

Met-AMC

Aminopeptidase substrate

Fluorogenic substrate for methionine aminopeptidase 2 (MetAP2) ($k_{\text{cat}}/K_{\text{m}} = 1.1 \times 10^5 \text{ M}^{-1}\text{min}^{-1}$; $K_{\text{m}} = 310 \text{ }\mu\text{M}$) or aminopeptidase N ($K_{\text{m}} = 377 \text{ }\mu\text{M}$). 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).

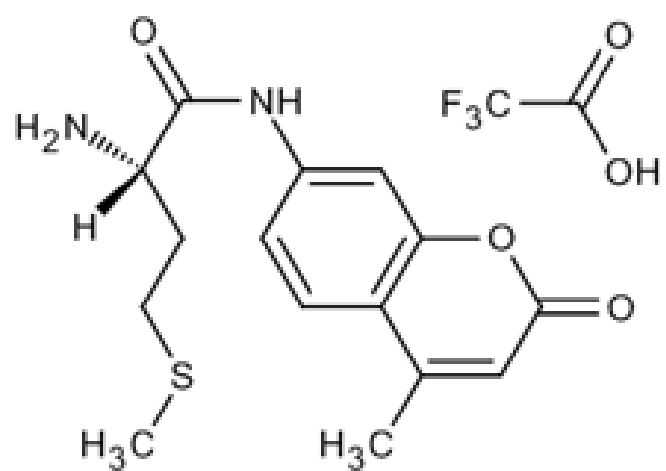
Ordering Information

[Order Online »](#)

BML-P236-0025	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

Appearance White to off white powder

CAS 94367-35-8

Formula $C_{17}H_{19}F_3N_2O_5S$

MW 420.4

Purity ≥98% (TLC)

Sequence H-Met-AMC . TFA [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)