

**Parvovirus B19 VP1**recombinant, *E. coli*

Cat. No.	Amount
PR-BA122VP1-S	250 µg
PR-BA122VP1-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**Molecular Weight:** 87 kDa**Form:** liquid**Concentration:** Lot specific, specified in mg/ml. Determined by Bradford assay (Biorad) using BSA as standard.**Description:**

Parvovirus B19 structural protein VP1.

**Purification:**

Chromatography

**Presentation:**

The protein is presented in buffer containing 50 mM Tris/HCl pH 8.0, 500 mM NaCl and 8 M Urea.

**Analysis:**

SDS-PAGE and ELISA