

SUMO3 Protein

Recombinant human protein expressed in E.coli cells

Catalog # S295-31H

Lot # F722-2

Product Description

Recombinant human SUMO3 (1-92) was expressed in E. coli cells using an N-terminal His tag. The gene accession number is [NM_006936](#).

Gene Aliases

SMT3A, SMT3H1, SUMO-3, Smt3B

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

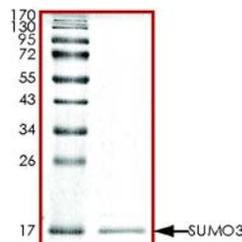
Scientific Background

SUMO3 is a member of the SUMO (small ubiquitin-like modifier) protein family which functions in a manner similar to ubiquitin. SUMO3 is bound to target proteins as part of a post-translational modification system and modifies numerous cellular proteins and affect their metabolism and function. SUMO3 is involved in a variety of cellular processes, such as nuclear transport, transcriptional regulation, apoptosis, and protein stability (1). SUMO3 predominantly detected in the cytoplasm (2).

References

1. Su, H.-L. et.al: Molecular features of human ubiquitin-like SUMO genes and their encoded proteins. *Gene* 296: 65-73, 2002.
2. Lapenta, V.et.al: SMT3A, a human homologue of the *S. cerevisiae* SMT3 gene, maps to chromosome 21qter and defines a novel gene family. *Genomics* 40: 362-366, 1997.

Purity



The purity of SUMO3 was determined to be >95% by densitometry. Approx. MW 17kDa.

SUMO3 Protein

Recombinant human protein expressed in E. coli cells

Catalog #	S295-31H
Lot #	F722-2
Purity	>95%
Concentration	0.5 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.