

Z-YVAD(OMe)-FMK

Inhibitor of caspase-1 and -4

Irreversible, cell permeable inhibitor of caspase-1. Inhibits also caspase-4.

Citations: 7

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions should be divided into aliquots and stored at -20°C.
Handling	Packaged under inert gas.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Caspase-1 inhibitor (fluoromethylketone)
Appearance	White to off-white powder.
Formula	$C_{31}H_{39}FN_4O_9$
MW	630.7
Purity	≥98% (TLC)
Sequence	Z-Tyr-Val-Ala-Asp(OMe)-fluoromethylketone
Solubility	Soluble in DMSO.

Technical Info / Product Notes

The carboxyl groups are OMe ester forms to enhance cell permeability. If purified enzyme is used, esterase should be added for the hydrolysis of the ester groups.

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