

Product Information

Description:	PI/PS (4:1, w/w) unilamellar vesicle, prepared with L- α phosphatidylinositol (Liver, Bovine) (sodium salt) and L- α -phosphatidylserine (Brain, Porcine) (sodium salt). Lipid monolayer was formed by sonication. MW PI= 909.12, MW PS= 812.05.
Size:	500 μ l at 0.625 mg/ml
Formulated In:	50 mM HEPES, pH 7.5 and 1 mM EGTA
Format:	Lipid
Stability:	2 to 3 months at +4°C. Do not freeze.
Storage:	+4°C
Assay Conditions:	Enzyme reaction was performed at 30°C for 15 min utilizing 100 μ M ATP and PI/PS unilamellar vesicles (0.2 mg/ml PI at final). Luminescence signal was generated using ADP-Glo Kinase Assay (Promega) reagents.
Applications:	Useful as a substrate for PI-3 kinase activity assays.

Quality Control Data

PI3K Substrate Activity

