

Z-Leu-Leu-Leu-B(OH)₂

Proteasome inhibitor

A potent cell-permeable proteasome inhibitor ($K_i=0.023$ nM).

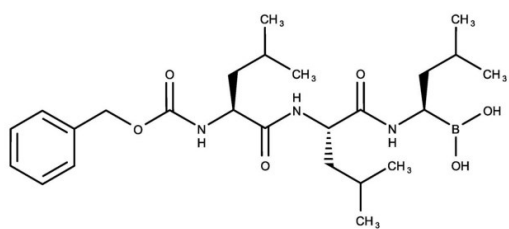
Ordering Information

[Order Online »](#)

BML-PI109-0100	100µg
----------------	-------

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 MG-262, Z-LLL-B(OH)₂, Proteasome inhibitor III

Appearance Solid.

Formula C₂₅H₄₂BN₃O₆

Formulation Lyophilized.

MW 491.4

Purity ≥99% (TLC)

Sequence Z-Leu-Leu-Leu-B(OH)₂

Solubility Soluble in 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)