

Proteasome substrate I (fluorogenic)

Fluorogenic proteasome substrate.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-088-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Z-LLL-AMC, Z-Leu-Leu-Leu-AMC

Appearance White lyophilized powder.

Formula $C_{36}H_{48}N_4O_7$

MW 648.8

Purity ≥98%

Sequence Z-Leu-Leu-Leu-AMC (AMC=7-Amino-4-methylcoumarin)

Solubility Soluble in DMSO.

Technical Info / Product Notes AMC has an excitation maximum of 340-360nm and an emission maximum of 440-460nm.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com