

Data Sheet

MEKK2

Human, recombinant, N-terminal GST-tag

Catalog #: 40122

Lot#: 120509-G12 **Conc.:** 0.062 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol and 3 mM DTT.

Stability: >6 months at -80 °C

References:

1. Blank, J. L. *et al. J. Biol. Chem.* **271:** 5361-5368, 1996.
2. Hammaker, D. R. *et al. J. Immun.* **172:** 1612-1618, 2004.

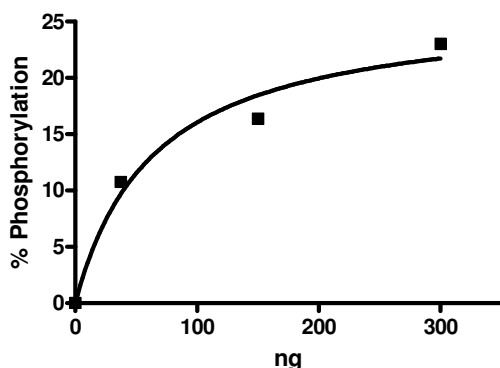
Description: Human MEKK2, also known as Mitogen-activated protein kinase kinase 2 (MAP3K2) and MEKK2B, GenBank Accession No. NM_006609, full length, N-terminal GST-tag, MW=96.5 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

Specific Activity: 18 pmol/min/μg
 Assay conditions: Enzyme reaction is conducted in a Kinase Buffer (50 mM Tris pH 7.4, 20 mM MgCl₂, 0.1 mg/ml BSA and 1 mM DTT) containing 20 μM ATP and 0.2 mg/ml poly-(Glu-Tyr) for 30 min at 30 °C. Enzyme activity is measured using Kinase-Glo Plus reagents (Promega Corporation, Madison, WI).

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

MEKK2 Specific Activity



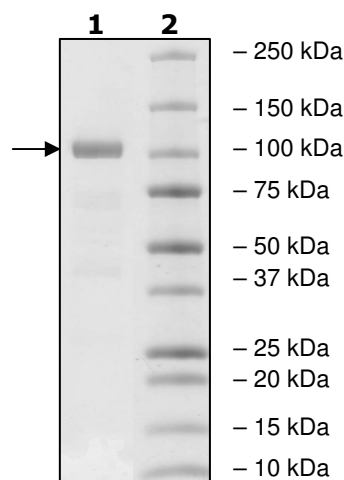
4-20% SDS-PAGE Coomassie staining

Lane 1:
0.7 μg MEKK2

Lane 2:
Protein Marker
Biorad (#161-0373)

MW: 96.5 kDa

Purity: ≥87%



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