

## Data Sheet

### TTBK1, GST-Tag, His-Tag

Human, Recombinant, N-terminal GST-Tag,  
C-terminal His-Tag  
**Catalog #:** 40309  
**Lot #:** 170119-HG                      **Conc:** 0.24 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 20% glycerol, 2 mM glutathione

**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. Sato, S., *et al.*, *Journal of Neurochemistry*. 2006; **98(5)**: 1573-1584.
2. Yu, N., *et al.*, *Neuroscience Letters*. 2011; **491(1)**: 83-86.

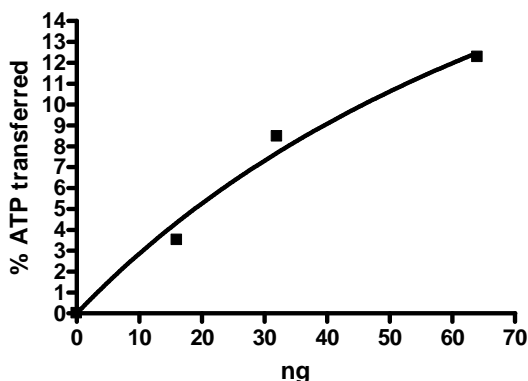
**Description:** Human TTBK1 protein, also known as Tau-tubulin kinase-1, GenBank Accession No. NM\_032538, a.a. 1-479, with N-terminal GST-tag and C-terminal His-tag. MW = 81 kDa, expressed in Sf9 cells via a baculovirus expression system.

**Assay Conditions:** Assay was done in a Kinase buffer using MBP (0.1 mg/ml) as a substrate with 20  $\mu\text{M}$  ATP at  $30^{\circ}\text{C}$  for 45 min. The 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

### Specific Activity

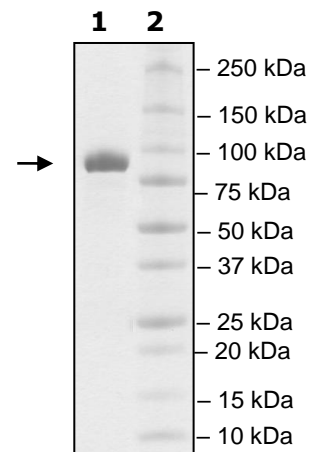


### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
TTBK, GST-Tag,  
His-Tag

**Lane 2:**  
Protein Marker

**MW:** 81 kDa  
**Purity:**  $\geq 90\%$



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)