

# CA-074 Me

## Cathepsin inhibitor

Potent and selective irreversible cell permeable cathepsin B inhibitor. Methyl ester (Me) is hydrolyzed by intracellular esterases releasing the active inhibitor. Prevents death of CA1 neurons after ischemia following IV delivery after the ischemic insult. In isolated rat osteoclasts it inhibited bone resorption with a maximal effect at 50  $\mu$ M. This product is only suitable for in vivo and whole cell experiments. The free acid form must be used for in vitro experiments. Can be used to discriminate between cathepsin B and L/S forms.

Citations: 16

[View Online »](#)

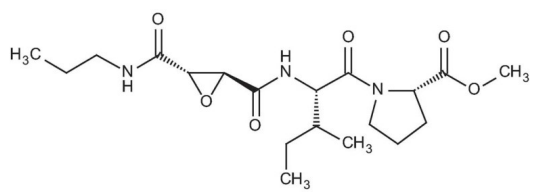
## Ordering Information

[Order Online »](#)

BML-PI126-0001	1mg
----------------	-----

## 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

**Alternative Name** N-(L-3-trans-Propylcarbonyl-oxirane-2-carbonyl)-L-isoleucyl-L-proline methyl ester

**Appearance** White solid.

**Formula**  $C_{19}H_{31}N_3O_6$

**MW** 397.5

**Purity** ≥98% (HPLC)

**Sequence** N-(L-3-trans-Propylcarbonyl-oxirane-2-carbonyl)-L-isoleucyl-L-proline methyl ester

**Solubility** Soluble in DMSO.



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)