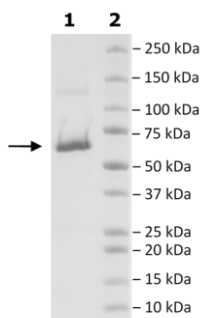


Product Information

Construct:	HPK1 (GST-Th-1-346-His-Avi)
Concentration:	0.57 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
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_007181
MW:	67 kDa
Purity:	≥90%
Specific Activity:	≥100 pmol/min/μg
Assay Conditions:	Assay was done according to the HPK1 Assay Kit (BPS Bioscience #79775) with HPK1 titrated from 120 ng/rxn – 1.875 ng/rxn. 20 μM ATP was used instead of 10 μM.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

