

TG MIC3 (residues 234-306) Toxoplasma Gondii Microneme Protein 3 *Toxoplasma gondii*, recombinant, *E. coli*

Cat. No.	Amount
PR-1243-1	1 mg

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.