

# clasto-Lactacystin $\beta$ -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  $\beta$ -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 »](#)

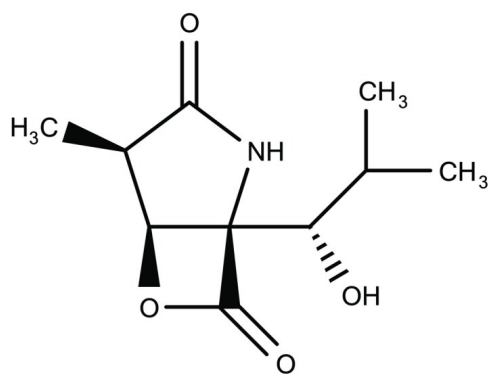
## Ordering Information

[Order Online »](#)

BML-PI108-0100	100 $\mu$ 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	154226-60-5
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.

Last modified: September 15, 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-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-fr@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)