

SUMO-2 (human) (1-93), (recombinant)

Active form of SUMO-2

Active form of human SUMO-2 for conjugation to protein substrates.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-089-C500	500µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability Stable for a few days at room temperature.

Long Term Storage -80°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

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

Application Notes Standard reaction for protein conjugation (10µl) includes the target protein, SUMO-1 activating enzyme (Prod. No. BML-UW9330) (100ng), UbcH9 (900ng) (Prod. No. ALX-201-046), SUMO-2 (1-93) (10µg) in 50mM TRIS, pH 7.6, 5mM magnesium chloride, 2mM ATP, 10mM creatine phosphate, 3.5U/ml of creatine kinase and 0.6U/ml of inorganic pyrophosphatase.

Formulation Liquid. In PBS.

Purity ≥95%

Source Produced in *E. coli*.

Technical Info / Product Notes Formerly this product was called SUMO-3 (1-93) (human) (recombinant), but the name was changed in accordance with UniProt.

UniProt ID P61956



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-
fr@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)