

# NEDP1 (human), (recombinant) (His- tag)

NEDP1, a human NEDD8-specific protease, is highly conserved throughout evolution and equivalent proteins are present in yeast, plants, insects, and mammals. NEDP1 appears to be specific for NEDD8 as neither ubiquitin nor SUMO bearing COOH- terminal extensions are utilized as substrates. Inhibition studies and mutagenesis indicate that NEDP1 is a cysteine protease with sequence similarities to SUMO-specific proteases and the class of viral proteases typified by the adenovirus protease. *in vivo* NEDP1 deconjugates NEDD8 from a wide variety of substrates including the cullin component of SCF-like complexes. Thus NEDP1 is likely to play an important role in ubiquitin-mediated proteolysis by controlling the activity of SCF complexes.

## Ordering Information

[Order Online »](#)

BML-UW9770-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -80°C

Shipping      Dry Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

**Alternative Name**      NEDD8-specific protease 1, SENP8, Sentrin-specific protease 8, Deneddylase 1

**Formulation**      Liquid. In 50mM TRIS-Cl, pH 7.7, containing 50% glycerol, 1mM DTT.

**Source**      Produced in *E. coli*.

**UniProt ID**      Q96LD8



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)