



11526 Sorrento Valley Rd SteA2
San Diego Ca 92121
Tel: 1.858.202.1401
1.858.829.3082
Fax: 1.858.259.5734
Email: info@bpsbioscience.com

Data Sheet

BRX, Active

human, recombinant, N-terminal GST tag
Catalog #: 40403

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

Stability: >6 months at -80°C

References:

1. Mitchell, PJ. et al. Oncogene. 1994 Aug;9(8):2383-90.
2. Mitchell, PJ. et al. Oncogene. 1997 Sep 18;15(12):1497-502. Erratum in: Oncogene 1998 Jul 9;17(1):129.

Description:

Recombinant full-length human BRK was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is NM_005975.

Specific Activity:

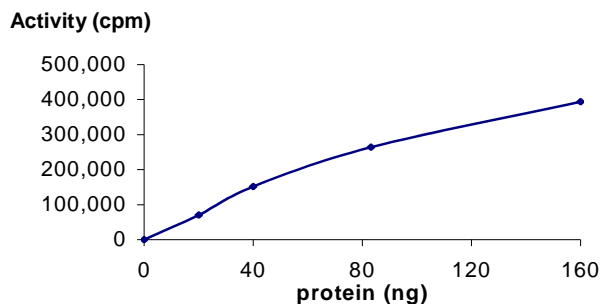
133 nmol/min/mg

Application:

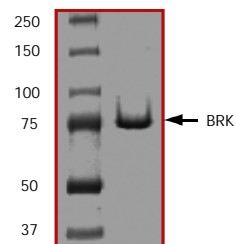
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

BRX Activity
133 nmol/min/mg



10% SDS-PAGE
Coomassie staining



MW: 80 kDa.

Purity: 90%

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** or **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