

# 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 »](#)

<b>LSI-SP-4110-0500</b>	500U
<b>LSI-SP-4110-1000</b>	1000U
<b>LSI-SP-4110-5000</b>	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)