

ProQinase™ CDK1 D146N/CycB1

cyclin Dependent Kinase 1

Recombinant Kinase Substrate (Human Protein Kinase, point mutant, inactive)

HGNC Symbol: CDK1

Synonyms: CDC2, CDC28A

Product No.: 1830-0135-1

Lot: 003

Description: Human CDK1, amino acids M₁-M₂₉₇ (as in [NCBI/Protein](#) entry NP_001777.1) with a point mutation D146N (resulting in loss of kinase activity) and human CycB1, amino acids M1-V433 (as in [NCBI/Protein](#) entry NP_114172.1), both N-terminal GST-HIS₆ fusion proteins with a 3C (CDK1) and Thrombin (CycB1) cleavage site, coexpressed in Sf9 insect cells

Theoretical MW_{GST-CDK1}: 62,997 Da

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

Expression host: Sf9 insect cells

Purification: GST Affinity Chromatography

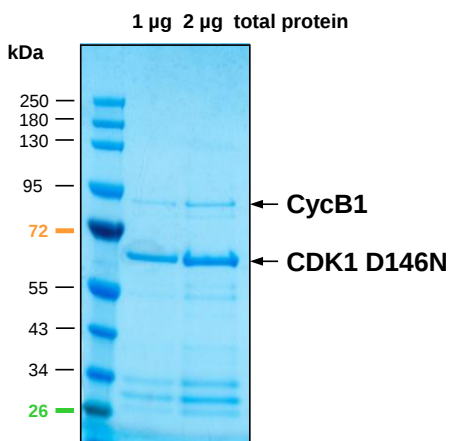
Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20 % glycerol

Storage temperature: -80°C

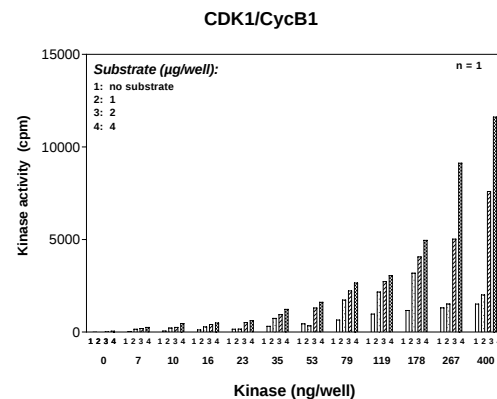
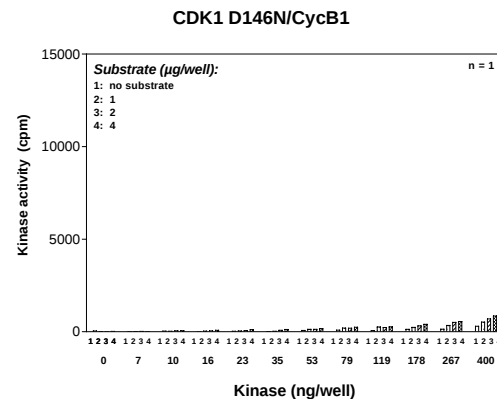
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10 µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 3.093 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

CDK1 D146N/CycB1 Lot 003:
Coomassie stain



Kinase activity CDK1 D146N/CycB1 vs. CDK1/CycB1



- 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-CHKtide (variable amounts)
 - Kinase: variable amounts
 - Assay: ³²PanQinase® Assay

Recombinant Proteins

CDK1 D146N/CycB1

Product No.: 1830-0135-1

GST-CDK1 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	RDSLEVLFGQ	240
241	PLAMGARGRM	EDYTKIEKIG	EGTYGVVYK	RHKTGQVVA	MKKIRLESEE	EGVPSTAI	300
301	ISLLKELRHP	NIVSLQDVLM	QDSRLYLIFE	FLSMDLKKYL	DSIPPGQYMD	SSLVKSILYQ	360
361	ILQGIVFCHS	RRVLHRDLKP	QNLLIDDKGT	IKLANFGLAR	AFGIPIRVYT	HEVVTWLWYRS	420
421	PEVLLGSARY	STPVDIWSIG	TIFAELATKK	PLFHGDSEID	QLFRIFRALG	TPNNEVWPEV	480
481	ESLQDYKNTF	PKWKPGSLAS	HVKNLDENGL	DLLSKMLIYD	PAKRISGKMA	LNHPYFNDLD	540
541	NQIKKM						600

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: CDK1 boxed: D146N

CDK1 wt ¹ Amino Acid Sequence							
1	MEDYTKIEKI	GEGTYGVVYK	GRHKTGQV	AMKKIRLESE	EEGVPSTAIR	EISLLKELRH	60
61	PNIVSLQDVL	MQDSRLYLIF	EFLSMDLKKY	LDSIPPGQYM	DSSLVKSILY	QILQGIVFCH	120
121	SRRLVLRDLK	PQNLLIDDKG	TIKLANFGLA	RAFGIPIRVY	THEVVTWLWYR	SPEVLLGSAR	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	DIRYGVSR	YSKDFETLKV	120
121	DFLSKLPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RRRASVAAGI	240
241	LVPRGSPGLD	GICSIEEFRP	PWELAMALRV	TRNSKINAEN	KAKINMAGAK	RVPTAPAATS	300
301	KPGLRPRAL	GDIGNKVSEQ	LQAKMPMKKE	AKPSATGKVI	DKKLKPLEK	VPMLVPVPS	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	QHYLSYTEES	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 wt ² Amino Acid Sequence							
1	MALRVTRNSK	INAENKAKIN	MAGAKRVPTA	PAATSKPGLR	PRTALGDIGN	KVSEQLQAKM	60
61	PMKKEAKPSA	TGKVIDKKLP	KPLEKVPMLV	PVPVSEPVPE	PEPEPEPEPV	KEEKLSPPEI	120
121	LVDTASPSPM	ETSGCAPAEE	DLCQAFSDVI	LAVNDVDAED	GADPNLCSEY	VKDIYAYLRQ	180
181	LEEEQAVRPK	YLLGREVTGN	MRAILIDWL	VQVMKFRLLQ	ETMYMTVSI	DRFMQNNCVP	240
241	KKMLQLVGV	AMFIASKYEE	MYPPEIGDFA	FVTDNTYTKH	QIRQMEMKIL	RALNFGGLRP	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