

CLK1

CDC-like kinase 1

Recombinant Human Active Protein Kinase

HGNC Symbol: CLK1

Synonyms: CLK, CLK4, STY

Product No.: 0447-0000-1

Lot: 001

Description: Human CLK1, full length, amino acids M₁-I₄₈₄ (as in [NCBI/Protein](#) entry NP_004062.2), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: CLK1 Lot 001, was confirmed as CLK1 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 86,695 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: This kinase was not activated by special procedures

Storage buffer: 50 mM TRIS-HCl pH 8.0, 100 mM NaCl, 5 mM DTT, 4 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: 0.139 µg/µl

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

Biochemical Parameters:

Specific kinase activity (P_i transfer): 0.4 pmol/µg × min

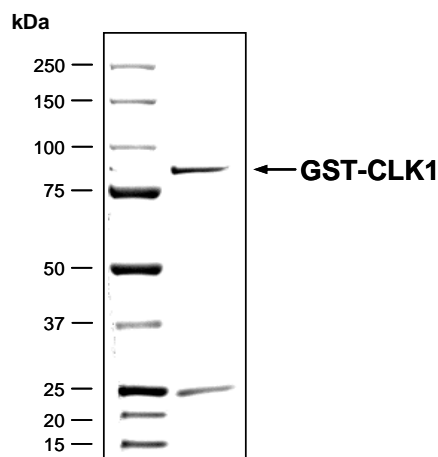
ATP-K_M: 0.2 µM

Additional assay technology:

CLK1 Lot 001 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

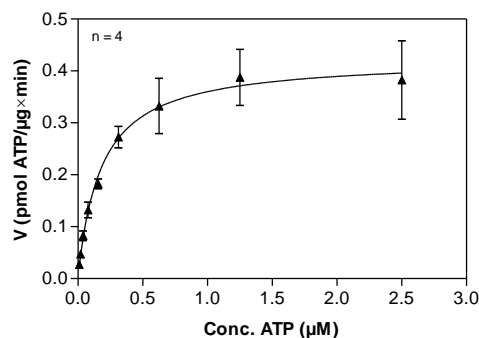
This product was manufactured at ProQinase in Freiburg, Germany, and is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.

CLK1 Lot 001: Coomassie stain



2.0 µg GST-CLK1

CLK1 Lot 001: 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: 1 µg/ml
- Filter binding assay
MSPH membrane (Millipore)

CLK1

Product No.: 0447-0000-1

GST-CLK1 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 HHHHHG	RRRASVAAGI	240
241	LVPRGS PGLD	GICSR MRHSK	RTYCPDWDDK	DWDYGKWRSS	SSHKRRKRSH	SSAHENKRCK	300
301	YNHSMCDSH	YLESRSINEK	DYHSRRYIDE	YRNDYTQCE	PGHRQRDHE	RYQNHSSKSS	360
361	GRSGRSSYKS	KHRIHHSTSH	RRSHGKSHRR	KRTRSVEDDE	EGHLICQSGD	VLSARYEIVD	420
421	TLGEGAFGKV	VECIDHKAGG	RHVAVKIVKN	VDRYCEAARS	EIQVLEHLNT	TDPNSTFRVCV	480
481	QMLEWFEHHG	HICIVFELLG	LSTYDFIKEN	GFLPFRLDHI	RKMAYQICKS	VNFLHSNKLT	540
541	HTDLKPENIL	FVQSDYTEAY	NPKIKRDERT	LINPDIKVD	FGSATYDDEH	HSTLVSTRHY	600
601	RAPEVILALG	WSQPCDVWSI	GCILIEYYLG	FTVFPHTDSK	EHLAMMERIL	GPLPKHMIQK	660
661	TRKRKYFHHD	RLDWDEHSSA	GRYVSRCKP	LKEFMLSQDV	HERLFDLIQ	KMLEYDPAKR	720
721	ITLREALKHP	FFDLLKCSI					780

1-218: GST **Red**: HIS6-tag **Pink**: Thrombin cleavage site **blue**: CLK1 **boxed**: variation from RefSeq

CLK1 wt ¹ Amino Acid Sequence							
1	MRHSKR TYCP	DWDDK DWDYG	KWRSSSS SHKR	RKRSHSSA QE	NKRCKY NHNSK	MCDSHY LESR	60
61	SINEKDY HSR	RYIDEY RNDY	TQCEP GHRO	RDHESRY QNH	SSKSSGR SGR	SSYKSK HRH	120
121	HSTSHRR SHG	KSHRRK RTRS	VEDDEE GHLI	CQSGD VLSAR	YEIVDT LGEG	AFGKV VECID	180
181	HKAGGR HVAV	KIVKNV DRYC	EAARSE IQVL	EHLNTT DPNS	TFRCVQ MLEW	FEHHG HICIV	240
241	FELLGL STYD	FIKENG FLLP	RLDHIR KMAY	QICKSV NFLH	SNKLT HSDLK	PENILF VQSD	300
301	YTEAYN PKIK	RDERTL INPD	IKVVD FGSAT	YDDEH HSTLV	STRHYR APEV	ILALG WSQPC	360
361	DVWSIG CILI	EYYLGF TVFP	THDSK EHLAM	MERILG PPLPK	HMIQK TRKRK	YFHHD RLDWD	420
421	EHSSAG RYVS	RRCKPL KEFM	LSQD VEHERL	FDLIQ KMLEY	DPAKRIT LRE	ALKHP FFDLL	480
481	KKSI						540

blue: CLK1 sequence expressed in recombinant protein **Red**: variant in recombinant protein

¹[NCBI/Protein](#) accession number NP_004062.2