

Ac-DEVD-CHO

Caspase-3 inhibitor

Highly specific, potent and reversible inhibitor of caspase-3. $K_i < 1\text{nM}$ for purified caspase-3. $\text{IC}_{50} = 0.2\text{nM}$ for PARP cleavage in cultured human osteosarcoma cell extracts. In contrast, the caspase-1 (ICE) specific inhibitor YVAD-CHO has a $K_i = 12\mu\text{M}$ for caspase-3 and $\text{IC}_{50} > 10\mu\text{M}$ for PARP cleavage in osteosarcoma cell extracts. The DEVD sequence is based on PARP cleavage site Asp-216 for caspase-3.

Citations: 25

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-030-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. Solutions are stable for at least 1 month when stored at -20°C in the dark and protected from moisture.
Handling	Keep under nitrogen. Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Caspase-3 inhibitor (aldehyde), Caspase-7 inhibitor (aldehyde)
CAS	169332-60-9
Formula	$C_{20}H_{30}N_4O_{11}$
Formulation	Lyophilized.
MW	502.5
Peptide Content	70-90%.
Purity	≥95% (HPLC)
Reconstitution	Add 0.2ml DMSO per 1mg peptide. Shake well until all the contents are dissolved. This furnishes a 10mM solution of this material.
Sequence	Ac-Asp-Glu-Val-Asp-CHO
Solubility	Soluble in distilled water or DMSO.



Last modified: May 29, 2024

ENZO LIFE SCIENCES
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-uk@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com