

USP8 (Ubiquitin Specific Protease 8)

Cat. # DB504

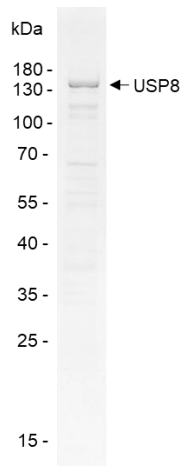
Background USP8 (Ubiquitin Specific Protease 8) is a deubiquitinase that converts both K48- and K63-linked ubiquitin chains. USP8 was shown to play a critical role in the stability of receptor tyrosine kinases.

Alternate names Deubiquitinating enzyme 8, hUBPy, HumORF8, KIAA0055, MGC129718, Ubiquitin carboxyl-terminal hydrolase 8, Ubiquitin-specific processing protease 8, Ubiquitin thioesterase 8, UBPY

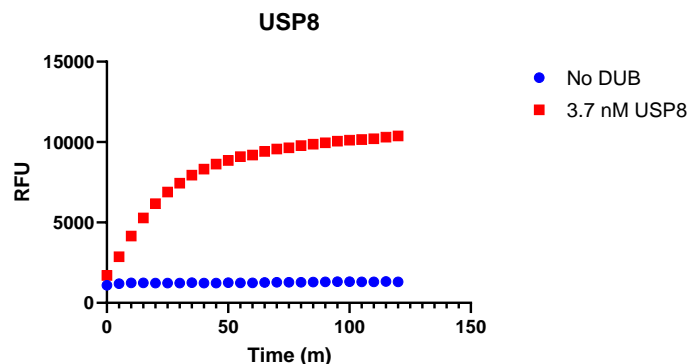
Product Information

Purity	≥ 80% by SDS-PAGE
Accession No	Q05DF5
Molecular Weight	138 kDa
Quantity	25 µg
Physical State	Liquid, 25 mM Tris-HCl, pH 7.4, 150 mM NaCl, 5 mM DTT, 10% glycerol
Source	Human Recombinant
Tag	His6-SUMO
Suggested Substrate	Ub-CHOP Reporter
Storage	-80° C. Avoid repeated freeze/thaw cycles

Product QC



SDS-Page Analysis of purified USP8. Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of USP8. 3.7 nM USP8 was tested in a CHOP assay showing robust DUB activity.

References

1. Xiong, W., et al., Nat Commun., 2022. 13(1):1700.
2. Zhu, Y., et al., Aging, 2021. 13(11):14999-15012.

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