

ProQinase™ TRK-C

Neurotrophic tyrosine kinase, receptor, type 3

Recombinant Human Active Protein Kinase

HGNC Symbol: NTRK3

Synonyms: Gp145-TrkC

Product No.: 0775-0000-1

Lot: 005

Description: Human TRK-C, C-terminal fragment, amino acids V₅₁₀-G₈₂₅ (as in NCBI/Protein entry NP_002521.2), activated, N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: TRK-C Lot 005, was confirmed as TRK-C by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 64,870 Da

Expression: Baculovirus infected Sf9 cells

Purification: GST-Affinity Chromatography

Activation: in vitro autoactivation

Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20% glycerol

Storage temperature: -80°C

For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.085 µg/µl

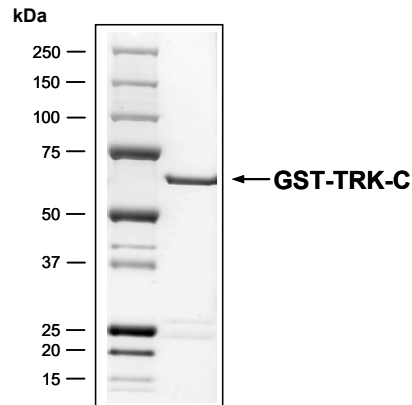
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

Specific kinase activity (P_i transfer): 35 pmol/µg×min

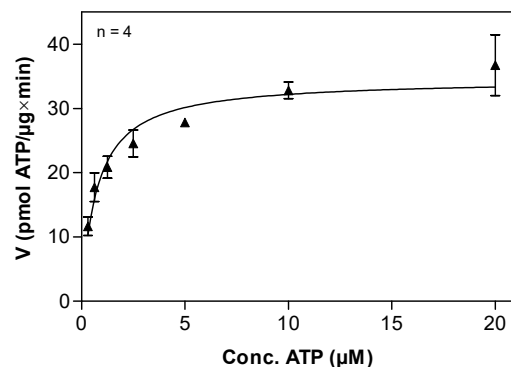
ATP-K_M: 0.7 µM

**TRK-C Lot 005:
Coomassie stain**



2.0 µg GST-TRK-C

**TRK-C Lot 005:
Determination of V_{max} and K_M value for ATP**



Determination of K_M value & Specific activity:

- Assay conditions:
 - 60 mM HEPES-NaOH, pH 7.5
 - 3 mM MgCl₂
 - 3 mM MnCl₂
 - 3 µM Na-orthovanadate
 - 1.2 mM DTT
 - 50 µg / ml PEG_{20,000}
 - ATP (variable)
 - Substrate: Poly(Glu:Tyr)_{4,1} (Sigma P-0275), 40 µg / ml
 - TRK-C: 0.4 µg / ml
- Filter binding assay
- MSFC membrane (Millipore)

Recombinant Proteins

Additional assay technology: TRK-C Lot 005

was also successfully tested by Reaction Biology for the use with the ADP-Glo™ Kinase assay from ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

ProQinase™ TRK-C

Product No.: 0775-0000-1

TRK-C Recombinant Fusion Protein Amino Acid Sequence

1	MSPILGYWKI	KGLVQPTLL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQ SMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQG WQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVLFGQ	240
241	PLAMFVIENP	QYFRQGHNCH	KPDTYVQH IK	RRDIVLKREL	GEGAFGKVFL	AECYNLSPTK	300
301	DKMLVAVKAL	KDPTLAARKD	FQREAE LLTN	LQHEHIVK FY	GVCGDG DPLI	MVFEYMKHGD	360
361	LNKFLRAHGP	DAMILVDGQP	RQAKGELGLS	QMLHIASQIA	SGMVYLA SQH	FVHRDLATRN	420
421	CLVGANLLVK	IGDFGMSRDV	YSTDYRVGG	HTMLPIRWMP	PESIMYRKFT	TESDVWSFGV	480
481	ILWEIFTY GK	QPWFQLS NTE	VIECITQGRV	LERPRVCPKE	VYDVMLGCWQ	REPQQR LNIK	540
541	EIYKILHALG	KATPIYLDIL	G				600

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: TRK-C fragment

TRK-C wt¹ amino acid sequence

1	MDVSLCPAKC	SFWRIFLLGS	VWLDYVGSVL	ACPANCVCSK	TEINCRRPDD	GNLFPLLEGQ	60
61	DSGNSNGNAS	INITDISRNI	TSIHENWRS	LHTLNAV DME	LYTGLQK LTI	KNSGLRSIQP	120
121	RAFAKNPHLR	YINLSSNRLT	TLSWQLFQTL	SLRELQLEQN	FFNCSCDIRW	MQLWQE QGEA	180
181	KLNSQNLYCI	NADGSQ LPLF	RMNISQCDLP	EISVSHVNL T	VREGDNAVIT	CNGSGSPLPD	240
241	VDWIVTGLQS	INTHQTNL NW	TNVHAINLTL	VNVTS EDNGF	TLTCIAENVV	GMSNASVALT	300
301	VYYP PVRVSL	EEPELRLEHC	IEFVVRGNPP	PTLHWHLNGQ	PLRESKI IHV	EYYQEGEISE	360
361	GCLLFNKPTH	YNNGN YTLIA	KNPLGTANQT	INGHFLKEPF	PESTDNFILF	DEVSP TPPIT	420
421	VTHKPEEDTF	GVSI AVGLAA	FACVLLV VLF	VMINKYGRRS	KFGMKGPVAV	ISGEEDSASP	480
481	LHHINH GITT	PSSLDAGPDT	VVIGMTRIPV	IENPQYFRQ	HNCHKPDTYV	QHIKRRDIVL	540
541	KRELGEGAFG	KVFLAECYNL	SPTKDKMLVA	VKALKDPTLA	ARKDFQREAE	LLTNLQHEHI	600
600	VKfygvcgdg	DPLIMVFEYM	KHGD LNKFLR	AHGPNAMILV	DGQPRQAKGE	LGLS QMLHIA	660
661	SQIASGMVYL	ASQH FVHRDL	ATRNCLVGAN	LLVKIGDFGM	SRDVYSTDY	RVGGHTMLPI	720
721	RWMPPESIMY	RKFTTESDVW	SFGVILWEIF	TYGKQ PWFQL	SNTEVIECIT	QGRVLERPRV	780
781	CPKEVYDVML	GCWQRE PQR	LNIKEIYKIL	HALGKATPIY	LDILG		840

blue: TRK-C sequence expressed in fusionprotein

¹NCBI/Protein accession number NP_002521.2

Recombinant Proteins