

Data Sheet

SYK, Catalytic Subunit, GST-Tag

Human, recombinant, N-terminal GST-tag
Catalog #: 40487
Lot #: 120806-G **Conc.:** 0.21 mg/ml

Formulated in: 40 mM Tris-HCl pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, and
 20% glycerol

Stability: >6 months at -80°C. *Avoid
 freeze/thaw cycle.*

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.

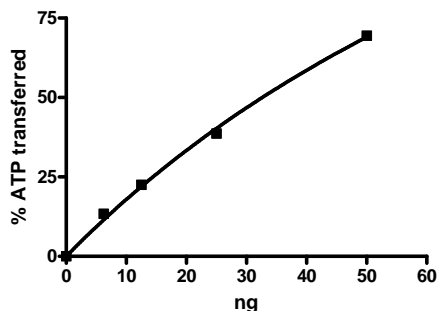
Description: Recombinant human SYK Catalytic Subunit, also known as SYK, and Spleen Tyrosine Kinase, GenBank Accession No. NM_003177, a.a. 359-635(end) with a N-terminal GST-tag, expressed in Sf9 insect cells using. MW=58 kDa.

Specific Activity: 451 pmol/min/μg
 Assay conditions: 40 mM Tris-HCl pH 7.4, 20 mM MgCl₂, 0.1 mg/mL BSA and 1 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



10% SDS-PAGE
Coomassie staining

Lane 1:
 1.8 μg SYK
Lane 2:
 Protein Marker
MW: 58 kDa
Purity: ≥90%

