

SUMO-GFP

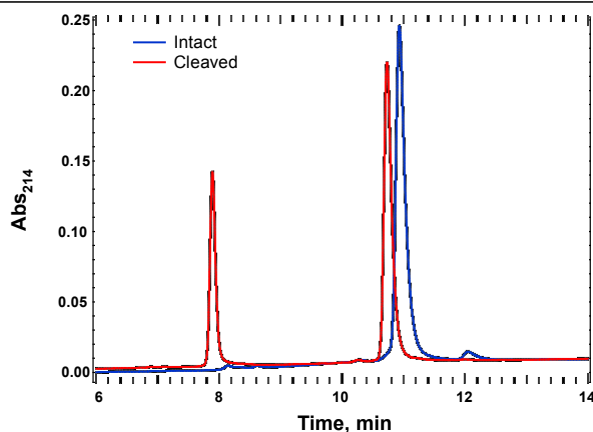
Cat. # 5000

Background: SUMO-GFP is a recombinant protein that contains yeast SUMO fused to a control polypeptide (GFP). This fusion protein is used as a positive control for SUMO Protease 1 activity. Uncleaved SUMO-GFP migrates on SDS-PAGE with an apparent MW of ~45 kDa while SUMO runs at 18-20 kDa and GFP runs at ~25 kDa following cleavage.

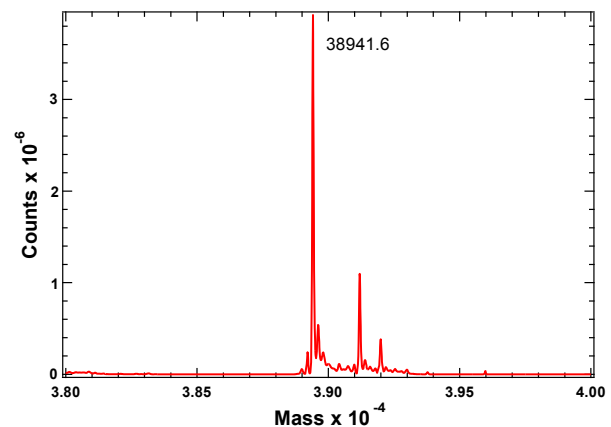
Application: Determination of the activity of SUMO Protease 1.

Product Information

Purity:	≥ 90% by RP-HPLC and SDS-PAGE
Molecular Weight:	38941.6 Da by MS (expected 38940.9 Da)
Physical State:	5mg/ml in TBS pH 7.5, 10% glycerol
Quantity:	100 µg
Solubility:	≥ 5 mg/mL
Storage:	-80° C. Avoid repeated freeze/thaw cycles



RP-HPLC



Deconvoluted mass spectrum

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**,1-9.
3. 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.

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