

Product Information

Construct:	NIK (GST-318-947(end))
Concentration:	0.17 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.4% Tween-20, 3 mM DTT, and 20% Glycerol.
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_003954
MW:	96 kDa
Purity:	≥75%

Quality Control Data

4-20% SDS-Page Coomassie Staining

