

Data Sheet

BRK, His-Tag

Human, Recombinant, N-terminal GST-tag

Catalog #: 40403

Lot #: 151203

Conc.: 0.28 mg/ml

Formulated in: 40mM Tris pH 8.0, 110mM NaCl, 2.2mM KCl, 0.04% Tween-20, 20% glycerol, 3mM DTT and variable amounts of glutathione.

Stability: >12 months at -80°C. Avoid freeze/thaw cycles.

References:

1. Mitchell, P.J., *et al. Oncogene* **9(8):** 2383-2390 (1994).
2. Mitchell, P.J., *et al. Oncogene* **15(12):** 1497-1502 (1997).

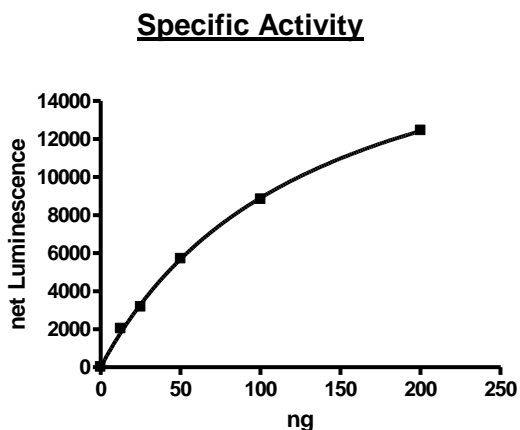
Description: Human BRK, also known as PTK6 (GenBank Accession No. NM_005975), 2-451 (end), with N-terminal GST-tag, MW=80 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

Specific Activity: 88 pmol/min/μg

Enzyme reaction was conducted in a buffer (25 μl) containing 40 mM Tris-HCl (pH 7.5), 20 mM MgCl₂, 100 μM ATP, 0.2 mg/ml poly-(Glu, Tyr), and BRK for 40 min at 30°C. ATP reduction is detected using ADP-Glo™ Kinase Assay Platform (Promega Corporation, Madison, WI).

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

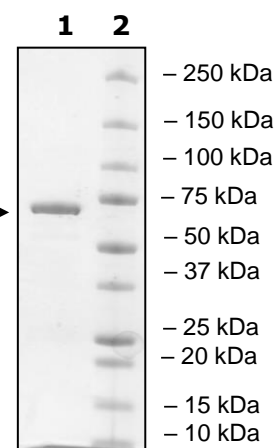
Quality Assurance



4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg BRK
Lane 2:
Protein Marker

MW: 80 kDa
Purity: ≥65%



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