

TAPI-1

Inhibitor of TNF- α processing

Inhibitor of TNF- α 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- α (IC_{50} =50-100 μ M), APP (1-8 μ M), ACE (10 μ M), IL6R (5-10 μ M), TNFRI (5-10 μ M), TNFRII (25-50 μ M), and ObRa (leptin receptor), but not CSF-1, from cells in culture.

Citations: 19

[View Online »](#)

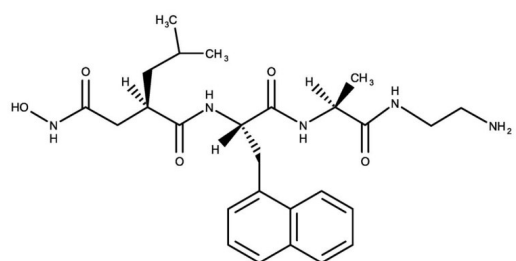
Ordering Information

[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



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)