

Data Sheet

TRKA(NRTK1), GST-Tag

Human, recombinant, N-terminal GST-tag

Catalog #: 40280

Lot#: 151113-G

Conc: 0.68 mg/mL

Formulated in: 40 mM Tris-HCl pH8.0, 110mM NaCl, 2.2 mM KCl, 0.04% Tween20, 3mM DTT, and 20% glycerol

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

References:

1. Koch, A., et al. (2000). *FEBS Lett.* **469** (1): 72–62.
2. Wiesmann, C., et al. (1999). *Nature* **401** (6749): 184–83.
3. Yamashita, H. et al. (1999). *J. Biol. Chem.* **274** (21): 15059–65

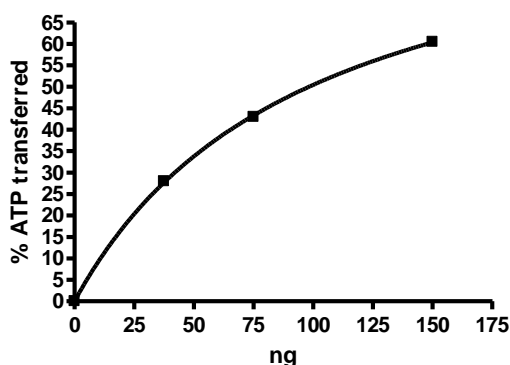
Description: Human TRKA, also known as NTKR1, and Neurotrophic Tyrosine Kinase Receptor, type 1, GenBank Accession No. NM_002529, a.a. 442-796(end), with N-terminal GST-tag, expressed in Sf9 cells via a baculovirus expression system. MW=65 kDa.

Specific Activity: 100 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. 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

Specific Activity



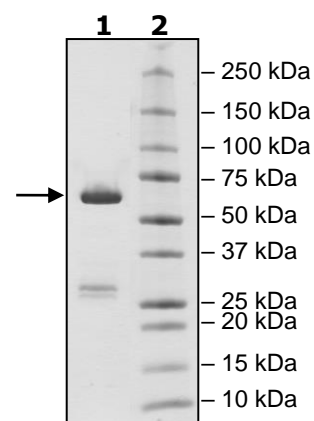
4-20% SDS-PAGE
Coomassie staining

Lane 1:
3 μg TrkA

Lane 2:
Protein Marker

MW: 67 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