

Product Information

Construct:	PDHK1 (GST-2-436(end))
Concentration:	0.13 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, 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:	NM_002610
MW:	76 kDa
Purity:	60%

Quality Control Data

4-20% SDS-Page Coomassie Staining

