

TG GRA7/p29 (residues 24-100) *Toxoplasma Gondii* Dense Granule Protein 7 *Toxoplasma gondii*, recombinant, *E. coli*

Cat. No.	Amount
PR-1245-1	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 p29 (GRA7) immunodominant regions, amino acids 24-100.

The protein is purified by proprietary chromatographic technique.

Background

Toxoplasma gondii is an obligate intracellular protozoan parasite that infects all warm-blooded animals, including humans, and causes toxoplasmosis.

GRA7 is a recently discovered 29 kDa protein. Like GRA1, it is secreted from the dense granules, and it reacts with sera from humans with acute and chronic infections.

Selected References:

- Hiszczynska-Sawicka *et al.* (2003) High yield expression and single-step purification of *Toxoplasma gondii* SAG1, GRA1, and GRA7 antigens in *Escherichia coli*. *Protein Expr. Purif.* **27**:150.
- Neudeck *et al.* (2002) Expression variance, biochemical and immunological properties of *Toxoplasma gondii* dense granule protein GRA7. *Microbes Infect.* **4**:581.
- Vercammen *et al.* (2000) DNA vaccination with genes encoding *Toxoplasma gondii* antigens GRA1, GRA7, and ROP2 induces partially protective immunity against lethal challenge in mice. *Infect. Immun.* **68**:38.
- Ferguson *et al.* (1999) *In vivo* expression and distribution of dense granule protein 7 (GRA7) in the exoenteric (tachyzoite, bradyzoite) and enteric (coccidian) forms of *Toxoplasma gondii*. *Parasitology.* **119**:259.
- Jacobs *et al.* (1999) Evaluation of recombinant dense granule antigen 7 (GRA7) of *Toxoplasma gondii* for detection of immunoglobulin G antibodies and analysis of a major antigenic domain. *Clin. Diagn. Lab. Immunol.* **6**:24.
- Bonhomme *et al.* (1998) Quantitative immunolocalization of a P29 protein (GRA7), a new antigen of *Toxoplasma gondii*. *J. Histochem. Cytochem.* **46**:1411.