

H-Arg-AMC . 2HCl

Substrate for cathepsin H and aminopeptidases

Fluorogenic substrate for Cathepsin H ($K_m = 150 \mu\text{M}$) and aminopeptidases. 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: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P135-0025	25mg
---------------	------

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

MW	404.3
Purity	≥99%
Sequence	H-Arg-AMC . 2HCl [AMC=7-amino-4-methylcoumarin]

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com