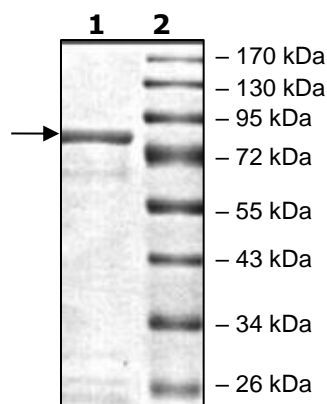


Product Information

Construct:	ROR1 (GST-429-end)
Concentration:	0.10 mg/ml
Species:	Human
Formulated In:	50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM glutathione, 0.1 mM EDTA, 0.25 mM DTT, 0.1 mM PMSF, and 25% 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_005012
MW:	82 kDa
Purity:	≥90%
Specific Activity:	0.2 pmol/min/μg
Assay Conditions:	ROR1 kinase activity was measured using substrate Myelin Basic Protein (MBP) diluted to 1 mg/ml in distilled water. ROR1 was diluted to 0.1 μg/ml in a buffer consisting of 5 mM MOPS, pH 7.2, 2.5 mM β-glycerol-phosphate, 5 mM MgCl ₂ , 1 mM EGTA, 0.4 mM EDTA, 50 μM fresh DTT and 50 ng/ml BSA (Bovine Serum Albumin). A serial dilution of ROR1 (10 μl of diluted kinase/tube) was incubated with 5 μl of 1 mg/ml stock MBP and 5 μl of distilled water. The "blank" had no substrate and was added water instead. The reaction was initiated by adding 5 μl of ATP cocktail: 250 μM [33P]-ATP (approximately 1 μCi/tube) diluted in 25mM MOPS, pH 7.2, 12.5mM β-glycerol-phosphate, 25 mM MgCl ₂ , 5 mM EGTA, 2 mM EDTA and 250 μM fresh DTT. The reaction was incubated for 15 min at 30°C and terminated by spotting 20 μl of the mixture onto phosphocellulose paper strips that were fixed in 1% phosphoric acid and washed three times. Radioactivity was determined using a scintillation counter.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Specific Activity

