

ProQinase™ STK32A

serine/threonine kinase 32A

Recombinant Human Active Protein Kinase

HGNC Symbol: STK32A

Synonyms: MGC22688, YANK1

Product No.: 2205-0000-1

Lot: 002

Description: Human STK32A, full length, amino acids M₁-L₃₉₆ (as in [NCBI/Protein](#) entry NP_001106195.1), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: STK32A Lot 002, has been verified by mass spectrometry LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 74459 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!

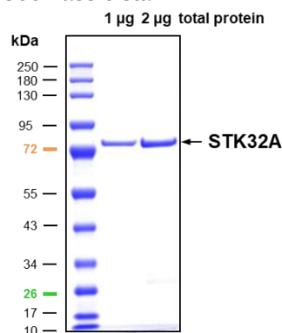
Protein concentration: 0.495 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

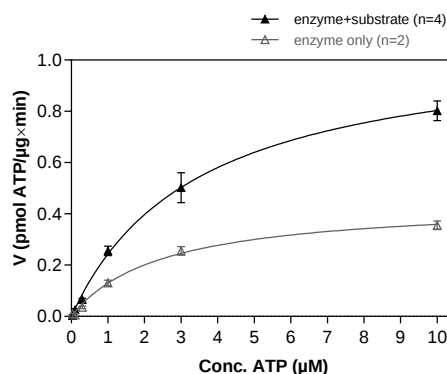
Specific kinase activity (P_i transfer): 1.1 pmol/µg*min

ATP-K_M: 3.4 µM

STK32A Lot 002: Coomassie stain



STK32A Lot 002: 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: 2 µg/ml

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

Sequence information

GST-STK32A 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	PMGHHHHHGG	RDSLEVLFGG	240
241	PMGANTSRKP	PVFDENEDVN	FDHFEILRAI	GKGSFGKVC	VQKNDTKKMY	AMKYMNKQKC	300
301	VERNEVRNVF	KELQIMQGLE	HPFLVNLWYS	FQDEEDMFMV	VDLLGGDLR	YHLQONVHFKE	360
361	EETVKLFICE	LVMALDYLQN	QRIIHRDMKP	DNILLDEHGH	VHITDFNIAA	MLPRETQITF	420
421	MAGTKPYMAP	EMFSSRKAG	YSFAVDWWSL	GVTAYELLRG	RRPYHIRSST	SSKEIVHTFE	480
481	TTVVTYPSAW	SQEMVSLKK	LLEPNPDQRF	QLSDVQNF	YMNDINWDAV	FQKRLIPGFI	540
541	PNKGRNLCDP	TFELEEMILE	SKPLHKKKKR	LAKKEKDMRK	CDSSQTCLLQ	EHLDSVQKEF	600
601	IIFNREKVNR	DFNKRQPNLA	LEQTKDPQGE	DGQNNNL			660

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: STK32A

STK32A wt¹ Amino Acid Sequence

1	MGANTSRKPP	VFDENEDVNF	DHFEILRAIG	KGSFGKVCIV	QKNDTKKMYA	MKYMNKQKCV	60
61	ERNEVRNVFK	ELQIMQGLEH	PFLVNLWYSF	QDEEDMFMV	DLLGGDLRY	HLQONVHFKE	120
121	ETVKLFICEL	VMALDYLQNQ	RIIHRDMKPD	NILLDEHGHV	HITDFNIAAM	LPRETQITTM	180
181	AGTKPYMAPE	MFSSRKAGY	SFAVDWWSLG	VTAYELLRGR	RPYHIRSSTS	SKEIVHTFET	240
241	TVVTYPSAWS	QEMVSLKKL	LEPNPDQRF	QLSDVQNF	YMNINWDAVF	QKRLIPGFIP	300
301	NKGRNLCDPT	FELEEMILES	KPLHKKKKRL	AKKEKDMRKC	DSSQTCLLQE	HLDSVQKEFI	360
361	IIFNREKVNR	DFNKRQPNLA	LEQTKDPQGE	DGQNNNL			420

blue: STK32A sequence expressed in recombinant protein

¹[NCBI/Protein](https://www.ncbi.nlm.nih.gov/Protein/NP_001106195.1) accession number NP_001106195.1