

Sivelestat sodium

Elastase inhibitor

Potent and selective leukocyte elastase inhibitor ($IC_{50}=44nM$). Displays various anti-inflammatory activities including attenuation of LTB_4 -induced neutrophil migration and serum cytokine levels. Reduces acute inflammatory lung injury *in vivo*.

Citations: 3

[View Online »](#)

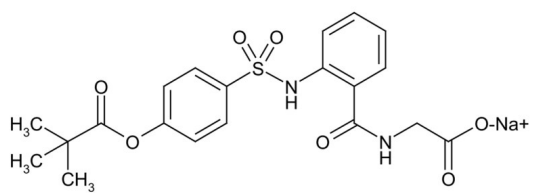
Ordering Information

[Order Online »](#)

BML-PI157-0010	10mg
BML-PI157-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at room temperature. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	ONO-5046, N-{2-[(4-[2,2-Dimethylpropanoyl]oxy)phenyl]sulfonyl}amino]benzoyl}glycine sodium salt
Appearance	White solid.
CAS	201677-61-4
Couple Target	Elastase
Couple Type	Inhibitor
Formula	$C_{20}H_{21}N_2NaO_7S \cdot 4H_2O$
Identity	Determined by NMR.
MW	528.5
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (>25mg/ml), ethanol (20mg/ml) or water (2mg/ml, warm).



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-fr@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