

Z-GPR-AMC

Cathepsin K substrate

Fluorogenic substrate specific for cathepsin K, among other cathepsin family members. This substrate is also cleaved by granzyme A, thrombin, tryptase, trypsin, and kallikreins. Ex.: 365 nm, Em.: 440 nm, although the following Ex/Em can also be used: 355,380/450,460. This substrate is useful for inhibitor screening and kinetic analysis. Also available: fluorogenic calibration standard AMC (BML-KI144).

Citations: 11

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P142-0020	20mg
---------------	------

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-Gly-Pro-Arg-AMC

Appearance White powder.

Formulation $C_{31}H_{37}N_7O_7$

MW 656.1

Purity ≥97% (HPLC)

Sequence Z-Gly-Pro-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)