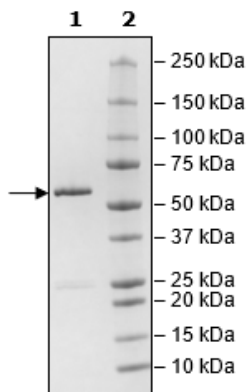


Product Information

Construct:	RIPK1 (GST-1-327) (Mouse)
Concentration:	0.38 mg/ml
Species:	Mouse
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
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_009068
MW:	64 kDa
Purity:	≥90%
Assay Conditions:	Assay was done according to the RIPK1 Kinase (Mouse) Assay Kit (BPS Bioscience #79897) with RIPK1 titrated from 400 ng/rxn to 0 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



RIPK1 (Mouse) Activity

