

# Z-VAD(OMe)-FMK

## Caspase inhibitor

Irreversible and cell permeable broad-spectrum caspase inhibitor. Inhibits Fas-mediated apoptosis in Jurkat T cells and apoptosis in THP.1 cells induced by diverse stimuli (including cycloheximide, thapsigargin, etoposide and staurosporine). For cell free assays use Prod. No. ALX-260-020.

Citations: 24

[View Online »](#)

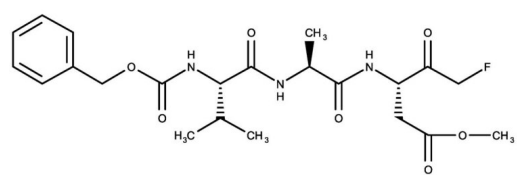
## Ordering Information

[Order Online »](#)

BML-P416-0001	1mg
BML-P416-0005	5mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for up to 1 year after receipt when stored at -20°C. Stock solutions should be divided into aliquots and stored at -20°C for up to 1 month.
Handling	Keep cool and dry.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Caspase family inhibitor (fluoromethylketone)
Appearance	White to off-white solid.
Application Notes	Irreversible, cell permeable, broad spectrum caspase inhibitor. Limited Fas-induced mouse liver damage and apoptosis to <1%, at a dose of 50µmol/kg. At 50µM, abrogated 7-hydroxystaurosporine and etoposide-induced apoptosis in HL60 cells.
CAS	187389-52-2
Formula	$C_{22}H_{30}N_3O_7F$
MW	467.5
Purity	≥95%
Sequence	Z-Val-Ala-Asp(OMe)-fluoromethylketone
Solubility	Soluble in DMSO (>10mg/ml), dimethyl formamide (7mg/ml) or 100% ethanol (1.3mg/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-fr@enzolifesciences.com](mailto:info-fr@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)