

Biotinyl-VAD(OMe)- FMK

Caspase inhibitor

Irreversible, cell permeable broad-spectrum caspase inhibitor.

Citations: 1

[View Online »](#)

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. The stock solution is stable at -20°C for 6-8 months.
Handling	Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Caspase family inhibitor (fluoromethylketone)
Appearance	Off-white to brown powder.
Formula	$C_{30}H_{49}FN_6O_8S$
MW	672.8
Purity	≥90% (TLC)
Sequence	Biotinyl-Ahx-Val-Ala-Asp(OMe)-fluoromethylketone
Solubility	Soluble in DMSO.

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)