

ProQinase™ PKC-zeta wt aa184-592 (PKM-zeta) protein kinase C zeta

Recombinant Human Active Protein Kinase

HGNC Symbol: PRKCZ

Synonyms: PKM-zeta, nPKC-zeta, PKC2

Product No.: 1240-0000-1

Lot: 001

Description: Human PKC-zeta, C-terminal fragment, amino acids M₁₈₄-V₅₉₂ (as in [NCBI/Protein](#) entry NP_002735.3), constitutive active fragment also denoted by PKM-zeta, N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: PKC-zeta wt aa184-592 Lot 001, was confirmed as PKC-zeta by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 74,896 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: 0.336 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

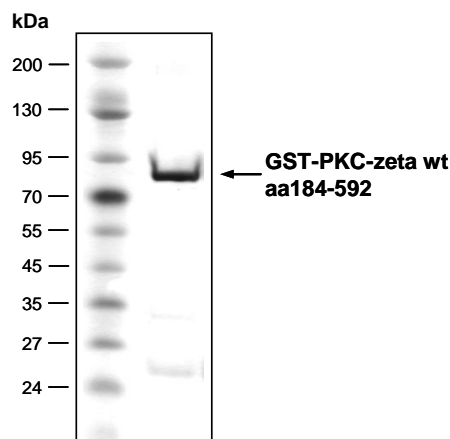
Biochemical Parameters:

Specific kinase activity (P_i transfer): 148 pmol/µg × min
ATP-K_M: 8.1 µM

Additional assay technology:

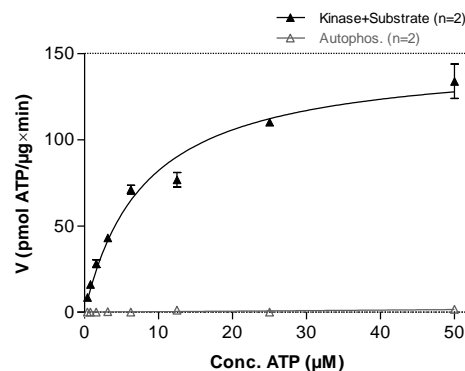
PKC-zeta wt aa184-592 Lot001 was also successfully tested by Reaction Biology for the use with the ADP-Glo™ Kinase assay from Promega ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

PKC-zeta wt aa184-592 Lot 001: Coomassie stain



2.0 µg GST-PKC-zeta wt aa184-592

PKC-zeta wt aa184-592 Lot 001: 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; 1 mM CaCl₂
 - 4 mM EDTA
 - 5 µg/ml Phosphatidylserine
 - 1 µg/ml 1,2 Dioleoyl-glycerol
 - 3 µM Na-orthovanadate
 - 1.2 mM DTT
 - 50 µg / ml PEG_{20,000}
 - ATP (variable)
 - Substrate: PKC-derived peptide 40 µg/ml
 - PKC-zeta wt aa184-592: 1 µg/ml
- Filter binding assay
MSPH membrane (Millipore)

ProQinase™ PKC-zeta wt aa184-592 (PKM-zeta)

Product No.: 1240-0000-1

GST-PKC-zeta wt aa184-592 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMG HHHHHG	RDS LEVLFQG	240
241	PLAMDSVMP	QEPPVDDKNE	DADLPSEETD	GIAYISSSRK	HDSIKDDSED	LKPVIDGMDG	300
301	IKISQGLGLQ	DFDLIRVIGR	GSYAKVLLVR	LKKNDQIYAM	KVVKKELVHD	DEDIDWVQTE	360
361	KHVFEQASSN	PFLVGLHSCF	QTTSRFLVI	EYVNGGDLMF	HMQRQRKLPE	EHARFYAAEI	420
421	CIALNFLHER	GIIYRDLKLD	NVLLDADGHI	KLTDYGMCKE	GLGPGD'TTST	FCGTPNYIAP	480
481	EILRGEYGF	SVDWWALGVL	MFEMMAGRSP	FDIITDNPDM	NTEDYLFQVI	LEKPIRIPRF	540
541	LSVKASHVLK	GFLNKDPKER	LGCRPQTGFS	DIKSHAFFRS	IDWDLLEKQ	ALPPFPQIT	600
601	DDYGLDNFDT	QFTSEPVQLT	PDEDAIKRI	DQSEFEGFEY	INPLLLSTEE	SV	660

1-218: GST **Red**: HIS6-tag **Green**: 3C cleavage site **blue**: PKC-zeta fragment

PKC-zeta wt ¹ Amino Acid Sequence							
1	MPSRTGPKME	GSGGRVRLKA	HYGGDIFITS	VDAATTFEEL	CEEVRDMCRL	HQQHPLTLKW	60
61	VDSEGDPCTV	SSQMELEEEAF	RLARQCRDEG	LIHVFPSTP	EQPGLPCPGE	DKSIYRRGAR	120
121	RWRKLYRANG	HLFQAKRFNR	RAYCGQCSE	IWGLARQGYR	CINCKLLVHK	RCHGLVPLTC	180
181	RKHMDSVMP	QEPPVDDKNE	DADLPSEETD	GIAYISSSRK	HDSIKDDSED	LKPVIDGMDG	240
241	IKISQGLGLQ	DFDLIRVIGR	GSYAKVLLVR	LKKNDQIYAM	KVVKKELVHD	DEDIDWVQTE	300
301	KHVFEQASSN	PFLVGLHSCF	QTTSRFLVI	EYVNGGDLMF	HMQRQRKLPE	EHARFYAAEI	360
361	CIALNFLHER	GIIYRDLKLD	NVLLDADGHI	KLTDYGMCKE	GLGPGD'TTST	FCGTPNYIAP	420
421	EILRGEYGF	SVDWWALGVL	MFEMMAGRSP	FDIITDNPDM	NTEDYLFQVI	LEKPIRIPRF	480
481	LSVKASHVLK	GFLNKDPKER	LGCRPQTGFS	DIKSHAFFRS	IDWDLLEKQ	ALPPFPQIT	540
541	DDYGLDNFDT	QFTSEPVQLT	PDEDAIKRI	DQSEFEGFEY	INPLLLSTEE	SV	600

blue: PKC-zeta sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_002735.3