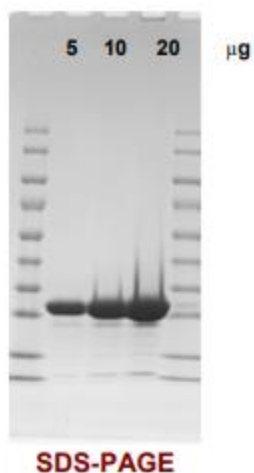


## Product Information

<b>Description:</b>	Active recombinant SENP2 (SUMO Protease 2). This construct contains an N-terminal His-tag. This protein was affinity purified.
<b>Background:</b>	SUMO Protease 2, 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 huSUMO3. As a result, SUMO Protease 2 will not cleave within the fused protein of interest.
<b>Species:</b>	Human
<b>Tag:</b>	His
<b>Concentration:</b>	10 U/μl
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	≥90% by SDS-PAGE and RP-HPLC
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	50 mM HEPES, pH 7.5, 150 mM NaCl, 10% glycerol
<b>MW:</b>	28 kDa
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Specific Activity</b>	> 1 x 10 <sup>5</sup> U/mg, 1 U will cleave >90 μg of hSUMO3-GFP in 1 hr at 37°C
<b>Assay Conditions:</b>	SUMO3-containing protein should be purified and dialyzed with PBS, pH 7.4 or 20 mM Tris Buffer and 150 mM NaCl pH 8. Incubation with Senp2 is then performed by adding 1 unit of Senp2 per each 10-100 μg of SUMO-3-containing protein and adding DTT to reach a 2 mM final concentration. Reactions can be incubated at 30°C for 1 hour with gentle agitation (no vortexing) or overnight at 4°C.
<b>Applications:</b>	Useful for removal of hSUMO3 from recombinant proteins over a range of temperature (30°C optimal) and ionic strength, and from pH 5.5 to 9.5.

## Quality Control Data

### SDS-PAGE Coomassie Staining



### RP-HPLC

