

Product Information

Construct:	CDK8 (GST-2-464(end)) / Cyclin C (GST-2-283(end))
Concentration:	0.12 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, 20% Glycerol.
Expression System:	Coexpressed in Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	CDK8: NM_001260; Cyclin C: NM_005190
MW:	CDK8: 79 kDa; Cyclin C: 59 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

