

Den1 (Deneddylase 1)

Cat. # DB801

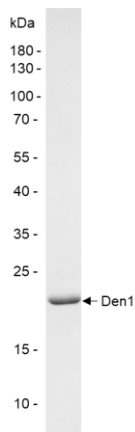
Background DEN1 (SENP8, NEDP1) is a deNEDDylase. It deconjugates the NEDD8 moiety from NEDDylated proteins. A recent study demonstrated that DEN1 deNEDDylated the tumor suppressor breast cancer-associated protein BCA3. Also, full length DEN1 was shown to cleave NEDD8-PLA₂ with high efficiency.

Alternate Names DEN1, Deneddylase-1, FKSG8, HsT17512, NEDD8-specific protease 1, NEDP1, Protease, cysteine 2, PRSC2, Sentrin/SUMO-specific protease SENP8, Sentrin-specific protease 8

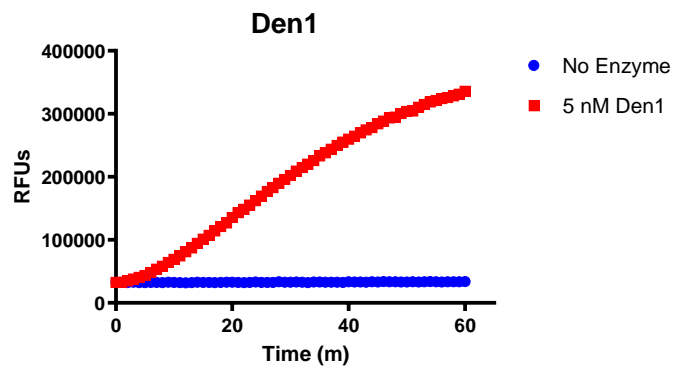
Product Information

Purity	≥ 95% by SDS-PAGE
Molecular Weight	51 kDa
Quantity	25 µg
Physical State	Liquid
Source	Human Recombinant
Tag	GST
Activity	This enzyme is active in the NEDD8-IsoPro Assay.
Storage	Store at -80°C, avoiding repeated freeze-thaw cycles.

Product QC



SDS-Page Analysis of purified Den1. Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of Den1. 5 nM Den1 was tested in a Nedd8-CHOP assay showing robust DUB activity.

References

- 1) Song, JM., et al., J Neurochem., 2023. 165(3):348-361.
- 2) Xie, P., et al., Cell Res., 2021. 31(3):291-311.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2007 LifeSensors, Inc. All Rights Reserved