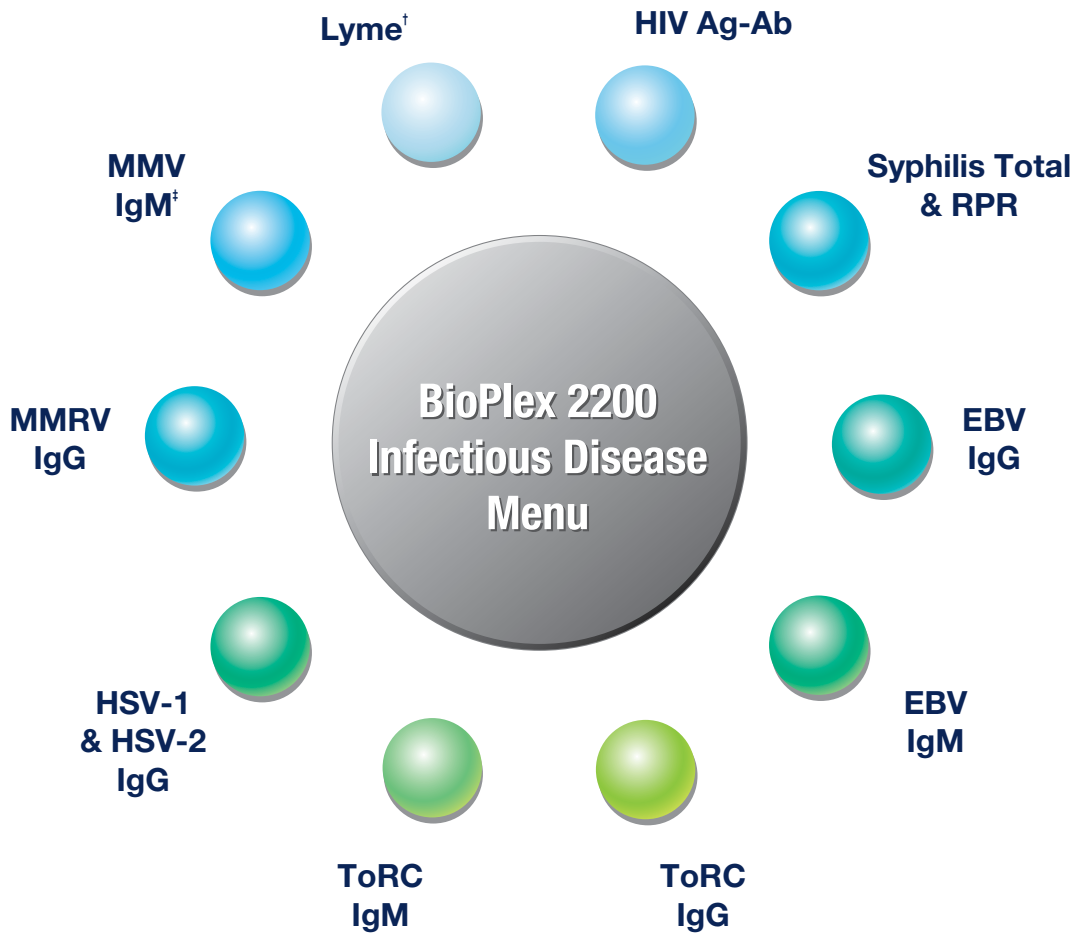


BioPlex 2200 Infectious Disease Panels

An Expanding Multiplexed Assay Menu





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 as little as 45 minutes
- Ready-to-use multiplexed calibrators and controls, with curve stability up to 30 days

‡ Available outside US only
† In development



BioPlex 2200 Infectious Disease Panels

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

- HIV Ag-Ab
- EBV IgG and IgM
- ToRC IgG and IgM
- HSV-1 & HSV-2 IgG
- MMRV IgG
- MMV IgM[†]
- Syphilis Total & RPR
- Lyme[‡]

[‡] Available outside US only
[†] In development

Human Immunodeficiency Virus

BioPlex 2200 HIV Ag-Ab Panel



The BioPlex 2200 HIV Ag-Ab kit offers enhanced multiplexed HIV screening by independently detecting and reporting antibodies and antigen, differentiating HIV-1 and HIV-2, and distinguishing early (acute/primary) from established infection.

BioPlex 2200 HIV Ag-Ab^Δ

Analytes

HIV Ag-Ab (composite) with

- HIV-1 p24 Ag
- HIV-1 Ab (Groups M & O)
- HIV-2 Ab

^Δ All analytes are reported in the US

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

Syphilis

BioPlex 2200 Syphilis Total & RPR Panel



The BioPlex 2200 Syphilis Total & RPR kit offers a one-step universal testing method by simultaneously detecting and differentiating treponemal and non-treponemal antibodies.

BioPlex 2200 Syphilis Total & RPR

Analytes

- Syphilis Total (IgG/IgM)
- Qualitative RPR (Rapid Plasma Reagin)
- Onboard RPR titer determination (1:4, 1:8, 1:16, 1:32, 1:64)

Measles, Mumps, Rubella and Varicella-zoster Virus

BioPlex 2200 MMRV IgG and MMV IgM[‡] Panels



The BioPlex 2200 MMRV IgG and MMV IgM kit provides a fast, accurate, and efficient method to simultaneously determine serological status (immunity).

BioPlex 2200 MMRV IgG

Analytes

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

BioPlex 2200 MMV IgM[‡]

Analytes

- Measles IgM
- Mumps IgM
- Varicella-zoster virus IgM

[‡] Available outside US only

Toxoplasma, Rubella and Cytomegalovirus

BioPlex 2200 ToRC IgG and IgM Panels



The BioPlex 2200 ToRC IgG and ToRC IgM kits offer a simple, centralized, and cost-effective method for the simultaneous identification of antibodies commonly tested in prenatal care.

BioPlex 2200 ToRC IgG

Analytes

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

BioPlex 2200 ToRC IgM

Analytes

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

Herpes Simplex Virus

BioPlex 2200 HSV-1 & HSV-2 IgG Panel



The BioPlex 2200 HSV-1 and HSV-2 IgG 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

‡ Available outside US only

Additional BioPlex 2200 Panels in Development

BioPlex 2200 Lyme

Ordering Information

Catalog No.	Description	Catalog No.	Description
BioPlex 2200 HIV Ag-Ab Panel			
665-3450	BioPlex 2200 HIV Ag-Ab Reagent Pack (OUS) 1 Pack	665-2450	BioPlex 2200 MMRV IgG Reagent Pack..... 1 Pack
665-3455	BioPlex 2200 HIV Ag-Ab Reagent Pack (US/Canada)..... 1 Pack	663-2400	BioPlex 2200 MMRV IgG Calibrator Set 1 Set
663-3400	BioPlex 2200 HIV Ag-Ab Calibrator Set (OUS)..... 3 Set	663-2430	BioPlex 2200 MMRV IgG Control Set 2 Set
663-3405	BioPlex 2200 HIV Ag-Ab Calibrator Set (US/Canada) 3 Set	BioPlex 2200 MMV IgM Panel†	
663-3430	BioPlex 2200 HIV Ag-Ab Control Set (OUS)..... 3 Set	12000930	BioPlex 2200 MMV IgM Reagent Pack..... 1 Pack
663-3435	BioPlex 2200 HIV Ag-Ab Control Set (US/Canada) 3 Set	12000933	BioPlex 2200 MMV IgM Calibrator Set..... 1 Set
		12000931	BioPlex 2200 MMV IgM Control Set 2 Set
BioPlex 2200 Syphilis Total & RPR Panel			
12000650	BioPlex 2200 Syphilis Total & RPR Reagent Pack... 1 Pack	BioPlex 2200 ToRC IgG Panel	
12000651	BioPlex 2200 Syphilis Total & RPR Calibrator Set 1 Set	665-1650	BioPlex 2200 ToRC IgG Reagent Pack..... 1 Pack
12000653	BioPlex 2200 Syphilis Total & RPR Control Set 2 Set	663-1600	BioPlex 2200 ToRC IgG Calibrator Set 1 Set
		663-1630	BioPlex 2200 ToRC IgG Control Set 2 Set
BioPlex 2200 EBV IgG Panel			
665-1250	BioPlex 2200 EBV IgG Reagent Pack..... 1 Pack	BioPlex 2200 ToRC IgM Panel	
663-1200	BioPlex 2200 EBV IgG Calibrator Set 1 Set	12000670	BioPlex 2200 ToRC IgM Reagent Pack..... 1 Pack
663-1201	BioPlex 2200 EBV IgG Calibrator Set (OUS) 1 Set	12000677	BioPlex 2200 ToRC IgM Calibrator Set 1 Set
663-1230	BioPlex 2200 EBV IgG Control Set 2 Set	12000678	BioPlex 2200 ToRC IgM Control Set..... 2 Set
663-1231	BioPlex 2200 EBV IgG Control Set (OUS)..... 2 Set	BioPlex 2200 HSV-1 & HSV-2 IgG Panel	
BioPlex 2200 EBV IgM Panel			
665-1350	BioPlex 2200 EBV IgM Reagent Pack 1 Pack	665-3350	BioPlex 2200 HSV-1 & HSV-2 IgG Reagent Pack 1 Pack
663-1300	BioPlex 2200 EBV IgM Calibrator Set 1 Set	663-3300	BioPlex 2200 HSV-1 & HSV-2 IgG Calibrator Set..... 1 Set
663-1330	BioPlex 2200 EBV IgM Control Set..... 2 Set	663-3330	BioPlex 2200 HSV-1 & HSV-2 IgG Control Set 2 Set

† Available outside US only



**Bio-Rad
Laboratories, Inc.**

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

Clinical
Diagnostics Group

Web site www.bio-rad.com/diagnostics **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