

Data Sheet

PI3 p110δ/p85α

Human, recombinant

Catalog #: 40628

Lot #: 121008 **Conc.** 0.2 mg/ml

Formulated in: 25 mM Tris, pH 8.0, 69 mM NaCl, 1.35 mM KCl, 0.03% Tween-20, 3 mM DTT, and 50% glycerol.

Stability: >6 months at -80 °C

References:

1. Vanhaesebroeck *et al.* (1997) *Proc. Natl. Acad. Sci. U.S.A.* **94**:4330.
2. Vanhaesebroeck *et al.* (1999) *EMBO J.* **18**:1292.

Description: Complex of N-terminal GST recombinant full-length human p110δ (GenBank Accession No. NM_005026.2), and recombinant full length, human p85α (GenBank Accession No. XM_043865). Coexpressed in a Baculovirus infected Sf9 cell expression system. p110δ MW=146 kDa, p85α MW=84 kDa.

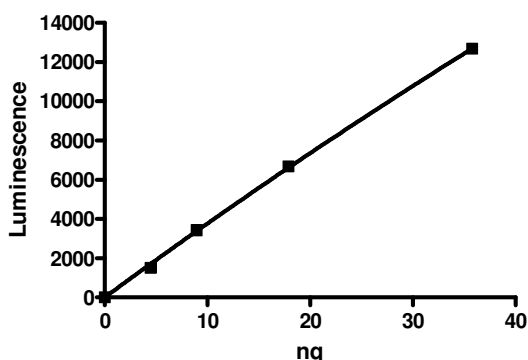
Specific Activity: 520 pmol/min/μg.

Assay conditions: 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 p110δ /p85α for 17 minutes 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

Specific Activity

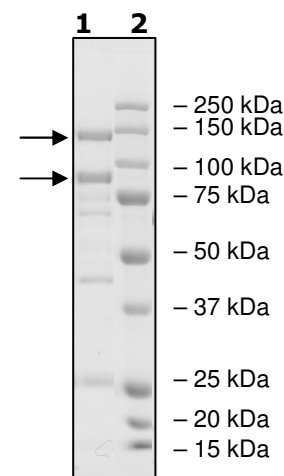


4-20% SDS-PAGE
Coomassie staining

Lane 1:
 1 μg p110δ /p85α

Lane 2:
 Protein Marker

MW:
 p110δ: 146 kDa
 p85α: 84kDa
Purity: ≥61%



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