

# Z-DEVD-pNA

## Caspase substrate

Chromogenic substrate for caspase-3 with increased cell permeability. Also substrate for caspase-1, -4 and -7.  $\lambda_{\text{max}}$  of pNA is 400nm. Similar to Ac-DEVD-pNA except the N-terminus is blocked by a benzyloxycarbonyl (Z) moiety which may increase cell permeability.

Citations: 5

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-260-048-M001	1mg
------------------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Caspase-3 substrate (chromogenic)
Appearance	White powder.
Formula	$C_{32}H_{38}N_6O_{14}$
Formulation	Lyophilized.
MW	730.7
Peptide Content	70-90%
Purity	≥99% (HPLC)
Sequence	Z-Asp-Glu-Val-Asp-pNa (pNA = p-Nitroaniline)
Solubility	Soluble in DMSO or acetonitrile.



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)