

MMP inhibitor

Potent inhibitor of human stromelysin-1 (MMP-3), $K_i=0.13 \mu\text{M}$. Also inhibits MMP-1, MMP-2, MMP-7, MMP-8, MMP-9, MMP-10, MMP-12 (metalloelastase), MMP-13, MMP-14, and MMP-20.

Citations: 7

[View Online »](#)

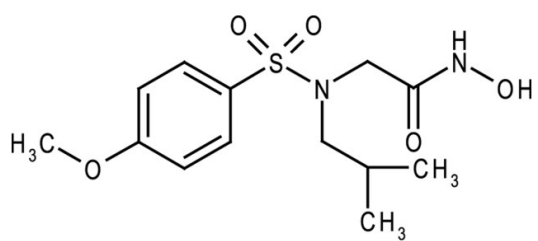
Ordering Information

[Order Online »](#)

BML-PI115-0005	5mg
BML-PI115-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N-Isobutyl-N-(4-methoxyphenylsulfonyl)glycyl hydroxamic acid

Appearance White solid.

CAS 161314-17-6

Couple Target MMP

Couple Type Inhibitor

Formula $C_{13}H_{20}N_2O_5S$

MW 316.4

Purity $\geq 99\%$ (TLC)

Solubility Soluble in DMSO (30mg/ml) or 100% ethanol (8mg/ml).

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-
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)