

## Certificate of Analysis

### MRCK $\gamma$ , active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 16-046, 16-046-K, 16-046M

Parent Lot # D17NP005N

The data presented in this document apply to the parent lot shown above and to all pack sizes derived from subsequent vialing runs of this parent lot. An alphabetical suffix after the parent lot number is used to denote each vialing run.

**Product Description:** *N*-terminal FLAG-tagged and *N*-terminal 6His-tagged, recombinant, human MRCK $\gamma$  amino acids 1-473 expressed by baculovirus in Sf21 insect cells. Purified using immunoaffinity chromatography.

Purity 99% by SDS-PAGE and Coomassie blue staining. MW = 56kDa.

**Specific Activity (Parent lot# D17NP005N):** 193U/mg, where one unit of MRCK $\gamma$  activity is defined as 1nmol phosphate incorporated into 50 $\mu$ M KKRNRTLTV per minute at 30°C with a final ATP concentration of 100 $\mu$ M.

**Formulation:** 0.58mg/ml of enzyme in 43mM Tris/HCl pH7.5, 129mM NaCl, 270mM sucrose, 129 $\mu$ g/ml FLAG peptide, 0.9mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution.

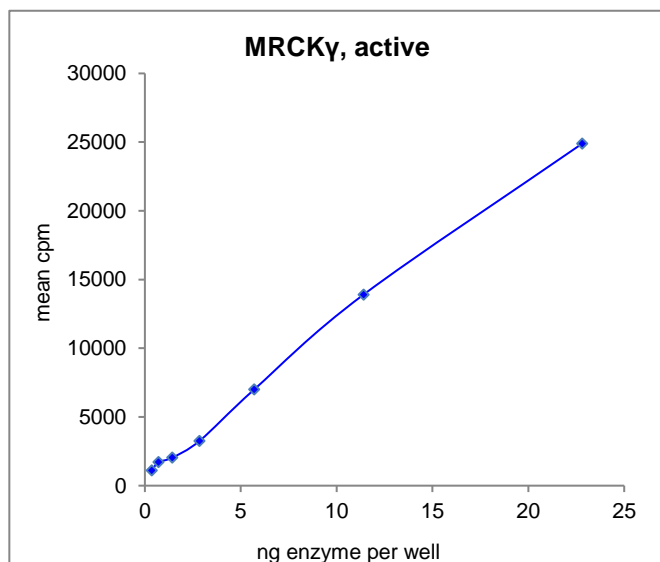
**Storage and Stability:** On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

**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.

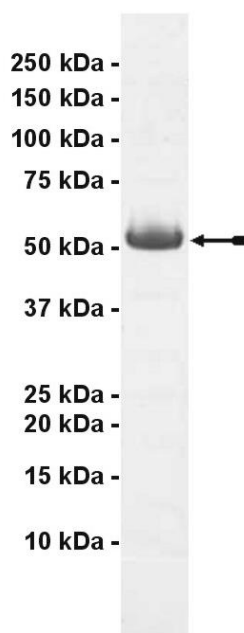
**FOR IN VITRO RESEARCH USE ONLY  
NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**Kinase Assay:** 0.4–22.8ng of this lot of enzyme phosphorylated 50 $\mu$ M KKRNRTLTV 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 MRCK $\gamma$  with the translated 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 MRCK $\gamma$ , active.

## Certificate of Analysis

### Kinase Assay Protocol

#### Stock Solutions:

- 1. 5 x Reaction Buffer:** 40mM MOPS/NaOH pH7.0, 1mM EDTA.
- 2. KKRNRTLTV:** Use at a final assay concentration of 50 $\mu$ M. Prepare a 1mM stock and add 1.25 $\mu$ l of stock per assay point.
- 3. MRCK $\gamma$ , active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 0.4–22.8ng per assay point.
- 4. [ $\gamma$ -<sup>33</sup>P]ATP:** 2.5 x MgAc/[ $\gamma$ -<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 1.25 $\mu$ l of KKRNRTLTV.
3. Add **2.5 $\mu$ l (0.4-22.8ng) MRCK $\gamma$ , active.**
4. Add 6.25 $\mu$ l of dH<sub>2</sub>O.
5. Add 10 $\mu$ l of diluted [ $\gamma$ -<sup>33</sup>P]ATP mixture.
6. Incubate for 30 minutes at 30°C.
7. Stop the reaction by adding 5 $\mu$ l of 3% phosphoric acid.
8. Transfer a 10 $\mu$ 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 dried 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 $\mu$ l of 30% phosphoric acid.

## Certificate of Analysis

### MRCKy, active Sequence Information

<b>Protein</b>	Human MRCKy
<b>Tags</b>	N-terminal FLAG and N-terminal 6His
<b>Native sequence</b>	M20 of the recombinant protein is equivalent to M1 of human MRCKy.
<b>Accession number</b>	GenPept NP_059995.1 (synthetic coding sequence)

#### Recombinant MRCKy amino acid sequence:

```

1  MGGDYKDDDD  KDPHHHHHHM  ERRRLALEQL  ARGEAGGCPG  LDGLLDL LLA  LHHELSSGPL
61  RRERSVAQFL  SWASPFVSKV  KELRLQRDDF  EILKVI GRGA  FGEVTVVRQR  DTGQIFAMKM
121 LHKWEMLKRA  ETACFREERD  VLVKGD SRWV  TTLHYAFQDE  EYLYLVMDYY  AGGDLLTLLS
181 RFEDRLPPEL  AQFYLAEMVL  AIHSLHQ LGY  VHRDVKPDNV  LLDVNGHIRL  ADFGSC LRLN
241 TNGMVDSSVA  VGTPDYISPE  ILQAMEEGKG  HYG PQCDWWS  LGVCAYELLF  GETPFYAESL
301 VETYGKIMNH  EDHLQFPPDV  PDVPASAQDL  IRQLLCRQEE  RLGRGGLDDF  RNHPFFEGVD
361 WERLASSTAP  YIPELRGPM  TSNFDVDDDT  LNHPGTLPPP  SHGAFSGHHL  PFVGFYTSG
421 SHSPESSSEA  WAALERKLQC  LEQEKVELSR  KHQEALHAPT  DHRELEQLRK  EVQTLRDRLP
481 EMLRDKASLS  QT

```

#### Recombinant MRCKy nucleotide sequence:

```

1  atgggaggtg  actacaaaga  cgatgacgac  aaggatccac  atcaccatca  ccatcacatg
61  gaaagaaggt  tgagggctct  ggaacaattg  gcaaggggag  aagccgggtg  ctgccagggc
121 ctggatggac  tcctggacct  gctgctggca  ctgcaccacg  aactgtcctc  tggacccttg
181 aggagggaga  ggtccgttgc  tcagttcctg  tcctgggcat  caccctttgt  ctctaagggt
241 aaggaactgc  gcctccagcg  cgatgatttc  gagatcctga  aagtgattgg  taggggtgct
301 tttggtgaag  ttactgtcgt  gcgccagcgc  gataccggcc  aaatcttcgc  catgaagatg
361 ctccacaaat  gggaaatgct  caagagggcc  gaaaccgcct  gcttccgcga  agagcgcgat
421 gttctcgtaa  aaggcgattc  aagggtgggt  acaaccttgc  attacgcctt  tcaggacgag
481 gagtacctgt  atctggtcat  ggactattac  gccggagggt  acctcctgac  tttgctcagc
541 cgcttcgaag  acaggctgcc  tcccagagtc  gctcaatctt  acctggctga  aatggttctg
601 gccattcata  gcctccacca  gctgggatat  gtccatcgcg  atgtgaaacc  agacaatgtc
661 ctggttgacg  tgaacggcca  cattagactc  gccgatttcg  gtagctgcct  gcgcttgaat
721 acaaacggaa  tggttgactc  cagcgttgc  gtgggtacct  ctgactacat  ctctcccag
781 attctgcaag  caatggagga  aggaaaagga  cactacggcc  cccaatgcga  ctggtggtct
841 ttgggagtg  gcgcctacga  gctgctcttt  ggcgaaacc  ctttctacgc  agaaagcctg
901 gtcgagacat  atggaaaaat  catgaaccac  gaggaccatc  tgcaatttcc  cctgacgtg
961 ccagacgtgc  cagctagcgc  tcaggatctg  attcgccagt  tgttgctgag  acaagaggaa
1021 cgcttgggcc  gcggcgccct  ggacgatttc  agaatcacc  ctttcttga  gggagtcgat
1081 tgggaaaggc  tcgcttcaag  caccgcccc  tacatcccag  agctgagggg  acccatggac
1141 acctccaatt  tcgatgtgga  tgacgacact  ctgaaccacc  caggcacctt  gccaccccc
1201 agccacggcg  cttctccgg  ccaccatctg  ccattcgttg  gctttactta  cacctccgga
1261 agccacagcc  cagaatcttc  ctctgaggct  tgggctgcct  tggagaggaa  attgcagtgc
1321 ctggaacaag  agaaggtcga  actgagccgc  aagcatcagg  aagctttgca  cgctcccact
1381 gaccataggg  aactggaaca  actgaggaag  gaagtgcaa  ctttgcgcga  tagactgcct
1441 gagatgttga  gggataaagc  ctcatgagc  caaacctaa

```

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.