

## ProQinase™ CDK9/CycT1 cyclin dependent kinase 9 / CyclinT1

### Recombinant Human Active Protein Kinase

HGNC Symbol: CDK9

Synonyms: C-2k; CDC2L4; PITALRE; TAK; CDC2-related kinase

Product No.: 0371-0345-1

Lot: 014

**Description:** Human CDK9, full length, amino acids M<sub>1</sub>-F<sub>372</sub> (as in [NCBI/Protein](#) entry NP\_001252.1), N-terminal GST-HIS<sub>6</sub> fusion protein with a Thrombin cleavage site, and human CycT1, full length, amino acids M<sub>1</sub>-K<sub>726</sub> (as in [NCBI/Protein](#) entry NP\_001231.2), N-terminal HIS<sub>6</sub> fusion protein with a Thrombin cleavage site, coexpressed in Sf9 insect cells

**Product identity:** CDK9/CycT1 Lot 014, has been verified by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW**<sub>GST-CDK9</sub>: 72,173 Da  
**Theoretical MW**<sub>HIS-CycT1</sub>: 85,441 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

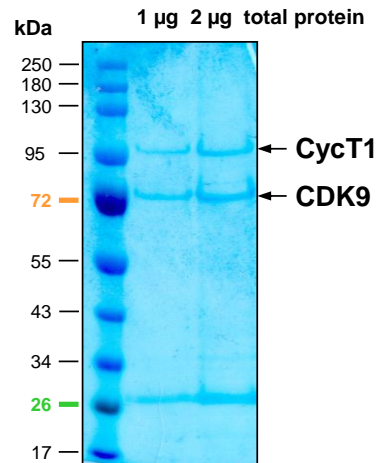
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

**Protein concentration:** 0.102 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

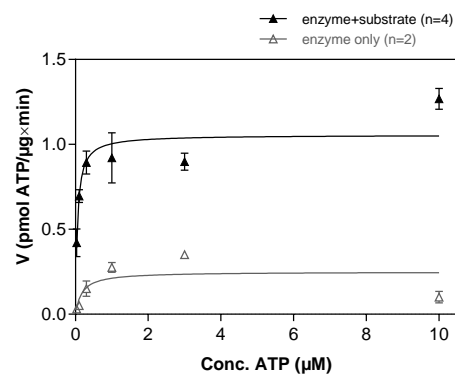
**Biochemical Parameters:**

Specific kinase activity (P<sub>i</sub> transfer): 1.1 pmol/µg x min  
ATP-K<sub>M</sub>: 0.05 µM

CDK9/CycT1 Lot 014:  
Coomassie stain



CDK9/CycT1 Lot 014:  
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: RBER-CHKtide 80 µg/ml  
Kinase: 0.5 µg/ml
- Filter binding assay  
MSFC membrane (Millipore)

Recombinant Proteins

Sequence information

GST-CDK9 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHGG	RRRASVAAGI	240
241	LVPRGSPGLD	GICSRMAKQY	DSVECPFCDE	VSKYEKLAKI	GQGTFGVEFK	ARHRKTGQKV	300
301	ALKKVLME	KEGFPITALR	EIKILQLLKH	ENVVNLIEIC	RTKASPYNRC	KGSIYLVDFD	360
361	CEHDLA	NVLVKFTLSE	IKRVMQMLLN	GLYYIHRNKI	LHRDMKAANV	LITRDGVLKL	420
421	ADFLARAFS	LAKNSQPNRY	TNRVTLWYR	PPELLLGERD	YGPPIDLWGA	GCIMAEMWTR	480
481	SPIMQGNTEQ	HQLALISQLC	GSITPEVWPN	VDNYELYEKL	ELVKGQKRKV	KDRLKAYVRD	540
541	PYALDLIDKL	LVLDPAQRID	SDDALNHDFD	WSDPMPSDLK	GMLSTHLTSM	FEYLAPPRRK	600
601	GSQITQOSTN	QSRNPATTNQ	TEFERVF				660

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue: CDK9

CDK9 wt <sup>1</sup> Amino Acid Sequence							
1	MAKQYDSVEC	PFCDEVSKYE	KLAKIGQGTG	GEVFKARHRK	TGQKVALKKV	LMENEKEGFP	60
61	ITALREIKIL	QLLKHENVVN	LIEICRTKAS	PYNRCKGSIY	LVDFCEHDL	AGLLSNVLVK	120
121	FTLSEIKRVM	QMLLNGLYYI	HRNKILHRDM	KAANVLITRD	GVLKLADFG	ARAFSLAKNS	180
181	QPNRYTNRVV	TLWYRPPPELL	IGERDYGPPI	DLWGAGCIMA	EMWTRSPIMQ	GNTEQHQLAL	240
241	ISQLCGSITP	EVWPNVDNYE	LYEKLELVKG	QKRKVKDRLK	AYVRDPYALD	LIDKLLVLDP	300
301	AQRIDSDDAL	NHDFFWSDPM	PSDLKGMMLST	HLTSMFEYLA	PPRRKGSQIT	QOSTNQSRNP	360
361	ATTNQTEFER	VF					420

blue: CDK9 sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](https://www.ncbi.nlm.nih.gov/protein/NP_001252.1) accession number NP\_001252.1

**HIS-CycT1 Recombinant Fusion Protein Amino Acid Sequence**

1	MSPIDPMGHH HHHHGRRRAS VAAGILVPRG SPGLDGIYAR GIQASMEGER KNNNKRWYFT	60
61	REQLNSPSR RFGVDPDKEL SYRQQAANLL QDMGQRLNVS QLTINTAIVY MHRFYMIQSF	120
121	TQFPGNVAP AALFLAAKVE EQPKKLEHVI KVAHTCLHPQ ESLPDTRSEA YLQQVQDLVI	180
181	LESIILOTLG FELTIDHPHT HVVKCTQLVR ASKDLAQTSY FMATNSLHLT TFSLYQTPPV	240
241	VACVCIHLAC KWSNWEIPVS TDGKHWWEYV DATVTLELLD ELTHEFLQIL EKTPNRLKRI	300
301	WNWRACEAAK KTKADDRGTD EKTSEQTILN MISQSSSDTT IAGLMSMSTS TTSAVPSLPV	360
361	SESSSNLTS VEMPLGKRWL SSQPSFKLEP TQGHRTSENL ALTGVDHSLP QDGSNAFISQ	420
421	KQNSKSVPSA KVSLKEYRAK HAEELAAQKR QLENMEANVK SQYAYAAQNL LSHHDSHSSV	480
481	ILKMPIEGSE NPERPFLEKA DKTALKMRIP VAGGDKAASS KP EEIKMRIK VHAAADKHNS	540
541	VEDSVTKSRE HKEKHKTHPS NHHHHHNNHS HKHSHSQLPV GTGNKRP GPD KHSSQTSNLA	600
601	HKTYLSSSF SSSSTRKRG PSEETGGAVF DHPAKIAKST KSSSLNFSFP SLPTMGQMPG	660
661	HSSDTSGLSF SQPSCKTRVP HSKLDKGP TG ANGHNTTQTI DYQDTVNMLH SLLSAQGVQP	720
721	TQPTAFEFVR PYS DYLNPRS GGISSRSNT KPRPPPLPS EPPPPLPLP K	780

Red: HIS6-tag Pink: Thrombin cleavage site blue: CycT1

**CycT1 wt<sup>2</sup> Amino Acid Sequence**

1	MEGERKNNNK RWFYFTREQLE NSPSRRFGVD PDKELSYRQQ AANLLQDMGQ RLNVSQLTIN	60
61	TAIVYMHRFY MIQSFTQFPG NSVAPAALFL AAKVEEQPKK LEHVIKVAHT CLHPQESLPD	120
121	TRSEAYLQQV QDLVILESII LQTLGFELTI DHPHSHVVKC TQLVRASKDL AQTSYFMATN	180
181	SLHLTTFSLQ YTPPVVACVC IHLACKWSNW EIPVSTDGKH WWEYVDATVT LELLDELTHE	240
241	FLQILEKTPN RLKRIWNWRA CEAAKTKAD DRGTDEKTSE QTILNMISQS SSDTTIAGLM	300
301	SMSTSTTSV PSLPVSEESS SNLTSVEMLP GKRWLSSQPS FKLEPTQGHR TSENALTGV	360
361	DHSLPQDGSN AFISQKQNSK SVPSAKVSLK EYRAKHAEL AAQRQLENM EANVKSQYAY	420
421	AAQNLLSHHD SHSSVILKMP IEGSENPERP FLEKADKTAL KMRI PVAGGD KAASSKPEEI	480
481	KMRIKVHAAA DKHNSVEDSV TKSREHKEKH KTHPSNHHHH HHHSHKHSH SQLPVG TG NK	540
541	RP GPKHSSQ TSNLAHKTY S LSSSFSSSS TRKRG PSEET GGAVFDHPAK IAKSTKSSSL	600
601	NFSFPSLPTM GQMPGHSSDT SGLSFSQPSC KTRVPHSKLD KGPTGANGHN TTQTIDYQDT	660
661	VNMLHLLSA QGVQPTQPTA FEFVRPYSY LNPRSGGISS RSGNTDKPRP PPLPSEPPPP	720
721	LPPLPK	780

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

<sup>2</sup>NCBI/Protein accession number NP\_001231.2  
 HGNC identifier CycT1: CCNT1