

# 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.

## Ordering Information

[Order Online »](#)

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

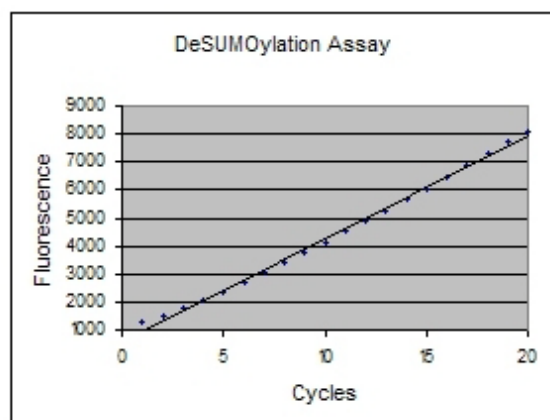
## Manuals, SDS & CofA

[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	<p>Useful fluorogenic substrate for deSUMOylating enzymes.</p> <p>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).</p> <p><b>Uses:</b></p> <ol style="list-style-type: none"><li>1. Substrate for deSUMOylating enzyme activity assays.</li><li>2. Identification/confirmation of enzyme deSUMOylation activity.</li><li>3. Investigation of deconjugating enzyme substrate specificity in comparison with alternative UBL-AMC/AMCA substrates (e.g. SUMO1-AMC, NEDD8-AMCA).</li></ol>
Formulation	Lyophilized solid.
MW	747
Purity	≥95% (HPLC)
Sequence	E <sup>93</sup> QTGG <sup>97</sup> -AMCA
Solubility	Soluble in DMSO. Can be subsequently diluted in aqueous buffers.



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-fr@enzolifesciences.com](mailto:info-fr@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)