

**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.