

## Data Sheet

### ITK

Human, recombinant, N-terminal GST tag

**Catalog #:** 40445

**Lot#:** 80610

**Conc.:** 1.02 mg/ml

**Formulated in:** 25 mM Tris-HCl, pH 8.0,  
100 mM NaCl, 1.37 mM KCl, 0.05% Tween-  
20, 30% glycerol, 10 mM glutathione, and 3  
mM DTT.

**Stability:** >6 months at -80°C

#### References:

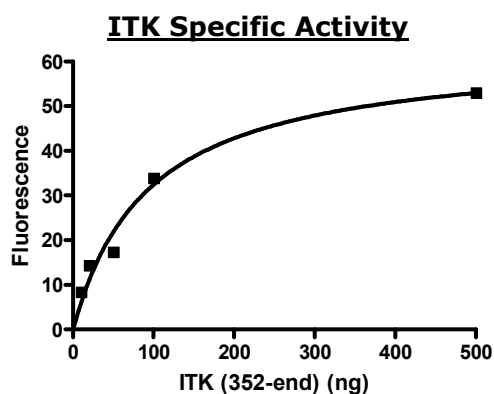
1. Joseph R.E., *et al.*, *J. Mol. Biol.* **373** (5), 1281-1292 (2007).
2. Hu J. *et al.*, *Eur. J. Immunol.* **37** (10), 2892-2899 (2007).

**Description:** Human ITK (GenBank Accession No. NM\_005546), a-a 352 to end with N-terminal GST tag, MW= 57 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

**Specific Activity:** 13 pmol/min/μg  
Assay condition: 50 mM HEPES (pH7.5), 10 mM MgCl<sub>2</sub>, 1 mM EGTA, 0.01% Brij-35, 200 μM ATP, 2 μM Tyr1 substrate (Z-lyte kinase assay kit from Invitrogen), 1 hour at room temperature.

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

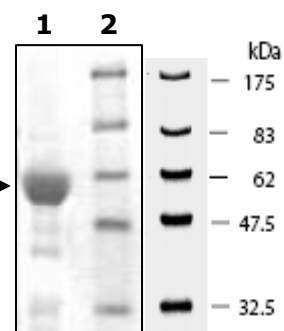
## Quality Assurance



### 10% SDS-PAGE Coomassie staining

**Lane 1:**  
6 μg ITK →

**Lane 2:**  
Protein Marker  
BioLabs (#P7708L)



**MW:** 57 kDa.  
**Purity:** >80%

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)