

## Data Sheet

### **PI3 Kinase [p110 $\alpha$ (E545K)/p85 $\alpha$ ], His-Tag**

Human, Recombinant, N-terminal His-Tag

**Catalog #:** 40644

**Lot #:** 161215      **Conc.:** 0.05 mg/ml

**Formulation:** 50 mM sodium phosphate, pH 7.0, 300 mM NaCl, 150mM imidazole, 0.1 mM PMSF, 0.25 mM DTT, 25% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

**References:**

1. Samuels, Y., *et al.* *Science* **304**: 554 (2008)
2. Graupera, M., *et al.*, *Nature* **453**: 662-666 (2008)

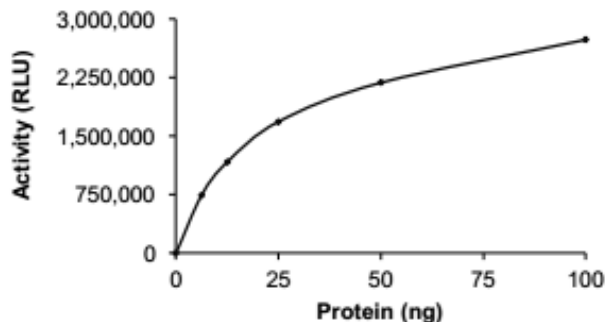
**Description:**

Complex of recombinant full-length human p110 $\alpha$  (GenBank Accession No. NM\_006218) with E545K mutation and N-terminal His-Tag, and recombinant full length, human p85 $\alpha$  (GenBank Accession No. NM\_181523), no tag. Co-expressed in Sf9 cells in a Baculovirus expression system. p110 $\alpha$  (E545K) MW=111 kDa, p85 $\alpha$  MW=86 kDa.

**Application:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## **Quality Assurance**

**Specific Activity**



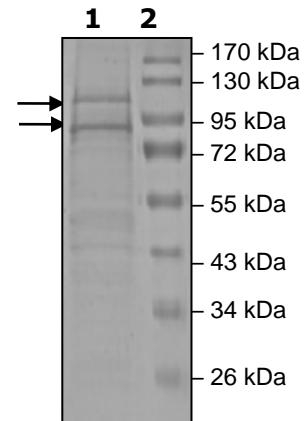
**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
p110a(E545K)  
and p85a

**Lane 2:**  
Protein Marker

**MW:**  
111 kDa, 110a(E545K)  
86 kDa, p85a

**Purity:**  $\geq 70\%$



OUR PRODUCTS ARE FOR RESEARCH 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)