

FYN wt

FYN proto-oncogene, Src family tyrosine kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: FYN

Synonyms: MGC45350, SLK, SYN, p59-Fyn

Product No.: 0352-0000-1

Lot: 002

Description: Human FYN (Isoform A), full length, amino acids M₁-L₅₃₇ (as in [NCBI/Protein](#) entry NP_002028.1), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

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

Theoretical MW_{Fusion Protein}: 90,358 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: 0.093 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

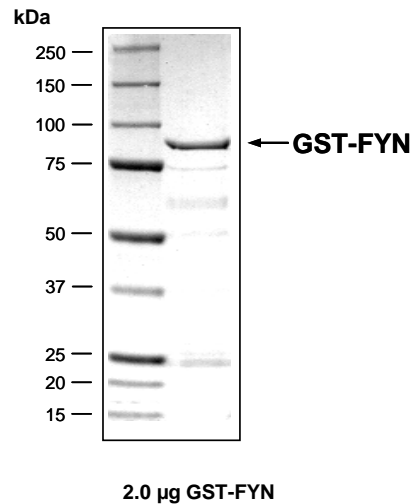
Biochemical Parameters:

Specific kinase activity (P_i transfer): 87 pmol/µg × min
ATP-K_M: 2.4 µM

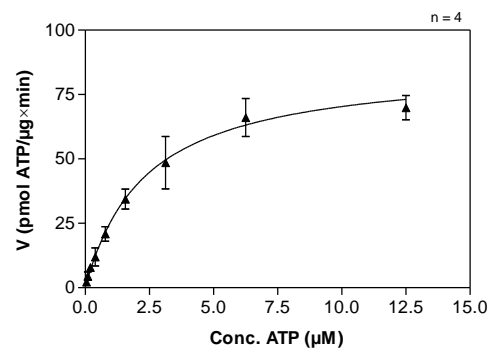
Additional assay technology:

FYN wt Lot 002 was also successfully tested by ProQinase 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

FYN wt Lot 002: Coomassie stain



FYN wt 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: Poly(Glu:Tyr)_{4:1}, 20 µg/ml
Kinase: 2 µg/ml
- Filter binding assay
MSFC membrane (Millipore)

FYN wt

Product No.: 0352-0000-1

GST-FYN wt Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDKVLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMG HHHHHG	RRRASVAAGI	240
241	LVPRGS PGLD	GICSRNS MGC	VQCKDKEATK	LTEERDGS LN	QSSGYRYGTD	PTPQHYP SFG	300
301	VTSIPNYN NF	HAAGQG LTV	FGGVN SSSHT	GTLRTRGGTG	VTLFVALYDY	EARTEDDL SF	360
361	HKGEKFQILN	SSEG DWWEAR	SLTTGETGYI	PSNYVAPVDS	IQAEEWYFGK	LGRKDAERQL	420
421	LSFGNPRGTF	LIRESETTKG	AYSLSIRDWD	DMKGDHV KHY	KIRKLDNGGY	YITTRAQFET	480
481	LQQLVQHYSE	RAAGLCCLV	VPCHKGMPRL	TDLSVKT KD V	WEIPRESLQL	IKRLGNGQFG	540
541	EVVMGTWNGN	TKVAIKTLKP	GTMSPE SFLE	EAQIMK LKH	DKLVQLYAVV	SEEPYIVTE	600
600	YMNGSLLDF	LKDGEGRALK	LPNLVDMAAQ	VAAGMAYIER	MNYIHRDLRS	ANILVGNGLI	660
661	CKIADFG LAR	LIEDNEYTAR	QGAKFPIKWT	APEAALYGRF	TIKSDVWSFG	ILLTELVT KG	720
721	RVPYPGMN NR	EVLEQVERGY	RMPCPQDCPI	SLHELM IHCW	KKDPEERPTF	EYLQSFLE DY	780
781	FTATEPQYQP	GENL					840

1-218: GST **Red**: HIS6-tag **Pink**: Thrombin cleavage site **blue**: FYN

FYN wt ¹ Amino Acid Sequence							
1	MGCVQCKDKE	ATKLTEERDG	SLNQSSGYRY	GTDPTPQHYP	SFGVTSIPNY	NNFHAAGGQG	60
61	LTVFGGVNSS	SHTGTLRTRG	GTGVTLFVAL	YDYEARTEDD	LSFHKGEK FQ	ILNSSEG DWW	120
121	EARSLTTGET	GYIPSNYVAP	VDSIQAE E WY	FGKLGRKDAE	RQLLSFGNPR	GTFLIRESET	180
181	TKGAYLSLIR	DWDDMKGDHV	KHYKIRKLDN	GGYITTRAQ	FETLQQLVQH	YSERAAGLCC	240
241	RLVVPCHKGM	PRLTDL SVKT	KDVWEIPRES	LQLIKRLGNG	QFGEVVMGTW	NGNTKVAIKT	300
301	LKPGTMSPE S	FLEEAQIMKK	LKHDKLVQLY	AVVSEEPYI	VTEYMNKGS L	LDFLKDGEGR	360
361	ALKLPNLVDM	AAQVAAGMAY	IERMNYIHRD	LRSANILVGN	GLICKIADFG	LARLIEDNEY	420
421	TARQGAKFPI	KWTAPEAALY	GRFTIKSDVW	SFGILLTEL V	TKGRVPYPM	NNREVLEQVE	480
481	RGYRMPCPQD	CPISLHELMI	HCWKKDPEER	PTFEYLQSF L	EDYFTATEPQ	YQPGENL	540

blue: kinase sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_002028.1

This product was manufactured at ProQinase in Freiburg, Germany, and is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.