

clasto-Lactacystin β -Lactone

Proteasome inhibitor

Highly specific, cell permeable, and irreversible 20S proteasome inhibitor. Inhibits ubiquitin proteasome pathway in cell culture ($IC_{50}=1\mu M$). It has been suggested that the natural product lactacystin acts as a precursor for *clasto*-lactacystin β -lactone and that the latter is the sole species that interacts with the proteasome. Inhibits cathepsin A. Leads to a reduction of adipogenesis. Induces neurite growth and inhibits cell cycle progression.

Citations: 14

[View Online »](#)

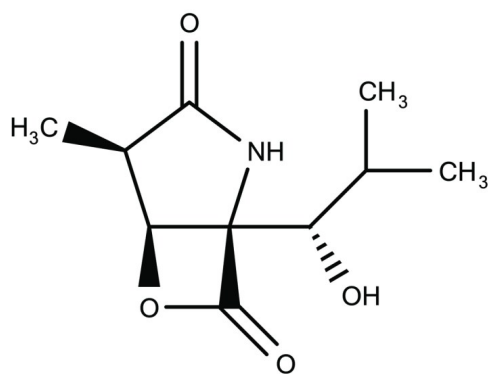
Ordering Information

[Order Online »](#)

BML-PI108-0100	100 μ g
----------------	-------------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 2 years after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 1 month at -80°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Omuralide
Appearance	White solid.
CAS	155975-72-7
Couple Target	Cathepsin, Proteasome 20S
Couple Type	Inhibitor
Formula	$C_{10}H_{15}NO_4$
MW	213.2
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (25mg/ml) or acetonitrile.



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)