

## Data Sheet

### **PI3 Kinase [p110α (N345K)/p85α], His-Taq**

Human, Recombinant, N-terminal His-Tag  
**Catalog #:** 40646  
**Lot #:** 161216      **Conc.:** 0.05 mg/ml

**Formulation:** 50 mM sodium phosphate, pH 7.0, 300 mM NaCl, 150 mM 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)

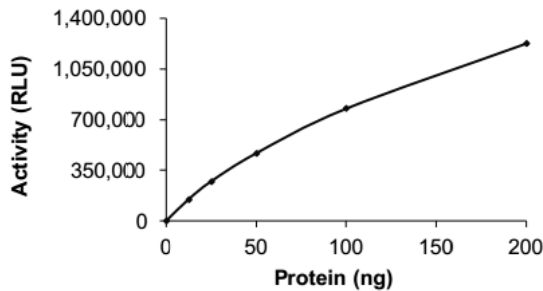
**Description:**

Complex of recombinant full-length human p110α (GenBank Accession No. NM\_006218.2) with N345K mutation and N-terminal His-Tag, and recombinant full length, human p85α (GenBank Accession No. NM\_181523), no tag. Co-expressed in Sf9 cells in a Baculovirus expression system. p110α (N345K) MW=111 kDa, p85α 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(N345K)  
 and p85a

**Lane 2:**  
 Protein Marker

**MW:**  
 111 kDa, p110a(N345K)  
 86 kDa, p85a

**Purity:** ≥80%

