

Ac-

# AAVALLPAVLLALLAPYVAD-CHO

## Caspase-1 inhibitor

The YVAD-CHO portion of this peptide is a highly-specific, potent ( $K_i=1$  nM) and reversible inhibitor of caspase-1 (interleukin-1 $\beta$  converting enzyme, ICE), whose sequence is based on the pre-interleukin-1 $\beta$  Asp-116 cleavage site for caspase-1. The sequence linked N-terminal to YVAD-CHO (residues 1-16) has been shown to enhance cell permeability. The efficacy of the resulting chimeric peptide, for caspase-1 inhibition, is unknown.

### Ordering Information

[Order Online »](#)

ALX-260-047-M001	1mg
------------------	-----

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Preincubate cultured cells 1-2 hrs. with 0.1-2.0µM of the peptide, before inducing for apoptosis. Measure indicator(s) of apoptosis at a number of time points, since the effect of the inhibitor may be a delay in the onset of apoptosis, rather than permanent inhibition. It may also be helpful to determine a minimum effective concentration for the apoptosis inducer and use that concentration, when attempting to observe the effects of this peptide.
<b>Handling</b>	Keep cool and dry. After reconstitution, prepare aliquots and store at -20°C.
<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-1 inhibitor (aldehyde)
<b>Appearance</b>	White to off-white powder.
<b>Formula</b>	$C_{97}H_{160}N_{20}O_{24}$
<b>MW</b>	1990.5
<b>Peptide Content</b>	78-98%
<b>Purity</b>	≥97% (HPLC)
<b>Sequence</b>	Ac-Ala-Ala-Val-Ala-Leu-Leu-Pro-Ala-Val-Leu-Leu-Ala-Leu-Leu-Ala-Pro-Tyr-Val-Ala-Asp-CHO
<b>Solubility</b>	Soluble in DMSO (0.25mg/ml).



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