

TG SAG1/p30 (residues 45-198)

Toxoplasma Gondii Surface Antigen 30

Toxoplasma gondii, recombinant, *E. coli*

Cat. No.	Amount
PR-1246	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.

Application

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

Specificity

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

Purity

>95% by SDS-PAGE

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.

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.