

Nafamostat . mesylate

Tryptase inhibitor

Competitive, reversible broad-spectrum serine proteinase inhibitor that can be used as a specific tryptase inhibitor at very low concentrations. Has been used *in vitro* and *in vivo*.

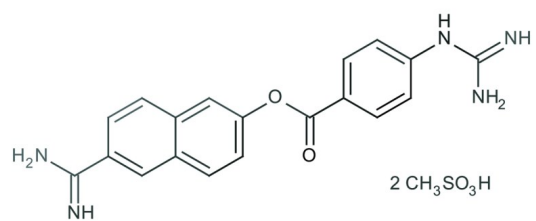
Ordering Information

[Order Online »](#)

BML-PI143-0010	10mg
BML-PI143-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Hygroscopic.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White crystalline solid.
CAS	82956-11-4
Couple Target	Serine protease, Tryptase
Couple Type	Inhibitor
Formula	$C_{19}H_{17}N_5O_2 \cdot 2CH_4O_3S$
MW	539.6
Purity	≥98% (HPLC)
Solubility	Soluble in water (25 mg/ml) Note: Phosphate-containing buffers may cause precipitation.

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)