

## Certificate of Analysis

### CaM Kinase II $\delta$ , active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-723, 14-723-K, 14-723M

Parent Lot # 1838058

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, full length CaM Kinase II $\delta$ , expressed by baculovirus in Sf21 insect cells. Purified using Ni<sup>2+</sup>/NTA agarose. Purity 90.3% by SDS-PAGE and Coomassie blue staining. MW = 58kDa.

**Specific Activity (Parent lot# 1838058):** 16527U/mg, where one unit of CaM Kinase II $\delta$  activity is defined as 1nmol phosphate incorporated into 250 $\mu$ M (KKLNRTL~~S~~FAEPG) per minute at 30°C with a final ATP concentration of 100 $\mu$ M.

**Formulation:** 0.38mg/ml of enzyme in 50mM Tris/HCl pH7.5, 300mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 1mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution.

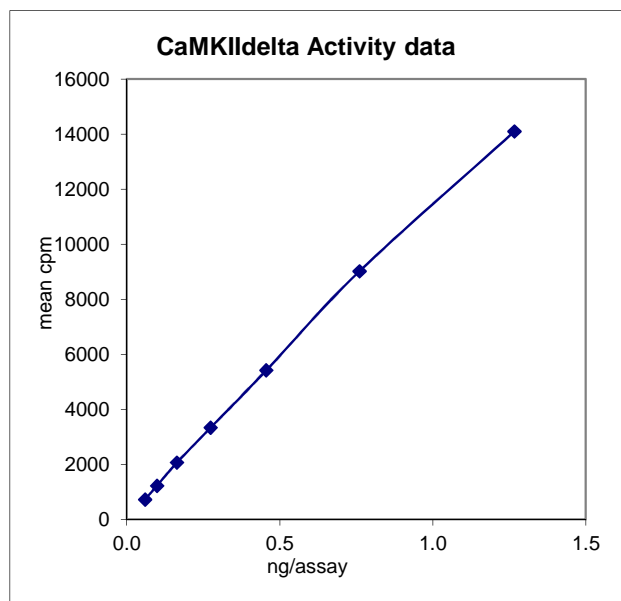
**Storage and Stability:** On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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

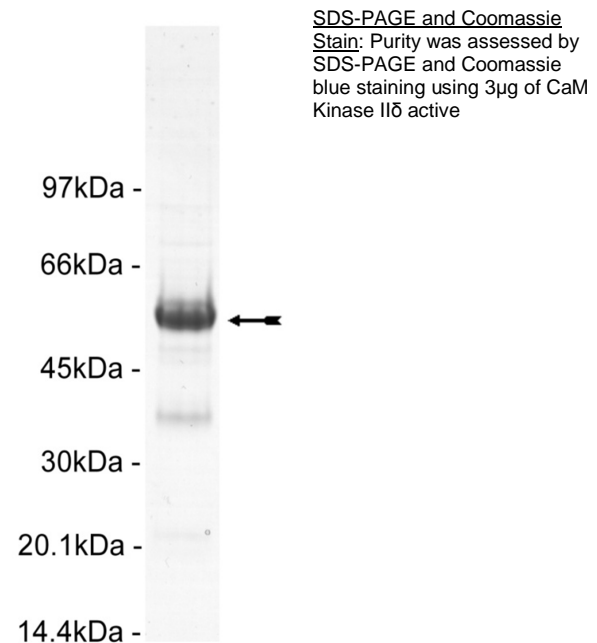
**FOR IN VITRO RESEARCH USE ONLY  
NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**Kinase Assay:** 0.1–1.3ng of this lot of enzyme phosphorylated 250 $\mu$ M (KKLNRTL~~S~~FAEPG) 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 CaM Kinase II $\delta$  with the translated native sequence listed on page three.



## Certificate of Analysis

### Kinase Assay Protocol

#### Stock Solutions:

1. **5 x Reaction Buffer:** 40mM MOPS-NaOH pH7.0, 1mM EDTA.
2. **(KCLNRTL SFAEPG):** Use at a final concentration of 250 $\mu$ M. Make up a 2.5mM stock. Add 2.5 $\mu$ l of stock per assay point.
3. **CaCl<sub>2</sub>:** Use at a final assay concentration of 500 $\mu$ M. Make up a 5mM stock. Add 2.5 $\mu$ l of stock per assay point.
4. **Calmodulin:** Use at a final assay concentration of 1 $\mu$ M. Make up a 0.3mg/ml stock. Add 1.33 $\mu$ l of stock per assay point.
5. **CaM Kinase II $\delta$ , active:** Dilute with 20mM MOPS-NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 0.1–1.3ng per assay point.
6. **[ $\gamma$ -<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 5 $\mu$ l of 5 x reaction buffer per assay to wells.
2. Add 2.5 $\mu$ l of **(KCLNRTL SFAEPG)**.
3. Add **2.5 $\mu$ l (0.1–1.3ng) CaM Kinase II $\delta$  active**.
4. Add 1.17 $\mu$ l of dH<sub>2</sub>O.
5. Add 2.5 $\mu$ l of CaCl<sub>2</sub>.
6. Add 1.33 $\mu$ l of Calmodulin.
7. Add 10 $\mu$ l of diluted [ $\gamma$ -<sup>33</sup>P]ATP mixture.
8. Incubate for 10 minutes at 30°C.
9. Stop the reaction by adding 5 $\mu$ l of 3% phosphoric acid.
10. Transfer a 10 $\mu$ l aliquot onto the appropriate area of a **P30 Filtermat**.
11. Wash the filtermat three times for 5 minutes with 75mM phosphoric acid.
12. Wash the filtermat once for 2 minutes with methanol.
13. Transfer the filtermat to a sealable plastic bag and add 4ml of scintillation cocktail.
14. 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

### CaM Kinase II $\delta$ Sequence Information

<b>Protein</b>	Human CaM Kinase II $\delta$
<b>Tags</b>	N-terminal 6His
<b>Native sequence</b>	M31 of recombinant protein is equivalent to M1 of human CaM Kinase II $\delta$
<b>Accession number</b>	GenBank NM_172115

#### Recombinant CaM Kinase II $\delta$ amino acid sequence:

```

1 MSYYHHHHHH DYDIPTTENL YFQGAMDPEF MASTTTCTRF TDEYQLFEEL GKGAFSVVRR
61 CMKIPTGQEY AAKIINTKKL SARDHQKLER EARICRLKHK PNIVRLHDSI SEEGFHLYLVF
121 DLVTGGELFE DIVAREYYS EADASHCIQQI LESVNHCHLN GIVHRDLKPE NLLLASKSKG
181 AAVKLADFGL AIEVQGDQQA WFGFAGTPGY LSPEVLRKDP YGKPVDMWAC GVILYILLVG
241 YPPFWDEDQH RLYQQIKAGA YDFPSPEWDT VTPEAKDLIN KMLTINPAKR ITASEALKHP
301 WICQRSTVAS MMHRQETVDC LKKFNARRKL KGAILTTMLA TRNFSAAKSL LKKPDGVKES
361 TESSNTTIED EDVKARKQEI IKVTEQLIEA INNGDFEAYT KICDPGLTAF EPEALGNLVE
421 GMDFHRFYFE NALSKSNKPI HTIILNPHVH LVGDDAACIA YIRLTQYMDG SGMPKTMQSE
481 ETRVWHRRDG KWQNVHFHRS GSPTVPIK
  
```

#### Recombinant CaM Kinase II $\delta$ nucleotide sequence:

```

1 atgtcgtact accatcacca tcaccatcac gattacgata tcccaacgac cgaaaacctg
61 tattttcagg gcgccatgga tccggaattc atggcttcga ccacaacctg caccaggttc
121 acggacgagt atcagctttt cggaggactt ggaaaggggg cattctcagt ggtgagaaga
181 tgtatgaaaa ttcctactgg acaagaatat gctgccaaaa ttatcaacac caaaaagctt
241 tctgctaggg atcatcagaa actagaaaga gaagctagaa tctgccgtct tttgaagcac
301 cctaataattg tgcgacttca tgatagcata tcagaagagg gctttcacta cttgggtgtt
361 gatttagtta ctggaggtga actgtttgaa gacatagtgg caagagaata ctacagtgaa
421 gctgatgcca gtcattgtat acagcagatt ctagaaagtg ttaatcattg tcacctaaat
481 ggcatagttc acagggacct gaagcctgag aatttgcttt tagctagcaa atccaagggg
541 cgacgtgtga aattggcaga ctttggctta gccatagaag ttcaagggga ccagcaggcg
601 tggtttggtt ttgctggcac acctggatat ctttctccag aagttttacg taaagatcct
661 tatggaaagc cagtggatat gtgggcattg ggtgtcattc tctatatct acttgtggg
721 tatccaccct tctgggatga agaccaacac agactctatc agcagatcaa ggctggagct
781 tatgattttc catcaccaga atgggacacg gtgactcctg aagccaaaga cctcatcaat
841 aaaatgctta ctatcaacc tgccaaacgc atcacagcct cagaggcact gaagcaccca
901 tggatctgtc aacgttctac tgttgcttcc atgatgcaca gacaggagac tgtagactgc
961 ttgaagaaat ttaatgctag aagaaaacta aagggtgcca tcttgacaac tatgctggct
1021 acaaggaatt tctcagcagc caagagtttg ttgaagaaac cagatggagt aaaggagtca
1081 actgagagtt caaatacaac aattgaggat gaagatgtga aagcacgaaa gcaagagatt
1141 atcaaagtca ctgaacaact gatcgaagct atcaacaatg gggactttga agcctacaca
1201 aaaatctgtg acccaggcct tactgctttt gaacctgaag ctttgggtaa tttagtgaa
1261 gggatggatt ttcaccgatt ctactttgaa aatgctttgt ccaaaagcaa taaaccaatc
1321 cacactatta ttctaaacc tcatgtacat ctggtagggg atgatgccgc ctgcatagca
1381 tatattaggc tcacacagta catggatggc agtggaatgc caaagacaat gcagtcagaa
1441 gagactcgtg tgtggcaccg ccgggatgga aagtggcaga atgttcattt tcatcgctcg
1501 ggtcaccaa cagtacccat caagtaa
  
```

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.