

Anti-SUMO1

Rabbit Polyclonal Antibody

Catalog # S293-63R

Lot # O2121-1

Cited Applications

WB, ELISA, IF, IHC

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the SUMO1 protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

SUMO1 antibody was raised against a 14 amino acid synthetic peptide from near the amino terminus of human SUMO1

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

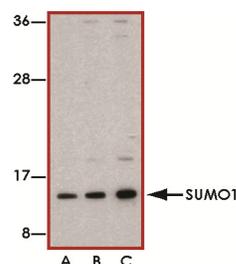
Scientific Background

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.

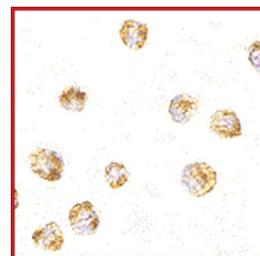
References

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.

Sample Data



Western blot analysis of SUMO1 in HL-60 cell lysate with SUMO1 antibody at (A) 0.5, (B) 1, and (C) 2 µg/ml.



Immunocytochemistry of SUMO1 in HL60 cells with SUMO1 antibody at 5 µg/ml.

Anti-SUMO1

Rabbit Polyclonal Antibody

Catalog Number

S293-63R

Specific Lot Number

O2121-1

Purification
Stability
Storage & Shipping

Affinity chromatography
1yr at -20°C from date of shipment
Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

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.