

Data Sheet

HPK1, GST-Tag, His-Avi-Tag, Biotin Labeled

Human, Recombinant, N-terminal GST tag, C-Terminal His-Avi-tag

Catalog #: 40399

Lot #: 161026

Conc.: 0.12 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 20% glycerol, 300 mM Imidazole

Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. Hu, M., *et al.*, *Genes Dev.* 1996; **10**: 2251-2264
2. Shui, J.R., *et al.*, *Nat. Immunol.* 2007; **8**: 84-91.

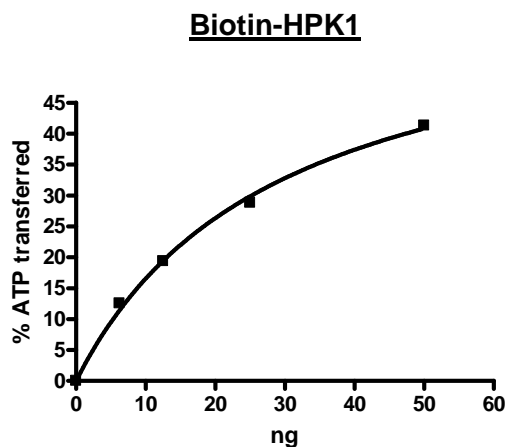
Description: Human hematopoietic progenitor kinase 1 (HPK1) also known as mitogen activated protein kinase 1 (MAP4K1), GenBank Accession No. NM_007181, a.a. 1-346, with N-terminal GST tag and C-terminal His-Avi-tag, MW = 67 kDa. HPK1 is expressed in Sf9 insect cells via a baculovirus expression system.

Specific Activity: ≥430 pmol/min/μg

Assay Description: Assay was done in Kinase buffer containing 1 mM DTT using MBP substrate (0.2 mg/ml) and 20 μM ATP. Reaction was done at 30°C for 45 min. Amount of ATP transferred was calculated using Kinase-Glo reagent (Promega)

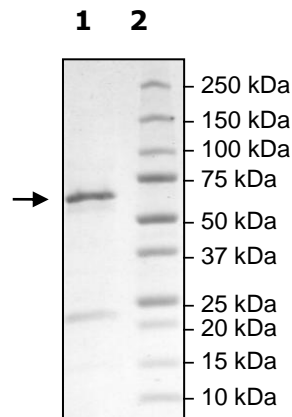
Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
HPK1
Lane 2:
Protein Marker
Purity: ≥72%
MW:
67 kDa



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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com