

SUMO Protease 2

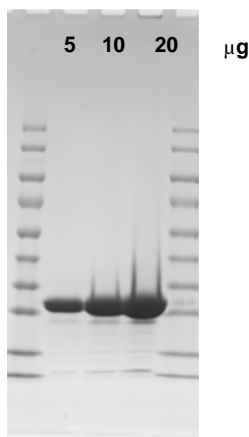
Cat. # 4020

Background: LifeSensors' industry leading SUMO Protease 2 (Senp2), a highly active and robust recombinant protease, cleaves hSUMO3 from recombinant fusion proteins. Unlike thrombin, EK, or TEV protease, which recognize short linear sequences, SUMO Protease 2 recognizes the tertiary structure of SUMO. As a result, SUMO Protease 2 will not cleave within the fused protein of interest. SUMO Protease 2 cleaves consistently over a broad range of temperature (30°C is optimal), pH [5.5 – 9.5], and ionic strength. SUMO Protease 2 contains a polyhistidine tag at the N-terminus, making it easy to remove from the cleavage reaction by affinity chromatography.

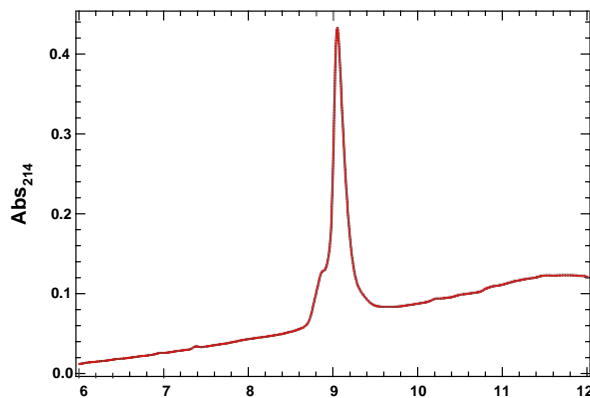
Application: Removal of hSUMO tags from recombinant, fusion proteins.

Product Information

Purity:	≥ 95% by SDS-PAGE and RP-HPLC
Molecular Weight:	27739.2 Da by MS (27739.1 expected)
Specific Activity:	≥ 1 x 10 ⁵ U/mg, 1 U will cleave >90 µg of hSUMO3-GFP in 1 hr at 37° C
Physical State:	10U/µl in 50 mM Hepes, pH 7.5, 150 mM NaCl, 10% glycerol
Quantity:	250, 500, 1000, or 5000 Units
Storage:	-80° C. Avoid repeated freeze/thaw cycles



SDS-PAGE



RP-HPLC

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 SUMO3-protein fusion is purified: dialyze sample against proper buffer (e.g. PBS, pH 7.4 or 20 mM TRIS buffer containing 150 mM NaCl, pH 8.0) at 4°C.
2. Add SUMO Protease 2 to substrate (1 unit enzyme to 10-100 µg substrate should suffice, depends on SUMO3 fusion protein); add DTT to 2 mM final.
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
 - i. (you can also perform a. followed by b.)
4. Check the cleavage using SDS-PAGE. If the SUMO3-fusion is not cleaved up to 95%, add more SUMO protease 2.

References

1. Marblestone JG, Edavettal SC, Lim Y, Lim P, Zuo X, Butt TR (2006). "Comparison of SUMO fusion technology with traditional gene fusion systems: Enhanced expression and solubility with SUMO," *Protein Science*, 15:182-9.
2. Butt TR, Edavettal SC, Hall JP, Mattern MR (2005). "SUMO fusion technology for difficult-to-express proteins," *Protein Expr Purif.*, 43(1):1-9.
3. Zuo X, Li S, Hall J, Mattern MR, Tran H, Shoo J, Tan R, Weiss SR, Butt TR (2005). "Enhanced expression and purification of membrane proteins by SUMO fusion in *Escherichia coli.*," *J Struct Funct Gen.*, 6:103-11.
4. Malakhov MP, Mattern MR, Malakhov OA, Drinker M, Weeks SD, Butt TR (2004). "SUMO fusions and SUMO-specific protease for efficient expression and purification of proteins," *J Struct Funct Gen*, 5:75-86.