

Data Sheet

p110β/p85α

Human, recombinant

Catalog #: 40622

Lot#: 110927

Conc.: 0.13 mg/ml

Formulated in: 25 mM Tris-HCl, pH 8.0, 69 mM NaCl, 1.35 mM KCl, 0.025 Tween-20, 0.05mg/ml Flag peptide, 3mMDTT, 50% Glycerol.

Stability: >6 months at -80°C

References:

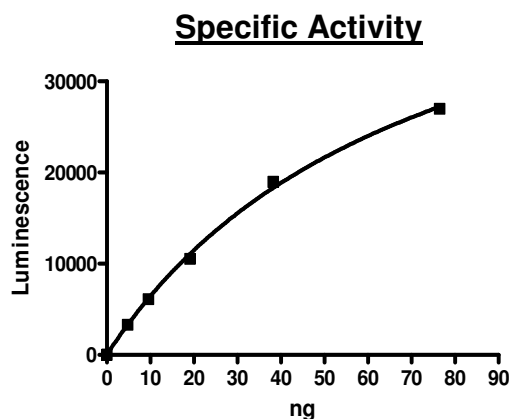
1. Marques M. *et al.*, (2008). *Mol. Cell. Biol.* **28 (8)**, 2803-2814.
2. Straub A. *et al.*, (2008). *Thromb. Haemost.* **99 (3)**, 609-615.

Description: Complex of N-terminal Flag-tagged recombinant full-length human p110β (GenBank Accession No. NM_006219), and untagged, recombinant full length, human p85α. (GenBank Accession No. NM_181523). Coexpressed in a Baculovirus infected Sf9 cell expression system. p110β MW=124 kDa, p85α MW=84 kDa.

Specific Activity: 1023 pmole/min/μg
 A 25 μl kinase reaction is conducted in a buffer containing 40 mM Tris-HCl (pH 7.5), 20 mM MgCl₂, 100 μM ATP, 200 μM 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

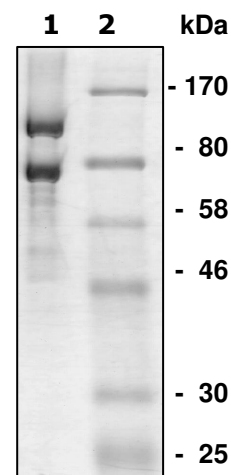


**10% SDS-PAGE
 Coomassie staining**

Lane 1:
 3.4 μg p110β/p85α →

Lane 2:
 Protein Marker →

MW:
 p110β: 124 kDa
 p85α: 84 kDa
Purity: ≥86%



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

Please visit our website at: www.bpsbioscience.com