

Data Sheet

MLK3

Human, recombinant, N-terminal GST-tag

Catalog #: 41104

Lot #: xxxxxx

Conc.: 0.1 mg/ml

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

Stability: >6 months at -80°C. Avoid freeze/thaw cycles.

References:

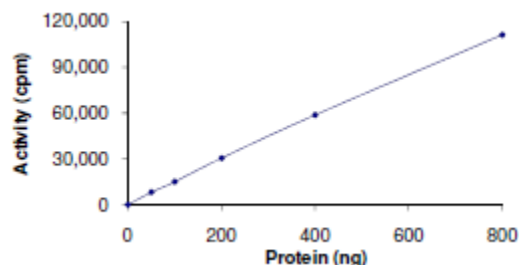
1. Ing, Y.L. *et al.*, *Oncogene*. 1994. **9**:1745-1750.
2. Chadee, D.N. *et al.*, *Nature Cell Biol.* 2004. **6**:770-776.

Description: Human MLK3, also known as Mixed Lineage Kinase 3, Mitogen-activated protein kinase kinase kinase 11, MAP3K11, PTK1, SPRK, and MGC17114, GenBank Accession No. NM_002419, a.a. 1-488, with N-terminal GST-tag, MW = 83 kDa, expressed in Sf9 cells via a baculovirus expression system.

Specific Activity: 9.5 pmole/min/μg
Assay Conditions: Varying concentrations of MLK3 were mixed with 5 μg of MBP protein substrate and 5 μl of 250 μM [³³P]-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.

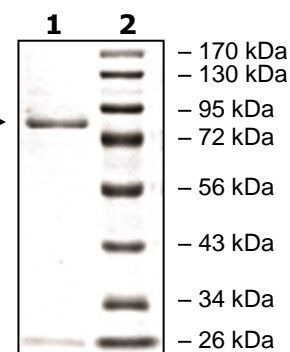
Quality Assurance



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

Lane 1: MLK3
Lane 2: Protein Marker

MW: 83 kDa
Purity: ≥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