

Bio-Rad Laboratories
Improving lives with every test.

Rubeola

Measles IgG EIA

Measles is a highly contagious, acute, exanthematous disease. It is generally self-limiting and without serious consequences, although complications such as bronchopneumonia and otitis media do occur.

The most serious consequence of measles is encephalomyelitis, which is fortunately rare (about 1 in 10,000 cases). Natural infection with measles virus confers permanent immunity.

The Bio-Rad Measles IgG test is an enzyme immunoassay test in a microwell format. Test results are obtained after a 90-minute incubation. The results are objective and normalized as index values, permitting uniformity of results.

The Bio-Rad Measles IgG test is intended for the qualitative and semi-quantitative detection of human IgG antibodies to measles virus in human serum by EIA. Individual serum specimens may be used for the determination of immune status. Paired (acute/convalescent) sera may be used to demonstrate seroconversion or significant rises in antibody level, as an aid in the diagnosis of primary infection.

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 Measles
- Monitor IgG levels to identify active infection



Measles IgG EIA Testing

Measles IgG 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)
- Incubate 30 minutes at room temperature
- 4 washes

3

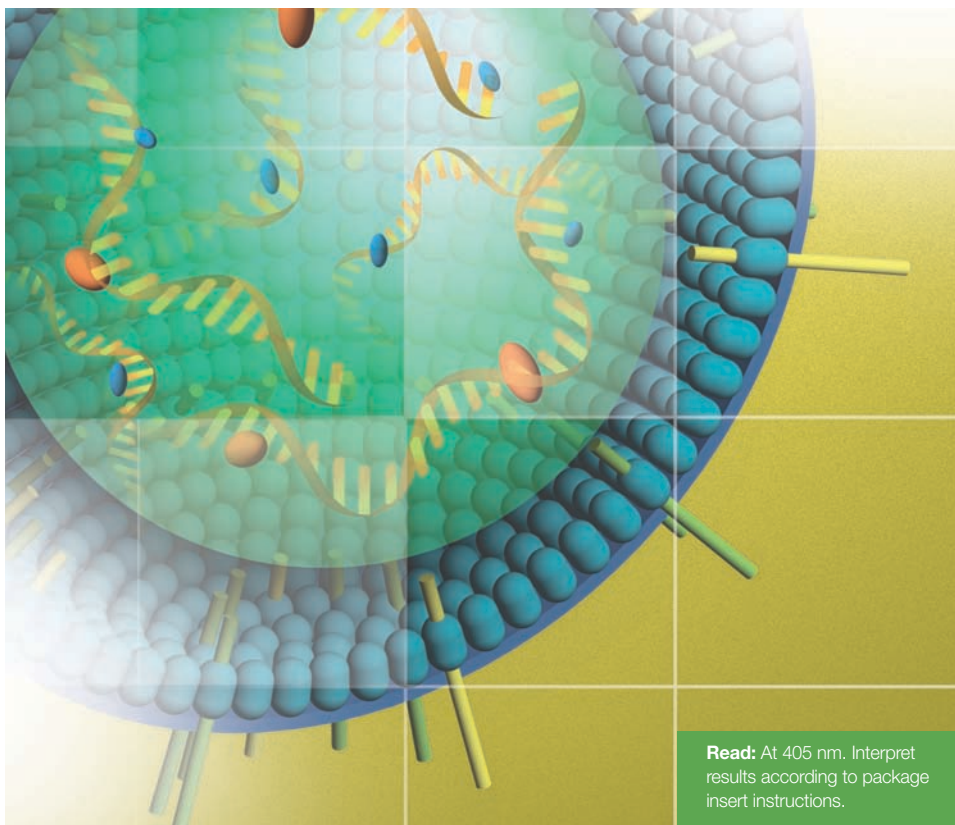


- Dispense 100 μ L of ready-to-use conjugate
- Incubate 30 minutes at room temperature
- 4 washes

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



Read: At 405 nm. Interpret results according to package insert instructions.

Ordering Information

Catalog No.	Description
25181	Measles IgG EIA Kit 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