

## ProQinase™ CDK8/CycC

cyclin-dependent kinase 8 / cyclin C

Recombinant Human Active Protein Kinase

HGNC Symbol: CDK8

Synonyms: K35

Product No.: 0376-0390-1

Lot: 004

**Description:** Human CDK8, full length, amino acids M<sub>1</sub>-Y<sub>463</sub> (as in [NCBI/Protein](#) entry XP\_005266272.1), N-terminal GST-HIS<sub>6</sub> fusion protein with a Thrombin cleavage site and human CycC, full length, amino acids M<sub>1</sub>-S<sub>283</sub> (as in [NCBI/Protein](#) entry NP\_005181.2), N-terminally fused to HIS<sub>6</sub>-Thrombin cleavage site, coexpressed in Sf9 insect cells

**Product identity:** CDK8/CycC Lot004 was confirmed as CDK8/CycC by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>GST-CDK8</sub>:** 83,051 Da

**Theoretical MW<sub>His-CycC</sub> :** 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:** 0.262 µg/µl

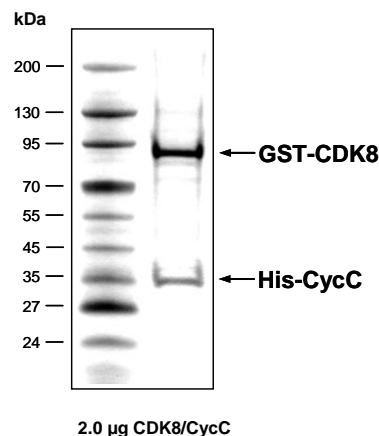
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Biochemical Parameters:**

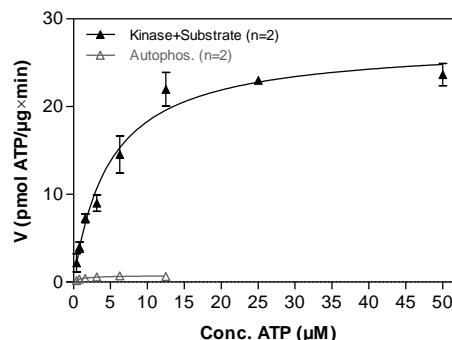
Specific kinase activity (P<sub>i</sub> transfer): 27 pmol/µg x min

ATP-K<sub>M</sub>: 4.9 µM

CDK8/CycC Lot 004:  
Coomassie stain



CDK8/CycC Lot 004:  
Determination of V<sub>max</sub> and K<sub>M</sub> value for ATP



**Determination of K<sub>M</sub> value & Specific activity:**

- 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, 100 µg/ml
  - CDK8/CycC: 2 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

## ProQinase™ CDK8/CycC

Product No.: 0376-0390-1

GST-CDK8 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	PMG <b>HHHHH</b> G	RRRASVAAGI	240
241	<b>LVPRGS</b> PGLD	GIYARGIQAS	<b>MDYDFKVKLS</b>	<b>SERERVEDLF</b>	<b>EYEGCKVGRG</b>	<b>TYGHVYKAKR</b>	300
301	<b>KDGKDDKDYA</b>	<b>LKQIEGTGIS</b>	<b>MSACREIAL</b>	<b>RELKHPNVIS</b>	<b>LQKVFLSHAD</b>	<b>RKVWLLFDYA</b>	360
361	<b>EHDLWHIIF</b>	<b>HRASKANKKP</b>	<b>VQLPRGMVKS</b>	<b>LLYQILDGIH</b>	<b>YLHANWVLR</b>	<b>DLKPANILVM</b>	420
421	<b>GEGPERGRVK</b>	<b>IADMGFARLF</b>	<b>NSPLKPLADL</b>	<b>DPVVVTFWYR</b>	<b>APELLLGARH</b>	<b>YTKAIDIWAI</b>	480
481	<b>GCIFAEELLS</b>	<b>EPIFHCROED</b>	<b>IKTSNPYHHD</b>	<b>QLDRIFNVMG</b>	<b>FPADKDWEDI</b>	<b>KKMPEHSTLM</b>	540
541	<b>KDFRRNTYTN</b>	<b>CSLIKMEKH</b>	<b>KVKPDSKAFH</b>	<b>LLQKLLTMDP</b>	<b>IKRITSEQAM</b>	<b>QDPYFLEDPL</b>	600
601	<b>PTSDVFAGCQ</b>	<b>IPYPKREFLT</b>	<b>EEEPDDKGDGK</b>	<b>NQQQQQGNH</b>	<b>TNGTGHPGNQ</b>	<b>DSSHTQGPPL</b>	660
661	<b>KKVRVVPPTT</b>	<b>TSGGLIMTSD</b>	<b>YQRSNPAAAY</b>	<b>PNPGPSTSQP</b>	<b>QSSMGYSATS</b>	<b>QPPQYSHQT</b>	720
721	<b>HRY</b>						780

1-218: GST **Red**: HIS6-tag **Pink**: Thrombin cleavage site **blue**: CDK8

CDK8 wt <sup>1</sup> Amino Acid Sequence							
1	<b>MDYDFKVKLS</b>	<b>SERERVEDLF</b>	<b>EYEGCKVGRG</b>	<b>TYGHVYKAKR</b>	<b>KDGKDDKDYA</b>	<b>LKQIEGTGIS</b>	60
61	<b>MSACREIAL</b>	<b>RELKHPNVIS</b>	<b>LQKVFLSHAD</b>	<b>RKVWLLFDYA</b>	<b>EHDLWHIIF</b>	<b>HRASKANKKP</b>	120
121	<b>VQLPRGMVKS</b>	<b>LLYQILDGIH</b>	<b>YLHANWVLR</b>	<b>DLKPANILVM</b>	<b>GEGPERGRVK</b>	<b>IADMGFARLF</b>	180
181	<b>NSPLKPLADL</b>	<b>DPVVVTFWYR</b>	<b>APELLLGARH</b>	<b>YTKAIDIWAI</b>	<b>GCIFAEELLS</b>	<b>EPIFHCROED</b>	240
241	<b>IKTSNPYHHD</b>	<b>QLDRIFNVMG</b>	<b>FPADKDWEDI</b>	<b>KKMPEHSTLM</b>	<b>KDFRRNTYTN</b>	<b>CSLIKMEKH</b>	300
301	<b>KVKPDSKAFH</b>	<b>LLQKLLTMDP</b>	<b>IKRITSEQAM</b>	<b>QDPYFLEDPL</b>	<b>PTSDVFAGCQ</b>	<b>IPYPKREFLT</b>	360
361	<b>EEEPDDKGDGK</b>	<b>NQQQQQGNH</b>	<b>TNGTGHPGNQ</b>	<b>DSSHTQGPPL</b>	<b>KKVRVVPPTT</b>	<b>TSGGLIMTSD</b>	420
421	<b>YQRSNPAAAY</b>	<b>PNPGPSTSQP</b>	<b>QSSMGYSATS</b>	<b>QPPQYSHQT</b>	<b>HRY</b>		480

**blue**: CDK8 sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](#) accession number XP\_005266272.1

HIS-CycC Recombinant Fusion Protein Amino Acid Sequence							
1	MSPIDPMG <b>HH</b>	<b>HHHH</b> GRRRAS	VAAGIL <b>LVPRG</b>	<b>SPGLDGIYAR</b>	GIQAS <b>MAGNF</b>	<b>WQSSHYLQWI</b>	60
61	<b>LDKQDLLKER</b>	<b>QKDLKFLSEE</b>	<b>EYWKLOIFFT</b>	<b>NVIQALGEHL</b>	<b>KLRQQVIATA</b>	<b>TVYFKRFYAR</b>	120
121	<b>YSLKSIDPVL</b>	<b>MAPTCVFLAS</b>	<b>KVEEFGVVS</b>	<b>TRLIAAATSV</b>	<b>LKTRFSYAFP</b>	<b>KEFPYRMNHI</b>	180
181	<b>LECEFYLLEL</b>	<b>MDCCLIVYHP</b>	<b>YRPLLQYVQD</b>	<b>MGQEDMLLPL</b>	<b>AWRIVNDTYR</b>	<b>TDLCLLYPPF</b>	240
241	<b>MIALACLHVA</b>	<b>CVVQKQDARQ</b>	<b>WFAELSDVME</b>	<b>KILEIIRVIL</b>	<b>KLYEQWKNFD</b>	<b>ERKEMATILS</b>	300
301	<b>KMPKPKPPPN</b>	<b>SEGEQGPNGS</b>	<b>QNSSYSQS</b>				360

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

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

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

<sup>1</sup>[NCBI/Protein](#) accession number NP\_005181.2

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