

SUMO-3 (human), (recombinant) (His- tag)

Mature protein for functional studies.

Conjugation of mature SUMO proteins to specific lysine residues on target proteins via isopeptide bonds, is involved in a range of processes including nuclear protein targeting, formation of sub-nuclear complexes, regulation of transcriptional activities, and control of protein stability.

Citations: 1

[View Online »](#)

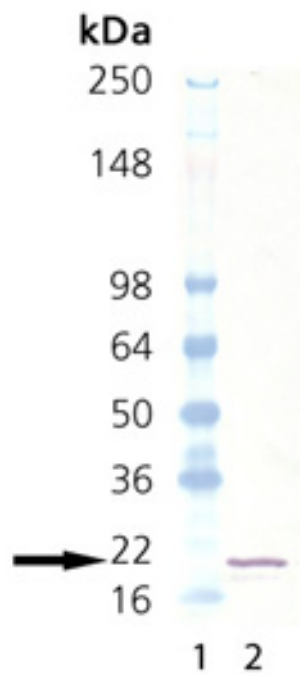
Ordering Information

[Order Online »](#)

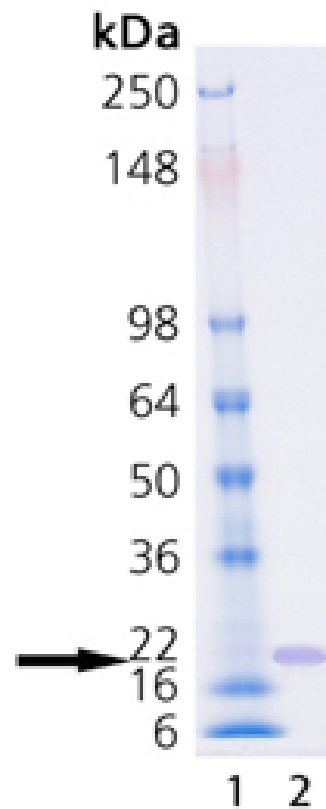
BML-UW9215-0500	500µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis: Lane 1: MW marker, Lane 2: SUMO-3 probed with SUMO-2/3 pAb.



SDS-PAGE analysis: Lane 1: MWM, Lane 2: 1 μ g SUMO-3 recombinant protein.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name

Sentrin-3, Small ubiquitin-related modifier 3, SUMO-2

Application Notes

Uses:

1. SUMO-3 modification of specific proteins in vitro.
2. Demonstrate novel proteins are potential targets for SUMOylation under in vitro conditions.
3. Generate substrates for deSUMOylating enzymes.

Formulation

Liquid. In 20mM HEPES, pH 8.0, containing 50mM sodium chloride and 1mM dithiothreitol.

MW

15.5 kDa

Purity

≥95% (SDS-PAGE)

Quality Control

Ability to SUMOylate RanGAP substrate (Prod. No. BML-UW9755) *in vitro*.

Source

Produced in *E. coli*. Human SUMO-3 is fused to an N-terminal His-tag.

UniProt ID

P55854

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-be@enzolifesciences.com)

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