

**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!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Purity: > 95 % (SDS-PAGE)

Form: liquid (Supplied in 50 mM Tris-HCl pH 8.0, 1.5 M urea and 50% glycerol)

Applications:

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

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.

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

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.