



Bio-Rad Laboratories
Improving lives with every test.

VZV IgG

Varicella Zoster Virus (VZV) IgG

Varicella-zoster virus is the etiological agent of varicella and herpes zoster.

Varicella (chicken pox) is a mild, highly infectious disease, which occurs in hosts (usually children) without immunity. Zoster (shingles) is the recurrent form of the disease, occurring in adults previously infected with VZV.

An important aspect of the control of varicella-zoster virus infection is the prevention of contact between susceptible individuals and those with active infection. Antibodies to varicella-zoster have been detected by CF, immunodiffusion, neutralizing antibody, or fluorescent antibody techniques, including the fluorescent-antibody-membrane-antigen (FAMA) assay. The FAMA assay has proven to be very sensitive in determining low levels of VZV antibody and immune status. However, it has the disadvantage of requiring tissue culture, and is performed only in specialized laboratories. The performance of EIA procedures is comparable to the FAMA assay and more convenient.

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

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 Varicella-Zoster Virus
- Monitor IgG levels to identify active infection

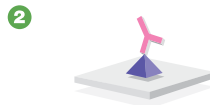
BIO-RAD

Varicella Zoster Virus IgG EIA Testing

Varicella Zoster Virus IgG Indirect EIA Procedure



- Microplate: breakable wells dry coated with Ag



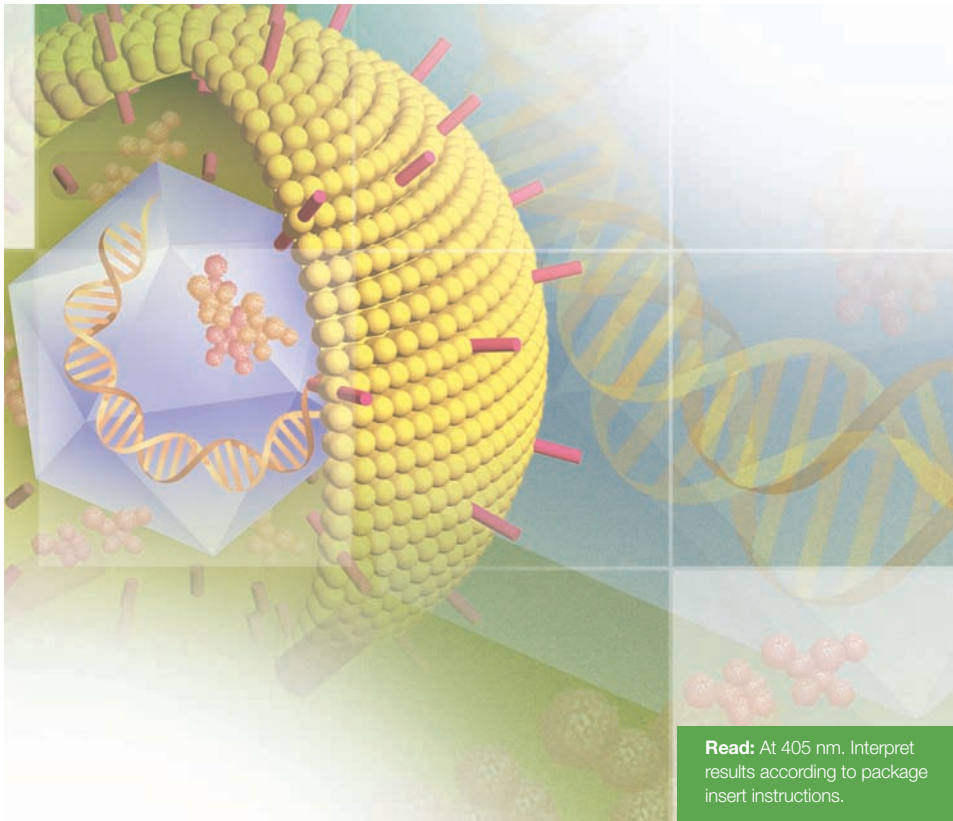
- Pipette 100 μ L of diluted test samples, calibrators or samples into the wells (1:51 for IgG)
- Incubate 30 minutes at room temperature
- 4 washes



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



- 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
25180	VZV IgG EIA..... 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 305-894-5960 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