

Data Sheet

p110α(E545K)/p85α

Human, recombinant, Complex

Catalog #: 40640

Lot #: 120409 **Conc.:** 0.56 mg/ml

Formulated in: 25 mM Tris-HCl, pH 8.0, 69 mM NaCl, 1.35 mM KCl, 0.025% Tween-20, 50% Glycerol and 3 mM DTT

Stability: At least 6 months at -80°C.
 Avoid freeze/thaw cycles.

References:

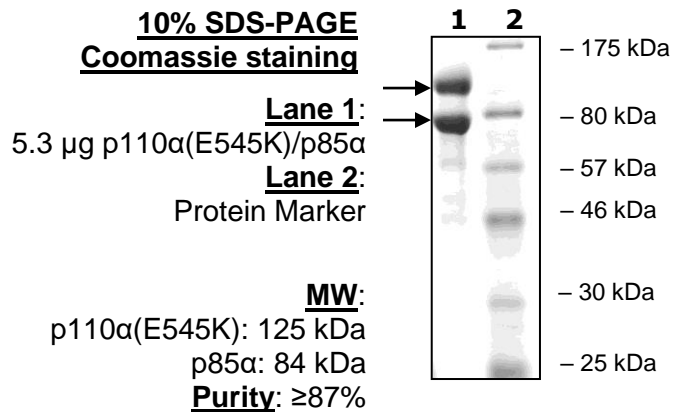
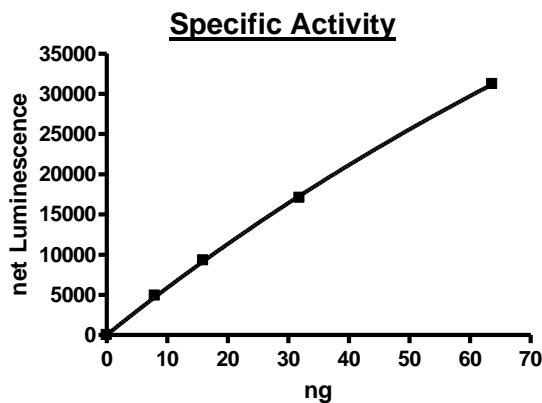
1. Engelman, J.A., *et al.* *Nature Med.* **14**: 1351-1356 (2008)
2. Hafner, C., *et al.*, *Proc. Natl. Acad. Sci. USA.* **104(33)**:13450-13454 (2007)
3. Serra, V., *et al.*, *Cancer Res.* **68(19)**:8022-8030 (2008)

Description: Complex of human p110α, also known as PIK3CA, GenBank Accession No. NM_006218, a.a. 2-1068(full length), with E545K mutation, with N-terminal FLAG-tag, MW=125 kDa, and human p85α, also known as PIK3R1, GenBank Accession No. NM_181523, a.a. 1-724(full length), MW=84 kDa. Coexpressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 1,200 pmol/min/μg.
 Assay Conditions: A unit of 110α(E545K)/p85α kinase activity is defined as the amount of enzyme required to produce 1 pmol of Phosphoinositol-3,4,5-Trisphosphate/min at 37°C. A 25 μl kinase reaction is conducted in 10 mM Tris-HCl (pH 7.5), 50 mM NaCl, 5 mM MgCl₂, 100 μM ATP, 10 μg mix of PS and Phosphoinositol-4,5-Bisphosphate, and 50 ng enzyme for 15 minutes at 30°C. Conversion of ATP to ADP is detected using PI3K ADP-Glo[®] Luminescent Kinase Assay (Promega Corp., Madison, WI).

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

Quality Assurance



OUR PRODUCTS ARE FOR RESEARCH ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
 To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com