

PKC-zeta aa184-592 K281M (PKM-zeta)

Protein kinase C, zeta

Recombinant Human inactivated Protein Kinase

HGNC Symbol: PRKCZ

Synonyms: PKM-zeta, nPKC-zeta, PKC2

Product No.: 1241-0000-1

Lot: 001

Description: Human PKC-zeta, C-terminal fragment, amino acids M₁₈₄-V₅₉₂ (as in NCBI/Protein entry NP_002735.3), fragment also termed PKM-zeta, mutationally inactivated by a K₂₈₁M mutation, N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Theoretical MW_{Protein}: 74,899 Da

Expression: Sf9 cells

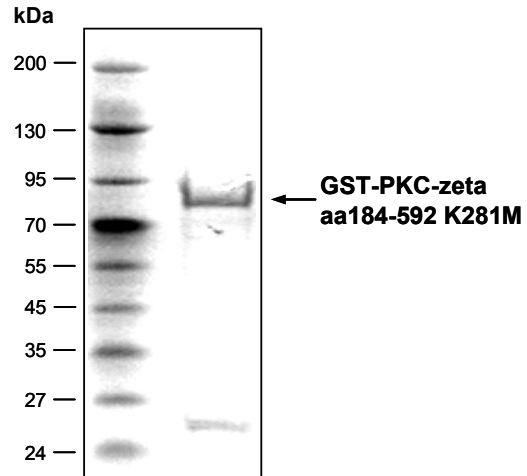
Purification: GST-Affinity Chromatography

Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 20% glycerol

Storage temperature: -80°C
Avoid repeated freeze-thaw cycles!

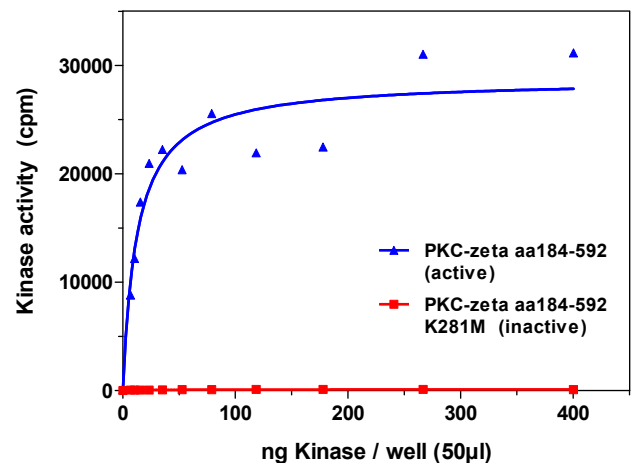
Protein concentration: 0.140 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Coomassie stain:



2.0 µg GST-PKC-zeta aa184-592 K281M

Kinase activity of PKC-zeta aa184-592 K281M (inactivated) vs PKC-zeta wt aa184-592:



Final assay concentrations:

- 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}
- 1 µM ATP
- Substrate: PKC-derived peptide (R₁₁-RFARKGSLRQKNV) 10 µg/ml
- Recombinant PKC-zeta aa184-592 K281M or PKC-zeta wt aa184-592 (variable concentrations)

Assay: ³³PanQinase® Assay

This product 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.

CoA V1.0 PKCzeta_aa184-592_K281M_inactivated_Lo001_V1.doc

PKC-zeta aa184-592 K281M (PKM-zeta)

Product No.: 1241-0000-1

PKC-zeta aa184-592 K281M Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQG WQATF	GGGDHPPKSD	PMGHHHHHGG	RDSLEVL FQG	240
241	PLAMDSVMPS	QEPPVDDKNE	DADLPSEETD	GIAYISSSRK	HDSIKD DSED	LKPVIDGMDG	300
301	IKISQGLGLQ	DFDLIRVIGR	GSYAKVLLVR	LKKNDQIYAM	MVVKKELVHD	DEDIDWVQTE	360
361	KHVFEQASSN	PFLVGLHSCF	QTTSRFLVI	EYVNGGDLMF	HMQRQRK LPE	EHARFYAAEI	420
421	CIALNFLHER	GIIYRDLKLD	NVLLDADGHI	KLTDYGMCKE	GLGPGDTTST	FCGTPNYIAP	480
481	EILRGEEYGF	SVDWWALGVL	MFEMMAGRSP	FDIITDNPDM	NTEDYLFQVI	LEKPIRIPRF	540
541	LSVKASHVLK	GFLNKDPKER	LGCRPQTGFS	DIKSHAFFRS	IDWDLLEKKQ	ALPPFPQIT	600
601	DDYGLDNFDT	QFTSEPVQLT	PDEDAIKRI	DQSEFEGFEY	INPLLLSTEE	SV	660

1-218: GST Red: HIS6-tag Green: 3C blue:PKC-zeta fragment M: K281M mutation

PKC-zeta wt ¹ amino acid sequence							
1	MPSRTGPKME	GSGGRVRLKA	HYGGDIFITS	VDAATTFEEL	CEEVRDMCRL	HQQHPLTLKW	60
61	VDSEGD PCTV	SSQMELEEEAF	RLARQCRDEG	LIIHVFPSTP	EQPGLPCPGE	DKSIYRRGAR	120
121	RWRKLYRANG	HLFQAKRFNR	RAYCGQC SER	IWGLARQGYR	CINCKLLVHK	RCHGLVPLTC	180
181	RKHMD SVMPS	QEPPVDDKNE	DADLPSEETD	GIAYISSSRK	HDSIKD DSED	LKPVIDGMDG	240
241	IKISQGLGLQ	DFDLIRVIGR	GSYAKVLLVR	LKKNDQIYAM	KVVKKELVHD	DEDIDWVQTE	300
301	KHVFEQASSN	PFLVGLHSCF	QTTSRFLVI	EYVNGGDLMF	HMQRQRK LPE	EHARFYAAEI	360
361	CIALNFLHER	GIIYRDLKLD	NVLLDADGHI	KLTDYGMCKE	GLGPGDTTST	FCGTPNYIAP	420
421	EILRGEEYGF	SVDWWALGVL	MFEMMAGRSP	FDIITDNPDM	NTEDYLFQVI	LEKPIRIPRF	480
481	LSVKASHVLK	GFLNKDPKER	LGCRPQTGFS	DIKSHAFFRS	IDWDLLEKKQ	ALPPFPQIT	540
541	DDYGLDNFDT	QFTSEPVQLT	PDEDAIKRI	DQSEFEGFEY	INPLLLSTEE	SV	600

blue: PKC-zeta sequence expressed in fusionprotein Red: K281M point mutation

¹NCBI/Protein accession number NP_002735.3