

Bio-Rad Laboratories
Improving lives with every test.

Rubella

Rubella IgG and IgM EIAs

Rubella, sometimes called German measles, is an acute viral infection that causes a mild illness in children and slightly more severe illness in adults.

The disease is spread person-to-person through airborne contact. Rubella infection in pregnant women, especially in the first trimester, may cause serious complications in the fetus, including a range of severe birth defects.

Demonstration of rubella IgG antibody in women prior to conception provides assurance of fetal protection from possible rubella viral infection during pregnancy. Aggressive immunization of pre-school children has greatly reduced the incidence of rubella epidemics, but a need still exists for accurate monitoring of immune status, especially for women of child-bearing age. Laboratory criteria for diagnosis of rubella include isolation of the rubella virus, or a significant rise in rubella antibody level by any standard serologic assay, or positive serologic test for rubella IgM antibody. Serology is the most common method of confirming a diagnosis.

Bio-Rad offers Rubella IgG and IgM Microplate EIA assays for qualitative and semi-quantitative detection.

Convenient

- Ready-to-use reagents
- Rapid turnaround

Objective

- Microplate format
- Qualitative and semi-quantitative detection
- Results normalized as index values

Simple

- 96-test kit
- Indirect EIA

IgG EIA Testing

- As an aid in the assessment of the patient's immunological response to Rubella
- Screening to determine immune status
- Identify active infection

IgM EIA Testing

- Estimate the time of infection by using IgM EIA in conjunction with the IgG EIA

BIO-RAD

Rubella EIA Testing



Rubella Indirect EIA Procedure

1



- Microplate: breakable wells dry coated with Ag

2



- Pipette 100 μ L of diluted test samples, calibrators or samples into the wells (1:51 for IgG/1:26 for IgM)
- Incubate 30 minutes at room temperature
- 4 washes

3

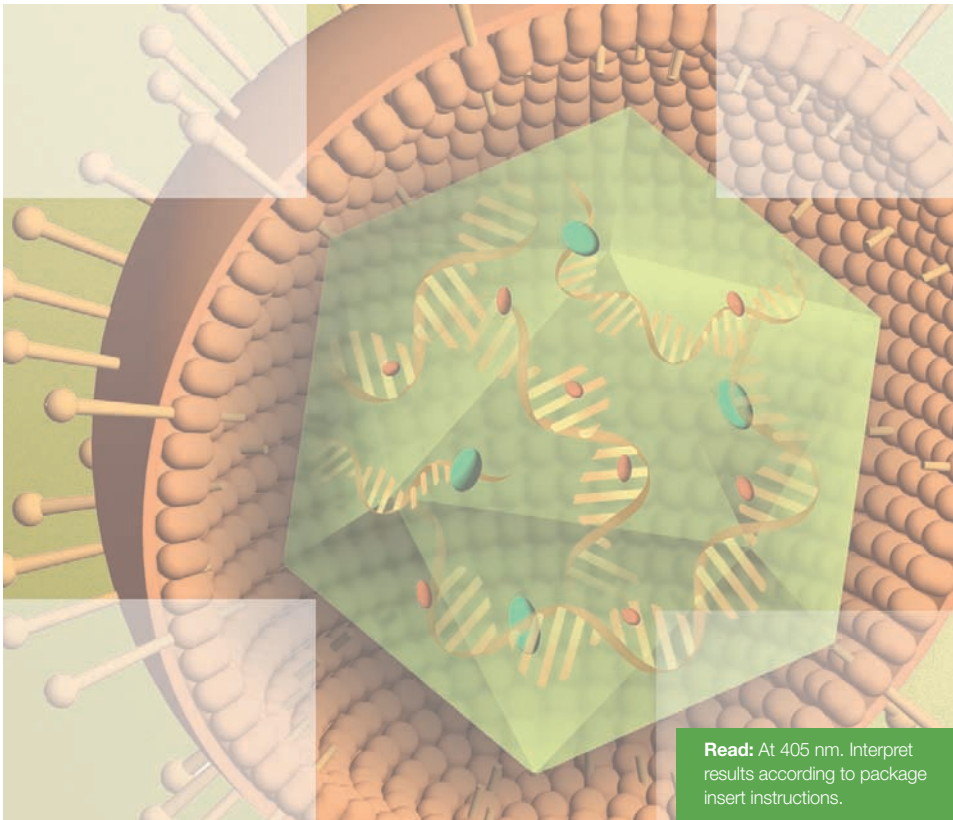


- Dispense 100 μ L of ready-to-use conjugate
- Incubate 30 minutes at room temperature
- 4 washes

4



- Dispense 100 μ L of ready-to-use chromogen/substrate
- Incubate 30 minutes at room temperature
- Add 100 μ L of ready-to-use Stop Reagent



Ordering Information

Catalog No.	Description	Tests
25173	Rubella IgG EIA, Microplate	96 tests
25174	Rubella IgM EIA, Microplate	96 tests



Bio-Rad Laboratories

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com.

Clinical Diagnostics Group

Website www.bio-rad.com U.S. 1-800-2-BIO-RAD Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium 32-9-385-5511 Brazil 5521-2507-6191 Canada 1-514-334-4372 Czech Republic 420-2-41430532 China 86-21-63052255 Denmark 45-4452-1000 Finland 358-9-804-22-00 France 33-1-4795-6000 Germany 49-89-31884-0 Hong Kong 852-2789-3300 India 91-124-6398112 Israel 972-3-9514127 Italy 39-02-216091 Japan 81-3-5811-6290 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 5255-5534-2552 The Netherlands 31-318-540666 New Zealand 64-9-415-2280 Norway 47-23-38-41-30 Poland 48-22-331-99-99 Portugal 351-21-472-7700 Russia 7-095-721-14-00 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 44-208-328-2000