

**Echo Virus Antigen**recombinant, *E. coli*

Cat. No.	Amount
PR-BA135VS-S	250 µg
PR-BA135VS-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**Form:** liquid**Protein concentration:** Lot specific. Determined by Bradford assay (Biorad) using BSA as standard.**Presentation:** The antigen is presented in buffer containing 39 mM Tris/HCl pH 8.0, 150 mM NaCl, 10 % glycerol.**Analysis:** SDS-Page**Purification:** Chromatography