

PKSI-527

Kallikrein inhibitor

Highly selective plasma kallikrein inhibitor ($K_i = 0.81 \mu\text{M}$). Can be used for affinity purification of kalikrein. Suppresses collagen-induced arthritis in a mouse model. Synergizes with NO donors producing antithrombotic effects *ex vivo*.

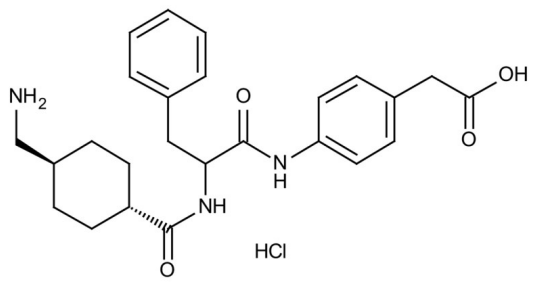
Ordering Information

[Order Online »](#)

BML-PI147-0005	5mg
BML-PI147-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store as supplied at -20°C for up to 1 year. Solutions in distilled water may be stored at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	trans-4-Aminomethylcyclohexanecarbonyl-phenylalanyl-4-carboxymethylanilide
Appearance	White solid.
CAS	128837-71-8
Couple Target	Kallikrein
Couple Type	Inhibitor
Formula	$C_{25}H_{31}N_3O_4 \cdot HCl$
Identity	Determined by NMR.
MW	474.0
Purity	≥95% (HPLC)
Solubility	Soluble in water (5 mg/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