

MeOSuc-AAPV-AMC

Elastase substrate

Highly sensitive substrate for neutrophil elastase (leukocyte elastase) ($K_m = 130 \mu M$), but not granzyme H. This substrate can be used for *in vitro* assays with purified enzyme, on live cells, or in conditioned medium. Ex: 380 nm, Em: 460 nm.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P224-0005	5mg
---------------	-----

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	Neutrophil elastase substrate (fluorogenic)
Appearance	White solid
CAS	72252-90-5
Formula	$C_{31}H_{41}N_5O_9$
MW	627.7
Purity	≥95% (HPLC)
Sequence	N-Methoxysuccinyl-Ala-Ala-Pro-Val-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)