

N-Acetyl-Leu-Leu-Methional

Calpain inhibitor

Cell-permeable, peptide aldehyde inhibitor of calpain and other neutral cysteine proteases. Potent inhibitor of cathepsin L ($K_i=0.6\text{nM}$) and cathepsin B ($K_i=100\text{nM}$).

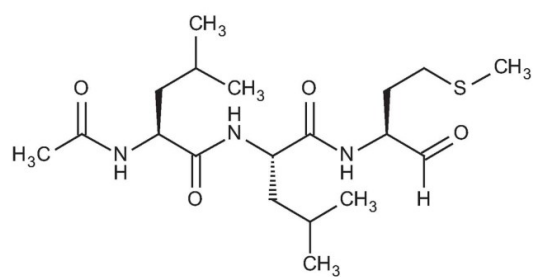
Ordering Information

[Order Online »](#)

BML-PI100-0005	5mg
BML-PI100-0025	25mg

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 ALLM, Ac-LLM-CHO, Calpain inhibitor II

Appearance White to off-white powder.

CAS 110115-07-6

Formula $C_{19}H_{35}N_3O_4S$

MW 401.6

Purity ≥95% (HPLC)

Sequence N-Acetyl-Leu-Leu-Met-H

Solubility 5mg/ml soluble in 100% ethanol, methanol or DMSO

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)