

Product Information

Construct:	CDK13 (GST-516-1512(end)) / Cyclin K (GST-2-580(end))
Concentration:	0.59 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 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:	CDK13: NM_003718; Cyclin K: NM_001099402
MW:	CDK13: 138 kDa; Cyclin K: 90 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

