

**Bio-Rad Laboratories**  
Improving lives with every test.

# Mumps

## Mumps IgG EIA

**Mumps is a generalized illness characterized by fever, and inflammation and swelling of the salivary glands, particularly the parotid glands.**

Mumps is usually not severe in children, but in adults the inflammation may involve the ovaries or testes (orchitis). Mumps is also one of the most common causes of aseptic meningitis and encephalitis. The etiological agent is a member of the paramyxovirus group.

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

The Bio-Rad Mumps IgG test is intended for the qualitative and semi-quantitative detection of human IgG antibodies to mumps virus in human serum by enzyme immunoassay. Individual serum specimens may be used for the determination of immune status. Paired (acute/convalescent) sera may be used to demonstrate sero-conversion 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 Mumps
- Monitor IgG levels to identify active infection

**BIO-RAD**



# Mumps IgG EIA Testing

## Mumps IgG Indirect EIA Procedure

1



- Microplate: breakable wells dry coated with Ag

2



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

3

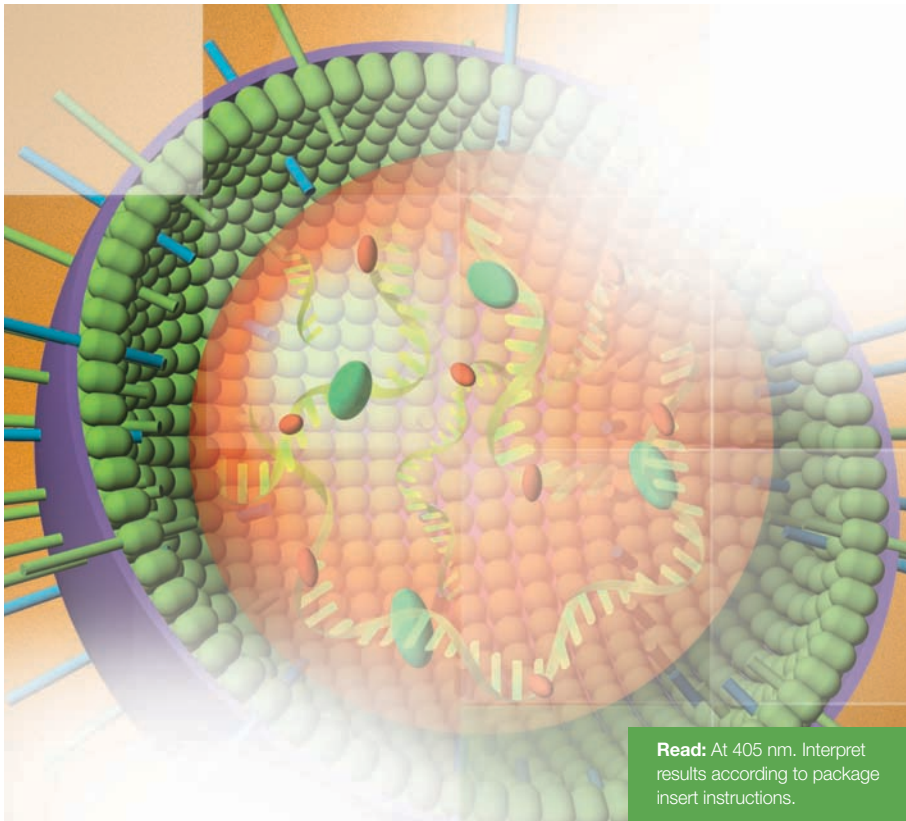


- Dispense 100  $\mu$ L of ready-to-use conjugate
- Incubate 30 minutes at room temperature
- 4 washes

4



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



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

## Ordering Information

Catalog No.	Description
25182	Mumps 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](http://www.bio-rad.com).

Clinical Diagnostics Group

Website [www.bio-rad.com](http://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 47-23-39-41-30 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