

Certificate of Analysis

CaM Kinase I γ , active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-719, 14-719-K, 14-719M

Parent Lot # D8NN026U

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 I γ , amino acids 1–330, expressed by baculovirus in Sf21 insect cells. Purified using Ni²⁺/NTA agarose. Purity 91.2% by SDS-PAGE and Coomassie blue staining. MW = 41.2kDa.

Specific Activity (Parent lot# D8NN026U): 20010U/mg, where one unit of CaM Kinase I γ activity is defined as 1nmol phosphate incorporated into 250 μ M (KKLNRTL γ FAEPG) per minute at 30°C with a final ATP concentration of 100 μ M.

Formulation: 3.885mg/ml of enzyme in 50mM Tris/HCl pH8.0, 300mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 1mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution.

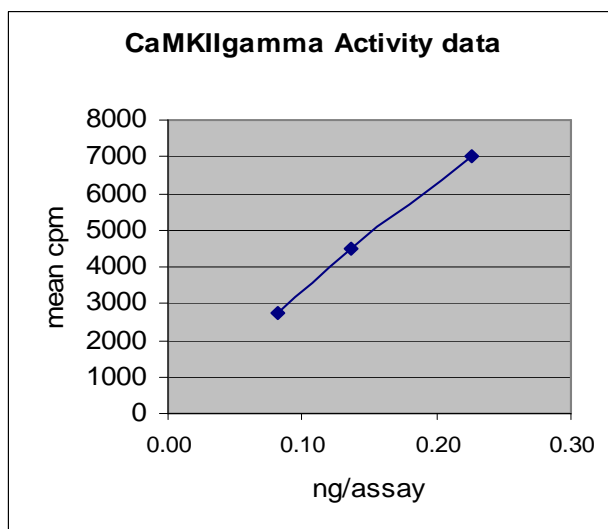
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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**

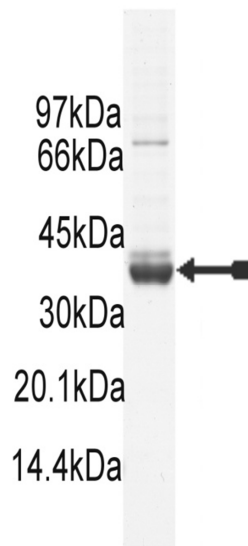
Quality Control Testing

Kinase Assay: 0.08–0.23ng of this lot of enzyme phosphorylated 250 μ M (KKLNRTL γ 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 I γ 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 μ g of CaM Kinase I γ , 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 250µM. Make up a 2.5mM stock. Use 2.5µl of stock per assay point.
3. **CaCl₂:** Use at a final assay concentration of 500µM. Make up a 5mM stock. Add 2.5µl of stock per assay point.
4. **Calmodulin:** Use at a final assay concentration of 1µM. Make up a 0.3 mg/ml Stock. Add 1.33µl of stock per assay point.
5. **CaM Kinase II γ , active:** Dilute with 20mM MOPS-NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 0.08–0.23ng per assay point.
6. **[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 5µl of 5 x reaction buffer per assay to wells.
2. Add 2.5µl of **(KKLNRTLSFAEPG)**.
3. Add **2.5µl (0.08–0.23ng) CaM Kinase II γ active**.
4. Add 1.17µl of dH₂O.
5. Add 2.5µl of CaCl₂.
6. Add 1.33µl of calmodulin.
7. Add 10µl of diluted [γ -³³P]ATP mixture.
8. Incubate for 10 minutes at 30°C.
9. Stop the reaction by adding 5µl of 3% phosphoric acid.
10. Transfer a 10µl 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 1µl of 30% phosphoric acid.

Certificate of Analysis

CaM Kinase Ily Sequence Information

<u>Protein</u>	Human CaM Kinase Ily
<u>Tags</u>	N-terminal 6His
<u>Native sequence</u>	M31 of the recombinant protein is equivalent to M1 of CaM Kinase Ily
<u>Accession number</u>	GenBank NM_172171

Recombinant CaM Kinase Ily amino acid sequence:

```

1 MSYYHHHHHH DYDIPTTENL YFQGAMDPEF MATTATCTRF TDDYQLFEEL GKGAFSVVR
61 CVKKTSTQEY AAKIINTKKL SARDHQKLER EARICRLLKH PNIVRLHDSI SEEGFHYLVF
121 DLVTGGELFE DIVAREYYSE ADASHCIHQI LESVNHIIHQH DIVHRDLKPE NLLLASKCKG
181 AAVKLADFGF AIEVQGEQQA WFGFAGTPGY LSPEVLRKDP YGKPVDIWAC GVILYILLVG
241 YPPFWDEDQH KLYQQIKAGA YDFPSPEWDT VTPEAKNLIN QMLTINPAKR ITADQALKHP
301 WVCQRSTVAS MMHRQETVEC LRKFNARRKL KGAILTTMLV SRNFSAAKSL LNKKSDGGVK
  
```

Recombinant CaM Kinase Ily nucleotide sequence:

```

1 atgtcgact accatcacca tcaccatcac gattacgata tcccaacgac cgaaaacctg
61 tattttcagg gcgccatgga tccggaattc atggccacca cgcaccactg caccggttcc
121 accgacgact accagctctt cgaggagcct ggcaaggggtg ctttctctgt ggtccgcagg
181 tgtgtgaaga aaacctccac gcaggagtag gcagcaaaaa tcatcaatac caagaagttg
241 tctgcccggg atcaccagaa actagaacgt gaggctcgga tatgtcgact tctgaaacat
301 ccaaaccatcg tgcgcctcca tgacagtatt tctgaagaag ggtttacta cctcgtgttt
361 gaccttgtaa ccggcgggga gctgtttgaa gacattgtgg ccagagagta ctacagtgaa
421 gcagatgcca gccactgtat acatcagatt ctggagagtg ttaaccacat ccaccagcat
481 gacatcgacc acagggacct gaagcctgag aacctgctgc tggcgagtaa atgcaagggg
541 gccgccgtca agctggctga ttttggccta gccatcgaag tacagggaga gcagcaggct
601 tggtttggtt ttgctggcac ccagggttac ttgtcccctg aggtcttgag gaaagatccc
661 tatggaaaac ctgtggatat ctgggcctgc ggggtcatcc tgtatatcct cctggggggc
721 taccctccct tctgggatga ggatcagcac aagctgtatc agcagatcaa ggctggagcc
781 tatgatattc catcaccaga atgggacacg gtaactcctg aagccaagaa cttgatcaac
841 cagatgctga ccataaacc agcaaagcgc atcacggctg accaggctct caagcaccgc
901 tgggtctgtc aacgatccac ggtggcatcc atgatgcatc gtcaggagac tgtggagtgt
961 ttgcgcaagt tcaatgcccg gagaaaactg aagggtgcca tctcacgac catgcttgtc
1021 tccaggaact tctcagctgc caaaagccta ttgaacaaga agtcggatgg cgggtgtcaag
1081 taa
  
```

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.