

TAPI-2

Inhibitor of TNF- α processing

Inhibitor of TACE, other ADAMs, ACE secretase and other metalloproteinases. Blocks release of TNF α , TGF- α , DCC, APP, L-selectin, ACE, IL6R, NOTCH, and erbB2/HER2. Has been used in culture (typical concentration 5-100 μ M).

Citations: 23

[View Online »](#)

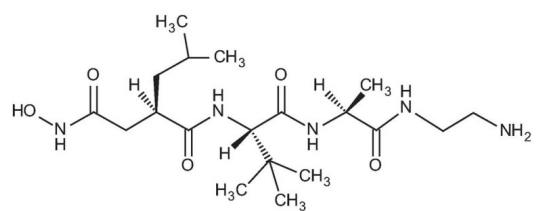
Ordering Information

[Order Online »](#)

BML-PI135-0001	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 IC3, N-(R)-(2-(Hydroxyaminocarbonyl)methyl)-4-methylpentanoyl-L-*t*-butyl-glycine-L-alanine 2-aminoethyl amide

Appearance Lyophilized solid.

Formula $C_{19}H_{37}N_5O_5$

MW 415.28

Purity ≥99% (HPLC)

Sequence (IC-3)
N-(R)-(2-(Hydroxyaminocarbonyl)methyl)-4-methylpentanoyl-L-*t*-butyl-glycine-L-alanine 2-aminoethyl amide

Solubility Soluble in water (at least 20mM) or 100% ethanol (at least 5mg/ml)



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)