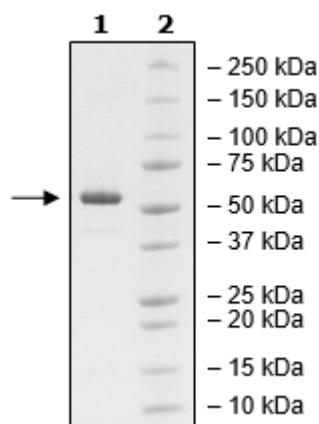


Product Information

Construct:	RIPK2 (GST-Pre-2-299)
Concentration:	0.29 mg/ml
Species:	Human
Formulated In:	45 mM Tris-HCl, pH 8.0, 500 mM NaCl, 2.5 mM KCl, 0.02% Tween-20, 3 mM DTT and 10% 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_003821
MW:	61 kDa
Purity:	≥90%
Specific Activity:	≥15 pmol/min/μg
Assay Conditions:	Assay was done in a Kinase buffer (BPS Bioscience #79334) using MBP (0.2 mg/ml) as a substrate with 20 μM ATP at 30°C for 45 minutes. The amount of ATP transferred was calculated using Kinase-Glo reagent (Promega).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Specific Activity

