

ProQinase™ BRK

protein tyrosine kinase 6

Recombinant Human Active Protein Kinase

HGNC Symbol: PTK6

Synonyms: n/a

Product No.: 0181-0000-1

Lot: 004

Description: Human BRK, full length, amino acids M₁-T₄₅₁ (as in [NCBI/Protein](#) entry NP_005966.1), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: BRK Lot 004, was confirmed as BRK by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 81,790 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: This kinase was not activated by special procedures

Storage buffer: 50 mM TRIS-HCl pH 8.0, 100 mM NaCl, 5 mM DTT, 4 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: 0.160 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

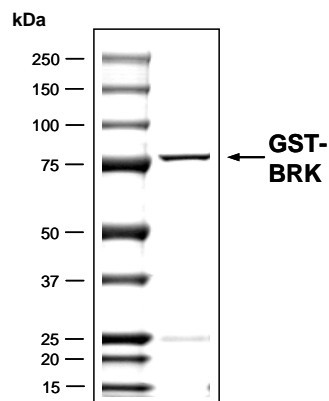
Biochemical Parameters:

Specific kinase activity (P_i transfer): 31 pmol/µg × min
ATP-K_M: 4.1 µM

Additional assay technology:

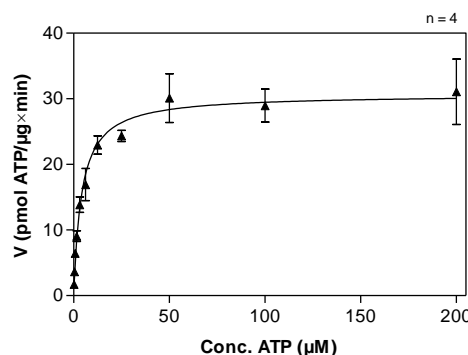
BRK Lot 004 was also successfully tested by Reaction Biology for the use with the ADP-Glo™ Kinase assay from Promega ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

BRK Lot 004: Coomassie stain



2.0 µg GST-BRK

BRK Lot 004: Determination of V_{max} and K_M value for ATP



- 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} 20 µg/ml
Kinase: 1 µg/ml
- Filter binding assay
MSFC membrane (Millipore)

ProQinase™ BRK

Product No.: 0181-0000-1

GST-BRK Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQG WQATF	GGGDHPPKSD	PMGHHHHHG	RRRASVAAGI	240
241	LVPRGSPGLD	GIYARGIQAS	MVSRDQAHLG	PKYVGLWDFK	SRTDEELSFR	AGDVFHVARK	300
301	EEQWWATLL	DEAGGAVAQG	YVPHNYLAER	ETVESEPWF	GYISRSEAVR	RLQAEGNATG	360
361	AFLIRVSEKP	SADYVLSVRD	TQAVRHYKI	RRAGGRLLHN	EAVSFLSLPE	LVNYHRAQSL	420
421	SHGLRLA APC	RKHEPEPLPH	WDDWERPREE	FTLCRKLGS	YFGEVFEGW	KDRVQVAIKV	480
481	ISRDNLLHQ	MLQSEIQAMK	KLRHKHILAL	YAVVSVGDPV	YIITELMAK	SLELLRDS	540
541	EKVLVSELL	DIAWQVAEGM	CYLESQNYIH	RDLAARNILV	GENTLCKVGD	FGLARLIKED	600
601	VYLSHDHNIP	YKWTAPEALS	RGHYSTKSDV	WSFGILLHEM	FSRGQVPYPG	MSNHEAFLRV	660
661	DAGYRMP CPL	ECPPSVHKLM	LTCWCRDPEQ	RPCFKALRER	LSSFTSYENP	T	720

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue: BRK

BRK wt ¹ Amino Acid Sequence							
1	MVSRDQAHLG	PKYVGLWDFK	SRTDEELSFR	AGDVFHVARK	EEQWWATLL	DEAGGAVAQG	60
61	YVPHNYLAER	ETVESEPWF	GISRSEAVR	RLQAEGNATG	AFLIRVSEKP	SADYVLSVRD	120
121	TQAVRHYKI	RRAGGRLLHN	EAVSFLSLPE	LVNYHRAQSL	SHGLRLA APC	RKHEPEPLPH	180
181	WDDWERPREE	FTLCRKLGS	YFGEVFEGW	KDRVQVAIKV	ISRDNLLHQ	MLQSEIQAMK	240
241	KLRHKHILAL	YAVVSVGDPV	YIITELMAK	SLELLRDS	EKVLVSELL	DIAWQVAEGM	300
301	CYLESQNYIH	RDLAARNILV	GENTLCKVGD	FGLARLIKED	VYLSHDHNIP	YKWTAPEALS	360
361	RGHYSTKSDV	WSFGILLHEM	FSRGQVPYPG	MSNHEAFLRV	DAGYRMP CPL	ECPPSVHKLM	420
421	LTCWCRDPEQ	RPCFKALRER	LSSFTSYENP	T			480

blue: kinase sequence expressed in recombinant protein Red: variant in recombinant protein

¹[NCBI/Protein](https://www.ncbi.nlm.nih.gov/protein/1005966.1) accession number NP_005966.1