

## CDK19/CycC

cyclin dependent kinase 19 / cyclin C

Recombinant Human Active Protein Kinase

HGNC Symbol: CDK19

Synonyms: CDC2L6, CDK11

Product No.: 1384-0390-1

Lot: 016

**Description:** Human CDK19, amino acids M<sub>1</sub>-Y<sub>502</sub> (as in [NCBI/Protein](#) entry NP\_055891.1), N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site and human CycC, amino acids M<sub>1</sub>-S<sub>283</sub> (as in [NCBI/Protein](#) entry NP\_005181.2), N-terminally fused to HIS<sub>6</sub>-Thrombin cleavage site, coexpressed in Sf9 insect cells

**Product identity:** CDK19/CycC Lot 016, was confirmed as CDK19/CycC by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>GST-CDK19</sub>** : 84,892 Da

**Theoretical MW<sub>CycC</sub>** : 37,999 Da

**Expression host:** Sf9 insect cells

**Purification:** 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, 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:** 0.186 µg/µl

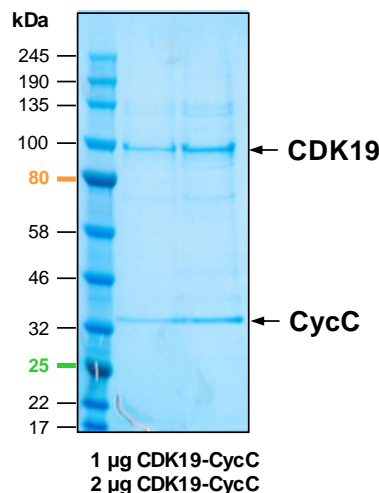
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Biochemical Parameters:**

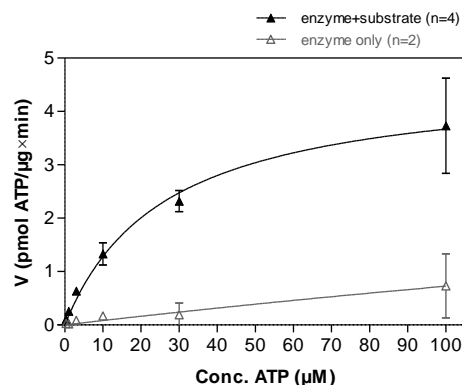
Specific kinase activity (P<sub>i</sub> transfer): 4.6 pmol/µg × min

ATP-K<sub>M</sub>: 26 µM

**CDK19/CycC Lot 016:**  
Coomassie stain



**CDK19/CycC Lot 016:**  
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: RBER-IRStide, 80 µg/ml
  - Kinase: 2 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

## CDK19/CycC

Product No.: 1384-0390-1

GST-CDK19 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI	PQID KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVLFCG	240
241	PMYDFKAKL	AAERERVEL	FEYEGCKVGR	GTYGHVYKAR	RKDGKDEKEY	ALKQIEGTGI	300
301	SMSACREIAL	LRELKHPNVI	ALQKVFLSHS	DRKVLLFDY	AEHDLWHIHK	FHRASKANKK	360
361	PMQLPRSMVK	SILLYQILDGI	HYLHANWVLH	RDLKPANILV	MGEGPERGRV	KIADMGFARL	420
421	FNSPLKPLAD	LDPVVVTFWY	RAPELLLGAR	HYTKAIDIWA	IGCIFAELLT	SEPIFHCROE	480
481	DIKTSNPFHH	DQLDRIFSVM	GFPADKWED	IRKMPEYPTL	QKDFRRTTYA	NSSLIKYMEK	540
541	HKVKPDSKVF	LLLQKLLTMD	PTKRITSEQA	LQDPYFQEDP	LPTLDVFAGC	QIPYPKREFL	600
601	NEDDPEEKGD	KNQQQQNQH	QQPTAPPQQA	AAPPQAPPQ	QNSTQTNGTA	GGAGAGVGGT	660
661	GAGLQHSQDS	SLNQVPPNKK	PRLGPSGANS	GGPVMPSDYQ	HSSRLNYQS	SVQSSSQS	720
721	TLGYSSSSQ	SSQYHPSHQA	HRY				780

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: CDK19

CDK19 wt <sup>1</sup> Amino Acid Sequence							
1	MDYDFKAKLA	AERERVELF	EYEGCKVGRG	TYGHVYKARR	KDGKDEKEYA	LKQIEGTGIS	60
61	MSACREIALL	RELKHPNVIA	LQKVFLSHSD	RKVLLFDYA	EHDLWHIIFK	HRASKANKKP	120
121	MQLPRSMVKS	LLYQILDGIH	YLHANWVLHR	DLKPANILVM	GEGPERGRVK	IADMGFARLF	180
181	NSPLKPLADL	DPVVVTFWYR	APELLLGARH	YTKAIDIWAI	GCIFAELLTS	EPIFHCROED	240
241	IKTSNPFHHD	QLDRIFSVMG	FPADKWEDI	RKMPEYPTLQ	KDFRRTTYAN	SSLIKYMEKH	300
301	KVKPDSKVFL	LLQKLLTMDP	TKRITSEQAL	QDPYFQEDPL	PTLDVFAGCQ	IPYPKREFLN	360
361	EDDPEEKGDK	NQQQQNQHQ	QPTAPPQQA	APPQAPPQ	NSTQTNGTAG	GAGAGVGGTG	420
421	AGLQHSQDSS	LNQVPPNKKP	RLGPSGANS	GPVMPSDYQH	SSRLNYQSS	VQSSSQSQST	480
481	LGYSSSSQS	SQYHPSHQA	RY				540

blue: CDK19 sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](#) accession number NP\_055891.1

HIS-CycC Recombinant Fusion Protein Amino Acid Sequence							
1	MSPIDPMGHH	HHHHGRRRAS	VAAGILVPRG	SPGLDGIYAR	GIQASMAGNF	WQSSHYLQWI	60
61	LDKQDLLKER	QKDLKFLSEE	EYWKLIFFT	NVIQALGEHL	KLRQQVIATA	TVYFKRFYAR	120
121	YSLKSIDPVL	MAPTCVFLAS	KVEEFGVSN	TRLIAAATSV	LKTRFSYAFP	KEFPYRMNHI	180
181	LECEFYLLEL	MDCCLIVYHP	YRPLLQYVQD	MGQEDMLLPL	AWRIVNDTYR	TDLCLLYPPF	240
241	MIALACLHVA	CVVQKQDARQ	WFAELSVDME	KILEIIRVIL	KLYEQWKNFD	ERKEMATILS	300
301	KMPKPKPPPN	SEGEQPNGS	QNSSYSQS				360

Red: HIS6-tag Pink: Thrombin cleavage site blue: CycC

CycC wt <sup>2</sup> Amino Acid Sequence							
1	MAGNFWQSSH	YLQWILDKQD	LLKERQKDLK	FLSEEEYWKL	QIFFTNVIAQ	LGEHLKLRQQ	60
61	VIATATVYFK	RFYARYSLKS	IDPVLMAPTC	VFLASKVEEF	GVSNTRLIA	AATSVLKTRF	120
121	SYAFPKEFPY	RMNHILECEF	YLLELMDCC	IVYHPYRPLL	QYVQDMGQED	MLLPLAWRIV	180
181	NDTYRTDLCL	LYPPFMIALA	CLHVACVVQQ	KDARQWFAEL	SVDMEKILEI	IRVILKLYEQ	240
241	WKNFDERKEM	ATILSKMPKP	KPPPSEGEQ	GPNGSQNSSY	SQS		300

blue: CycC sequence expressed in recombinant protein

<sup>2</sup>[NCBI/Protein](#) accession number NP\_005181.2

[HGNC](#) symbol CycC: CCNC

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.