

# Biotinyl-DEVD-CHO

Inhibitor of caspase-3 and -7

Reversible inhibitor of caspase-3 and -7. Useful for the purification or extract depletion of caspase-3 in conjunction with streptavidin conjugates.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-034-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Packaged under inert gas.
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), CPP32 affinity reagent
Appearance	White powder.
Formula	$C_{28}H_{42}N_6O_{12}S$
Formulation	Lyophilized.
MW	686.7
Peptide Content	65-90%.
Purity	≥99% (HPLC)
Sequence	Biotinyl-Asp-Glu-Val-Asp-CHO
Solubility	Soluble in 50% acetic acid with agitation (1mg/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)