

## Data Sheet

### p110 $\alpha$ (N345K)/p85 $\alpha$

Human, Recombinant, Complex

**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.

**References:**

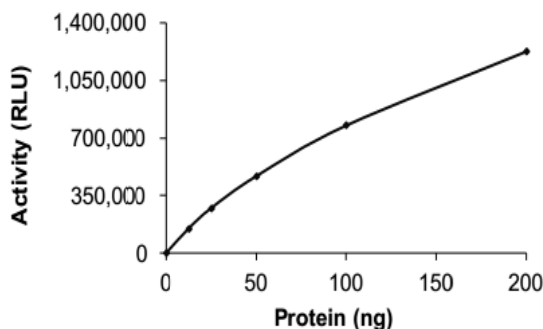
1. Samuels, Y., *et al.* *Science* **304**: 554 (2008)
2. Graupera, M., *et al.*, *Nature* **453**: 662-666 (2008)

**Description:** Complex of human p110 $\alpha$ , also known as PIK3CA, GenBank Accession No. NM\_006218, a.a. 2-1068(full length), with N345K mutation, with N-terminal His-tag, MW=111 kDa, and human p85 $\alpha$ , also known as PIK3R1, GenBank Accession No. NM\_181523, a.a. 1-724(full length), MW=85 kDa. Coexpressed in a Baculovirus infected Sf9 cell expression system.

**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)/p85a

**Lane 2:**  
Protein Marker

**MW:**  
p110a(N345K): 111 kDa  
p85a: 85 kDa

**Purity:**  $\geq 80\%$

