

Data Sheet

MAPKAPK2, GST-tag

Human, Recombinant, N-terminal GST-tag

Catalog #: 40088

Lot#: 130531-G2

Conc.: 0.44 mg/ml

Formulated in: 40 mM Tris pH 8, 110 mM NaCl, 2.2mM KCl, 20% glycerol, 3mM DTT.

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

References:

1. Ter Haar, E. et al., J. Biol. Chem. 282 (13), 9733-9739 (2007).
2. Hillig, RC. et al., J Mol Biol. 369(3):735-45 (2007).

Description:

Human MAPKAPK2 also known as MK2, MAPK Activated Protein Kinase 2, and MAPKAP Kinase 2, GenBank Accession No. NM_032960, a.a 1-400 (full length) with N-terminal GST-tag, expressed in a Baculovirus infected Sf9 cell expression system. MW=72 kDa.

Quality Assurance

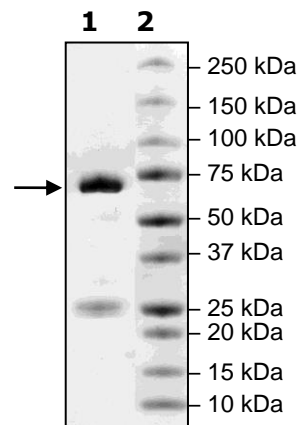
4-20% SDS-PAGE Coomassie staining

Lane 1:
2.5 μg MAPKAPK2

Lane 2:
Protein Marker

MW: 72 kDa

Purity: $\geq 74\%$



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