

Z-VVR-AMC

Substrate for cathepsin S and cathepsin L

Sensitive fluorogenic substrate for cathepsin S and cathepsin L. Also cleaved by cathepsin B, cathepsin V, kallikrein 8, and factor VIIa/TF but not cathepsin K nor kallikrein 2. 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: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P199-0010	10mg
---------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Cathepsin S & L substrate (fluorogenic)
Appearance	Lyophilized solid.
CAS	124485-41-2
MW	663.8
Purity	≥99% (TLC)
Sequence	Z-Val-Val-Arg-AMC . HCl [Z=Cbz=Benzyloxycarbonyl; AMC=7-amino-4-methylcoumarin]
Solubility	Soluble in DMSO.

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)