

Bortezomib

Proteasome inhibitor

Potent and selective proteasome inhibitor. Inhibits proliferation of a number of tumor cell lines. Inhibits TNF synthesis and FGF-induced angiogenesis. Reversible. Cell permeable.

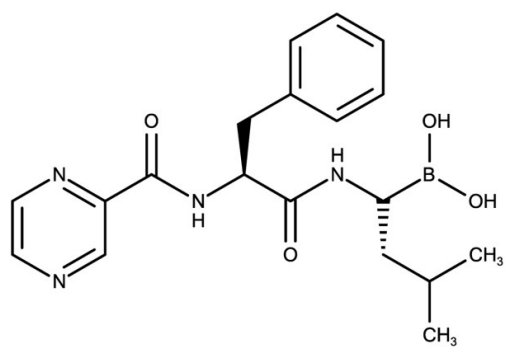
Ordering Information

[Order Online »](#)

ENZ-CHM488-0025	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 Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N-(2-Pyrazine)carbonyl-L-phenylalanine-L-leucine boronic acid

Appearance White solid.

CAS 179324-69-7

Couple Target Proteasome

Couple Type Inhibitor

Formula $C_{19}H_{25}BN_4O_4$

Identity Determined by NMR.

MW 384.25

Purity ≥98% (TLC)

Solubility Soluble in DMSO (50 mg/ml) or ethanol (35 mg/ml).

Last modified: February 27, 2025



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)