

Gabexate . mesylate

Protease inhibitor

A competitive, reversible inhibitor of serine proteases tryptase ($K_i=3.4\text{nM}$), thrombin ($K_i=500\text{nM}$), uPA ($K_i=1.3\mu\text{M}$), plasmin ($K_i=1.5\mu\text{M}$), trypsin ($K_i=2\mu\text{M}$), factor Xa ($K_i=2\mu\text{M}$). At higher concentrations it inhibits kallikrein-1 ($K_i=2.6\text{mM}$), C_1 esterase ($\text{IC}_{50}=520\mu\text{M}$), as well as phospholipase A_2 ($\text{IC}_{50}=500\mu\text{M}$), rat iNOS ($K_i=5\text{mM}$) and cNOS ($K_i=150\mu\text{M}$), and swine kidney copper amine oxidase ($K_i=27\mu\text{M}$), but not chymotrypsin. Has been used in tissue culture and *in vivo*. In clinical use for pancreatitis.

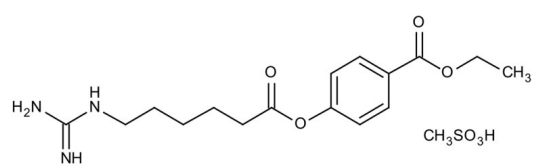
Ordering Information

[Order Online »](#)

BML-PI144-0010	10mg
BML-PI144-0050	50mg

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 Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White solid.

CAS 56974-61-9

Couple Target C1 esterase, Clotting factor, Kallikrein, NOS, Phospholipase, PLA, Plasmin, Protease, Serine protease, Thrombin, Trypsin, Tryptase, uPA

Couple Type Inhibitor

Formula $C_{16}H_{23}N_3O_4 \cdot CH_3SO_3H$

MW 417.5

Purity ≥98% (HPLC)

Solubility Soluble in water (25 mg/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)