

USP28 (Ubiquitin Specific Protease 28)

Cat. # DB525

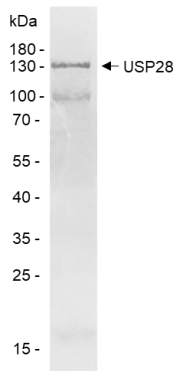
Background USP28 (Ubiquitin Specific Protease 28) is a deubiquitinase involved in DNA damage response checkpoint and MYC proto-oncogene stability.

Alternate names Deubiquitinating enzyme 28, KIAA1515, Ubiquitin carboxyl-terminal hydrolase 28, Ubiquitin-specific-processing protease 28, Ubiquitin thioesterase 28

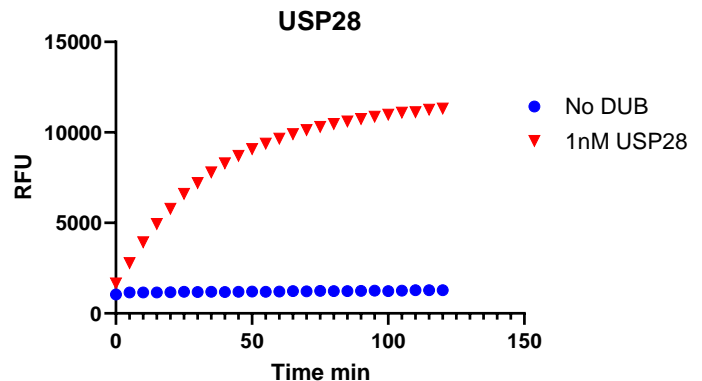
Product Information

Purity	≥ 85% by SDS-PAGE
Molecular Weight	123 kDa
Quantity	25 µg
Physical State	Liquid
Source	Human Recombinant
Tag	His6
Activity	This enzyme is active with ubiquitin rhodamine 110 (Cat. # SI230) and in the Ub-CHOP assay (Cat. # PR1001).
Storage	-80° C. Avoid repeated freeze/thaw cycles

Product QC



SDS-Page Analysis of purified USP28. Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of USP28. 1.0 nM USP28 was tested in a Ub-Rh110 assay showing robust DUB activity.

References

1. Chen, L., et al., Cell Death Dis., 2021. 12(10):887.
2. Zhao, LJ., et al., J Cell Biochem., 2019. 120(5):7657-7666.

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