

## Product Information

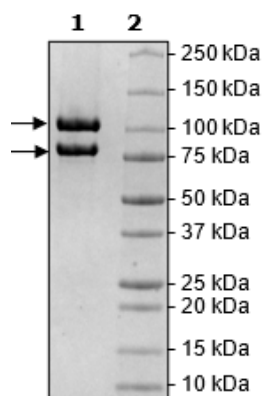
<b>Description:</b>	Recombinant human PI3 Kinase complex of p110 $\alpha$ , with an E542K mutation, and p85 $\alpha$ variant 1. Subunit p110 $\alpha$ encompasses amino acids 2-1068(end), and an N-terminal FLAG tag. Subunit p85 $\alpha$ encompasses amino acids 1-724(end). The two constructs were expressed together, and the complex was affinity purified.
<b>Background:</b>	PI3 (phosphoinositide 3) kinases, or phosphatidylinositol 3 kinases, are a family of proteins that can be subdivided into four classes: I, II, III and IV. Class I is involved in converting PI (4, 5) P2 (phosphatidylinositol (4, 5)-biphosphate) into PI (3, 4, 5) P3 (phosphatidylinositol (3, 4, 5)-triphosphate) when activated by tyrosine kinase receptors and G-protein coupled receptors. They are heterodimeric proteins with a regulatory and a catalytic subunit. The heterodimer between p110 (catalytic subunit) and p85 (regulatory subunit) belongs to class IA. P110 and p85 have three variants each. P110 $\alpha$ is ubiquitously expressed, and p85 $\alpha$ is the most abundant variant of p85. Class I PI3K participates in cell signaling, mostly via the activation of PKB (protein kinase B) and the PI3K/AKT/mTOR pathway. Dysfunction of these kinases impacts cell growth and differentiation, and mutations in p110 $\alpha$ have been linked to cancer. At least three isoform-specific inhibitors are approved by FDA for the treatment of lymphoma and leukemia. Further studies will help identify more selective inhibitors with a good tolerance that can bypass the development of drug resistance.
<b>Species:</b>	Human
<b>Construct:</b>	p110 $\alpha$ (E542K) (FLAG-2-1068(end)) / p85 $\alpha$ (1-724(end))
<b>Mutation:</b>	p110 $\alpha$ : E542K
<b>Concentration:</b>	1.54 mg/ml
<b>Expression System:</b>	Sf9
<b>Purity:</b>	$\geq 90\%$
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	25 mM Tris-HCl, pH 8.0, 69 mM NaCl, 1.3 mM KCl, 0.025% Tween-20, 50% glycerol, and 3 mM DTT
<b>MW:</b>	p110 $\alpha$ : 125 kDa; p85 $\alpha$ : 84 kDa
<b>Genbank Accession:</b>	p110 $\alpha$ : NM_006218; p85 $\alpha$ : NM_181523
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Assay Conditions:</b>	Assay was done according to Chemi-Verse™ PI3 Kinase p110 $\alpha$ (E542K)/p85 $\alpha$ Kinase Assay Kit (BPS Bioscience #82113) with various concentrations of PI3 Kinase p110 $\alpha$ (E542K)/p85 $\alpha$ .
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

# PI3 Kinase p110 $\alpha$ (E542K)/p85 $\alpha$ , FLAG-Tag Recombinant

Catalog: 101913  
Lot: 230808

## Quality Control Data

### 4-20% SDS-PAGE Coomassie Staining



### p110 $\alpha$ (E542K)/p85 $\alpha$ Activity

