

# Z-IE(OMe)TD(OMe)-FMK

## Caspase and granzyme inhibitor

Irreversible, cell permeable inhibitor of caspase-8 and granzyme B. Inhibits also caspase-10.

Citations: 6

[View Online »](#)

## Ordering Information

[Order Online »](#)

|                  |     |
|------------------|-----|
| ALX-260-073-M001 | 1mg |
|------------------|-----|

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

|                   |                                                                                                              |
|-------------------|--------------------------------------------------------------------------------------------------------------|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. Store stock solutions in aliquots 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-8 inhibitor (fluoromethylketone), Granzyme B inhibitor (fluoromethylketone) |
| Appearance       | White to off-white solid.                                                           |
| Formula          | $C_{30}H_{43}FN_4O_{11}$                                                            |
| MW               | 654.7                                                                               |
| Purity           | ≥95% (TLC)                                                                          |
| Sequence         | Z-Ile-Glu(OMe)-Thr-Asp(OMe)-fluoromethylketone                                      |
| Solubility       | Soluble in DMSO.                                                                    |

## 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](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@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)