

ProQinase™ B-RAF wt

B-Raf proto-oncogene, serine/threonine kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: BRAF

Synonyms: B-Raf, BRAF1, RAFB1, B-raf 1

Product No.: 0403-0000-1

Lot: 001

Description: Human B-RAF, C-terminal fragment, amino acids Q₄₁₇-H₇₆₆ (as in [NCBI/Protein](#) entry NP_004324.2), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: B-RAF wt Lot 001, was confirmed as B-RAF by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 69,127 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 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: 0.294 µg/µl

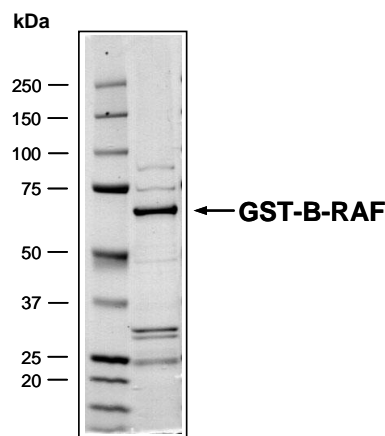
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

Specific kinase activity (P_i transfer): 139 pmol/µg × min

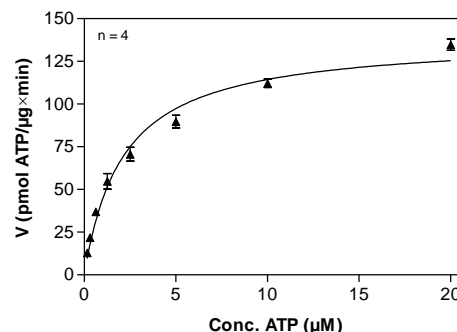
ATP-K_M: 2.1 µM

B-RAF wt Lot 001: Coomassie stain



2.0 µg GST-B-RAF

B-RAF wt Lot 001: 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: MEK1 K97M, 40 µg/ml
- Kinase: 0.5 µg/ml

- Filter binding assay
- MSFC membrane (Millipore)

Additional assay technology:

B-RAF wt Lot 001 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

ProQinase™ B-RAF wt

Product No.: 0403-0000-1

GST-B-RAF wt Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWAQTF	GGGDHPPKSD	PMG HHHHHG	RRRASVAAGI	240
241	LVPRGSPGLD	GICSRNS QKS	PGPQERKSS	SSSEDNRMK	TLGRRDSSDD	WEIPDGQITV	300
301	GQRIGSGSFG	TVYKKGWHGD	VAVKMLNVTA	PTPQQLQAFK	NEVGVLKTR	HVNILLFMGY	360
361	STKPQLAIVT	QWCEGSSLYH	HLHIETKFE	MIKLIDIARQ	TAQGM DYLHA	KSIHRDLKS	420
421	NNIFLHEDLT	VKIGDFGLAT	VKSRWSGSHQ	FEQLSGSILW	MAPEVIRMQD	KNPYSFQSDV	480
481	YAFGIVLYEL	MTGQLPYSNI	NNRDQIIFMV	GRGYLSPDLS	KVRSNCPKAM	KRLMAECLKK	540
541	KRDERPLFPQ	ILASIELLAR	SLPKIHRAS	EPSLNRAGFQ	TEDFSLYACA	SPKTPIQAGG	600
601	YGAFPVH						660

1-218: GST **Red**: HIS6-tag **Pink**: Thrombin cleavage site **blue**: B-RAF wt fragment **boxed**: variation from RefSeq

B-RAF wt ¹ Amino Acid Sequence							
1	MAALSGGGGG	GAEPGQALFN	GDMEPEAGAG	AGAAASSAAD	PAIPEEVWNI	KQMIKLTQEH	60
61	IEALLDKFGG	EHNPPSIYLE	AYEEYTSKLD	ALQOREQQLL	ESLGNGTDFS	VSSSASMDTV	120
121	TSSSSSSLSV	LPSSLSVFQN	PTDVARSNPK	SPQKPIRVVF	LPNKQRTVVP	ARCGVTVRDS	180
181	LKKALMMRGL	IPECCAVYRI	QDGEKKPIGW	DTDISWLTGE	ELHVEVLENV	PLTTHNFVRK	240
241	TFFTLAFCDF	CRKLLFQGF	CQTCGYKFHQ	RCSTEVPLMC	VNYDQLDLLE	VSKFFEHHP	300
301	PQEEASLAET	ALTSGSSPSA	PASDSIGPQI	LTSPSPSKSI	PIPQPPRPAD	EDHRNQFGQR	360
361	DRSSAPNVH	INTIEPVNID	DLIRDQGFRG	DGGSTTGLSA	TPPASLPGSL	TNVKAL QKSP	420
421	GPQERKSSS	SSEDNRMKT	LGRRDSSDDW	EIPDGQITVG	QRIGSGSFGT	VYKKGWHGDV	480
481	AVKMLNVTAP	TPQQLQAFKN	EVGVLKTRH	VNILLFMGYS	TKPQLAIVTQ	WCEGSSLYHH	540
541	LHIETKFEM	IKLIDIARQT	AQGM DYLHAK	SIIHRDLKSN	NIFLHEDLTV	KIGDFGLATV	600
601	KSRWSGSHQF	EQLSGSILWM	APEVIRMQDK	NPYSFQSDVY	AFGIVLYELM	TGQLPYSNIN	660
661	NRDQIIFMVG	RGYLSPLSK	VRSNCPKAMK	RLMAECLKKK	RDERPLFPQI	LASIELLARS	720
721	LPKIHRSASE	PSLNRAGFQT	EDFSLYACAS	PKTPIQAGGY	GAFPVH		780

blue: B-RAF sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_004324.2