

## Data Sheet

### MAPKAPK2

Human recombinant, with N-terminal His-tag

**Catalog #:** 40116

**Lot #:** 111111      **Conc.:** 0.1 mg/ml

**Formulated in:** 50 mM sodium phosphate, pH 7.5, 300 mM NaCl, 150 mM imidazole, 0.1 mM PMSF, 0.25 mM DTT, 25% glycerol

**Stability:** >6 months at -80°C

**References:**

1. Hagood JS, Olman MA. *Am. J. Respir. Cell Mol. Biol.* **37(5)**:503-6 (2007).
2. Kotlyarov A, Gaestel M. *Biochem. Soc. Trans.* **30(Pt 6)**:959-63 (2002).

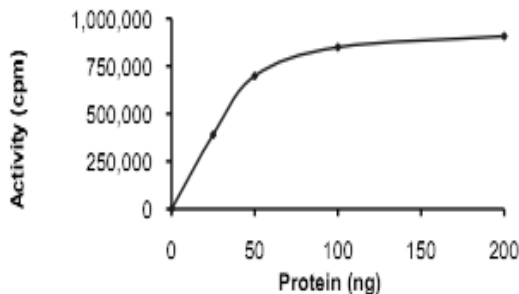
**Description:** Recombinant human MAPKAPK2 (MAPK-activated protein kinase 2), also known as MK2, Genbank Accession #NM\_032960, a.a. 46-end, with an N-terminal His-tag, MW = 41 kDa, expressed in Sf9 cells via a baculovirus expression system.

**Specific Activity:** 450 pmol/min/μg.  
**Assay conditions:** 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, 50 μg/ml BSA, 2 mM ATP. Incubate MEK1 with 0.2 mg/ml HSP27tide peptide as 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 studying enzyme kinetics, screening inhibitors, and selectivity profiling.

### Quality Assurance

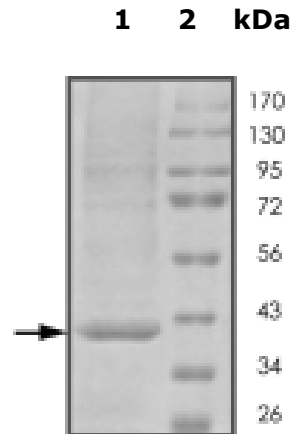
**Specific Activity**



**10% SDS-PAGE**  
**Coomassie staining**

**Lane 1:** MAPKAPK2  
**Lane 2:** Protein Marker

**MW:** 41 kDa  
**Purity:** ≥85%



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)