

NEDD8 (human), (recombinant) (biotin conjugate)

Biotinylated NEDD8 is a useful reagent for studying NEDDylation *in vitro*, where it represents a readily detectable alternative to radio-iodinated NEDD8. Biotinylated NEDD8 is activated by NEDD8 E1, is a competent substrate for the NEDD8 E2 conjugating enzyme UbcH12 and can be conjugated to target proteins in the presence of a suitable NEDD8 E3 ligase. The biotinyl moiety also provides a potential route to affinity purification of conjugated substrates (or thiol ester-linked enzymes) on avidin/streptavidin matrices. NEDD8 contains 9 lysine residues, any/many of which are likely to be biotinylated in this product. These modification sites have not been mapped. Biotinylated NEDD8 is prepared from recombinant human His-tagged NEDD8 (Prod. No. BML-UW9225).

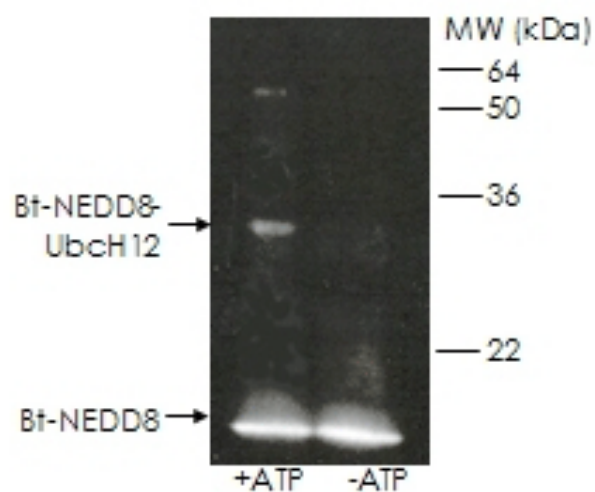
Ordering Information

[Order Online »](#)

BML-UW0560-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot of biotinylated-NEDD8 (Prod. No. BML-UW0560) thioester assay. Bt-NEDD8 conjugated to UbcH12 (Prod. No. BML-UW9145) in the presence of NEDD8 E1 (Prod. No. ENZ-PRT112). Bt-NEDD8 conjugated species detected with streptavidin-HRP.

Handling & Storage

Use/Stability	Stable for at least 12 months after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Formulation	Lyophilized. Contains PBS.
MW	~11kDa (based on Bt3-NEDD8 as predominant form present).
Reconstitution	Reconstitute with sterile distilled water.
Source	Produced in <i>E. coli</i> .
Technical Info / Product Notes	<p>Biotinylated mature protein for NEDD8 conjugation and NEDDylation pathway studies.</p> <p>Suggested uses:</p> <ol style="list-style-type: none">1. Ready detection of NEDD8 modification of specific proteins <i>in vitro</i>.2. Demonstration that novel proteins are potential targets for NEDDylation under <i>in vitro</i> conditions.3. Purification of NEDD8-conjugates via streptavidin affinity chromatography.
UniProt ID	Q15843



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)