

Ac-YVAD-pNA

Caspase-1 substrate

Sequence is a modification of pro-interleukin-1 β Asp-116 cleavage site for caspase-1. Similar to Ac-YVAD-AMC but cleavage is monitored colorimetrically at 405nm.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-026-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Caspase-1 substrate (chromogenic)
Appearance	White to off-white powder.
CAS	149231-66-3
Formula	$C_{29}H_{36}N_6O_{10}$
Formulation	Lyophilized.
MW	628.6
Peptide Content	75-95%
Purity	≥97% (HPLC)
Sequence	Ac-Tyr-Val-Ala-Asp-pNA (pNA=p-Nitroaniline)
Solubility	Soluble in DMSO. Dilute with buffer pH 7.5.
Technical Info / Product Notes	λ_{max} of pNA is 400nm.

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-
fr@enzolifesciences.com](mailto:info-fr@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)