

## Certificate of Analysis

### PKC mu, active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-508, 14-508-K, 14-508M

Parent Lot # 25876U

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, full length, human PKC mu expressed in Sf21 insect cells. Purified using Ni<sup>2+</sup>/NTA agarose. Purity 78.8% by SDS-PAGE and Coomassie blue staining. MW = 105kDa.

**Specific Activity (Parent lot# 25876U):** 414U/mg, where one unit of PKC mu, active activity is defined as 1nmol phosphate incorporated into 30µM (KKLNRTLVA) per minute at 30°C with a final ATP concentration of 100µM.

**Formulation:** 0.222mg/ml of enzyme in 20mM Tris/HCl pH7.5, 0.02% Triton X-100, 5% glycerol, 1mM EGTA, 1mM EDTA, 10mM benzamidine, 1mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution.

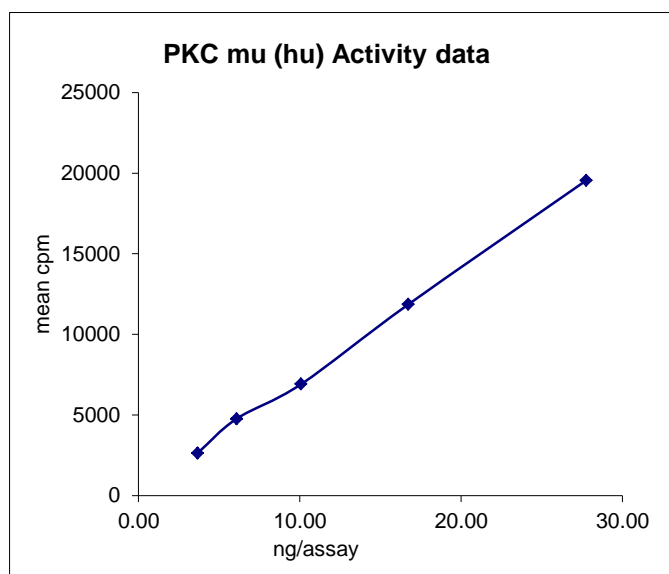
**Storage and Stability:** On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

**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.

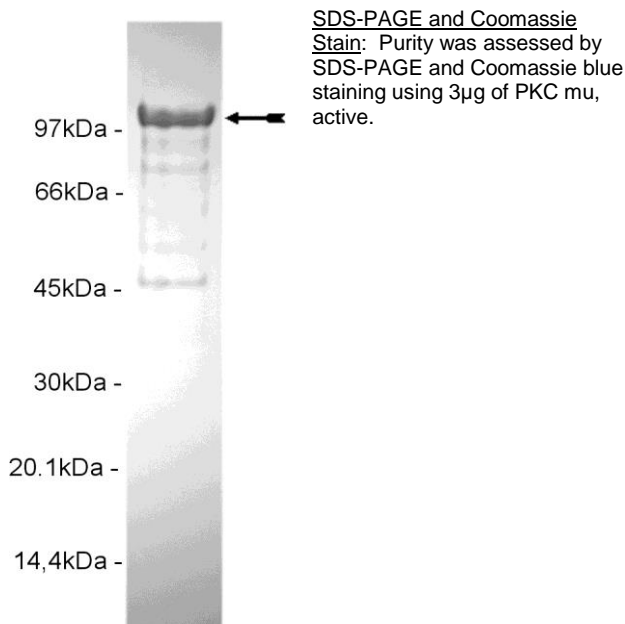
**FOR IN VITRO RESEARCH USE ONLY  
NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**Kinase Assay:** 3.7–27.8ng of this lot of enzyme phosphorylated 30µM (KKLNRTLVA) 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 mu 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, 0.3% Triton X-100.
- 2. (KKLNRTLVA):** Use at a final assay concentration of 30 $\mu$ M. Make up a 30 $\mu$ M stock. Add 2.5 $\mu$ l of stock per assay point.
- 3. PKC mu, active:** Dilute with 20mM HEPES/NaOH pH7.4. Use 3.7–27.8ng per assay point.
- 4. [ $\gamma$ -<sup>33</sup>P]ATP:** 2.5 x magnesium acetate/[ $\gamma$ -<sup>33</sup>P]ATP cocktail: 25mM MgAc and 0.25mM ATP to which is added [ $\gamma$ -<sup>33</sup>P]ATP (specific activity approximately 500 - 800cpm/pmol as required.)

#### Assay Procedure (96 well plate format):

1. Add 2.5 $\mu$ l of 10 x reaction buffer per assay to wells.
2. Add 2.5 $\mu$ l of **(KKLNRTLVA)**.
3. Add **2.5 $\mu$ l (3.7–27.8ng) PKC mu, active**.
4. Add 7.5 $\mu$ l of dH<sub>2</sub>O.
5. Add 10 $\mu$ l of diluted [ $\gamma$ -<sup>33</sup>P]ATP mixture.
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5 $\mu$ l of 3% phosphoric acid.
8. Transfer a 10 $\mu$ l aliquot onto the appropriate area of a **P30 Filtermat**.
9. Wash the filtermat three times for 5minutes 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 1 $\mu$ l of 30% phosphoric acid.

## Certificate of Analysis

### PKC mu Sequence Information

<b><u>Protein</u></b>	Human PKC mu
<b><u>Tags</u></b>	N-terminal 6His
<b><u>Native sequence</u></b>	M29 of the recombinant protein is equivalent to M1 of human PKC mu
<b><u>Accession number</u></b>	GenBank NM_002742

#### Recombinant PKC mu amino acid sequence:

```

1 MSYYHHHHHH DYDIPTTENL YFQGAMGSMS APPVLRPPSP LLPVAAAAAA AAAALVPGSG
61 PGPAPFLAPV AAPVGGISFH LQIGLSREPV LLLQDSSGDY SLAHVREMAC SIVDQKFPEC
121 GFYGMYSKIL LFRHDPTSEN ILQLVKAASD IQEGDLIEVV LSASATFEDF QIRPHALFVH
181 SYRAPAFCDH CGEMLWGLVR QGLKCEGCGL NYHKRCAFKI PNNCSGVRRR RLSNVSLTGV
241 STIRTSSAEL STSAPDEPLL QKSPSEFIG REKRSNSQSY IGRPIHLDKI LMSKVVPHT
301 FVIHSYTRPT VCQYCKKLLK GLFRQGLQCK DCRFNCHKRC APKVPNNCLG EVTINGDLLS
361 PGAESDVVME EGSDDNDSER NSGLMDDMEE AMVQDAEMAM AECQNDSGEM QDPDPDHEDA
421 NRTISPSTSN NIPLMRVVQS VKHTKRKSSST VMKEGWMVHY TSKDTLRKRH YWRLDSKCIT
481 LFQNDTGSRY YKEIPLSEIL SLEPVKTSAL IPNGANPHCF EITTANVVYY VGENVVNPSS
541 PSPNNSVLTS GVGADVARMW EIAIQHALMP VIPKGSSVGT GTNLHRDISV SISVSNQCIO
601 ENVDISTVYQ IFPDEVLGSG QFGIVYGGKH RKTGRDVAIK IIDKLRFPFK QESQLRNEVA
661 ILQNLHHPGV VNLECMFETP ERVFFVMEKL HGDMLEMILS SEKGRLEPHI TKFLITQILV
721 ALRHLHFKNL VHCDLKPENV LLASADFPFQ VKLCDFGFAR IIGEKSFRRS VVGTPAYLAP
781 EVLRNKGYNR SLDMWSVGVI IYVSLSGTFP FNEDEDIHDQ IQNAAFMYPP NPWKEISHEA
841 IDLINLLQV KMRKRYSDVK TLSHPWLQDY QTWLDLRELE CKIGERYITH ESDDLREWEKY
901 AGEQRLQYPT HLINPSASHS DTPETEETEM KALGERVSIL

```

#### Recombinant PKC mu nucleotide sequence:

```

1 atgtcgtact accatcacca tcaccatcac gattacgata tccaacgac cgaaaacctg
61 tattttcagg gcgccatggg atccatgagc gcccctccgg tcttgcggcc gccagtcctg
121 ctgctgcccg tggcggcggc agctgccgca gcgcccgccg cactgggtccc aggggtccggg
181 cccgggcccg cgccgttctt ggctcctgtc gcgcccccgg tcgggggcat ctcgttccat
241 ctgcagatcg gctgagccg tgagccggtg ctgctgctgc aggactcgtc cggggactac
301 agcctggcgc acgtccgcga gatggcttgc tccattgtcg accagaagt cctgaatgt
361 ggtttctacg gaatgtatga taagatcctg ctttttcgcc atgaccctac ctctgaaaac
421 atccttcagc tggtgaaagc ggccagtgat atccaggaag gcatcttat tgaagtggtc
481 ttgtcagctt ccgccacctt tgaagacttt cagattcgtc cccacgctct ctttgttcat
541 tcatacagag ctccagcttt ctgtgatcac tgtggagaaa tgctgtgggg gctggtactg
601 caaggtctta aatgtgaagg gtgtggtctg aattaccata agagatgtct atttaaaata
661 cccaacaatt gcagcgggtg gaggcggaga aggtctcaa acgtttccct cactggggtc
721 agcaccatcc gcacatcatc tgctgaactc tctacaagtg cccctgatga gcccttctg
781 caaaaatcac catcagagtc gtttattggt cgagagaaga ggtcaaattc tcaatcatac
841 attggacgac caattcacct tgacaagatt ttgatgtcta aagttaaagt gccgcacaca
901 tttgtcatcc actcctacac ccggcccaca gtgtgccagt actgcaagaa gcttctgaag
961 gggcttttca ggcagggctt gcagtgcaa gattgcagat tcaactgcca taaacgttgt
1021 gcaccgaaag taccaaaca ctgccttggc gaagtgacca ttaatggaga tttgcttagc
1081 cctggggcag agtctgatgt ggtcatggaa gaaggagtg atgacaatga tagtgaaagg
1141 aacagtgggc tcatggatga tatggaagaa gcaatggtcc aagatgcaga gatggcaatg
1201 gcagagtgcc agaacgacag tggcgagatg caagatccag acccagacca cgaggacgcc
1261 aacagaacca tcagtccatc aacaagcaac aatatcccac tcatgagggt agtgcagtct
1321 gtcaaacaca cgaagaggaa aagcagcaca gtcatgaaag aaggatggat ggtccactac
1381 accagcaagg acacgctgcg gaaacggcac tattggagat tggatagcaa atgtattacc
1441 ctctttcaga atgacacagg aagcaggtac tacaaggaaa ttcctttatc tgaaattttg
1501 tctctggaac cagtaaaaaa ttcagcttta attcctaatt gggccaatcc tcattgtttc

```

## Certificate of Analysis

```
1561 gaaatcacta cggcaaagt agtgtattat gtgggagaaa atgtgggtcaa tccttccagc
1621 ccatcaccaa ataacagtgt tctcaccagt ggcgttggtg cagatgtggc caggatgtgg
1681 gagatagcca tccagcatgc ctttatgcc gtcattccca agggctcctc cgtgggtaca
1741 ggaaccaact tgcacagaga tatctctgtg agtatttcag tatcaaattg ccagattcaa
1801 gaaaatgtgg acatcagcac agtatatcag atttttcctg atgaagtact gggttctgga
1861 cagtttgaa ttgtttatgg aggaaaacat cgtaaaacag gaagagatgt agctattaaa
1921 atcattgaca aattacgatt tccaacaaaa caagaaagcc agcttcgtaa tgaggttgca
1981 attctacaga accttcatca ccctggtggt gtaaatttgg agtgtatggt tgagacgcct
2041 gaaagagtgt ttgttggtat ggaaaaactc catggagaca tgctggaaat gatcttgtca
2101 agtgaaaagg gcaggttgcc agagcacata acgaagtttt taattactca gatactcgtg
2161 gctttgcggc accttcattt taaaaatc gttcactgtg acctcaaacc agaaaatgtg
2221 ttgctagcct cagctgatcc ttttctcag gtgaaacttt gtgattttgg ttttgcccgg
2281 atcattggag agaagtcttt ccggagggtca gtgggtgggta cccccgctta cctggctcct
2341 gaggtcctaa ggaacaagg ctacaatcgc tctctagaca tgtggtctgt tggggctcct
2401 atctatgtaa gcctaagcgg cacattccca tttaatgaag atgaagacat acacgaccaa
2461 attcagaatg cagctttcat gtatccacca aatccctgga aggaaatc tcatgaagcc
2521 attgatctta tcaacaattt gctgcaagta aaaatgagaa agcgctacag tgtggataag
2581 accttgagcc accttggct acaggactat cagacctggt tagatttgcg agagctggaa
2641 tgcaaaatcg gggagcgcata catcacccat gaaagtgatg acctgaggtg ggagaagtat
2701 gcaggcgagc agcggctgca gtaccccaca cacctgatca atccaagtgc tagccacagt
2761 gacactcctg agactgaaga aacagaaatg aaagccctcg gtgagcgtgt cagcatcctc
2821 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.