

USP2 Core (Ubiquitin specific protease 2)

Cat. # DB501

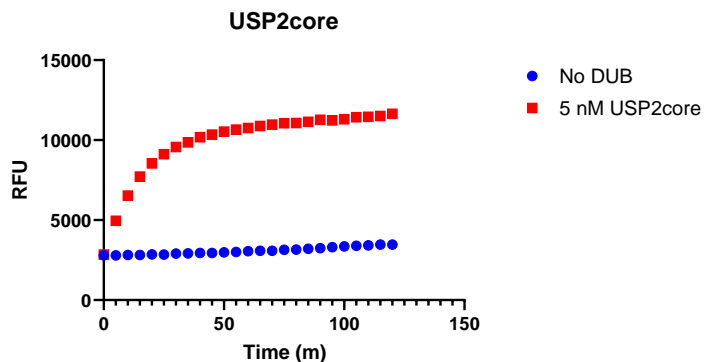
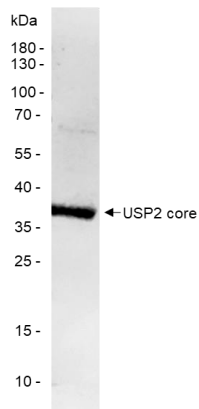
Background USP2 core represents the common catalytic core domain of the two isoforms of USP2 (Ubiquitin specific protease 2), USP2a and USP2b. It deubiquitinates polyubiquitinated target proteins such as MDM2, MDM4 and CCND1. USP2 plays a role in the G1/S cell-cycle progression in normal and cancer cells.

Alternate names 41 kDa ubiquitin-specific protease, Deubiquitinating enzyme 2, Ubiquitin carboxyl-terminal hydrolase 2, Ubiquitin-specific-processing protease 2, Ubiquitin thioesterase 2, UBP41, USP9

Product Information

Purity	≥ 90% by SDS-PAGE
Molecular Weight	40 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 USP2 core. Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.

Activity Assay of USP2core. 5 nM USP2core was tested in a CHOP assay showing robust DUB activity.

References

1. Nadolny, C., et al., Am J Cancer Res., 2021. 11(10):4746-4767.
2. Zhu, H., et al., Exp Cell Res., 2023. 427(1):113597.

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