

ProQinase™ SRC (GST-HIS-tag)

SRC proto-oncogene, non-receptor tyrosine kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: SRC

Synonyms: ASV, SRC1

Product No.: 0200-0000-1

Lot: 005

Description: Human SRC, full length, amino acids M₁-L₅₃₆ (as in [NCBI/Protein](#) entry NP_005408.1), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: SRC Lot 005, was confirmed as SRC by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 89,744 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, 4 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: 0.239 µg/µl

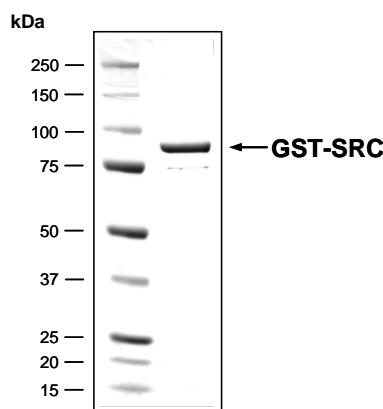
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

Specific kinase activity (P_i transfer): 254 pmol/µg × min

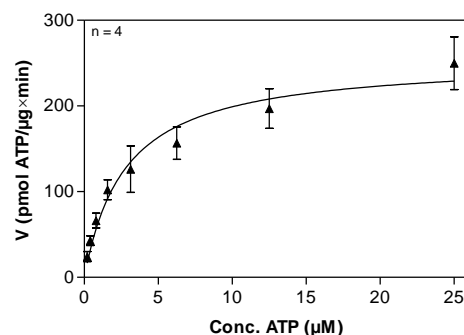
ATP-K_M: 2.8 µM

SRC Lot 005:
Coomassie stain



2.0 µg GST-SRC

SRC Lot 005:
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: Poly(Glu:Tyr)_{4:1}, 20 µg/ml
 - Kinase: 0.4 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

Additional assay technology:

SRC Lot 005 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™ SRC (GST-HIS-tag)

Product No.: 0200-0000-1

GST-SRC 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	WPLQGWQATF	GGGDHPPKSD	PMG HHHHHG	RRRASVAAGI	240
241	LVPRG SPGLD	GIYARGIQAS	M SNKSKPKD	ASQRRRSLEP	AENVHGAGGG	AFPASQTPSK	300
301	PASADGHRGP	SAAFAPAAAE	PKLFGGFNSS	DTVTSPQRAG	PLAGGVTFV	ALYDYESRTE	360
361	TDLSFKKGER	LQIVNNTGEG	WWLAHSLSTG	QTGYIPSNYV	APSDSIQAE	WYFGKITRRE	420
421	SERLLLNAEN	PRGTFLVRES	ETTKGAYCLS	VSDFDNAKGL	NVKHYKIRKL	DSGGFYITSR	480
481	TQFNSLQQLV	AYYSKHADGL	CHRLTVCPT	SKPQTQGLAK	DAWEIPRESL	RLEVKLGQGC	540
541	FGEVWMTWN	GTTRVAIKTL	KPGTMSPEAF	LQEAQVMKKL	RHEKLVQLYA	VVSEPIYIV	600
601	TEYMSKGSLL	DFLKGETGKY	LRLPQLVDMA	AQIASGMAYV	ERMNVHRDL	RAANILVGEN	660
661	LVCKVADFGL	ARLIEDNEYT	ARQGAKFPIK	WTAPEAALYG	RFTIKSDVWS	FGILLTELTT	720
721	KGRVPYPGMV	NREVLDQVER	GYRMPCPPEC	PESLHDLMCQ	CWRKEPEERP	TFEYLQAFLE	780
781	DYFTSTEPQY	QPGENL					840

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

SRC wt ¹ Amino Acid Sequence							
1	M SNKSKPKD	ASQRRRSLEP	AENVHGAGGG	AFPASQTPSK	PASADGHRGP	SAAFAPAAAE	60
61	PKLFGGFNSS	DTVTSPQRAG	PLAGGVTFV	ALYDYESRTE	TDLSFKKGER	LQIVNNTGEG	120
121	WWLAHSLSTG	QTGYIPSNYV	APSDSIQAE	WYFGKITRRE	SERLLLNAEN	PRGTFLVRES	180
181	ETTKGAYCLS	VSDFDNAKGL	NVKHYKIRKL	DSGGFYITSR	TQFNSLQQLV	AYYSKHADGL	240
241	CHRLTVCPT	SKPQTQGLAK	DAWEIPRESL	RLEVKLGQGC	FGEVWMTWN	GTTRVAIKTL	300
301	KPGTMSPEAF	LQEAQVMKKL	RHEKLVQLYA	VVSEPIYIV	TEYMSKGSLL	DFLKGETGKY	360
361	LRLPQLVDMA	AQIASGMAYV	ERMNVHRDL	RAANILVGEN	LVCKVADFGL	ARLIEDNEYT	420
421	ARQGAKFPIK	WTAPEAALYG	RFTIKSDVWS	FGILLTELTT	KGRVPYPGMV	NREVLDQVER	480
481	GYRMPCPPEC	PESLHDLMCQ	CWRKEPEERP	TFEYLQAFLE	DYFTSTEPQY	QPGENL	540

blue: SRC sequence expressed in recombinant protein **Red**: variant in recombinant protein

¹[NCBI/Protein](#) accession number NP_005408.1

G2A: legacy from cloning procedure