

ProQinase™ STK17B

serine/threonine kinase 17b

Recombinant Human Active Protein Kinase

HGNC Symbol: STK17B

Synonyms: DRAK2

Product No.: 1504-0000-1

Lot: 001

Description: Human Kinase, full length, amino acids M₁-C₃₇₂ (as in [NCBI/Protein](#) entry NP_004217.1), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: STK17B Lot 001, was confirmed as STK17B by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 70,434 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: This kinase was not activated by special procedures

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 10 µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.634 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

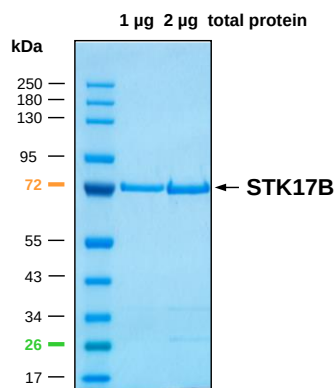
Biochemical Parameters:

Specific kinase activity (P_i transfer): 1.1 pmol/µg x min

ATP-K_M: 0.2 µM

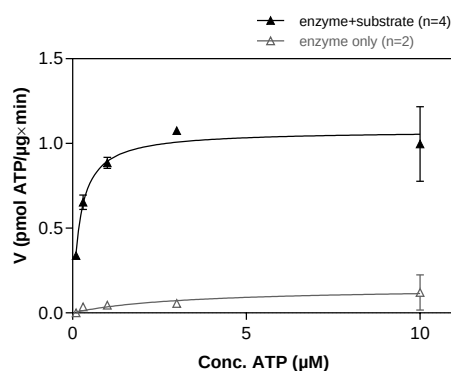
STK17B Lot 001:

Coomassie stain



STK17B 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: MRCL3 40 µg/ml
Kinase: 3 µg/ml
- Filter binding assay
MSFC membrane (Millipore)

Sequence information

GST-STK17B Recombinant Fusion Protein Amino Acid Sequence

1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLIERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHGG	RDSLEVLFGG	240
241	MSRRRFDCR	SISGLLTTP	QPIKMENFN	NFYILTSKEL	GRGKFAVVRQ	CISKSTGQEY	300
301	AAKFLKKRRR	QDCRAEILH	EIAVLELAKS	CPRVINLHEV	YENTSEIILI	LEYAAGGEIF	360
361	SLCLPELAEM	VSENDVIRLI	KQILEGVYYL	HQNNIVHLDL	KPQNILLSSI	YPLGDIKIVD	420
421	FGMSRKIGHA	CELREIMGTP	EYLAPEILNY	DPITTATDMW	NIGIIAYMLL	THTSPFVGED	480
481	NQETYNISQ	VNVDYSEETF	SSVSQLATDF	IQSLLVKNPE	KRPTAEICLS	HSWLQQWDFE	540
541	NLFHPEETSS	SSQTQDHSVR	SSDKTSKSS	CNGTCGDRED	KENIPEDSSM	VSKRFRFDDS	600
601	LPNPHELVD	LLC					660

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: STK17B

STK17B wt¹ Amino Acid Sequence

1	MSRRRFDCRS	ISGLLTTPQ	IPIKMENFNN	FYILTSKELG	RGKFAVVRQC	ISKSTGQEYA	60
61	AKFLKKRRRG	QDCRAEILHE	IAVLELAKSC	PRVINLHEVY	ENTSEIILIL	EYAAGGEIFS	120
121	LCLPELAEMV	SENDVIRLIK	QILEGVYYLH	QNNIVHLDL	PQNILLSSY	PLGDIKIVDF	180
181	GMSRKIGHAC	ELREIMGTPE	YLAPEILNYD	PITTATDMWN	IGIIAYMLLT	HTSPFVGEDN	240
241	QETYNISQV	NVDYSEETF	SVSSQLATDFI	QSLLVKNPEK	RPTAEICLSH	SWLQQWDFEN	300
301	LHFPEETSSS	SQTQDHSVRS	SEDKTSKSSC	NGTCGDREDK	ENIPEDSSMV	SKRFRFDDSL	360
361	PNPHELVDL	LC					420

blue: STK17B sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_004217.1