

USP15 (Ubiquitin Specific Protease 15)

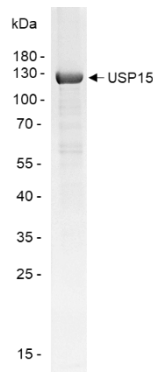
Cat. # DB513

Background	USP15 (Ubiquitin Specific Protease 15, UNPH4) is a deubiquitinase that is associated with the COP9 signalosome (CSN) complex. It regulates various pathways such as the TGF-beta receptor signaling, NFkB and RNF41/NRDP1-PRKN pathways.
Alternate names	Deubiquitinating enzyme 15, KIAA0529, MGC131982, MGC74854, Ubiquitin carboxyl-terminal hydrolase 15, Ubiquitin-specific-processing protease 15, Ubiquitin thioesterase 15, Unph-2, Unph4, UNPH4

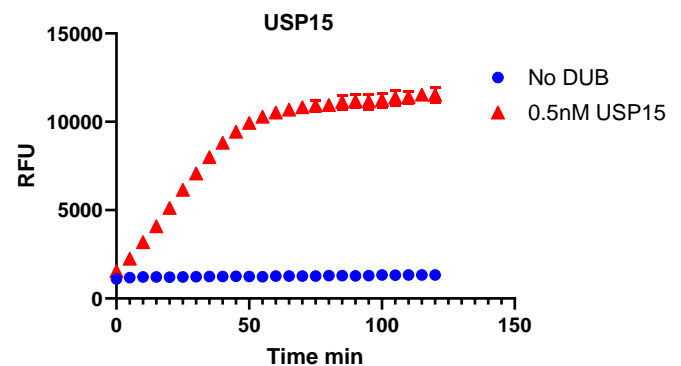
Product Information

Purity	≥ 90% by SDS-PAGE
Accession No.	Q08AL5
Molecular Weight	121 kDa
Quantity	25 µg
Physical State:	Liquid, 50 mM sodium phosphate pH 7.5, 50 mM NaCl, 1 mM DTT
Source:	Human recombinant
Tag:	His6-SUMO
Activity:	This enzyme is active with IQF-Diubiquitin DU4802 and in the Ub-CHOP assay.
Storage:	-80° C. Avoid repeated freeze/thaw cycles

Product QC



SDS-Page Analysis of purified USP15. Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of USP15. 0.5 nM USP15 was tested in Ub-Rh110 assay showing robust DUB activity.

References

1. Kim, MJ., et al., Cell Death Dis., 2022. 13(4):348.
2. Timani, K., et al., Front Oncol., 2022. 12:843157.

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