

ProQinase™ TSK2

testis specific serine kinase 2

Recombinant Human Active Protein Kinase

HGNC Symbol: TSSK2

Synonyms: STK22B; SPOGA2

Product No.: 0220-0000-1

Lot: 002

Description: Human TSK2, full length, amino acids M₁-T₃₅₈ (as in [NCBI/Protein](#) entry NP_443732.3), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: TSK2 Lot 002, was confirmed as TSK2 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 73,916 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, 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: 0.378 µg/µl

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

Biochemical Parameters:

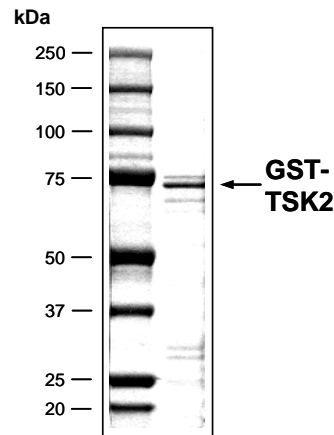
Specific kinase activity (P_i transfer): 5.9 pmol/µg × min

ATP-K_M: 0.2 µM

Additional assay technology:

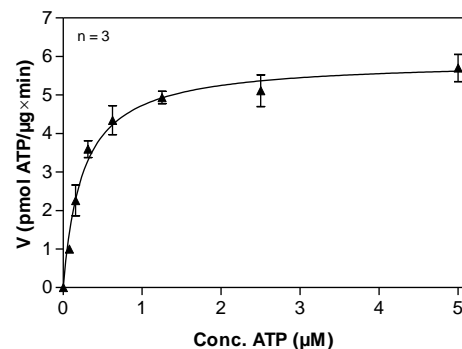
TSK2 Lot 002 was also successfully tested by Reaction Biology for the use with the ADP-Glo™ Kinase assay from Promega ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

TSK2 Lot 002:
Coomassie stain



SP002, 2.0 µg

TSK2 Lot 002:
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: Histone H2A, 50 µg/ml
 - TSK2: 4 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

ProQinase™ TSK2

Product No.: 0220-0000-1

GST-TSK2 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	PMGHHHHHG	RRRASVAAGI	240
241	LVPKSGPGLD	GIYARGIQAS	MGARGRQCDG	YLQNSPLMDD	ATVLRKKGVI	VGINLGKGSY	300
301	AKVKSAYSER	LKFNVAVKII	DRKKTPTDFV	ERFLPREMDI	LATVNHGSI	KTYEIFETSD	360
361	GRIYIIMELG	AQGDLLLEFIK	CQCALHEDVA	RKMFRQLSSA	VKYCHDLDIV	HRDLKCNLL	420
421	LDKDFNIKLS	DFGFSKRCLR	DSNGRIILSK	TFCGSAAYAA	PEVLQSIPIQ	PKVYDIWSLG	480
481	VILYIMVCGS	MPYDDSDIRK	MLRIQKEHRV	DFPRSKNLTC	ECKDLIYRML	QPDVSQRLHI	540
541	DEILSHSWLQ	PPKPKATSSA	SFKREGEKGY	RAECKLDTKT	GLRPDHRPDH	KLGAKTQHRL	600
601	LVPVENENRM	EDRLAETSRA	KDHHISGAEV	GKASTKGEFQ	HTGGRY		660

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue: TSK2 boxed: variation from RefSeq

TSK2 wt ¹ Amino Acid Sequence							
1	MDDATVLRKK	GYIVGINLGK	GSYAKVKSAY	SERLKFNVAV	KIIDRKKTPT	DFVERFLPRE	60
61	MDILATVNHG	SIIKTYEIFE	TSDGRIYIIM	ELGVQGDLE	FIKQOGALHE	DVARKMFRQL	120
121	SSAVKYCHDL	DIVHRDLKCE	NLLLDKDFNI	KLSDFGFSKR	CLRDSNGRI	LSKTFGGSAA	180
181	YAAPEVLQSI	PYQPKVYDIW	SLGVILYIMV	CGSMPYDDSD	IRKMLRIQKE	HRVDFPRSKN	240
241	LTCECKDLIY	RMLQPDVSR	LHIDEILSHS	WLQPPKPKAT	SSASFKREGE	GKYRAECKLD	300
301	TKTGLRPDHR	PDHKLGAQTQ	HRLLVVPENE	NRMEDRLAET	SRAKDHHISG	AEVGKAST	360

blue: TSK2 sequence expressed in recombinant protein Red: variant in recombinant protein

¹[NCBI/Protein](#) accession number NP_443732.3