



# BioPlex™ 2200 EBV IgG and IgM

The first and only fully-automated, multiplexed solution for Epstein-Barr Virus Antibody testing





# Like no other

**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:

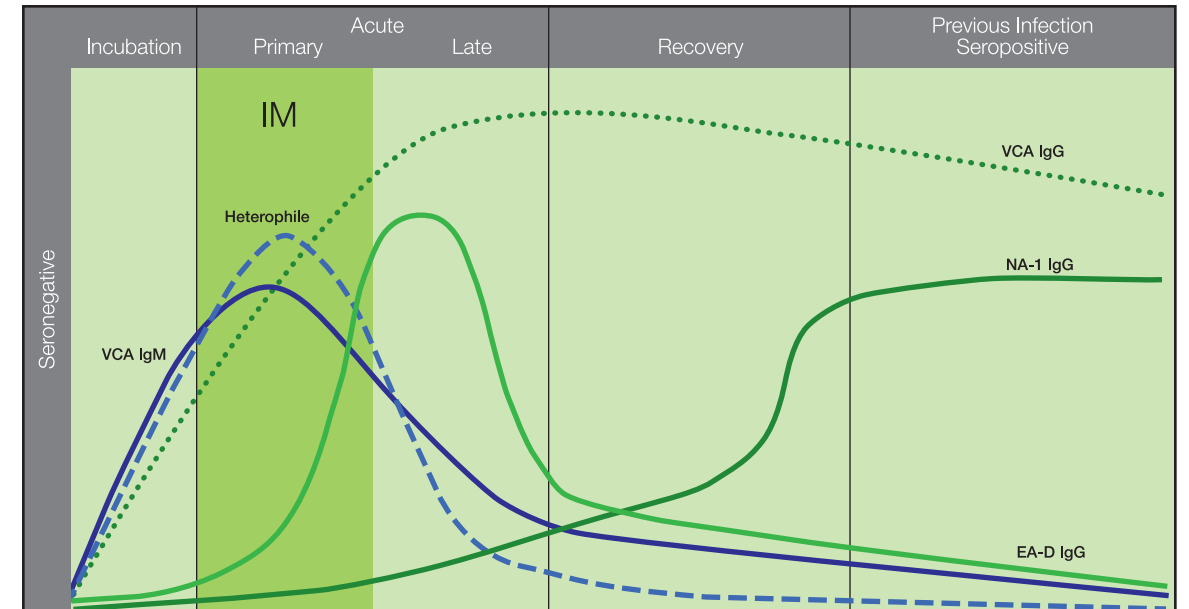
- EBV VCA IgM
- Heterophile IgM

## EBV Infection Diagnosis

Infectious Mononucleosis (IM) diagnosis is typically based on the detection of Heterophile antibodies and antibodies against specific Epstein-Barr Virus antigens:

- Viral Capsid Antigen: EBV VCA IgG and EBV VCA IgM
- Nuclear Antigen: EBV NA-1 IgG
- Early Antigen: EBV EA-D IgG

Combining Heterophile IgM and VCA IgM detection enhances the detection of acute IM. Determining the complete Patient EBV Serological profile is essential to define the stage of the infection.



	Incubation	Primary Acute	Late	Recovery	Previous Infection Seropositive
<b>Heterophile</b>		<b>+/-</b> Normally present in acute IM	<b>+/-</b>	<b>-</b> Typically absent during recovery phase	<b>-</b>
<b>VCA IgM</b>		<b>+/-</b> Normally present in acute IM	<b>+/-</b>	<b>-</b> Generally absent during recovery phase	<b>-</b> A rise can be seen in reactivation phase
<b>VCA IgG</b>		<b>+/-</b> Normally present in acute IM	<b>+</b>	<b>+</b> Generally present during recovery phase	<b>+</b> Normally persists for lifetime
<b>EA-D IgG</b>		<b>+/-</b> Frequently present in acute IM	<b>+/-</b>	<b>-/+</b> Generally absent during recovery phase	<b>-</b> A rise can be seen in reactivation phase
<b>NA-1 IgG</b>		<b>-</b> Rarely present in acute IM	<b>+/-</b>	<b>+/-</b> Typically rise during recovery phase	<b>+/-</b> Normally persists for lifetime

Adapted from: J.M. Seigneurin, Apport du laboratoire dans l'infection à virus Epstein-Barr Laboratory diagnosis of Epstein-Barr virus infections. Immuno-analyse & Biologie Spécialisée 1 (2002) 33-39.

## Quality



### Automated EIA-like Heterophile Antibody Detection

All the benefits of full automation and EIA technology:

- Minimizes subjectivity linked to non-EIA methods
- Improves reproducibility

### Enhanced Acute Infection Diagnosis

Combining heterophile with VCA IgM enhances the clinical diagnosis of acute Infectious Mononucleosis.

## Confidence



### Full Automation

From primary tube to final result, Bioplex™ 2200 is the only fully-automated, multiplexed EBV testing system:

- Minimizes sample and reagent handling errors
- Standardizes the assay procedure for improved performance

### Excellent Precision

Heterogeneous assay format and approximately 200 readings per analyte dramatically enhance assay precision.

### Antibodies of Clinical Utility

By associating the five essential EBV antibodies, BioPlex™ 2200 EBV IgG and IgM kits provide complete information to precisely determine EBV patient status.

### Internal Quality Controls

For every sample processed, three internal quality controls are also processed, dramatically improving confidence in the results:

- Serum Verification Bead (SVB) confirms presence of serum
- Internal Standard Bead (ISB) standardizes detector performance
- Reagent Blank Bead (RBB) identifies sample problems associated with non-specific binding

## Efficiency



### High Throughput Capability

Up to 190 IgG and IgM results per hour for improved workflow and reduced turnaround time.

### Optimal Ease-of-Use

All reagents, calibrators and controls are ready-to-use and all the procedure steps are fully-automated. No operator intervention is required from sample loading to result release, significantly reducing hands-on-time, labor costs and cost per patient result.

### eFlex™ Software

The system features a powerful package of software tools that provides optimum flexibility:

- “Add-on” of new test orders from previously run patient samples
- Creation of custom EBV panels
- Medical Decision Support Software (MDSS)\* that correlates patient antibody results with specific EBV profiles for improved information to clinicians about EBV status
- Bi-directional LIS connectivity

\* Under development for future software versions.

## Comparison Using Serological Status

618 results were obtained using commercially available ELISA microplate and agglutination tests and the BioPlex™ 2200 EBV IgG and IgM kits. Samples were then assigned a serological status for both methods using a generally accepted EBV Serology Interpretation table. The BioPlex™ 2200 EBV IgG and IgM kits show excellent correlation with commercially available assays for each disease status, with fewer samples categorized as Inconclusive.

EBV Serological Status	BioPlex 2200 EBV IgG & IgM Profile									
	Primary Acute	Late Acute	Recovering	Previous Infection	Susceptible	Inconclusive	Total	% Serological Agreement	95% Confidence Interval	
Primary Acute	30	0	0	0	0	1	31	96.8%	83.8 - 99.4%	
Late Acute	5	90	1	13	0	1	110	81.8%	73.6 - 87.9%	
Recovering	1	0	3	0	0	0	4	75.0%	30.0 - 95.4%	
Previous Infection	0	31	2	263	4	5	305	86.2%	81.9 - 89.7%	
Susceptible	4	0	0	0	122	1	127	96.1%	91.1 - 98.3%	
Inconclusive	6	10	0	7	11	7	41	17.1%	8.5 - 31.3%	
Overall	46	131	6	283	137	15	618	83.3%	80.2 - 86.1%	

EBV Serological Status	BioPlex 2200 EBV IgG & IgM Profile					
	Primary Acute	Non-Acute	Inconclusive	Total	% Serological Agreement	95% Confidence Interval
Acute	125	14	2	141	88.7%	82.4 - 92.9%
Non-Acute	36	394	6	436	90.4%	87.2 - 92.8%
Inconclusive	16	18	7	41	17.1%	8.5 - 31.3%
Overall	177	426	15	618	85.1%	82.1 - 87.7%

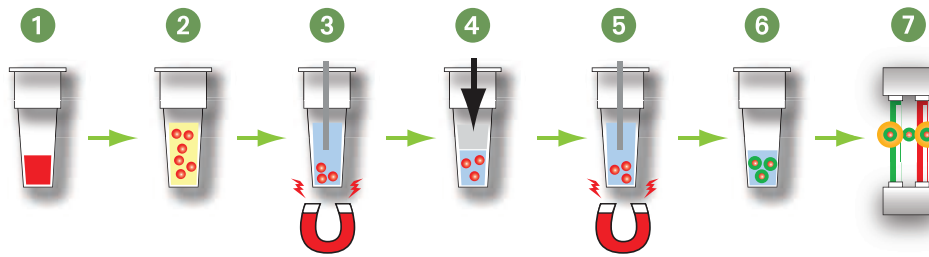
## Reproducibility

Following CLSI EP5-A guidelines, the BioPlex™ 2200 EBV IgG and IgM kit positive controls were run at 3 clinical sites for 3 days, 4 times a day, for a total of 36 replicates. The BioPlex™ 2200 EBV IgG and IgM kits offer excellent reproducibility for both Within-Run and Between-Run calculations. Outstanding reproducibility can be attributed to: a) full automation, b) magnetic bead, heterogenous format of the assay and c) approximately 200 measurements per analyte (up to 4 times other multiplex methods).

Positive Control	Sample N	Grand Mean AI	Within-Run		Between-Run	
			SD	%CV	SD	%CV
EBV VCA IgG	36	2.3	0.1	2.8%	0.1	2.7%
EBV NA-1 IgG	36	2.9	0.1	1.9%	0.1	2.4%
EBV EA-D IgG	36	3.0	0.1	3.1%	0.1	2.5%

Positive Control	Sample N	Grand Mean AI	Within-Run		Between-Run	
			SD	%CV	SD	%CV
EBV VCA IgM	36	2.0	0.1	5.5%	0.1	3.0%
Heterophile IgM	36	2.5	0.1	5.3%	0.0	0.0%

## EBV IgG and IgM Assays Procedure



1. Patient sample (5  $\mu$ L) automatically added to Reaction Vessel
2. Sample diluent & beads (dilution 1:60) added and incubated at 37°C
3. Wash step
4. Conjugate added and incubated at 37°C
5. Wash step
6. Bead resuspended in wash buffer
7. Flow-based, dual laser detection



### EBV IgG and IgM Reagent Packs

Each contains all the necessary reagents (sample diluent, beads and conjugate) to process 100 samples per pack and up to 3 (IgG) and 2 (IgM) results per sample.



### EBV IgG and IgM Calibrator Sets

The EBV IgG and IgM kit calibrators are multiplexed and ready-to-use.



### EBV IgG and IgM Control Sets

The EBV IgG and IgM kit controls are multiplexed and ready-to-use.

## Ordering Information

**Catalog No.**      **Description**

### BioPlex 2200 EBV IgG Panel

665-1250	EBV IgG Panel .....	1 pack
663-1200	EBV IgG Calibrator Set .....	1 set
663-1230	EBV IgG Control Set .....	1 set

### BioPlex 2200 EBV IgM Panel

665-1350	EBV IgM Panel .....	1 pack
663-1300	EBV IgM Calibrator Set .....	1 set
663-1330	EBV IgM Control Set .....	1 set

### BioPlex 2200 System and Bulk Reagents

660-0000	BioPlex™ 2200 .....	1 system
660-0800	Sample Racks (30 per box) .....	1 box
660-0801	Sample Tray .....	1 tray
660-0817	Instrument Sheath Fluid (2 per box) .....	1 box
660-0818	Instrument Wash Buffer .....	1 box
660-2003	Reaction Vessels (1000 per bag) .....	1 bag
666-0001	Instrument Detector Calibration Pack .....	1 pack
666-0002	Instrument Detector Clean Pack .....	1 pack
666-0003	Instrument Probe Cleaning Solution (4 per box) .....	1 box



**Bio-Rad  
Laboratories**

*For further information, please contact the Bio-Rad office nearest you or visit our website at [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics).*

*Clinical  
Diagnostics Group*

**Website** [www.bio-rad.com/diagnostics](http://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** 36-1-455-8800    **India** 91-124-4029300    **Israel** 972-3-9636050    **Italy** +39-02-216091  
**Japan** 81-3-5811-6290    **Korea** 82-2-3473-4460    **Mexico** +52-(555)-200-0520    **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