

SUMO-1 (93-97)-AMCA

SUMO1-AMCA is a synthetic AMCA-peptide substrate based on the C-terminal residues of SUMO1 and is a useful fluorogenic substrate for deSUMOylating enzymes, including SENP1 (BML-UW9760). It has particular utility for the study of deSUMOylating activity where continuous monitoring of activity is essential. It provides an alternative substrate to SUMO1-AMC.

Citations: 1

[View Online »](#)

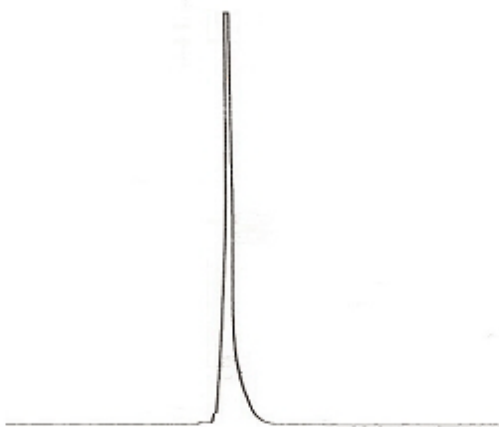
Ordering Information

[Order Online »](#)

BML-UW0500-1000	1mg
-----------------	-----

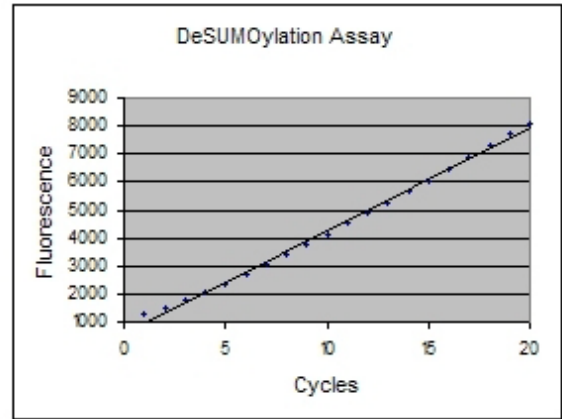
Manuals, SDS & CofA

[View Online »](#)



HPLC TRACE

HPLC trace



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solid at -20°C for up to twelve months. Solutions can be stored for up to three months at -20°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application Notes

Useful fluorogenic substrate for deSUMOylating enzymes.

Typical assay conditions:

Assay substrate concentration: 12.5µM-100µM.

Enzyme concentrations: 0.5µM-4µM.

Release of AMCA fluorescence by deSUMOylating enzymes can be monitored using fluorimetry (380nm excitation and 460nm emission wavelengths).

Uses:

1. Substrate for deSUMOylating enzyme activity assays.
2. Identification/confirmation of enzyme deSUMOylation activity.
3. Investigation of deconjugating enzyme substrate specificity in comparison with alternative UBL-AMC/AMCA substrates (e.g. SUMO1-AMC, NEDD8-AMCA).

Formulation

Lyophilized solid.

MW

747

Purity

≥95% (HPLC)

Sequence

E⁹³QTGG⁹⁷-AMCA

Solubility

Soluble in DMSO. Can be subsequently diluted in aqueous buffers.

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)