

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 β converting enzyme, ICE), whose sequence is based on the pre-interleukin-1 β 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.

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	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.
Handling	Keep cool and dry. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Caspase-1 inhibitor (aldehyde)
Appearance	White to off-white powder.
Formula	$C_{97}H_{160}N_{20}O_{24}$
MW	1990.5
Peptide Content	78-98%
Purity	≥97% (HPLC)
Sequence	Ac-Ala-Ala-Val-Ala-Leu-Leu-Pro-Ala-Val-Leu-Leu-Ala-Leu-Leu-Ala-Pro-Tyr-Val-Ala-Asp-CHO
Solubility	Soluble in DMSO (0.25mg/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-
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)