

## CDK6/CycD1

cyclin dependent kinase 6

Recombinant Human Active Protein Kinase

HGNC Symbol: CDK6

Synonyms: PLSTIRE; cell division protein kinase 6

Product No.: 0051-0143-2

Lot: 006

**Description:** Human CDK6, full length, amino acids M<sub>1</sub>-A<sub>326</sub> ([NCBI/Protein](#) entry NP\_001250.1), N-terminally fused to GST-Thrombin cleavage site and human CycD1, full length, amino acids Q<sub>4</sub>-I<sub>295</sub> ([NCBI/Protein](#) entry NP\_444284.1), N-terminally fused to GST-Thrombin cleavage site, expressed in Sf9 insect cells

**Product identity:** CDK6/CycD1 Lot 006, was confirmed as CDK6/CycD1 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>GST-CDK6</sub>:** 63,279 Da

**Theoretical MW<sub>GST-CycD1</sub>:** 59,995 Da

**Expression host:** Sf9 insect cells

**Purification:** GST-Affinity Chromatography

**Activation:** in vitro by CAK1

**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:** 0.388 µg/µl

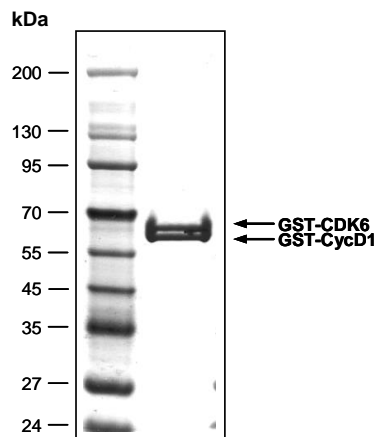
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Biochemical Parameters:**

Specific kinase activity (P<sub>i</sub> transfer): 104 pmol/µg × min

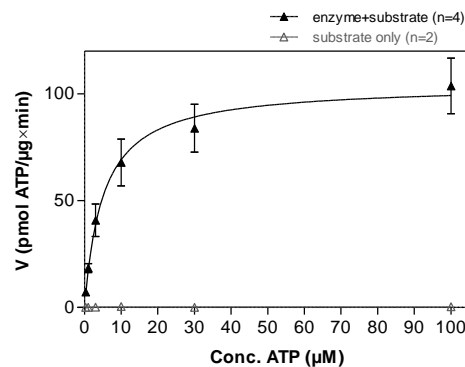
ATP-K<sub>M</sub>: 8.1 µM

### CDK6/CycD1 Lot 006: Coomassie stain



2.0 µg GST-CDK6/GST-CycD1

### CDK6/CycD1 Lot 006: Determination of V<sub>max</sub> and K<sub>M</sub> value for ATP



#### Determination of K<sub>M</sub> value & Specific activity:

- 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: RB-CTF, 100 µg/ml
  - Kinase: 4 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

#### Additional assay technology:

CDK6/CycD1 Lot 006 was also successfully tested by ProQinase for the use with the ADP-Glo™ Kinase assay from Promega. ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

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.

## CDK6/CycD1

Product No.: 0051-0143-2

GST-CDK6 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	LVPRGSM	GLCRADQQYE	240
241	CVAEIGEGAY	GKVF	NGGRFV	VRVQTG	PLSTIREVAV	LRHLETFEHP	300
301	NVRLFDVCT	VSRTDRETKL	TLVFEHVDQD	LTTYLDKVPE	PGVPTETIKD	MMFQLLRGLD	360
361	FLSHR	DLKPQ	SSGQIKLADF	GLARIYSFQM	ALTSVVVTLW	YRAPEVLLQS	420
421	SYATPVDLWS	VGCIFAEMFR	RKPLFRGSSY	VDQLGKILDV	IGLPGEEDWP	RDVALPRQAF	480
481	HSKSAQPIEK	FVTDIDELGK	DLLLKCLTFN	PAKRISAYS	LSPYFQDLE	RCKENLDSHL	540
541	PPSQNTSELN	TA					600

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue: CDK6 boxed: variation from RefSeq

CDK6 wt <sup>1</sup> Amino Acid Sequence								
1	MEKDGLCRAD	QQYECVAEIG	EGAYGKVFKA	RDLKNGGRFV	ALKRVRVQTG	EEGMPLSTIR	60	
61	EVAVLRHLET	FEHPNVVRLF	DVCTVSR	ETKLT	LVFEH	VDQDLTTYLD	KVPEPGVPTE	120
121	TIKDMMFQ	RLDFLHSHR	VVHRDLKPQ	ILVTSSGQIK	LADFLGARIY	SFQMALTSVV	180	
181	VTLWYRAPEV	LLQSSYATPV	DLWSVGCIFA	EMFRRKPLFR	GSSVDQLGK	ILDVIGLPE	240	
241	EDWPRDVALP	RQAFHSKSAQ	PIEKFVTDID	ELGKDLLLKC	LTFNPAKRIS	AYSALSHPHYF	300	
301	QDLRCKENL	DSHLP	PPSQNT	SELNTA			360	

blue: CDK6 sequence expressed in recombinant protein Red: variant in recombinant protein

<sup>1</sup>NCBI/Protein accession number NP\_001250.1

GST-CycD1 Recombinant Fusion Protein Amino Acid Sequence								
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60	
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120	
121	DFLSKLP	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180	
181	KRIEAI	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	LVPRGSAWI	LLCCEVETIR	240	
241	RAYPDANLLN	DRVLRAMLKA	EETCAPSVSY	FKCVQKEVLP	SMRKIVATWM	LEVCEEQKCE	300	
301	EEVFPLAMNY	LDRFLSLEPV	KKSRLQLLGA	TCMFVASKMK	ETIPLTAEKL	CIYTDNSIRP	360	
361	EELLQ	MELLV	VNKLKWNLAA	MTPHDFIEHF	LSKMPEAEEN	KQIRKHAQT	FVALCATDVK	420
421	FISNPPSMVA	AGSVAAVQ	LNLRSPNNFL	SYRRLTRFLS	RVIKCDPDCL	RACQEIQEAL	480	
481	LESSLRQAQ	NMDPKAAEEE	EEEEEEVDLA	CTPTDVRD	V		540	

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue: CycD1

CycD1 wt <sup>2</sup> Amino Acid Sequence									
1	MEHQ	LLCCEV	ETIRRAYPDA	NLLNDRVLR	MLKAAETCAP	SVSYFKCVQK	EVLPSMRKIV	60	
61	ATWML	EVCEE	QKCEEEVFPL	AMNYLDRFLS	LEPVKKSRLQ	LLGATCMFVA	SKMKETIPLT	120	
121	AEKLCIYTDN	SIRPEELLQ	ELLLVNKLKW	NLAAMTPHDF	IEHFLSKMPE	AEENKQIRK		180	
181	HAQTFVALCA	TDVKFISNPP	SMVAAGSVVA	AVOGLNLRSP	NNFLSYRRLT	RFLSRVIKCD		240	
241	PDCLRACQE	Q	IEALLESLR	QAQQNMDPKA	AE	EEEEEE	VDLACTPTDV	RDVDI	300

blue: CycD1 sequence expressed in recombinant protein

<sup>2</sup>NCBI/Protein accession number NP\_444284.1  
HGNC identifier CycD1: CCND1

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.