

Certificate of Analysis

CK1 ϵ , active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 16-045, 16-045-K, 16-045M

Parent Lot # D17NP008N

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 GST-tagged, recombinant, human CK1 ϵ amino acids 1-348 expressed by baculovirus in Sf21 insect cells. Purified using glutathione agarose followed by size exclusion chromatography.

Purity 85% by SDS-PAGE and Coomassie blue staining. MW = 67kDa.

Specific Activity (Parent lot# D17NP008N): 1462U/mg, where one unit of CK1 ϵ activity is defined as 1nmol phosphate incorporated into 200 μ M KRRRALS(p)VASLPGL per minute at 30°C with a final ATP concentration of 100 μ M.

Formulation: 0.17mg/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.

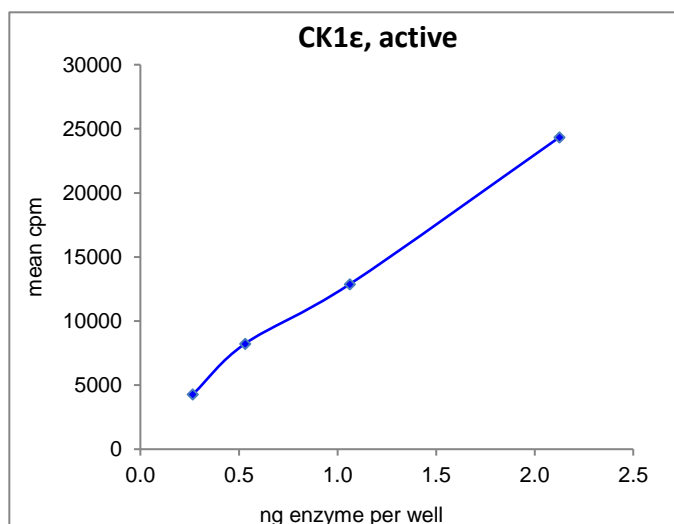
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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

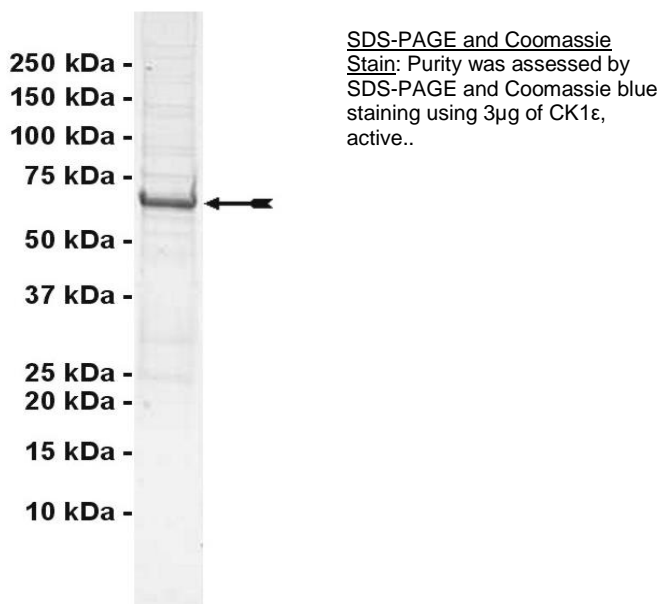
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 0.3–2.1ng of this lot of enzyme phosphorylated 200 μ M KRRRALS(p)VASLPGL 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 CK1 ϵ with the translated 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. KRRRALS(p)VASLPGL:** Use at a final assay concentration of 200 μ M. Prepare a 2mM stock and add 2.5 μ l of stock per assay point.
- 3. CK1 ϵ , active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 0.3–2.1ng per assay point.
- 4. [γ -³³P]ATP:** 2.5 x MgAc/[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 5 μ l of 5 x reaction buffer per assay to wells.
2. Add 2.5 μ l of KRRRALS(p)VASLPGL.
3. Add **2.5 μ l (0.3–2.1ng) CK1 ϵ , active.**
4. Add 5 μ l of dH₂O.
5. Add 10 μ l of diluted [γ -³³P]ATP mixture.
6. Incubate for 30 minutes at 30°C.
7. Stop the reaction by adding 5 μ l of 3% phosphoric acid.
8. Transfer a 10 μ l 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 1 μ l of 30% phosphoric acid.

Certificate of Analysis

CK1ε, active Sequence Information

Protein	Human CK1ε
Tags	N-terminal GST
Native sequence	M230 of the recombinant protein is equivalent to M1 of human CK1ε
Accession number	GenBank NM_001894.4

Recombinant CK1ε amino acid sequence:

```
1  MSPILGYWKI  KGLVQPTRL  LEYLEEKYEE  HLYERDEGDK  WRNKKFELGL  EFPNLPYYID
61  GDVKLTQSM  IIRYIADKHN  MLGGCPKERA  EISMLEGAVL  DIRYGVSRIA  YSKDFETLKV
121 DFLSKLPEML  KMFEDRLCHK  TYLNGDHVTH  PDFMLYDALD  VVLYMDPMCL  DAFPKLVCFK
181 KRIEAIPOID  KYLKSSKYIA  WPLQGWQATF  GGGDHPPKSD  LVPRGSKELM  ELRVGNKYRL
241 GRKIGSGSFG  DIYLGANIAS  GEEVAIKLEC  VKTKHPQLHI  ESKFYKMMQG  GVGIPSIKWC
301 GAEGDYNVMV  MELLGPSLED  LFNFCSRKFS  LKTVLLLDQ  MISRIEYIHS  KNFIHRDVKP
361 DNFLMGLGKK  GNLVYIIDFG  LAKKYRDART  HQHIPYRENK  NLTGTARYAS  INTHLGIEQS
421 RRDDLESGLY  VLMYFNLGSL  PWQGLKAATK  RQKYERISEK  KMSTPIEVL  KGYPSEFSTY
481 LNFCSRSLRF  DKPDYSYLRQ  LFRNLFHRQG  FSYDYVFDWN  MLKFGAARNP  EDVDRERREH
541 EREERMGQLR  GSATRALPPG  PPTGATANRL  RSAAEPV
```

Recombinant CK1ε nucleotide sequence:

```
1  atgtccccta  tactaggtta  ttgaaaatt  aagggccttg  tgcaaccac  tcgacttctt
61  ttggaatc  ttgaagaaa  atatgaagag  catttgatg  agcgcgatga  aggtgataaa
121  tggcgaaaca  aaaagtttga  attgggttg  gagtttcca  atcttcctta  ttatattgat
181  ggtgatgtta  aattaacaca  gtctatggcc  atcatacggt  atatagctga  caagcacaac
241  atgttgggtg  gttgtccaaa  agagcgtgca  gagatttcaa  tgcttgaagg  agcggttttg
301  gatattagat  acggtgtttc  gagaattgca  tatagtaaag  actttgaaac  tctcaaagtt
361  gattttctta  gcaagctacc  tgaatgctg  aaaatgttcg  aagatcgttt  atgtcataaa
421  acatatttaa  atggtgatca  tgtaaccat  cctgacttca  tgttgatga  cgctcttgat
481  gttgttttat  acatggacc  aatgtgcctg  gatgcgttcc  caaaattagt  ttgttttaa
541  aaacgtattg  aagctatccc  acaaattgat  aagtacttga  aatccagcaa  gtatatagca
601  tggcctttgc  agggctggca  agccacgttt  ggtggtggcg  accatcctcc  aaaatcggat
661  ctggttccgc  gtggatccaa  ggaattgatg  gagctacgtg  tggggaacaa  gtaccgcctg
721  ggacggaaga  tcgggagcgg  gtccttcgga  gatattctac  tgggtgccaa  catcgcctct
781  ggtgaggaag  tcgccatcaa  gctggagtgt  gtgaagacaa  agcaccacca  gctgcacatc
841  gagagcaagt  tctacaagat  gatgcagggt  ggcgtgggga  tcccgtccat  caagtgggtg
901  ggagctgagg  gcgactacaa  cgtgatggtc  atggagctgc  tggggcctag  cctcgaggac
961  ctggtcaact  tctgttccc  caaattcagc  ctcaagacgg  tgctgctctt  ggccgaccag
1021 atgatcagcc  gcatcgagta  tatccactcc  aagaacttca  tccaccggga  cgtcaagccc
1081 gacaacttcc  tcatggggtc  ggggaagaag  gggcaactgg  tctacatcat  cgactctggc
1141 ctggccaaga  agtaccggga  cgcccgcacc  caccagcaca  tccctaccg  ggaaaacaag
1201 aacctgaccg  gcacggcccg  ctacgcttcc  atcaacacgc  acctgggcat  tgagcaaagc
1261 cgtcgagatg  acctggagag  cctgggctac  gtgctcatgt  acttcaacct  gggctccctg
1321 ccctggcagg  ggctcaaagc  agccaccaag  cgccagaagt  atgaacggat  cagcgagaag
1381 aagatgtcaa  cgcccatcga  ggtcctctgc  aaaggctatc  cctccgaatt  ctcaacatac
```

Certificate of Analysis

```
1441 ctcaacttct gccgctccct gcggtttgac gacaagcccg actactetta cctacgtcag
1501 ctcttccgca acctcttcca ccggcagggc ttctcctatg actacgtctt tgactggaac
1561 atgctgaaat tcggtgcagc ccggaatccc gaggatgtgg accgggagcg gcgagaacac
1621 gaacgcgagg agaggatggg gcagctacgg gggtcgcgca cccgagccct gccccctggc
1681 ccaccacagg gggccactgc caaccggctc cgcagtgccg ccgagcccg gtaa
```

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.