

**Rubella Virus Antigen**

Strain Judith  
native, Vero cell culture

Cat. No.	Amount
PR-BA129G2-S	250 µg
PR-BA129G2-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

**Culture system:** Vero cell culture

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

**Preservation:** None

**Analysis:** ELISA

**Preparation:** Infected cells and culture supernatant are harvested, concentrated, purified by sucrose cushion centrifugation and Tween-ether treated.

**Inactivation:** This product has been inactivated by Tween-ether 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.