

## ProQinase™ CDK2/CycE2 cyclin dependent kinase 2

### Recombinant Human Active Protein Kinase

HGNC Symbol: CDK2

Synonyms: p33

Product No.: 0050-2123-1

Lot: 001

**Description:** Human CDK2, full length, amino acids M<sub>1</sub>-L<sub>298</sub> (as in [NCBI/Protein](#) entry NP\_001789.2), N-terminal GST-HIS<sub>6</sub> fusion protein with a Thrombin cleavage site, and human CycE2, amino acids M<sub>1</sub>-H<sub>404</sub> (as in [NCBI/Protein](#) entry NP\_477097.1), N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site, expressed in Sf9 insect cells.

**Product identity:** CDK2/CycE2 Lot 001, has been verified by mass spectrometry LC-ESI-MS/MS

**Theoretical MW**<sub>GST-CDK2</sub>: 60222 Da  
**Theoretical MW**<sub>GST-CycE2</sub>: 74847 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

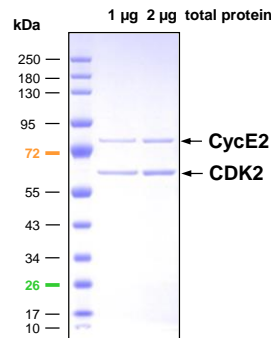
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

**Protein concentration:** 0.583 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

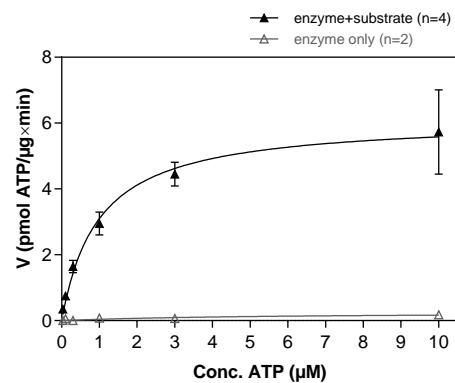
**Biochemical Parameters:**

Specific kinase activity (P<sub>i</sub> transfer): 6.1 pmol/µg\*min  
ATP-K<sub>M</sub>: 1 µM

**CDK2/CycE2 Lot001:  
Coomassie stain**



**CDK2/CycE2 Lot001:  
Determination of V<sub>max</sub> and K<sub>M</sub> value for ATP**



- Assay conditions:
  - 60 mM HEPES-NaOH, pH 7.5
  - 3 mM MgCl<sub>2</sub>
  - 3 mM MnCl<sub>2</sub>
  - 3 µM Na-orthovanadate
  - 1.2 mM DTT
  - 50 µg/ml PEG<sub>20,000</sub>
  - ATP (variable)
  - Substrate: Histone H1 10 µg/ml
  - Kinase: 1 µg/ml

Assay technology:  
Radiometric filter binding assay  
MSFC membrane (96 well plate, Millipore)

Recombinant Proteins

Sequence information

GST-CDK2 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	LVPRGSMENF	QKVEKIGEGT	240
241	YGVVYKARNK	LTGEVVALKK	IRLDTETEGV	PSTAIREISL	LKELNHPNIV	KLLDVIHTEN	300
301	KLYLVFEFLH	QDLKKFMDAS	ALTGIPLPLI	KSYLFQLLQG	LAFCHSHRVL	HRDLKPQNLL	360
361	INTEGAIKLA	DFGLARAFGV	PVRTYTHEVV	TLWYRAPEIL	LGCKYYSTAV	DIWSLGCIFA	420
421	EMVTRRALFP	GDSEIDQLFR	IFRTLGTPE	VVWPGVTSMP	DYKPSFPKWA	RQDFSKVVP	480
481	LDEDGRSLLS	QMLHYDPNKR	ISAKAALAH	FFQDVTKPVP	HLRL		540

1-218: GST **Pink**: Thrombin cleavage site **blue**: CDK2

CDK2 wt <sup>1</sup> Amino Acid Sequence							
1	MENFQVKEKI	GEGTYGVVYK	ARNKLTGEVV	ALKKIRLDT	TEGVPSTAIR	EISLLKELNH	60
61	PNIVKLLDVI	HTENKLYLVF	EFLHQDLKKE	MDASALTGIP	LPLIKSYLFQ	LLQGLAFCHS	120
121	HRVLHRDLKP	QNLLINTEGA	IKLADFLGAR	AFGVPVRTYT	HEVVTLWYRA	PEILLGCKYY	180
181	STAVDIWSLG	CIFAEMVTRR	ALFPGDSEID	QLFRIFRITG	TPDEVVWPGV	TSMPDYKPSF	240
241	PKWARQDFSK	VVPLDEDGR	SLLSQMLHYD	PNKRISAKAA	LAHPFFQDVT	KVPVHLRL	300

**blue**: CDK2 sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](#) accession number NP\_001789.2

GST-CycE2 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	RDSLEVLFGG	240
241	MSRRSSRLQ	AQQPQPSQT	ESQEAQIIQ	AKRKTQDV	KRREEVTKK	HOYEIRNCWP	300
301	PVLSGGISPC	IIETPHKEI	GTSDFSRTN	YRFKNLFINP	SPLPDLWGCS	SKEVWLNMLK	360
361	KESRYVHDKH	FEVLHSDLEP	QMSILLDWL	LEVCEVYTLH	RETFYLAQDF	FDRFMLTQKD	420
421	INKNMLQLIG	ITSLFIASKL	EIYAPKLQE	FAYVTDGACS	EEDILRMELI	ILKALKWELC	480
481	PVTIISWLN	FLQVDALKDA	PKVLLPQYSQ	ETFIQIAQLL	LCILAIKSL	EFQYRILTAA	540
541	ALCHFTSIEV	VKASGLEWD	SISECVDMV	PVNVVKSTS	PVKLKTFFKI	PMEDRHNIQT	600
601	HTNYLAMLEE	VNYINTFRKG	QLSPVCNGG	IMTPPKSTEK	PPGKH		660

1-218: GST **Red**: HIS6-tag **Green**: 3C cleavage site **blue**: CycE2

CycE2 wt <sup>2</sup> Amino Acid Sequence							
1	MSRRSSRLQA	KQQPQPSQTE	SPQEAQIIQA	KRKTQDVK	KRREEVTKKH	QYEIRNCWPP	60
61	VLSGGISPCI	IIETPHKEIG	TSDFSRFTNY	RFKNLFINP	PLPDLWGCS	KEVWLNMLK	120
121	ESRYVHDKHF	EVLHSDLEPQ	MRSILLDWL	EVCEVYTLHR	ETFYLAQDF	DRFMLTQKDI	180
181	NKNMLQLIGI	TSLFIASKLE	EIYAPKLQEF	AYVTDGACSE	EDILRMELII	LKALKWELCP	240
241	VTIISWLNLF	LQVDALKDAP	KVLLPQYSQE	TFIQIAQLLD	LCILAIKSL	FQYRILTAAA	300
301	LCHFTSIEVV	KKASGLEWDS	ISECVDWVMP	FVNVVKSTS	VKLKTFFKI	MEDRHNIQTH	360
361	TNYLAMLEEV	NYINTFRKGG	QLSPVCNGGI	MTPPKSTEK	PGKH		420

**blue**: CycE2 sequence expressed in recombinant protein

<sup>2</sup>[NCBI/Protein](#) accession number NP\_477097.1

[HGNC](#) identifier CycE2: CCNE2