

Ac-YVAD-CHO

Inhibitor of caspase-1 and -4

Potent and reversible inhibitor of caspase-1. Also inhibits caspase-4. Strongly inhibits anti-Fas induced apoptosis in L929-Fas cells.

Citations: 18

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-027-M001	1mg
ALX-260-027-M005	5mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions should be divided into aliquots and stored at -20°C.
Handling	Keep under inert gas. Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Caspase-1 inhibitor (aldehyde)
Appearance	White powder.
Formula	$C_{23}H_{32}N_4O_8$
Formulation	Lyophilized.
MW	492.5
Peptide Content	70-90%
Purity	≥96% (HPLC)
Reconstitution	Add 0.2ml DMSO per 1.0mg peptide. Shake well until all the contents are dissolved. This furnishes a 10mM solution of this material.
Sequence	Ac-Tyr-Val-Ala-Asp-CHO
Solubility	Soluble in DMSO or water (both 50mg/ml).

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)