
SAP18 Protein

(Catalog No. E20000)

Description

Human SAP18 recombinant protein containing 153 amino acids (aa 20- 172) with His tag at N-terminus

Molecular Weight (kDa)

19.7 kDa

Formulation

20mM Tris-HCl buffer (pH8.0), 0.1M NaCl, 30% glycerol, 1mM DTT

Source

Escherichia Coli.

Size

20 µg

Storage Temp & Conditions

6 months at -80°C. Please prevent freeze-thaw cycles. Best used within three months from date of receipt.

Purification Notes

Purity : 90.0%

Application Notes

Enzyme-linked Immunoabsorbent Assay; Western Blot (Recombinant protein); Antibody Production; Protein Array; Preparation Method: in vitro wheat germ expression system.

Ordering Information:

Products	Size	Cat. No.
SAP18 Protein	20 µg	E20000-1

This product is for research purposes only. Not intended for use in diagnostic procedures.