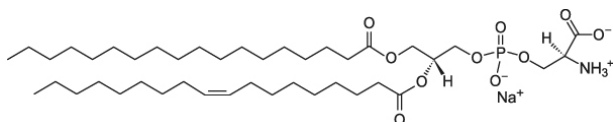


PS

L- α -Phosphatidylserine porcine brain, sodium salt

Cat. No.	Amount
LI-006	5 mg



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

Form

Lyophilized, sodium salt.

Solubility

Soluble in Chloroform, Toluene.
Insoluble in Ethanol, Methanol, Water.

Molecular Formula

C₄₂H₇₉NO₁₀PNa.

Molecular Weight

812.05 g/mol

Purity

≥99%

Description

Phosphatidylserine (PS) is involved in cell signaling and apoptosis.

In higher eukaryotes, PS is synthesized from phosphatidylethanolamine or phosphatidylcholine through the exchange of the nitrogen base with free serine.

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

Selected References:

Wymann (2003) Phosphoinositide 3-kinase signalling – which way to target? *Trend Pharmacol. Sci.* **24**:323.

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.

Suzuki *et al.* (2009) Biochemical characterization of tau protein and its associated syndapin 1 and protein kinase C for their functional regulation in rat brain. *Biochimica et Biophysica Acta (BBA) - General Subjects* **1790(3)**:188-197.