

TG SAG1/p30 (residues 45-198) (Toxoplasma Gondii Surface Antigen 30) *Toxoplasma gondii*, Recombinant, *E. coli*

| Cat. No. | Amount |
|----------|--------|
| PR-1246 | 100 µg |

Liquid. Supplied as 1 mg/ml solution containing 50 mM Tris-HCl, pH 8.0, 1.5 M urea, and 50% glycerol.

Description: The protein contains the p30 (SAG1) immunodominant regions, amino acids 45-198.

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. SAG1, the major surface molecule of *Toxoplasma gondii*, is an important attachment ligand for the host cell. It is a highly immunogenic protein which has generated great interest as a diagnostic reagent, as a potential subunit vaccine, and for its role in invasion.

AVOID FREEZE/THAW CYCLES!

For in vitro use only!

Purity: >95% by SDS-PAGE.

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.

Store: -20 °C

Selected References:

Hiszczynska-Sawicka *et al.* (2003) High yield expression and singlestep purification of *Toxoplasma gondii* SAG1, GRA1, and GRA7 antigens in *Escherichia coli*. *Protein Expr. Purif.* **27**:150.

Robinson *et al.* (2004) *Toxoplasma gondii* major surface antigen (SAG1): in vitro analysis of host cell binding. *Parasitology.* **128**:391.

Seng *et al.* (2004) SAG1 is a host-targeted antigen for protection against *toxoplasma gondii* infection. *Pathobiology.* **71**:144.

Chen *et al.* (2003) Induction of immune responses in mice by vaccination with Liposome-entrapped DNA complexes encoding *Toxoplasma gondii* SAG1 and ROP1 genes. *Chin. Med. J. (Engl)* **116**:1561.

Sager *et al.* (2003) Immunodiagnosis of primary *Toxoplasma gondii* infection in sheep by the use of a P30 IgG avidity ELISA. *Parasitol. Res.* **91**:171.

Mohamed *et al.* (2003) Induction of protective immunity by DNA vaccination with *Toxoplasma gondii* HSP70, HSP30 and SAG1 genes. *Vaccine.* **21**:2852.