

USP5 (Ubiquitin specific protease 5)-(isoT)

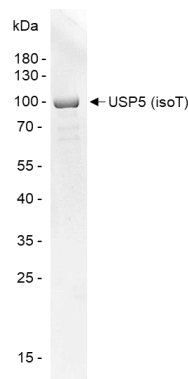
Cat. # DB516

Background	USP5 (Isopeptidase T; IsoT) is a deubiquitinase that is responsible for the deconjugation of the majority of unanchored polyubiquitin chains. USP5 is a Zinc-finger containing deubiquitinating enzyme that releases one ubiquitin at a time from the proximal end of the chain.
Alternate names	Deubiquitinating enzyme 5, Isopeptidase T, IsoT, ISOT, Ubiquitin carboxyl-terminal hydrolase 5, Ubiquitin-specific-processing protease 5, Ubiquitin thioesterase 5

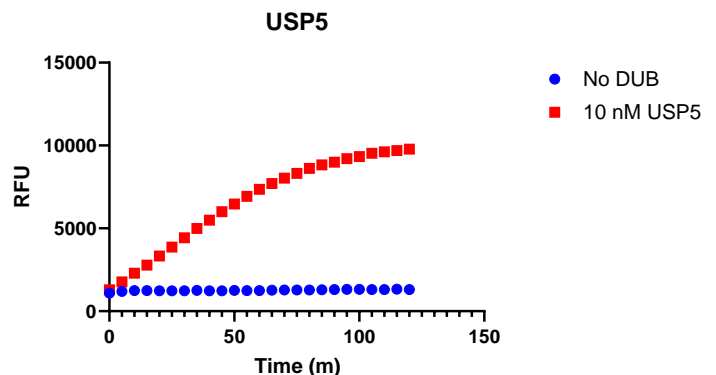
Product Information

Purity	≥ 90% by SDS-PAGE
Molecular Weight	94 kDa
Quantity	25 µg
Physical State	Liquid
Source	Human Recombinant
Tag	His6
Activity	This enzyme is active in the Ub-CHOP assay.
Storage	-80° C. Avoid repeated freeze/thaw cycles

Product QC



SDS-Page Analysis of purified USP5. Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of USP5. 10 nM USP5 was tested in a Ub-Rh110 assay showing robust DUB activity.

References

1. Cai, B., et al., Autophagy, 2022. 18(5):990-1004.
2. Zhang, Z., et al., Transl Lung Cancer Res., 2021. 10(10):3995-4011.

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