

## USP30 (Ubiquitin Specific Protease 30)

Cat. # DB530

### Background

USP30 (Ubiquitin Specific Protease 30) is a deubiquitinase involved in the regulation of mitophagy. It is attached to the outer mitochondrial membrane by insertion of a transmembrane domain (residues 36-56). It removes ubiquitin from target proteins ubiquitinated by the E3 Ligase Parkin (PARK2). This form of USP30 has been engineered to increase its stability while retaining the cleavage preference for K6-linked ubiquitin chains over other chains. USP30 shows reduced efficiency to cleave phosphorylated ubiquitin chains.

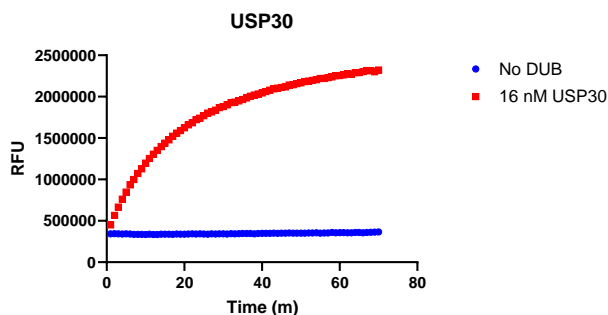
### Alternate names

Deubiquitinating enzyme 30, Ubiquitin thioesterase 30, Ubiquitin-specific-processing protease 30

### Product Information

<b>Purity</b>	≥ 90% by SDS-PAGE
<b>Molecular Weight</b>	51 kDa
<b>Quantity</b>	25, 50 µg
<b>Physical State:</b>	Liquid
<b>Tag:</b>	His6-SUMO
<b>Storage:</b>	-80° C. Avoid repeated freeze/thaw cycles

### Product QC



**Activity Assay of USP30.** 16 nM USP30 was tested in a CHOP assay showing robust DUB activity.

### References

1. Gu, L., et al., Hepatology, 2021. 73(1):160-174.
2. Qin, X., et al., Autophagy, 2022. 18(9):2178-2197.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2012 LifeSensors, Inc. All Rights Reserved