

Z-WEHD-FMK (Ready-to-Use)

Inhibitor of caspase-1, -4, -5 and -14

Irreversible, cell permeable inhibitor of caspase-1 and -5. Inhibits also caspase-4 and-14. Designed as a methyl ester to facilitate cell permeability.

Ordering Information

[Order Online »](#)

ALX-260-140-R020	10mM - 20µl
ALX-260-140-R100	2mM - 100µl

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Caspase-1 inhibitor (fluoromethylketone), Caspase-5 inhibitor (fluoromethylketone)
Formula	$C_{37}H_{42}FN_7O_{11}$
Formulation	Liquid. In DMSO.
MW	763.8
Purity	≥99% (HPLC)
Sequence	Z-Trp-Glu(OMe)-His-Asp(OMe)-fluoromethylketone

Technical Info / Product Notes

We recommend a final concentration of 2µM for inhibiting caspase activity in Jurkat 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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com