

## 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 M<sub>1</sub>-M<sub>460</sub> (as in [NCBI/Protein](#) entry NP\_002604.1), N-terminal HIS<sub>6</sub> 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**<sub>Fusion Protein</sub>: 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 10µl 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)

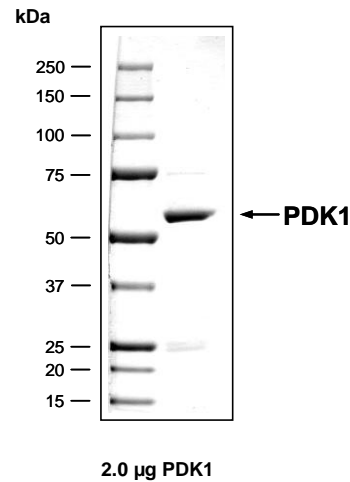
### Biochemical Parameters:

Specific kinase activity (P<sub>i</sub> transfer): 8 pmol/µg × min  
ATP-K<sub>M</sub>: 0.7 µM

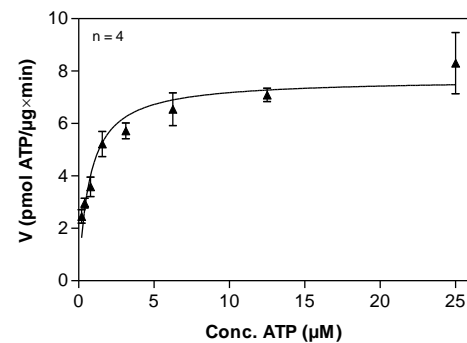
### Additional assay technology:

PDK1 Lot 002 was also successfully tested by ProQinase 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

### PDK1 Lot 002: Coomassie stain



### PDK1 Lot 002: Determination of V<sub>max</sub> and K<sub>M</sub> value for ATP



### Determination of K<sub>M</sub> value & Specific activity:

- Assay conditions:
  - 60 mM HEPES-NaOH, pH 7.5
  - 3 mM MgCl<sub>2</sub>
  - 3 mM MnCl<sub>2</sub>
  - 3 µM Na-orthovanadate
  - 1.2 mM DTT
  - 50 µg / ml PEG<sub>20,000</sub>
  - ATP (variable)
  - Substrate: tetra(LRRWSLG), 40 µg/ml
  - PDK1: 0.8 µg/ml
- Filter binding assay  
MSFC membrane (Millipore)

# PDK1

Product No.: 0367-0000-1

HIS-PDK1 Recombinant Fusion Protein Amino Acid Sequence											
1	MSPIDPMG	HH	HHH	GRRRAS	VAAGI	LVPRG	SPGLDGICSR	NSMARTTSQL	YDAVPIQSSV	60	
61	VLCS	CPSPSM	VRTQ	TESSTP	PGIP	GGSRQG	PAMDGTAAEP	RPGAGSLQHA	QPPPQPRKKR	120	
121	PEDFK	FGKIL	GEGS	FSTVVL	AREL	ATSREY	AIKILEKRHI	IKENKVPYVT	RERDVMSRLD	180	
181	HPFF	VKLYFT	FQD	DEKLYFG	LSYAK	NGELL	KYIRKIGSFD	ETCTRFYTAE	IVSALEYLHG	240	
241	KGI	IHRDLKP	ENILL	NEDMH	IQIT	DFGTAK	VLSPESKQAR	ANSFVGTQY	VSPELLTEKS	300	
301	ACK	SSLWAL	GCI	IYQLVAG	LPP	FRAGNEY	LIFQKI	IKLE	YDFPEKFFPK	ARDLVEKLLV	360
361	LDAT	KRLGCE	EMEG	YGPLKA	HPFF	FESVTWE	NLHQQT	PPKL	TAYLPAMSED	DEDCYGNYDN	420
421	LLS	QFGCMQV	SSSS	SSHSLS	ASDT	TGLPQRS	GSNIEQYIHD	LDSNSFELD	QFSEDEKRL	480	
481	LEK	QAGGNPW	HQF	VENNLIL	KM					540	

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

PDK1 wt <sup>1</sup> Amino Acid Sequence															
1	MARTT	SQLYD	AVPI	QSSVVL	CSC	SPSMVR	TQTES	STPPG	IPG	SRQGPA	MDGTAAEPRP	60			
61	GAGS	LQHAQP	PPQP	RKRPE	DFK	FGKILGE	GSF	STVVLAR	ELAT	SREYAI	KILEKRHIK	120			
121	ENK	VPYVTRE	RDV	MSRLDHP	FFV	KLYFTFQ	DDE	KLYFGLS	YAK	NGELLKY	IRKIGSFDET	180			
181	CTRF	YTAEIV	SALE	YLGK	IIHR	DLKPEN	ILL	NEDMHIQ	ITD	FGTAKVL	SPESKQARAN	240			
241	SFV	GTAQYVS	PELL	TEKSAC	KSS	DLWALGC	IIY	QLVAGLP	PFR	AGNEYLI	FQKI	IKLEYD	300		
301	FPE	KFFPKAR	DLVE	KLLVLD	ATK	RLGCEEM	EGY	GPLKAHP	FF	FVTWENL	HQOT	PPKLTA	360		
361	YLP	AMSEDE	DCY	GNYNLL	SQF	GCMQVSS	SSSS	SHLSAS	DTGL	PQRS	GS	NIEQYIHDLD	420		
421	SNS	FELD	LQF	SEDE	KRLLE	KQAG	GNPW	HQF	VENNLIL	KM	GPV	DKRGLF	ARRR	QLLLE	480
481	GPH	LYYVDPV	NKVL	KGEIPW	SQEL	RPEAKN	FKT	FFVHTPN	RTY	YLM	DPSG	NAHK	WCR	KIQ	540
541	EVWR	QRYQSH	PDA	AVQ											600

blue: PDK1 sequence expressed in recombinant protein

<sup>1</sup>NCBI/Protein accession number NP\_002604.1

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.