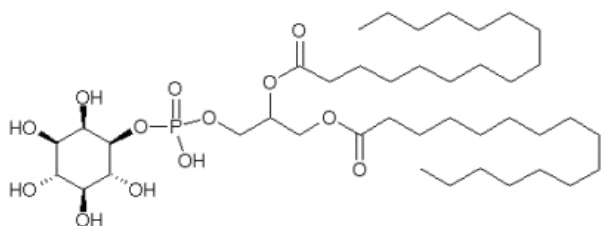


## PI synthetic L- $\alpha$ -Phosphatidylinositol synthetic

Cat. No.	Amount
LI-008	100 $\mu$ g



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

**Form**  
 Lyophilized.

**Solubility**  
 Soluble in water.

**Molecular Formula**  
 $C_{41}H_{79}O_{13}P$ .

**Molecular Weight**  
 811.03 g/mol

**Purity**  
 $\geq 98\%$

### Description

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

PI can be used as a substrate for PI3K lipid kinase activity assay together with a mix of PIP, PIP<sub>2</sub>, PC, PE, PS and SM.

Di-C18PtdIns contains two chains of the 18:0 saturated fatty acid Octadecanoic Acid (common name: stearic acid).

### 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.