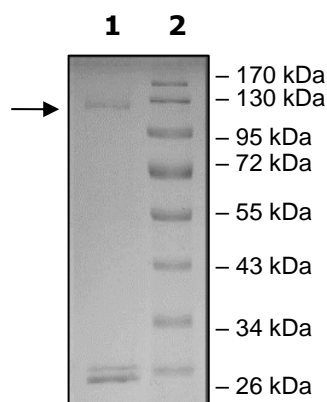


Product Information

Construct:	(GST-TPM3 (1-258)_TRKA (399-end))
Concentration:	0.05 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, 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:	TPM3: NM_152263; TRKA: NM_002529
MW:	105 kDa
Purity:	70%
Specific Activity:	13 pmol/min/μg
Assay Conditions:	<p>Kinase activity was measured using a Poly (Glu₄, Tyr₁) synthetic peptide substrate diluted in distilled water to a final concentration of 1 mg/ml.</p> <p>Increasing amounts of kinase were mixed with Poly (Glu₄,Tyr₁) synthetic peptide substrate with a final concentration of 200 μg/ml in a buffer containing 5 mM MOPS, pH 7.2, 2.5 mM β-glycerol-phosphate, 5 mM MgCl₂, 1 mM EGTA, 0.4 mM EDTA and 0.05 mM fresh DTT to a final volume of 20 uL. The reaction was initiated by addition of 5 μl of [33P]-ATP diluted in kinase buffer: 6 ml kinase buffer containing 1 mCi [33P]-ATP, 0.25 mM ATP, 25 mM MOPS, pH 7.2, 12.5 mM β-glycero-phosphate, 25 mM MgCl₂, 5 mM EGTA, 2 mM EDTA, and 0.25 mM fresh DTT.</p> <p>After incubating for 30°C for 15 minutes, the reaction was 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. The blank was determined from a “no substrate” sample.</p>
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

