

CA-074

Cathepsin B inhibitor

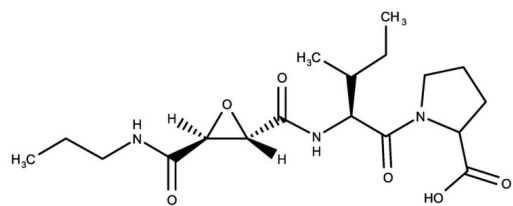
Potent and specific inhibitor of cathepsin B *in vitro* and *in vivo*. Free acid form used *in vitro* or in isolated enzyme studies but not in whole cells. Can discriminate between cathepsin B and L/S forms.

Citations: 4

[View Online »](#)

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-[L-3-trans-Propylcarbamoyloxirane-2-carbonyl]-Ile-Pro-OH
Appearance	White to off-white powder.
CAS	134448-10-5
Formula	$C_{18}H_{29}N_3O_6$
MW	383.5
Purity	≥98% (HPLC)
Solubility	Soluble in ethyl acetate (50mg/ml) or DMSO.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com