

## USP4 (Ubiquitin specific protease 4)

Cat. # DB511

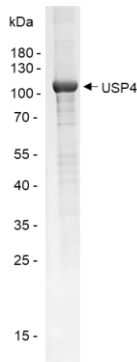
**Background** USP4 (Ubiquitin specific protease 4, UNP) has been shown cleave the ubiquitin precursor UBICEP52 and the isopeptide (isopeptide linked) ubiquitin conjugate ubiquitin-(εN)-lysine *in vitro*. It deubiquitinates PDPK 1 and TRIM21. USP4 may play a role in regulating mRNA splicing and quality control in the ER.

**Alternate names** Deubiquitinating enzyme 4, Ubiquitin carboxyl-terminal hydrolase 4, Ubiquitin-specific-processing protease 4, Ubiquitin thioesterase 4, Ubiquitous nuclear protein homolog, UNP, Unph, UNPH

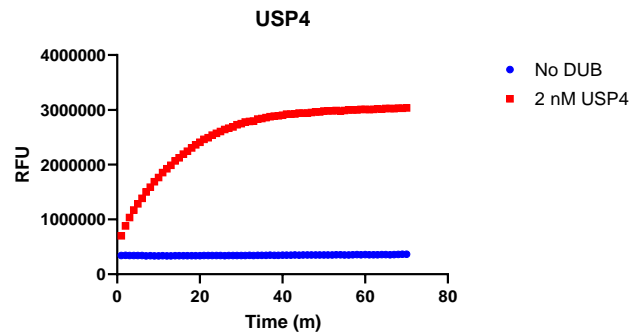
### Product Information

<b>Purity</b>	≥ 85% by SDS-PAGE
<b>Accession No.</b>	Q13107
<b>Molecular Weight</b>	120 kDa
<b>Quantity</b>	25 µg
<b>Physical State</b>	Liquid, 25 mM Tris-HCl, pH 7.5, 150 mM NaCl, 5 mM DTT, 10% glycerol
<b>Source</b>	Human Recombinant
<b>Tag</b>	None
<b>Suggested substrate</b>	Ub-CHOP reporter. Ubiquitin-rhodamine
<b>Storage</b>	-80° C. Avoid repeated freeze/thaw cycles

### Product QC



**SDS-Page Analysis of purified USP4.** Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



**Activity Assay of USP4.** 2 nM USP4 Protein was tested in a Ub-CHOP assay showing robust DUB activity.

### References

1. Chen, Y., et al., PLoS One., 2023. 18(8)e0290688.
2. Wang, Y., et al., Cancer Cell Int., 2020. 20:298.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2009 LifeSensors, Inc. All Rights Reserved