

## Certificate of Analysis

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

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-718, 14-718-K, 14-718M

Parent Lot # 31119U

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, CaM Kinase II $\beta$ , amino acids 1–315, expressed by baculovirus in Sf21 insect cells. Purified using Ni<sup>2+</sup>/NTA agarose. Purity 86% by SDS-PAGE and Coomassie blue staining. MW = 39.6kDa.

**Specific Activity (Parent lot# 31119U):** 14744U/mg, where one unit of CaM Kinase II $\beta$  activity is defined as 1nmol phosphate incorporated into 250 $\mu$ M (KKLNRTL $\beta$ FAEPG) per minute at 30°C with a final ATP concentration of 100 $\mu$ M.

**Formulation:** 3.6mg/ml of enzyme in 50mM Tris/HCl pH7.5, 300mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 1mM benzimidazole, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution.

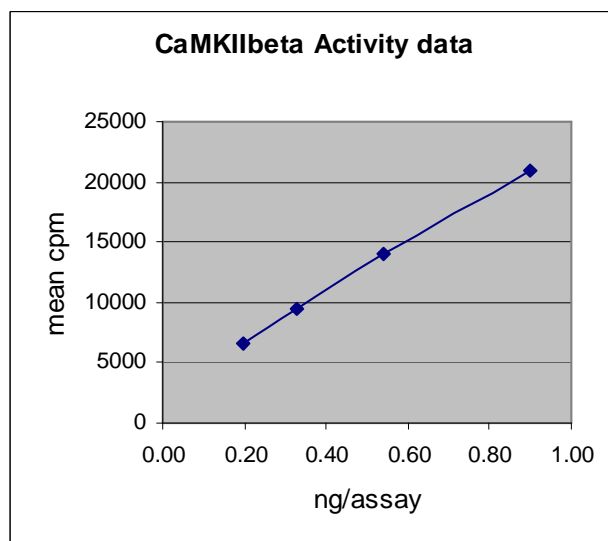
**Storage and Stability:** On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

**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.

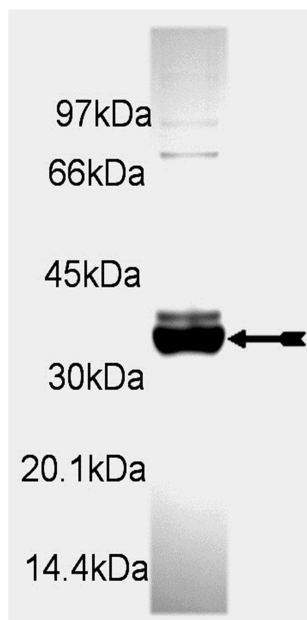
**FOR IN VITRO RESEARCH USE ONLY  
NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**Kinase Assay:** 0.2–0.9ng of this lot of enzyme phosphorylated 250 $\mu$ M (KKLNRTL $\beta$ 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 $\beta$  with the translated native sequence listed on page three.



**SDS-PAGE and Coomassie Stain:** Purity was assessed by SDS-PAGE and Coomassie blue staining using 3 $\mu$ g of CaM Kinase II $\beta$  , active

## Certificate of Analysis

### Kinase Assay Protocol

#### Stock Solutions:

1. **5 x Reaction Buffer:** 40mM MOPS-NaOH pH7.0, 1mM EDTA.
2. **(KKLNRTLSFAEPG):** Use at a final concentration of 250 $\mu$ M. Make up a 2.5mM stock. Use 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 $\beta$ , active:** Dilute with 20mM MOPS-NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 0.2–0.9ng 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 **(KKLNRTLSFAEPG)**.
3. Add **2.5 $\mu$ l (0.2–0.9ng) CaM Kinase II $\beta$  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 $\beta$ Sequence Information

<b><u>Protein</u></b>	Human CaM Kinase II $\beta$
<b><u>Tags</u></b>	N-terminal 6His
<b><u>Native sequence</u></b>	M31 of the recombinant protein is equivalent to M1 of human CaM Kinase II $\beta$
<b><u>Accession number</u></b>	GenBank AF081572

#### ***Recombinant CaM Kinase II $\beta$ amino acid sequence:***

```

1 MSYYHHHHHH DYDIPTTENL YFQGAMDPEF MATTVTCTRF TDEYQLYEDI GKGAFSVVR
61 CVKLCTGHEY AAKIINTKKL SARDHQKLER EARICRLLKH SNIVRLHDSI SEEGFHYLVF
121 DLVTGGELFE DIVAREYSE ADASHCIQQI LEAVLHCHQM GVVHRDLKPE NLLLASKCKG
181 AAVKLADDFGL AIEVQGDQQA WFGFAGTPGY LSPEVLRKEA YGKPVDIWAC GVILYILLVG
241 YPPFWDEDQH KLYQQIKAGA YDFPSPewDT VTPEAKNLIN QMLTINPAKR ITAHEALKHP
301 WVCQRSTVAS MMHRQETVEC LKKFNARRKL KGAILTTMLA TRNFS

```

#### ***Recombinant CaM Kinase II $\beta$ nucleotide sequence:***

```

1 atgtcgtact accatcacca tcaccatcac gattacgata tccaacgac cgaaaacctc
61 tattttcagg ggcctatgga tccggaattc atggccacca cggtagactg caccgccttc
121 accgacgagt accagctcta cgaggatatt ggcaaggggg ctttctctgt ggtccgacgc
181 tgtgtcaagc tctgcaccgg ccatgagtat gcagccaaga tcatcaacac caagaagctg
241 tcagccagag atcaccagaa gctggagaga gaggctcgga tctgccgcct tctgaagcat
301 tccaacatcg tgcgtctcca cgacagcatc tccgaggagg gcttccacta cctggctctc
361 gatctgggtca ctgggtgggga gctctttgaa gacattgtgg cgagagagta ctacagcgag
421 gctgatgcca gtcactgtat ccagcagatc ctggaggccg ttctccattg tcaccaaatg
481 ggggtcgtcc acagagacct caagccggag aacctgcttc tggccagcaa gtgcaaaggg
541 gctgcagtga agctggcaga cttcggccta gctatcgagg tgcaggggga ccagcaggca
601 tggtttggtt tcgctggcac accaggctac ctgtcccctg aggtccttcg caaagaggcg
661 tatggcaagc ctgtggacat ctgggcatgt ggggtgatcc tgtacatcct gctcgtgggc
721 taccaccctt tctgggacga ggaccagcac aagctgtacc agcagatcaa ggctgggtgc
781 tatgacttcc cgtcccctga gtgggacacc gtcactcctg aagccaaaaa cctcatcaac
841 cagatgctga ccatcaacc tgccaagcgc atcacagccc atgaggcctt gaagcaccgc
901 tgggtctgcc aacgctccac ggtagcatcc atgatgcaca gacaggagac tgtggagtgt
961 ctgaaaaggt tcaatgccag gagaaagctc aaggaggcca tcctcaccac catgctggcc
1021 acacggaatt tctcataa

```

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.