

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)

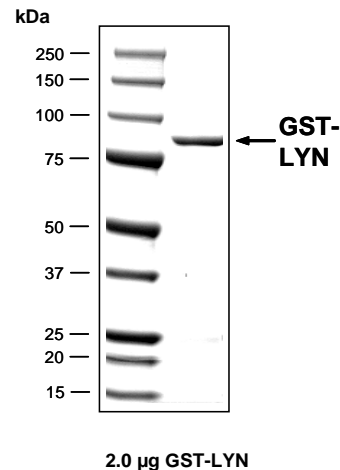
Biochemical Parameters:

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

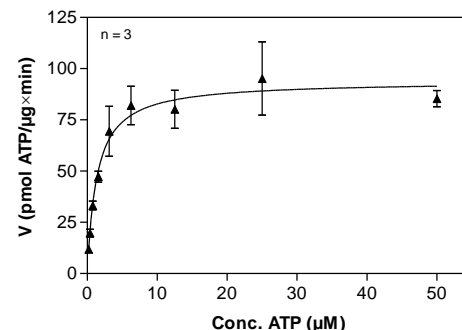
Additional assay technology:

LYN Lot 001 was also successfully tested by 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

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)

LYN

Product No.: 0358-0000-1

GST-LYN Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDKVLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMG HHHHHG	RRRASVAAGI	240
241	LVPRGS PGLD	GIYARGIQAS	MGCIKSKGKD	SLSDDGVDLK	TQPVRNTER	IYVRDPTS SNK	300
301	QQRVPV ESQL	LPGQRFQ TKD	PEEQD IVVA	LYPYD GIHPD	DLSFKK GEKM	KVLEE HGEWW	360
361	KAKSLL TKKE	GFIPS NYVAK	LNTLE TEEF	FKDIT RKDAE	RQLLA PGNSA	GAFLI RESET	420
421	LKGSF SLSVR	DFDPV HGDVI	KHYKIR SLDN	GGYI SPRIT	FPCIS DMIKH	YQQA DGLCR	480
481	RLEKAC ISPK	PQKP WDKAW	EIPRES IKLV	KRLGAG QFGE	VWGM YYNNST	KVAVK TLKPG	540
541	TMSVQA FLEE	ANLMK TLOHD	KLVR LYAVVT	REEPI YIITE	YMAK GSLDF	LKSDE GGKVL	600
601	LPKLID FSAQ	IAEGM AYIER	KNYI HRDLRA	ANVL VSESLM	CKIAD FGLAR	VIEDN EY TAR	660
661	EGAKF PIKWT	APEAI NFGCF	TIKSD VWSFG	ILLYE IVTYG	KIPYP GR TNA	DVMTA LSQGY	720
721	RMPRV ENCPD	ELYDI MKMCW	KEKAE ERTF	DYLSV LDDF	YTATE GQYQQ	QP	780

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

LYN wt ¹ Amino Acid Sequence								
1	MGCIK SKGKD	SLSDD GVDLK	TQPVR NTER	IYVRD PTS	SNK	QQRVP VESQL	LPGQR FQTKD	60
61	PEEQD IVVA	LYPYD GIHPD	DLSFK KGEKM	KVLEE HGEWW	KAKSLL TKKE	GFIPS NYVAK		120
121	LNTLE TEEF	FKDIT RKDAE	RQLLA PGNSA	GAFLI RESET	LKGSF SLSVR	DFDPV HGDVI		180
181	KHYKIR SLDN	GGYI SPRIT	FPCIS DMIKH	YQQA DGLCR	RLEKAC ISPK	PQKP WDKAW		240
241	EIPRES IKLV	KRLGAG QFGE	VWGM YYNNST	KVAVK TLKPG	TMSVQA FLEE	ANLMK TLOHD		300
301	KLVR LYAVVT	REEPI YIITE	YMAK GSLDF	LKSDE GGKVL	LPKLID FSAQ	IAEGM AYIER		360
361	KNYI HRDLRA	ANVL VSESLM	CKIAD FGLAR	VIEDN EY TAR	EGAKF PIKWT	APEAI NFGCF		420
421	TIKSD VWSFG	ILLYE IVTYG	KIPYP GR TNA	DVMTA LSQGY	RMPRV ENCPD	ELYDI MKMCW		480
481	KEKAE ERTF	DYLSV LDDF	YTATE GQYQQ	QP				540

blue: LYN sequence expressed in recombinant protein

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

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.