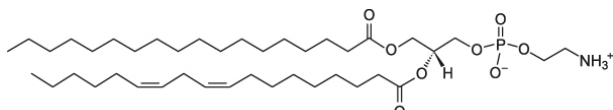


PE

L- α -Phosphatidylethanolamine chicken egg

Cat. No.	Amount
LI-005	5 mg



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

Form

Lyophilized.

Solubility

Soluble in Chloroform.
Insoluble in Acetone, Water.

Molecular Formula

C₄₁H₇₈NO₈P.

Molecular Weight

744.05 g/mol

Purity

≥99%

Description

Phosphatidylethanolamine together with phosphatidylcholine is the substrate for PS synthase 1 and 2 to synthesize phosphatidylserine by a base-exchange reaction. PE can be used in a lipid mix for PI3K lipid kinase activity assay together with PI, PIP, PIP₂, PS, PC and SM.

Selected References:

Wymann (2003) Phosphoinositide 3-kinase signalling – which way to target? *Trend Pharmacol. Sci.* **24**:323.
Vance (2003) Molecular and cell biology of phosphatidylserine and phosphatidylethanolamine metabolism. *Prog. Nucleic Acid Res. Mol. Biol.* **75**:69.