

Data Sheet

TRKC(NTKR3)

Human, recombinant, N-terminal GST -tag
Catalog #: 40282
Lot #: 101007 **Conc.:** 0.73 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0,
 124 mM NaCl, 2.4 mM KCl, 20 mM
 glutathione, 10% glycerol
 and 3 mM DTT.

Stability: >6 months at -80°C

References:

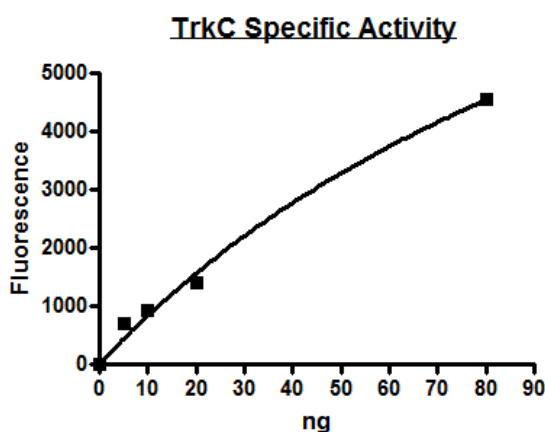
1. Lamballe, F., *et al.* (1991). *Cell* **66** (5): 967-979.
2. Tessarollo, L., *et al.* (1993). *Development* **118** (2): 463-75.
3. McGregor, L.M., *et al.* (1994). *Genomics* **22** (2): 267-72.

Description: Human TRKC, also known as NTKR3 (Neurotrophic Tyrosine Kinase Receptor, type 3) or NT-3 Receptor, GenBank Accession No. NM_002530, a.a. 456-825(end), with N-terminal GST-tag, MW=68kDa, expressed in Sf9 cells via a baculovirus expression system.

Specific Activity: 1.0 pmole/min/μg
 Activity determined using a DELFIA-assay with a poly-(Glu:Tyr 4:1)-biotin-coated plate, for 30 min. at room temperature.

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

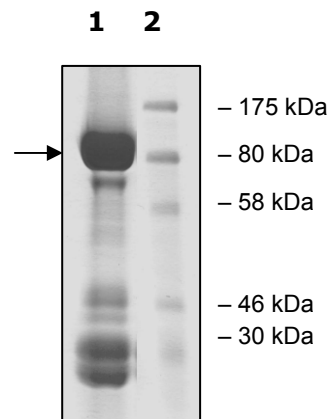
Quality Assurance



**10% SDS-PAGE
 Coomassie staining**

Lane 1:
 12 μg TRKC
Lane 2:
 Protein Marker
 BioLabs (#P7708L)

MW: 68 kDa
Purity: ≥38%



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