

Certificate of Analysis

PKC ϵ , active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-518, 14-518-K, 14-518M

Parent Lot # D8PN019U

The data presented in this document apply to the parent lot shown above and to all pack sizes derived from subsequent vialling runs of this parent lot. An alphabetical suffix after the parent lot number is used to denote each vialling run.

Product Description: N-terminal 6His-tagged, recombinant PKC epsilon, amino acids 2–end, expressed in Sf21 insect cells. Purified using Ni²⁺/NTA agarose. Purity 96.3% by SDS-PAGE and Coomassie blue staining. MW = 84.6kDa.

Specific Activity (Parent lot# D8PN019U): 2420U/mg, where one unit of PKC epsilon, active activity is defined as 1nmol phosphate incorporated into 50 μ M PKCtide (ERM^RPRKRQGSVRRRV) in the presence of lipid activator (cat# 20-133) per minute at 30°C with a final ATP concentration of 100 μ M.

Formulation: 0.126mg/ml of enzyme in 20mM Tris/HCl pH 7.5, 1mM EGTA, 1mM EDTA, 0.02% Triton X-100, 5% glycerol, 1mM PMSF, 10mM benzamidine, 1mM DTT. Frozen solution.

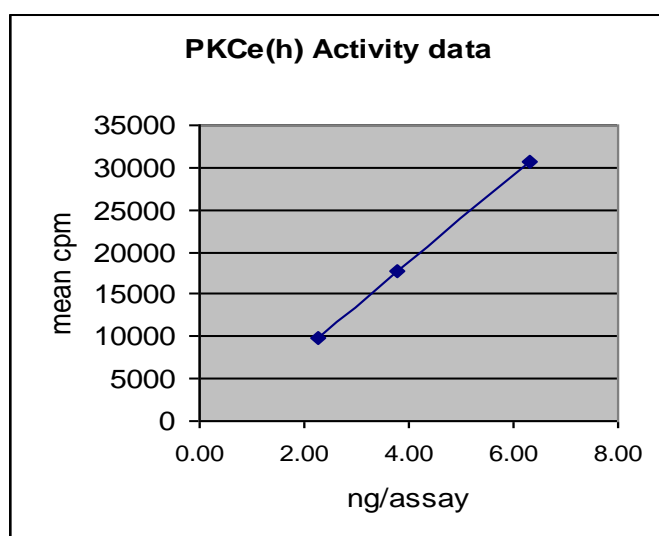
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

Handling Recommendations: Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled micro-centrifuge tubes and immediately snap-freeze the vials in liquid nitrogen prior to re-storage at -70°C.

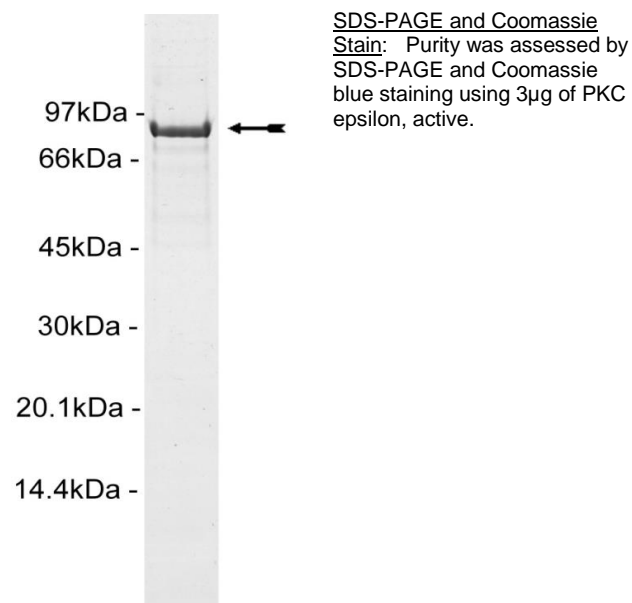
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 2.3–6.3ng of this lot of enzyme phosphorylated 50 μ M PKC tide in the assay described on page two. Assay background was subtracted from the actual counts to yield the results shown below.



MS Tryptic Fingerprint: Confirmed identity as human PKC epsilon with the translated sequence listed on page three.



Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

- 1. 10 x Reaction Buffer:** 200mM HEPES/NaOH pH7.4.
- 2. PKCtide(ERM₂PRKRQGSVRRRV):** Use at a final assay concentration of 50µM. Prepare a 500µM stock. Use 2.5µl of stock per assay point.
- 3. 10 x Lipid Activator:** 0.3% Triton X-100, 1mg/ml phosphatidylserine, 0.1mg/ml diacylglycerol). Use 2.5µl of stock per assay point.
- 4. PKC epsilon, active:** Dilute with 20mM HEPES/NaOH pH7.4, 0.03% Triton X-100. Use 2.3–6.3ng per assay point.
- 5. [γ -³³P]ATP:** 2.5 x magnesium acetate/[γ -³³P]ATP cocktail: 25mM MgAc and 0.25mM ATP to which is added [γ -³³P]ATP (specific activity approximately 500 - 800cpm/pmol as required.)

Assay Procedure (96 well plate format):

1. Add 2.5µl of 10 x reaction buffer per assay to wells.
2. Add 2.5µl of **PKCtide**.
3. Add 2.5µl of lipid activator.
4. Add **2.5µl (2.3–6.3ng) PKC epsilon, active**.
5. Make up to 15µl with dH₂O.
6. Add 10µl of diluted [γ -³³P]ATP mixture.
7. Incubate for 10 minutes at 30°C.
8. Transfer a 10µl aliquot onto the appropriate area of a **P30 Filtermat**.
9. Wash the filtermat three times for 5 minutes with 75mM phosphoric acid.
10. Wash the filtermat once for 2 minutes with methanol.
11. Transfer the filtermat to a sealable plastic bag and add 4ml of scintillation cocktail.
12. Read in a scintillation counter. Compare cpm of enzyme samples with cpm of control samples that contain all assay components plus 1µl of 30% phosphoric acid.

Certificate of Analysis

PKC epsilon Sequence Information

<u>Protein</u>	human PKC epsilon
<u>Tags</u>	N-terminal 6His
<u>Native sequence</u>	V8 of the recombinant protein is equivalent to V2 of human PKC epsilon
<u>Accession number</u>	GenBank X65293

Recombinant PKC epsilon amino acid sequence:

```

1  MHHHHHHVVF  NGLLKIKICE  AVSLKPTAWS  LRHAVGPRPQ  TFLLDPYIAL  NVDDSRIGQT
61  ATKQKTNSPA  WHDEFVTDVC  NGRKIELAVF  HDAPIGYDDF  VANCTIQFEE  LLQNGSRHFE
121 DWIDLEPEGR  VYVIDLSGS  SGEAPKDNEE  RVFRERMRPR  KRQGAVRRRV  HQVNGHKFMA
181 TYLRQPTYCS  HCRDFIWGVI  GKQGYQCQVC  TCVVHKRCHE  LIITKCAGLK  KQETPDQVGS
241 QRFVSNMPHK  FGIHNYKVPT  FCDHCGSLLW  GLLRQGLQCK  VCKMNVHRRC  ETNVAPNCGV
301 DARGIAKVLA  DLGVTDPKIT  NSGQRRKKLI  AGAESPQPAS  GSSPSEEDRS  KSAPTSPCDQ
361 EIKELENNIR  KALSFDNRGE  EHRAASSPDG  QLMSPGENGE  VRQGQAKRLG  LDEFNFIKVL
421 KGKSFQKVM  AELKKGDEVY  AVKVLKKDVI  LQDDVDCTM  TEKRLALAR  KHPYLTQLYC
481 CFQTKDRLFF  VMEYVNGGDL  MFQIQRSRKF  DEPRSRFYAA  EVTSALMFLH  QHGVIYRDLK
541 LDNILLDAEG  HCKLADFGMC  KEGILNGVTT  TFCGTPDYI  APEILQELEY  GPSVDWWALG
601 VLMYEMMAGQ  PPFEADNEDD  LFESILHDDV  LYPVWLSKEA  VSILKAFMTK  NPHKRLGCVA
661 SQNGEDAIAK  HPFFKEIDWV  LLEQKKIKPP  FKPRIKTKRD  VNNFDQDFTR  EEPVLTLDVE
721 AIVKQINQEE  FKGFSYFGED  LMP
    
```

Recombinant PKC epsilon nucleotide sequence:

```

1  atgcatcacc  atcatcacca  tgtagtgttc  aatggccttc  ttaagatcaa  aatctgcgag
61  gccgtgagct  tgaagccac  agcctggctg  ctgcccattg  cgggtgggacc  cgggccgcag
121 accttccttc  tcgacccta  cattgccttc  aatgtggacg  actcgcgcac  cggccaaacg
181 gccaccaagc  agaagaccaa  cagcccggcc  tggcacgacg  agttcgtcac  cgatgtgtgc
241 aacggacgca  agatcgagct  ggctgtcttt  cacgatgccc  ccataggcta  cgacgacttc
301 gtggccaact  gcaccatcca  gtttgaggag  ctgctgcaga  acgggagccg  ccacttcgag
361 gactggattg  atctggagcc  agaaggaaga  gtgtatgtga  tcatcgatct  ctcagggtcg
421 tcgggtgaag  ccctaaaga  caatgaagag  cgtgtgttca  gggaaacgat  cggcccgagg
481 aagcggcagg  gggcgtcag  gcgcagggtc  catcaggta  acggccaca  gttcatggcc
541 acatatcttc  ggcagccac  ctactgtcc  cattgcagag  acttcatctg  ggggtgcata
601 gaaagcagg  gataccagtg  tcaagtctgc  acctgcgtgg  tccacaagcg  gtggccagag
661 ctataatca  caaagtgtgc  tgggttaaag  aagcaggaga  cccccgacca  ggtgggctcc
721 cagcgtttca  gcgtcaacat  gccccacaag  ttcggtatcc  acaactaca  ggtccctacc
781 tctgcgatc  actgtgggtc  cctgctctgg  ggactcttgc  ggagggttt  gcagtgtaaa
841 gtctgcaaaa  tgaatgttca  ccgtcgatgt  gagaccaacg  tggctcccaa  ctgtggagtg
901 gatccagag  gaatcgcaa  agtactggcc  gacctgggcg  ttacccaga  caaatcacc
961 aacagcggcc  agagaaggaa  aaagctcatt  gctggtgccg  agtccccga  gcctgcttct
1021 ggaagctcac  catctgagga  agatcgatcc  aagtcagcac  ccacctccc  ttgtgaccag
1081 gaaataaaag  aacttgagaa  caacattcgg  aaagccttgt  catttgacaa  ccgaggagag
1141 gagcaccggg  cagcatcgtc  tcctgatggc  cagctgatga  gccccggtga  gaatggcgaa
1201 gtccggcaag  gccaggccaa  gcgcctgggc  ctggatgagt  tcaacttcat  caaggtgttg
1261 ggcaaaggca  gctttggcaa  ggtcatgttg  gcagaactca  agggcaaaga  tgaagtatat
1321 gctgtgaagg  tcttaaagaa  ggacgtcatc  cttcaggatg  atgacgtgga  ctgcacaatg
1381 acagagaaga  ggattttggc  tctggcacgg  aacaccctgt  accttacc  actctactgc
1441 tgcttcaga  ccaaggaccg  cctcttttc  gtcattggaat  atgtaaattg  tggagacctc
1501 atgtttcaga  ttcagcgtc  ccgaaaattc  gacgagcctc  gttcacggtt  ctatgctgca
1561 gaggtcacat  cggccctcat  gttcctccac  cagcatggag  tcatctacag  ggatttga
    
```

Certificate of Analysis

```
1621 ctggacaaca tccttctgga tgcagaaggt cactgcaagc tggctgactt cgggatgtgc
1681 aaggaaggga ttctgaatgg tgtgacgacc accacgttct gtgggactcc tgactacata
1741 gtccttgaga tcctgcagga gttggagtat ggcccctccg tggactgggtg ggccctgggg
1801 gtgctgatgt acgagatgat ggctggacag cctccccttg aggccgacaa tgaggacgac
1861 ctatctgagt ccatcctcca tgacgacgtg ctgtaccag tctggctcag caaggaggct
1921 gtcagcatct tgaagcttt catgacgaag aatccccaca agcgcctggg ctgtgtggca
1981 tgcagaatg gcgaggacgc catcaagcag cacccattct tcaaagagat tgactgggtg
2041 ctcttgagc agaagaagat caagccacc ttcaaaccac gcattaaaac caaaagagac
2101 gtcaataatt ttgaccaaga ctttaccgg gaagagccgg tactcaccct tgtggacgaa
2161 gcaattgtaa agcagatcaa ccaggaggaa ttcaaaggtt tctcctactt tggatgaagac
2221 ctgatgccct ga
```

Reviewed and approved by site quality representative.

Unless otherwise stated in our catalogue or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

© 2014 Eurofins Pharma Discovery Services UK Limited is an independent member of Eurofins Discovery Services.