

Certificate of Analysis

PKC ζ , active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-525, 14-525-K, 14-525M

Parent Lot # 1718700

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 human PKC ζ , amino acids 2–end. Expressed by baculovirus in Sf21 insect cells. Purified using Ni²⁺/NTA agarose. Purity 91.8% by SDS-PAGE and Coomassie blue staining. MW = 68.9kDa.

Specific Activity (Parent lot# 1718700): 1377U/mg, where one unit of PKC ζ activity is defined as 1nmol phosphate incorporated into 50 μ M PKCtide (ERM_RPRKRQGSVRRRV) per minute at 30°C with a final ATP concentration of 100 μ M.

Formulation: 0.538mg/ml of enzyme in 20mM Tris/HCl pH7.5, 0.02% Triton X-100, 5% glycerol, 1mM EDTA, 1mM EGTA, 0.1% 2-mercaptoethanol, 1mM PMSF, 10mM benzamidine. Frozen solution.

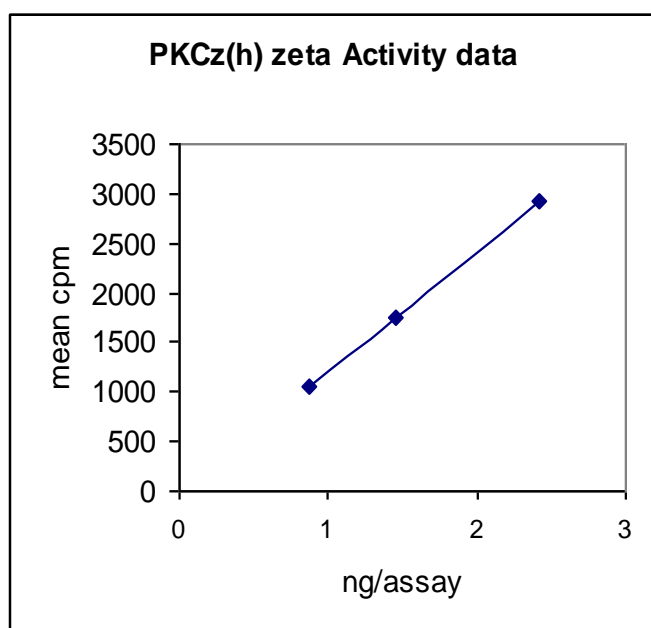
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 6 months from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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

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

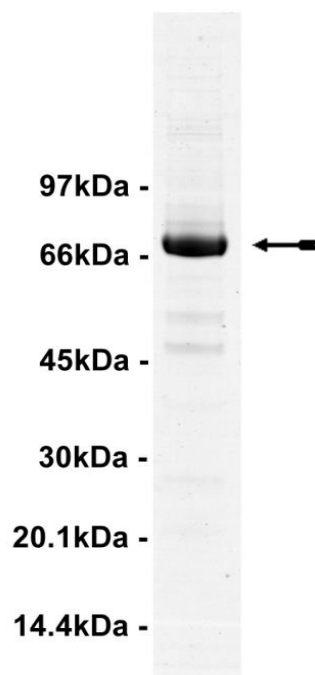
Quality Control Testing

Kinase Assay: 0.9–2.4ng of this lot of enzyme phosphorylated 50 μ M PKCtide 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 PKC ζ with the translated sequence listed on page three.

SDS-PAGE and Coomassie Stain: Purity was assessed by SDS-PAGE and Coomassie blue staining using 3 μ g of active PKC ζ .



Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

- 1. 5 x Reaction Buffer:** 40mM MOPS/NaOH pH7.0, 1mM EDTA.
- 2. PKCtide (ERM_RPRKRQGSVRRRV):** Use at a final assay concentration of 50µM. Prepare a 500µM stock and add 2.5µl of stock per assay point.
- 3. PKCζ, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 5% glycerol, 0.01% Brij-35, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 0.9–2.4ng per assay point.
- 4. [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 assay):

1. Add 5µl of 5 x reaction buffer per assay to wells.
2. Add 2.5µl of **PKCtide**.
3. Add **2.5µl (0.9–2.4ng) PKCζ, active**.
4. Add 5µl of dH₂O.
5. Add 10µl of diluted [γ-³³P]ATP mixture.
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5µl of 3% phosphoric acid.
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 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 ζ Sequence Information

| | |
|-------------------------|---|
| Protein | Human PKC ζ |
| Tags | N-terminal 6His |
| Native sequence | P10 of the recombinant protein is equivalent to P2 of human PKC ζ |
| Accession number | GenBank BC014270 |

Recombinant PKC ζ amino acid sequence:

```

1 MHHHHHHEFP SRTGPKMEGS GGRVRLKAHY GGDIFITSVD AATTFEELCE EVRDMCRLHQ
61 QHPLTLKWVD SEGDPCTVSS QMELEEAFLR ARQCRDEGLI IHVFPSTPEQ PGLPCPGEDK
121 SIYRRGARRW RKLRYRANGHL FQAKRFNRRA YCGQCSEIRW GLARQGYRCI NCKLLVHKRC
181 HGLVPLTCRK HMDVMPSQE PPVDDKNEDA DLPSEETDGI AYISSSRKHD SIKDDSEDLK
241 PVIDGMDGIK ISQGLGLQDF DLIRVIGRGS YAKVLLVRLK KNDQIYAMKV VKKELVHDDE
301 DIDWVQTEKH VFEQASSNPF LVGLHSCFQT TSRLFLVIEY VNGGDLMFHM QRQRKLPEEH
361 ARFYAAEICI ALNFLHERGI IYRDLKLDNV LLDADGHIKL TDYGMCKEGL GPGDTTSTFC
421 GTPNYIAPEI LRGEYGFVS DWWALGVLMF EMMAGRSPFD IITDNPDMNT EDYLFQVILE
481 KPIRIPRFLS VKASHVLKGF LNKDPKERLG CRPQTGFSDI KSHAFFRSID WDLLEKKQAL
541 PPFQPQITDD YGLDNFDTQF TSEPVLTPD DEDAIKRIDQ SEFEGFEYIN PLLLSTEEVS
  
```

Recombinant PKC ζ nucleotide sequence:

```

1 atgcatcatc accatcacca tgaattcccc agcaggaccg gccccaagat ggaagggagc
61 ggcggccgcg tccgcctcaa ggcgcattac ggggggggaca tcttcatcac cagcgtggac
121 gccgccacga ccttcgagga gctctgtgag gaagtgagag acatgtgtcg tctgcaccag
181 cagcaccgca tcaccctcaa gtgggtggac agcgaaggatg acccttgacg ggtgtcctcc
241 cagatggagc tggaaaggc tttccgctg gcccgtcagt gcagggatga aggcctcatc
301 attcatgttt tcccagcac cctgagcag cctggcctgc catgtccggg agaagacaaa
361 tctatctacc gccggggagc cagaagatgg aggaagctgt accgtgcca cggccacctc
421 ttccaagcca agcgccttaa caggagagcg tactgcggtc agtgcagcga gaggatatgg
481 ggcttcgca ggcaaggcta caggtgcatc aactgcaaac tgctggtcca taagcgtgctc
541 caggcctcg tcccgtgac ctgcaggaag catatggatt ctgtcatgcc ttccaagag
601 cctccagtag acgacaagaa cgaggacgcc gaccttcctt ccgaggagac agatggaatt
661 gcttacatct cctcatccc gaagcatgac agcattaaag acgactcggg ggaccttaag
721 ccagttatcg atgggatgga tggaaatcaa atctctcagg ggcttgggct gcaggacttt
781 gacctaatca gactcatcgg gcgaggagc tacgccaagg ttctcctggt gcggttgaag
841 aagaatgacc aaatttacgc catgaaagtg gtgaagaaag agctggtgca tgatgacgag
901 gatattgact ggttacagac agagaagcac gtgtttgagc aggcattccag caacccttc
961 ctggtcggat tacactcctg cttccagacg acaagtcggt tgttcctggt cattgagtac
1021 gtcaacggcg gggacctgat gttccacatg cagagcaga ggaagctccc tgaggagcac
1081 gccaggttct acgaggcga gatctgcatc gccctcaact tcctgcacga gagggggac
1141 atctacaggg acctgaagct ggacaacgct ctcctggatg cggacgggca catcaagctc
1201 acagactacg gcatgtgcaa ggaaggctg ggccctggtg acacaacgag cactttctgc
1261 ggaacccgga attacatcgc ccccgaatc ctgcggggag aggagtacgg gttcagcgtg
1321 gactggtggg cgctgggagt cctcatgttt gagatgatgg ccgggcgctc cccgttcgac
1381 atcatcaccg acaaccgga catgaacaca gaggactacc tttccaagt gatcctggag
1441 aagccatcc ggatccccg gttcctgtcc gtcaaagcct cccatgtttt aaaaggattt
1501 ttaaataagg accccaaaga gaggctcggc tgccggccac agactggatt ttctgacatc
1561 aagttccacg cgttcttcg cagcatagac tgggacttgc tggagaagaa gcaggcgtc
1621 cctccattcc agccacagat cacagacgac tacggtctgg acaactttga cacacagttc
  
```

Certificate of Analysis

```
1681 accagcgagc ccgtgcagct gaccccagac gatgaggatg ccataaagag gatcgaccag
1741 tcagagttcg aaggctttga gtatatcaac ccattattgc tgtccaccga ggagtcggtg
1801 tga
```

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.