

Ubiquitin-AMC

High purity industry standard substrate for deubiquitinating enzymes (DUBs).

Ubiquitin-AMC is a fluorogenic substrate for a wide range of deubiquitinating enzymes (DUBs), including ubiquitin C-terminal hydrolases (UCHs) and ubiquitin specific proteases (USPs). It is a particularly useful reagent for the study of deubiquitinating activity where detection sensitivity or continuous monitoring of activity is essential. Ubiquitin-AMC has been shown to be a sensitive substrate for UCH-L3 (Prod. No. BML-UW9745) and for Isopeptidase-T (Prod. No. BML-UW9690), and is particularly useful for studying deubiquitinating activity where detection sensitivity or continuous monitoring of activity is essential.

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE211-0025	25µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)

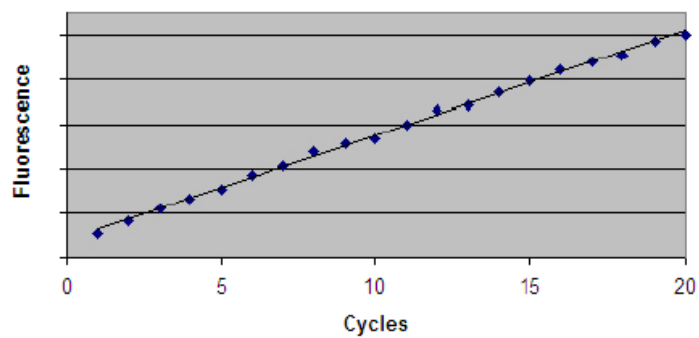


Figure: DUB Assay: 50nM UCLH1, His6-tagged (UW9740); 500nM Ubiquitin-AMC; 50mM HEPES pH7.8, 0.5mM EDTA, 1mM DTT.

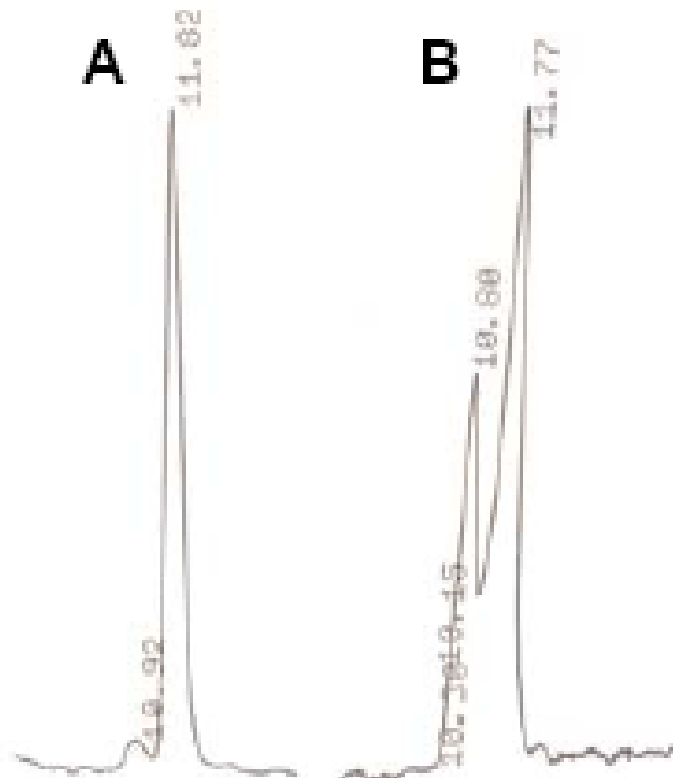


Figure: HPLC analysis. Comparison of Prod. No. BML-SE-211 (A) with contaminated alternatively sourced material (B).

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

Application Notes

Can be used for the following applications:

1. Substrate for deubiquitinating enzyme activity assays.
2. Identification/confirmation of enzyme deubiquitinylation activity.
3. Investigation of deconjugating enzyme substrate specificity in comparison with alternative UBL substrates.

Formulation

In DMSO

Purity

≥95% (HPLC)

Purity Detail

Please note: considerable efforts are taken to ensure that ubiquitin-AMC is homogeneous and not contaminated with ubiquitin and ubiquitin deletions, which are common in material obtained from other sources. The chromatogram below demonstrates the homogeneity of BML-SE211.

Source

Produced in *E.coli*.

UniProt ID

P0CG47 (UBB), P0CG48 (UBC), P62979 (RPS27A), P62987 (UBA52)

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)