

Data Sheet

HIPK4

Human, recombinant, N-terminal GST tag

Catalog #: 40111

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:

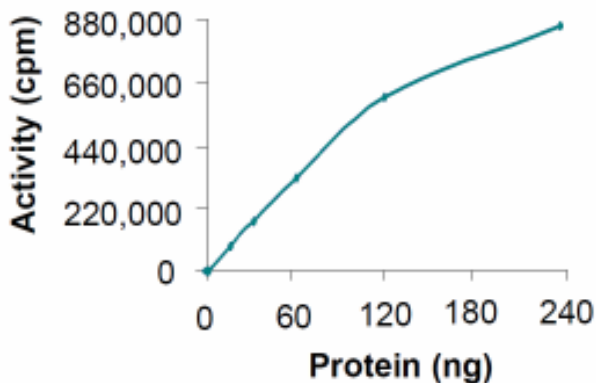
1. Arai, S., *et al.* *FEBS Lett.* **581(29):** 5649-5657 (2007).
2. Di Stefano, V., *et al.* *Exp. Cell Res.* **293(2):** 311-320 (2004).

Description: Human HIPK4 (homeodomain-interacting protein kinase 4), GenBank Accession No. NM_144685, full length, with N-terminal GST tag, MW ~110 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

Specific Activity: 202 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 [³²P]-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.

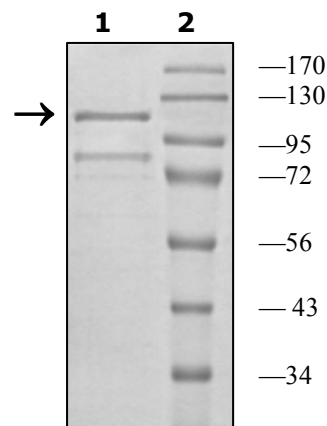
Quality Assurance



10% SDS-PAGE
Coomassie staining

Lane 1: HIPK4
Lane 2: Protein Marker

MW: 110 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