

CLK3

CDC-like kinase 3

Recombinant Human Active Protein Kinase

HGNC Symbol: CLK3

Synonyms: PHCLK3

Product No.: 0921-0000-1

Lot: 002

Description: Human CLK3, full length, amino acids M₁-R₄₉₀ (as in [NCBI/Protein](#) entry NP_003983.2), untagged, expressed in Sf9 insect cells

Product identity: CLK3 Lot 002, was confirmed as CLK3 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 59,157 Da

Expression host: Sf9 insect cells

Purification: GST-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: 0.138 µg/µl

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

Biochemical Parameters:

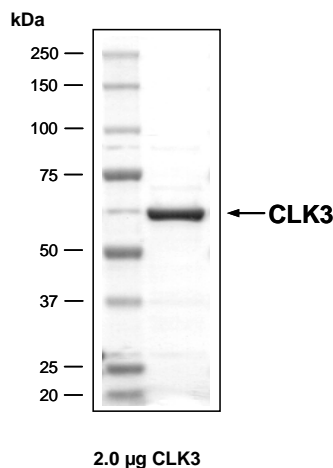
Specific kinase activity (P_i transfer): 76 pmol/µg × min

ATP-K_M: 0.6 µM

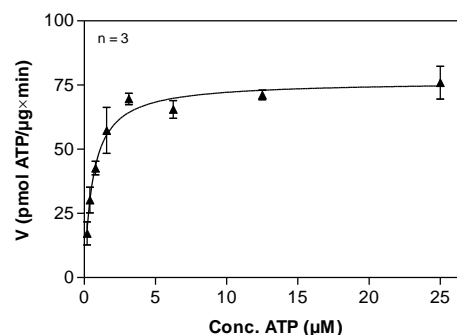
Additional assay technology:

CLK3 Lot 002 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

**CLK3 Lot 002:
Coomassie stain**



**CLK3 Lot 002:
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: RS-peptide 40 µg/ml
 - Kinase: 0.2 µg/ml
- Filter binding assay
MSPH membrane (Millipore)

CLK3

Product No.: 0921-0000-1

| CLK3 Recombinant Fusion Protein Amino Acid Sequence | | | | | | | | |
|---|------------|------------|------------|------------|------------|-------------|------------|-----|
| 1 | GPLAMV | MHHC | KRYRSPEPDP | YLSYRWKRRR | SYSREHEGRL | RYPSRREPPP | RRSRSRSHDR | 60 |
| 61 | LPYQRRYRER | RSDTYRCEE | RSPSFGEDYY | GPSRSRHRRR | SRERGPYRTR | KHAHHCHKRR | | 120 |
| 121 | TRSCSSASSR | SQQSSKRSSR | SVEDDKEGHL | VCRIGDWLQE | RYEIVGNLGE | GTFGKVVECL | | 180 |
| 181 | DHARGKSQVA | LKIIRNVGKY | REARLEINV | LKKIKEKDKE | NKFLCVLMSD | WFNFHGHMCI | | 240 |
| 241 | AFELLGKNTF | EFLKENNFQP | YPLPHVRHMA | YQLCHALRFL | HENQLTHTDL | KPENILFVNS | | 300 |
| 301 | EFETLYNEHK | SCEEKSVKNT | SIRVADFGSA | TFDHEHHTTI | VATRHYRPE | VILELGWAQP | | 360 |
| 361 | CDVWSIGCIL | FEYYRGFTLF | QTHENREHLV | MMEKILGPIP | SHMIHRTRKQ | KYFYKGGVLVW | | 420 |
| 421 | DENSSDGRYV | KENCKPLKSY | MLQDSLEHVQ | LFDLMRRMLE | FDPAQRITLA | EALLHPFFAG | | 480 |
| 481 | LTPEERSFHT | SRNPSR | | | | | | 540 |

1-6: legacy from 3C cleavage **blue**: CLK3

| CLK3 wt ¹ Amino Acid Sequence | | | | | | | |
|--|------------|------------|------------|------------|------------|------------|-----|
| 1 | MHCKRYRSP | EPDPYLSYRW | KRRRSYSREH | EGRLRYPSRR | EPPRRRSR | SHDRLPYQRR | 60 |
| 61 | YRERRSDTY | RCEERSPSFG | EDYYGPSRSR | HRRRSRERGP | YRTRKHAHHC | HKRRTRSCSS | 120 |
| 121 | ASSRSQSSK | RSSRSVEDDK | EGHLVCRIGD | WLQERYEIVG | NLGGFTFGKV | VECLDHARGK | 180 |
| 181 | SQVALKIIRN | VGKYREARL | EINVLKKIKE | KDKENKFLCV | LMSDWFNFHG | HMCIAFELLG | 240 |
| 241 | KNTFEFLKEN | NFQPYPLPHV | RHMAYQLCHA | LRFLHENQLT | HTDLKPENIL | FVNSEFETLY | 300 |
| 301 | NEHKSCEEKS | VKNTSIRVAD | FGSATFDHEH | HTTIVATRHY | RPPEVILELG | WAQPCDVWSI | 360 |
| 361 | GCILFEYYRG | FTLFQTHENR | EHLVMMEKIL | GPIPSHMIHR | TRKQYFYKG | GLVWDENSSD | 420 |
| 421 | GRYVKENCKP | LKSYMLQDSL | EHVQLFDLMR | RMLEFDPAQR | ITLAEALLHP | FFAGLTPEER | 480 |
| 481 | SFHTSRNPSR | | | | | | 540 |

blue: CLK3 sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_003983.2

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.