

Data Sheet

VRK1, GST-tag

Human, recombinant, N-terminal GST-tag

Catalog #: 40397

Lot#: 130816

Conc.: 0.1 mg/ml

Formulated in: 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM glutathione, 0.1 mM EDTA, 0.1 mM PMSF, 25% glycerol, and 0.25 mM DTT.

Stability: 12 months at -70°C . Avoid freeze/thaw cycles.

References:

1. Nezu, J., *et al.*, *Genomics* 1997;**45**:327-331.
2. Renbaum, P., *et al.*, *Am. J. Hum. Genet.* 2009;**85**:281-289.

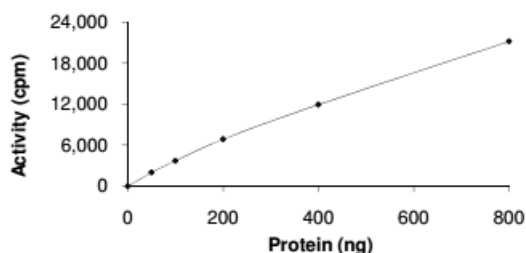
Description: Human vaccinia related kinase 1 (VRK1), GenBank Accession No. NM_003384), full length with N-terminal GST tag, MW=71 kDa, expressed in Sf9 insect cells.

Specific Activity: 2 pmol/min/ μg
 Assay buffer: 5 mM MOPS, 2.5 mM β -glycerophosphate, 5 mM MgCl_2 , 1 mM EGTA, 0.4 mM EDTA, 0.05 mM DTT, 2 mM ATP. Incubate VRK1 with 0.25 mg/ml dephosphorylated casein as substrate and $[^{33}\text{P}]\text{-ATP}$ 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

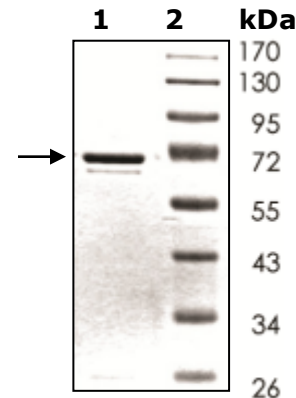
Specific Activity



10% SDS-PAGE
Coomassie staining

Lane 1: VRK1
Lane 2: Protein Marker

MW: 71 kDa
Purity: >95%



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 **1.858.259.5734**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com