

## CDC7/DBF4

cell division cycle 7

Recombinant Human Active Protein Kinase

HGNC Symbol: CDC7

**Synonyms:** CDC7L1, HsCdc7, HsCDC7, Hsk1, huCdc7

**Product No.:** 1513-1516-1

**Lot:** 006

**Description:** Human CDC7 amino acids M<sub>1</sub>-L<sub>574</sub> (as in [NCBI/Protein](#) entry NP\_003494.1) N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site and DBF4, amino acids M<sub>1</sub>-F<sub>674</sub> (as in [NCBI/Protein](#) entry NP\_006707.1), N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site, coexpressed in Sf9 insect cells

**Product identity:** CDC7/DBF4 Lot 006, was confirmed as CDC7/DBF4 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>GST-CDC7</sub>:** 91,978 Da  
**Theoretical MW<sub>GST-DBF4</sub>:** 104,938 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, 15 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:** 0.197 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

### Biochemical Parameters:

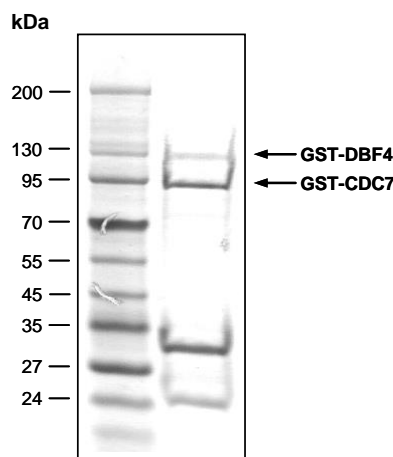
Specific kinase activity (P<sub>i</sub> transfer): 1.8 pmol/µg × min  
ATP-K<sub>M</sub>: 0.04 µM

### Additional assay technology:

CDC7/DBF4 Lot 006 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

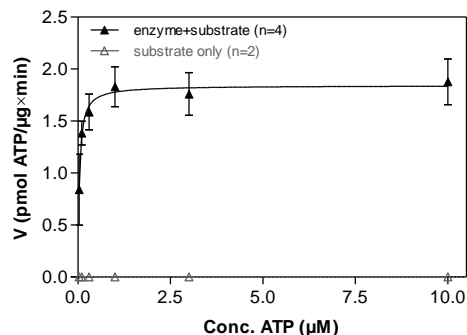
This product was manufactured at ProQinase in Freiburg, Germany, and is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless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

### CDC7/DBF4 Lot 006: Coomassie stain



2.0 µg GST-CDC7/DBF4

### CDC7/DBF4 Lot 006: 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: Histone H1, 20 µg/ml
  - CDC7/DBF4: 2 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

## CDC7/DBF4

Product No.: 1513-1516-1

GST-CDC7 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDKVLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVLFCG	240
241	MEASLGIQM	DEPMAFSPQR	DRFOAEGSLK	KNEQNFKLAG	VKKDIEKLYE	AVPQLSNVFK	300
301	IEDKIGEGTF	SSVYLATAQL	QVGPEEKIAL	KHLIPTSHPI	RIAAELQCLT	VAGGQDNVMG	360
361	VKYCFRKNNDH	VVIAMPYLEH	ESFLDILNSL	SFQEVREYML	NLFKALKRIH	QFGIVHRDVK	420
421	PSNFLYNRRL	KKYALVDFGL	AQGTHDTKIE	LLKFVQSEAO	QERCSONKSH	IITGNKIPLS	480
481	GPVPKELDQQ	STTKASVKRP	YTNAQIQIKQ	GKDGKEGSVG	LSVQRSVFGE	RNFNIHSSIS	540
541	HESPAVKLMK	QSKTVDVLSR	KLATKKKAIS	TKVMNSAVMR	KTASSCPASL	TCDCYATDKV	600
601	CSICLSRRQQ	VAPRAGTPGF	RAPEVLTKCP	NQTTAIDMWS	AGVIFLSLLS	GRYPFYKASD	660
661	DLTALAQIMT	IRGSRETIQA	AKTFGKSILC	SKEVPAQDLR	KLCERLRGMD	SSTPKLTSDI	720
721	QGHASHQPAI	SEKTDHKASC	LVQTPPGQYS	GNSFKKGDSN	SCEHCFFEYN	TNLEGWNEVP	780
781	DEAYDLLDKL	LDLNPASRIT	AEALLHPFF	KDMSL			840

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: CDC7

CDC7 wt <sup>1</sup> Amino Acid Sequence							
1	MEASLGIQMD	EPMAFSPQRD	RFQAEGLSK	NEQNFKLAGV	KKDIEKLYEA	VPQLSNVFKI	60
61	EDKIGEGTFS	SVYLATAQLQ	VGPEEKIALK	HLIPTSHPIR	IAAELQCLTV	AGGQDNVMGV	120
121	KYCFRKNNDHV	VIAMPYLEHE	SFLDILNSLS	FQEVREYMLN	LFKALKRIHQ	FGIVHRDVKP	180
181	SNFLYNRRLK	KYALVDFGLA	QGTHDTKIEL	LKFVQSEAOQ	ERCSONKSHI	ITGNKIPLSG	240
241	PVPKELDQQS	TTKASVKRPY	TNAQIQIKQG	KDGKEGSGVL	SVQRSVFGER	NFNIHSSISH	300
301	ESPAVKLMKQ	SKTVDVLSRK	LATKKKAIST	KVMNSAVMRK	TASSCPASLT	CDCYATDKVC	360
361	SICLSRRQQV	APRAGTPGFR	APEVLTKCPN	QTTAIDMWSA	GVIFLSLLSG	RYPFYKASDD	420
421	LTALAQIMTI	RGSRETIQAA	KTFGKSILCS	KEVPAQDLRK	LCERLRGMDS	STPKLTSDIQ	480
481	GHASHQPAIS	EKTDHKASCL	VQTPPGQYSG	NSFKKGDSNS	CEHCFFEYNT	NLEGWNEVPD	540
541	EAYDLLDKLL	DLNPASRITA	EEALLHPFFK	DMSL			600

blue: CDC7 sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](https://www.ncbi.nlm.nih.gov/protein/NP_003494.1) accession number NP\_003494.1

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.

**GST-DBF4 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	KRIEAIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMG <b>HHHHHG</b>	RDS <b>LEVLFQG</b>	240
241	<b>MNSGAMRIH</b>	<b>SKGHFQGGIQ</b>	<b>VKNEKNRPSL</b>	<b>KSLKTDNRPE</b>	<b>KSKCKPLWGK</b>	<b>VFYLDLPSVT</b>	300
301	<b>ISEKLQKDIK</b>	<b>DLGGRVEEFL</b>	<b>SKDISY LISN</b>	<b>KKEAKFAQTL</b>	<b>GRISPVPSPE</b>	<b>SAYTAETTSP</b>	360
361	<b>HPSHDGSSFK</b>	<b>SPDTVCLSRG</b>	<b>KLLVEKAIKD</b>	<b>HDFIPNSIL</b>	<b>SNALSWG VKI</b>	<b>LHIDDIRYI</b>	420
421	<b>EQKKKELYLL</b>	<b>KKSSTSVRDG</b>	<b>GKRVGSGAQK</b>	<b>TRTGRLKPF</b>	<b>VKVEDMSQLY</b>	<b>RPFYLQLTNM</b>	480
481	<b>PFINYSIQKP</b>	<b>CSPFDVDPKS</b>	<b>SMQKQTQVKL</b>	<b>RIQTDGDKYG</b>	<b>GTSIQLQLKE</b>	<b>KKKKGYCECC</b>	540
541	<b>LQKYEDLETH</b>	<b>LLSEQHRNFA</b>	<b>QSNQYQVDD</b>	<b>IVSKLVDFV</b>	<b>EYEKDTPKKK</b>	<b>RIKYSVGSLS</b>	600
601	<b>PVSASVLKKT</b>	<b>ISQKDCQEDD</b>	<b>TVKQNFY</b>	<b>KETQETKLL</b>	<b>LFISEPI<b>S</b>HP</b>		660
661	<b>SNELRGLNEK</b>	<b>MSNKCSMLST</b>	<b>AEDDIRQNTQ</b>	<b>QLPLHKNQEC</b>	<b>CILDISEHTL</b>	<b>SENDLEELRV</b>	720
721	<b>DHYKCNIQAS</b>	<b>VHVSDFSTDN</b>	<b>SGSQPKQKSD</b>	<b>TVLFPAKDLK</b>	<b>EKDLHSIFTH</b>	<b>DSGLITINSS</b>	780
781	<b>QEHLTVQAKA</b>	<b>PFHTPPEEPN</b>	<b>ECDFKNMDSL</b>	<b>PSGKIHRKVK</b>	<b>IILGRNRKEN</b>	<b>LEPNAEFDKR</b>	840
841	<b>TEFITQEENR</b>	<b>ICSSPVQSLD</b>	<b>DLFQTSEEKS</b>	<b>EFLGFTSYTE</b>	<b>KSGICNVLDI</b>	<b>WEEENSNDLL</b>	900
901	<b>TAFFSSPSTS</b>	<b>TFTGF</b>					960

1-218: GST **Red**: HIS6-tag **Green**: 3C cleavage site **blue**: DBF4 **boxed**: variation from RefSeq

**DBF4 wt<sup>2</sup> Amino Acid Sequence**

1	<b>MNSGAMRIHS</b>	<b>KGHFQGGIQV</b>	<b>KNEKNRPSLK</b>	<b>SLKTDNRPEK</b>	<b>SKCKPLWGKV</b>	<b>FYLDLPSVTI</b>	60
61	<b>SEKLQKDIKD</b>	<b>LGGRVEEFLS</b>	<b>KDISY LISNK</b>	<b>KEAKFAQTLG</b>	<b>RISPVPSPE</b>	<b>AYTAETTSPH</b>	120
121	<b>PSHDGSSFKS</b>	<b>PDTVCLSRGK</b>	<b>LLVEKAIKDH</b>	<b>DFIPNSILS</b>	<b>NALSWG VKIL</b>	<b>HIDDIRYIE</b>	180
181	<b>QKKKELYLLK</b>	<b>KSSTSVRDGG</b>	<b>KRVGSGAQKT</b>	<b>RTGRLKPFV</b>	<b>KVEDMSQLYR</b>	<b>PFY LQLTNMP</b>	240
241	<b>FINYSIQKPC</b>	<b>SPFDVDPKSS</b>	<b>MQKQTQVKLR</b>	<b>IQTDGDKYGG</b>	<b>TSIQLQLKEK</b>	<b>KKKGYCECCL</b>	300
301	<b>QKYEDLETHL</b>	<b>LSEQHRNFAQ</b>	<b>SNQYQVDDI</b>	<b>VSKLVDFVE</b>	<b>YEKDTPKKKR</b>	<b>IKYSVGSLSP</b>	360
361	<b>VSASVLK KTE</b>	<b>QKEKVELQHI</b>	<b>SQKDCQEDDT</b>	<b>TVKQNFYK</b>	<b>ETQETEKLL</b>	<b>FISEPI<b>P</b>HPS</b>	420
421	<b>NELRGLNEKM</b>	<b>SNKCSMLSTA</b>	<b>EDDIRQNTQ</b>	<b>LPLHKNQEC</b>	<b>ILDISEHTLS</b>	<b>ENDLEELRVD</b>	480
481	<b>HYKCNIQASV</b>	<b>HVSDFSTDNS</b>	<b>GSQPKQKSDT</b>	<b>VLFPKDLKE</b>	<b>KDLHSIFTHD</b>	<b>SGLITINSSQ</b>	540
541	<b>EHLTVQAKAP</b>	<b>FHTPPEEPNE</b>	<b>CDFKNMDSL</b>	<b>SGKIHRKVKI</b>	<b>IILGRNRKENL</b>	<b>EPNAEFDKRT</b>	600
601	<b>EFITQEENRI</b>	<b>CSSPVQSLLD</b>	<b>LFQTSEEKSE</b>	<b>FLGFTSYTEK</b>	<b>SGICNVLDIW</b>	<b>EEENSNDLLT</b>	660
661	<b>AFFSSPSTST</b>	<b>FTGF</b>					720

**blue**: DBF4 sequence expressed in recombinant protein **Red**: variant in recombinant protein

<sup>2</sup>[NCBI/Protein](https://www.ncbi.nlm.nih.gov/protein/NP_006707.1) accession number NP\_006707.1