

ProQinase™ CDK11B 357-795/CycK

cyclin dependent kinase 11B

Recombinant Human Active Protein Kinase

HGNC Symbol: CDK11B

Synonyms: CDC2L1, CDK11-p110, CDK11-p58, CDK11-p46

Product No.: 1779-1484-1

Lot: 002

Description: Human CDK11B/CycK complex, CDK11B C-terminal fragment, amino acids M₃₅₇-F₇₉₅ (as in [NCBI/Protein](#) entry NP_001778.2), CycK N-terminal fragment, amino acids M1-S300 (as in [NCBI/Protein](#) entry NP_001092872.1) both N-terminal GST-HIS₆ fusion proteins with a 3C cleavage site, co-expressed in Sf9 insect cells

Product identity: CDK11B 357-795/CycK Lot 002, was confirmed as CDK11B/CycK by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{CDK11B 357-795}: 78,059 Da

Theoretical MW_{CycK}: 63,031 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: This kinase was not activated by special procedures

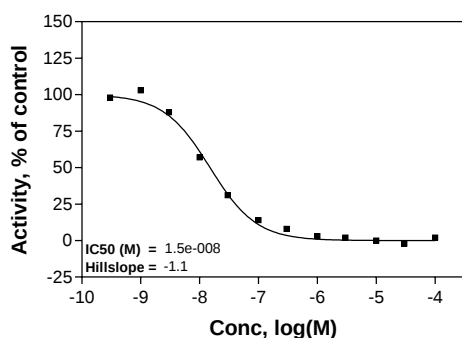
Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20 % glycerol

Storage temperature: -80°C

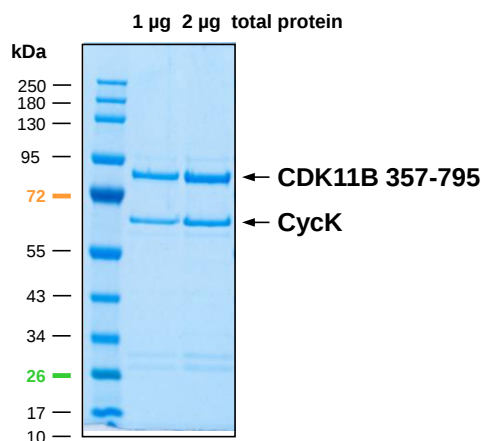
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.524 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

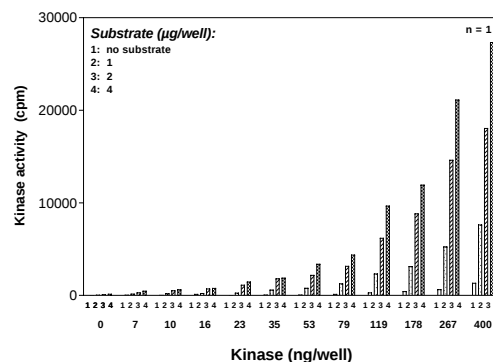
Inhibition of CDK11B 357-795/CycK by AT-7519



CDK11B 357-795/CycK Lot 002: Coomassie stain



CDK11B 357-795/CycK Lot 002: in-vitro 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}
0.1 µM ATP
Substrate: SUPT5H
Kinase: variable
³³PanQinase assay

CDK11B 357-795/CycK

Product No.: 1779-1484-1

GST-CDK11B 357-795 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLIERDEGDK	WRNKKFELGL	EFNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHGG	RDSL LEVLFQ	240
241	PLAMVMS EDE	ERENENHLLV	VPESRFD RDS	GESEEA EEV	GEGTPQSSAL	TEGDYVPD SP	300
301	ALSPIELKQ E	LPKYLPA LQ	CRSVEEFQ CL	NRIEEGTYG V	VYRAKD KKT	EIVALKRL KM	360
361	EKEKEGFPIT	SLREINTILK	AQHPNIVTVR	EIVVGSNMDK	IYIVMNYVEH	DLKSLMETMK	420
421	QPFPLGEVKT	LMIQLLRGVK	HLHDNWILHR	DLKTSNLLLS	HAGILKVGDF	GLAREYGSPL	480
481	KAYTPVVVTL	WYRAPELLLG	AKEYSTAVDM	WSVGCIFGEL	LTQKPLFP GK	SEIDQINKV F	540
541	KDLGTPSEKI	WPGYSELPV	KKMTFSEHPY	NNLRKRFGAL	LSDQGF DL	MNKF LT	600
601	ISAEDGLKHE	YFRETPLPID	PSMFPTWPAK	SEQQRV KRG	TSPRPPEGGLG	YSQLGDDDLK	660
661	ETGFHLTTTN	QGASAAGPGF	SLKF				720

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: CDK11B fragment

CDK11B wt ¹ Amino Acid Sequence							
1	MGDEKDSWKV	KTLDEILQEK	KRRKEQEEKA	EIKRLKNSDD	RDSKRDSLEE	GELRDHRMEI	60
61	TIRNSPYRRE	DSMEDRGEED	DSLAIKPPQQ	MSRKEKAHHR	KDEKRKEKRR	HRSHSAEGGK	120
121	HARVKEKERE	HERRKRHREE	QDKARREW	QKRREMAH	SRRERDRLEQ	LERKRERERK	180
181	MREQQKEQRE	QKERERRAEE	RRKEREARE	VSAHRTMRE	DYSDKVKASH	WSRSPRP	240
241	ERFELGDGRK	PGEARPAPAQ	KPAQLKEEKM	EERDLLSDLQ	DISDSERKTS	SAESSAESG	300
301	SGSEEEEEEE	EEEEEEGSTS	EESEEEEEEE	EEEEETGSN	SEEASEQSAE	EVSEEE MSED	360
361	EERENENHLL	VVPESRFD RD	SGSEEA EEE	VGEGTPQSSA	LTEGDYVPDS	PALSPIELKQ	420
421	ELPKYLPALQ	GCRSVEEFQC	LNRIEETG	VYRAKD KKT	DEIVALKRLK	MEKEKEGFP	480
481	TSLREINTIL	KAQHPNIVTV	REIVVGSNMD	KIYIVMNYVE	HDLKSLMETM	KQPFPLGEVK	540
541	TLMIQLLRGV	KHLHDNWILH	RDLKTSNLLS	SHAGILKVG	FGLAREYGS	LKAYTPVVVT	600
601	LWYRAPELLL	GAKEYSTAVD	MWSVGCIFGE	LLTQKPLFP	KSEIDQINKV	FKDLGTPSEK	660
661	IWPGYSELPA	VKKMTFSEHP	YNNLRKRFGA	LLSDQGF DL	MNKF LT	RISAEDGLKH	720
721	EYFRETPLPI	DPSMFPTWPA	KSEQQRV KRG	TSPRPPEGGL	GYSQLGDDDL	KETGFHLTTT	780
781	NQGASAAGPG	FSLKF					840

blue: CDK11B sequence expressed in recombinant protein

¹NCBI/Protein accession number NP_001778.2

GST-CycK Recombinant Fusion Protein Amino Acid Sequence

1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLIERDEGDK	WRNKKFELGL	EFNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSR	YSKDFETLKV	120
121	DFLSKLPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVLFG	240
241	PMKENKENS	PSVTSANLDH	TKPCWYWDKK	DLAHTPSQLE	GLDPATEARY	RREGARFIFD	300
301	VGTRLGLHYD	TLATGIIYFH	RFYMFHSFKQ	FPRYVTGACC	LFLAGKVEET	PKCKDIIKT	360
361	ARSLNDVQF	QFGDDPKEE	VMVLERILLQ	TIKFDLQVEH	PYQFLLKYAK	QLKGDKNKIQ	420
421	KLVQMAWTFV	NDSLCTTSL	QWEPEIIAVA	VMYLAGRLCK	FEIQEWTSKP	MYRRWWEQFV	480
481	QDVPVDVLED	ICHQILDLYS	QKGQMPHHT	PHQLQPPSL	QPTPQVPQVQ	QSQPSQSSEP	540
541	S						600

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

CycK wt² Amino Acid Sequence

1	MKENKENS	SVTSANLDHT	KPCWYWDKKD	LAHTPSQLEG	LDPATEARYR	REGARFIFDV	60
61	GTRLGLHYD	LATGIIYFHR	FYMFHSFKQF	PRYVTGACCL	FLAGKVEETP	KKCKDIIKTA	120
121	RSLNDVQFG	QFGDDPKEEV	MVLERILLQT	IKFDLQVEHP	YQFLLKYAQ	LKGDKNKIQK	180
181	LVQMAWTFVN	DSLCTTSLQ	WEPEIIAVAV	MYLAGRLCKF	EIQEWTSKPM	YRRWWEQFVQ	240
241	DVPVDVLEDI	CHQILDLYSQ	GKQMPHHTP	HQLQPPSLQ	PTPQVPQVQ	SQPSQSSEPS	300
301	QPQKDPQQP	AQQQPAQQP	KKPSPQSSP	RQVKRAVVVS	PKEENKAAEP	PPPKIPKIET	360
361	THPPLPPAHP	PPDRKPPLAA	ALGEAEPPGP	VDATDLPKVQ	IPPAHPAPV	HQPPPLPHRP	420
421	PPPPSSSYMT	GMSTTSSYMS	GEGYQSLQSM	MKTEGPSYGA	LPPAYGPPAH	LPYHPHYVPP	480
481	NPPPPPVPPP	PASFPPPAIP	PPTPGYPPP	PTYNPNFPPP	PPRLPPTHAV	PPHPPPGI GL	540
541	PPASYPPPAV	PPGGQPPVPP	PIPPPGMPPV	GGLGRAAWMR			600

blue: CycK sequence expressed in recombinant protein

²[NCBI/Protein](#) accession number NP_001092872.1[HGNC](#) identifier CycK: CCNK