
SUMO-2 Protein

(Catalog No. E28002)

Description

All SUMO isoforms are translated with additional C-terminal residues that have to be removed to generate the active protein. Pro-SUMO-2 (95 amino acids) is the inactive precursor of SUMO-2 (93 amino acids) and is processed at the C-terminus by SUMO-2 specific proteases (SENPs). The resulting SUMO-2 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)

10.9 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.

Purification Notes

Purity: >95% by SDS-PAGE

Application Notes

Typical in vitro concentrations are 10-50 µM depending on conditions

Ordering Information:

Products	Size	Cat. No.
SUMO-2 Protein	100 µg	E28002-1

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