

Data Sheet

PDK1, GST-tag

Human, recombinant, N-terminal GST-tag
Catalog #: 40080
Lot #: 151111-G2 **Conc.:** 0.13 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT and 20% Glycerol.

Stability: At least 6 months from date of receipt at -80°C.

References:

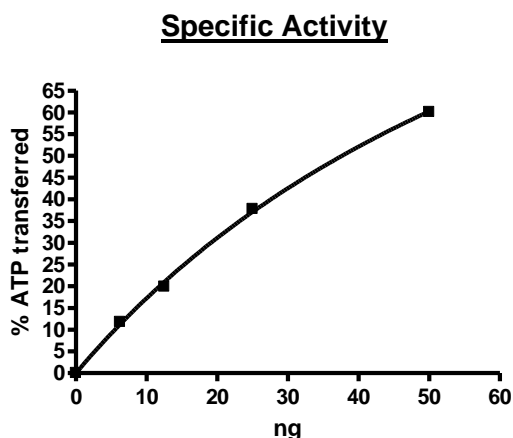
1. Yang, KJ., *et al. J Biol Chem.* 2008 Jan 18;**283(3)**:1480-91.
2. Kikani, CK., *et al. J Cell Biochem.* 2005 Dec 15;**96(6)**:1157-62.

Description: Human PDK1, GenBank Accession No. NM002613, a.a 1-556(full length) with N-terminal GST-tag, expressed in a Baculovirus infected Sf9 cell expression system MW=89 kDa.

Specific Activity: 310 pmol/min/μg
Assay Conditions: Assay was done in Kinase buffer containing 0.2 mM DTT using PDKtide 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

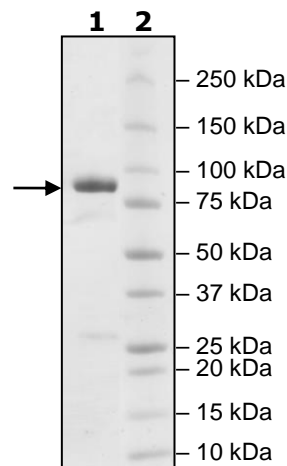


4-20% SDS-PAGE
Coomassie staining

Lane 1:
 2 μg PDK1
Lane 2:
 Protein Marker

MW: 89 kDa

Purity: ≥84%



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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com