

**Parvovirus B19 VP2 C-term**recombinant, *E. coli*

Cat. No.	Amount
PR-BA122VP1B-S	250 µg
PR-BA122VP1B-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**Molecular Weight:** 35 kDa**Form:** liquid**Concentration:** Lot specific, specified in mg/ml. Determined by Bradford assay (Biorad) using BSA as standard.**Description:**

The protein contains the C-terminal fragment of parvovirus B19 structural protein VP2.

Purification:

Chromatography

Presentation:

The protein is presented in buffer containing PBS pH 7.3 and 4 M Urea.

Analysis:

SDS-PAGE and ELISA