

## Certificate of Analysis

### MAPKAP Kinase 2, unactive (Recombinant enzyme expressed in *E.coli* cells)

Item # 14-349, 14-349-K, 14-349M

Parent Lot # 1691879

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 MAPKAP Kinase 2, amino acids 46–end, expressed in *E.coli* cells. Purified using glutathione agarose. Purity 78.7% by SDS-PAGE and Coomassie blue staining. MW=70.2kDa.

**Formulation:** 0.914mg/ml of enzyme in 50mM Tris/HCl pH7.5, 150mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 50% glycerol, 1mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol, 5mM DTT. Liquid at -20°C.

**Specific Activity (Parent lot# 1691879):** As provided, this lot demonstrated <1% of maximum activity. Activated by phosphorylation with MAP Kinase 2 (cat# 14-173.)

**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.

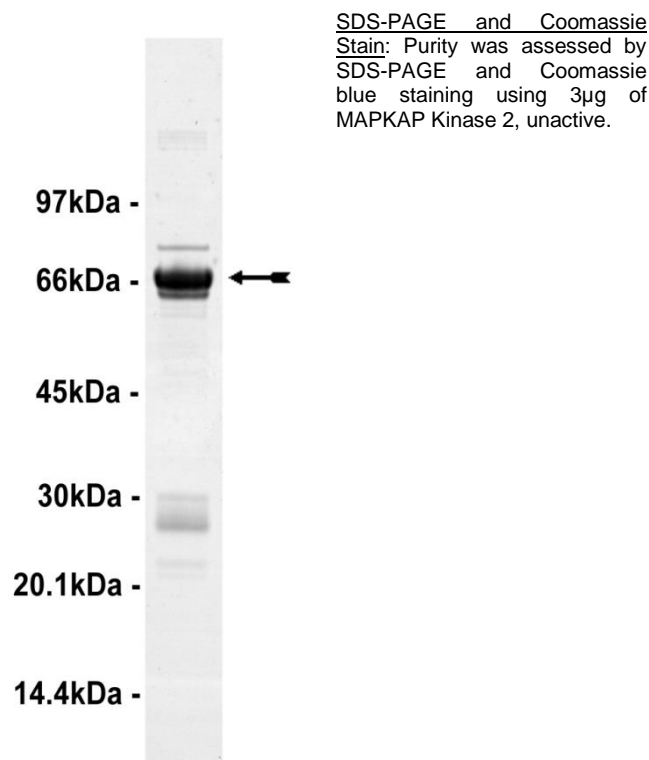
**FOR IN VITRO RESEARCH USE ONLY  
NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**Activation Assay:** 4µM unactive MAPKAP Kinase 2 was activated using 0.2µM active MAP Kinase 2 (cat# 14-173) diluted 500–1000 fold, and the increased activity against (KKLNRTLVA, cat# 12-240) determined. The activation and assay are described on page two. Results of this assay are shown below.

**MS Tryptic Fingerprint:** Confirmed identity as MAPKAP Kinase 2 with the translated sequence listed on page three.

Active MAP Kinase 2	Unactive MAPKAP Kinase 2	Mean cpm	Comments
0.34µg	7µg	7480	Kinase activity
none	7µg	640	Background



## Certificate of Analysis

### Kinase Assay Protocol

#### Stock Solutions:

- 10 x Activation Buffer:** 50mM Tris/HCl pH7.5, 1mM EGTA, 1% 2-mercaptoethanol.
- 20 x Reaction Buffer:** 1M Na- $\beta$ -glycerophosphate, 2mM EGTA.
- Enzyme Dilution Buffer:** 20mM MOPS/NaOH pH7.0, 1mM EDTA, 5% glycerol, 0.01% Brij-35, 1% 2-mercaptoethanol, 1mg/ml BSA.
- MAPKAP Kinase 2, inactive:** Use at a final assay concentration of 4 $\mu$ M (0.28mg/ml). Prepare a 1.4mg/ml stock and add 5 $\mu$ l of stock per assay point.
- MAP Kinase 2, active (Catalogue # 14-173):** Use at a final assay concentration of 0.2 $\mu$ M (0.014mg/ml). Prepare a 0.14mg/ml stock and add 2.5 $\mu$ l of stock per assay point.
- Stage One 5 x Mg/ATP:** 50mM magnesium acetate, 0.5mM ATP.
- [ $\gamma$ -<sup>33</sup>P]ATP:** 2.5 x magnesium acetate/[ $\gamma$ -<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.)
- MAPKAP Kinase 2 substrate peptide (KKLNRTLVA) (Catalogue# 12-240):** Use at a final assay concentration of 30 $\mu$ M. Prepare a 300 $\mu$ M stock. Add 2.5 $\mu$ l of stock per assay point.

#### Assay Protocol:

##### **Stage One:** *Activation of MAPKAP Kinase 2, by MAP Kinase 2:*

- Add 2.5 $\mu$ l of 10 x activation buffer to a microcentrifuge tube.
- Add 5 $\mu$ l (7 $\mu$ g) of **MAPKAP Kinase 2, inactive**.
- Add 2.5 $\mu$ l (0.34 $\mu$ g) of **MAP Kinase 2, active**.
- Add 10 $\mu$ l of dH<sub>2</sub>O.
- Add 5 $\mu$ l of stage one Mg/ATP mixture.
- Incubate for 30 minutes at 30°C.
- Stop reaction by diluting the tubes 500–1000 fold and incubating on ice.

##### **Stage Two:** *Phosphorylation of KKLNRTLVA substrate peptide with MAPKAP Kinase2 (96 well plate format):*

- Add 1.25 $\mu$ l of reaction buffer per assay to wells.
- Add 2.5 $\mu$ l of substrate peptide (**KKLNRTLVA**).
- Add 2.5 $\mu$ l of diluted MAPKAP Kinase 2 from **Stage One**.
- Add 8.7 $\mu$ l of dH<sub>2</sub>O.
- Add 10 $\mu$ l of the diluted [ $\gamma$ -<sup>33</sup>P] ATP.
- Incubate for 10 minutes at 30°C.
- Stop the reaction by adding 5 $\mu$ l of 3% phosphoric acid.
- Slowly transfer 10 $\mu$ l onto appropriate area of a **P30 Filtermat**.
- Wash filtermat three times for 5 minutes with 75mM phosphoric acid.
- Wash filtermat once for 2 minutes with methanol.
- Transfer filtermat to a sealable plastic bag and add 4ml scintillation cocktail.
- Read in scintillation counter. Compare cpm of enzyme samples with cpm of control samples that contain all appropriate assay components plus 1 $\mu$ l of 30% phosphoric acid enzyme (background control).

## Certificate of Analysis

## Certificate of Analysis

### MAPKAP Kinase 2, Sequence Information

<b><u>Protein</u></b>	human MAPKAP Kinase 2
<b><u>Tags</u></b>	N-Terminal GST
<b><u>Native sequence</u></b>	F243 of the recombinant protein is equivalent to F46 of human MAPKAP Kinase 2
<b><u>Accession number</u></b>	GenBank NM_032960. The recombinant sequence is D116H (native co-ordinates) with respect to GenBank NM_032960. This conflict is reported in SWISSPROT P49137. The cDNA also contains the translational conflict A399G, and a c-myc epitope at the C-terminus.

#### **Recombinant MAPKAP Kinase 2 amino acid sequence:**

```

1  MSPILGYWKI  KGLVQPTRLL  LEYLEEKYEE  HLYERDEGDK  WRNKKFELGL  EFPNLPYYID
61  GDVKLTQ SMA  IIRYIADKHN  MLGGCPKERA  EISMLEGAVL  DIRYGVSRIA  YSKDFETLKV
121  DFLSKLP EML  KMFEDRLCHK  TYLNGDHSVH  PDFMLYDALD  VVLYMDPMCL  DAFPKLVCFK
181  KRIEAI P QID  KYLKSSKYIA  WPLQGWAQTF  GGGDHPPKSD  LVPRGSPGIS  GGGGGILEAT
241  MEFHVK SGLQ  IKKNAIIDDY  KVTSQVLGLG  INGKVLQIFN  KRTQEKFALK  MLQDCPKARR
301  EVELHW RASQ  CPHIVRIVDV  YENLYAGRKC  LLIVMECLDG  GELFSRIQDR  GDQAFTEREA
361  SEIMKS IGEA  IQYLHSINIA  HRDVKPENLL  YTSKRPNAIL  KLTDFGFAKE  TTSHNSLTTP
421  CYTPYV VAPE  VLGPEKYDKS  CDMWSLGVIM  YILLCGYPPF  YSNHGLAISP  GMKTRIRMGQ
481  YEFNP E WSE  VSEEVKMLIR  NLLKTEPTQR  MTITEFMNHP  WIMQSTKVPQ  TPLHTSRVLK
541  EDKERW EDVK  EEMTSALATM  RVDYEQIKIK  KIEDASNPLL  LKRRKKARAL  EAAALGHMEQ
601  KLISEEDLK

```

#### **Recombinant MAPKAP Kinase 2 nucleotide sequence:**

```

1  atgtccccta  tactaggtta  ttggaaaatt  aagggccttg  tgcaaccac  tcgacttctt
61  ttggaatata  ttgaagaaaa  atatgaagag  catttgtatg  agcgcgatga  aggtgataaaa
121  tggcgaaaaca  aaaagtttga  attgggtttg  gagtttccca  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  tgaaatgctg  aaaatgttcg  aagatcgttt  atgtcataaaa
421  acataattta  atggtgatca  tgtaacccat  cctgacttca  tgttgtatga  cgctcttgat
481  gttgttttat  acatggacc  aatgtgctg  gatgcttcc  caaaattagt  ttgttttaaa
541  aaacgtattg  aagctatccc  acaaattgat  aagtacttga  aatccagcaa  gtatatagca
601  tggcctttg  agggctggca  agccacgttt  ggtggtggcg  accatctcc  aaaatcggat
661  ctggttccgc  gtggatcccc  gggaatttcc  ggtggtggtg  gtggaattct  agaggccacc
721  atggagtcc  acgtcaagtc  cggcctgcag  atcaagaaga  acgcatcat  cgatgactac
781  aaggtcacca  gccaggtcct  gggctgggg  atcaacggca  aagttttgca  gatcttcaac
841  aagaggacc  aggagaaatt  cgccctcaa  atgcttcagg  actgccccaa  ggcccgcagg
901  gaggtggagc  tgcactggcg  ggctcccag  tgcccgcaca  tcgtacggat  cgtggatgtg
961  tacgagaatc  tgtacgcagg  gaggaagtgc  ctgctgattg  tcatggaatg  tttggacggt
1021  ggagaactct  ttagccgaat  ccaggatcga  ggagaccagg  cattcacaga  aagagaagca
1081  tccgaaatca  tgaagagcat  cggtgaggcc  atccagtatc  tgcattcaat  caacattgcc
1141  catcgggatg  tcaagcctga  gaatctctta  tacacctcca  aaaggcccaa  cgccatcctg
1201  aaactcactg  actttggctt  tgccaaggaa  accaccagcc  acaactcttt  gaccactcct
1261  tgttatacac  cgtactatgt  ggctccagaa  gtgctgggtc  cagagaagta  tgacaagtcc
1321  tgtgacatgt  ggtccctggg  tgtcatcatg  tacatcctgc  tgtgtgggta  tcccccttc
1381  tactccaacc  acggccttgc  catctctccg  ggcatgaaga  ctgcacccg  aatgggccag
1441  tatgaatttc  ccaaccaga  atggtcagaa  gtatcagagg  aagtgaagat  gctcattcgg
1501  aatctgctga  aaacagagcc  caccagaga  atgaccatca  ccgagtttat  gaaccaccct
1561  tggatcatgc  aatcaacaaa  ggtccctcaa  accccactgc  acaccagccg  ggtcctgaag
1621  gaggacaagg  agcgggtggga  ggatgtcaag  gaggagatga  ccagtgcctt  ggccacaatg

```

## Certificate of Analysis

```
1681 cgcgttgact acgagcagat caagataaaa aagattgaag atgcatccaa ccctctgctg
1741 ctgaagaggc ggaagaaagc tcgggccctg gaggctgcgg ctctgggcca catggagcag
1801 aagctgatca gcgaggagga cctgaagtga
```

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.