

ProQinase™ CDK9/CycK

Cyclin dependent kinase 9 / CyclinK

Recombinant Human Active Protein Kinase

HGNC Symbol: CDK9

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

Product No.: 0371-1484-1

Lot: 004

Description: Coexpression of human CDK9, full length, amino acids M₁-F₃₇₂ (as in NCBI/Protein entry NP_001252.1), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, and human human CycK, amino acids M₁-S₃₀₀ (as in NCBI/Protein entry NP_001092872.1), N-terminally fused to GST-HIS₆ 3C cleavage site, both expressed in Sf9 insect cells

Product identity: CDK9/CycK Lot 004, was confirmed as CDK9/CycK by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{GST-CDK9}: 72,173 Da

Theoretical MW_{GST-CycK}: 63,031 Da

Expression: Baculovirus infected Sf9 cells

Purification: GST-Affinity Chromatography

Activation: This kinase was activated by coexpression with its physiological cofactor CycK

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: 0.369 µg/µl

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

Biochemical Parameters:

Specific kinase activity (P_i transfer): 13 pmol/µg×min

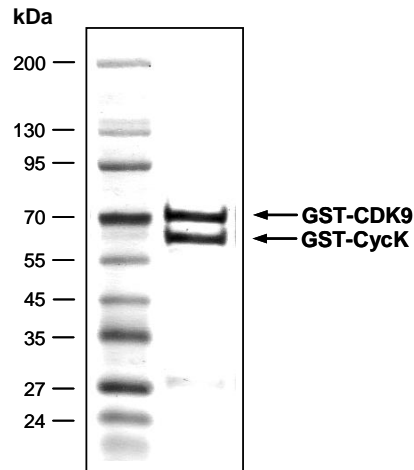
ATP-K_M: 1.3 µM

Additional assay technology: CDK9/CycK Lot 004

was also successfully tested by Reaction Biology for the use with the ADP-Glo™ Kinase assay from ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details



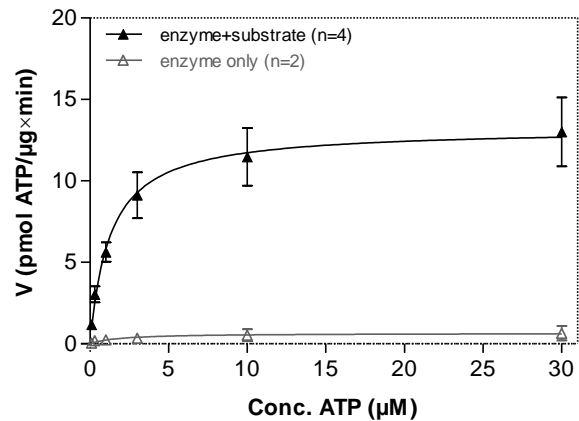
**CDK9/CycK Lot 004:
Coomassie stain**



2.0 µg CDK9/CycK

CDK9/CycK Lot 004:

Determination of V_{max} and K_M value for ATP



Determination of K_M value & Specific activity:

- Assay conditions:
 - 60 mM HEPES-NaOH, pH 7.5
 - 3 mM MgCl₂
 - 3 mM MnCl₂
 - 3 µM Na-orthovanadate
 - 1.2 mM DTT
 - 50 µg / ml PEG_{20,000}
 - ATP (variable)
 - Substrate: RBER-CHKtide, 80 µg/ml
 - CDK9/CycK: 2.0 µg/ml
- Filter binding assay
 - MSFC membrane (Millipore)

Recombinant Proteins

ProQinase™ CDK9/CycK

Product No.: 0371-1484-1

Recombinant Proteins

CDK9 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDKVLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEM	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHGG	RRRASVAAGI	240
241	LVPRGSPGLD	GICSRMAQOY	DSVECPFCDE	VSKYEKLAKI	GQGTFFGEVFK	ARHRKTGQKV	300
301	ALKKVLME	KEGFPITALR	EIKILQLLKH	ENVVNLIEIC	RTKASPYNRC	KGSIIYLVDF	360
361	CEHDLA	NVLVKFTLSE	IKRVMQMLLN	GLYYIHRNKI	LHRDMKAANV	LITRDGVLKL	420
421	ADFLARAFS	LAKNSQPNRY	TNRVVTWYR	PELLLGERD	YGPPIDLWGA	GCIMAEMWTR	480
481	SPIMQGNTEQ	HQLALISQLC	GSITPEVWPN	VDNYELYEKL	ELVKGQKRKV	KDRLKAYVRD	540
541	PYALDLIDKL	LVLDPQRID	SDDALNHDFD	WSDPMPSDLK	GMLSTHLSM	FEYLAPPRK	600
601	GSQITQOSTN	QSRNPATTNQ	TEFERVF				660

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

CDK9 wt ¹ amino acid sequence							
1	MAQYDSVEC	PFCDEVSKYE	KLAKIGQGT	GEVFKARHRK	TGQKVALKKV	LMENEKEGFP	60
61	ITALREIKIL	QLLKHENVVN	LIEICRTKAS	PYNRCKGSIY	LVDFCEHDL	AGLLSNVLVK	120
121	FTLSEIKRVM	QMLLNGLYYI	HRNKILHRDM	KAANVLITRD	GVLKLADFGL	ARAFSLAKNS	180
181	QPNRYTNRV	TLWYRPELL	LGERDYGPPI	DLWGAGCIMA	EMWTRSPIMQ	GNTAQHQLAL	240
241	ISQLCGSITP	EVWPNVDNYE	LYEKLELVKG	QKRKVKDRLK	AYVRDPYALD	LIDKLLVLDP	300
301	AQRIDSDDAL	NHDFFWSDPM	PSDLKGLMST	HLTSMFEYLA	PARRKGSQIT	QQSTNQSRNP	360
361	ATTNQTEFER	VF					420

blue: CDK9 sequence expressed in fusionprotein

¹NCBI/Protein accession number NP_001252.1

CycK Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDKVLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEM	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHGG	RDSLEVLFGG	240
241	PMKENKENS	PSVTSANLDH	TKPCWYWDKK	DLAHTPSQLE	GLDPATEARY	RREGARFIFD	300
301	VGTRLGLHYD	TLATGIIYFH	RFYMFHSFKQ	FPRYVTGACC	LFLAGKVEET	PKCKDKIIT	360
361	ARSLNDVQF	GQFGDDPKEE	VMVLERILLQ	TIKFDLQVEH	PYQFLLKYAK	QLKGDKNKIQ	420
421	KLVQMAWTFV	NDSLCTTSL	QWEPEIIAVA	VMYLAGRLCK	FEIQEWTSKP	MYRRWWEQFV	480
481	QDVPVDVLED	ICHQILDLYS	QKQQMPHHT	PHQLQQPPSL	QPTPQVPQVQ	QSQPSQSSEP	540
541	S						600

Red: HIS6-tag Green: 3C protease cleavage site blue: CycK fragment

CycK wt ² amino acid sequence							
1	MKENKENS	SVTSANLDHT	KPCWYWDKDD	LAHTPSQLEG	LDPATEARYR	REGARFIFDV	60
61	GTRLGLHYDT	LATGIIYFHR	FYMFHSFKQF	PRYVTGACCL	FLAGKVEETP	KKCKDKIKTA	120
121	RSLNDVQFG	QFGDDPKEEV	MVLERILLQT	IKFDLQVEHP	YQFLLKYAKQ	LKGDKNKIQK	180
181	LVQMAWTFVN	DSLCTTSLQ	WEPEIIAVAV	MYLAGRLCKF	EIQEWTSKPM	YRRWWEQFVQ	240
241	DVPVDVLEDI	CHQILDLYSQ	GKQQMPHHTP	HQLQQPPSLQ	PTPQVPQVQQ	SQPSQSSEPS	300
301	QPQQKDPQQP	AQQQQPAQQP	KKPSPQSSP	RQVKRAVVVS	PKEENKAAEP	PPPPIPKIET	360
361	THPPLPAHP	PPDRKPLAA	ALGEAEPGP	VDATDLPKVQ	IPPPAHPAPV	HQPPPLPHRP	420
421	PPPPSSYMT	GMSTTSSYMS	GEGYQSLQSM	MKTEGPSYGA	LPPAYGPPAH	LPYHPHYVPP	480
481	NPPPPVPPP	PASFPPAIP	PPTPGYPPP	PTYNPNFPPP	PPRLPPTHAV	PPHPPGLGL	540
541	PPASYPPPAV	PPGQPPVPP	PIPPPGMPPV	GGLGRAAWMR			600

blue: CycK sequence expressed in fusionprotein

²NCBI/Protein accession number NP_001092872.1

This product was manufactured at Reaction Biology in Freiburg, Germany, and is for in vitro research use only, not for use in humans or animals.
© European Union, 2020. Material may not be reproduced or distributed without written permission from Reaction Biology Europe GmbH