

Z-Leu-Arg-Gly-Gly-AMC

Substrate for ubiquitin C-terminal hydrolases (UCHs)

Fluorogenic substrate for ubiquitin C-terminal hydrolases (UCHs), including UCH-L3 and Isopeptidase T. Sequence is based on the C-terminal residues of ubiquitin. Ex.: 380 nm Em.: 460 nm.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P801-0005	5mg
---------------	-----

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-LRGG-AMC, Ubiquitin C-terminal hydrolase substrate (fluorogenic), UCH substrate (fluorogenic)

Appearance Lyophilized solid.

MW 692.8

Purity ≥95%

Sequence Z-Leu-Arg-Gly-Gly-AMC (7-amino-4-methylcoumarin)

Solubility Soluble in DMSO.

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)