

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

# Toxoplasmosis

## Toxo IgG & IgM EIAs

**Toxoplasmosis is a zoonotic infection (disease transmitted from animals to humans) by the protozoan *Toxoplasma gondii*. Cats are the predominant source of infection.**

Most *T. gondii* infections are asymptomatic or present with a transient febrile illness, often with lymphadenopathy. Primary infections acquired during pregnancy can cause stillbirth or congenital abnormalities, including blindness, encephalomyelitis, impaired brain functions and convulsions in the newborn. Infections can also be more severe in individuals who are infected with HIV or who are otherwise immunocompromised. Toxoplasmosis can be treated with antibiotics but early detection of the disease is crucial for effective therapy.

Diagnostic serology for Toxoplasmosis has progressed from the use of a specialized dye test to immunofluorescence, latex and enzyme immunoassay (EIA) techniques. In addition to Toxo IgG indirect EIA's, Bio-Rad offers both Toxo IgM immunocapture and indirect EIA tests.

### Convenient and Simple

- Rapid turnaround, 96-test kit
- Ready-to-use reagents

### Objective

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

### IgG EIA Testing

- As an aid in the assessment of the patient's immunological response to *Toxoplasma*
- Monitor IgG levels to identify active infection
- Indirect EIA testing

### IgM EIA Testing


- Estimate the time of infection by using IgM EIA in conjunction with the IgG EIA
- Platelia® Toxo IgM is an immunocapture EIA
- *Toxoplasma* IgM EIA is an indirect EIA

**BIO-RAD**

# Toxoplasmosis EIA Testing


**Platelia® Toxo IgG Indirect EIA Procedure**

**1**




- Microplate: breakable wells dry coated with Ag
- Wash microplate strips once

**2**




- Pipette 200 µL of diluted test samples and calibrators into the wells (1:101 for IgG)
- Incubate for 1 hour at 37°C
- 3 washes

**3**



- Dispense 200 µL of working conjugate
- Incubate for 1 hour at 37°C
- 4 washes

**4**




- Dispense 200 µL of chromogen/substrate
- Incubate 30 minutes at room temperature in the dark
- Add 100 µL of stopping solution

**Read:** At 450/620 nm. Interpret results according to package insert instructions.


**Platelia® Toxo IgM Immunoenzymatic Capture Method**

**1**




- Microplate: breakable wells dry coated with anti human mu-chain antibodies
- Wash microplate strips once

**2**




- Pipette 200 µL of diluted test samples and calibrators into the wells (1:101 for IgM)
- Incubate for 1 hour at 37°C
- 3 washes

**3**



- Dispense 200 µL of working conjugate
- Incubate for 1 hour at 37°C
- 4 washes

**4**




- Dispense 200 µL of chromogen/substrate
- Incubate 30 minutes at room temperature in the dark
- Add 100 µL of stopping solution

**Read:** At 450/620 nm. Interpret results according to package insert instructions.


**Toxoplasma IgG/IgM 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
25096	Platelia® Toxo IgG TMB, Microplate .....	96 tests
25095	Platelia® Toxo IgM TMB, Microplate.....	96 tests
25175	Toxoplasma IgG EIA, Microplate .....	96 tests
25176	Toxoplasma 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