

## TG SAG1/p30 (residues 45-198)

### Toxoplasma Gondii Surface Antigen 30

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

| Cat. No.  | Amount |
|-----------|--------|
| PR-1246-1 | 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.

#### **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.