

Ac-VEID-CHO

Inhibitor of caspase-6, -8 and -10

Reversible aldehyde inhibitor of caspase-6. Also inhibits caspase-8 and -10. Sequence is based on lamin A site of cleavage.

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-062-M001	1mg
ALX-260-062-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. The solution is stable for weeks when stored at -20°C and protected from light. Do not dissolve in aqueous buffer due to possible racemization.
Handling	Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Caspase-6 inhibitor (aldehyde)
Appearance	White powder.
Formula	$C_{22}H_{36}N_4O_9$
Formulation	Lyophilized.
MW	500.6
Peptide Content	70-90%.
Purity	≥95% (HPLC)
Sequence	Ac-Val-Glu-Ile-Asp-CHO
Solubility	Soluble in DMSO or water.

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)