

Ac-VDVAD-pNA

Caspase-2 substrate

Chromogenic substrate for caspase-2. Preferred substrate for caspase-2 ($K_m=53\mu\text{M}$)

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-059-M001	1mg
ALX-260-059-M005	5mg

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-2 substrate (chromogenic)

Appearance White to off-white powder.

CAS 189684-53-5

Formula $C_{29}H_{41}N_7O_{12}$

Formulation Lyophilized.

MW 679.7

Peptide Content 70-90%

Purity ≥98% (HPLC)

Sequence Ac-Val-Asp-Val-Ala-Asp-pNA (pNA=p-Nitroanilin)

Solubility Soluble in DMSO (100mM). Dilute stock solution with distilled water or buffer.

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)