

## ProQinase™ CDK12/CycK cyclin dependent kinase 12

### Recombinant Human Active Protein Kinase

HGNC Symbol: CDK12

Synonyms: CRK7, CRKR, CrkRS, CRKRS, hCDK12

Product No.: 1483-1484-1

Lot: 013

**Description:** Human CDK12, amino acids Q696-S1082 (as in [NCBI/Protein](#) entry NP\_057591.2), N-terminal GST-HIS6 fusion protein with a 3C cleavage site and human CycK, amino acids M1-S300 (as in [NCBI/Protein](#) entry NP\_001092872.1), N-terminally fused to GST-HIS6 3C cleavage site, co-expressed in Sf9 insect cells

**Product identity:** CDK12/CycK Lot 013, was confirmed as CDK12/CycK by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW**<sub>GST-CDK12</sub>: 73,066 Da

**Theoretical MW**<sub>GST-CycK</sub>: 63,031 Da

**Expression host:** Sf9 insect cells

**Purification:** GST-Affinity Chromatography

**Activation:** With CAK

**Storage buffer:** 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20 % glycerol

**Storage temperature:** -80°C

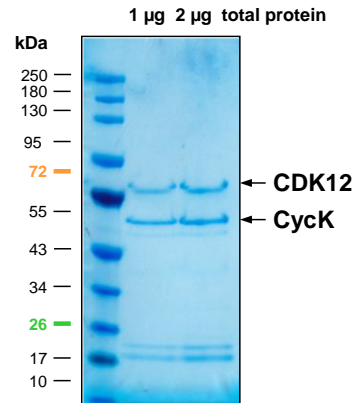
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

**Protein concentration:** 0.244 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

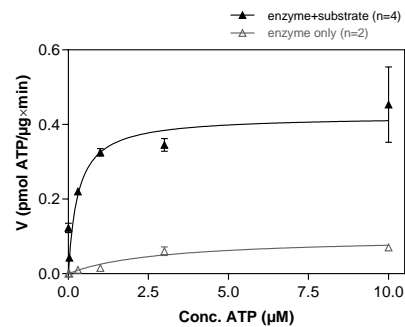
**Biochemical Parameters:**

Specific kinase activity (P<sub>i</sub> transfer): 0.4 pmol/µg x min  
ATP-K<sub>M</sub>: 0.3 µM

**CDK12/CycK Lot 013:  
Coomassie stain**



**CDK12/CycK Lot 013:  
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-IRStide 80 µg/ml  
Kinase: 4 µg/ml
- Filter binding assay  
MSFC membrane (Millipore)

Recombinant Proteins

Sequence information

GST-CDK12 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLIERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEM	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHGG	RDSLEVLFGG	240
241	PQQPYKKRPK	ICCPRYGERR	QTESDWGKRC	VDKFDIIGII	GEGTYGQVYK	AKDKDTGELV	300
301	ALKKVRDNE	KEGFPIAIR	EIKILRQLIH	RSVVMKEIV	TDKQDALDFK	KDKGAFYLVF	360
361	EYMDHDLMLG	LESGLVHFSE	DHIKSFMKQL	MEGLEYPCHK	NFLHRDIKCS	NILLNNSGQI	420
421	KLADFGRLAR	YNSESRPYT	NKVITLWYRP	PELLLGEERY	TPAIDVWSCG	CILGELFTKK	480
481	PIFQANLELA	QLELISRLCG	SPCPAVWPDV	IKLPYFNTMK	PKKQYRRRLR	EEFSFIPSA	540
541	LDLLDHMLTL	DPSKRCTAEQ	TLQSDFLKDV	ELSKMAPDDL	PHWQDCHELW	SKKRRRQRQS	600
601	GVVVEEPPPS	KTSRKETTSG	TSTEPVKNS				660

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: CDK12 fragment

CDK12 wt <sup>1</sup> Amino Acid Sequence							
1	MPNSERHGGK	KDGSGGASGT	LQPSSGGGSS	NSRERHRLVS	KHKRHKSKHS	KDMGLVTPEA	60
61	ASLGTVIKPL	VEYDDISSDS	DTFSDDMAFK	LDRRENDERR	GSDRSDDLHK	HRHHQHRRSR	120
121	DLLKAKQTEK	EKSQEVSSKS	GSMKDRISGS	SKRSNEETDD	YGKAQVAKSS	SKESRSSKLH	180
181	KEKTRKEREL	KSGHKDRSKS	HRKRETPKSY	KTVDSPKRRS	RSPHRKWSDS	SKQDDSPSGA	240
241	SYGQDYDLSP	SRSHTSSNYD	SYKKSPPGTS	RRQSVSPYK	EPSAYQSSTR	SPSPYSRRQR	300
301	SVSPYSRRRS	SSYERSGSYS	GRSPSPYGRR	RSSSPFLSKR	SLSRSPPLSR	KSMKSRSRSP	360
361	AYSRHSSSHS	KKKRSSSRSR	HSSISPVRLP	LNSSLGAELS	RKKKRAAAA	AAAKMDGKES	420
421	KGSPVFLPRK	ENSSVEAKDS	GLESKKLPKS	VKLEKSAPDT	ELVNVTHLNT	EVKNSSDTGK	480
481	VKLDENSEKH	LVKDLKAQGT	RDSKPIALKE	EIVTPKETET	SEKETPPPLP	TIASPPPLP	540
541	TTTPPPQTPP	LPPLPPIPAL	PQQPPLPPSQ	PAFSQVPASS	TSTLPPSTHS	KTSAVSSQAN	600
601	SQPPVQVSVK	TQVSVTAAIP	HLKTSTLPLP	PLPPLLPDGD	DMDSPKETLP	SKPVKKEKEQ	660
661	RTRHLLTDLP	LPPELPGGDL	SPPDSPEPKA	ITPPQPYKK	RPKICCPRYG	ERRQTESDWG	720
721	KRCVDKFDII	GIIGEGTYGQ	VYKAKDKDTG	ELVALKKVRL	DNEKEGFPI	AIREIKILRQ	780
781	LIHRSVVNMK	EIVTDKQDAL	DFKDKGAFY	LVFEYMDHDL	MGLLESLVH	FSEDHIKFSM	840
841	KQLEMEGYC	HKKNFLHRDI	KCSNILLNNS	GQIKLADFLG	ARLYNSEESR	PYTNKVITLW	900
901	YRPELLLGE	ERYTPAIDVW	SCGCILGELF	TKKPIFQANL	ELAQLELISR	LCGSPCPAVW	960
961	PDVIKLPYFN	TMKPKKQYRR	RLREEFSPFIP	SAALDLLDHM	LTLDPKRCRT	AEQTLQSDFL	1020
1021	KDVELSKMAP	PDLPHWQDCH	ELWSKRRRQ	RQSGVVVEEP	PPSKTSRKET	TSGTSTEPVK	1080
1081	NSSPAPPQPA	PGKVESGAGD	AIGLADITQQ	LNQSELAVLL	NLLQSQTDL	IPQMAQLLNI	1140
1141	HSNPEMQQL	EALNQSISAL	TEATSQQQDS	ETMAPEESLK	EAPSAPVILP	SAEQTTLEAS	1200
1201	STPADMQNIL	AVLLSQLMKT	QEPAGSLEEN	NSDKNSGPQG	PRRTPTMPQE	EAAACPPHIL	1260
1261	PPEKRPPEPP	GPPPPPPPPP	LVEGDLSSAP	QELNPAVTAA	LLQLLSQPEA	EPPGHLPEH	1320
1321	QALRPMEYST	RPRPNRTYGN	TDGPETGFSA	IDTDERNNGP	ALTESLVQTL	VKNRTFSGSL	1380
1381	SHLGESSSYQ	GTGSVQFPGD	QDLRFARVPL	ALHPVVGQPF	LKAEGSSNSV	VHAETKLQNY	1440
1441	GELGPGTTGA	SSSGAGLHWG	GPTQSSAYGK	LYRGPTRVPP	RGGRGRGVPY		1500

blue: CDK12 sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](https://www.ncbi.nlm.nih.gov/Protein) accession number NP\_057591.2

**GST-CycK 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	KRIEAIPIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVLFGG	240
241	PMKENKENS	PSVTSANLDH	TKPCWYWDKK	DLAHTPSQLE	GLDPATEARY	RREGARFIFD	300
301	VGTRLGLHYD	TLATGIIYFH	RFYMFHSFKQ	FPRYVTGACC	LFLAGKVEET	PKKCKDIKT	360
361	ARSLNDVQF	QQFGDDPKEE	VMVLERILLQ	TIKFDLQVEH	PYQFLLKYAK	QLKGDKNKIQ	420
421	KLVMAWTFV	NDSLCTTSL	QWEPEIIAVA	VMYLAGRLCK	FEIQEWTSKP	MYRRWWEQFV	480
481	QDVPVDVLED	ICHQILDLYS	QKQQMPHHT	PHLQQPPSL	QPTPQVPQVQ	QSQPSQSSEP	540
541	S						600

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

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

1	MKENKENS	SVTSANLDHT	KPCWYWDK	LAHTPSQLE	LDPATEARY	REGARFIFDV	60
61	GTRLGLHYD	LATGIIYFHR	FYMFHSFKQ	PRYVTGACCL	FLAGKVEET	KKCKDIKTA	120
121	RSLNDVQFG	QFGDDPKEE	MVLERILLQ	IKFDLQVEH	YQFLLKYAK	LKGDKNKIQ	180
181	LVQMAWTFV	DSLCTTSLQ	WEPEIIAVAV	MYLAGRLCKF	EIQEWTSKPM	YRRWWEQFVQ	240
241	DVPVDVLEDI	CHQILDLYS	GKQQMPHHT	HQLQQPPSLQ	PTPQVPQVQ	SQPSQSSEPS	300
301	QFQQKDPQQP	AQQQQPAQQP	KKPSPQPSSP	RQVKRAVVVS	PKEENKAAEP	PPPKIPKIET	360
361	THPPLPPAHP	PPDRKPPLAA	ALGEAEPGP	VDATDLPKVQ	IPPPAPPAV	HQPPPLPHRP	420
421	PPPPSSYMT	GMSTTSSYMS	GEGYQSLQSM	MKTEGPSYGA	LPPAYGPPAH	LPYHPHYVPP	480
481	NPPPPVPPP	PASFPPPAIP	PPTPGYPPP	PTYNPNFPPP	PPRLPPTHAV	PPHPPGLGL	540
541	PPASYPPPAV	PPGGQPPVPP	PIPPGMPPV	GGLGRAAWMR			600

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

<sup>2</sup>[NCBI/Protein](#) accession number NP\_001092872.1  
[HGNC](#) identifier CycK: CCNK