

Bio-Rad Laboratories
Improving lives with every test.

CMV

Cytomegalovirus IgG and IgM EIAs

Approximately 0.5 to 2.5% of newborn infants are identified each year with congenital Cytomegalovirus (CMV) infection. Fewer than 5% of congenitally-infected infants develop severe disease. The immune response to CMV involves synthesis of IgM antibodies several weeks after primary infection and one week later, IgG antibodies appear.

CMV infection remains difficult to identify by symptoms alone, and laboratory diagnosis is needed to identify acute infections as well as to establish the immune status of high risk individuals. Primary CMV infection in adults is often manifested by a mononucleosis-like syndrome and is usually self-limiting. However, in neonates and immunocompromised individuals, CMV can be debilitating and life-threatening.

Bio-Rad CMV IgG and CMV IgM EIA's

The CMV IgG and CMV IgM enzyme immunoassays are for the qualitative detection of antibodies to cytomegalovirus.

The IgG assay is used with serum for diagnostic assessment of prior infection with CMV. The CMV IgM assay is primarily indicated for the detection of early or recent infection.

Convenient

- Ready-to-use reagents
- Rapid turn-around

Objective

- Microplate format
- Qualitative 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 Cytomegalovirus
- Monitor IgG levels to identify active infection

IgM EIA Testing


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


BIO-RAD





Cytomegalovirus EIA Testing

Cytomegalovirus Indirect EIA Procedure

1 
 • Microplate: breakable wells dry coated with Ag

2 
 • Pipette 100 µL 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 µL of ready-to-use conjugate
 • Incubate 30 minutes at room temperature
 • 4 washes

4 
 • Dispense 100 µL of ready-to-use chromogen/substrate
 • Incubate 30 minutes at room temperature
 • Add 100 µL of ready-to-use Stop Reagent



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

Ordering Information

Catalog No.	Description	Tests
25177	CMV IgG EIA, Microplate	96 tests
25178	CMV 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