

Data Sheet

TTK

Human, recombinant, N-terminal GST-tag

Catalog #: 40291

Lot #: 140701

Conc.: 0.1 mg/ml

Formulated in: 50 mM Tris-HCl, 7.5, 150 mM NaCl, 10 mM glutathione, 0.25 mM, 0.1 mM PMSF, 25% glycerol.

Stability: >6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. Fisk, H.A., *et al. Proc. Nat. Acad. Sci.* 2003; **100**:14875-14880
2. Huang, Y.F., *et al. Mol. Cell Biol.* 2009; **11**:2935-2944.

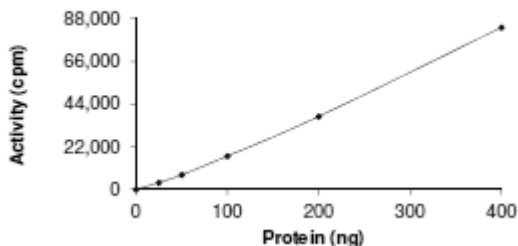
Description: Human Dual specificity protein kinase TTK, also known as MPS1, MPS1L1 GenBank Accession No. NM_003318, amino-acids 1-857(end) with an N-terminal GST-tag, MW=130 kDa, expressed in a baculovirus-infected Sf9 cell expression system.

Specific Activity: 18 pmole/min/ μg

Assay Conditions: Varying concentrations of MLK1 were mixed with 5 μg of MBP protein substrate and 5 μl of 250 μM [^{33}P]-ATP was added for a total volume of 25 μl . Mixture was incubated for 15 min at 30°C . Reaction was stopped by placing 20 μl of reaction onto phosphocellulose P81 paper. Paper was washed 3 times with 1% phosphoric acid solution for 10 min each and air dried. Radioactivity was measured from the P81 paper in the presence of scintillation fluid using a scintillation counter.

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

Quality Assurance

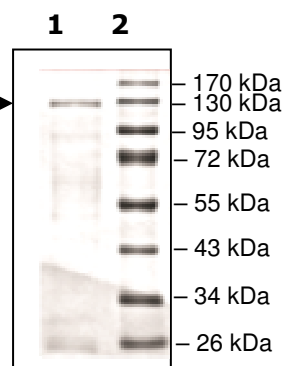


4-20% SDS-PAGE Coomassie staining

Lane 1:
TTK

Lane 2:
Protein Marker

MW: 130 kDa
Purity: $\geq 75\%$



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