

MG-115

Proteasome inhibitor

Potent, reversible proteasome inhibitor. Specifically inhibits the chymotrypsin-like activity of the proteasome. Has been shown to induce apoptosis in Rat-1 and PC12 cells via a p53-dependent pathway. The ubiquitin-proteasome system (UPS) and autophagy serve as two complementary, reciprocally regulated protein degradation systems, thus blockade of UPS by MG-115 activates autophagy.

Citations: 10

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-091-M005	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 Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Z-LL-Nva-CHO, Z-Leu-Leu-Nva-CHO

Appearance White to off-white powder.

CAS 133407-86-0

Formula $C_{25}H_{39}N_3O_5$

MW 461.6

Purity ≥95% (HPLC)

Sequence Z-Leu-Leu-Nva-CHO

Solubility Soluble in DMSO.



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)