

Boc-D(OMe)-FMK

Caspase inhibitor

Irreversible, cell permeable broad-spectrum caspase inhibitor.

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|------------------|-----|
| ALX-260-071-M001 | 1mg |
| ALX-260-071-M005 | 5mg |

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. For long term storage keep aliquots at -70°C. |
| Handling | Keep cool and dry. After reconstitution, prepare aliquots and store at -80°C. |
| Long Term Storage | -20°C |
| Shipping | Ambient Temperature |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|------------------|-----------------------------------------------|
| Alternative Name | Caspase family inhibitor (fluoromethylketone) |
| Appearance | White to off-white powder. |
| CAS | 187389-53-3 |
| Formula | $C_{11}H_{18}FNO_5$ |
| MW | 263.3 |
| Purity | ≥95% (TLC) |
| Sequence | Boc-Asp(OMe)-fluoromethylketone |
| Solubility | Soluble in DMSO or dimethyl formamide. |

Technical Info / Product Notes

The carboxyl groups are OMe ester form 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](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)