

USP8core (Ubiquitin Specific Protease 8 core domain)

Cat. # DB503

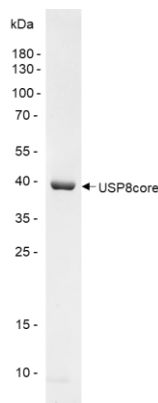
Background USP8 (Ubiquitin Specific Protease 8) is a deubiquitinase that converts both K48- and K63-linked ubiquitin chains. USP8 was shown to play a critical role in the stability of receptor tyrosine kinases. The USP8core contains only the core domain of the USP8 protein.

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

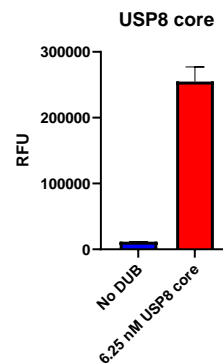
Product Information

Purity	≥ 95% by SDS-PAGE
Molecular Weight	46 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 USP8core. Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of USP8core. 6.25 nM USP8core was tested in a Ub-Rh110 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