

PRODUCT DATA SHEET



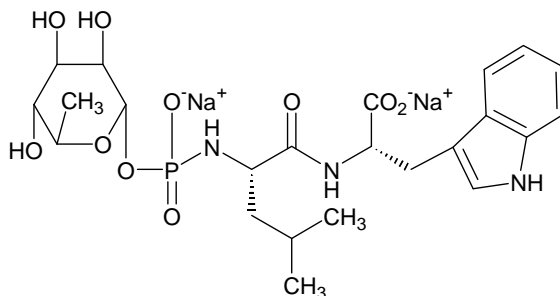
PRODUCT: Phosphoramidon
N-(α -Rhamnopyranosyloxyhydroxy-phosphinyl)-Leu-Trp \cdot 2Na

CAS NO. 119942-99-3

CATALOG NO.: PI-111

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: White solid

MOLECULAR FORMULA: $C_{23}H_{34}N_3O_{10}P \cdot 2Na$

MOLECULAR WEIGHT: 587.5

PURITY: 97%

SOLUBILITY: Soluble in water, DMSO and methanol

STORAGE: Store, as supplied, at $-20^{\circ}C$ for up to 1 year. Store solutions at $-20^{\circ}C$ for up to 3 months.

APPLICATION NOTES: Inhibits neutral endopeptidase 24.11 (NEP; Neprilysin; EC 3.4.24.11) ($K_i=2$ nM)¹ - a cell surface zinc metalloprotease that hydrolyzes bioactive peptides such as bradykinin², neurotensin³ and enkephalin⁴. Also inhibits neprilysin and neprilysin 2 ($K_i \sim 2$ nM)^{1,5}. Potent inhibitor of endothelin-converting enzyme ($IC_{50} = 0.68$ μ M)⁶.

- REFERENCES:
1. C. Rose *et al.* *Biochem. J.* 2002 **363** 697
 2. X. Zhang *et al.* *J. Cardiovasc. Pharmacol.* 1998 **31** 623
 3. T.W. Moody *et al.* *Peptides* 1998 **19** 253
 4. T. Hiranuma *et al.* *J. Pharmacol. Exp. Therap.* 1997 **281** 769
 5. A.R. Whiteside and A.J. Turner *FEBS Lett.* 2008 **582** 2382
 6. K. Ahn *et al.* *Arch. Biochem. Biophys.* 1998 **359** 258

The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.