

# 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

<b>Use/Stability</b>	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.
<b>Handling</b>	Keep cool and dry.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

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



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)