

Boc-Leu-Arg-Arg-AMC

A substrate for the Kex2 endopeptidase from *S. cerevisiae*. This peptide is also a substrate for kallikrein 5 (KLK5), and a candidate substrate for the investigation of proteasome subunit function.

Citations: 21

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-BW8515-0005	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 Boc-LRR-AMC

Appearance White powder.

Formula $C_{33}H_{52}N_{10}O_7 \cdot 2 HCl$

MW 773.8

Purity ≥97% (HPLC)

Sequence Boc-LRR-AMC, Fluorogenic Substrate

Solubility Soluble in water (50mg/ml).

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)