

## Data Sheet

### MLK4

Human, recombinant, N-terminal GST-tag

**Catalog #:** 41105

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:**

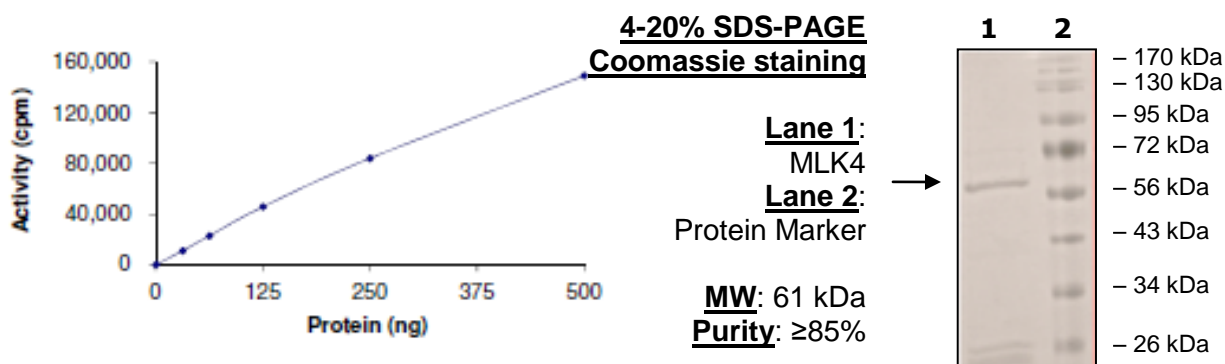
1. Nagase, T., *et al.*, *DNA Res.* 2001. **8**:85-95.
2. Seit-Nebi, A., *et al.*, *Cell. Molec. Immun.* 2012. **9**:27-33.

**Description:** Human MLK4, also known as Mixed Lineage Kinase 4, and KIAA1804, GenBank Accession No. NM\_032435, a.a. 1-557, with N-terminal GST-tag, MW = 61 kDa, expressed in Sf9 cells via a baculovirus expression system.

**Specific Activity:** 21 pmole/min/μg  
 Assay Conditions: Varying concentrations of MLK4 were mixed with 5 μg of MBP protein substrate and 5 μl of 250 μM [<sup>33</sup>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



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact: 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)