

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: 166

[View Online »](#)

Ordering Information

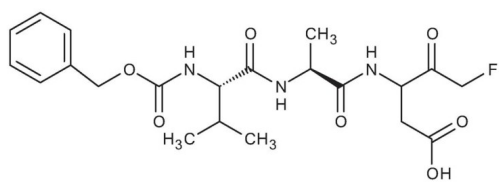
[Order Online »](#)

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

Use/Stability	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.
Handling	After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Caspase family inhibitor (fluoromethylketone), Z-Val-Ala-DL-Asp-FMK
Appearance	Solid
Application Notes	Can be used as a broad spectrum caspase inhibitor in conjunction with apoptosis detection assays.
CAS	220644-02-0
Formula	$C_{21}H_{28}FN_3O_7$
MW	453.5
Peptide Content	77-97%
Purity	≥95% (HPLC)
Sequence	Z-Val-Ala-Asp-fluoromethylketone
Solubility	Soluble in DMSO or acetonitrile (10mg/ml).