

Data Sheet

p110α(K111N)/p85α

Human, recombinant, FLAG-tag (p110α)
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. Storing diluted
 enzyme is not recommended, if
 necessary, use carrier protein (BSA 0.1 –
 0.5%).

References:

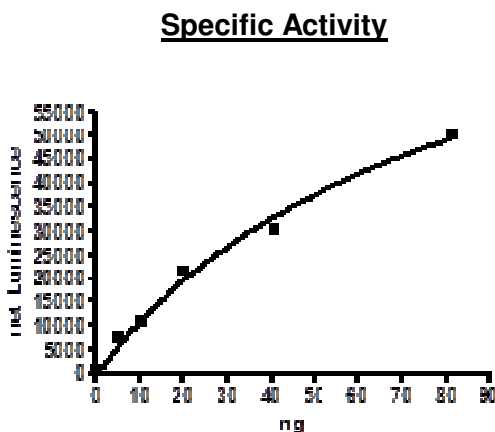
1. Rudd, M.L., *et al.*, *Clin. Canc. Res.* 2011; **17(6)**: 1331-1340.
2. Brady, S.W., *et al.*, *Molec. Canc. Ther.* 2014; **13(1)**: 60-70.

Description: Complex of N-terminal FLAG-tagged recombinant full-length human p110α (GenBank Accession No. U79143), with K111N mutation, and recombinant full length, human p85α (no tag, GenBank Accession No. XM_043865). Co-expressed in Sf9 cells in a Baculovirus expression system. p110α(K111N) MW=124 kDa, p85α MW=84 kDa.

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.

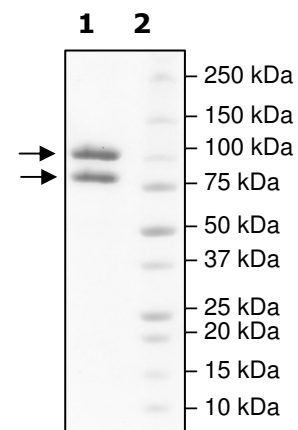
Quality Assurance



**4-20% SDS-PAGE
 Coomassie staining**

Lane 1: 4.2 µg P110α(K111N)/p85α
Lane 2: Protein Marker

Purity: ≥95%
MW:
 P110α(K111N):124 kDa
 P85α:84 kDa



OUR PRODUCTS ARE FOR RESEARCH ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
 To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com