



EBV

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

## EB NA IgG EIA and EB VCA IgG and IgM EIAs

**Epstein-Barr Virus (EBV) is a common human pathogen, that infects the majority of individuals sometime during their lifetime.**

Since its discovery in 1964, EBV has been etiologically implicated in an increasing number of human diseases, such as Infectious Mononucleosis (IM). EBV is classified as a member of the herpes virus family. All herpes viruses share the ability to establish a latent infection in their hosts.

The EBV laboratory diagnosis is generally confirmed by a positive heterophile antibody test result. Severity of the disease however, is not indicated by the heterophile antibody test. Heterophile negative IM occurs in 10–20 percent of adults and in an even greater percentage of children. Detection and identification of antibodies to EBV are necessary to fully ascertain the stage of EBV infection. Bio-Rad offers microplate EIA kits for EB NA IgG, EB VCA IgG and EB VCA IgM testing. Results are obtained after a 90-minute incubation.

### Convenient

- Ready-to-use reagents
- Rapid turnaround

### Objective

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

### Simple

- 96-test kit
- Indirect EIA

### IgG EIA Testing

- As an aid in the diagnosis of infectious mononucleosis

### IgM EIA Testing

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

**BIO-RAD**

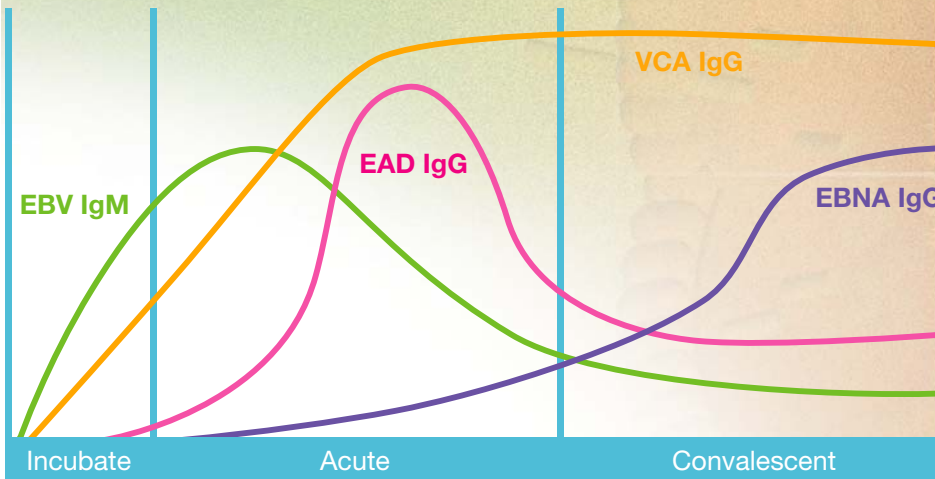


# Epstein-Barr Virus Testing





## Interpretation of Results – Same Criteria for EB NA IgG, EB VCA IgG or IgM

Index Value	Interpretation
<0.9	Negative for EBV antibody
≥1.1	Positive for EBV antibody
≥0.9 <1.1	Equivocal

## EBV Typical Antibody Response



## Epstein-Barr Virus Indirect EIA Procedure

- 
  - Microplate: breakable wells dry coated with Ag
- 
  - 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
- 
  - Dispense 100 µL of ready-to-use conjugate
  - Incubate 30 minutes at room temperature
  - 4 washes
- 
  - 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
25183	EB NA IgG EIA, Microplate..... 96 tests
25184	EB VCA IgG EIA, Microplate..... 96 tests
25185	EB VCA 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](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 47-23-38-41-30 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