

**Toxoplasma gondii Antigen**

Strain RH  
native, Vero cell culture

Cat. No.	Amount
PR-BA110-S	250 µg
PR-BA110-L	1 mg

**For general laboratory use.**

**Shipping:** shipped on dry ice

**Storage Conditions:** store at -80 °C

**Additional Storage Conditions:** avoid freeze/thaw cycles

**Shelf Life:** 12 months

**Form:** liquid

**Culture system:** Vero cell culture

**Protein concentration:** Lot specific. Determined by Bradford assay (Biorad) using BSA as standard.

**Preservation:** None

**Analysis:** ELISA

**Preparation:** Disrupted Toxoplasma 'ghosts' were treated with detergent and further purified. The antigen is presented in cell culture medium containing n-octylglucopyranoside.

**Inactivation:** This product has been inactivated by disruption and detergent treatment. Confirmed inactive by attempted growth under original culture conditions. Since no procedure can guarantee absolute sterility, the reagent should be handled with appropriate safety precautions.