

Data Sheet

TRKB(NTKR2)

Human, recombinant, N-terminal GST-tag

Catalog #: 40281

Lot #: 130826

Conc.: 1.0 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0. 124 mM NaCl, 2.4 mM KCl, 3 mM DTT, 5 mg/ml Glutathione and 10% glycerol.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

References:

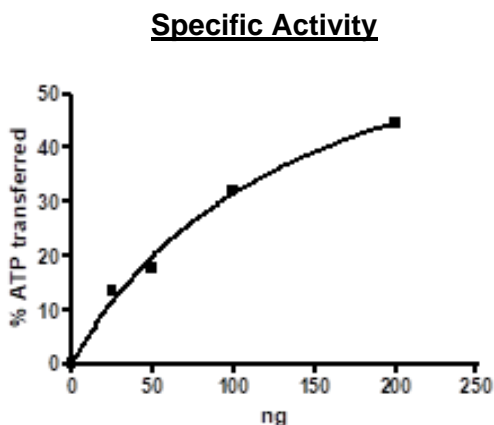
1. Naylor, R.L., *et al.* (2002). *Biochem. Biophys. Res. Commun.* **291 (3)**: 501-7.
2. Soppet, D., *et al.* (1991). *Cell* **65 (5)**: 895-903.
3. Squinto, S.P., *et al.* (1991). *Cell* **65 (5)**: 885-893.

Description: Human TRKB, also known as NTKR2 (Neurotrophic Tyrosine Kinase Receptor, type 2) or BDNF/NT-3 Receptor, GenBank Accession No. NM_001018064, a.a. 456-822(end), with N-terminal GST-tag, MW=67.5 kDa, expressed in Sf9 cells via a baculovirus expression system.

Specific Activity: ≥ 76 pmole/min/ μg
Assay Conditions: 40 mM Tris-HCl pH 7.4, 20 mM MgCl_2 , 0.1 mg/mL BSA and 1 mM DTT using 0.2 mg/mL Poly (Glu, Tyr) substrate and 20 μM ATP. Reaction was done at 30°C for 45 min. The amount of ATP transferred was calculated using Kinase-Glo reagent from Promega.

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

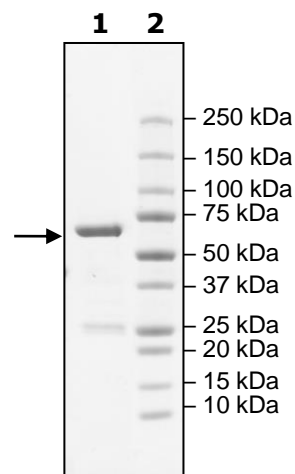
Quality Assurance



**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2.1 μg TrkB
Lane 2:
Protein Marker

MW: 67.5 kDa
Purity: $\geq 80\%$



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