

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

Caspase-6 inhibitor

Irreversible, cell permeable inhibitor of caspase-6. The inhibitor is designed as a methyl ester to facilitate cell permeability.

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-6 inhibitor (fluoromethylketone)

Formulation Liquid. In DMSO.

MW 652.0

Purity ≥98% (HPLC)

Sequence Z-Val-Glu-Ile-Asp-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-be@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