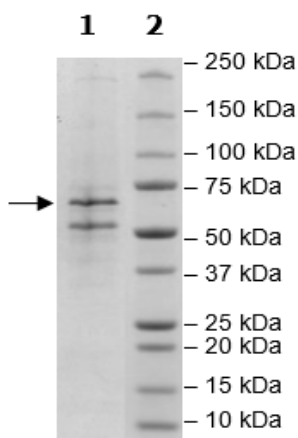


## Product Information

<b>Description:</b>	Recombinant human USP2 (ubiquitin specific peptidase 2), full length, encompassing amino acids 2-605(end). This construct contains an N-terminal Flag-tag. The recombinant protein was affinity purified.
<b>Species</b>	Human
<b>Construct:</b>	USP2 (Sf9-derived) (FLAG-2-605(end))
<b>Concentration:</b>	0.13 mg/ml
<b>Expression System:</b>	Sf9
<b>Purity:</b>	55%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 3 mM DTT
<b>MW:</b>	69 kDa
<b>Genbank Accession:</b>	NM_004205
<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>	2.9 pmol/min/μg
<b>Assay Conditions:</b>	Assay was done according to USP7 Inhibitor Screening Assay Kit (BPS Bioscience #79256) with USP2 (BPS Bioscience #80352) titrated at various concentrations.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



USP2 Activity

