

## TG MIC3 (residues 234-306)

### Toxoplasma Gondii Microneme Protein 3

#### *Toxoplasma gondii*, recombinant, *E. coli*

Cat. No.	Amount
PR-1243	100 µg

For *in vitro* use only  
Quality guaranteed for 12 months  
Store at -20°C

#### **Avoid freeze / thaw cycles**

#### **Form**

Liquid. Supplied in 50 mM Tris-HCl, pH 8.0, 1.5 M urea, and 50% glycerol.

#### **Specificity**

Immunoreactive with sera of *T. gondii*-infected individuals.

#### **Application**

Antigen in ELISA and Western blots, excellent antigen for detection of *Toxoplasma gondii* with minimal specificity problems.

#### **Purity**

>95% by SDS-PAGE

#### **Description**

The protein contains the *Toxoplasma gondii* MIC3 immunodominant regions, amino acids 234-306.

The protein is purified by proprietary chromatographic technique.

#### **Background**

*Toxoplasma gondii* is an obligate intracellular protozoan parasite that infects all warmblooded animals, including humans, and causes toxoplasmosis.

The homodimeric micronemal protein MIC3, which is a potent adhesin molecule of *T. gondii*, could be a significant candidate vaccine against toxoplasmosis.

#### **Selected References:**

Ismael et al. (2003) The MIC3 gene of *Toxoplasma gondii* is a novel potent vaccine candidate against toxoplasmosis. *Infect. Immun.*

**71**:6222.

Cerede et al. (2002) The *Toxoplasma gondii* protein MIC3 requires pro-peptide cleavage and dimerization to function as adhesin. *EMBO J.*

**21**:2526.

Garcia-Reguet et al. (2000) The microneme protein MIC3 of *Toxoplasma gondii* is a secretory adhesin that binds to both the surface of the host cells and the surface of the parasite. *Cell. Microbiol.* **2**:353.