

Certificate of Analysis



CDC42BPB

CDC42-binding protein kinase beta

Recombinant Human Active Protein Kinase

HGNC Symbol: CDC42BPB

Synonyms: MRCKB, DMPK-like

Product No.: 0382-0000-1

Lot: 001

Description: Human CDC42BPB, N-terminal fragment, amino acids M₁-H₄₇₂ (as in NCBI/Protein entry NP_006026.2), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: CDC42BPB Lot 001, was confirmed as CDC42BPB by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 84,444 Da

Expression: Baculovirus infected Sf9 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: 0.205 µg/µl

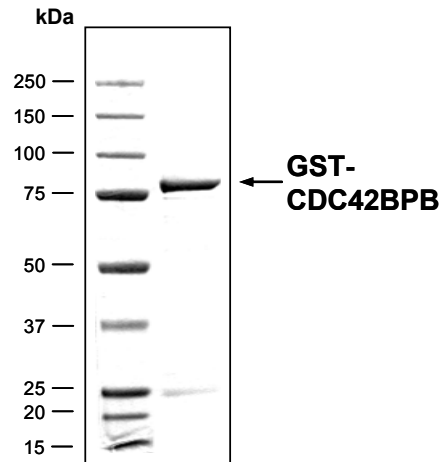
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

Specific kinase activity (P_i transfer): 14 pmol/µg×min

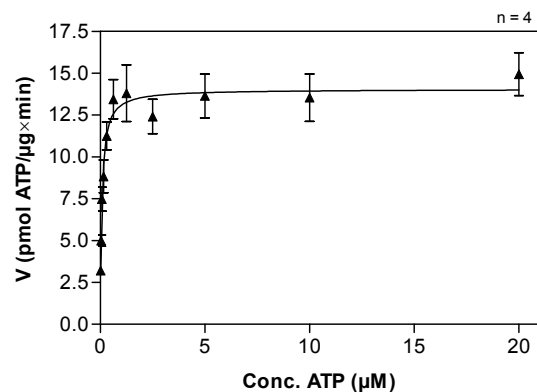
ATP-K_M: 0.1 µM

CDC42BPB Lot 001: Coomassie stain



2.0 µg GST-CDC42BPB

CDC42BPB 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: INSRR-derived peptide 20 µg/ml
 - CDC42BPB: 1.0 µg/ml
- Filter binding assay
 - MSFC membrane (Millipore)

CDC42BPB

Product No.: 0382-0000-1

CDC42BPB Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRLL	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 HHHHHH HG	RRRASVAAGI	240
241	LVPRGS PGLD	GIYARGIQAS	MSAKVRLKKL	EQLLLDGPWR	NESALSVETL	LDVLVCLYTE	300
301	CSHSALRRDK	YVAEFLEWAK	PFTQLVKEMQ	LHREDFEIIK	VIGRGAFGEV	AVVKMKNTER	360
361	IYAMKILNKW	EMLKRAETAC	FREERDVLVN	GDCQWITALH	YAFQDENHLY	LVMDYYVGGD	420
421	LLTLLSKFED	KLPEDMARFY	IGEMVLAIDS	IHQHLYVHRD	IKPDNVLLDV	NGHIRLADFG	480
481	SCLKMNDGT	VQSSVAVGTP	DYISPEILQA	MEDGMGKYGP	ECDWWSLQVC	MYEMLYGETP	540
541	FYAESLVETY	GKIMNHEERF	QFSPSHVTDVS	EEAKDLIQLR	ICSRERRLGQ	NGIEDFKKHA	600
601	FFEGLNWENI	RNLEAPYIPD	VSSPDSNFS	DVDDDLRNT	EILPPGSHTG	FSGLHLPFIG	660
661	FTFTTESCFS	DRGSLKSIMQ	SNTLTKDEDV	QRDLEHSLQM	EAYERRIRRL	EQEKLELSRK	720
721	LQESTQTVQS	LH					780

1-218: GST **Red:** HIS6-tag **Pink:** Thrombin cleavage site **blue:** CDC42BPB fragment

CDC42BPB wt ¹ Amino Acid Sequence							
1	MSAKVRLKKL	EQLLLDGPWR	NESALSVETL	LDVLVCLYTE	CSHSALRRDK	YVAEFLEWAK	60
61	PFTQLVKEMQ	LHREDFEIIK	VIGRGAFGEV	AVVKMKNTER	IYAMKILNKW	EMLKRAETAC	120
121	FREERDVLVN	GDCQWITALH	YAFQDENHLY	LVMDYYVGGD	LLTLLSKFED	KLPEDMARFY	180
181	IGEMVLAIDS	IHQHLYVHRD	IKPDNVLLDV	NGHIRLADFG	SCLKMNDGT	VQSSVAVGTP	240
241	DYISPEILQA	MEDGMGKYGP	ECDWWSLQVC	MYEMLYGETP	FYAESLVETY	GKIMNHEERF	300
301	QFSPSHVTDVS	EEAKDLIQLR	ICSRERRLGQ	NGIEDFKKHA	FFEGLNWENI	RNLEAPYIPD	360
361	VSSPDSNFS	DVDDDLRNT	EILPPGSHTG	FSGLHLPFIG	FTFTTESCFS	DRGSLKSIMQ	420
421	SNTLTKDEDV	QRDLEHSLQM	EAYERRIRRL	EQEKLELSRK	LQESTQTVQS	LHGSSRALSN	480
481	SNRDKEIKKL	NEEIERLKNK	IADSNRLERQ	LEDTVALRQE	REDSTQRLRG	LEKQHRVVRQ	540
541	EKEELHKQLV	EASERLKSQA	KELKDAHQQR	KLALQEFSEL	NERMAELRAQ	KQKVSRLRD	600
601	KEEEME VATQ	KVDAMRQEMR	RAEKLRKELE	AQLDDAVAEA	SKERKLREHS	ENFCKQMESE	660
661	LEALKVKQGG	RGAGATLEHQ	QEISKIKSEL	EKKVLFYEEE	LVRREASHVL	EVKNVKEEVH	720
721	DSESHQLALQ	KEILMLKDKL	EKSKRERHNE	MEEAVGTIKD	KYERERAMLF	DENKKLTAEN	780
781	EKLCSFVDKL	TAQNRQLEDE	LQDLAAKKE	VAHWEAQIAE	IIQVWSDEKD	ARGYLQALAS	840
841	KMTEELEALR	SSSLGSRTLD	PLWKVRRSQK	LDMSARLELQ	SALEAEIRAK	QLVQEELRKV	900
901	KDANLTLESK	LKDSEAKNRE	LLEEMEILKK	KMEEKFRADT	GLKLPDFQDS	IFEYFNTAPL	960
961	AHDLTFRFSS	ASEQETQAPK	PEASPSMSVA	ASEQQEDMAR	PPQRPSAVPL	PTTQALALAG	1020
1021	PKPKAHQFSI	KSFSSPTQCS	HCTSLMVGLI	RQGYACEVCS	FACHVCKDG	APQVCPIPE	1080
1081	QSKRPLGVDV	QRGIGTAYKG	HVKVPKPTGV	KKGWQRAYAV	VCECKLFLYD	LPEGKSTQPG	1140
1141	VIASQVLDLR	DDEFSVSSVL	ASDVIHATRR	DIPCIFRVTA	SLLGAPSKTS	SLLILTENEN	1200
1201	EKRKSWGILE	GLQSILHKNR	LRNQVVHVPL	EAYDSSLPLI	KAILTAAIVD	ADRIAVGLEE	1260
1261	GLYVIEVTRD	VIVRAADCKK	VHQIELAPRE	KIVILLCGRN	HHVHLYPWSS	LDGAEGSFDI	1320
1321	KLPETKGCQL	MATATLKRNS	GTCLFVAVKR	LILCYEIQRT	KPFHRKFNEI	VAPGSVQCLA	1380
1381	VLRDRLCVGY	PSGFCLLSIQ	GDGQPLNLVN	PNDPSLAFLS	QQSFDALCAV	ELESEEYLLC	1440
1441	FSHMGLYVDP	QGRRARAQEL	MWPAAPVACS	CSPTHVTVYS	EYGVDFDVR	TMEWVQTIGL	1500
1501	RRIRPLNSEG	TLNLLNCEPP	RLIYFKSKFS	GAVLNPPTS	DNSKKQMLRT	RSKRRFVFKV	1560
1561	PEEERLQQR	EMLRDPELRS	KMISNPTNFN	HVAHMGPGDG	MQVLMDLPLS	AVPPSQEER	1620
1621	GPAPTNLARQ	PPSRNKPYIS	WPSSGGSEPS	VTVPLRMSD	PDQDFDKEPD	SDSTKHSTPS	1680
1681	NSSNPSGPPS	PNSPHRSQLP	LEGLEQPACD	T			1740

blue: CDC42BPB sequence expressed in fusionprotein

¹NCBI/Protein accession number NP_006026.2