

Data Sheet

HIPK1

Human, recombinant, N-terminal GST tag

Catalog #: 40179

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. Ecsedy, J.A., *et al. Mol. Cell Biol.* **23**: 950-960 (2003).
2. Kondo, S., *et al. Proc. Natl. Acad. Sci. USA* **100**: 5431-5436 (2003).

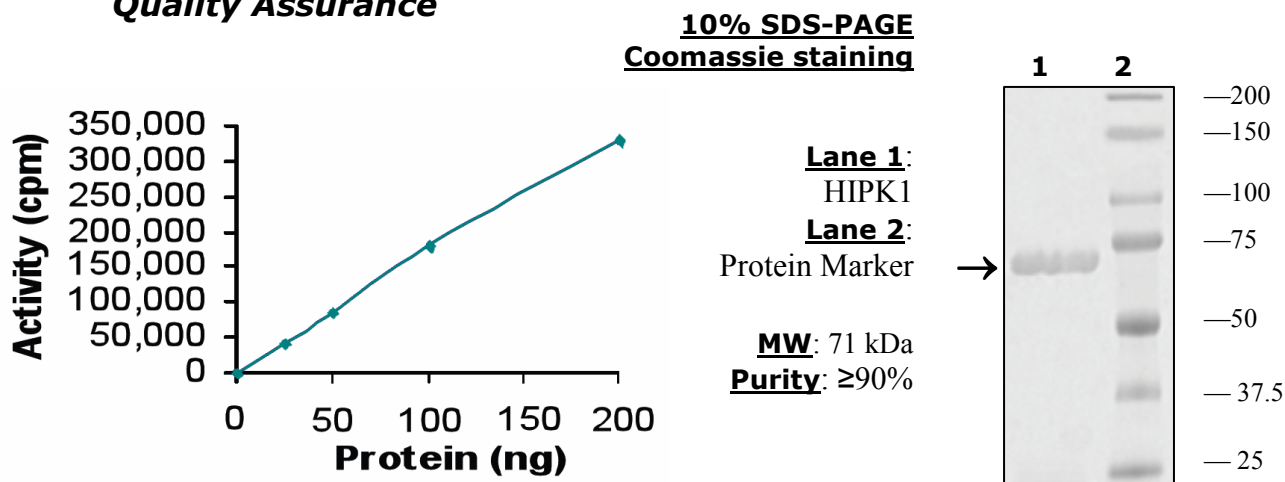
Description: Human HIPK1 (homeodomain-interacting protein kinase 1), also known as Myak and Nbak2 (GenBank Accession No. NM_152696), full length, with N-terminal GST tag, MW ~71 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

Specific Activity: 79 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 HIPK1 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

Quality Assurance



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