

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

Inhibitor of caspase-3 and -7

Irreversible, cell permeable inhibitor of caspase-3 and -7.

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-1410	2mM - 100µl
--------------	-------------

Manuals, SDS & CofA

[View Online »](#)

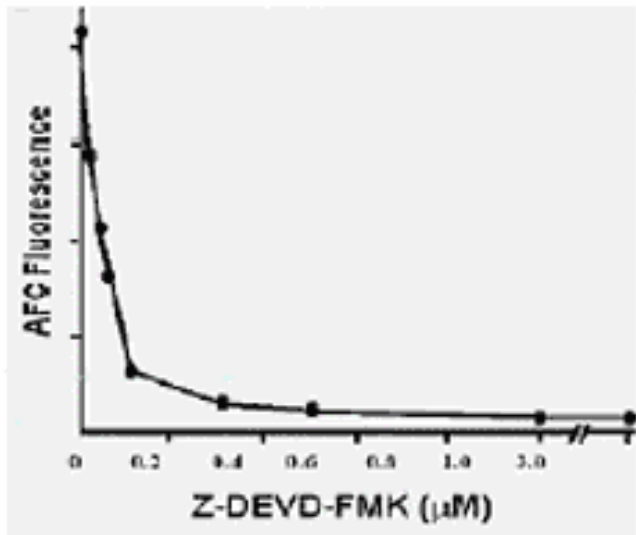


Figure: Apoptosis was induced in Jurkat cells with camptothecin in the presence of various concentrations (0-3µM) of Z-DEVD-FMK (Ready-to-Use) (Prod. No. [ALX-260-141](#)) as indicated. Caspase activity was analyzed using Caspase-3 Fluorometric Assay Kit (Prod. No. ALX-850-216). """"""

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-3 inhibitor (fluoromethylketone), Caspase-7 inhibitor (fluoromethylketone)
Formulation	Liquid. In DMSO.
MW	668.0
Purity	≥98% (HPLC)
Sequence	Z-Asp-Glu-Val-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