

PIP5K1C

phosphatidylinositol-4-phosphate 5-kinase type 1 gamma

Recombinant Human Active Lipid Kinase

HGNC Symbol: PIP5K1C

Synonyms: PIP5K1-gamma, PIP5Kgamma, PIP5K-GAMMA, PIPK1g_v4

Lipid Kinase Family: PIP5K

(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.: 1222-0000-1

Lot: 003

Description: Human PIP5K1C, full length, amino acids M₁-T₆₆₈ (as in [NCBI/Protein](#) entry NP_036530.1), untagged, expressed in Sf9 insect cells

Product identity: PIP5K1C Lot 003, was confirmed as PIP5K1C by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 73,260 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, 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: 0.159 µg/µl

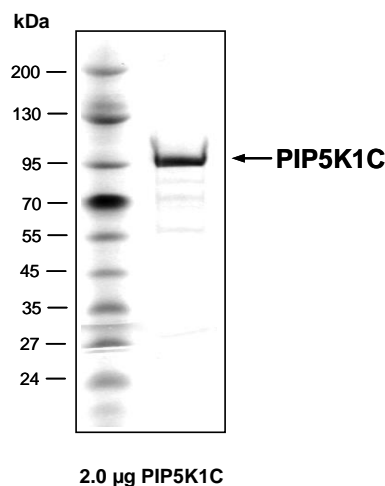
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

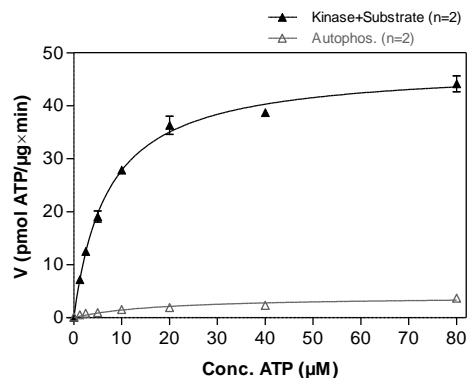
Specific kinase activity (P_i transfer): 47 pmol/µg × min

ATP-K_M: 7 µM

PIP5K1C Lot 003:
Coomassie stain



PIP5K1C Lot 003:
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 MnCl₂
- 3 µM Na-orthovanadate
- 1.2 mM DTT
- 50 µg/ml PEG_{20,000}
- ATP (variable)
- Substrate: PI: 25 µM / PS: 225 µM
- PI: L-alpha-phosphatidylinositol
- PS: 1-Palmitoyl-2-Oleoyl-sn-Glycero-3-[Phospho-L-Serine]
- PIP5K1C: 1 µg/ml

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.

PIP5K1C

Product No.: 1222-0000-1

PIP5K1C Recombinant Fusion Protein Amino Acid Sequence							
1	GPLAML	MELE	VPDEAESAEA	GAVPSEAAWA	AESGAAAGLA	QKKAAPTEVL SMTAQP	60
61	GKKLGHRGVD	ASGETTYKKT	TSSTLKGAIQ	LGIGYTVGHL	SSKPERDVL	MQDFYVVESIF	120
121	FPSEGSNLTP	AHHFQDFRFK	TYAPVAFRYF	RELFGIRPDD	YLYSLCNEPL	IELSNPGASG	180
181	SLFYVTSDD	FIIKTVMHKE	AEFLQKLLPG	YYMNLNQNPR	TLLPKFYGLY	CVQSGGKNIR	240
241	VVVMNINLPR	VVKMHLKFDL	KGSTYKRRAS	KKEKEKSFPT	YKDLDFMQDM	PEGLLLDADT	300
301	FSALVKTLQR	DCLVLESFKI	MDYSLLLGVH	NIDQHERERQ	AQGAQSTSDE	KRPVGQKALY	360
361	STAMESIQGG	AARGEAIESD	DTMGGIPAVN	GRGERLLLHI	GIIDILQSYR	FIKLEHTWK	420
421	ALVHDGDTVS	VHRPSFYAER	FFKFMSTVVF	RKNSSLKSSP	SKKGRGGALL	AVKPLGPTAA	480
481	FSASQIPSER	EEAQYDLRGA	RSYPTLEDEG	RPDLLPCTPP	SFEEATTASI	ATTLSSTLS	540
541	IPERSPSETS	EQPRYRRRTQ	SSGQDGRPQE	EPPAEEDLQQ	ITVQVEPACS	VEIVVPKEED	600
601	AGVEASPAGA	SAAVEVETAS	QASDEEGAPA	SQASDEEDAP	ATDIYFPTDE	RSWVYSPLHY	660
661	SAQAPPASDG	ESDT					720

1-6: legacy of 3C cleavage blue: PIP5K1C

PIP5K1C wt ¹ Amino Acid Sequence							
1	MELEVPDEAE	SAEAGAVPSE	AAWAAESGAA	AGLAQKKAAP	TEVLSMTAQP	GPGHGKKLGH	60
61	RGVDASGETT	YKKTTSSTLK	GAIQLGIGYT	VGHLSSKPER	DVLMQDFYVV	ESIFFPSEGS	120
121	NLTPAHHFQD	FRFKTYAPVA	FRYFRELFGI	RPDDYLYSLC	NEPLIELSNP	GASGSLFYVT	180
181	SDDEFIIKTV	MHKEAEFLQK	LLPGYYMNLN	QNPRTLLPKF	YGLYCVQSGG	KNIRVVMMN	240
241	ILPRVVKMHL	KFDLKGSTYK	RRASKKEKEK	SFPTYKDLDF	MQDMPEGLL	DADTFVSALVK	300
301	TLQRDCLVLE	SFKIMDYSL	LGVHNIDQHE	RERQAQGAQS	TSDEKRPVGG	KALYSTAMES	360
361	IQGAARGEAE	IESDDTMGGI	PAVNGRGERL	LLHIGIIDIL	QSYRFIKKLE	HTWKALVHDG	420
421	DTVSVHRPSF	YAERFFKFMS	NTVFRKNSSL	KSSPSKGRG	GALLAVKPLG	PTAAFSASQI	480
481	PSEREEAQYD	LRGARSYPTL	EDEGRPDLLP	CTPPSFEEAT	TASIATTLSS	TLSIPERSP	540
541	SETSEQPRYR	RRTQSSGQDG	RPQEEPPAEE	DLQITVQVE	PACSV EIVVP	KEEDAGVEAS	600
600	PAGASAAVEV	ETASQASDEE	GAPASQASDE	EDAPATDIYF	PTDERSWVYS	PLHYSAQAPP	660
661	ASDGESDT						720

blue: PIP5K1C sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_036530.1

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.