

Z-RR-AMC

Cathepsin substrate

Fluorogenic substrate specific for cathepsin B ($k_{\text{cat}}/K_m = 10^5 \text{ M}^{-1}\text{s}^{-1}$). The following proteases do *not* cleave, or inefficiently cleave, this substrate: papain, cathepsins C/DPP-I, F, H, K/O2, L, L2/V, O, S, X/Z. Ex.: 365 nm, Em.: 440 nm, although the following Ex/Em can also be used: 355,380/450,460. This substrate is useful for analysis of cathepsin B activity in cell extracts, inhibitor screening, and kinetic analysis. Also available: fluorogenic calibration standard AMC (BML-KI144).

Citations: 11

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P137-0010	10mg
BML-P137-0050	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-Arg-Arg-AMC
CAS	88937-61-5
MW	621.7
Purity	≥98%
Sequence	Z-Arg-Arg-AMC . HCl [Z=Cbz=Benzyloxycarbonyl; AMC=7-amino-4-methylcoumarin]

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)