

CDK13/CycK

cyclin-dependent kinase 13

Recombinant Human Active Protein Kinase

HGNC Symbol: CDK13

Synonyms: CDC2L, CDC2L5, CHED

Product No.: 1530-1484-1

Lot: 004

Description: Human CDK13, internal fragment amino acids P₆₅₆-G₁₀₅₀ (as in [NCBI/Protein](#) entry NP_003709.3), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site and human CycK, M₁-S₃₀₀ (as in [NCBI/Protein](#) entry NP_001092872.1), N-terminally fused to GST-HIS₆ 3C cleavage site, coexpressed in Sf9 insect cells

Product identity: CDK13/CycK Lot 004, was confirmed as CDK13/CycK by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{GST-CDK13}: 73,831 Da

Theoretical MW_{GST-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: 0.239 µg/µl

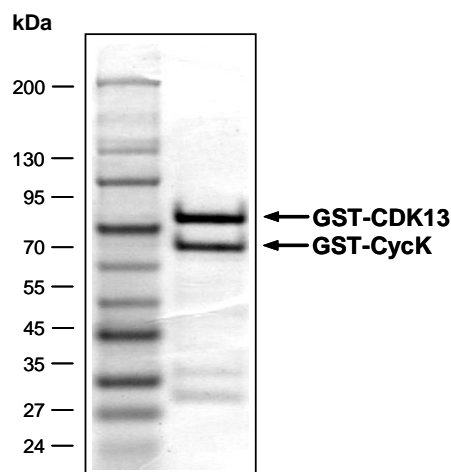
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

Specific kinase activity (P_i transfer): 3 pmol/µg x min

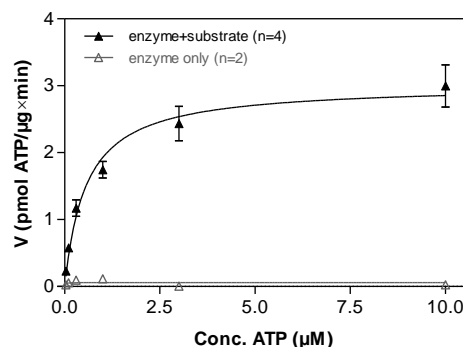
ATP-K_M: 0.56 µM

CDK13/CycK Lot 004:
Coomassie stain



2.0 µg CDK13/CycK

CDK13/CycK Lot 004:
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: RBER-CDC25tide, 40 µg/ml
 - CDK13/CycK: 4 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

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.

CDK13/CycK

Product No.: 1530-1484-1

GST-CDK13 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPE	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVLFCG	240
241	PPGGDDLSKS	PEEKKTATQL	HSKRPRKICG	PRYGETKEKD	IDWGKRCVDK	FDIIGIIGEG	300
301	TYGQVYKARD	KDTGEMVALK	KVRLDNEKEG	FPITAIREIK	ILRQLTHQSI	INMKEIVTDK	360
361	EDALDFKKDK	GAFYLVFEYM	DHDLMLGLES	GLVHFNENHI	KSFMRQLEMG	LDYCHKKNFL	420
421	HRDIKCSNIL	LNNRGQIKLA	DFGLARLYSS	EESRPYTNKV	ITLWYRPPPEL	LLGEERYTPA	480
481	IDVWSCGCIL	GELFTKKPIF	QANQELAQLE	LISRICGSPC	PAVWPDVIKL	PYFNTMKPKK	540
541	QYRRKLREEF	VFIPAAALDL	FDYMLALDPS	KRCTAEQALQ	CEFLRDVEPS	KMPPDPLPLW	600
601	QDCHELWSKK	RRRQKQMGMT	DDVSTIKAPR	KDLSLG			660

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: CDK13 fragment

CDK13 wt ¹ Amino Acid Sequence							
1	MPSSSDTALG	GGGGLSWAEK	KLEERRKRRR	FLSPQQPPLL	LPLLQPQLLQ	PPPPPPPLLF	60
61	LAAPGTAAAA	AAAAAASSSC	FSPGPPLEVK	RLARGKRRAG	GRQKRRRGPR	AGQEAEKRRV	120
121	FSLPQPQQDG	GGGASSGGGV	TPLVEYEDVS	SQSEQGLLLG	GASAATAATA	AGGTGGSGGS	180
181	PASSSGTQRR	GEGSERRPRR	DRRSSSGRSK	ERHREHRRRD	GQRGGSEASK	SRSRHSHSGE	240
241	ERAEVAKSGS	SSSSGGRKRS	ASATSSSSSS	RKDRDSKAHR	SRTKSSKEPP	SAYKEPPKAY	300
301	REDKTEPKAY	RRRSLSPICG	GRDDSPVSHR	ASQSLRSRKS	PSPAGGGSSP	YSRRLPRSPS	360
361	PYSRRRSPSY	SRHSSYERGG	DVSPSPYSSS	SWRRSRSPYS	PVLRRSQKSR	SRSPYSSRHS	420
421	RSRHRHLSLR	SRSRHSSISP	STLTLKSSLA	AELNKNKKAR	AAEAARAAEA	AKAAEATKAA	480
481	EAAAKAAKAS	NTSTPTKGNT	ETSASASQTN	HVKDVKKIKI	EHAPSPSSGG	TLKNDKAKTK	540
541	PPLQVTKVEN	NLIVDKATKK	AVIVGKESKS	AATKEESVSL	KEKTKPLTPS	IGAKEKEQHV	600
601	ALVTSTLPPL	PLPPMLPEDK	EADSLRGNIS	VKAVKKEVEK	KLRCLLADLP	LPPELPGGDD	660
661	LSKSPEEKT	ATQLHSKRRP	KICGPRYGET	KEKDIDWGKR	CVDKFDIIGI	IGEGTYGQVY	720
721	KARDKDTGEM	VALKKVRLDN	EKEGFPITAI	REIKILRQLT	HQSIINMKEI	VTDKEDALDF	780
781	KKDKGAFYLV	FEYMDHDLMG	LLESGLVHFN	ENHIKSFMRQ	LMEGLDYCHK	KNFLHRDIKC	840
841	SNILLNNRGQ	IKLADFGLAR	LYSSEESRPY	TNKVITLWYR	PELLLGEER	YTPAIDVWSC	900
901	GCILGELFTK	KPIFQANQEL	AQLELISRIC	GSPCPAVWPD	VIKLPYFNTM	KPKKQYRRKL	960
961	REEFVFIPAA	ALDLFDYMLA	LDPSKRCTAE	QALQCEFLRD	VEPSKMPPPD	LPLWQDCHEL	1020
1021	WSKRRRQKQ	MGMTDDVSTI	KAPRKDLSLG	LDDSRNTNPQ	GVLPSQLKLS	QGSSNVAPVK	1080
1081	TGPGQHLNHS	ELAILLNLQ	SKTSVNMAF	VQVLNIKVNS	ETQQQLNKIN	LPAGILATGE	1140
1141	KQTDPTPQQ	ESSKPLGGIQ	PSSQTIQPKV	ETDAAQAAVQ	SAFVLLTQL	IKAQQSKQKD	1200
1201	VLLEERENG	GHEASLQLRP	PPEPSTPVSG	QDDLIQHQM	RILELTPEPD	RPRILPPDQR	1260
1261	PPEPPEPPV	TEEDLDYRTE	NQHVPTTSSS	LTDPHAGVKA	ALLQLLAHQH	PQDDPKREGG	1320
1321	IDYQAGDTYV	STSDYKDNFG	SSSFSSAPYV	SNDGLGSSSA	PPLERRSFIG	NSDIQSLDNY	1380
1381	STASSHSGGP	PQPSAFSESF	PSSVAGYGDI	YLNAGPMLFS	GDKDHRFEYS	HGPIAVLANS	1440
1441	SDPSTGPEST	HPLPAKMHNY	NYGGNLQENP	SGPSLMHGQT	WTSPAQGGPY	SQGYRGHIST	1500
1501	STGRGRGRL	PY					1560

blue: CDK13 sequence expressed in recombinant protein

¹NCBI/Protein accession number NP_003709.3

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.

GST-CycK Recombinant Fusion Protein Amino Acid Sequence

1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVLFG	240
241	MKENKENS	PSVTSANLDH	TKPCWYWDKK	DLAHTPSQLE	GLDPATEARY	RREGARFIFD	300
301	VGTRLGLHYD	TLATGIIYFH	RFYMFHSFKQ	FPRYVTGACC	LFLAGKVEET	PKCKDIIKT	360
361	ARSLNDVQF	QFGDDPKEE	VMVLERILLQ	TIKFDLQVEH	PYQFLLYAK	QLKGDKNKIQ	420
421	KLVQMAWTFV	NDSLCTTSL	QWEPEIIAVA	VMYLAGRLCK	FEIQEWTSKP	MYRRWWEQFV	480
481	QDVPVDVLED	ICHQILDLYS	QGKQMPHHT	PHLQPPSL	QPTPQVPQVQ	QSQPSQSEEP	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	LAHTPSQLE	LDPATEARYR	REGARFIFDV	60
61	GTRLGLHYDT	LATGIIYFHR	FYMFHSFKQF	PRYVTGACCL	FLAGKVEETP	KKCKDIIKTA	120
121	RSLNDVQFG	QFGDDPKEEV	MVLERILLQT	IKFDLQVEHP	YQFLLYAKQ	LKGDKNKIQK	180
181	LVQMAWTFVN	DSLCTTSLQ	WEPEIIAVAV	MYLAGRLCKF	EIQEWTSKPM	YRRWWEQFVQ	240
241	DVPVDVLEDI	CHQILDLYSQ	GKQMPHHTP	HQLQPPSLQ	PTPQVPQVQ	SQPSQSSEPS	300
301	QPQQKDPQQP	AQQQQPAQQP	KKPSPQSSP	RQVKRAVVVS	PKEENKAAEP	PPPKIPKIET	360
361	THPPLPPAHP	PPDRKPPLAA	ALGEAEPGP	VDATDLPKVQ	IPPPAHPAPV	HQPPPLPHRP	420
421	PPPPSSYMT	GMSTTSSYMS	GEGYQSLQSM	MKTEGPSYGA	LPPAYGPPAH	LPYHPHYVPP	480
481	NPPPPVPPP	PASFPPPAIP	PPTPGYPPP	PTYNPNFPPP	PPRLPPTHAV	PPHPPGLGL	540
541	PPASYPPPAV	PPGQPPVPP	PIPPGMPPV	GGLGRAAWMR			600

blue: CycK sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_001092872.1

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.