
SUMO-1 Protein

(Catalog No. E28000)

Description

All SUMO isoforms are translated with additional C-terminal residues that have to be removed to generate the active protein. Pro-SUMO-1 (101 amino acids) is the inactive precursor of SUMO-1 (97 amino acids) and is processed at the C-terminus by SUMO-1 specific proteases (SENPs). The resulting SUMO-1 protein has the conserved C-terminal Gly-Gly residues that function in activation and conjugation reactions. This protein can be used as a negative control in sumoylation reactions or as a substrate for SENPs.

Molecular Weight (kDa)

11.6 kDa

Formulation

50 mM HEPES pH 8.0, 150 mM NaCl, 1mM DTT, 10% glycerol

Source

Expressed in E.coli expression system

Size

100 µg

Storage Temp & Conditions

Store at -80°C. Avoid multiple freeze/thaw cycles. Best use within three months from date of receipt.

Purification Notes

Purity: >95% by HPLC

Application Notes

Enzyme-linked Immunoabsorbent Assay; Western Blot (Recombinant protein); Antibody Production; Protein Array.

Ordering Information:

Products	Size	Cat. No.
SUMO-1 Protein	100 µg	E28000 -1

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