

BioPlex 2200 MMRV IgG and MMV IgM Kits

Simultaneous detection of multiple antibodies to highly contagious, vaccine-preventable diseases

BIO-RAD



BioPlex 2200 MMRV IgG and MMV IgM Assays

The BioPlex 2200 MMRV IgG and MMV IgM assays provide a fast, accurate and efficient method to determine serological status or acute infection for measles, mumps, rubella, and varicella- zoster virus. The BioPlex 2200 System also offers aid in diagnosis of current or recent infection of rubella IgM in the BioPlex 2200 ToRC IgM assay. The BioPlex 2200 System provides laboratories with efficient workflow and faster turnaround time for these commonly ordered infectious disease assays.

The BioPlex 2200 MMRV IgG assay reports qualitative results for the following

- Measles IgG
- Mumps IgG
- Rubella IgG
- Varicella-Zoster Virus IgG

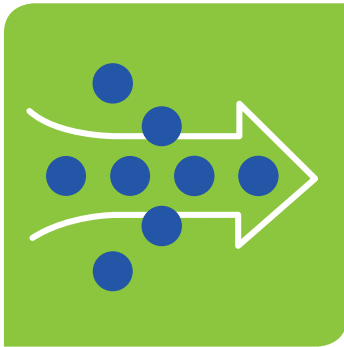
The BioPlex 2200 MMV IgM assay reports qualitative results for the following

- Measles IgM
- Mumps IgM
- Varicella-Zoster Virus IgM



Efficiency

- Multiplex your workflow – offering four tests simultaneously for MMRV IgG and three tests simultaneously for MMV IgM
- Multiplex your inventory – dramatically reducing reagent and calibrator storage requirements
- Multiplex your samples – significantly reducing your sample volume requirements
- Multiplex your tubes – accommodating serum, heparinized plasma, and EDTA



Throughput

- Market-leading throughput with up to 400 multiplexed test results per hour to meet heavy workload demands
- Random access processing ensures results are available when you need them
- Time to first result of 45 minutes maximizes speed and flexibility in result reporting



Confidence

- Full automation from primary tube to final result, minimizing the chance of human error
- Internal quality control beads included with each sample to verify serum addition and standardize detector performance
- One system with an expanding menu to meet your autoimmune and serology testing needs



Performance

Performance of the BioPlex 2200 MMRV IgG and MMV IgM assays were evaluated against corresponding commercially available assays. Results from these studies are shown in the tables below.

Clinical studies demonstrate excellent positive agreement for the BioPlex 2200 MMRV IgG assay

Population	Pregnant Women (N=396)	Pre-Employment Screening (N=393)	Test Ordered (N=790)
Measles IgG	97.6% (359/368)	99.7% (355/356)	98.2% (638/650)
Mumps IgG	99.7% (346/347)	100% (341/341)	95.8% (646/674)
Rubella IgG	99.5% (362/364)	98.9% (344/348)	97.6% (702/719)
VZV IgG	98.9% (348/352)	97.8% (359/367)	96.1% (688/716)

Note: % agreement calculations exclude equivocal results

Clinical studies demonstrate excellent negative agreement for the BioPlex 2200 MMRV IgG assay

Assay	Negative % Agreement
Measles IgG	100% (93/93)
Mumps IgG	100% (83/83)
Rubella IgG	97.6% (123/126)
VZV IgG	100% (143/143)

Note: % agreement calculations exclude equivocal results

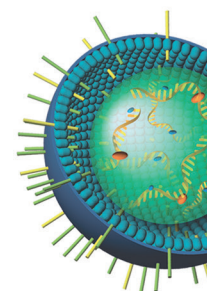
Clinical studies demonstrate excellent positive agreement for the BioPlex 2200 MMV IgM assay

Assay	Positive % Agreement
Measles IgM	96.0% (95/99)
Mumps IgM	96.6% (86/89)
VZV IgM	98.8% (84/85)

Note: % agreement calculations exclude equivocal results

Clinical studies demonstrate excellent negative agreement for the BioPlex 2200 MMV IgM assay

Population	Healthy Donor (n=300)	Test Ordered (n=300)
Measles IgM	100% (299/299)	99.3% (294/296)
Mumps IgM	97.9% (285/291)	96.9% (251/259)
VZV IgM	100% (297/297)	97.9% (279/285)





Precision

Precision was assessed using serum, EDTA plasma and heparinized plasma samples according to the CLSI EP5-A3 guideline "Evaluation of Precision Performance of Clinical Chemistry Devices." In this study, 40 runs were performed on one BioPlex 2200 System over 20 working days using a reproducibility panel of samples.

MMRV

	Within-Run % CV	Between-Run % CV	Total Precision % CV
Measles IgG	3.1-9.8	3.4-7.4	9.4-16.0
Mumps IgG	2.4-5.6	2.1-5.3	5.3-10.2
Rubella IgG	3.3-6.2	2.7-7.4	8.0-15.2
VZV IgG	2.4-7.3	1.9-10.6	6.1-10.0

MMV IgM

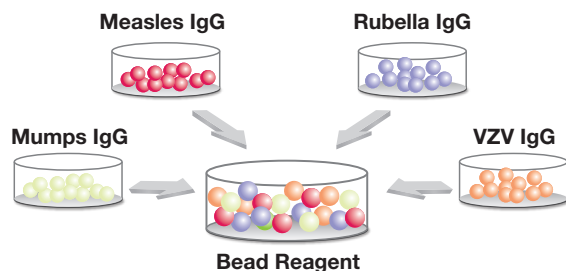
	Within-Run % CV	Between-Run % CV	Total Precision % CV
Measles IgM	0.0-4.2	0.0-2.2	0.0-4.2
Mumps IgM	2.7-6.2	0.0-2.1	3.0-6.2
VZV IgM	2.5-9.5	0.0-2.3	2.5-9.8



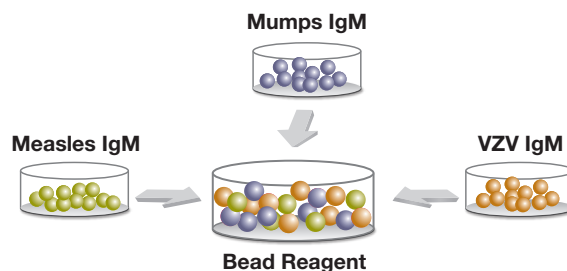
Assay Principle

The BioPlex 2200 MMRV IgG and MMV IgM assays are multiplexed bead based immunoassays designed to detect and differentiate antibodies to Measles, Mumps, Rubella, and Varicella-zoster Virus. Each bead reagent is a mixture of disease specific dyed microparticle beads. The beads are individually manufactured and combined with internal QC beads into a bead reagent for multiplex analysis.

MMRV IgG Multiplex Testing



MMV IgM Multiplex Testing



MMRV IgG and MMV IgM Reagent Packs

The MMRV IgG and MMV IgM Reagent Packs contain all the necessary reagents (sample diluent, beads and conjugate) to process 100 samples per pack for MMRV IgG and 150 samples per pack for MMV IgM.



MMRV IgG and MMV IgM Calibrator Sets

The MMRV IgG and MMV IgM calibrators are multiplexed and ready to use offering 30 day calibration curve stability.



MMRV IgG and MMV IgM Control Sets

The MMRV IgG and MMV IgM controls are multiplexed and ready to use.

Ordering Information

Catalog No. Description

665-2450 BioPlex 2200 MMRV IgG Reagent Pack, 100 samples..... 1 pack
663-2400 BioPlex 2200 MMRV IgG Calibrator Set..... 1 set
663-2430 BioPlex 2200 MMRV IgG Control Set2 sets

12000930 BioPlex 2200 MMV IgM Reagent Pack, 150 samples* 1 pack
12000933 BioPlex 2200 MMV IgM Calibrator Set* 1 set
12000931 BioPlex 2200 MMV IgM Control Set*2 sets

* Not Available in the U.S.

BIO-RAD

**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.** 1-800-224-6723 **Australia** 61-2-9914-2800 **Austria** 43-1-877-8901 **Belgium** +32 (3)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