

ProQinase™ CDK17/p35NCK

Cyclin-dependent kinase 17

Recombinant Human Active Protein Kinase

HGNC Symbol: CDK17

Synonyms: PCTAIRE2, PCTK2

Product No.: 1526-0355-1

Lot: 003

Description: Co-expression of human CDK17, amino acids M₁-F₅₂₃ (as in NCBI/Protein entry NP_002586.2), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site and human p35NCK, amino acids M₁-R₃₀₇ (as in NCBI/Protein entry NP_003876.1), N-terminally fused to GST-Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: CDK17/p35NCK Lot 003, was confirmed as CDK17/p35NCK by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{GST-CDK17}: 87,672 Da

Theoretical MW_{GST-p35NCK}: 63,956 Da

Expression: Baculovirus infected Sf9 cells

Purification: Immobilized Metal Affinity Chromatography

Activation: This kinase was activated by coexpression with its physiological cofactor p35NCK

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

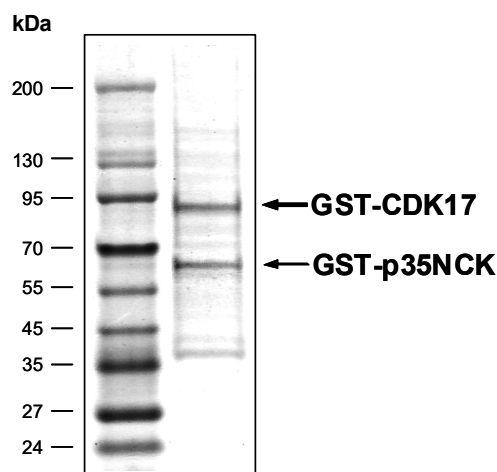
Storage temperature: -80°C
Avoid repeated freeze-thaw cycles!

Protein concentration: 0.286 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:
Specific kinase activity (P_i transfer): 22 pmol/µg×min
ATP-K_M: 15.0 µM

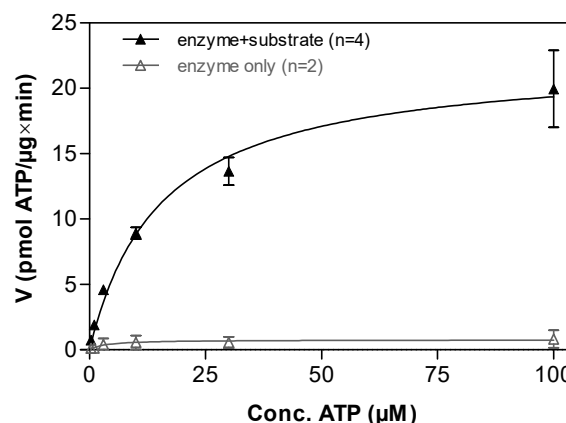
Additional assay technology: CDK17/p35NCK Lot 003 was also successfully tested by Reaction Biology for the use with the ADP-Glo™ Kinase assay from ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

CDK17/p35NCK Lot 003:
Coomassie stain



2.0 µg CDK17/p35NCK

CDK17/p35NCK Lot 003:
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: GSK3-derived peptide (R11-SGRARTSSFAEPGGK), 100 µg / ml
 - CDK17/p35NCK: 4.0 µg / ml
- Filter binding assay
 - MSFC membrane (Millipore)

ProQinase™ CDK17/p35NCK

Product No.: 1526-0355-1

| CDK17 Recombinant Fusion Protein Amino Acid Sequence | | | | | | | |
|--|--------------|------------|-------------|------------|-------------|-------------|-----|
| 1 | MSPILGYWKI | KGLVQPTRL | LEYLEEKYEE | HLYERDEGDK | WRNKKFELGL | EFPNLPYYID | 60 |
| 61 | GDVKLTQ SMA | IIRYIADKHN | MLGGCPKERA | EISMLEGAVL | DIRYGVSRIA | YSKDFETLKV | 120 |
| 121 | DFLSKLP EML | KMFKDRLCHK | TYLNGDHVTH | PDFMLYDALD | VVLYMDPMCL | DAFPKLVCFK | 180 |
| 181 | KRIEAI P QID | KYLKSSKYIA | WPLQGWQATF | GGGDHPPKSD | PMGHHHHHHG | RDSLEVL FQG | 240 |
| 241 | PMKKFKRRLS | LTLRGSQTID | ESLSELAEQM | TIEENSSKDN | EPIVKNGRPP | TSHSMHSFLH | 300 |
| 301 | QYTG SFKKPP | LRRPHSVIGG | SLGSFMAMPR | NGSRLDIVHE | NLKMGS DGES | DQASGTSSDE | 360 |
| 361 | VQSPTGVCLR | NRIHRRISME | DLNKRLSLPA | DIRIPDGYLE | KLQINSPPFD | QPMSRRSRA | 420 |
| 421 | SLSEIGFGKM | ETYIKLEKLG | EGTYATVYKG | RSKLTENLVA | LKEIRLEHEE | GAPCTAIREV | 480 |
| 481 | SLKDLKHAN | IVTLHDIVHT | DKSLTLVFEY | LDKDLKQYMD | DCGNIMSMHN | VKFLYQILR | 540 |
| 541 | GLAYCHRRKV | LHRDLKPQNL | LINEK GELKL | ADFGLARAKS | VPTKTY SNEV | VTLWYRPPDV | 600 |
| 601 | LLGSSEYSTQ | IDMWGVGCIF | FEMASGRPLF | PGSTVEDELH | LIFRLLGTPS | QETWPGISSN | 660 |
| 661 | EEFKNYNFPK | YKPQPLINHA | PRLDSEGIEL | ITKFLQYESK | KRVSAEEAMK | HVYFRSLGPR | 720 |
| 721 | IHALPESVSI | FSLKEIQLQK | DPGFRNSSYP | ETGHGKNRRQ | SMLF | | 780 |

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: CDK17

| CDK17 wt ¹ Amino Acid Sequence | | | | | | | |
|---|-------------|------------|-------------|------------|-------------|------------|-----|
| 1 | MKKFKRRLSL | TLRGSQTIDE | SLSELAEQMT | IEENSSKDNE | PIVKNGRPPT | SHSMHSFLHQ | 60 |
| 61 | YTGSFKK PPL | RRPHSVIGGS | LGSFMAMPRN | GSRLDIVHEN | LKMGS DGEDS | QASGTSSDEV | 120 |
| 121 | QSPTGVCLRN | RIHRRISMED | LNKRLSLPAD | IRIPDGYLEK | LQINSPPFDQ | PMSRRSRRAS | 180 |
| 181 | LSEIGFGKME | TYIKLEKLGE | GTATVYKGR | SKLTENLVAL | KEIRLEHEEG | APCTAIREVS | 240 |
| 241 | LLKDLKHANI | VTLHDIVHTD | KSLTLVFEYL | DKDLKQYMD | CGNIMSMHNV | KLFLYQILRG | 300 |
| 301 | LAYCHRRKVL | HRDLKPQNLL | INEK GELKLA | DFGLARAKSV | PTKTY SNEVV | TLWYRPPDVL | 360 |
| 361 | LGSSEYSTQI | DMWGVGCIFF | EMASGRPLFP | GSTVEDELHL | IFRLLGTPSQ | ETWPGISSNE | 420 |
| 421 | EFKNYNFPKY | KPQPLINHAP | RLDSEGIELI | TKFLQYESKK | RVSAAEEAMKH | VYFRSLGPRI | 480 |
| 481 | HALPESVSIF | SLKEIQLQKD | PGFRNSSYPE | TGHGKNRRQS | MLF | | 540 |

blue: CDK17 sequence expressed in fusionprotein

¹NCBI/Protein accession number NP_002586.2

| p35NCK Recombinant Fusion Protein Amino Acid Sequence | | | | | | | |
|---|--------------|-------------|-------------|------------|------------|-------------|-----|
| 1 | MSPILGYWKI | KGLVQPTRL | LEYLEEKYEE | HLYERDEGDK | WRNKKFELGL | EFPNLPYYID | 60 |
| 61 | GDVKLTQ SMA | IIRYIADKHN | MLGGCPKERA | EISMLEGAVL | DIRYGVSRIA | YSKDFETLKV | 120 |
| 121 | DFLSKLP EML | KMFEDRLCHK | TYLNGDHVTH | PDFMLYDALD | VVLYMDPMCL | DAFPKLVCFK | 180 |
| 181 | KRIEAI P QID | KYLKSSKYIA | WPLQGWQATF | GGGDHPPKSD | PMGHHHHHHG | RRRASVAAGI | 240 |
| 241 | LVPRGSPGLD | GIYARGIQAS | MGTVLSLSPS | YRKATLFEDG | AATVGHYTAV | QNSKNAKDKN | 300 |
| 301 | LKRHSIISVL | PWKRIVAVSA | KKKNSKQVQP | NSSYQNNITH | LNNENLKKSL | SCANLSTFAQ | 360 |
| 361 | PPPAQPPAPP | ASQLSGSQTG | GSSSVK KAPH | PAVTSAGTPK | RVIVQASTSE | LLRCLGEFLC | 420 |
| 421 | RRCYRLK HLS | PTDPVLW LRS | VDRSLLLQGW | QDQGFITPAN | VVFLYMLCRD | VISSEVGS DH | 480 |
| 481 | ELQAVLLTCL | YLSYSYMGNE | ISYPLK PFLV | ESCKEAFWDR | CLSVINLMSS | KMLQINADPH | 540 |
| 541 | YFTQVFS DLK | NESGQEDK KR | LLLGLDR | | | | 600 |

1-218: GST Red: HIS6-tag Pink: Thrombin blue: p35NCK boxed: variation from RefSeq

| p35NCK wt ² Amino Acid Sequence | | | | | | | |
|--|-------------|------------|------------|-------------|-------------|-------------|-----|
| 1 | MGTVLSLSPS | YRKATLFEDG | AATVGHYTAV | QNSKNAKDKN | LKRHSIISVL | PWKRIVAVSA | 60 |
| 61 | KKKNSKQVQP | NSSYQNNITH | LNNENLKKSL | SCANLSTFAQ | PPPAQPPAPP | ASQLSGSQTG | 120 |
| 121 | GSSSVK KAPH | PAVTSAGTPK | RVIVQASTSE | LLRCLGEFLC | RRCYRLK HLS | PTDPVLW LRS | 180 |
| 181 | VDRSLLLQGW | QDQGFITPAN | VVFLYMLCRD | VISSEVGS DH | ELQAVLLTCL | YLSYSYMGNE | 240 |
| 241 | ISYPLK PFLV | ESCKEAFWDR | CLSVINLMSS | KMLQINADPH | YFTQVFS DLK | NESGQEDK KR | 300 |
| 301 | LLLGLDR | | | | | | 360 |

blue: p35NCK sequence expressed in fusionprotein Red: variant in fusionprotein

²NCBI/Protein accession number NP_003876.1
Q193K: SNP variation see NCBI/dbSNP ID: rs17852832