

## Data Sheet

### MST4

Human, recombinant, N-terminal GST tag

**Catalog #:** 40134

**Lot #:** 110419      **Conc.:** 0.1 mg/ml

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

**Stability:** >12 months at -80°C

**References:**

1. Dan, I., *et al. J. Biol. Chem.* **277**: 5929-5939 (2002).
2. Sung, V., *et al. Cancer Res.* **63**: 3356-3363 (2003).

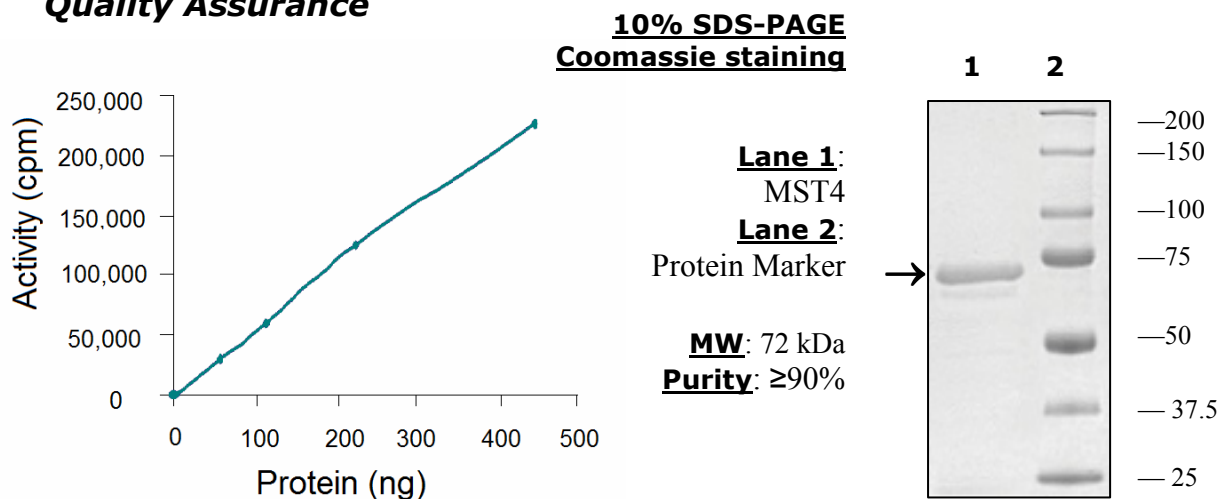
**Description:** Human MST4, also known as MASK (GenBank Accession No. NM\_016542), full length, with N-terminal GST tag, MW ~72 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

**Specific Activity:** 23 pmol/min/μg

Assay Buffer: 5 mM MOPS, 2.5 mM β-glycerophosphate, 5 mM MgCl<sub>2</sub>, 1 mM EGTA, 0.4 mM EDTA, 0.05 mM DTT, 2 mM ATP. Incubate MST4 with 0.1 mg/ml Myelin basic protein (substrate) and [<sup>33</sup>P]-ATP at 30°C for 15 minutes, then spot reaction on phosphocellulose paper, fix in 1% phosphoric acid, and assay with 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)