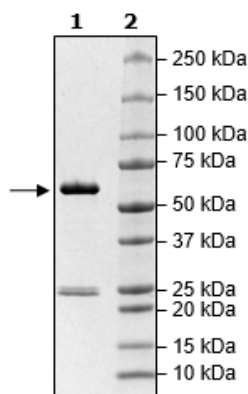


Product Information

Construct:	TNIK (GST-2-325)
Concentration:	0.49 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, and 2 mM glutathione
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_015028
MW:	63 kDa
Purity:	75%
Specific Activity:	301 pmol/min/μg
Assay Conditions:	Assay was done according to TNIK Kinase Assay Kit (BPS Bioscience #78370) with various amounts of TNIK.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



TNIK Activity

