

ProQinase™ STK32B

serine/threonine kinase 32B

Recombinant Human Active Protein Kinase

HGNC Symbol: STK32B

Synonyms: STKG6, YANK2, STK32, HSA250839

Product No.: 2213-0000-1

Lot: 002

Description: Human STK32B, amino acids M₁-S₄₁₇ (as in [NCBI/Protein](#) entry NP_060871.1), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: STK32B Lot 002, has been verified by mass spectrometry LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 75972 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: 0.314 µg/µl

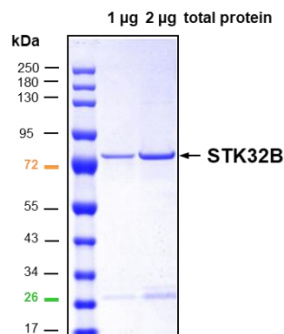
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

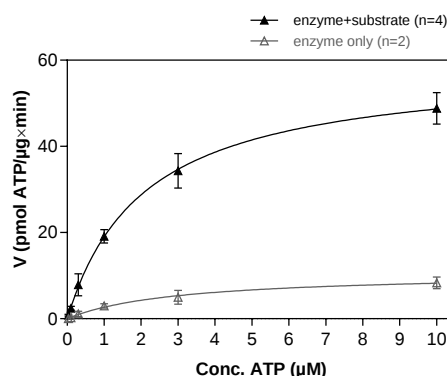
Specific kinase activity (P_i transfer): 59 pmol/µg*min

ATP-K_M: 2.1 µM

STK32B LOT002: Coomassie stain



STK32B LOT002: 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: Casein 20 µg/ml
- Kinase: 1 µg/ml

Assay technology:

Radiometric filter binding assay
MSFC membrane (96 well plate, Millipore)

Sequence information

GST-STK32B Recombinant Fusion Protein Amino Acid Sequence

1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLIERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHIG	RDSLEVLFGG	240
241	PMGGNHSKPP	PVFDENEEVN	FDHFQILRAI	GKGSFGKVCV	VQKRDTKKMY	AMKYMNKQKC	300
301	IERDEVNRVF	RELQIMQGLE	HPFLVNLWYS	FQDEEDMFMV	VDLLGGDLR	YHLQONVHFT	360
361	EGTVKLYICE	LALALEYLQR	YHIIHRDIKP	DNILLDEHGH	VHITDFNIAT	VVKAERASS	420
421	MAGTKPYMAP	EVFQVYMDRG	PGSYFVDWW	SLGITAYELL	RGWRPYEIH	VTPIDEILNM	480
481	FKVERVHYSS	TWCKGMVALL	RKLLTKDPES	RVSSLHDIQS	VPYLADMNWD	AVFKKALMPG	540
541	FVPNKGRLLC	DPTFELEEMI	LESKPLHKKK	KRLAKNRSRD	GTKDSCPLNG	HLQHCLETVR	600
601	EEFIIFNREK	LRRQQGQGSQ	LLDTSRGGG	QAQSKLQDGC	NNLLTHTCT	RGCSS	660

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: STK32B

STK32B wt¹ Amino Acid Sequence

1	MGGNHSKPP	VFDENEEVN	DHFQILRAIG	KGSFGKVCIV	QKRDTKKMYA	MKYMNKQKCI	60
61	ERDEVNRVFR	ELQIMQGLEH	PFLVNLWYSF	QDEEDMFMV	DLLGGDLRY	HLQONVHFTE	120
121	GTVKLYICEL	ALALEYLQRY	HIIHRDIKPD	NILLDEHGHV	HITDFNIATV	VKAERASSM	180
181	AGTKPYMAPE	VFQVYMDRGP	GYSYPVDWWS	LGITAYELLR	GWRPYEIHVS	TPIDEILNMF	240
241	KVERVHYSS	WCKGMVALLR	KLLTKDPESR	VSSLHDIQSV	PYLADMNWD	VFKKALMPGF	300
301	VPNKGRLLCD	PTFELEEMIL	ESKPLHKKKK	RLAKNRSRDG	TKDSCPLNGH	LQHCLETVRE	360
361	EFIIFNREKL	RRQQGQGSQ	LDTSRGGGQ	AQSKLQDGCN	NNLLTHTCTR	GCSS	420

blue: STK32B sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_060871.1