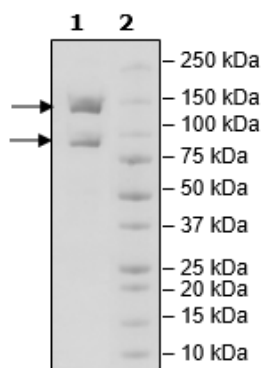


Product Information

Construct:	PI3 kinase (p110 α (GST-2-1068(end)) / p85 α (1-724(end)))
Concentration:	0.24 mg/ml
Species:	Human
Formulated In:	25 mM Tris-HCl, pH 8.0, 69 mM NaCl, 1.3 mM KCl, 0.03% Tween-20, 50% glycerol, and 3 mM DTT
Expression System:	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 : 151 kDa; p85 α : 84 kDa
Purity:	≥90%
Specific Activity:	880 pmol/min/μg
Assay Conditions:	Assay was done according to PI3K β (p110 β /p85 α) Assay Kit (BPS Bioscience #79802) with PI3K (BPS Bioscience #40621) titrated from 50 ng/rxn- 0 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining

PI3K (p110 α /p85 α) Activity