

Data Sheet

MNK2

Human, recombinant, N-terminal GST tag

Catalog #: 40128

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. Waskiewicz, A.J., *et al. EMBO J.* **277:** 5929-5939 (1997).
2. Slentz-Kesler, K., *et al. Genomics.* **69:** 63-71 (2000).

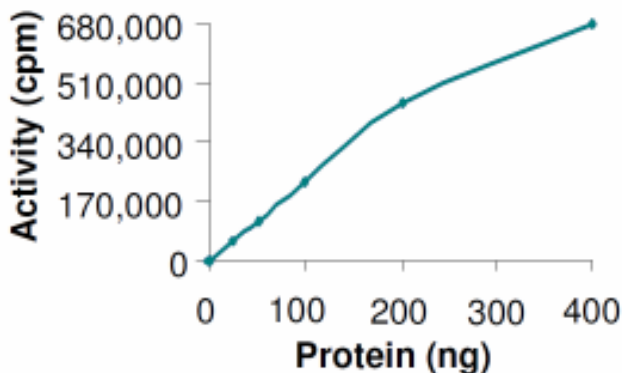
Description: Human MNK2 (MAP kinase interacting serine/threonine kinase 2), also known as GPRK7 (G protein-coupled receptor kinase 7), GenBank Accession No. NM_017572, full length, with N-terminal GST tag, MW ~74 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

Specific Activity: 93 pmol/min/μg

Assay Buffer: 5 mM MOPS, 2.5 mM β-glycerophosphate, 5 mM MgCl₂, 1 mM EGTA, 0.4 mM EDTA, 0.05 mM DTT, 2 mM ATP. Incubate MNK2 with 0.25 mg/ml Myelin basic Protein (substrate) and [³³P]-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.

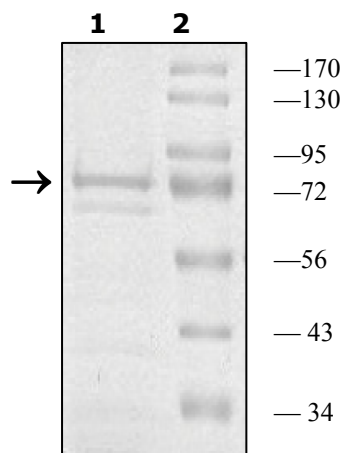
Quality Assurance



**10% SDS-PAGE
 Coomassie staining**

Lane 1: MNK2
Lane 2: Protein Marker

MW: 74 kDa
Purity: ≥70%



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