

## Anti-SUMO3

Rabbit Polyclonal Antibody

**Catalog # S295-63R**

Lot # O2121-32

### Cited Applications

WB, ELISA, IF, IHC

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

### Specificity

Recognizes the SUMO3 protein

### Cross Reactivity

Human, Mouse and Rat

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

SUMO3 antibody was raised against a 13 amino acid synthetic peptide near the carboxy terminus of human SUMO3.

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

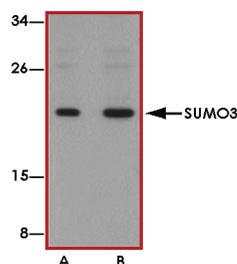
### Scientific Background

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).

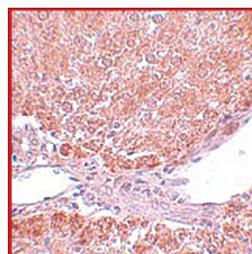
### References

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.

### Sample Data



Western blot analysis of SUMO3 in mouse liver tissue lysate with SUMO3 antibody at (A) 1 µg/ml and (B) 2 µg/ml.



Immunohistochemistry of SUMO2/3 in rat liver tissue with SUMO2/3 antibody at 5 µg/ml.

## Anti-SUMO3

Rabbit Polyclonal Antibody

Catalog Number

**S295-63R**

Specific Lot Number

**O2121-32**

Purification Affinity chromatography

Stability 1yr at -20°C from date of shipment

Storage & Shipping

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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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