

ProQinase™ CDK15/CycB1 cyclin dependent kinase 15

Recombinant Human Active Protein Kinase

HGNC Symbol: CDK15

Synonyms: PFTAIRE2; ALS2CR7; PFTK2

Product No.: 1889-0135-1

Lot: 002

Description: Human CDK15, internal fragment, amino acids Y₁₀₃-F₃₈₇ (as in [NCBI/Protein](#) entry NP_001248364.1), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site and human cyclin B1, full length, amino acids M₁-V₄₃₃ (as in [NCBI/Protein](#) entry NP_114172.4), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, co-expressed in Sf9 insect cells

Product identity: CDK15/CycB1 Lot 002, was confirmed as CDK15/CycB1 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW _{CDK15} :	60,389	Da
Theoretical MW _{CycB1} :	78,919	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

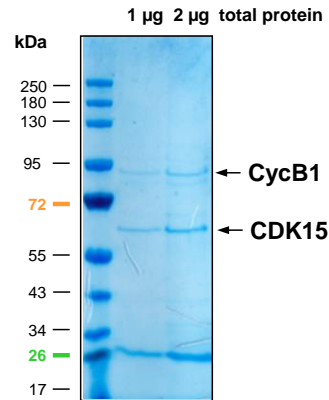
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.156 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

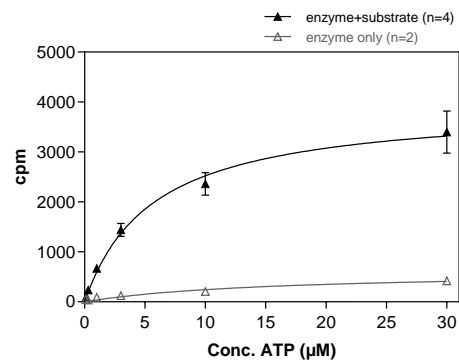
Biochemical Parameters:

Specific kinase activity (P_i transfer): 1.9 pmol/µg x min
ATP-K_M: 2.8 µM

CDK15/CycB1 Lot 002:
Coomassie stain



CDK15/CycB1 Lot 002:
Determination of V_{max} and K_M value for ATP



- 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: RBER-NTRK3tide 80 µg/ml
 - Kinase: 4 µg/ml
- Filter binding assay
MSFC membrane (Millipore)

Recombinant Proteins

Sequence information

GST-CDK15 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEM	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHGG	RDSLEVLFGG	240
241	PYLNLEKLGE	GSYATVYKGI	SRINGQLVAL	KVISMNAEEG	VPFTAIREAS	LLKGLKHANI	300
301	VLLHDIHTK	ETLTFVFEYM	HTDLAQYMSQ	HPGGLHPHNV	RLFMFQLLRG	LAYIHHQHVL	360
361	HRDLKPQNL	ISHLGELKLA	DFGLARAKSI	PSQYTSSEVV	TLWYRPPDAL	LGATEYSSEL	420
421	DIWGAGCIFI	EMFQGQPLFP	GVSNIHQLE	KIWEVLGVP	EDTWPVSKL	PNYNPEWFPL	480
481	PTPRSLHVW	NRLGRVPEAE	DLASQMLKGF	PRDRVSAQEA	LVHDYF		540

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: CDK15 fragment

CDK15 wt ¹ Amino Acid Sequence							
1	MGQELCAKTV	QPGCSCYHCS	EGGEAHSCRR	SQPETTEAAF	KLTDLKEASC	SMTSFHPRGL	60
61	QAARAQKFKS	KRPRSNSDCF	QEEDLRQGFQ	WRKSLPFGAA	SSYLNLEKLG	EGSYATVYKG	120
121	ISRINGQLVA	LKVISMNAEE	GVPFTAIREA	SLLKGLKHAN	IVLLHDIHT	KETLTFVFEY	180
181	MHTDLAQYMS	QHPGGLHPHN	VRLFMFQLLR	GLAYIHHQHV	LHRDLKPQNL	LISHLGELKL	240
241	ADFGLARAKS	IPSQYTSSEV	VTLWYRPPDA	LLGATEYSSE	LDIWGAGCIF	IEMFQGQPLF	300
301	PGVSNILEQL	EKIWEVLGVP	TEDTWPVSK	LPNYPWFPP	LPTPRSLHV	WNRLGRVPEA	360
361	EDLASQMLKG	FPRDRVSAQE	ALVHDYFSAL	PSQLYQLPDV	SGVRLKPEMC	DLLASYQKGH	420
421	HPAQFSKCW						480

blue: CDK15 sequence expressed in recombinant protein

¹NCBI/Protein accession number NP_001248364.1

GST-CycB1 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEM	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHGG	RRRASVAAGI	240
241	LVPRGSPGLD	GICSIEEFRP	PWELAMALRV	TRNSKINAEN	KAKINMAGAK	RVPTAPAATS	300
301	KPGLRPRTAL	GDIGNKVSEQ	LQAKMPMKKE	AKPSATGKVI	DKKLPKPLEK	VPMLVPVPS	360
361	EPVPEPEPEP	EPEPVKEEKL	SPEPILVDTA	SPSPMETSGC	APAEEDLCOA	FSDVILAVND	420
421	VDAEDGADPN	LCSEYVKDIY	AYLRQLEEEQ	AVRPKYLLGR	EVTGNMRAIL	IDWLQVQMK	480
481	FRLLOETMYM	TVSIIIDRFMQ	NNCVPKMLQ	LVGVTAMFIA	SKYEEMYPPE	IGDFAFVTDN	540
541	TYTKHQIRQM	EMKILRALNF	GLGRPLPLHF	LRRASKIGEV	DVEQHTLAKY	LMELTMSDYD	600
601	MVHFPPSQIA	AGAFCLALKI	LDNGEWTPTL	QHLYSYTEES	LLPVMQHLAK	NVVMVNOGLT	660
661	KHMTVKNKYA	TSKHAKISTL	PQLNSALVQD	LAKAVAKV			720

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue: CycB1 boxed: variation from RefSeq

GST-CycB1 wt ² Amino Acid Sequence							
1	MALRVTRNSK	INAENKAKIN	MAGAKRVPTA	PAATSKPGLR	PRTALGDIGN	KVSEQLQAKM	60
61	PMKKEAKPSA	TGKVIDKKLP	KPLEKVPMLV	PVPVSEPVPE	PEPEPEPEPV	KEEKLSPEPI	120
121	LVDTASPSPM	ETSGCAPAEE	DLCQAFSDVI	LAVNDVDAED	GADPNLCSEY	VKDIYAYLRQ	180
181	LEEEQAVRPK	YLLGREVTGN	MRAILIDLWV	QVQMKFRLLQ	ETMYMTVSI	DRFMQNNCVP	240
241	KKMLQLVGVT	AMFIASKYEE	MYPPEIGDFA	FVTDNTYTKH	QIRQMEMKIL	RALNFGLRP	300
301	LPLHFLRRAS	KIGEVDVEQH	TLAKYLMELT	MLDYDMVHFP	PSQIAAG AFC	LALKILDNGE	360
361	WTPTLQHYLS	YTESLLPVM	QHLAKNVVMV	NOGLTKHMTV	KNKYATSKHA	KISTLPQLNS	420
421	ALVQDLAKAV	AKV					480

blue: CycB1 sequence expressed in recombinant protein Red: variant in recombinant protein

²NCBI/Protein accession number NP_114172.1
 HGNC identifier CycB1: CCNB1