

Z-Leu-Leu-Glu-AMC

Fluorogenic substrate for the peptidylglutamyl-peptide hydrolysing (caspase-like) activity of the proteasome which may be stimulated in the presence of Mg^{2+} ions and rapidly inactivated by N-acetylimidazole.

Citations: 41

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-ZW9345-0005	5mg
BML-ZW9345-0050	50mg

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Z-LLE-AMC, Proteasome substrate II (fluorogenic)

Appearance Lyophilized. Off white powder.

Formula $C_{35}H_{44}N_4O_9$

MW 664.75 Da

Purity ≥95% (HPLC)

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

Solubility Soluble in dimethylformamide.

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)