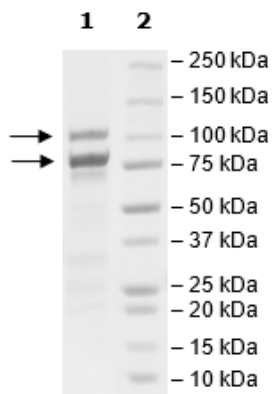


Product Information

Construct:	p110 α (H1047R) (FLAG-2-1068(end)) / p85 α (1-724(end))
Mutation:	p110 α : H1047R
Concentration:	1.80 mg/ml
Species:	Human
Formulated In:	25 mM Tris-HCl, pH 8.0, 69 mM NaCl, 1.35 mM KCl, 0.025% Tween-20, 3 mM DTT and 50% Glycerol
Expression System:	Coexpressed in Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	p110 α : NM_006218; p85 α : NM_181523
MW:	p110 α : 125 kDa; p85 α : 84 kDa
Purity:	80%
Specific Activity:	1,770 pmol/min/ μ g
Assay Conditions:	25 μ l kinase reaction is conducted in 40 mM Tris-HCl (pH 7.5), 10 mM MgCl ₂ , 100 μ M ATP, 200 μ M of PI3K substrate (BPS Bioscience #40560) and the PI3K enzyme for 18 minutes at 30°C. Conversion of ATP to ADP is detected using ADP-Glo [®] Luminescent Kinase Assay reagents (Promega Corp., Madison, WI).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

