

Data Sheet

MAP4K3 (GLK), GST-tag

Human, Recombinant, N-terminal GST-tag
Catalog #: 100624
Lot #: 200130 **Conc.:** 0.1 mg/ml

Formulated in: 50 mM Tris-HCl, pH 7.5,
150 mM NaCl, 10 mM glutathione, 0.25 mM
DTT, 0.1 mM EDTA, 0.1 mM PMSF, 25%
glycerol.

Stability: At least 6 months at -80°C . Avoid
freeze/thaw cycles.

References:

1. Lijun, Y. *et al*, *Molec. Cell.* **37(5)**:
633-642, 2010.
2. Zhao, B. *et al*, *Cancer Letters.*
342(1): 43-51, 2014

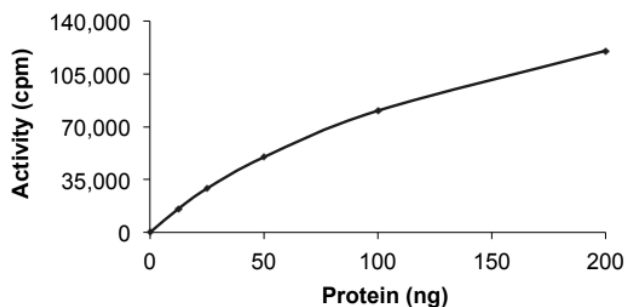
Description: Human MAP4K3, also
known as GLK, MAPKKK3, MEKKK3
and RAB8IPL1, GenBank Accession No.
NM_003618, a.a. 1-380, with N-terminal
GST-tag, expressed in a Baculovirus
infected Sf9 cell expression system.
MW= 70 kDa.

Specific Activity: ≥ 78 nmol/min/mg

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

Quality Assurance

Specific Activity

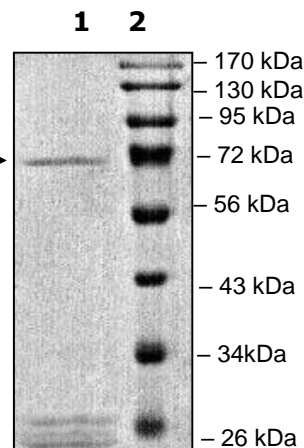


4-20% SDS-PAGE Coomassie staining

Lane 1:
MAP4K3(GLK)
Lane 2: →
Protein Marker

Purity: $\geq 70\%$

MW: 70 kDa



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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com