

Certificate of Analysis

MAPKAP Kinase 3, unactive

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

Item # 14-586, 14-586-K, 14-586M

Parent Lot # 1952459

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 3, amino acids 2-end, expressed in *E.coli* cells. Purified using glutathione agarose. Purity 81.1% by SDS-PAGE and Coomassie blue staining. MW=69.8kDa.

Specific Activity (Parent lot# 1952459): As provided, this lot demonstrated <1% of maximum activity. Activated by phosphorylation with SAPK2a (cat# 14-587).

Formulation: 3.177mg/ml of enzyme in 50mM Tris/HCl pH7.5, 300mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 1mM benzamide, 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**

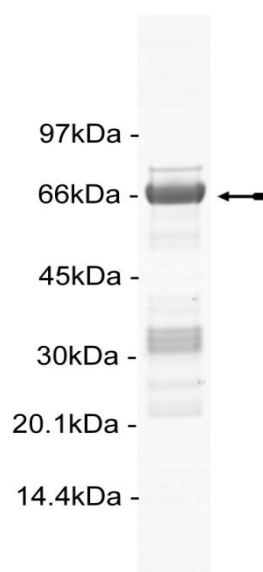
Quality Control Testing

Activation Assay: 4µM unactive MAPKAP Kinase 3 was activated using 0.2µM SAPK2a (cat# 14-587) diluted 200–2500 fold, and the increased activity against (KKLNRTLSVA) (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 3 with the translated sequence listed on page three

SDS-PAGE and Coomassie Stain: Purity was assessed by SDS-PAGE and Coomassie blue staining using 3µg of MAPKAP Kinase 3, unactive.

Active SAPK2a	Unactive MAPKAP Kinase 3	Mean cpm	Comments
0.218µg	None	74	Background
none	3.5µg	115	Background
0.218µg	3.5µg	4834	Kinase activity



Certificate of Analysis

Kinase Activation Assay Protocol

Stock Solutions:

1. **10 x Activation Buffer:** 50mM Tris/HCl pH7.5, 1mM EGTA, 1% 2-mercaptoethanol.
2. **20 x Assay Buffer:** 1M Na- β -glycerophosphate, 2mM EGTA, pH7.5.
3. **Enzyme Dilution Buffer:** 20mM MOPS/NaOH pH7.0, 1mM EDTA, 5% glycerol, 0.01% Brij-35, 1% 2-mercaptoethanol, 1mg/ml BSA.
4. **MAPKAP Kinase 3, inactive:** Use at a final assay concentration of 4 μ M (0.28mg/ml). Prepare a 1.4mg/ml stock and add 5 μ l of stock per assay point.
5. **SAPK2a, active (Catalogue# 14-587):** Use at a final assay concentration of 0.2 μ M (0.009mg/ml). Prepare a 0.09mg/ml stock and add 2.5 μ l of stock per assay point.
6. **Stage One 5 x Mg/ATP:** 50mM magnesium acetate, 0.5mM ATP.
7. **(KKLNRTLVA):** Use at a final concentration of 30 μ M. Make up a 300 μ M stock. Use 2.5 μ l of stock per assay point.
8. **[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 268cpm/pmol as required.)

Assay Procedure:

Stage One: *Activation of MAPKAP Kinase 3 by SAPK2a*

1. Add 2.5 μ l of 10 x activation buffer to a microcentrifuge tube.
2. Add **5 μ l of MAPKAP Kinase 3, inactive.**
3. Add 2.5 μ l of **SAPK2a, active.**
4. Add 10 μ l of dH₂O.
5. Add 5 μ l of stage one Mg/ATP mixture.
6. Incubate for 15 minutes at 30°C.
7. Stop reaction by diluting the tubes 200–2500 fold and incubating on ice.

Stage Two: *Phosphorylation of KKLNRTLVA by MAPKAP Kinase 3*

1. Add 1.25 μ l of 20 x reaction buffer to a microcentrifuge tube.
2. Add 2.5 μ l of **(KKLNRTLVA).**
3. Add **2.5 μ l (0.2–1.4ng) MAPKAP Kinase3, active** from **Stage One.**
4. Add 8.75 μ l dH₂O
5. Add 10 μ l of [γ -³³P]ATP mix.
6. Incubate for 10 minutes at 30°C.
7. Slowly transfer 20 μ l onto the centre of a 2 cm x 2 cm **P81** paper.
8. Wash assay squares twice for 5 minutes with 75mM phosphoric acid.
9. Wash assay squares once for 2 minutes with acetone.
10. Transfer assay squares to scintillation vials and add 1ml scintillation cocktail.
11. Read in scintillation counter. Compare cpm of enzyme samples with cpm of control samples that contain all assay components and 1 μ l of 30% phosphoric acid (background control).

Certificate of Analysis

MAPKAP Kinase 3 Sequence Information

<u>Protein</u>	Human MAPKAP Kinase 3
<u>Tags</u>	N-terminal GST
<u>Native sequence</u>	D232 of the recombinant protein is equivalent to D2 of human MAPKAP Kinase 3
<u>Accession number</u>	GenBank NM_004635

Recombinant MAPKAP Kinase 3 amino acid sequence:

```

1  MSPILGYWKI  KGLVQPTRL  LEYLEEKYE  HLYERDEGD  WRNKKFELG  EFPNLPYYI
61  GDVKLTQSM  IIRYIADKH  MLGGCPKER  EISMLEGAV  DIRYGVSRI  YSKDFETLK
121  DFLSKLP  KMFEDRLCH  TYLNGDHV  PDFMLYDAL  VVLYMDPM  DAFPKLVC
181  KRIEAI  KYLKSSKY  WPLQGWA  GGDHPPK  LEVLFQGP  SDGETAEE
241  GPVPPP  GPGLGGAP  RREPKKY  DDYQLSK  GLGVNGKV  CFHRRTGQ
301  ALKLLY  ARQEVDDH  ASGGPHIV  LDVYENMH  KRCLLIIM  MEGGELFS
361  QERGDQ  REAAEIMR  GTAIQFL  NIAHRDV  NLLYTSKE  AVLKLTDF
421  AKETTQ  TPCYTPY  PEVLGPE  KSCDMWS  IMYILLCG  PFYSNTGQ
481  SPMKRR  GQYGFNP  SEVSEDA  IRLLLKT  ERLTITQ  HPWINQSM
541  PQTPLH  LQEDKDH  VKEEMTS  TMRVDYD  IKDLKTS  LLNKRRKQ
601  GSSSAS  NQ

```

Recombinant MAPKAP Kinase 3 nucleotide sequence:

```

1  atgtccccta  tactaggtta  ttggaaaatt  aagggccttg  tgcaaccac  tcgacttctt
61  ttggaatata  ttgaagaaaa  atatgaagag  catttgatg  agcgcgatg  aggtgataaa
121  tggcgaaa  aaaagtttga  attgggttg  gagtttcca  atcttcctta  ttatattgat
181  ggtgatgta  aattaacaca  gtctatggc  atcatacg  atatagctga  caagcacaac
241  atgttgggtg  gttgtccaaa  agagcgtgca  gagatttcaa  tgcttgaagg  agcggttttg
301  gatattagat  acggtgtttc  gagaattgca  tataagtaa  actttgaaac  tctcaaagtt
361  gattttctta  gcaagctacc  tgaaatgctg  aaaatgttcg  aagatcgtt  atgtcataaa
421  acatatttaa  atggtgatca  tgtaaccat  cctgacttca  tgttgatga  cgctcttgat
481  gttgttttat  acatggacc  aatgtgctg  gatgcttcc  caaaattagt  ttgttttaa
541  aaacgtattg  aagctatccc  acaaattgat  aagtacttga  aatccagcaa  gtatatagca
601  tggcctttgc  agggctggca  agccactgtt  ggtggtggcg  accatcctcc  aaaatcggat
661  ctggaagtct  tgttccagg  gccctggga  tccgatggtg  aaacagcaga  ggagcagggg
721  ggcctgtg  ccccgcagt  tgcaccggc  ggaccggct  tggcggtgc  tccggggggg
781  cggcgggagc  ccaagaagta  cgcagtgacc  gacgactacc  agttgtccaa  gcaggtgctg
841  ggcctgggtg  tgaacggcaa  agtctggag  tgcttccatc  ggcgactgg  acagaagtgt
901  gccctgaagc  tctgtatga  cagcccaag  gcccggcagg  aggtagacca  tctctggcag
961  gcttctggcg  gccccatat  tgtctgcatc  ctggatgtgt  atgagaacat  gcaccatggc
1021  aagcgtgtc  tcctcatcat  catggaatgc  atggaaggtg  gtgagttgtt  cagcaggatt
1081  caggagcgtg  gcgaccaggc  tttcactgag  agagaagctg  cagagataat  gcgggatatt
1141  ggcactgcc  tccagtttct  gcacagccat  aacattgcc  accgagatgt  caagcctgaa
1201  aacactactc  acacatctaa  ggagaaagac  gcagtgtta  agctcaccga  ttttggcttt
1261  gtaaggaga  ccacccaaa  tgccctgcag  acaccctgct  atactcccta  ttatgtggcc
1321  cctgagggtc  tgggtccaga  gaagtatgac  aagtcagtgt  acatgtggtc  cctgggtgtc
1381  atcatgtaca  tctctctttg  tggcttccca  cccttctact  ccaacacggg  ccaggccatc
1441  tccccgggga  tgaagaggag  gattegcctg  ggccagtacg  gcttcccaa  tctctgagtgg
1501  tcagaagtct  ctgaggatgc  caagcagctg  atccgcctcc  tgttgaagac  agacccaca
1561  gagaggctga  ccatcactca  gttcatgaac  caccctgga  tcaaccaatc  gatggtagtg
1621  ccacagacc  cactccacac  ggcccagtg  ctgcaggagg  acaagacca  ctgggacgaa
1681  gtcaaggagg  agatgaccag  tgccttgcc  actatgcggg  tagactacga  ccaggtgaa

```

Certificate of Analysis

1741 atcaaggacc tgaagacctc taacaaccgg ctctcaaca agaggagaaa aaagcaggca
1801 ggcagctcct ctgcctcaca gggctgcaac aaccagtag

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.