

ProQinase™ CDK1/CycB1

cyclin Dependent Kinase 1

Recombinant Human Active Protein Kinase

HGNC Symbol: CDK1

Synonyms: CDC2, CDC28A

Product No.: 0134-0135-1

Lot: 034

Description: Human CDK1, amino acids M₁-M₂₉₇ (as in [NCBI/Protein](#) entry NP_001777.1) and human CycB1, amino acids M1-V433 (as in [NCBI/Protein](#) entry NP_114172.1), both N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, coexpressed in Sf9 insect cells

Product identity: CDK1/CycB1 Lot 034, was confirmed as CDK1/CycB1 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{GST-CDK1}: 63,882 Da

Theoretical MW_{GST-CycB1}: 78,919 Da

Expression host: Sf9 insect cells

Purification: GST/Immobilized Metal-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

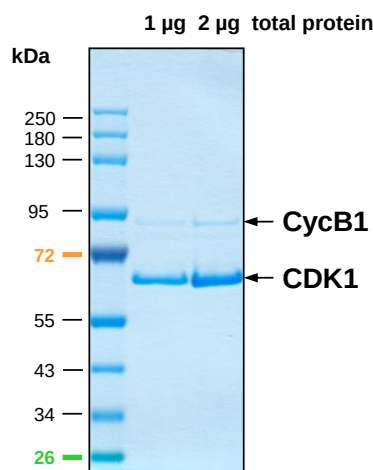
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10 µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.762 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

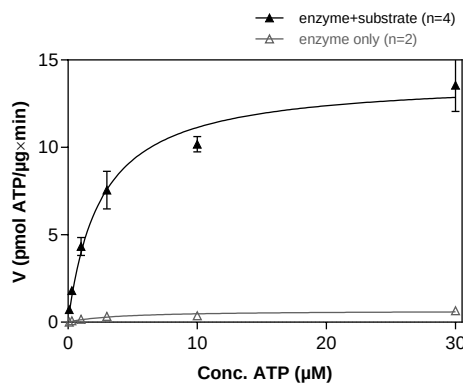
Biochemical Parameters:

Specific kinase activity (P_i transfer): 14 pmol/µg x min
ATP-K_M: 2.5 µM

CDK1/CycB1 Lot 034: Coomassie stain



CDK1/CycB1 Lot 034: 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: RB-CTF 80 µg/ml
Kinase: 4 µg/ml
- Filter binding assay
MSFC membrane (Corning)

CDK1/CycB1

Product No.: 0134-0135-1

GST-CDK1 Recombinant Fusion Protein Amino Acid Sequence

1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLIERDEGDK	WRNKKFELGL	EFNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHGG	RRRASVAAGI	240
241	LVPRGSPGLD	GICSIEEFTM	EDYTKIEKIG	EGTYGVVYKG	RHKTGQVVA	MKKIRLESE	300
301	EGVPSTAI	ISLLKELRHP	NIVSLQDVLM	QDSRLYLIFE	FLSMDLKKYL	DSIPPGQYMD	360
361	SSLVKS	SYLYQ	ILQGI	VFC	HS	RRVLHRDLKP	420
421	HEVVT	LWYRS	PEVLLGSARY	STPVDIWSIG	TIFAELATKK	PLFHGDSEID	480
481	TPNNEVWPEV	ESLQDYKNTF	PKWKPGSLAS	HVKNLDENGL	DLLSKMLIYD	PAKRISGKMA	540
541	LNHPYFNDLD	NQIKKM					600

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue: CDK1 boxed: variation from RefSeq

CDK1 wt¹ Amino Acid Sequence

1	MEDYTKIEKI	GEGTYGVVYK	GRHKTGQV	AMKKIRLESE	EEGVPSTAIR	EISLLKELRH	60
61	PNIVSLQDVL	MQDSRLYLIF	EFLSMDLKKY	LDSIPPGQYM	DSSLVKS	SYLYQ	120
121	SRRLVHRDLK	PQNLLIDDKG	TIKLADFLA	RAFGIPIRVY	THEVVT	LWYR	180
181	YSTPVDIWSI	GTIFAELATK	KPLFHGDSEI	DQLFRIFRAL	GTPNNEVWPE	VESLQDYKNT	240
241	FPKWKPGSLA	SHVKNLDENG	LDLLSKMLIY	DPAKRISGKM	ALNHPYFNDL	DNQIKKM	300

blue: CDK1 sequence expressed in recombinant protein Red: variant in recombinant protein

¹NCBI/Protein accession number NP_001777.1

GST-CycB1 Recombinant Fusion Protein Amino Acid Sequence

1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLIERDEGDK	WRNKKFELGL	EFNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHGG	RRRASVAAGI	240
241	LVPRGSPGLD	GICSIEEFRP	PWELAMALRV	TRNSKINAEN	KAKINMAGAK	RVPTAPAATS	300
301	KPGLRPR	TAL	GDIGNKVSEQ	LQAKMPMKKE	AKPSATGKVI	DKKLKPLEK	360
361	EPVPEPEPEP	EPEPVKEEKL	SPEPILVDTA	SPSPMETSGC	APAEEDLCQA	FSDVILAVND	420
421	VDAEDGADPN	LCSEYVKDIY	AYLRQLEEEQ	AVRPKYLLGR	EVTGNMRAIL	IDWLQVQVMK	480
481	FRLQETMYM	TVSIIDRFMQ	NNCVPKKMLQ	LVGVTAMFIA	SKYEEMYPPE	IGDFAFVTDN	540
541	TYTKHQIRQM	EMKILRALNF	GLGRPLPLHF	LRRASKIGEV	DVEQHTLAKY	LMELTMSDYD	600
601	MVHFPPSQIA	AGAFCLALKI	LDNGEWTPTL	QHYLSYTES	LLPVMQHLAK	NVVMVNQGLT	660
661	KHMTVKNKYA	TSKHAKISTL	PQLNSALVQD	LAKAVAKV			720

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue: CycB1 boxed: variation from RefSeq

CycB1 wt² Amino Acid Sequence

1	MALRVTRNSK	INAENKAKIN	MAGAKRVPTA	PAATSKPGLR	PRTALGDIGN	KVSEQLOAKM	60
61	PMKKEAKPSA	TGKVIDKKLP	KPLEKVPMLV	PVPVSEPVPE	PEPEPEPEPV	KEEKLSP	120
121	LVDTASPSM	ETSGCAPAEE	DLCQAFSDVI	LAVNDVDAED	GADPNLCSEY	VKDIYAYLRQ	180
181	LEEEQAVRPK	YLLGREVTGN	MRAILIDWL	VQVMKFRLLQ	ETMYMTVSI	DRFMQNNCVP	240
241	KKMLQLVGV	T	AMFIASKYEE	MYPPEIGDFA	FVTDNTYTKH	QIRQMEMKIL	300
301	LPLHFLRRAS	KIGEVDVEQH	TLAKYLMELT	MLDYDMVHFP	PSQIAAGAF	LALKILDNGE	360
361	WTPTLQHYLS	YTEESLLPVM	QHLAKNVVMV	NQGLTKHMTV	KNKYATSKHA	KISTLPQLNS	420
421	ALVQDLAKAV	AKV					480

blue: CycB1 sequence expressed in recombinant protein Red: variant in recombinant protein

²NCBI/Protein accession number NP_114172.1