

Z-LE(OMe)HD(OMe)-FMK

Caspase-9 inhibitor

Irreversible, cell permeable inhibitor of caspase-9. The carboxyl groups are OMe ester form to enhance cell permeability. If purified enzyme is used, esterase should be added for the hydrolysis of the ester groups.

Ordering Information

[Order Online »](#)

ALX-260-076-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions should be stored in aliquots at -20°C.
Handling	Very hygroscopic.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Caspase-9 inhibitor (fluoromethylketone)
Appearance	Off-white to tan solid.
Formula	$C_{32}H_{43}FN_6O_{10}$
Formulation	TFA salt.
MW	690.7
Purity	≥95% (HPLC)
Sequence	Z-Leu-Glu(OMe)-His-Asp(OMe)-fluoromethylketone
Solubility	Soluble in DMSO (20mM).

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)