

**Rubella Virus Antigen**

Strain Judith
native, Vero cell culture

Cat. No.	Amount
PR-BA129G2-S	250 µg
PR-BA129G2-L	1 mg

For *in vitro* use only!

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.