

SUMOstar Protease

Recombinant SUMOstar Protease 1 specifically recognizes and cleaves the SUMOstar tag from recombinant fusion proteins made with SUMOstar expression systems (Cat.#s: 1100A, 1106, 1100K, 1101, 2100, 2101, 2160, 2161, 3100, 3101, 7110, 7111, 7120, 7121). Unlike thrombin, EK, or TEV proteases, which recognize short, linear sequences, SUMOstar Protease 1 recognizes the tertiary structure of SUMOstar. As a result, SUMOstar Protease 1 will not cleave within the protein of interest. SUMOstar Protease 1 will not cleave SUMO-tags from proteins expressed with the SUMOpro or SUMOpro3 systems. To cleave the SUMOpro tag (original yeast SUMO tag), please use SUMO Protease 1 and to cleave SUMOpro3 tag please use SUMO Protease 2.

Ordering Information

[Order Online »](#)

LSI-SP-4110-0250	250U
LSI-SP-4110-0500	500U
LSI-SP-4110-1000	1000U
LSI-SP-4110-5000	5000U

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping

Dry Ice

Regulatory Status RUO - Research Use Only

Product Details



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)