

Data Sheet

p110α(K111N)/p85α

Human, recombinant, Complex

Catalog #: 40639

Lot #: 150716-1 **Conc.:** 0.13 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 0.04%
 Tween-20, 20% glycerol, 0.2 mM TCEP,
 80 µg/ml FLAG peptide

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

References:

- Rudd, M.L., *et al.*, *Clin. Canc. Res.* 2011; **17(6)**: 1331-1340.
- Brady, S.W., *et al.*, *Molec. Canc. Ther.* 2014; **13(1)**: 60-70.

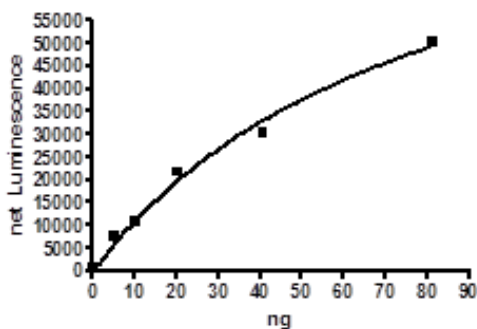
Description: Complex of human p110α, also known as PIK3CA, GenBank Accession No. NM_006218, a.a. 2-1068(full length), with K111N mutation, with N-terminal FLAG-tag, MW=124 kDa, and human p85α, also known as PIK3R1, GenBank Accession No. NM_181523, a.a. 1-724(full length), MW=84 kDa. Coexpressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 1326 pmol/min/µg.
 Assay Conditions: Reactions (25 µl) were done at 30°C for 15 min using BPS PI3K substrate, and the samples were incubated with the ADP-Glo reagent (25 µl) for 45 min followed by incubation with the Kinase Detection reagent (50 µl) for 55 min.

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

Quality Assurance

Specific Activity



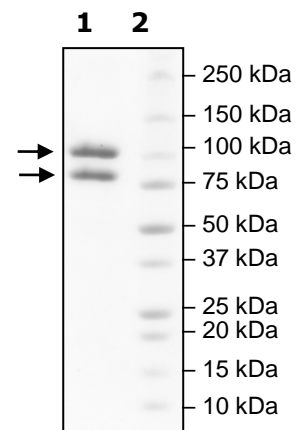
4-20% SDS-PAGE Coomassie staining

Lane 1:
4.2 µg P110a(K111N)/p85a

Lane 2:
Protein Marker

Purity: ≥95%

MW:
P110a(K111N):124 kDa
P85a:84 kDa



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