

OMNICATHEPSIN[®]

fluorogenic substrate

Fluorogenic substrate cleaved by plasma kallikrein and kallikrein 8, but not plasmin; by papain; and by most cathepsins (cathepsins B, C/DPP-I, F, K/O2, L, L2/V, O, S, and X/Z). Ex.: 380 nm, Em.: 460 nm, although the following Ex/Em can also be used: 355,375/440,450. This substrate is useful for inhibitor screening and kinetic analysis. Also available: fluorogenic calibration standard AMC (BML-KI144).

Citations: 15

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P139-0010	10mg
BML-P139-0050	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-FR-AMC, Z-Phe-Arg-AMC

CAS 65147-22-0

Formula $C_{33}H_{36}N_6O_6 \cdot HCl$

MW 649.2

Purity ≥99%

Sequence Z-FR-AMC, Z-Phe-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-
fr@enzolifesciences.com](mailto:info-fr@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)