

SUMO Protease 1

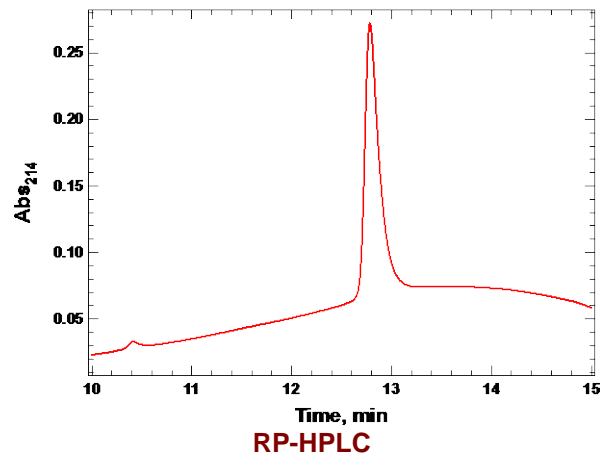
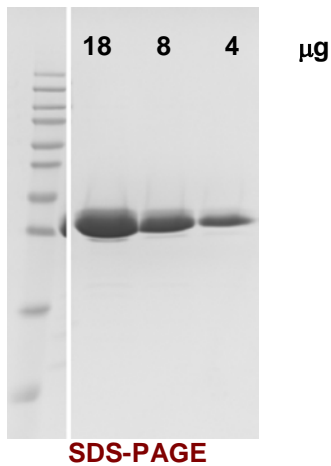
Cat. # 4010

Background: LifeSensors' industry leading SUMO Protease 1 (Ulp1), a highly active and robust recombinant protease, cleaves SUMO from recombinant fusion proteins. Unlike thrombin, EK, or TEV protease, which recognize short linear sequences, SUMO Protease 1 recognizes the tertiary structure of SUMO. As a result, SUMO Protease 1 never cleaves within the fused protein of interest. SUMO Protease 1 cleaves consistently over a broad range of temperature (30°C is optimal), pH [5.5 – 9.5], and ionic strength. SUMO Protease 1 contains a polyhistidine tag at the N-terminus, making it easy to remove from the cleavage reaction by affinity chromatography.

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

Product Information

Purity:	≥ 95% by SDS-PAGE and RP-HPLC
Molecular Weight:	26595.9Da by MS (26595.4 expected)
Specific Activity:	≥ 1 x 10 ⁵ U/mg, 1 U will cleave >90 µg of SUMO-GFP in 1 hr at 37° C
Physical State:	10U/µl in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 5mM DTT, 10% Glycerol
Quantity:	250, 500, 1000, or 5000 Units
Storage:	-80° C. Avoid repeated freeze/thaw cycles



All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
 Copyright © 2013 LifeSensors, Inc. All Rights Reserved

Protocol

1. After SUMO-protein fusion is purified: dialyze sample against proper buffer (e.g. PBS, pH 7.4 or 20mM TRIS, pH 8.0 containing 150 mM NaCl) at 4°C
2. Add SUMO Protease 1 to substrate (1 unit enzyme to 10-100 µg substrate should suffice, depends on SUMO fusion protein); add DTT to final 2 mM
3. Either:
 - a. incubate the mixture at 30°C for 1 h (mix gently do not vortex), or
 - b. incubate the mixture at 4°C overnight
 - c. (you can also perform a. followed by b.)
4. Check the cleavage using SDS-PAGE. If the SUMO-fusion is not cleaved up to 95% add more SUMO protease 1.

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. Mossessova, E. and C.D. Lima. 2000. Ulp1-SUMO crystal structure and genetic analysis reveal conserved interactions and a regulatory element essential for cell growth in yeast. *Mol Cell* **5**,865-
4. 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.