

SUMOstar-GFP

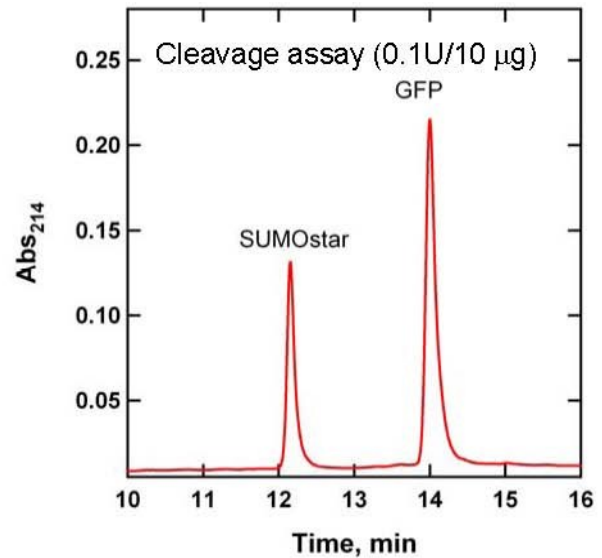
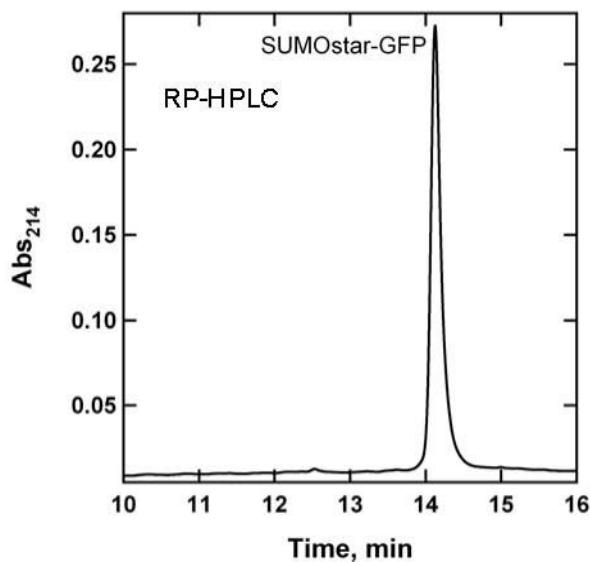
Cat. # 5010

Background: SUMOstar-GFP is a recombinant fusion protein that contains a modified yeast SUMO fused to a control polypeptide (GFP).

Application: Determination of the activity of SUMOstar Protease 1.

Product Information

Purity:	≥ 90% by RP-HPLC and SDS-PAGE
Molecular Weight:	38881.7 Da by MS (expected 38888.8 Da)
Physical State:	liquid; 5mg/ml in TBS pH 7.5, 10% glycerol
Quantity:	100 µg
Solubility:	≥ 5 mg/mL in aqueous buffer
Storage:	-80° C. Avoid repeated freeze/thaw cycles



References

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