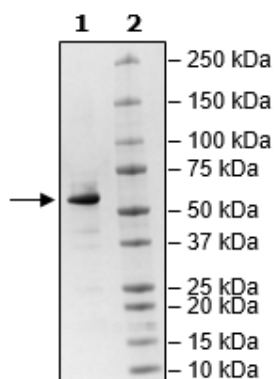


Product Information

Construct:	RIPK1 (GST-Th-1-327)
Concentration:	0.33 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 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_003804
MW:	64 kDa
Purity:	≥90%
Assay Conditions:	Assay was done according to RIPK1 Kinase Assay Kit (BPS Bioscience #79560) with RIPK1 (BPS Bioscience #40371) titrated at various concentrations.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



RIPK1 Activity

