

# Z-

# D(OMe)E(OMe)VD(OMe)- FMK

## Inhibitor of caspase-3 and -7

Irreversible, cell permeable inhibitor of caspase-3 and -7.

Citations: 11

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-260-072-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stable for 1 year after receipt when stored at -20°C.
<b>Handling</b>	Packaged under inert gas. After reconstitution, prepare aliquots and store at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Caspase-3 inhibitor (fluoromethylketone), Caspase-7 inhibitor (fluoromethylketone)
<b>Formula</b>	$C_{30}H_{41}FN_4O_{12}$
<b>Formulation</b>	Lyophilized.
<b>MW</b>	668.7
<b>Purity</b>	≥95% (TLC)
<b>Sequence</b>	Z-Asp(OMe)-Glu(OMe)-Val-Asp(OMe)-fluoromethylketone
<b>Solubility</b>	Soluble in DMSO (20mM) or dimethyl formamide.

### Technical Info / Product Notes

We recommend a final concentration of 2-10µM for inhibiting caspase activity in cell culture. The optimal doses may vary for different cells and culture conditions.

**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)