

CDK3/CycC

cyclin dependent kinase 3

Recombinant Human Active Protein Kinase

HGNC Symbol: CDK3

Synonyms: CDKN3

Product No.: 0212-0390-1

Lot: 004

Description: Co-expression of human CDK3, full length, amino acids M₁-H₃₀₅ (as in [NCBI/Protein](#) entry NP_001249.1), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site and human CycC, full length, amino acids M₁-S₂₈₃ (as in [NCBI/Protein](#) entry NP_005181.2), N-terminally fused to HIS₆-Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: CDK3/CycC Lot 004, was confirmed as CDK3/CycC by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{GST-CDK3}: 68,193 Da

Theoretical MW_{HIS-CycC} : 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: 0.182 µg/µl

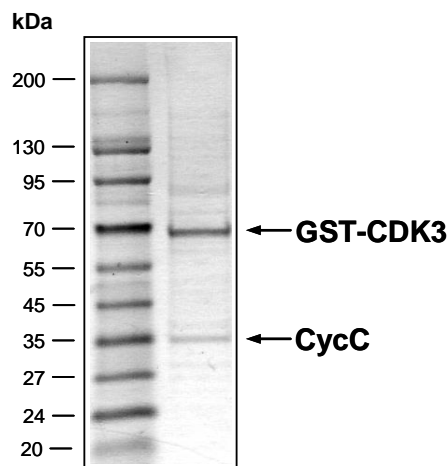
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

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

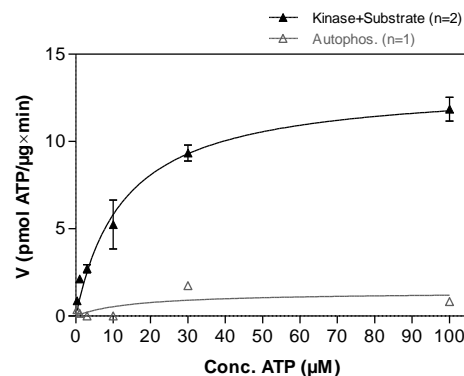
ATP-K_M: 13 µM

CDK3/CycC Lot 004: Coomassie stain



2.0 µg CDK3/CycC

CDK3/CycC Lot 004: Determination of V_{max} and K_M value for ATP



- 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, 40 µg/ml
Kinase: 2 µg/ml
- Filter binding assay
MSFC membrane (Millipore)

Additional assay technology:

CDK3/CycC Lot 004 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.



CDK3/CycC

Product No.: 0212-0390-1

GST-CDK3 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	PMGHHHHHG	RRRASVAAGI	240
241	LVPRGSPGLD	GIYARGIQAS	MGARGRQCDG	YLQNSPLAMD	MFQKVEKIGE	GTYGVVYKAK	300
301	NRETQQLVAL	KKIRLDLEME	GVPSTAIRES	SLLKELKHPN	IVRLLDVVHN	ERKLYLVFEF	360
361	LSQDLKKYMD	STPGSELPLH	LIKSYLFQLL	QGVSFCHSHR	VIHRDLKPQN	LLINELGAIK	420
421	LADFGLARAF	GVPLRITYTHE	VVTLWYRAPE	ILLGSKFYTT	AVDIWSIGCI	FAEMVTRKAL	480
481	FPGDSEIDQL	FRIFRMLGTP	SEDTWPGVTQ	LPDYKGSFPK	WTRKGLEEIV	PNLEPEGRDL	540
541	LMQLLQYDPS	QRITAKTALA	HPYFSSPEPS	PAARQYVLQR	FRHGKGEFQH	TGARY	600

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue: CDK3

CDK3 wt ¹ Amino Acid Sequence							
1	MDMFQVEKI	GEGTYGVVYK	AKNRETQQLV	ALKKIRLDLE	MEGVPSTAIR	EISLLKELKH	60
61	PNIVRLLDV	HNERKLYLVF	EFLSQDLKKY	MDSTPGSELP	LHLIKSYLFQ	LLQGVSFCHS	120
121	HRVIHRDLKP	QNLLINELGA	IKLADFGLAR	AFGVPLRITY	HEVVTWYRA	PEILLGSKFY	180
181	TTAVDIWSIG	CIFAEMVTRK	ALFPGDSEID	QLFRIFRMLG	TPSEDTWPGV	TQLPDYKGSF	240
241	PKWTRKLEE	IVPNLEPEGR	DLLMQLLQYD	PSQRITAKTA	LAHPYFSSPE	PSPAARQYVL	300
301	QRFRRH						360

blue: CDK3 sequence expressed in recombinant protein

¹NCBI/Protein accession number NP_001249.1

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

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

CycC wt ² Amino Acid Sequence							
1	MAGNFWQSSH	YLQWILDKQD	LLKERQKDLK	FLSEEEYWKL	QIFFTNVIQA	LGEHLKLRQQ	60
61	VIATATVYFK	RFYARYSLKS	IDPVLMAPTC	VFLASKVEEF	GVVSNTRLIA	AATSVLKTRF	120
121	SYAFPKEFPY	RMNHILECEF	YLLELMDCCL	IVYHPYRPLL	QYVQDMGQED	MLLPLAWRIV	180
181	NDTYRTDLCL	LYPPFMIALA	CLHVACVVQQ	KDARQWFAEL	SVDMEKILEI	IRVILKLYEQ	240
241	WKNFDERKEM	ATILSKMPKP	KPPPNSEGEQ	GPNGSQNSSY	SQS		300

blue: CycC sequence expressed in recombinant protein

²NCBI/Protein accession number NP_005181.2

HGNC identifier 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.