



BioPlex™ 2200 EBV IgG and IgM

For a complete EBV serological diagnostic profile

The BioPlex™ 2200 EBV IgG and IgM kits simultaneously detect the five most clinically relevant antibodies.

By detecting the five antibodies of highest clinical significance, including heterophile antibody, 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:

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

BioPlex™ 2200 EBV IgM:

- EBV VCA IgM
- Heterophile IgM



BioPlex™ 2200 EBV IgG and EBV IgM tests allow physicians to receive a complete EBV serological diagnostic profile of their patient.

1 sample



5 Tests

VCA IgG
EAD IgG
EBNA IgG
VCA IgM
Heterophile



Result:
Complete Serological Profile for Disease Interpretation

Primary Acute
Late Acute
Recovering
Previous Infection
Susceptible

Efficiency

- Fully-automated, random access multiplex assay procedure for EBV IgG & EBV IgM testing from primary tube to final result
- Consolidates 3 traditional single-analyte tests into 1 EBV IgG kit; 2 traditional single-analyte tests into 1 EBV IgM kit
- Up to 190 results per hour and up to 5 different reportable test results from 1 small patient sample
- Time to first result in approximately 45 minutes
- Ready-to-use multiplexed calibrators and controls
- Curve stability up to 30 days

Quality

- Good serological status agreement with commercially available ELISA microplate and agglutination tests methods resulting in 88.7% for Acute status and 90.4% in Non-Acute status*
- Excellent lot-to-lot reproducibility with %CV ≤ 3.1% for EBV IgG tests, and ≤ 5.5% for EBV IgM tests
- For every sample processed, three internal quality controls are also processed, improving confidence in results

* Please refer to Instructions for Use (IFU)

Ordering Information

Catalog No. Description

BioPlex™ 2200 EBV IgG

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

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



**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 **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