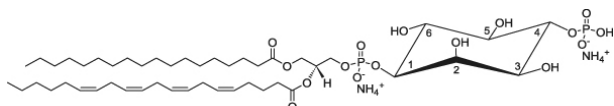


PI-4-P

L- α -Phosphatidylinositol-4-phosphate porcine brain

Cat. No.	Amount
LI-002	1 mg



For *in vitro* use only
 Quality guaranteed for 12 months
 Store at -20°C

Form

Lyophilized, Diammonium Salt.

Molecular Formula

$C_{47}H_{90}N_2O_{16}P_2$

Molecular Weight

1001.18 g/mol

Purity

≥99%

Description

Phosphatidylinositol-4-phosphate is one of the *in vitro* substrates for PI3-kinases. Phosphatidylinositol 3-kinase is a lipid kinase that phosphorylates the inositol ring of phosphatidylinositol-4-phosphate and related compounds at the 3'-OH position.

PIP can be used as a substrate for PI3K lipid kinase activity assay together with a mix of PI, PIP₂, PC, PE, PS and SM.

Selected References:

- Vanhaesebroeck *et al.* (2001) Synthesis and function of 3-phosphorylated inositol lipids. *Ann. Rev. Biochem.* **70**:535.
 Balla (2001) Pharmacology of phosphoinositides, regulators of multiple cellular functions. *Curr. Pharm. Des.* **7**:475.
 Wymann (2003) Phosphoinositide 3-kinase signalling – which way to target? *Trend Pharmacol. Sci.* **24**:323.
 Foukas *et al.* (2002) Direct effects of caffeine and theophylline on p110 delta and other phosphoinositide 3-kinases. Differential effects on lipid kinase and protein kinase activities. *J. Biol. Chem.* **277**:37124.