

BioPlex 2200 ToRC IgM Assay

Simultaneous detection of multiple antibodies for prenatal screening

The BioPlex 2200 ToRC IgM Assay targets IgM antibodies required for determining serological status for *T. gondii*, Rubella and Cytomegalovirus (CMV). The assay utilizes the power of multiplex testing to simultaneously identify and differentiate antibodies commonly tested in prenatal care.

BioPlex is a registered trademark of Bio-Rad Laboratories, Inc. in various jurisdictions.

Consolidated into one test panel

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

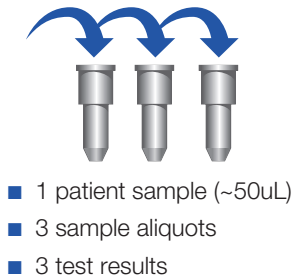
- *Toxoplasma gondii*
- Rubella
- Cytomegalovirus

BIO-RAD

The Power of Multiplex Testing

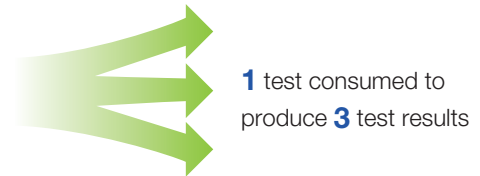
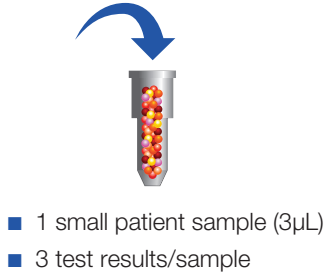
Traditional Testing:

Physician orders 3 individual tests
T.gondii IgM, Rubella IgM, and
 CMV IgM



Multiplex Testing:

Physician orders 3 individual tests
T.gondii IgM, Rubella IgM, and
 CMV IgM



Efficiency

- Fully-automated multiplex assay procedure for *T.gondii* IgM, Rubella IgM, and Cytomegalovirus IgM testing from primary tube to final result
- Takes one patient sample and consolidates 3 traditional single analyte tests
- Time to first result in 45 minutes, 180 results per hour thereafter
- Ready-to-use calibrators and controls
- Curve stability up to 30 days
- Kit open vial stability up to 60 days

Quality and Performance

- Two internal quality beads are processed with every sample ensuring system and result reliability
- Excellent negative agreement with commercially available methods resulting in 97.4%, 99.6%, and 97.9% for *T.gondii*, Rubella, and Cytomegalovirus respectively
- Excellent positive agreement with commercially available methods resulting in 97.1%, 98.0%, and 98.6% for *T.gondii*, Rubella, and Cytomegalovirus, respectively

Ordering Information

Catalog No.	Description	
12000670	BioPlex 2200 ToRC IgM Reagent Pack, 150 samples.....	1 pack
12000677	BioPlex 2200 ToRC IgM Calibrator Set	1 set
12000678	BioPlex 2200 ToRC IgM Control Set	2 sets



**Bio-Rad
 Laboratories, Inc.**

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. 1-800-224-6783 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