

Ac-Gly-Pro-Leu-Asp-AMC

Highly specific substrate for the caspase-like peptidic activity of the 20S proteasome.

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-AW9560-0005	5mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at or below -20°C for up to 3 months
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Ac-GPLD-AMC
Appearance	White solid.
Formula	$C_{29}H_{37}N_5O_9$
MW	599.64
Purity	≥98% (HPLC)
Sequence	Ac-GPLD-AMC, Fluorogenic Substrate
Solubility	Soluble in 20% aqueous acetic acid (1mg/ml).
Source	Synthetic.

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)