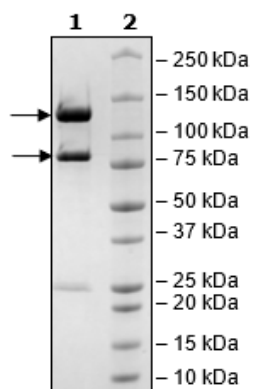


## Product Information

<b>Construct:</b>	PI3 Kinase p110δ (GST-1-1044(end)) / p85α (1-724(end))
<b>Concentration:</b>	0.90 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	25 mM Tris-HCl pH 8.0, 69 mM NaCl, 1.3 mM KCl, 0.03% Tween-20, 50% glycerol, and 3 mM DTT
<b>Expression System:</b>	Sf9
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	p110δ: NM_005026.2; p85α: NM_181523
<b>MW:</b>	p110δ: 145 kDa; p85α: 84 kDa
<b>Purity:</b>	≥90%
<b>Specific Activity:</b>	520 pmol/min/μg
<b>Assay Conditions:</b>	Assay was done according to PI3Kδ (p110δ/p85α) Assay Kit (96-well) (BPS Bioscience #79799) with various amounts of PI3 kinase (p110δ/p85α).
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-Page Coomassie Staining



PI3 Kinase (p110δ/p85α) Activity

