

Data Sheet

SRC

Human, recombinant, N-terminal HIS tag

Catalog #: 40484

Lot #: 130909 **Conc.:** 1.2 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 106 mM imidazole and 20% glycerol.

Stability: >6 months at -80°C . *Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

References:

1. Finlay, D. and Vuori, K.. Cancer Res. **67(24)**: 11704-11711 (2007).
2. Vielreicher, M. et al. J. Biol. Chem. **282(46)**: 33623-33631 (2007).

Description:

Human SRC (GenBank Accession No. NM_005417), full length with N-terminal His-tag, MW = 61 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

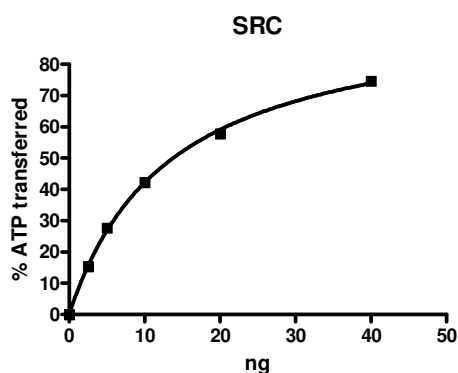
Specific Activity: 1007 pmol/min/ μg
 Assay conditions: Kinase buffer containing 1 mM DTT using Poly-(Glu4:Tyr) substrate (0.2 mg/ml) and 20 mM ATP. 35 min at room temperature.

Application:

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

Quality Assurance

SRC Specific Activity



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

Lane 1:
3.5 μg His-Src
Lane 2:
Protein Marker

Purity: $\geq 85\%$
MW: 61 kDa

