

Epoxomicin

Key inhibitor for use in proteasome research.

Cell permeable, potent and selective proteasome inhibitor originally isolated from *Actinomycetes* strain based on its potent *in vivo* antitumor activity. More potent inhibitor of the chymotrypsin-like activity of the proteasome than lactacystin (Prod. No. BML-PI104). Blocks also trypsin-like and PGPH activities of the proteasome. Regulates antigen presentation at non-toxic doses. Effectively inhibits NF- κ B activation *in vitro* and potently blocks inflammation *in vivo* in the mouse ear edema assay. The ubiquitin-proteasome system (UPS) and autophagy serve as two complementary, reciprocally regulated protein degradation systems, thus blockade of UPS by Epoxomicin activates autophagy.

Citations: 38

[View Online »](#)

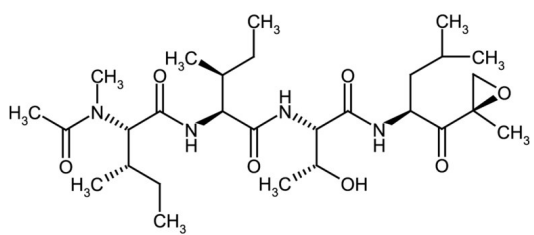
Ordering Information

[Order Online »](#)

BML-PI127-0100	100µg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Lyophilized. Very thin transparent film.
CAS	134381-21-8
Couple Target	Proteasome
Couple Type	Inhibitor
Formula	$C_{28}H_{50}N_4O_7$
MI	14: 3630
MW	554.7
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (15mg/ml) or dichloromethane:methanol (9:1); insoluble in water.

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-
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)