

USP7 (Ubiquitin Specific Protease 7)

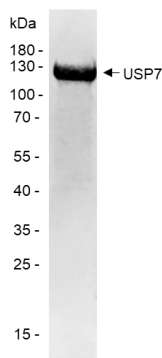
Cat. # DB502

Background	Ubiquitin-specific protease 7 (USP7), also known as herpesvirus-associated ubiquitin-specific protease (HAUSP) was originally identified by its binding activity to a herpes viral protein, ICPO. This DUB is known for its wide-ranging influence on various cellular processes, including the control of p53, MDM2, and other key regulatory proteins involved in cell cycle regulation and DNA damage response. USP7 plays a key role in regulating the ubiquitination and the stability of the RING-finger E3 ligase MDM2 (and its human homolog HDM2). Additional targets of USP7 have been identified including the forkhead transcription factor, FOXO44.
Alternate names	Deubiquitinating enzyme 7, HAUSP, Herpesvirus-associated ubiquitin-specific protease, TEF1, Ubiquitin carboxyl-terminal hydrolase 7, Ubiquitin-specific-processing protease 7, Ubiquitin thioesterase 7

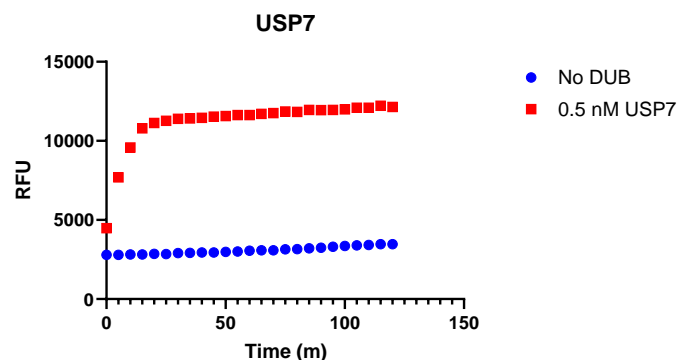
Product Information

Purity	≥ 95% by SDS-PAGE
Molecular Weight	135 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 USP7. Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of USP7. 0.5 nM USP7 was tested in a CHOP assay showing robust DUB activity.

References

1. Harakandi, C., et al., Bioorg Chem., 2021. 116:105273.
2. Qi, SM., et al., Front Cell Dev Biol., 2020. 8:233.

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