

Z-Leu-Leu-Leu-vinyl sulfone

Inhibits the trypsin-like, chymotrypsin-like and caspase-like (PGPH) activity in isolated proteasomes and living cells by covalent modification of the amino-terminal threonine of the catalytically active β subunits. The vinyl sulfone acts as a Michael acceptor for soft nucleophiles such as thiols, leading to the formation of a covalent bond, while being resistant to attack by free thiols, which can cause inactivation of other classes of protease inhibitors.

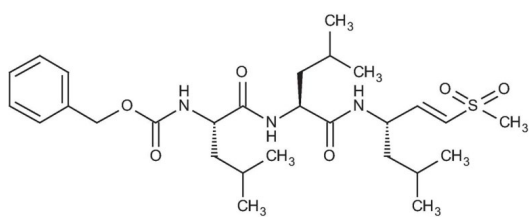
Ordering Information

[Order Online »](#)

BML-ZW9170-0500	500 μ g
-----------------	-------------

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

MW ~551.3 Da

Purity ≥99%

Sequence Z-(Leu)₃-vinyl sulfone



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)