

**Respiratory Syncytial Virus (RSV) Antigen**

Strain Long  
native, Vero cell culture

Cat. No.	Amount
PR-BA113-S	250 µg
PR-BA113-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

**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:** RSV infected cells are harvested and virus is extracted by alkaline treatment and sonication. The suspension is purified by centrifugation. The resulting antigen preparation consists of virus and viral components and residual host cell material. The antigen is presented in 100 mM Glycine buffer at pH 9.5.

**Inactivation:** This product has been treated with Beta-propiolactone. 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.