

Mu-Phe-hPhe-FMK

Cathepsin inhibitor

Irreversible, non-toxic, cell-permeable inhibitor of cysteine proteases calpain, cathepsins B and L, and *Trypanosome trypanopain*. Can be used *in vitro* (typically 20 μ M) or *in vivo*. Inhibits tumor formation and growth.

Citations: 7

[View Online »](#)

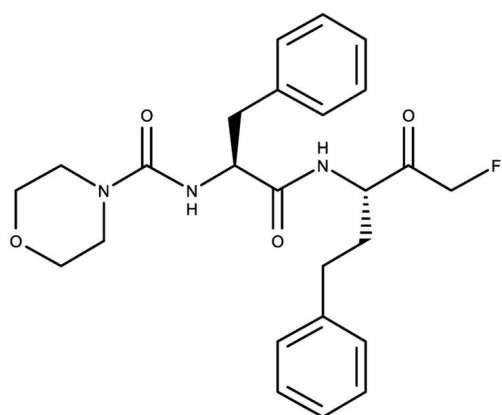
Ordering Information

[Order Online »](#)

BML-EI323-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at 4°C for up to 1 year. Store DMSO concentrates at -20°C for up to 1 year. Do not store dilute solutions.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-Morpholineurea-phenylalanyl-homophenylalanylfluoromethyl ketone
Appearance	White solid.
Formula	$C_{25}H_{30}FN_3O_4$
MW	455.5
Purity	≥98% (TLC)
Sequence	Mu-Phe-hPhe-FMK
Solubility	Soluble in DMSO (>25 mg/ml).



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-
be@enzolifesciences.com](mailto:info-be@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)