

**Coxsackievirus B6 Antigen**

Strain Schmitt  
native, Vero cell culture

Cat. No.	Amount
PR-BA134B6-S	250 µg
PR-BA134B6-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:** Infected cells are disrupted and clarified to remove cell debris. Virus is concentrated by ultracentrifugation and presented in cell culture medium free of fetal calf serum.

**Inactivation:** This product has been inactivated by heat treatment at 56°C. 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.