

ProQinase™ PIK3CG

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

Recombinant Human Active Lipid Kinase

HGNC Symbol: PIK3CG

Synonyms: p120-PI3K, PI3CG, PI3K, PI3K-gamma, PIK3

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.: 1163-0000-1

Lot: 002

Description: Human PIK3CG, full length, amino acids M₁-A₁₁₀₂ (as in [NCBI/Protein](#) entry NP_002640.2), N-terminal HIS₆ fusion protein with a Thrombin and TEV cleavage site, expressed in Sf9 insect cells

Product identity: PIK3CG Lot 002, was confirmed as PIK3CG by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 132,406 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

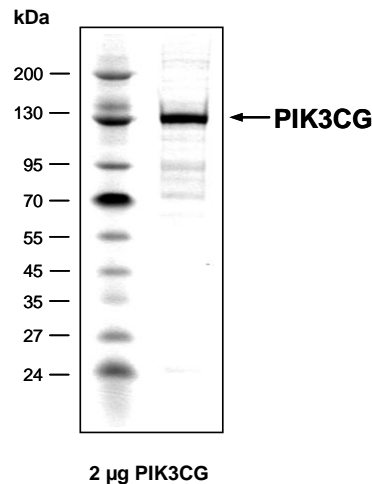
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.577 µg/µl (Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

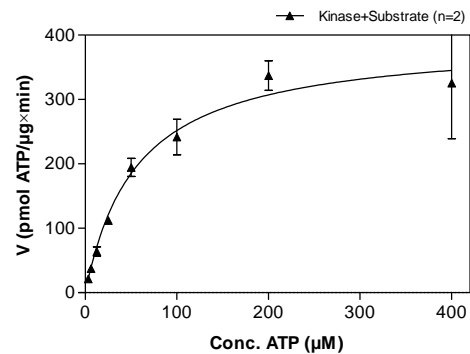
Biochemical Parameters:

Specific kinase activity (P_i transfer): 394 pmol/µg x min
ATP-K_M: 56 µM

**PIK3CG Lot 002:
Coomassie stain**



**PIK3CG Lot 002:
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 concentration
 - 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]
- PIK3CG: 1 µg/ml
- Assay technology: ADP-Glo (Promega)

ProQinase™ PIK3CG

Product No.: 1163-0000-1

HIS-PIK3CG Recombinant Fusion Protein Amino Acid Sequence												
1	MSPIDPMG	HH	HHHH	GRRRAS	VAAGI	LVPRG	SPGLDGIYAR	TENLYFQ	GAM	GARGR	MELEN	60
61	YKQPVV	LRED	NCRRRR	RMKP	RSAAAS	LSM	ELIPIEFVLP	TSQRKCK	SPE	TALLH	VAGHG	120
121	NVEQMK	QVW	LRALET	SVAA	DFYHRL	GPHH	FLLLYQKK	GQ	WYEIYDK	YQV	VQTLDC	180
181	KATHRSP	QI	HLVQR	HPPSE	ESQAFQ	RQLT	ALIGYD	VTDV	SNVHDE	LEF	TRRGLV	240
241	AEVASR	DPKL	YAMHP	PWVTSK	PLPEYL	WKKI	ANNCIF	FIVIH	RSTTSQ	TIKV	SPDDTP	300
301	QSFFT	KMAK	KSLMD	IPE	SEQDF	VRVC	GRDEYL	VGET	PIKNFQ	VRH	CLKNGE	360
361	VLDTPP	DPAL	DEV	RKEEWPL	VDDCTG	VGTGY	HEQLT	IHGKD	HESVFT	TVSLW	DCDRK	420
421	RGIDIP	VLP	NTDLT	VFVEA	NIQHGO	QVLC	QRRTSP	KPFT	EEVLWN	VWLE	FSIKIK	480
481	GALLNL	QIYC	GKAPAL	SSKA	SAESPS	SESK	GKVQLL	YYVN	LLLIDH	RFL	RRGEYV	540
541	QISGK	GEDQG	SFNAD	KLTS	TNPDK	ENSMS	ISILLD	NYCH	PIALPK	HQPT	PDPEGD	600
600	EMPNQ	LRKQL	EAI	IATDPLN	PLTAED	KELL	WHFRY	ESLKH	PKAYPK	LFS	VKWGQ	660
661	KTYQL	LARRE	VWDQ	SALDVG	LTMQLL	CNF	SDENV	RAIAV	QKLES	LEDD	VLHYLL	720
721	AVKFE	PYHDS	ALARF	LLKRG	LRNKR	IGHFL	FWFLR	SEIAQ	SRHYQ	RFV	ILEAYL	780
781	TAMLH	DFTQ	VQVI	EMLQKV	TLDIK	SLSAE	KYDVSS	QVIS	QLKQ	KLENLQ	NSQLP	840
841	PYDP	PGLK	GAGA	LAIEK	CKVMA	SKKK	PWLWF	KCADPT	ALSN	ETIGI	IFKHG	900
901	QILR	IMESI	W	ETESL	DLCLL	PYGC	ISTGDK	IGMIE	IVKDA	TTIAKI	QOST	960
961	VLNH	WLKE	KS	PTEEK	FQAAV	ERFV	YSCAGY	CVATF	VLGIG	DRHND	NIMIT	1020
1021	GHIL	GFYS	M	LGINK	ERV	VPF	VLTPD	FLFVM	GTSGK	KTS	PH	1080
1081	LLIIL	FSM	ML	MTGMP	QLTSK	EDIEY	IRDAL	TVGK	NEDAK	KYFLD	QIEV	1140
1141	WFLH	LVLG	IK	QGEK	SA							1200

Red: HIS6-tag Pink: Thrombin cleavage site Green: TEV cleavage site blue: PIK3CG

PIK3CG wt ¹ Amino Acid Sequence												
1	MELENY	KQPV	VLRED	NCRRR	RRMKPR	SAAA	SLSSMELIPI	EFVLPTS	QRK	CKSPET	ALLH	60
61	VAGHGN	VEQM	KAQV	WLRALE	TSVAAD	FYHR	LGP	HHFL	LLLY	QKKGQ	WYEIY	120
121	CLRYWK	KATHR	SPGQ	IHLVQR	HPPSE	ESQAF	QRQLT	ALIGY	DVTDV	SNVH	DELEF	180
181	VTPRMA	EVAS	RDPK	LYAMHP	WVTSK	PLPEY	LWKKI	ANNCI	FIVIH	RSTTS	QTIK	240
241	PGAILQ	SFFT	KMAK	KSLMD	IPESQ	SEQDF	VLRVC	GRDEY	LVGET	PIKNF	QWVRH	300
301	EEIHV	VLDTP	PDPAL	DEV	REEW	PLVDDCT	GVTGY	HEQLT	IHGKD	HESV	TVSLW	360
361	FRVKIR	GIDI	PVLPR	NTDLT	VFVEA	NIQHGO	QVLC	QRRTS	PKPFT	EEVLW	NVWLE	420
421	KDLPK	GALLN	LQIY	CGKAPA	LSSK	ASAESP	SSESK	GKVQL	LYYVN	LLLID	HRFLL	480
481	VLHMV	QISGW	GEDQ	SFNAD	KLTS	SATN	PDK	ENSMS	ISILL	NYCH	PIALP	540
541	DRVRA	EMP	NQ	LRKQ	LEAIIA	TDPLN	PLTAE	DKELL	WHFRY	ESLKH	PKAY	600
600	QEI	VAKTY	QL	LARRE	VWDQ	SALDVG	LTMQ	LDCNF	SDENV	RAIAV	QKLES	660
661	LQLV	QAVK	FE	PYHDS	SALARF	LLKRG	LRNKR	IGHFL	FWFLR	SEIAQ	SRHYQ	720
721	LRGCG	TAM	LH	DFTQ	QVQVIE	MLQK	VTL	DIK	SLSAE	KYDVS	SQVIS	780
781	ESFRV	PYDP	G	LKAG	ALAIEK	CKVM	SKKP	LWLEF	KCADP	TALS	NETIGI	840
841	DMLIL	QILRI	MESI	WETESL	DLCLL	PYGC	I	STGDK	IGMIE	IVKDA	TTIAK	900
901	AFKDE	VLNH	W	LKEK	SPT	EEK	FQAAV	ERFV	YSCAGY	CVATF	VLGIG	960
961	FHID	FGHILG	NYKS	F	LGIN	K	ERV	VPF	VLTPD	FLFVM	GTSGK	1020
1021	RHHTN	LLIIL	FSM	ML	MTGMP	QLTSK	EDIEY	IRDAL	TVGK	NEDAK	KYFLD	1080
1081	TVQFN	WFLHL	L	VLG	IKQGEK	SA						1140

blue: PIK3CG sequence expressed in recombinant protein

¹NCBI/Protein accession number NP_002640.2