

Phosphoramidon

Neprilysin inhibitor

Inhibits neutral endopeptidase 24.11 (NEP, neprilysin; EC 3.4.24.11) ($K_i = 2\text{nM}$), a cell surface zinc metalloprotease which cleaves bioactive peptides such as bradykinin, neurotensin, and enkephalin. Also inhibits neprilysin 2 ($K_i \sim 2\text{nM}$). Potent inhibitor of endothelin-converting enzyme (ECE) ($IC_{50} = 0.68\mu\text{M}$).

Citations: 10

[View Online »](#)

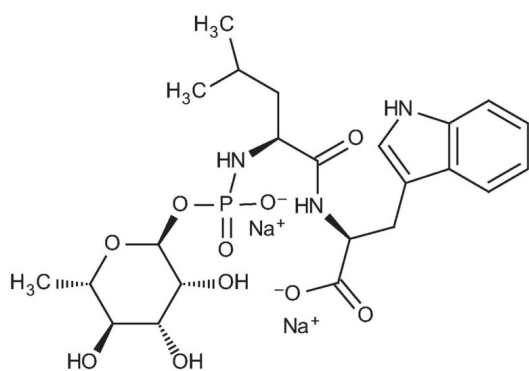
Ordering Information

[Order Online »](#)

BML-PI111-0005	5mg
BML-PI111-0025	25mg

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 at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-(α-Rhamnopyranosyloxyhydroxyphosphinyl)-L-leucyl-L-tryptophan
Appearance	White to slightly yellow solid.
CAS	119942-99-3
Formula	$C_{23}H_{34}N_3O_{10}P \cdot 2Na$
MW	541.5 . 46.0
Purity	≥97% (HPLC)
Sequence	N-(α-Rhamnopyranosyloxyhydroxyphosphinyl)-Leu-Trp . 2Na
Solubility	Soluble in water, DMSO or methanol.

Technical Info / Product Notes	Note: Product is not sterile.
---------------------------------------	-------------------------------



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)