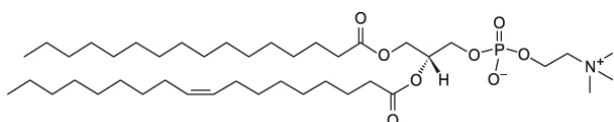


PC

L- α -Phosphatidylcholine

chicken egg

Cat. No.	Amount
LI-004	5 mg



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

Form

Lyophilized, sodium salt.

Solubility

Soluble in Chloroform, Hexane-ethanol, Methanol, Ethanol, Toluene, Ether.
 Insoluble in Water (CMC < 0.001 nM), cold Acetone.

Molecular Formula

C₄₂H₈₂NO₈P.

Molecular Weight

760.09 g/mol

Purity

≥99%

Description

Phosphatidylcholine is a major phospholipid in the outer plasma membrane of mammalian cells.

PC together with phosphatidylethanolamine is the substrate for PS synthase-1 and 2 to synthesize phosphatidylserine by a base-exchange reaction.

It is used by Phospholipase D to generate the second messenger phosphatidic acid (PA).

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

Selected References:

- Wymann (2003) Phosphoinositide 3-kinase signalling – which way to target? *Trend Pharmacol. Sci.* **24**:323.
 Ramoni *et al.* (2004) Phosphatidylcholine-specific phospholipase C in mitogen-stimulated fibroblasts. *Exp. Cell Res.* **299**:370.
 Farese (1996) Insulin-sensitive phospholipid signaling systems and glucose transport: an update. *Proc. Soc. Exp. Biol. Med.* **213**:1.
 Martelli *et al.* (2004) Metabolism and signaling activities of nuclear lipids. *Cell Mol. Life Sci.* **61**:1143.
 Mozzi *et al.* (2003) Metabolism and functions of phosphatidylserine in mammalian brain. *Neurochem Res.* **28**:195.