

## Data Sheet

### MLK1

Human, recombinant, N-terminal GST-tag

**Catalog #:** 41102

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  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

#### References:

1. Dorrow, D.S., *et al.*, *Europ. J. Biochem.* 1993. **213**:701-710.
2. Durkin, J.T., *et al.*, *Biochemistry.* 2004. **43(51)**:16348-16355.

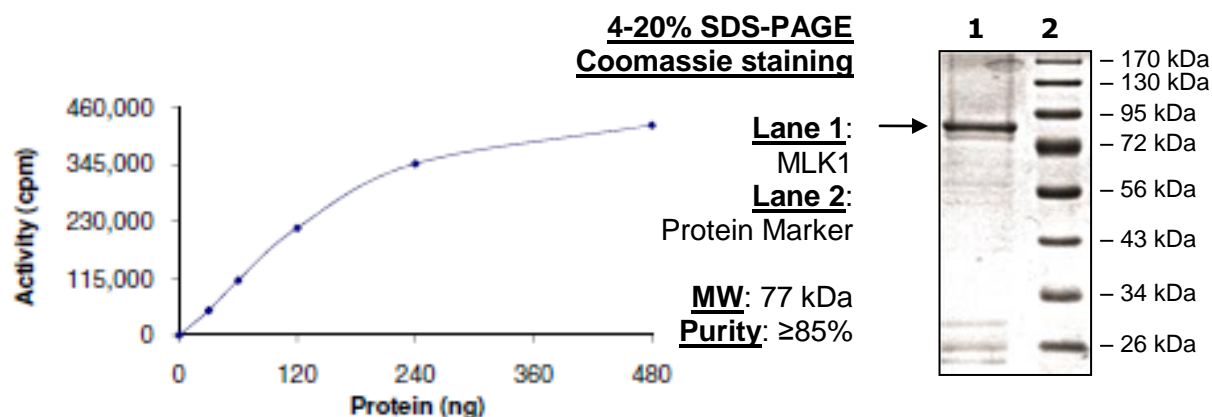
**Description:** Human MLK1, also known as Mixed Lineage Kinase 1, Mitogen-activated protein kinase kinase kinase 9, MAP3K9, MEKK9, and PRKE1, GenBank Accession No. NM\_033141, a.a. 1-433, with N-terminal GST-tag, MW = 77 kDa, expressed in Sf9 cells via a baculovirus expression system.

**Specific Activity:** 116 pmole/min/ $\mu\text{g}$

**Assay Conditions:** Varying concentrations of MLK1 were mixed with 5  $\mu\text{g}$  of MBP protein substrate and 5  $\mu\text{l}$  of 250  $\mu\text{M}$  [ $^{33}\text{P}$ ]-ATP was added for a total volume of 25  $\mu\text{l}$ . Mixture was incubated for 15 min at  $30^{\circ}\text{C}$ . Reaction was stopped by placing 20  $\mu\text{l}$  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



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)