

Data Sheet

PDK1

Human, recombinant, N-terminal GST tag

Catalog #: 40080

Lot#: 90610

Conc.: 0.5 mg/ml

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

Stability: >6 months at -80°C

References:

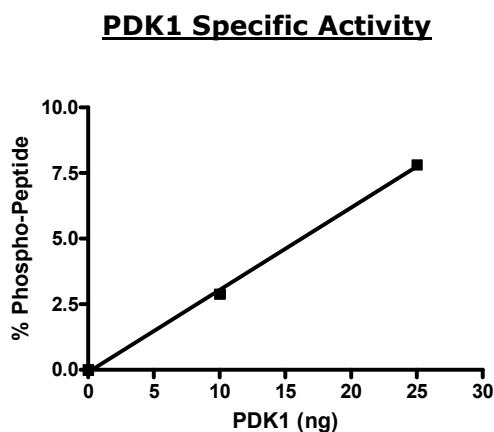
1. Yang, KJ., et al. J Biol Chem. 2008 Jan 18;**283(3)**:1480-91.
2. Kikani, CK., et al. J Cell Biochem. 2005 Dec 15;**96(6)**:1157-62.

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

Specific Activity: 5.02 pmol/min/μg
Enzyme reaction is conducted in a buffer containing 50 mM HEPES (pH7.5), 10 mM MgCl₂, 1 mM EGTA, 200 μM ATP, 0.01% Brij-35 and 2 μM substrate (SER/THR 6, Invitrogen) at room temperature for 1 hour.

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

Quality Assurance

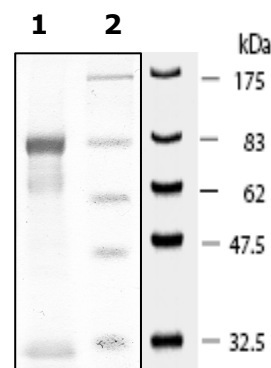


10% SDS-PAGE Coomassie staining

Lane 1:
6 μg PDK1

Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 89 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