

Data Sheet

ITK(Lyk), Rat

Rat recombinant, with N-terminal GST-tag

Catalog #: 40447

Lot #: 100623

Conc.: 1 mg/ml

Formulated in: 25 mM Tris-HCl, pH 8.0,
 100 mM NaCl, 0.05% Tween-20, 3 mM
 DTT, and 10% glycerol.

Stability: >6 months at -80°C

References:

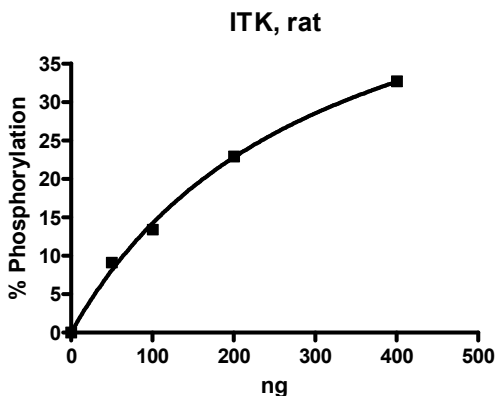
1. Hawkins, J. and Marcy, A. (2001). *Protein Expr. Purif.* **22 (2)**: 211-9.
2. Kosaka, Y., et al. (2006). *Trends Immunol.* **27 (10)**: 453-60
3. Tanaka, N., et al. (1993). *FEBS Lett.* **324 (1)**: 1-5.

Description: Rat recombinant IL2-inducible T-cell kinase (ITK), also known as Lyk and TSK (GenBank Accession #NM_001108825), a.a. 352-620(end) with N-terminal GST-tag, MW= 57 kDa, expressed in Sf9 cells via a baculovirus expression system.

Specific activity: 2.1 pmole/min/μg
 Assay conditions: Z-Lyte® (Invitrogen) screening protocol and condition were used. Enzyme reaction as conducted at 30°C for 1 hours with 2 μM Tyr 01 Peptide in 50 mM HEPES pH 7.5, 0.01% BRIJ-35, 10 mM MgCl₂, 1 mM EGTA. After the 1 hour incubation, 1:128 dilution of Development Reagent B is added.

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

Quality Assurance

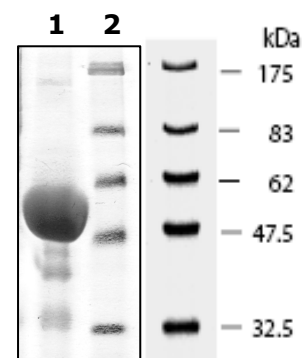


% Phosphorylation was calculated based on the emission ratio of 445 nm/520 nm after excitation

10% SDS-PAGE Coomassie staining

Lane 1:
20 μg ITK
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: ITK: 57 kDa
Purity: ≥58%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com