

Data Sheet

SYK, Full Length, GST-Tag

Human, recombinant, full length
 N-terminal GST tag
Catalog #: 40485
Lot# 100310 **Conc.:** 1 mg/ml

Fomulated in: 45mM Tris-HCl, pH 8,
 124mM NaCl, 2.4mM KCl, 18mM
 Glutathione, 3 mM DTT and 10% glycerol.

Stability: At least 6 months at -80°C. *Avoid freeze/thaw cycles.*

References:

1. Takano, T. et al. Antioxid Redox Signal. 2002 Jun;**4(3)**:533-41.
2. Navara, C S. Curr Pharm Des. 2004;**10(15)**:1739-44.

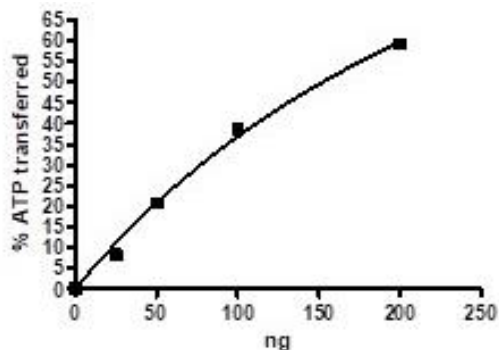
Description: Recombinant full-length human SYK also known as Spleen Associated Tyrosine Kinase and Tyrosine-Protein Kinase SYK, GenBank Accession No. NM_003177, a.a. full length with N-terminal GST-tag, expressed in a Sf9 expression system. MW = 100 kDa.

Specific Activity: 108 pmol/min/μg
 Assay conditions: 40 mM Tris-HCl pH 7.4, 20 mM MgCl₂, 0.1 mg/mL BSA and 0.2 mM DTT using 0.2 mg/ml Poly(Glu:Tyr) substrate and 20 μM ATP. Reaction was done at 30°C for 35 min. Amount of ATP transferred was calculated using Kinase-Glo reagent from Promega.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

Specific Activity



4-20 % SDS-PAGE
Coomassie staining

