

Certificate of Analysis

Casein Kinase 1 δ , active

(Recombinant enzyme expressed in *E.coli* cells)

Item # 14-520, 14-520-K, 14-520M

Parent Lot # D8KN032U

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 casein kinase 1 delta, amino acids 1–294, expressed in *E.coli* cells. Purified using glutathione-agarose. Purity 56.4% by SDS-PAGE and Coomassie blue staining. MW = 61kDa.

Specific Activity (Parent lot# D8KN032U): 2407U/mg, where one unit of CK1delta activity is defined as 1nmol phosphate incorporated into 200 μ M (KRRRALS(p)VASLPGL) (CK1tide) per minute at 30°C with a final ATP concentration of 100 μ M.

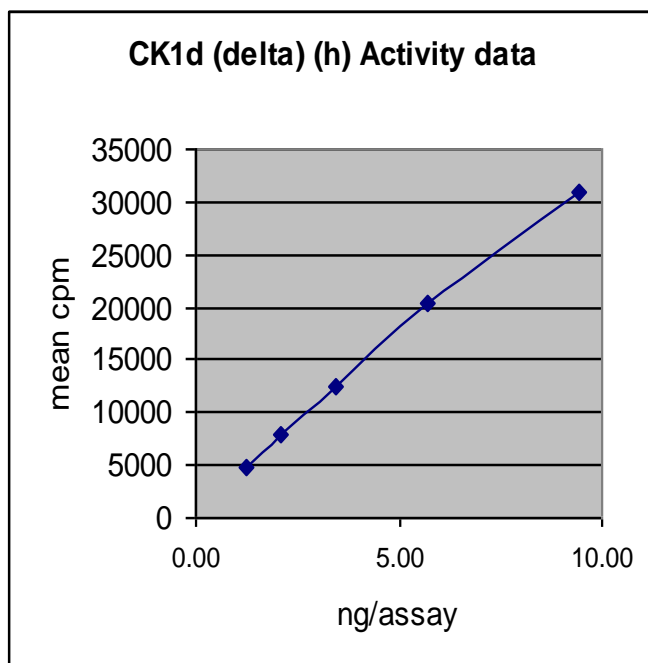
Formulation: 2.265mg/ml of enzyme in 50mM Tris/HCl, pH7.5, 150mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 50% glycerol, 0.2mM PMSF, 1mM benzamidine, 0.1 % 2-mercaptoethanol. Liquid at -20°C.

Storage and Stability: On receipt of material store at -20°C. Unopened reagent is stable for a minimum of 6 months 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.

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

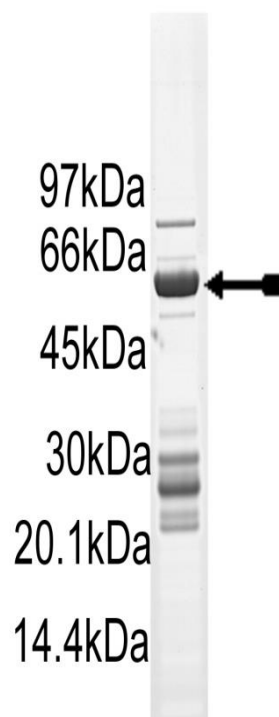
Quality Control Testing

Kinase Assay: 1.2–9.4ng 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 product identity as Casein Kinase 1 delta with the translated sequence listed on page three.

SDS-PAGE and Coomassie Stain: Representative gel from this lot. Purity was assessed by SDS-PAGE and Coomassie blue staining using 3 μ g of casein kinase 1 delta, active.



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. Use 2.5 μ l of stock per assay point.
3. **Casein Kinase 1 δ active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 1.2–9.4ng 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 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 (1.2–9.4ng) casein kinase 1 δ , active.**
4. Add 5 μ l of dH₂O.
5. Add 10 μ l of diluted [γ -³³P] ATP mixture.
6. Incubate for 10 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

Casein kinase 1 delta Sequence Information

Protein	human casein kinase 1 delta (1–294)
Tags	N-terminal GST
Native sequence	M232 of the recombinant protein is equivalent to M1 of human casein kinase 1 delta
Accession number	GenBank AB063114. This is a rat cDNA that is 100% identical to human casein kinase 1 delta over 1-294 of the predicted amino acid sequence.

Recombinant casein kinase 1 delta amino acid sequence:

```

1  MSPILGYWKI  KGLVQPTRLL  LEYLEEKYEE  HLYERDEGDK  WRNKKFELGL  EFPNLPYYID
61  GDVKLTQ SMA  IIRYIADKHN  MLGGCPKERA  EISMLEGAVL  DIRYGVSRIA  YSKDFETLKV
121 DFLSKLPEML  KMFEDRLCHK  TYLNGDHVTH  PDFMLYDALD  VVLYMDPMCL  DAFPKLVCFK
181 KRIEAI PQID  KYLKSSKYIA  WPLQGWQATF  GGGDHPPKSD  LVPRGSPGIR  LMELRVGNRY
241 RLGRKIGSGS  FGDIYLGTDI  AAGEEVAIKL  ECVKTKHPQL  HIESKIYKMM  QGGVGIPTIR
301 WCGAEGDYNV  VMPELLGPSL  EDLNFCSRK  FSLKTVLLLA  DQMISRIEYI  HSKNFIHRDV
361 KPDNFLMGLG  KKGNLVYIID  FGLAKKYRDA  RTHQHIPPYR  E  NKNLTGTARY  ASINTHLGIE
421 QSRRDDLESL  GYVLMYFNLG  SLPWQGLKAA  TKRQKYERIS  EKKMSTPIEV  LCKGYPSEFA
481 TYLNFCSRSL  FDDKPDYSYL  RQLFRNLFHR  QGFSYDYVFD  WNMLK
  
```

Recombinant casein kinase 1 delta nucleotide sequence:

```

1  atgtccccta  tactaggtta  ttggaaaatt  aagggccttg  tgcaaccac  tcgacttctt
61  ttggaatata  ttgaagaaaa  atatgaagag  catttgatg  agcgcgatga  aggtgataaa
121  tggcgaaaca  aaaagtttga  attgggtttg  gagtttccca  atcttcctta  ttatatgtat
181  ggtgatgtta  aattaacaca  gtctatggcc  atcatacggt  atatactgta  caagcacaac
241  atgttgggtg  gttgtccaaa  agagcgtgca  gagatttcaa  tgcttgaagg  agcggttttg
301  gatattagat  acggtgtttc  gagaattgca  tatagtaaag  actttgaaac  tctcaaagtt
361  gattttctta  gcaagctacc  tgaaatgctg  aaaatgttcg  aagatcgttt  atgtcataaa
421  acataattaa  atggtgatca  tgtaaccat  cctgacttca  tgttgatgta  cgctcttgat
481  gttgttttat  acatggacc  aatgtgcctg  gatgcgttcc  caaaattagt  ttgttttaaa
541  aaacgtattg  aagctatccc  acaaattgat  aagtacttga  aatccagcaa  gtatatagca
601  tggcctttgc  agggctggca  agccacgttt  ggtgggtggc  accatcctcc  aaaatcggat
661  ctggttccgc  gtggatcccc  gggaattcgg  cttatggagc  tgagggtcgg  gaataggtac
721  cgactgggcc  gcaagatcgg  cagcggctcc  ttcggagaca  tctatctcgg  tacggacatt
781  gctgcaggag  aagaagttgc  catcaagctt  gaatgtgtca  aaaccaaaaca  tcctcagctc
841  cacattgaga  gcaagatcta  caaatgatg  cagggaggag  tgggcatccc  taccatcaga
901  tgggtgtggg  ctgaggggga  ctacaatgtc  atggtgatgg  agctactggg  accagcctg
961  gaagacctat  tcaacttctg  ttcaaggaag  tttagtctca  aaactgttct  gttgcttgct
1021 gaccaaata  taagtcgtat  tgagtacatt  cattcgaaga  attttatcca  ccgagacgtg
1081 aaaccagata  acttctcat  ggggctggga  aagaaaggca  acctggtcta  catcattgac
1141 tttgggctgg  ccaagaagta  tcgggatgcc  cgcaccacc  agcatatccc  ctatcgagag
1201 aacaagaacc  tcacaggac  agcacgctat  gcctccatca  acacacacct  tggcattgaa
1261 caatctcgaa  gggatgactt  ggagtctctg  gggtagctgc  tgatgtactt  caacctgggc
1321 tctctcccct  ggcaggggct  gaaggccgcc  accaaaaggc  agaagtatga  gaggatcagt
1381 gagaagaaga  tgtccactcc  aattgaagtg  ctgtgcaagg  gctatccttc  tgaatttgcc
1441 acatacctga  atttctgccc  ttccttacgt  tttgatgaca  aacctgacta  ctctacctg
1501 agacagctct  tcagaaatct  gttccatcgc  cagggcttct  cctacgacta  tgtgttcgac
1561  ttggaacatg  tcaaatga
  
```

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.