

ProQinase™ LYN

LYN proto-oncogene, Src family tyrosine kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: LYN

Synonyms: JTK8

Product No.: 0358-0000-1

Lot: 001

Description: Human LYN, full length, amino acids M₁-P₅₁₂ (as in [NCBI/Protein](#) entry NP_002341.1), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: LYN Lot 001, was confirmed as LYN by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 88,470 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, 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!

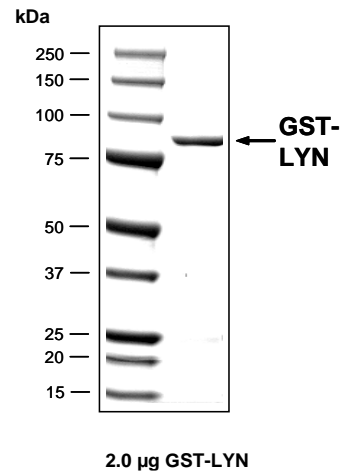
Protein concentration: 0.222 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:
Specific kinase activity (P_i transfer): 94 pmol/µg × min
ATP-K_M: 1.4 µM

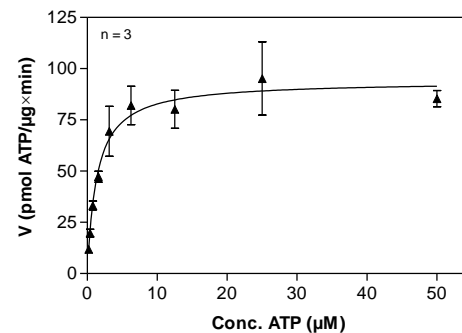
Additional assay technology:

LYN Lot 001 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

**LYN Lot 001:
Coomassie stain**



**LYN Lot 001:
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}, 40 µg/ml
 - LYN: 1 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

ProQinase™ LYN

Product No.: 0358-0000-1

GST-LYN 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	MGCIKSKGKD	SLSDDGVDLK	TQPVRNTER	IYVRDPTS SNK	300
301	QQRVPV ESQL	LPGQRFQTKD	PEEQDIVVA	LYPYDGIHPD	DLSFKKGEKM	KVLEEHGEWW	360
361	KAKSLLTKKE	GFIPSNYVAK	LNTLETEEWF	FKDITRKDAE	RQLLAPGNSA	GAFLIRESET	420
421	LKGSFSLSVR	DFDPVHGDVI	KHYKIRSLDN	GGYIISPRIT	FPCISDMIKH	YQKQADGLCR	480
481	RLEKACISPK	PQKPWDKDAW	EIPRESIKLV	KRLGAGQFGE	VWMGYNNST	KVAVKTLKPG	540
541	TMSVQAFLEE	ANLMKTLQHD	KLVRLYAVVT	REEPIYIITE	YMAKGSLLDF	LKSDEGGKVL	600
601	LPKLIDFSAQ	IAEGMAYIER	KNYIHRDLRA	ANVLVSESLM	CKIADFGLAR	VIEDNEYTAR	660
661	EGAKFPIKWT	APEAINFGCF	TIKSDVWSFG	ILLYEIVTYG	KIPYPGRNTA	DVMTALSQGY	720
721	RMPRVENC PD	ELYDIMKMCW	KEKAERPTF	DYLQSVLDDF	YTATEGQYQQ	QP	780

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

LYN wt ¹ Amino Acid Sequence							
1	MGCIKSKGKD	SLSDDGVDLK	TQPVRNTER	IYVRDPTS SNK	QQRVPV ESQL	LPGQRFQTKD	60
61	PEEQDIVVA	LYPYDGIHPD	DLSFKKGEKM	KVLEEHGEWW	KAKSLLTKKE	GFIPSNYVAK	120
121	LNTLETEEWF	FKDITRKDAE	RQLLAPGNSA	GAFLIRESET	LKGSFSLSVR	DFDPVHGDVI	180
181	KHYKIRSLDN	GGYIISPRIT	FPCISDMIKH	YQKQADGLCR	RLEKACISPK	PQKPWDKDAW	240
241	EIPRESIKLV	KRLGAGQFGE	VWMGYNNST	KVAVKTLKPG	TMSVQAFLEE	ANLMKTLQHD	300
301	KLVRLYAVVT	REEPIYIITE	YMAKGSLLDF	LKSDEGGKVL	LPKLIDFSAQ	IAEGMAYIER	360
361	KNYIHRDLRA	ANVLVSESLM	CKIADFGLAR	VIEDNEYTAR	EGAKFPIKWT	APEAINFGCF	420
421	TIKSDVWSFG	ILLYEIVTYG	KIPYPGRNTA	DVMTALSQGY	RMPRVENC PD	ELYDIMKMCW	480
481	KEKAERPTF	DYLQSVLDDF	YTATEGQYQQ	QP			540

blue: LYN sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_002341.1