

PKC-eta

protein kinase C eta

Recombinant Human Active Protein Kinase

HGNC Symbol: PRKCH

Synonyms: nPKC-eta, PKC-L

Product No.: 0420-0000-1

Lot: 006

Description: Human PKC-eta, full length, amino acids M₁-P₆₈₃ (as in [NCBI/Protein](#) entry NP_006246.2), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: PKC-eta Lot 006, was confirmed as PKC-eta by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 107,725 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, 15 mM reduced glutathione, 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: 0.117 µg/µl

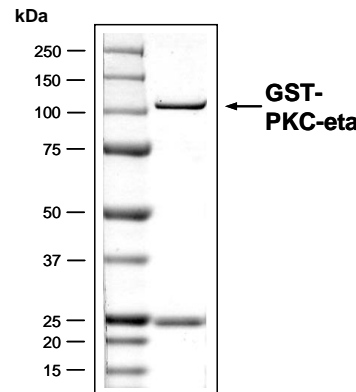
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

Specific kinase activity (P_i transfer): 71 pmol/µg × min

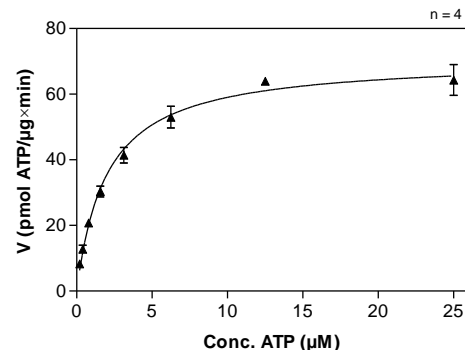
ATP-K_M: 2 µM

PKC-eta Lot 006: Coomassie stain



2.0 µg GST-PKC-eta

PKC-eta Lot 006: 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: Histone H2B 30 µg/ml
 - Kinase: 0.5 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

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.

PKC-eta

Product No.: 0420-0000-1

GST-PKC-eta Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RRRASVAAGI	240
241	LVPRGSPGLD	GICSRNSMSS	GTMKFNGLR	VRIGEAVGLQ	PTRWSLRHSL	FKKGHQLLDP	300
301	YLTVSV DQVR	VGQTSTKQKT	NKPTYNEEFC	ANVTDGGHLE	LAVFHETPLG	YDHFVANCTL	360
361	QFQELLRTTG	ASDTFEGWVD	LEPEGKVFV	ITLTGSFTEA	TLQRDRIFKH	FTRKRQRAMR	420
421	RRVHQINGHK	FMATYLRQPT	YCSHCREFIW	GVEGKQGYQC	QVCTCVVHKR	CHHLIVTACT	480
481	CQNNINKVDS	KIAEQRFGIN	IPHKFSIHNY	KVPTFCDHCG	SLLWGIMRQG	LQCKICKMNV	540
541	HIRCQANVAP	NCGVNAVELA	KTLAGMGLQP	GNISPTSCLV	SRSTLRRQ GK	ESSKEGNGIG	600
600	VNSSNRLGID	NFEFIRVLGK	GSFGKVM LAR	VKETGDLYAV	KVLK KDVILO	DDDVECTMTE	660
661	KRILSLARNH	PFLTQLFCCF	QTPDRLFFVM	EFVNGGDLMF	HIQSRRFDE	ARARFYAAEI	720
721	ISALMFLHDK	GIIYRDLKLD	NVLLDHEGHC	KLADFGMCKE	GICNGVTTAT	FCGTPDYIAP	780
781	EILQEMLYGP	AVDWWAMGVL	LYEMLCGHAP	FEAENEDDLF	EAILNDEVVY	PTWLHEDATG	840
841	ILKSFMTKNP	TMRLGSLTQG	GEHAILRHPF	FKEIDWAQLN	HRQIEPPFRP	RIKSREDVSN	900
901	FDPDFIKEEP	VLTPIDEGHL	PMINQDEFRN	FSYVSP ELQP			960

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue: PKC-eta

PKC-eta wt ¹ Amino Acid Sequence							
1	MSSGTMKFNG	YLRVRIGEAV	GLQPTRWSLR	HSLFKKGHQL	LDPYLTVSVD	QVRVGQTSTK	60
61	QKTNKPTYNE	EFCANVTDGG	HLELAVFHET	PLGYDHFVAN	CTLQFQELLR	TTGASDTFEG	120
121	WVDLEPEGKV	FVVITLTGSF	TEATLQRDRI	FKHFTRKRQR	AMRRRVHQIN	GHKFMATYLR	180
181	QPTYCSHCRE	FIWGVFGKQG	YQCQVCTCVV	HKRCHHLIVT	ACTCQNNINK	VDSKIAEQRF	240
241	GINIPHKFSI	HNYKVPTFCD	HCGSLLWGIM	RQGLQCKICK	MNVHIRCQAN	VAPNCGVNAV	300
301	ELAKTLAGMG	LQPGNISPTS	KLVSRSTLRR	QKESKEGN	GIGVNSSNRL	GIDNFEFIRV	360
361	LGKGSFGKVM	LARVKETGDL	YAVKVLKDV	ILQDDDV ECT	MTEKRILSLA	RNH PFLTQLF	420
421	CCFQTPDRLF	FVMEFVNGGD	LMFHIOKSRR	FDEARARFYA	AEIISALMFL	HDKGIIYRDL	480
481	KLDNVLLDHE	GHCKLADFGM	CKEGICNGVT	TATFCGTPDY	IAP EILQ EML	YGP AVDWWAM	540
541	GVLLYEMLCG	HAPFEAENED	DLFEAILNDE	VVYPTWLHED	ATGILKS FMT	KNPTMRLGSL	600
601	TQGG EHA ILR	HPFFKEIDWA	QLNHRQIEPP	FRPRIKSRED	VSNFDPDFIK	EEPVLTPIDE	660
661	GHLPMINQDE	FRNFSYVSPE	LQP				

blue: PKC-eta sequence expressed in recombinant protein

¹[NCBI/Protein](https://www.ncbi.nlm.nih.gov/Protein) accession number NP_006246.2

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.