

Certificate of Analysis

PKC iota, active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-505, 14-505-K, 14-505M

Parent Lot # 1590317

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 iota expressed in Sf21 insect cells. Purified using Ni²⁺/NTA-agarose. Purity 98.3% by SDS-PAGE and Coomassie blue staining. MW = 68.6kDa.

Specific Activity (Parent lot# 1590317): 828U/mg, where one unit of PKC iota, active activity is defined as 1nmol phosphate incorporated into 50µM PKCtide (ERM₂PRKRQGSVRRRV) per minute at 30°C with a final ATP concentration of 100µM.

Formulation: 1.453mg/ml of enzyme in 20mM HEPES/NaOH pH7.4, 250mM NaCl, 50% glycerol, 5mM DTT, 2mM EDTA, 2mM EGTA, 0.05% Triton X-100. Frozen solution.

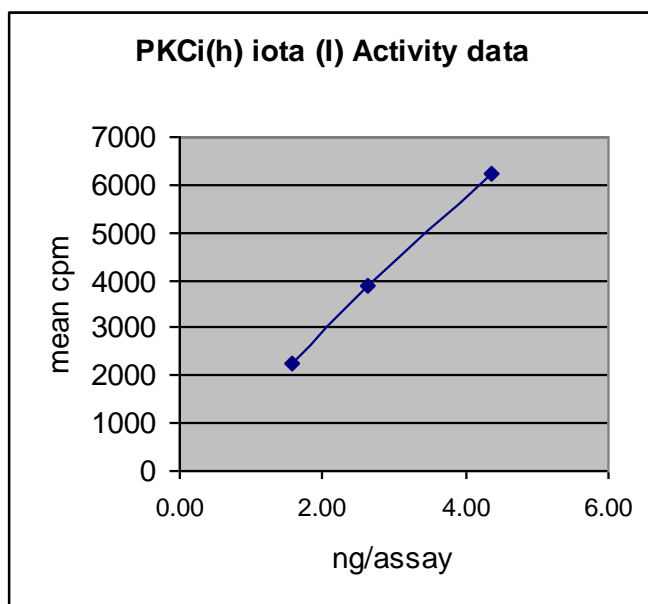
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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.

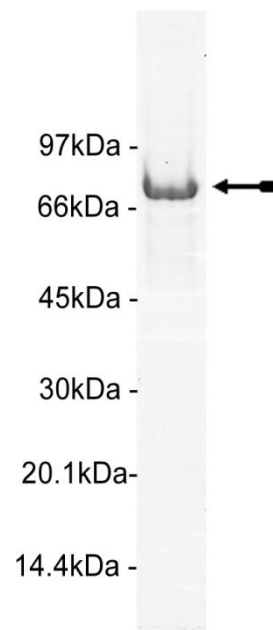
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 1.6–4.4ng of this lot of enzyme phosphorylated 50µM 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 product identity as PKC iota with the translated sequence listed on page three.



SDS-PAGE and Coomassie Stain: Purity was assessed by SDS-PAGE and Coomassie blue staining using 3µg of PKC iota, active.

Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

- 1. 10 x Reaction Buffer:** 200mM HEPES/NaOH pH7.4, 0.3% Triton X-100.
- 2. PKCtide (ERM_RPRKRQGSVRRRV):** Use at a final assay concentration of 50µM. Make up a 500µM stock. Add 2.5µl of stock per assay point.
- 3. PKC iota, active:** Dilute with 20mM HEPES/NaOH pH7.4, 0.03% Triton X-100. Use 1.6–4.4ng 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 format):

1. Add 2.5µl of 10 x reaction buffer per assay to wells.
2. Add 2.5µl of **PKCtide (ERM_RPRKRQGSVRRRV)**.
3. Add **2.5µl (1.6–4.4ng) PKC iota, active**.
4. Add 7.5µl of dH₂O.
5. Add 10µl of diluted [γ -³³P]ATP mixture.
6. Incubate for 10minutes 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 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 iota Sequence Information

<u>Protein</u>	Human PKC iota
<u>Tags</u>	N-terminal 6His
<u>Native sequence</u>	M10of the recombinant protein is equivalent to M1 of human PKC iota
<u>Accession number</u>	GenBank NM_002740

Recombinant PKC iota amino acid sequence:

```

1 MHHHHHHEFM SHTVAGGGSG DSHQVRVKA YYRGDIMITH FEPSISFEGL CNEVRDMCSF
61 DNEQLFTMKW IDEEGDPCTV SSQLELEEF RLYELNKDSE LLIHVFPVCP ERPGMPCPGE
121 DKSIYRRGAR RWRKLYCANG HTFQAKRFNR RAHCAICTDR IWGLGRQGYK CINCKLLVHK
181 KCHKLVITIEC GRHSLPQEPV MPMDQSSMHS DHAQTVIPYN PSSHESLDQV GEEKEAMNTR
241 ESGKASSSLG LQDFDLLRVI GRGSYAKVLL VRLKKTDRIV AMKVVKKELV NDDDEDIDWVQ
301 TEKHVFEQAS NHPFLVGLHS CFQTESRLFF VIEYVNGGDL MFHMQRQRKL PEEHARFYSA
361 EISLALNYLH ERGIYRDLK LDNVLLDSEG HIKLTDYGMC KEGLRPGDIT STFCGTPNYI
421 APEILRGEDY GFSVDWWALG VLMFEMMAGR SPFDIVGSSD NPDQNTEDYL FQVILEKQIR
481 IPRSLSVKAA SVLKSFLNKD PKERLGCHPQ TGFADIQGHP FFRNVDWDMM EQKQVPPFK
541 PNISGEFGLD NFDQSFTNEP VQLTPDDDDI VRKIDQSEFE GFEYINPLLM SAEECV
  
```

Recombinant PKC iota nucleotide sequence:

```

1 atgcatcatc accatcacca tgaattcatg tcccacacgg tcgcaggcgg cggcagcggg
61 gaccattccc accaggtccg ggtgaaagcc tactaccgcg gggatatcat gataacacat
121 tttgaacctt ccattctcct tgaggccctt tgcaatgagg ttcgagacat gtgttctttt
181 gacaacgaac agctcttcac catgaaatgg atagatgagg aaggagaccg gtgtacagta
241 tcatctcagt tggagttaga agaagccttt agactttatg agctaaacaa ggattctgaa
301 ctcttgattc atgtgttccc ttgtgtacca gaacgtcctg ggatgccttg tccaggagaa
361 gataaatcca tctaccgtag aggtgcacgc cgctggagaa agctttattg tgccaatggc
421 cacactttcc aagccaagcg tttcaacagg cgtgctcact gtgccatctg cacagaccga
481 atatggggac ttggacgcca aggatataag tgcatacaact gcaaactctt ggttcataag
541 aagtgcata aactcgtcac aattgaatgt gggcggcatt ctttgccaca ggaaccagtg
601 atgccatgg atcagtcac atcgattct gaccatgcac agacagtaat tccatataat
661 cttcaagtc atgagagttt ggatcaagtt ggtgaagaaa aagaggcaat gaacaccagg
721 gaaagtggca aagcttcac cagtctaggt cttcaggatt ttgatttgct ccgggtaata
781 ggaagaggaa gttatgccaa agtactgttg gttcgattaa aaaaaacaga tcgtatttat
841 gcaatgaaag ttgtgaaaaa agagcttggt aatgatgatg aggatattga ttgggtacag
901 acagagaagc atgtgtttga gcaggcatcc aatcatcctt tccttgttgg gctgcattct
961 tgctttcaga cagaaagcag attgttcttt gttatagagt atgtaaattg aggagacctt
1021 atgtttcata tgcagcgaca aagaaaactt cctgaagaac atgccagatt ttactctgca
1081 gaaatcagtc tagcattaaa ttatcttcat gagcgaggga taatttatag agatttgaaa
1141 ctggacaatg tattactgga ctctgaaggc cacattaaac tctactgacta cggcatgtgt
1201 aaggaaggat tacggccagg agatacaacc agcactttct gtggtactcc taattacatt
1261 gctcctgaaa ttttaagagg agaagattat ggtttcagtg ttgactggtg ggctcttgga
1321 gtgctcatgt ttgagatgat ggcaggaagg tctccatttg atattggttg gagctccgat
1381 aacctgacc agaacacaga ggattatctc ttccaagtta ttttgaaaaa acaaattcgc
1441 ataccagtt ctctgtctgt aaaagctgca agtgttctga agagttttct taataaggac
1501 cctaaggaac gattgggttg tcatcctcaa acaggatttg ctgatattca gggacaccgg
1561 ttcttccgaa atgttgattg ggatgatgat gagcaaaaac aggtggtacc tccctttaa
1621 ccaaatatth ctggggaatt tggtttggac aactttgatt ctcagtttac taatgaacct
  
```

Certificate of Analysis

1681 gtccagctca ctccagatga cgatgacatt gtgaggaaga ttgatcagtc tgaattttaa
1741 ggttttgagt atatcaatcc tcttttgatg tctgcagaag aatgtgtctg a

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