

Data Sheet

LCK, GST-Tag

Human, recombinant, N-terminal GST tag

Catalog #: 40470

Lot #: 170801-B2

Conc.: 0.53 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 20% glycerol mM DTT, 1.6 mM glutathione.

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

References:

1. Fischer, S. *et al. Biochem Biophys Res Commun.* 1987 Mar 30;**143(3)**:819-26.
2. Veillette, A. *et al. Oncogene Res.* 1987 Sep-Oct;**1(4)**:357-74.

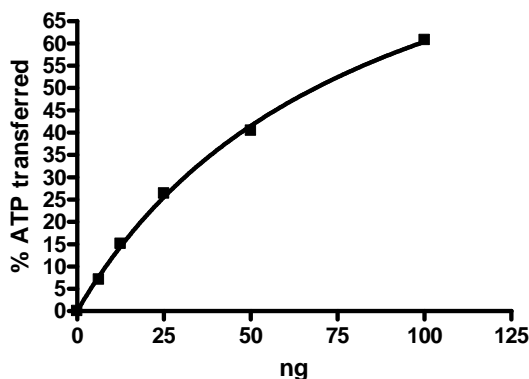
Description: Human LCK, also known as Leukocyte C-terminal Src kinase, Lymphocyte cell-specific protein-tyrosine kinase, IMD22, and LSK, GenBank Accession No. NM_005356, a.a. 2-509(end), with N-terminal GST-tag, expressed in Sf9 insect cells using a baculovirus expression system. MW=85 kDa.

Specific Activity: ≥200 pmol/min/μg.
 Assay Conditions: Assay was done in Kinase buffer containing 5 mM MnCl₂, 1 mM DTT using Poly-(Glu:Tyr) substrate (0.2 mg/ml) and 20 μM ATP. Reaction was done at 30°C for 45 min. Amount of ATP transferred was calculated using Kinase-Glo reagent (Promega).

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

Quality Assurance

Specific Activity



**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
LCK(2-e)

Lane 2:
Protein Marker

Purity: ≥90%

MW: 85 kDa

