

BioPlex™ 2200 ToRC IgG

Simultaneous detection of multiple antibodies for antenatal monitoring

Determining the serological status of potentially infected individuals, especially pregnant women, is critical in the prevention of life-threatening infections that may lead to congenital abnormalities. The BioPlex™ 2200 ToRC IgG offers a simple, centralized, cost-effective method for the simultaneous identification of antibodies associated with the most commonly tested infectious diseases such as:

- Toxoplasmosis
- Rubella (German Measles)
- Cytomegalovirus Infection

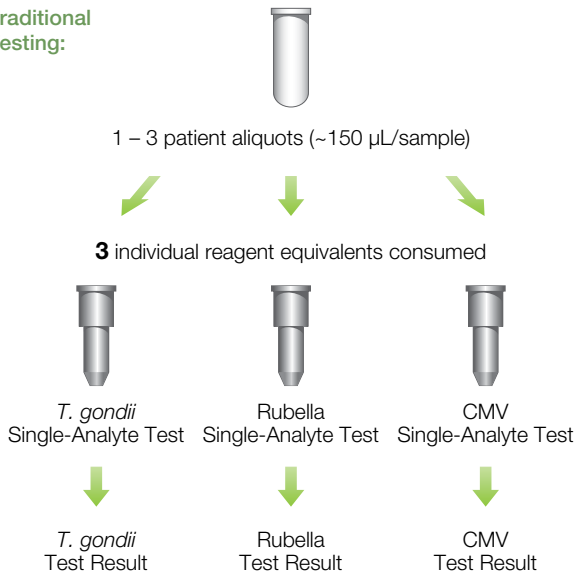
Consolidated into one test

A multiplex flow immunoassay for the detection in human serum or plasma of IgG antibodies to:

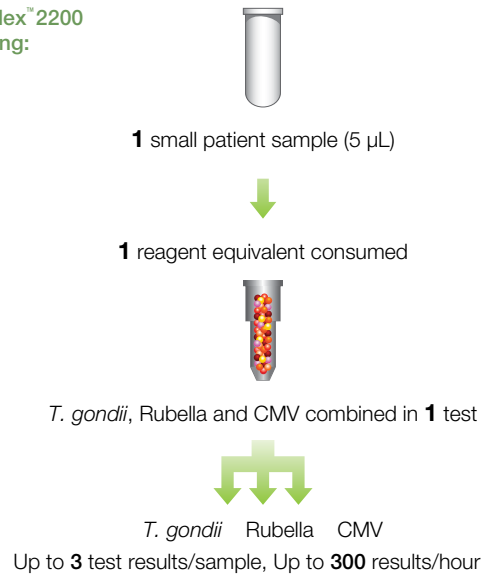
- *Toxoplasma gondii* (quantitative)
- Rubella (quantitative)
- Cytomegalovirus (qualitative)

BioPlex™ 2200 ToRC IgG

Traditional Testing:



BioPlex™ 2200 Testing:



Benefits of Multiplex ToRC Testing

- Consolidates 3 traditional single-analyte tests into 1 infectious disease panel
- Reduces reagent storage requirements
- Fewer hands-on steps reduces the chance of errors
- Dramatically reduces sample volume required to run multiple tests

Quality and Performance

- Excellent correlation with CDC evaluation serum panel resulting with 100% positive and 100% negative agreement for each *T. gondii* IgG, Rubella IgG and Cytomegalovirus IgG*
- Excellent lot-to-lot reproducibility near the cutoff (breakpoint) with %CV ≤ 13.6% for *T. gondii* IgG, ≤ 8.7% for Rubella IgG and ≤ 11.1% for Cytomegalovirus IgG*
- Assays calibrated to reference standards – 3rd International Standard, TOXM for *T. gondii* IgG and the 1st International Standard, RUBI-1-94 for Rubella IgG
- Three internal quality beads are processed with every sample ensures system and results reliability

* Please refer to Instructions for Use (IFU)

Ordering Information

Catalog No.	Description	
665-1650	BioPlex™ 2200 ToRC IgG Reagent Pack	1 pack
663-1600	BioPlex™ 2200 ToRC IgG Calibrator Set	1 set
663-1630	BioPlex™ 2200 ToRC 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

Clinical Diagnostics Group

Website 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 +31-318-540686 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