

## Data Sheet

### TBK1, GST-Tag

Human, Recombinant, N-terminal GST-Tag

**Catalog #:** 40286

**Lot#:** 170614-G      **Conc:** 0.12 mg/mL

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
 110 mM NaCl, 2.2 mM KCl, 0.04%  
 Tween20, 8.8mM glutathione, 20% glycerol

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid  
 freeze/thaw cycles. Storing diluted enzyme  
 is not recommended, if necessary, use  
 carrier protein (BSA 0.1 – 0.5%).

**References:**

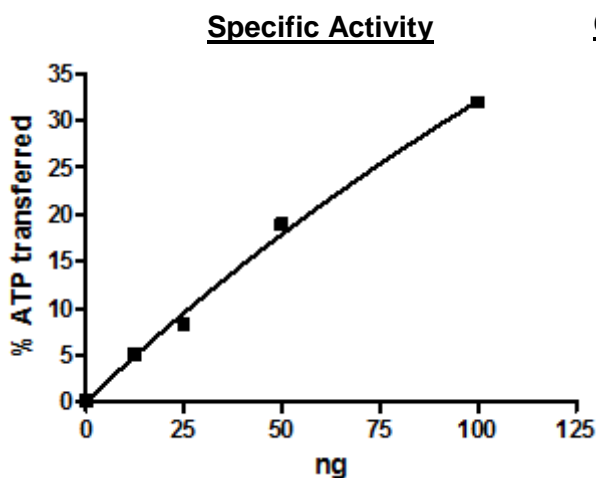
1. Tojima, Y. *et al. Nature* **404**: 778-782, 2000.
2. Sharma, S. *et al. Science* **300**: 1148-1151, 2003.

**Description:** Recombinant human TBK1, also known as NF-kappa-B-activating kinase, T2K, and TANK-binding kinase 1, with N-terminal GST-tag, 2-729 (end), Genbank accession number NM\_013254, expressed in Sf9 insect cells via a baculovirus expression system. MW=110 kDa.

**Specific Activity:** 80 pmol/min/ $\mu\text{g}$   
 Assay Conditions: Assay was done in Kinase buffer containing 1 mM DTT using MBP as a substrate (0.2 mg/ml) and 20  $\mu\text{M}$  ATP. Reaction was done at 30 $^{\circ}\text{C}$  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



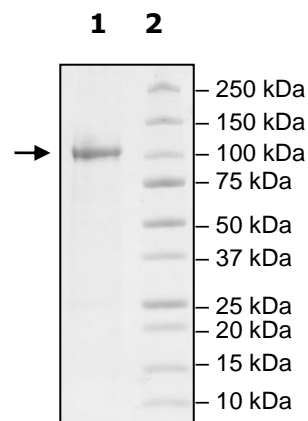
**4-20% SDS-PAGE  
 Coomassie staining**

**Lane 1:**  
 TBK1(2-e)

**Lane 2:**  
 Protein Marker

**Purity:**  $\geq 90\%$

**MW:** 110 kDa



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)