

Boc-Ala-Ala-Nva-SBzl

Proteinase substrate

Much more sensitive substrate for leukocyte proteinase 3 (PR-3, myeloblastin) ($K_{cat}/k_m = 1.06 \times 10^6 \text{ M}^{-1}\text{s}^{-1}$, $K_m = 63 \text{ }\mu\text{M}$) than MeOSuc-AAPV-pNA (Chromogenic Substrate) (Prod. No. BML-P213) ($K_{cat}/k_m = 1.06 \times 10^3 \text{ M}^{-1}\text{s}^{-1}$, $K_m = 470 \text{ }\mu\text{M}$). Also cleaved by neutrophil elastase ($K_{cat}/k_m = 9.2 \times 10^6 \text{ M}^{-1}\text{s}^{-1}$, $K_m = 2.2 \text{ }\mu\text{M}$), granzyme H, cathepsin G, chymotrypsin, and rat MCPI and II. Typical assay conditions: substrate concentration 88-250 μM ; enzyme concentration 0.2-20 nM; buffer [100 mM HEPES pH 7.5, 500 mM NaCl, 10% DMSO]; with 170 μM either 5,5'-dithiobis(2-nitrobenzoic acid) (DTNB) or 4,4'-dithiobis(pyridine) (from 15 mM DMSO stocks); 25°C. Hydrolysis of substrate can be detected at 412 nm for DTNB (

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P303-0005	5mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions in DMSO are stable for up to 3 months when stored at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white powder.
Formula	$C_{23}H_{35}N_3O_5S$
Formulation	Lyophilized.
MW	465.6
Purity	≥95% (HPLC)
Sequence	<i>t</i> -butyloxycarbonyl-Ala-Ala-Nva-thiobenzyl ester
Solubility	Soluble in DMSO (1mg/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-
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)