

Leu-pNA

Aminopeptidase substrate

Substrate for aminopeptidases such as aminopeptidase N (APN) cystinyl aminopeptidase, and leucine aminopeptidase (LAP). This substrate can be used for *in vitro* assays with purified enzyme, or with cell extracts. Also available: calibration standard pNA (BML-KI145).

Ordering Information

[Order Online »](#)

BML-P230-0100	100mg
---------------	-------

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	Aminopeptidase substrate (chromogenic)
Appearance	Light yellow powder.
CAS	4178-93-2
Formula	$C_{12}H_{17}N_3O_3$
MW	251.3
Purity	≥98% (TLC)
Sequence	H-L-Leu-pNA [pNA=p-nitroanilide]

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)