

## USP20 (Ubiquitin Specific Protease 20)

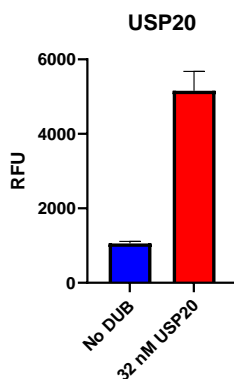
Cat. # DB520

<b>Background</b>	USP20 (Ubiquitin Specific Protease 20, VDU2) was identified by its homology to USP33 and binds competitively to the substrate recognition beta-domain of pVHL, a region commonly mutated in patients with pVHL disease. Furthermore, USP20 has been shown to be ubiquitinated by the pVHL E3 ligase complex <i>in vitro</i> and in cells. The enzyme acts as a regulator of G-protein coupled receptor (GPCR) signaling by mediating the deubiquitination beta-2 adrenergic receptor (ADRB2).
<b>Alternate names</b>	Deubiquitinating enzyme 20, KIAA1003, LSFR3A, Ubiquitin carboxyl-terminal hydrolase 20, Ubiquitin-specific-processing protease 20, Ubiquitin thioesterase 20, VDU2

### Product Information

<b>Molecular Weight</b>	105 kDa
<b>Quantity</b>	25 µg
<b>Physical State</b>	Liquid
<b>Source</b>	Human Recombinant
<b>Tag</b>	His6
<b>Activity</b>	This enzyme is active in the Ub-CHOP assay.
<b>Storage</b>	-80° C. Avoid repeated freeze/thaw cycles

### Product QC



**Activity Assay of USP20.** 32 nM USP20 was tested in a CHOP assay showing robust DUB activity.

### References

1. Li, Q., et al., Cell Death Dis., 2022. 13(5):434.
2. Li, W., et al., Genes Dev., 2020. 34(19-20):1310-1315.

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