

## Data Sheet

### p110 $\alpha$ /p85 $\alpha$

Human, recombinant, GST-tag  
**Catalog #:** 40621  
**Lot#:** 100426 **Conc.:** 0.3 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0,  
 124 mM NaCl, 2.4 mM KCl, 18 mM  
 glutathione, 10% glycerol, and 3 mM DTT.

**Stability:** >6 months at -80°C

#### References:

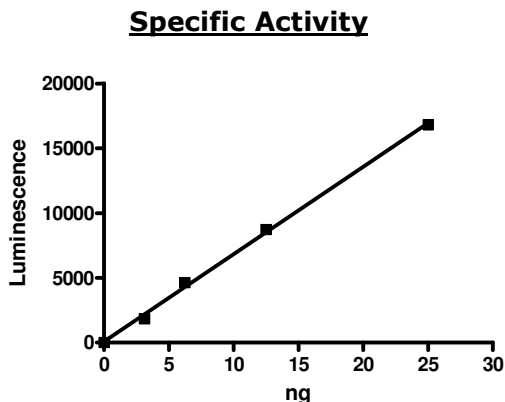
1. Vogt, PK *et al.*, *Virology* **344(1)**:131-  
 (2006)
2. Bader, AG *et al.*, *Nat Rev Cancer*  
**5(12)**:921-9 (2005)

**Description:** Complex of N-terminal GST-tagged recombinant full-length human p110 $\alpha$  (GenBank Accession No. U79143), and recombinant full length, human p85 $\alpha$  (no tag) (GenBank Accession No. XM\_043865). Coexpressed in a Baculovirus infected Sf9 cell expression system. p110 $\alpha$  MW=151 kDa, p85 $\alpha$  MW=88 kDa.

**Specific Activity:** 756 pmol/min/ $\mu$ g.  
 A 25  $\mu$ l kinase reaction is conducted in a buffer containing 40 mM Tris-HCl (pH 7.4), 20 mM MgCl<sub>2</sub>, 100  $\mu$ M ATP, 200  $\mu$ M PI:PS substrate (BPS #40560) and enzyme for 15 min at 30°C. ATP reduction is detected using ADP-Glo™ Kinase Assay Platform (Promega Corporation, Madison, WI).

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

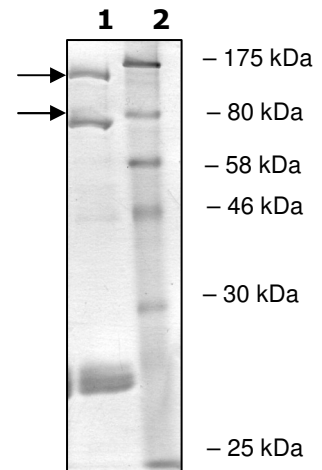
### Quality Assurance



#### 10% SDS-PAGE Coomassie staining

**Lane 1:**  
p110 $\alpha$ /p85 $\alpha$   
**Lane 2:**  
Protein Marker  
BioLabs (#P7708L)

**MW:**  
p110 $\alpha$ : 151 kDa  
p85 $\alpha$ : 88 kDa  
**Purity:**  $\geq$ 38%



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)