

# Z-VAD-FMK

**Pan-Caspase Inhibitor.** Z-VAD-FMK acts as an effective irreversible caspase inhibitor with no cytotoxic effects and, therefore, is a useful tool for studying caspase activity.

Irreversible, cell permeable broad-spectrum caspase inhibitor.

Citations: 162

[View Online »](#)

## Ordering Information

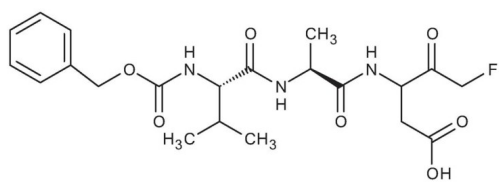
[Order Online »](#)

ALX-260-020-M001	1mg
ALX-260-020-M005	5mg

Manuals, SDS & CofA

[View Online »](#)

- Irreversible inhibitor
- Cell-permeable
- High purity peptide
- Non-cytotoxic



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Keep sealed after removing from the freezer until its temperature equilibrates with room temperature. Dissolve only the amount needed at one time because of the possibility of decomposition in solution.
<b>Handling</b>	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 family inhibitor (fluoromethylketone), Z-Val-Ala-DL-Asp-FMK
<b>Appearance</b>	Solid
<b>Application Notes</b>	Can be used as a broad spectrum caspase inhibitor in conjunction with apoptosis detection assays.
<b>CAS</b>	220644-02-0
<b>Formula</b>	$C_{21}H_{28}FN_3O_7$
<b>MW</b>	453.5
<b>Peptide Content</b>	77-97%
<b>Purity</b>	≥95% (HPLC)
<b>Sequence</b>	Z-Val-Ala-Asp-fluoromethylketone
<b>Solubility</b>	Soluble in DMSO or acetonitrile (10mg/ml).



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)