

Certificate of Analysis

MOK, active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-960, 14-960-K, 14-960M

Parent Lot # D15DP013N

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 and C-terminal 6His tagged, recombinant, human MOK amino acids 1-343 expressed by baculovirus in Sf21 insect cells. Purified using immobilized metal affinity chromatography followed by glutathione agarose. Purity 88% by SDS-PAGE and Coomassie blue staining. MW = 70kDa.

Specific Activity (Parent lot# D15DP013N): 410U/mg, where one unit of MOK activity is defined as 1nmol phosphate incorporated into 300µM RSRSRSRSRSRSR per minute at 30°C with a final ATP concentration of 100µM.

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.

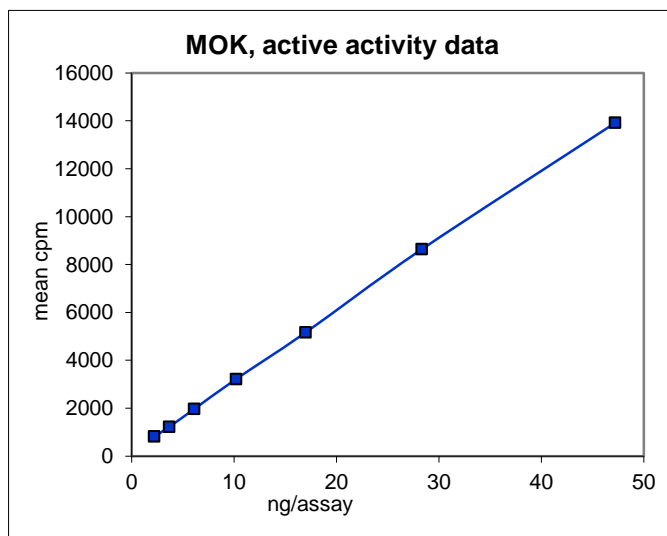
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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..

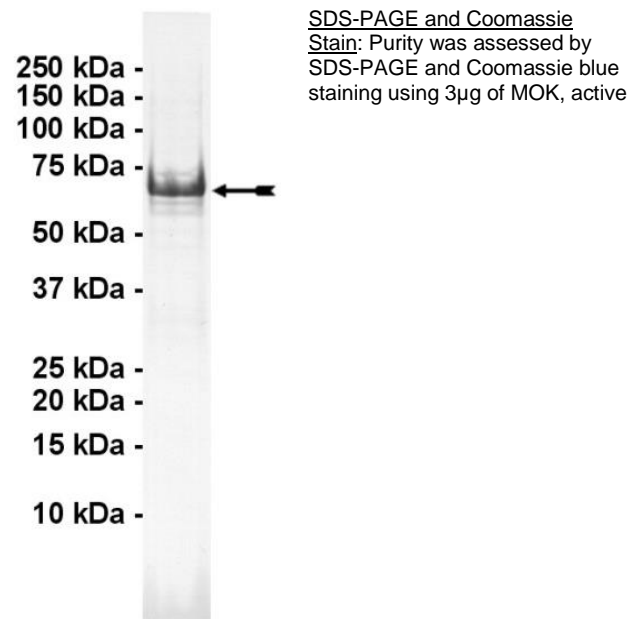
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 2–47ng of this lot of enzyme phosphorylated 300µM RSRSRSRSRSRSR 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 MOK 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. RSRSRSRSRSRSR:** Use at a final assay concentration of 300 μ M. Prepare a 3mM stock and add 2.5 μ l of stock per assay point.
- 3. MOK, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 2–47ng 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 RSRSRSRSRSRSR.
3. Add **2.5 μ l (2–47ng) MOK, 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

MOK, active Sequence Information

Protein	human MOK
Tags	N-terminal GST and C-terminal 6His
Native sequence	M230 of the recombinant protein is equivalent to M1 of human MOK
Accession number	GenBank NM_014226.2

Recombinant MOK amino acid sequence:

```

1  MSPILGYWKI  KGLVQPTRL  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  LVPRGSKEFM  KNYKAIGKIG
241  EGTFFSEVMKM  QSLRDGNYYA  CKQMKQRFES  IEQVNNLREI  QALRRLNPHP  NILMLHEVVF
301  DRKSGSLALI  CELMDMNIYE  LIRGRRYPLS  EKKIMHYMYQ  LCKSLDHIHR  NGIFHRDVKP
361  ENILIKQDVL  KLGDFGSCRS  VYSKQPYTEY  ISTRWYRAPE  CLLTDGFYTY  KMDLWSAGCV
421  FYEIASLQPL  FPGVNELDQI  SKIHDVIGTP  AQKILT KFKQ  SRAMNFDFFP  KKGSGIPLLT
481  TNLSFQCLSL  LHAMVAYDPD  ERIAAHQALQ  HPYFQEQRKT  EKRALGSHRK  AGFPEHPVAP
541  EPLSNSCQIS  KEGRKQKQSL  KQEEDRPKRR  GPENLYFQGV  EACQLGTDDY  DIPTTHHHHH
601  H

```

Recombinant MOK nucleotide sequence:

```

1  atgtccccta  tactaggtta  ttggaaaatt  aagggccttg  tgcaaccac  tcgacttctt
61  ttggaatata  ttgaagaaaa  atatgaagag  catttगतग  agcgcgatga  aggtgataaa
121  tggcgaaaaca  aaaagtttga  attgggtttg  gagtttccca  atcttcctta  ttatattgat
181  ggtगतगtta  aattaacaca  gtctatggcc  atcatacgtt  atatagctga  caagcacaac
241  atgttgggtg  gttgtccaaa  agagcगतगca  gagatttcaa  tgcttgaagg  agcggttttg
301  gatattगतग  acggtगतगt  gagaattgca  tatagtaaag  actttgaaac  tctcaaagtt
361  gattttctta  gcaagctacc  tgaaatगतग  aaaatगतग  aagatगतग  atgtcataaa
421  acatatttaa  atggtगतगca  tgtaaccat  cctगतगtca  tgttगतगta  cgctctगतग
481  gttgttttat  acatggacc  aatगतगcctg  gatगतगtcc  caaaattगतग  ttgttttaaa
541  aaacगतगt  aagctatccc  acaaattगतग  aagtactगतग  aatccगतगca  gatatगतगca
601  tggccttttg  agggcगतगca  agccacगतगt  ggtगतगggcg  accatcctcc  aaaatcगतग
661  ctgggtcccgc  gtggatccaa  ggaattगतग  aagaactata  aagcaattगतग  caaaatगतगga
721  gagggaaacgt  tttctgaagt  tatgaगतगt  caaagcctga  gagatगतगaaa  ctactगतगca
781  tgtaaaacaaa  tgaagcगतगcg  ctttgaaagt  attगतगcaag  tcaacaacct  acgगतगatc
841  caagcactga  ggcगतगcctgaa  tccगतगcacca  aacattctta  tgttगतगcatga  agtगतगtttt
901  gacagaaaat  ctggttctct  tgcactaata  tgtgaaactta  tggगतगatgaa  tatttatगतगg
961  ctaatacगतग  ggagaगतगata  cccattगतगca  gaaaaaaaa  ttatगतगcacta  tatगतगtaccag
1021  ttatगतगtaagt  ccctगतगgatca  tattcगतगcaga  aatगतगgaatat  ttcगतगcagaga  tgtaaaacca
1081  gaaaatatac  taataaगतगca  ggatगतगtccctg  aaattगतगggg  actttगतगgctc  ctgcccगतगagt
1141  gtctattcca  agcगतगccgta  cacगतगgaatac  atctccacct  gctगतगgtaccg  ggccccगतगgag
1201  tgtctcctca  ctगतगatgggtt  ctacगतगcgtac  aगतगatggacc  tगतगtgगतगgcg  cggctगतगgtg
1261  ttctacगतगga  tcgccगतगagtct  gcगतगccccctc  tttcctगतगgag  taaatगतगact  ggaccaaaatc
1321  tcaaaaatcc  acगतगatगतगcat  cggcगतगcacacc  gctcगतगagaaga  tcctcगतगaccaa  gttcaaacगतग
1381  tcgगतगagctta  tgaattttg  ttttctttt  aaaaगतगggat  caggगतगatacc  tctactaaca
1441  accaatttgt  ccccacगतगatg  cctctccctc  ctगतगcacgcaa  tggtगतगgccta  tgatcccगतग
1501  gagagaatcg  ccgcccगतगca  ggccctगतगcg  caccctगतगtact  tccaगतगagaaca  gaggaaaaca
1561  gagaagcगतगg  ctctगतगggcag  ccacगतगaaaa  gctगतगgctttc  cगतगगतगcacc  tगतगtगतगcaccg
1621  gaaccactca  gtaacगतगctg  ccगतगatttcc  aगतगगतगggca  gaaगतगcagaa  acagtccccta
1681  aagcaगतगag  agगतगaccगतगtcc  caगतगगतगcga  ggaccगतगgaaa  acctगतगtatt  tcaggगतगcggtg

```

Certificate of Analysis

1741 gaagcctgtc aattgggcac ggacgattac gacatcccga ctaccacca tcaccatcac
1801 cattaa

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