



# BioPlex™ 2200 MMRV IgG

**For comprehensive determination of serological status that meets evolving clinical practice**

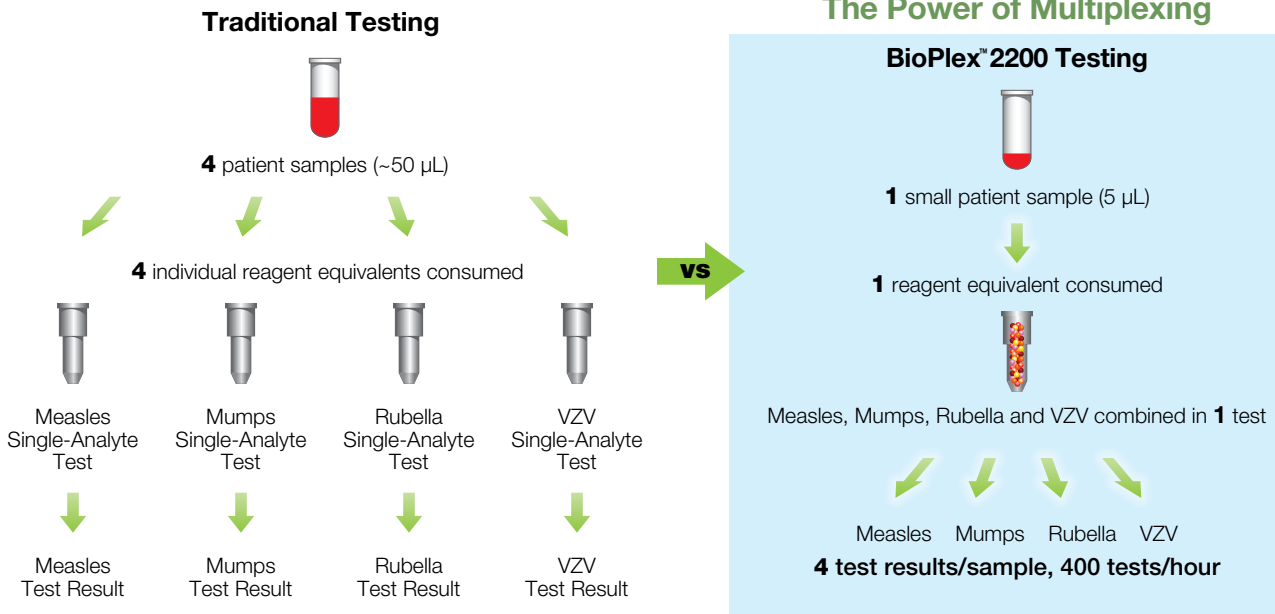
Measles, Mumps, Rubella and Varicella-zoster virus (VZV) are infectious diseases that are preventable through the practice of vaccination. Testing for serological status is an alternative to vaccination in cases where serological status is uncertain. Ideally, healthcare workers should be immune to these infectious diseases. The BioPlex™ 2200 MMRV IgG kit provides a fast, accurate and efficient method to determine serological status for all four infectious agents, simultaneously.

Consolidated into one test panel

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

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

**BioPlex™ 2200 MMRV IgG**



**Benefits of Multiplex MMRV Testing**

- Consolidates 4 traditional single-analyte tests in 1 infectious disease panel
- Reduces reagent storage requirements
- Reduces the chance of errors with fewer hands-on steps
- Reduces sample volume required to run multiple tests

**Precision**

- Excellent assay reproducibility with within-run and between-run %CV less than 10% for Measles, Mumps, Rubella and VZV

**Efficiency**

- High throughput with 100 samples per hour and up to 400 multiplexed test results per hour
- Rapid turnaround time with first result available in 45 minutes

**A Leader in Serology Testing**

- Offering the only fully-automated, random access MMRV kit
- Offering an expanding menu of BioPlex™ 2200 assays including EBV, Syphilis, ToRC and HSV

**Ordering Information**

Catalog No.	Description	
665-2450	BioPlex™ 2200 MMRV IgG Reagent Pack.....	1 pack
663-2400	BioPlex™ 2200 MMRV IgG Calibrator Set .....	1 set
663-2430	BioPlex™ 2200 MMRV IgG 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](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 +36-1-459-6100 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

Printed on recycled paper with soy-based inks

© 2009 Bio-Rad Laboratories, Inc. Printed in USA 10/09 DG09-0282 T-160