

Z-Phe-Ala-FMK

Caspase inhibitor

Irreversible inhibitor of cathepsins B, L, and S, and likely cathepsin H, as well as cruzain and papain. Has been shown to inhibit effector, but not initiator, caspases *in vitro*, and suppress some forms of apoptosis. Has been used *in vivo*.

Citations: 1

[View Online »](#)

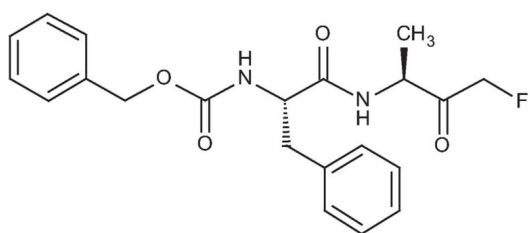
Ordering Information

[Order Online »](#)

BML-PI138-0001	1mg
BML-PI138-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid
CAS	197855-65-5
Formulation	Lyophilized
MW	386.4
Peptide Content	≥95%
Purity	≥95% (HPLC)
Sequence	Benzyloxycarbonyl-Phe-Ala-fluoromethylketone
Solubility	Soluble in DMSO (10mM)

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-eu@enzolifesciences.com)

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