

ProQinase™ PDK1

3-phosphoinositide dependent protein kinase-1

Recombinant Human Active Protein Kinase

HGNC Symbol: PDPK1

Synonyms: PKB-like

Product No.: 0367-0000-1

Lot: 002

Description: Human PDK1, N-terminal fragment, amino acids M₁-M₄₆₀ (as in [NCBI/Protein](#) entry NP_002604.1), N-terminal HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: PDK1 Lot 002, was confirmed as PDK1 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 56,192 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 10 µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 1.223 µg/µl

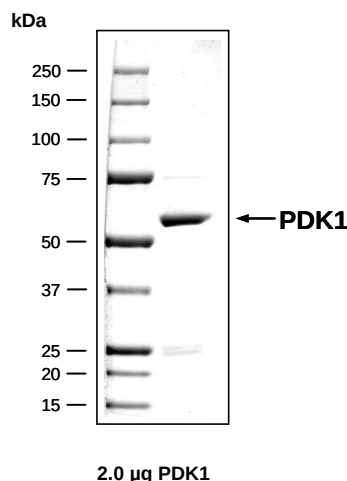
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

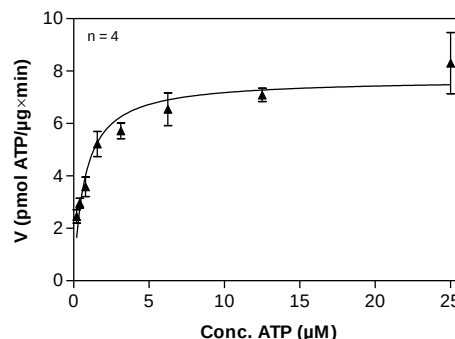
Specific kinase activity (P_i transfer): 8 pmol/µg × min

ATP-K_M: 0.7 µM

PDK1 Lot 002: Coomassie stain



PDK1 Lot 002: 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: tetra(LRRWSLG), 40 µg/ml
 - PDK1: 0.8 µg/ml
- Filter binding assay
 - MSFC membrane (Millipore)

Additional assay technology:

PDK1 Lot 002 was also successfully tested by Reaction Biology for the use with the ADP-Glo™ Kinase assay from Promega ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

ProQinase™ PDK1

Product No.: 0367-0000-1

HIS-PDK1 Recombinant Fusion Protein Amino Acid Sequence

1	MSPIDPMGHH HHHHGRRRAS	VAAGILVPRG SPGLDGICSR	NSMARTTSQL YDAVPIQSSV	60
61	VLCSPPSPSM VRTQTESSTP	PGIPGGSROG PAMDGTAAP	RPGAGSLQHA QPPPQPRKKR	120
121	PEDFKFGKIL GEGSFSTVVL	ARELATSREY AIKILEKRHI	IKENKVPYVT RERDVMSRLD	180
181	HPFFVKLYFT FQDDEKLYFG	LSYAKNGELL KYIRKIGSFD	ETCTRFYTAE IVSALEYLHG	240
241	KGIIHRDLKP ENILLNEDMH	IQITDFGTAK VLSPEKQAR	ANSFVGTAQY VSPELLTEKS	300
301	ACKSSDLWAL GCIIYQLVAG	LPPFRAGNEY LIFQKIIKLE	YDFPEKFFPK ARDLVEKLLV	360
361	LDATKRLGCE EMEGYGPLKA	HPFFESVTWE NLHQQTTPKL	TAYLPAMSED DEDCYGNYDN	420
421	LLSQFGCMQV SSSSSSHSLS	ASDTGLPQRS GSNIEQYIHD	LDSNSFELDL QFSEDEKRL	480
481	LEKQAGGNPW HQFVENNLIL	KM		540

Red: HIS6-tag Pink: Thrombin cleavage site blue: PDK1 fragment

PDK1 wt¹ Amino Acid Sequence

1	MARTTSQLYD AVPIQSSVVL	CSCPPSPSMVR TQTESSTPPG	IPGGSROGPA MDGTAAPRP	60
61	GAGSLQHAQP PPQPRKKRPE	DFKFGKILGE GSFSTVVLAR	ELATSREYAI KILEKRHIK	120
121	ENKVPYVTRE RDVMSRLDHP	FFVKLYFTFQ DDEKLYFGLS	YAKNGELLYK IRKIGSFDET	180
181	CTRFYTAEIV SALEYLHGKG	IIHRDLKPEN ILLNEDMHIQ	ITDFGTAKVL SPESQARAN	240
241	SFVGTAQYVS PELLTEKSAC	KSSDLWALGC ILYQLVAGLP	PFRAGNEYLI FQKIIKLEYD	300
301	FPEKFFPKAR DLVEKLLVLD	ATKRLGCEEM EGYGPLKAHP	FFESVTWENL HQQTTPKLTA	360
361	YLPAMSEDE DCYGNYNLL	SQFGCMQVSS SSSSHLSAS	DTGLPQRS GS NIEQYIHDLD	420
421	SNSFELDLOF SEDEKRLLE	KQAGGNPWHQ FVENNLILKM	GPVDRKGLF ARRRQLLLTE	480
481	GPHLYYVDPV NKVLKGEIPW	SQELRPEAKN FKTFVHTPN	RTYYLMDPSG NAHKWCRKIQ	540
541	EVWRQRYQSH PDAAVQ			600

blue: PDK1 sequence expressed in recombinant protein

¹NCBI/Protein accession number NP_002604.1