

# 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  $\mu$ 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.



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)