

GM-6001

MMP inhibitor

Potent broad-spectrum hydroxamate inhibitor of matrix metalloproteinases (MMPs). K_i values have been reported for the following human MMPs: MMP-1 (fibroblast collagenase): 0.4nM; MMP-2 (gelatinase A): 0.5nM; MMP-3 (stromelysin-1): 27nM; MMP-7 (matrilysin): 3.7nM; MMP-8 (neutrophil collagenase): 0.1nM; MMP-9 (gelatinase B): 0.2nM; MMP-12 (metalloelastase): 3.6nM; MMP-14 (MT1-MMP): 13.4nM; MMP-26 (endometase): 0.36nM. Also inhibits MMP-10, MMP-13, MMP-15, MMP-17, MMP-20, MMP-21, TACE, ADAM19, other ADAMs, anthrax lethal factor, neprilysin, leucine aminopeptidase, and DPPIII. Typical concentration range for use in tissue culture is 10-25 μ M. This inhibitor has also been used *in vivo*.

Citations: 41

[View Online »](#)

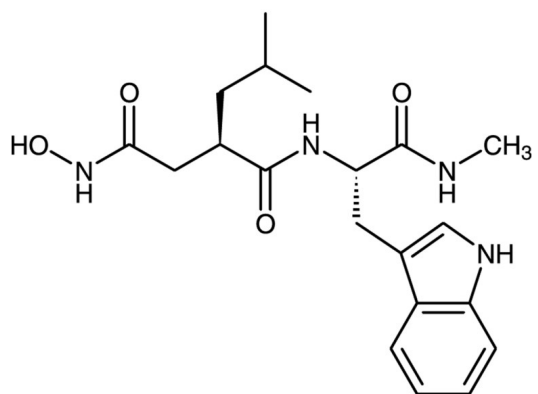
Ordering Information

[Order Online »](#)

BML-EI300-0001	1mg
BML-EI300-0005	5mg

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 Galardin, Ilomastat, N-[(2R)-2-(Hydroxamidocarbonylmethyl)-4-methylpentanoyl]-L-tryptophan methylamide

Appearance White to off-white powder.

CAS 142880-36-2

Couple Target ADAM, DPP, MMP

Couple Type Inhibitor

Formula $C_{20}H_{28}N_4O_4$

Formulation Lyophilized solid.

MW 388.5

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (at least 25mM) or 100% ethanol (10mg/ml)

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-be@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