

Z-RR-pNA

Cathepsin substrate

Chromogenic substrate specific for cathepsin B. Release of pNA in monitored at 405-410 nm. This substrate is useful for inhibitor screening and kinetic analysis. Also available: calibration standard pNA (BML-KI145).

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P138-0025	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

Alternative Name Z-Arg-Arg-pNA

CAS 201807-90-1

Formula $C_{26}H_{36}N_{10}O_6 \cdot 2HCl$

MW 657.6

Purity ≥96%

Sequence Z-Arg-Arg-pNA . 2HCl [Z=Cbz=Benzyloxycarbonyl; pNA=p-nitroanilide]

Last modified: December 1, 2025



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)