

NEDD8-AMC

NEDD8-AMC is a fluorogenic substrate for deNEDDylating enzymes, including NEDP1 and UCH-L3. It is a particularly useful reagent for the study of deNEDDylating activity where detection sensitivity or continuous monitoring of activity is essential.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-UW0050-0025	25µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Liquid.
Application Notes	Uses: 1. Substrate for deNEDDylating enzyme activity assays. 2. Identification/confirmation of enzyme deNEDDylation activity. 3. Investigation of deconjugating enzyme substrate specificity in comparison with alternative UBL-AMC substrates (e.g. Ubiquitin-AMC, Prod. No. BML-SE211).
Concentration	1 mg/ml
Formulation	Solution in DMSO.
MW	8.735 kDa
Purity	≥95% (HPLC)
Source	Produced in <i>E. coli</i> .
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)