

MG-132

A potent proteasome inhibitor that blocks protein degradation, crucial for investigating protein turnover and the ubiquitin-proteasome pathway.

MG-132 is a potent, cell permeable and selective proteasome inhibitor ($K_i = 4\text{nM}$).¹ It inhibits NF- κ B activation by preventing I κ B degradation ($IC_{50} = 3\mu\text{M}$). The peptide blocks degradation of short-lived proteins, which in turn induces HSP and ER chaperone expression, leading to thermotolerance ($1\mu\text{M}$ MG-132, 2 h.). It also stimulates neurite outgrowth in PC12 cells (20nM optimal). The peptide has also been reported to increase the survival rate of mesenchymal stem cells following their transplantation. IC_{50} 's for inhibition of Suc-LLVY-AMC and Z-LLL-AMC cleaving activities of proteasome were 0.85 and $0.1\mu\text{M}$ respectively. The ubiquitin-proteasome system (UPS) and autophagy serve as two complementary, reciprocally regulated protein degradation systems, thus blockade of UPS by MG-132 activates autophagy.

Citations: 145

[View Online »](#)

Ordering Information

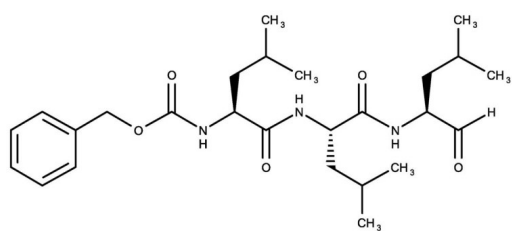
[Order Online »](#)

BML-PI102-0005	5mg
BML-PI102-0025	25mg

Manuals, SDS & CofA

[View Online »](#)

- High purity peptide
- Most economical MG-132 on the market
- Widely cited for more than two decades
- Well-characterized with a variety of applications



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions are stable for up to one week if stored at -20°C. Solutions are stable for up to two months if stored at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Z-LLL-CHO
Appearance	White solid.
CAS	133407-82-6
Formula	$C_{26}H_{41}N_3O_5$
MW	475.6
Purity	≥98%
Sequence	Z-Leu-Leu-Leu-CHO
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (25mg/ml).

Technical Info / Product Notes Replacement for **ADI-HPK-116**

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-be@enzolifesciences.com)

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