



## BioPlex 2200 MMV IgM Assay

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

Measles, mumps, and varicella-zoster virus (VZV) are highly contagious diseases that are endemic in many regions. Although these diseases are preventable through vaccination, outbreaks still occur worldwide due to lower immunization rates, waning immunity, and lack of antibody response to the vaccines. The BioPlex 2200 MMV IgM assay allows laboratories to simultaneously detect and differentiate IgM class antibodies to all three diseases.

### **Consolidated into one test kit**

A multiplex flow immunoassay for the qualitative detection of IgM antibodies to:

- Measles (Rubeola)
- Mumps
- Varicella-Zoster Virus

**BIO-RAD**

# The Power of Multiplex Testing

## Traditional Testing:

Physician orders 3 individual tests  
Measles IgM, Mumps IgM, and  
VZV IgM



## Multiplex Testing:

Physician orders 3 individual  
tests Measles IgM, Mumps IgM,  
and VZV IgM



## Efficiency

- Fully automated multiplex assay for Measles, Mumps, and VZV IgM testing
- Consolidates 3 separate traditional analyte tests into a single test
- Time to first result in 45 minutes, up to 189 results per hour
- Ready-to-use calibrators and controls
- Calibrator curve stability up to 30 days
- Reagent pack open vial stability up to 90 days

## Quality and Performance

- Excellent assay reproducibility with within-run and between-run % CV of less than 10% for Measles, Mumps, and VZV
- Two internal quality beads are processed with every sample ensuring system and result reliability
- Excellent positive agreement with commercially available methods of 96.0%, 96.6%, and 100% for Measles IgM, Mumps IgM, and VZV IgM, respectively
- Excellent negative agreement with commercially available methods of 99.3%, 96.9%, and 97.9% for Measles IgM, Mumps IgM, and VZV IgM, respectively

## Ordering Information

Catalog No.	Description	
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  
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.** 1-800-224-6723 **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