

USP21 (Ubiquitin Specific Protease 21)

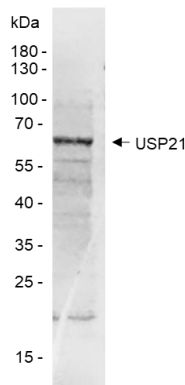
Cat. # DB509

Background	USP21 (Ubiquitin Specific Protease 21) regulates transcription initiation via deubiquitination of histone H2A and exhibits both deubiquitinase and deNEDDylase activities.
Alternate names	Deubiquitinating enzyme 21, MGC3394, NEDD8-specific protease, Ubiquitin carboxyl-terminal hydrolase 21, Ubiquitin-specific-processing protease 21, Ubiquitin thioesterase 21, USP16, USP23

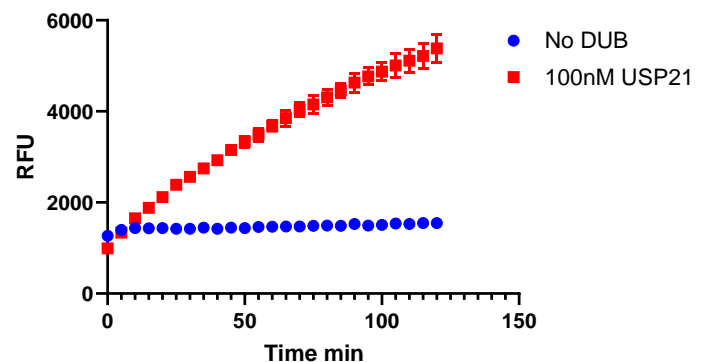
Product Information

Purity	≥ 85% by SDS-PAGE
Accession No.	Q9UK80
Molecular Weight	74.7 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	His ₆ -SUMO
Suggested Substrate	UB-CHOP Reporter, Ubiquitin-Rhodamine
Storage	-80°C. Avoid repeated freeze/thaw cycles

Product QC



SDS-Page Analysis of purified USP21. Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of USP21. 100 nM USP21 was tested in Ub-Rh110 assay showing robust DUB activity.

References

- Hou, P., et al., Genes Dev., 2019. 33(19029):1361-1366.
- Zhang, Q., et al., Cell Death Dis., 2022. 13(8):712.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2012 LifeSensors, Inc. All Rights Reserved