

## USP33core (Ubiquitin specific protease 33 catalytic domain)

Cat. # DB508

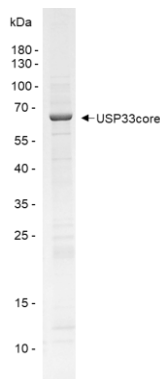
**Background** USP33 (VDU1) was originally identified as a protein that interacts with the E3 ubiquitin ligase pVHL. It is involved in various processes such as centrosome duplication, cellular migration and beta-2 adrenergic receptor/ADRB2 recycling. The USP33 catalytic core spans amino acids 181-712.

**Alternate names** Deubiquitinating enzyme 33, KIAA1097, MGC16868, Ubiquitin carboxyl-terminal hydrolase 33, Ubiquitin-specific-processing protease 33, Ubiquitin thioesterase 33, VDU1, VHL-interacting deubiquitinating enzyme 1

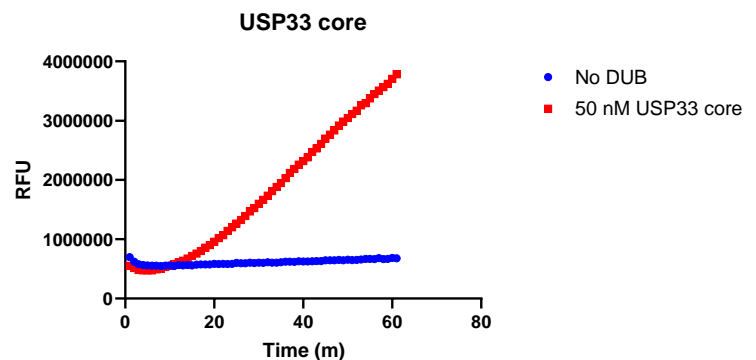
### Product Information

<b>Purity</b>	≥ 85% by SDS-PAGE
<b>Accession No</b>	Q8TEY7
<b>Molecular Weight</b>	61 kDa
<b>Quantity</b>	25 µg
<b>Physical State</b>	Liquid, 25 mM Tris-HCl, pH 7.5, 150 mM NaCl, 5 mM DTT, 10% glycerol
<b>Source</b>	Human Recombinant
<b>Tag</b>	His6-SUMO
<b>Activity</b>	This enzyme is active in the Ub-CHOP assay.
<b>Storage</b>	-80° C. Avoid repeated freeze/thaw cycles

### Product QC



**SDS-Page Analysis of purified USP33core.** Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



**Activity Assay of USP33core.** 50 nM USP33core was tested in a Ub-CHOP assay showing robust DUB activity.

### References

1. Niu, K., et al., Autophagy, 2020. 16(4):724-734/
2. Niu, Y., et al., Diabetes, 2022. 71(5):921-933.

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