

## Data Sheet

### ***PI-3 Kinase Lipid Substrate***

Unilamellar vesicle of phosphatidylinositol and phosphatidylserine

**Catalog #:** 40560

**Lot#:** 180731

**Size:** 1 mg @ 0.625 mg/ml

---

**Formulation:** 50 mM HEPES, pH 7.5 and 1 mM EGTA

**Stability:** 2 to 3 months at +4°C. **Do not freeze.**

**References:**

Vidugiriene J *et al.*, *Assay and Drug Development Technologies*, 2009 Dec; **7(6)**; 585-597.

**Description:**

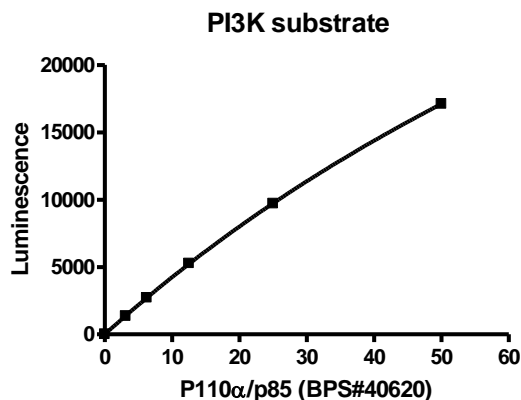
PI/PS (4:1, w/w) unilamellar vesicle, prepared with L- $\alpha$ -phosphatidylinositol (Liver, Bovine) (sodium salt) and L- $\alpha$ -phosphatidylserine (Brain, Porcine) (sodium salt). Lipid monolayer was formed by sonication. MW PI= 909.12, MW PS= 812.05.

**Application:**

Useful as a substrate for PI-3 kinase activity assays.

---

### **Quality Assurance**



Enzyme reaction was performed at 30°C for 15 min utilizing 100  $\mu$ 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.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082**, Fax **1.858.481.8694**  
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)