

## Data Sheet

### ITK(Lyk), Murine

Murine recombinant, with N-terminal GST-tag

**Catalog #:** 40446

**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:**

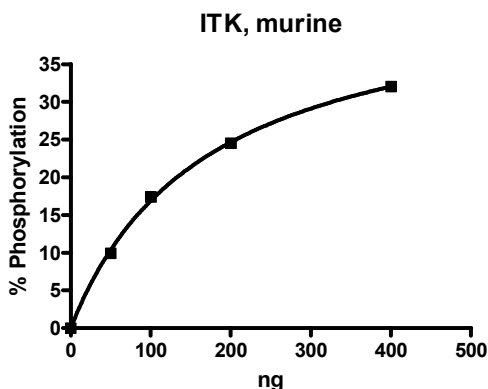
1. Hawkins, J. and Marcy, A. (2001). *Protein Expr. Purif.* **22 (2)**: 211-9.
2. Kosaka, Y., et al. (2006). *Trends Immunol.* **27 (10)**: 453-60
3. Tanaka, N., et al. (1993). *FEBS Lett.* **324 (1)**: 1-5.

**Description:** Murine recombinant IL2-inducible T-cell kinase (ITK), also known as Lyk and TSK (GenBank Accession #NM\_010583), a.a. 352-620(end) with N-terminal GST-tag, MW= 57 kDa, expressed in Sf9 cells via a baculovirus expression system.

**Specific activity:** 2.4 pmole/min/μg  
 Assay conditions: Z-Lyte<sup>®</sup> (Invitrogen) screening protocol and condition were used. Enzyme reaction was 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 MgCl<sub>2</sub>, 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

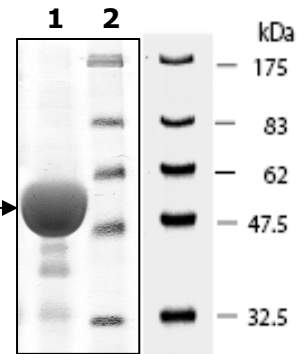


% Phosphorylation was calculated based on the emission ratio of 445 nm/520 nm after excitation at 400 nm.

**10% SDS-PAGE  
 Coomassie staining**

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

**MW:** ITK: 57 kDa  
**Purity:** ≥62%



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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)