

SUMOstar Protease

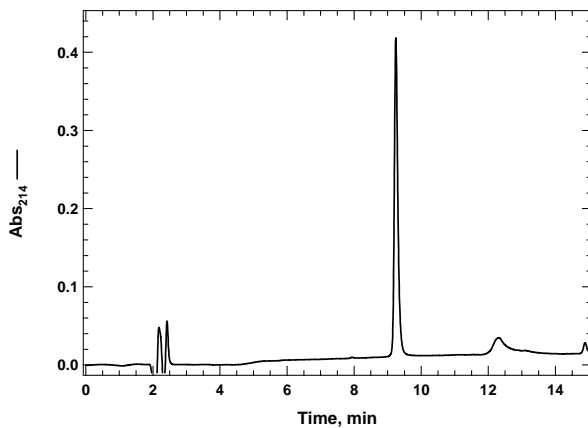
Cat. # 4110

Background: LifeSensors' SUMOstar Protease is a highly active and robust enzyme genetically engineered from yeast Ulp1 (see Cat. # 4010L). SUMOstar, derived from yeast SUMO, was engineered to be resistant to cleavage by endogenous desumoylases to allow expression of SUMOstar fusion proteins in eukaryotic expression systems. SUMOstar Protease was generated to specifically remove SUMOstar. Like SUMO Protease 1, SUMOstar Protease recognizes the tertiary structure of SUMOstar and as a result, never cleaves within the protein of interest. SUMOstar Protease acts over a broad range of temperature (30°C is optimal), pH [5.5 – 9.5], and ionic strength. SUMOstar Protease contains a His₆ tag at the N-terminus, making it easy to remove by affinity chromatography.

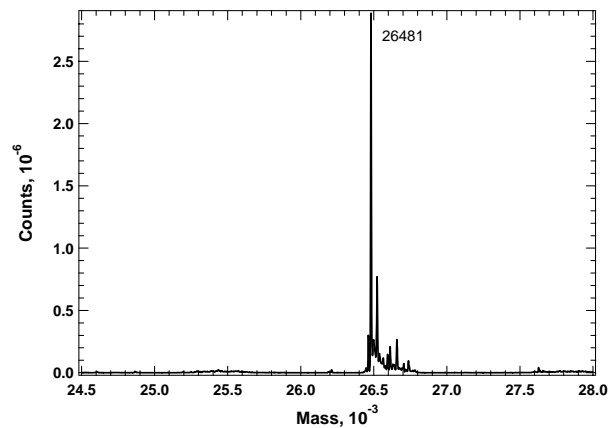
Application: Removal of SUMOstar tags from recombinant, fusion proteins.

Product Information

Purity:	≥ 95% by SDS-PAGE and RP-HPLC
Molecular Weight:	26,481.3Da
Specific Activity:	1 U will cleave >90 µg of SUMOstar-GFP in 1 hr at 37° C
Physical State:	Liquid, 10U/µL
Quantity:	500, 1000, 5000, or 10,000 Units
Storage:	-80° C. Avoid repeated freeze/thaw cycles



RP-HPLC



Deconvoluted mass spectrum

References

1. Li, S.-J. and M. Hochstrasser. 1999. A new protease required for cell cycle progression in yeast. *Nature* **398**,246-251.
2. Li, S.-J., W. Hankey, and M. Hochstrasser. 2005. Preparation and characterization of yeast and human desumoylating enzymes. *Meth Enzymol* **398**,457-467.
3. Malakhov, M.P., M.R. Mattern, O.A. Malakhova, M. Drinker, S.D. Weeks, and T.R. Butt. 2004. SUMO fusions and SUMO-specific protease for efficient expression and purification of proteins. *J Struct Funct Genomics* **5**,75-86.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2009 LifeSensors, Inc. All Rights Reserved