

# pro-SUMO-1 (human), (recombinant) (His- tag)

SUMO family members are expressed in their pro-forms with short C-terminal extensions, which are processed to expose the C-terminal glycine residue that is essential for conjugation to target proteins.

**Manuals, SDS & CofA**

[View Online »](#)

## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

### Application Notes

### Uses:

1. Study of SUMO-1 regulation and processing.
2. Substrate for deSUMOylating enzymes (C-terminal hydrolase activity).

### Formulation

Liquid. In 20mM HEPES, pH 8.0, 50mM NaCl, 1mM DTT.

### MW

16.4 kDa

### Purity

≥95% (SDS-PAGE)

### Source

Produced in *E. coli*. Human pro-SUMO-1 is fused at the N-terminus to a His-tag.

### UniProt ID

P63165

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)