

MeOSuc-AAPV-pNA

Substrate for neutrophil elastase and neutrophil proteinase 3

Highly sensitive substrate for human and mouse neutrophil elastase (leukocyte elastase) and neutrophil proteinase 3 (PR-3, myeloblastin), but not cathepsin G nor chymotrypsin. This substrate can be used for *in vitro* assays with purified enzyme, or with biological fluids or conditioned medium. Hydrolysis of the substrate can be detected at 405-410 nm.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P213-0005	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

CAS 70967-90-7

MW 590.6

Purity ≥98%

Sequence N-Methoxysuccinyl-Ala-Ala-Pro-Val-*p*-nitroanilide

Solubility Soluble in DMSO (>20 mM) or methanol (1 mg/mL).

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)