

Product Information

Construct:	PDK3 (GST-Th-2-406(end))
Concentration:	0.38 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 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:	NM_005391.5
MW:	73 kDa
Purity:	81%

Quality Control Data

4-20% SDS-Page Coomassie Staining

