

## CDK18/CycY

cyclin dependent kinase 18

Recombinant Human Active Protein Kinase

HGNC Symbol: CDK18

Synonyms: PCTAIRE, PCTAIRE3, PCTK3

Product No.: 1546-1619-1

Lot: 008

**Description:** Human CDK18, amino acids M<sub>1</sub>-F<sub>474</sub> (as in [NCBI/Protein](#) entry NP\_997667.1) and human CyclinY, amino acids M<sub>1</sub>-S<sub>341</sub> (as in [NCBI/Protein](#) entry NP\_659449.3), both N-terminal GST-HIS<sub>6</sub> fusion proteins with a 3C cleavage site, coexpressed in Sf9 insect cells.

**Product identity:** CDK18/CycY Lot 008, was confirmed as CDK18/CycY by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>CDK18</sub>** : 82,514 Da

**Theoretical MW<sub>CycY</sub>** : 67,841 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:** 0.335 µg/µl

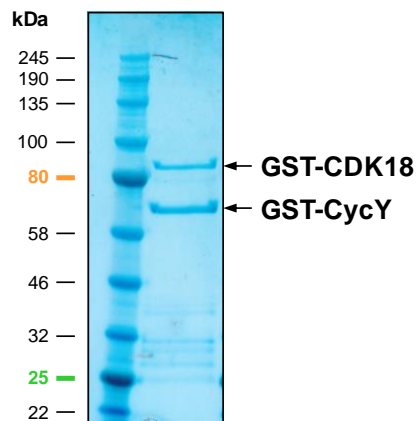
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Biochemical Parameters:**

Specific kinase activity (P<sub>i</sub> transfer): 30 pmol/µg × min

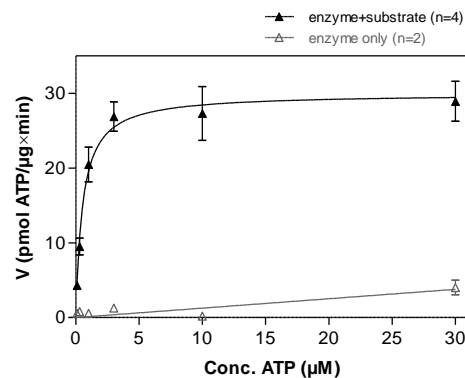
ATP-K<sub>M</sub>: 0.53 µM

**CDK18/CycY Lot 008:**  
Coomassie stain



1 µg GST-CDK18/CycY

**CDK18/CycY Lot 008:**  
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-NTRK3tide 4 µg/ml
  - Kinase: 0.5 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

**Additional assay technology:**

CDK18/CycY Lot 008 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.

## CDK18/CycY

Product No.: 1546-1619-1

GST-CDK18 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMG <b>HHHHHG</b>	RDS <b>LEVLFCG</b>	240
241	<b>PMIMNMKNF</b>	<b>KRRFSLSVPR</b>	<b>TETIEESLAE</b>	<b>FTEQFNQLHN</b>	<b>RRNENLQLGP</b>	<b>LGRDPPQEC</b>	300
301	<b>TFSPDTSGEE</b>	<b>PGQLSPGVQF</b>	<b>QRRQNORRFS</b>	<b>MEDVSKRLSL</b>	<b>PMDIRLPQEF</b>	<b>LQKLQMESPD</b>	360
361	<b>LPKPLSRMSR</b>	<b>RASLSDIGFG</b>	<b>KLETYVKLDK</b>	<b>LGEGTYATVF</b>	<b>KGRSKLTENL</b>	<b>VALKEIRLEH</b>	420
421	<b>EEGAPCTAIR</b>	<b>EVSLKLNKLNK</b>	<b>ANIVTLHDLI</b>	<b>HTDRSLTLVF</b>	<b>EYLDSDLKQY</b>	<b>LDHCGNLMMS</b>	480
481	<b>HNVKIFMFQL</b>	<b>LRGLAYCHHR</b>	<b>KILHRDLKPQ</b>	<b>NLLINERDEL</b>	<b>KLADFLGARA</b>	<b>KSVPTKTYSN</b>	540
541	<b>EVVTLWYRPP</b>	<b>DVLLGSTEYS</b>	<b>TPIDMWGVC</b>	<b>IHYEMATGRP</b>	<b>LFPGSTVKEE</b>	<b>LHLIFRLLGT</b>	600
601	<b>PTEETWPGVT</b>	<b>AFSEFRYTSF</b>	<b>PCYLPQPLIN</b>	<b>HAPRLDTDGI</b>	<b>HLLSSLLLYE</b>	<b>SKSRMSAEAA</b>	660
661	<b>LSHSYFRSLG</b>	<b>ERVHQLEDTA</b>	<b>SIFSLKEIQL</b>	<b>QKDPGYRGLA</b>	<b>FQPGRGKNR</b>	<b>RQSIF</b>	720

1-218: GST   **Red**: HIS6-tag   **Green**: 3C cleavage site   **blue**: CDK18

CDK18 wt <sup>1</sup> Amino Acid Sequence							
1	MIMNMKNFK	RRFSLSVPR	ETIEESLAEF	TEQFNQLHNR	RNENLQLGPL	GRDPPQECST	60
61	FSPTDSGEEP	QQLSPGVQFQ	RRQNORRFSM	EDVSKRLSLP	MDIRLPQEF	QKLQMESPD	120
121	PKPLSRMSR	ASLSDIGFGK	LETYVKLDKL	GEGTYATVFK	GRSKLTENLV	ALKEIRLEHE	180
181	EGAPCTAIR	VSLKLNKLNK	NIVTLHDLIH	HTDRSLTLVF	YLDSDLKQYL	DHCGNLMMSH	240
241	NVKIFMFQQL	RGLAYCHHRK	ILHRDLKPQN	LLINERDELK	LADFLGARA	SVPTKTYSNE	300
301	VVTLWYRPPD	VLLGSTEYST	PIDMWGVC	IHYEMATGRPL	FPGSTVKEEL	HLIFRLLGTP	360
361	TEETWPGVTA	FSEFRYTSFP	CYLPQPLINH	APRLDTDGIH	LLSSLLLYES	KSRMSAEAA	420
421	SHSYFRSLGE	RVHQLEDTAS	IFSLKEIQLQ	KDPGYRGLAF	QPGRGKNR	QSIF	480

**blue**: CDK18 sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](#) accession number NP\_997667.1

GST-CycY Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMG <b>HHHHHG</b>	RDS <b>LEVLFCG</b>	240
241	<b>PLAMVMGNTT</b>	<b>SCCVSSPKL</b>	<b>RRNAHSRLES</b>	<b>YRPDSDLRE</b>	<b>DTGCNLQHS</b>	<b>DRENIDDLNM</b>	300
301	<b>EFNPSDHPRA</b>	<b>STIFLSKSQT</b>	<b>DVREKRKSLF</b>	<b>INHHPPGQIA</b>	<b>RKYSSCSTIF</b>	<b>LDDSTVSQPN</b>	360
361	<b>LKYTIKCVL</b>	<b>AIYYHIKNRD</b>	<b>PDGRMLLDIF</b>	<b>DENLHPLSKS</b>	<b>EVPPDYDKHN</b>	<b>PEQKQIYRFV</b>	420
421	<b>RTLFSAAQLT</b>	<b>AECAIVTLVY</b>	<b>LERLLTYAEI</b>	<b>DICPANWKRI</b>	<b>VLGAILLASK</b>	<b>VWDDQAVWNV</b>	480
481	<b>DYQCILKDI</b>	<b>VEDMNELE</b>	<b>RLLELLFNIN</b>	<b>VPSSVYAKYY</b>	<b>FDLRSLAEAN</b>	<b>NLSFPLEPLS</b>	540
541	<b>RERAKLEAT</b>	<b>SRLCEDKYKD</b>	<b>LRRSARKRSA</b>	<b>SADNLTLP</b>	<b>SPAIIS</b>		600

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

CycY wt <sup>2</sup> Amino Acid Sequence							
1	MGNTTSCCVS	SSPKLRRNAH	SRLESYRPDT	DLSREDTGCN	LQHISDRENI	DDLNMEFNPS	60
61	DHPRASTIFL	SKSQTDVREK	RKSLFINHHP	PGQIARKYSS	CSTIFLDDST	VSQPNLKYTI	120
121	KCVALAIYH	IKNRDPDGRM	LLDIFDENLH	PLSKSEVPPD	YDKHNPEQKQ	IYRFVTRLES	180
181	AAQLTAECAI	VTLYLERLL	TYAEIDICPA	NWKRIVLGAI	LLASKVWDDQ	AVWNVDYQCI	240
241	LKDITVEDMN	ELERQFLELL	QFNINVPSSV	YAKYFDLRS	LAEANLSFP	LEPLSRERAH	300
301	<b>KLEAISRLCE</b>	<b>DKYKDLRSA</b>	<b>RKRSASADNL</b>	<b>TLPRWSPAIIS</b>			360

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

<sup>2</sup>[NCBI/Protein](#) accession number NP\_659449.3  
HGNC identifier CycY: CCNY

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.