

ProQinase™ PIK3CA/PIK3R1

phosphatidylinositol-4,5-bisphosphate 3-kinase catalytic subunit alpha

Recombinant Human Active Lipid Kinase

HGNC Symbol: PIK3CA

Synonyms PIK3CA: p110-alpha, PI3K, PI3K-alpha
Synonyms PIK3R1: GRB1, p85, p85-ALPHA

Lipid Kinase Family: PI3K Class I

(according to: Phylogenomics of phosphoinositide lipid kinases: perspectives on the evolution of second messenger signaling and drug discovery: James R Brown & Kurt R Auger; BMC Evolutionary Biology 11, 4-14 (2011))

Product No.: 1161-1165-1

Lot: 003

Description: Human PIK3CA, full length, amino acids M₁-N₁₀₆₈ (as in [NCBI/Protein](#) entry NP_006209.2), N-terminal HIS₆ fusion protein with a Thrombin/TEV cleavage site and PIK3R1 full length, amino acids M₁-R₇₂₄ (as in [NCBI/Protein](#) entry NP_852664.1), N-terminal fused to a MYC-tag, coexpressed in Sf9 insect cells

Product identity: PIK3CA/PIK3R1-SV1 Lot 003, was confirmed as PIK3CA/PIK3R1 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{PIK3CA} : 130,237 Da

Theoretical MW_{PIK3R1} : 85,371 Da

Expression host: Sf9 insect cells

Purification: Immobilized Metal 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: 0.380 µg/µl

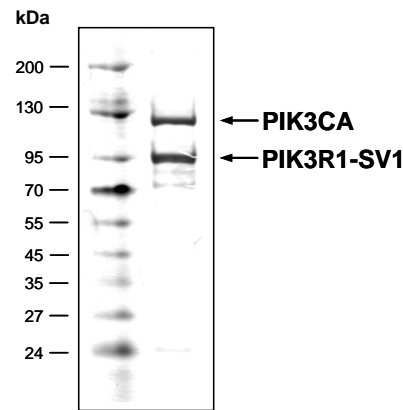
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

Specific kinase activity (P_i transfer): 15189 pmol/µg × min

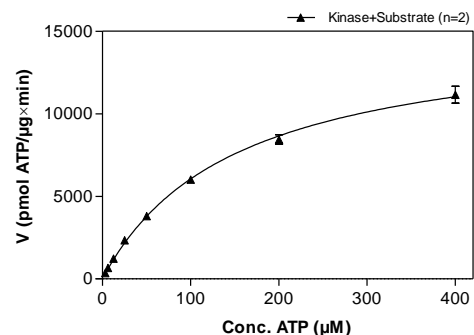
ATP-K_M: 151 µM

**PIK3CA/PIK3R1 Lot 003:
Coomassie stain**



4 µg PIK3CA/PIK3R1-SV1

**PIK3CA/PIK3R1 Lot 003:
Determination of V_{max} and K_M value for ATP**



Determination of K_M value & Specific activity:

• Assay conditions:

50 mM HEPES-NaOH, pH 7.5

3 mM MgCl₂

1 mM EGTA

100 mM NaCl

0,03% CHAPS

2 mM DTT

ATP (variable)

1 % (v/v) DMSO

Substrate: PIP₂: 50 µM / PS: 950 µM

PIP₂: 08:0 PI(4,5)P₂ (1,2-Dioctanoyl-sn-Glycero-3-(Phosphoinositol-4,5-Bisphosphate))

PS: 1-Palmitoyl-2-Oleoyl-sn-Glycero-3-(Phospho-L-Serine)

PIK3CA/PIK3R1: 1 µg/ml

• Assay technology:

ADP-Glo (Promega)

ProQinase™ PIK3CA/PIK3R1

Product No.: 1161-1165-1

HIS-PIK3CA Recombinant Fusion Protein Amino Acid Sequence												
1	MSPIDPMG	HH	HHH	GRRRAS	VAAGI	LVPRG	SPGLDGIYAR	TENLYFQ	GAM	GARGR	MPPRP	60
61	SSGELWGIHL	MPPRILVECL	LPNGMIVTLE	CLREATLITI	KHELFEARK	YPLHQLLQDE						120
121	SSYIFVSVTQ	EAEREFFDE	TRRLCDLRLF	QPFLKVIEPV	GNREEKILNR	EIGFAIGMPV						180
181	CEFDMVKDPE	VQDFRRNILN	VCKEAVDLRD	LNSPHSRAMY	VYPPNVESP	ELPKHIYNKL						240
241	DKGQIIVVIW	VIVSPNNDKQ	KYTLKINHDC	VPEQVIAEAI	RKKTRSMLLS	SEQLKLCVLE						300
301	YQGYILKVC	GCDEYFLEKY	PLSQYKIRS	CIMLGRMPNL	MLMAKESLYS	QLPMDCFTMP						360
361	SYSRRISTAT	PYMNGETSTK	SLWVINSALR	IKILCATYVN	VNIRDIDKIY	VRTGIYHGGE						420
421	PLCDNVNTQR	VPCSNPRWNE	WLNYDIYPD	LPRAARLCLS	ICSVKGRKGA	KEEHCPLAWG						480
481	NINLFDYTD	LVSQKMALNL	WPVPHGLEDL	LNPIGVTGSN	PNKETPCLEL	EFDWFSSVVK						540
541	FPDMSVIEEH	ANWSVSREAG	FSYSHAGLSN	RLARDNELRE	NDKEQLKATS	TRDPLSEITE						600
601	QEKDFLWSHR	HCVTPEIL	PKLLSVKWN	SRDEVAQMYC	LVKDWPPIKP	EQAMELLDCN						660
661	YPDPMVRGFA	VRCKEYLT	DKLSQYLQ	VQVLKYEYQL	DNLLVRFLLK	KALTNQRIGH						720
721	FFFWHLKSEM	HNKTYSQRF	LLLESYCRAC	GMYLKHLNRQ	VEAMEKLINL	TDILKQEKD						780
781	ETQKVQMKFL	VEQMRPDM	DALQGFSLP	NPAHQGLNLR	LEECRIMSSA	KRPLWLNWEN						840
841	PDIMSELLFQ	NNEIIFKNGD	DLRQDMLTLQ	IIRIMENIWQ	NOGLDLRMLP	YGCLSIGDCV						900
901	GLIEVVRNSH	TIMQIQCKGG	LKGALQFNH	TLHQWLKDKN	KGEIYDAAD	LFTRSCAGYC						960
961	VATFILGIGD	RHNSNIMVKD	DGQLFHIDFG	HFLDHKKKKF	GKREVRVPFV	LTQDFLIVIS						1020
1021	KGAQECTKTR	EFERFQEMCY	KAYLAIRQHA	NLFINLFSMM	LGSMPPELQS	FDDIAYIRKT						1080
1081	LALDKTEQEA	LEYFMQMND	AHGGWTTKM	DWIFHTIKQH	ALN							1140

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site Green: TEV cleavage site blue: PIK3CA

PIK3CA wt ¹ Amino Acid Sequence							
1	MPPRPSSGEL	WGIHLMPPRI	LVECLLPNGM	IVTLECLREA	TLITIKHELF	KEARKYPLHQ	60
61	LLQDESSYIF	VSVTQEAERE	EFFDETRRLC	DLRLFQPFK	VIEPVGNREE	KILNREIGFA	120
121	IGMPVCEFD	VKDPEVQDFR	RNILNVCKEA	VDLRDLNSPH	SRAMYVYPPN	VESSPELPKH	180
181	IYNKLDKQI	IVVIWVIVSP	NNDKQKYLK	INHDCVPEQV	IAEAIRKTR	SMLLSSEQLK	240
241	LCVLEYQGY	ILKVCGCDEY	FLEKYPLSQY	KYIRSCIMLG	RMPNMLMAK	ESLYSQLPMD	300
301	CFTMPSYSRR	ISTATPYMNG	ETSTKSLWVI	NSALRIKILC	ATYVNVNIRD	IDKIYVRTGI	360
361	YHGGEPLCDN	VNTQRVPCSN	PRWNEWLNVD	IYIPDLPRAA	RLCLSICSVK	GRKGAKKEHC	420
421	PLAWGNINLF	DYTDTLVSGK	MALNLWPVPH	GLEDLLNPIG	VTGSNPNET	PCLELEFDWF	480
481	SSVVKFPDMS	VIEEHANWSV	SREAGFSYSH	AGLSNRLARD	NELRENDKEQ	LKAISTRDPL	540
541	SEITEQEKDF	LWSHRHYCVT	IPEILPKLL	SVKWSRDEV	AQMYCLVKDW	PPIKPEQAME	600
601	LLDCNYPPDM	VRGFAVRCL	KYLTDKLSQ	YLIQLVQVLK	YEQYLDNLLV	RFLKALTN	660
661	QRIGHFFFWH	LKSEMHNKTV	SQRFGLLES	YCRACGMYLK	HLNRQVEAME	KLINLTDILK	720
721	QEKKDETQKV	QMKFLVEQMR	RPDFMDALQ	FLSPLNPAHQ	LGNLRLLECR	IMSSAKRPLW	780
781	LNWENPDIMS	ELLFQNEII	FKNGDDLQD	MLTLQIIRIM	ENIQNQGLD	LRMLPYGCLS	840
841	IGDCVGLIEV	VRNSHTIMQI	QCKGGLKCAL	QFNSHTLHQW	LKDKNKGIEY	DAAIDLFRS	900
901	CAGYCVATFI	LGIGDRHNSN	IMVKDDGQLF	HIDFGHFLDH	KKKFGYKRE	RVPFVLTQDF	960
961	LIVISKGAQE	CTKTREFERF	QEMCYKAYLA	IRQHANLFIN	LFSMMLGSGM	PELQSFDDIA	1020
1021	YIRKTLALDK	TEQEALEYFM	QOMNDAHGG	WTTKMDWIFH	TIQHALN		1080

blue: PIK3CA sequence expressed in recombinant protein

¹NCBI/Protein accession number NP_006209.2

MYC-PIK3R1 Recombinant Fusion Protein Amino Acid Sequence

1	MEEQKLISEE	DL	PMVMSAEG	YQYRALYDYK	KEREEDIDLH	LGDILTVMKG	SLVALGFSDG	60
61	QEARPEEIGW		LNGYNETTGE	RGDFPGTYVE	YIGRKKISPP	TPKPRPPRPL	PVAPGSSKTE	120
121	ADVEQQALTL		PDLAEQFAPP	DIAPPLLIK	VEAIEKKGLE	CSTLYRTQSS	SNLAELRQLL	180
181	DCDTPSVGLE		MIDVHVLADA	FKRYLLDLPN	PVIPAAYVSE	MISLAPEVQS	SEEIYQLLKK	240
241	LIRSPSIPHQ		YWLTLQYLLK	HFFKLSQTSS	KNLLNARVLS	EIFSPMLFRF	SAASSDNTEN	300
301	LIKVIEILIS		TEWNERQPAP	ALPPKPPKPT	TVANNGMNNN	MSLQDAEWYW	GDISREEVNE	360
361	KLRDTADGTF		LVRDASTKMH	GDYTLTLRKG	GNNKLIKIFH	RDGKYGFSDP	LTFSVVVELI	420
421	NHYRNESLAQ		YNPKLDVKLL	YPVSKYQQDQ	VVKEDNIEAV	GKKLHKYNTQ	FQEKSSREYDR	480
481	LYEEYTRTSQ		EIQMKRTAIE	AFNETIKIFE	EQCQTQERYS	KEYIEKFKRE	GNEKEIQRIM	540
541	HNYDKLKSRI		SEIIDSRRRL	EEDLKKQAAE	YREIDKRMNS	IKPDLIQLRK	TRDQYLMWLT	600
601	QKQVVRQKLN		EYLNQNTED	QYSLVEDDED	LPHHDEKTNW	VGSSNRNKAE	NLLRQKRDGT	660
661	FLVRESSKQG		CYACSVVDG	EVKHCVINKT	ATGYGFAEPE	NLYSSLKELV	LHYQHTSLVQ	720
721	HNDLNLVTLA		YPVYAQQRR					780

1-218: GST Red: MYC-tag blue: PIK3R1 boxed: variation from RefSeq

PIK3R1 wt² Amino Acid Sequence

1	MSAEGYQYRA	LYDYKKEREE	DIDLHLGDIL	TVNKGSLVAL	GFSQGQEARP	EEIGWLNQYN	60
61	ETTGERGDFP	GTIVEYIGRK	KISPPTPKPR	PPRPLPVAPG	SSKTEADVEQ	QALTLPLDLAE	120
121	QFAPPDIAPP	LLIKLVEAIE	KGLEKSTLY	RTQSSSNLAE	LRQLLDCDTP	SVDLEMIDVH	180
181	VLADAFKRYL	LDLNPVIPA	AVYSEMISLA	PEVQSSEEIY	QLLKKLIRSP	SIPHQYWLTL	240
241	QYLLKHFFKL	SQTSSKNLLN	ARVLSEIFSP	MLFRFSAASS	DNTENLIKVI	EILISTEWNE	300
301	RQPAPALPPK	PPKPTTVANN	GMNNMSLQD	AEWYWGDISR	EEVNEKLRDT	ADGTFVLRDA	360
361	STKMHGDYTL	TLRKGGNKLN	IKIFHRDGKY	GFSQDPLTFSS	VVELINHYRN	ESLAQYNPKL	420
421	DVKLLYPVSK	YQQDQVVKED	NIEAVGKKLH	EYNTQFQEK	REYDRLYEEY	TRTSQEIOMK	480
481	RTAIEAFNET	IKIFEEQCQT	QERYKEYIE	KFKREGNEKE	IQRIMHNYDK	LKSRISEIID	540
541	SRRLEEDLK	KQAAEYREID	KRMNSIKPDL	IQLRKTRDQY	LMWLTQKQV	QKKLNEWLGN	600
601	ENTEDQYSLV	EDDEDLPHHD	EKTWNVGSN	RNKAENLLRG	KRDGTFVRE	SSKQGCYACS	660
661	VVDGGEVVKH	VINKTATGYG	FAEPYNYLSS	LKELVLHYQH	TSLVQHNDL	NVTLAYPVYA	720
721	QQRR						780

blue: PIK3R1 sequence expressed in recombinant protein

²NCBI/Protein accession number NP_852664.1
E451K: SNP variation see NCBI/dbSNP ID: rs17852841