Phosphoramidon

Neprilysin inhibitor

Inhibits neutral endopeptidase 24.11 (NEP, neprilysin; EC 3.4.24.11) (K $_{\rm i}$ =2nM), a cell surface zinc metalloprotease which cleaves bioactive peptides such as bradykinin, neurotensin, and enkephalin. Also inhibits neprilysin 2 (K $_{\rm i}$ ~2nM). Potent inhibitor of endothelin-converting enzyme (ECE) (IC $_{\rm 50}$ =0.68µM).

Citations: 10

View Online »

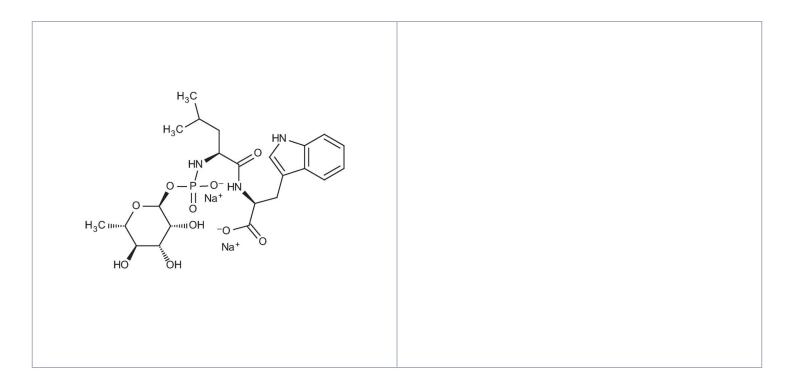
Ordering Information

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

White to slightly yellow solid. **Appearance**

CAS 119942-99-3

Formula $C_{23}H_{34}N_3O_{10}P$. 2Na

MW 541.5 . 46.0

Purity ≥97% (HPLC)

 $N-(\alpha-Rhamnopyranosyloxyhydroxyphosphinyl)-Leu-Trp$. Sequence

2Na

Solubility Soluble in water, DMSO or methanol.

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

Last modified: May 29, 2024

info-

