

Ac-IEPD-pNA

Caspase substrate

Colorimetric granzyme B substrate. Combinatorial and kinetic studies have found the IEPD sequence to be optimal for cleavage of a tetrapeptide substrate by recombinant granzyme B, with a K_m for Ac-IEPD-pNA of 57 μ M. Cleavage of the substrate by granzyme B releases pNA, which can be detected by its absorbance at 405 nm.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P133-0005	5mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solution in DMSO (e.g. 0.1M) is highly stable at -20°C or -70°C.
Handling	Keep cool and dry.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Caspase-8 substrate (chromogenic), Granzyme B substrate (chromogenic)
Appearance	Lyophilized solid.
Formula	$C_{28}H_{38}N_6O_{11}$
MW	634.7
Purity	≥95%
Sequence	N-acetyl-Ile-Glu-Pro-Asp-pNA (<i>p</i> -nitroanilide)
Solubility	Soluble to at least 100mM in DMSO.

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)