

# TAPI-1

## Inhibitor of TNF- $\alpha$ processing

Inhibitor of TNF- $\alpha$  processing ( $IC_{50}$  is approximately 100 nM using *in vitro* TNF peptide cleavage assay), similar to TAPI-0, but more stable in tissue culture and in vivo. Inhibits TACE and other metalloproteases. Blocks release of TNF- $\alpha$  ( $IC_{50}$ =50-100  $\mu$ M), APP (1-8  $\mu$ M), ACE (10  $\mu$ M), IL6R (5-10  $\mu$ M), TNFRI (5-10  $\mu$ M), TNFRII (25-50  $\mu$ M), and ObRa (leptin receptor), but not CSF-1, from cells in culture.

Citations: 19

[View Online »](#)

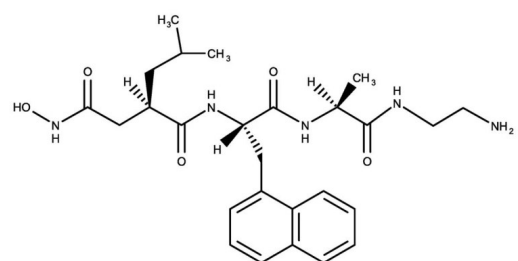
## Ordering Information

[Order Online »](#)

BML-PI134-0001	1mg
----------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Aliquot stock solutions and freeze at -20°C for up to one month. Do not store diluted solutions.
Handling	Hygroscopic. Protect from light.
Long Term Storage	-20°C
Shipping	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	IC2, N-(R)-(2-(Hydroxyaminocarbonyl)methyl)-4-methylpentanoyl-L-naphthylalanyl-L-alanine 2-aminoethyl amide
Appearance	Lyophilized solid.
Formula	$C_{26}H_{37}N_5O_5$
MW	499.3
Purity	≥98% (HPLC)
Sequence	N-(R)-(2-(Hydroxyaminocarbonyl)methyl)-4-methylpentanoyl-L-naphthylalanyl-L-alanine 2-aminoethyl amide
Solubility	Soluble in DMSO or methanol

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)