

SUMO1 (UBL1) Protein

Recombinant human protein expressed in E.coli cells

Catalog # S293-31H

Lot # F722-1

Product Description

Recombinant human SUMO1 (UBL1) (2-97) was expressed in E. coli cells using an N-terminal His tag. The gene accession number is [NM_003352](#).

Gene Aliases

UBL1; DAP1; GMP1; SMT3; OFC10; SENP2; SMT3C; SMT3H3; PIC1

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.

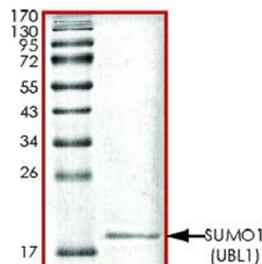
Scientific Background

SUMO1 (also known as UBL1) is a member of the SUMO (small ubiquitin-like modifier) protein family which functions in a manner similar to ubiquitin. SUMO1 is bound to target proteins as part of a post-translational modification system and it is involved in a variety of cellular processes, such as nuclear transport, transcriptional regulation, apoptosis, and protein stability (1). SUMO1 may be involved in the repair of TOP1-mediated DNA damage (2). The sumoylation pathway plays a significant role in mammalian DNA damage response.

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. Mao, Y. et.al: SUMO-1 conjugation to topoisomerase I: a possible repair response to topoisomerase-mediated DNA damage. *Proc. Nat. Acad. Sci.* 97: 4046-4051, 2000.

Purity



The purity of SUMO1 (UBL1) was determined to be >90% by densitometry. Approx. MW **19kDa**.

SUMO1 (UBL1) Protein

Recombinant human protein expressed in E. coli cells

Catalog #	S293-31H
Lot #	F722-1
Purity	>90%
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.