

Certificate of Analysis

MAP Kinase 2/Erk2, unactive (Recombinant enzyme expressed in *E.coli* cells.)

Item # 14-536, 14-536-K, 14-536M

Parent Lot # 1606930

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, full length MAP Kinase 2, expressed in *E.coli* cells. Purified using glutathione agarose. Purity 97.6% by SDS-PAGE and Coomassie blue staining. MW = 67.8kDa.

Formulation: 1.463mg/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. Liquid at -20°C.

Specific Activity (Parent lot# 1606930): As provided, this lot demonstrated <1% of maximum activity. Activated by phosphorylation with MEK1(h) (cat# 14-429.)

Storage and Stability: On receipt of material store at -20°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.

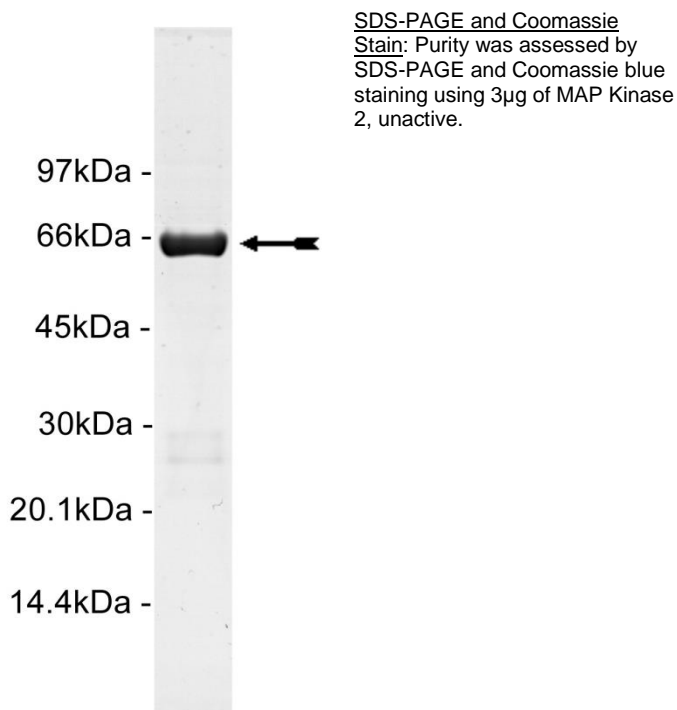
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 2µM unactive MAP Kinase 2 was activated using 0.4µM MEK1(h) (cat# 14-429) diluted 250-fold, and the increased activity against myelin basic protein determined. The activation and assay are described on page two. Results of this assay are shown below.

MS Tryptic Fingerprint: Confirmed product identity as MAP Kinase 2, unactive with the translated sequence listed on pages three and four.

Active MEK1(h)	Unactive MAP Kinase 2	Mean cpm	Comments
0.7µg	3.4µg	17243	Kinase Activity
None	3.4µg	44	Background



Certificate of Analysis

Kinase Assay Procedure

Stock Solutions:

1. **10 x Activation Buffer:** 500mM Tris/HCl pH7.5, 1mM EGTA, 1mM Na₃VO₄, 1% 2-mercaptoethanol, 0.3% Brij-35.
2. **5 x Reaction buffer:** 125mM Tris/HCl pH7.5, 0.1mM EGTA.
3. **Dilution Buffer:** 50mM Tris/HCl pH7.5, 0.1mM EGTA, 0.1% 2-mercaptoethanol, 0.1mM Na₃VO₄, 1mg/ml BSA.
4. **MAP Kinase 2, unactive:** Use at a final assay concentration of 2µM (0.136 mg/ml). Dilute in Dilution Buffer to 1.36mg/ml. Use 2.5µl per assay point.
5. **MEK1, active (Catalogue# 14-429):** Use at a final assay concentration of 0.4µM (0.028mg/ml). Dilute with Dilution Buffer to 0.28mg/ml. Use 2.5µl per assay point.
6. **Stage One Mg/ATP:** 50mM MgAc, 0.5mM ATP.
7. **[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.)
8. **Myelin Basic Protein (MBP) substrate:** Use a final assay concentration of 330µg/ml. Make up a 3.3mg/ml stock. Add 2.5µl of stock per assay point.

Assay Protocol:

Stage One: *Activation of MAP kinase 2 by MEK1*

1. Add 2.5µl of 10 x activation buffer to a microcentrifuge tube.
2. Add 12.5µl distilled H₂O.
3. Add 2.5µl (3.4µg) **MAP kinase 2, unactive.**
4. Add 2.5µl (0.7µg) **MEK1, active.**
5. Add 5µl of stage one 5 x Mg/ATP.
6. In appropriate controls, add dilution buffer to a final volume of 25µl.
7. Incubate at 30°C for 30 minutes.
8. Stop the reaction by diluting 100-fold in dilution buffer and storing on ice.

Stage Two: *Phosphorylation of MBP by MAP kinase 2 (96 well plate format)*

1. Add 5µl of 5 x reaction buffer per well.
2. Add 2.5µl of MBP substrate.
3. Add 2.5µl of diluted **MAP kinase 2 (3.4ng) from Stage One.**
4. Make up to 15µl with dH₂O.
5. Add 10µl of the diluted [γ-³³P]ATP mixture.
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5µl of 3% phosphoric acid.
8. Spot 10µl 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 scintillation cocktail.
12. Read in scintillation counter. Compare cpm of enzyme samples with cpm of control samples containing all components and 1µl of 30% phosphoric acid.

Certificate of Analysis

MAP Kinase 2 Sequence Information

<u>Protein</u>	human MAP Kinase 2
<u>Tags</u>	N-terminal GST
<u>Native sequence</u>	M227 of the recombinant protein is equivalent to M1 of human MAP Kinase 2
<u>Accession number</u>	GenBank NM_002745. The cDNA sequence corresponds to a humanised version of the mouse MAP Kinase 2 coding sequence described in GenBank D10939. At the protein level the sequence is identical to GenBank NM_002745.

Recombinant MAP Kinase 2 amino acid sequence:

```

1  MSPIILGYWKI  KGLVQPTRL  LEYLEEKYEE  HLYERDEGDK  WRNKKFELGL  EFPNLPYYID
61  GDVKLTQSMA  IIRYIADKHN  MLGGCPKERA  EISMLEGAVL  DIRYGVSRIA  YSKDFETLKV
121  DFLSKLP EML  KMFEDRLCHK  TYLNGDHVTH  PDFMLYDALD  VVLYMDPMCL  DAFPKLVCFK
181  KRIEAI PQID  KYLKSSKYIA  WPLQGWQATF  GGGDHPPKSD  LVPRGSMAAA  AAAGAGPEMV
241  RGQVFDV GPR  YTNLSYIGEG  AYGMVCSAYD  NVNKVRVAIK  KISPFEHQTY  CQRTLREIKI
301  LLRFRHENII  GINDIIRAPT  IEQMKDVYIV  QDLMETDLYK  LLKTQHLSND  HICYFLYQIL
361  RGLKYIHSAN  VLHRDLKPSN  LLLNTTCDLK  ICDFGLARVA  DPDHDHTGFL  TEYVATRWR  YR
421  APEIMLNSKG  YTKSIDIWSV  GCILAEMLSN  RPIFPGKHYL  DQLNHILGIL  GSPSQEDLNC
481  IINLKARNYL  LSLPHKNKVP  WNRLFNPADS  KALDLLDKML  TFNPHKRIEV  EQALAHPYLE
541  QYYDPSDEPI  AEAPFKFDME  LDDLPEKELK  ELIFEETARF  QPGYRS

```

Recombinant MAP Kinase 2 nucleotide sequence:

```

1  atgtccccta  tactaggtta  ttggaaaatt  aagggccttg  tgcaaccac  tcgacttctt
61  ttggaatata  ttgaagaaaa  atatgaagag  catttgatg  agcgcgatga  aggtgataaa
121  tggcgaaaca  aaaagtttga  attgggtttg  gagtttccca  atcttcctta  ttatattgat
181  ggtgatgta  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  atgtcataaa
421  acatatttaa  atggtgatca  tgtaaccat  cctgacttca  tgttgatga  cgctcttgat
481  gttgttttat  acatggacc  aatgtgectg  gatgcttcc  caaaattagt  ttgttttaa
541  aaacgtattg  aagctatccc  acaaattgat  aagtacttga  aatccagcaa  gtatatagca
601  tggcctttgc  agggctggca  agccacgttt  ggtggtggcg  accatcctcc  aaaatcggat
661  ctggttccgc  gtggatccat  ggcggcgcga  gcagcggccg  gcgctggacc  ggagatggtc
721  cgcgggcagg  tgtttgactg  agggcgcgcg  tacaccaacc  tctcgtacat  cggagaaggc
781  gcctacggca  tggtttgctc  tgcttatgat  aatgtcaaca  aagttcgagt  tgctatcaag
841  aaaatcagtc  cttttgagca  ccagacctac  tgtcaaagaa  ccctaagaga  gataaaaatc
901  ttactgcgct  tcagacatga  gaacatcatt  ggcatcaatg  acatcatccg  ggcaccaacc
961  attgagcaaa  tgaaagatgt  atatatagta  caggacctca  tggagacgga  cctttacaag
1021  ctcttgaaga  cacagcacct  cagcaatgac  cacatctgct  attttcttta  tcagatcctg
1081  agagggctaa  agtatatcca  ttcagctaac  gttctgcacc  gtgacctcaa  gccttccaac
1141  ctctgctga  acaccacttg  tgatctcaag  atctgtgact  ttggccttgc  ccgtgttgca
1201  gatccagatc  atgatcacac  agggttcttg  acagagtacg  tagccacacg  ttggtacaga
1261  gctccagaaa  ttatggtgaa  ttccaagggt  tataccaagt  ccattgatat  ttggtctgtg
1321  ggctgcatcc  tggcagagat  gctatccaac  aggcctatct  tcccaggaaa  gcattacctt
1381  gaccagctga  atcacatcct  gggatttctt  ggatctccat  cacaggaaga  tctgaattgt
1441  ataataaatt  taaaagctag  aaactatttg  ctttctctcc  cgcacaaaaa  taaggtgcca
1501  tggaacaggt  tgttcccaa  tgctgactcc  aaagctctgg  atttactgga  taaaatggtg
1561  acatttaacc  ctcaaaagag  gattgaagtt  gaacaggctc  tggcccacc  atacctggag
1621  cagtattatg  acccaagtga  tgagcccatt  gctgaagcgc  cattcaagtt  tgacatggag

```

Certificate of Analysis

1681 ttggacgact tacctaagga gaagctcaaa gaactcattt ttgaagagac tgctagattc
1741 cagccaggat acagatctta a

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.