

E-64-C

Calpain and cathepsin inhibitor

Cell-impermeable epoxysuccinyl peptide inhibitor of calpain and other cysteine proteases. Inhibits calpain and cathepsins B, H, and L.

Citations: 1

[View Online »](#)

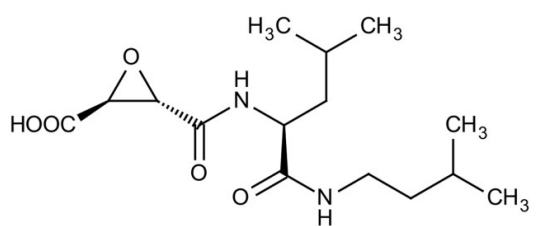
Ordering Information

[Order Online »](#)

BML-PI106-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



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 2S,3S-trans-(Carboxyoxirane-2-carbonyl) -L-leucine-(3-methylbutyl) amide

Appearance White to off-white solid

CAS 76684-89-4

Couple Target Calpain, Cathepsin, Protease

Couple Type Inhibitor

MW 314.4

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (5mg/ml).

Source Synthetic.

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)