

NEDD8 activating enzyme E1 (human), (recombinant) (His-tag)

Conjugation of mature NEDD8 to specific lysine residues on target proteins via isopeptide bonds allows NEDD8 to play a critical regulatory role in cell proliferation and development, a process closely associated to the ubiquitin-proteasome pathway. An increasing number of substrates for NEDDylation are being identified including the tumor suppressors p53 and VHL and members of the well-characterized cullin ubiquitin E3 ligase family, underlining the vital role NEDD8 modification plays in cellular processes.

Citations: 3

[View Online »](#)

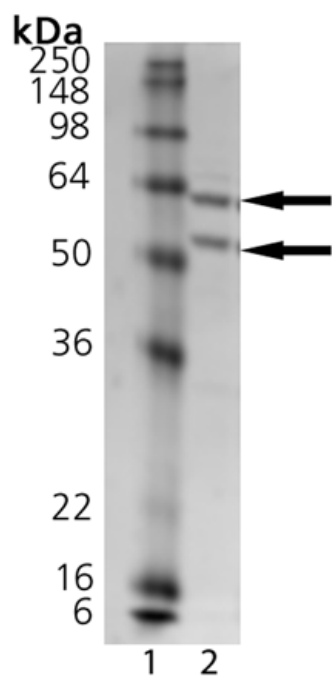
Ordering Information

[Order Online »](#)

ENZ-PRT112-0050	50µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis. Lane 1: MW Marker; Lane 2: 1 ug NEDD8 E1.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name

NEDD8-activating enzyme E1 catalytic subunit, NEDD8-activating enzyme E1C, Ubiquitin-activating enzyme E1C, Ubiquitin-like modifier-activating enzyme 3, Ubiquitin-activating enzyme 3

Application Notes

For use in NEDD8 conjugation reactions and NEDDylation pathway studies.

Formulation

Liquid. In 25mM HEPES, pH 7.5, containing 150mM NaCl, 1mM DTT.

MW

62kDa + 53kDa (heterodimer: 112.15kDa)

Purity

≥95% (SDS-PAGE)

Purity Detail

Copurified in His-tagged form with Uba3; copurified with APP-BP1.

Source

Produced in *E. coli*.

UniProt ID

Q8TBC4

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)