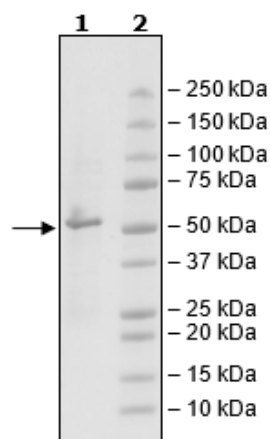


Product Information

Description:	Recombinant human SYK (spleen associated tyrosine kinase), encompassing amino acids 359-635(end). This construct contains an N-terminal GST-tag. A short peptide linker placed between GST and the protein includes a Thrombin Cleavage Site to facilitate removal of the GST portion. The recombinant protein was affinity purified.
Species	Human
Construct:	SYK Catalytic (GST-359-635(end))
Concentration:	0.13 mg/ml
Expression System:	Sf9
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 3 mM DTT
MW:	59 kDa
Genbank Accession:	NM_003177
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Specific Activity:	≥1000 pmole/min/μg
Assay Conditions:	Assay was done in Kinase Buffer 1 (BPS Bioscience #79334) using, 0.2 mg/ml Poly(Glu:Tyr) as a substrate with 20 μM ATP incubated at 30°C for 35 min. SYK, Catalytic, GST-Tag Recombinant (BPS Bioscience #40487) was titrated at various concentrations. The amount of ATP transferred was calculated using Kinase Glo max reagent (Promega).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



SYK Activity

