

Certificate of Analysis

MEK1, unactive

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

Item # 14-420, 14-420-K, 14-420M

Parent Lot # 2350988

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, and C-terminal 6His-tagged, recombinant full-length human MEK1 expressed in *E.coli* cells. Purified using glutathione-agarose followed by Ni²⁺/NTA-agarose. Purity 76.8% using SDS-PAGE and Coomassie blue staining. MW = 71kDa.

Formulation: 2.957mg/ml of enzyme in 50mM Tris/HCl pH7.5, 150mM NaCl, 5% glycerol, 1mM benzamidine, 0.2mM PMSF, 0.1mM EGTA, 0.1% 2-mercaptoethanol, 0.03% Brij-35. 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.

Specific Activity (Parent lot# 2350988): Less than 1% of the value obtained with MEK1 that has been activated with c-Raf. MEK1 activity determined by activation of 1µM MAP kinase 2.

Handling Recommendations: Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled microcentrifuge 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**

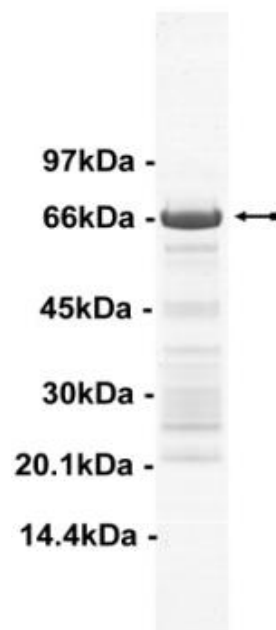
Quality Control Testing

Activation Assay: 5µM unactive MEK1 was activated using 0.1µM cRaf, then diluted 100-fold, used to activate 1µM unactive MAP kinase 2, and the increased activity of the MAP kinase 2 against MBP determined. The activation of MEK1, subsequent activation of MAP kinase 2, and assay are described on page two. Results of this assay are shown below.

MS Tryptic Fingerprint: Confirmed product identity as MEK1 with the translated sequence listed on page four.

Active c-Raf	Unactive MEK1	Mean cpm	Comment
None	8.875µg	85	Inactive MEK1
110ng	8.875µg	13653	Activated MEK1

SDS-PAGE and Coomassie Stain: Representative gel from this lot. Purity was assessed by SDS-PAGE and Coomassie blue staining using 3µg of MEK1.



Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

1. **10 x Activation Buffer:** 500mM Tris/HCl pH7.5, 1mM EGTA, 1mM Na₃VO₄, 1% 2-mercaptoethanol, 0.3% Brij-35.
2. **Dilution Buffer:** 50mM Tris/HCl pH7.5, 0.1mM EGTA, 0.1% 2-mercaptoethanol, 0.1mM Na₃VO₄, 1mg/ml BSA.
3. **Stage One 5 x Mg/ATP:** 50mM magnesium acetate, 1mM ATP.
4. **MEK1, unactive (Catalogue# 14-420):** Add to a final assay concentration of 5µM (0.355mg/ml). The volume is calculated by the formula, 0.355mg/ml divided by the concentration **2.957mg/ml** times 25µl.
5. **c-Raf-1, active (Catalogue# 14-352):** Add to a final assay concentration of 0.1µM (4.4µg/ml). Dilute with Dilution Buffer to 44µg/ml. Use 2.5µl per assay point.
6. **MAP kinase 2, unactive (Catalogue# 14-198):** Add to a final assay concentration of 1µM (67µg/ml). Dilute with Dilution Buffer to 0.67mg/ml. Use 2.5µl per assay point.
7. **Stage Two 5 x Mg/ATP:** 50mM magnesium acetate, 0.5mM ATP.
8. **10 x Assay Buffer:** 250mM Tris/HCl pH7.5, 2mM EGTA.
9. **MBP substrate (Catalogue# 13-104):** Dilute with dH₂O to prepare a 3.33mg/ml stock. Use 2.5µl stock per assay point.
10. **[γ-³³P]ATP:** 2.5 x magnesium acetate/[γ-³³P]ATP cocktail: 25mM magnesium acetate and 0.25mM ATP to which is added [γ-³³P]ATP (specific activity approximately 500cpm/pmol as required.)

Assay Procedure:

Stage One: *Activation of MEK1*

1. Add 2.5µl of 10 x activation buffer to a microcentrifuge tube.
2. Add 12µl dH₂O.
3. Add **3.0µl MEK1, unactive.**
4. Add 2.5µl **c-Raf, active.**
5. Add 5µl stage one 5 x Mg/ATP.
6. In appropriate controls, add dilution buffer to a final volume of 25µl.
7. Incubate for 30 minutes at 30°C.
8. Stop the reaction by diluting the MEK-1 250-fold in dilution buffer. Store on ice.

Stage Two: *Activation of MAPK2*

1. Add 2.5µl 10 x activation buffer to a microcentrifuge tube.
2. Add 12.5µl of dH₂O.
3. Add 2.5µl of **MAP kinase 2, unactive.**
4. Add 2.5µl of diluted **MEK1** from **Stage One.**
5. Add 5µl of stage two 5 x Mg/ATP.
6. Incubate for 15 minutes at 30°C.
7. Immediately transfer 1µl of the mixture to the stage three component mixture.

Certificate of Analysis

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

1. Add 2.5µl 10 x assay buffer per assay to wells.
2. Add 9µl dH₂O.
3. Add 2.5µl **MBP**.
4. Add 1µl of the **Stage Two** reaction product.
5. Add 10µl of the 2.5 x magnesium acetate/ [γ-³³P]ATP cocktail.
6. Incubate for 15 minutes at 30°C.
7. Stop the reaction by adding 5µl of 3% phosphoric acid.
8. Transfer a 10µl 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.
9. Transfer the filtermat to a sealable plastic bag and add 4ml of scintillation cocktail.
10. Read in a scintillation counter. Compare cpm of enzyme samples with cpm of control samples that contain all assay components plus 1µl 30% phosphoric acid and control samples with all assay components except c-Raf.

Certificate of Analysis

MEK1 Sequence Information

<u>Protein</u>	Human MEK1
<u>Tags</u>	N-terminal GST and C-terminal 6His tags
<u>Native sequence</u>	M231 of the recombinant protein is equivalent to M1 of human MEK1
<u>Accession number</u>	EMBL Z30163, possessing amino acid substitution A327G. Please note, introduction of the A327G mutation to the rabbit sequence (Z30163) results in expression of recombinant MEK1 protein that is 100% identical to wild-type human MEK1 described in SwissProt Q02750, note this SwissProt entry fails to record M1 and starts at P2.

Recombinant MEK1 amino acid sequence:

```

1 MSPILGYWKI KGLVQPTRL L LEYLEEKYEE HLYERDEGDK WRNKKFELGL EFPNLPYYID
61 GDVKLTQ SMA IIRYIADKHN MLGGCPKERA EISMLEGAVL DIRYGVSR IA YSKDFETLKV
121 DFLSKLPEML KMFEDRLCHK TYLNGDHVTH PDFMLYDALD VVLYMDPMCL DAFPKLVCFK
181 KRIEAI PQID KYLKSSKYIA WPLQGWQATF GGGDHPPKSD IEGRGIPRSA MPKKKPTPIQ
241 LNPAPDGS AV NGTSSAETNL EALQKKLEEL ELDEQQRKRL EAFLTQKQKV GELKDDDFEK
301 ISELGAGNGG VVFKVSHKPS GLVMARKLIH LEIKPAIRNQ IIRELQVLHE CNSPYIVGFY
361 GAFYS DGEIS ICMEHMDGGS LDQVLK KAGR IPEQILGKVS IAVIKGLTYL REKHKIMHRD
421 VKPSN ILVNS RGEIKLCDFG VSGQLIDSMA NSFVGTRSYM SPERLQGHY SVQSDIWSMG
481 LSLVEMAVGR YPIPPDAKE LELMFGCQVE GDA AETPPRP RTPGRPLSSY GMDSRPPMAI
541 FELLDIYVNE PPPKLPSGVF SLEFQDFV NK CLIKNPAERA DLKQLMVHAF IKRSDAEEVD
601 FAGWLCSTIG LNQPSTPTHA AGVHHHHHH
  
```

Recombinant MEK1 nucleotide sequence:

```

1 atgtccccta tactaggtta ttgga a aatt aagggccttg tgcaaccac tcgacttctt
61 ttggaat atc ttgaagaaaa atatgaagag cttttgtatg agcgcgatga aggtgataaa
121 tggcga aaca aaaagtttga attgggtttg gagtttccca atcttcctta ttatattgat
181 ggtgatg tta aattaacaca gtctatggcc atcatacgtt atatagctga caagcacaac
241 atgttgg gttgtccaaa agagcgtgca gagatttcaa tgcttgaagg agcgggtttg
301 gatattag at acggtgtttc gagaattgca tatagtaaag actttgaaac tctcaaagtt
361 gattttct ta gcaagctacc tgaaatgctg aaaatgctcg aagatcgttt atgtcataaa
421 acatattt aa atggtgatca tgtaaccat cctgacttca tgttgtatga cgctcttgat
481 gttgtttt at acatggacc c aatgtgctg gatgcttcc caaaattagt ttgttttaa
541 aaacgtat tg aagctatccc acaaattgat aagtacttga aatccagcaa gtatatagca
601 tggccttt gc agggctggca agccacgttt ggtggtggcg accatcctcc aaaatcggat
661 atcgaagg tc gtgggatccc ccgatccgcc atgccaaaga agaagcccac ccccatccag
721 ctgaatc ctg cccctgacgg ctcggcggtg aacggcacca gctcggcgga gaccaacctg
781 gaggcctt gc agaagaagct ggaggagctg gagcttgacg agcagcagcg gaagcgcctg
841 gaggcctt cc tcaccagaa gcagaaagtg ggagagctga aggacgatga cttcagagaag
901 atcagtga gc tgggagccgg caacggcggc gtggtgttca aggtctccca caagccgtcc
961 ggcctggt ca tggccagaaa gcttattcac ctggagatca aacctgctat ccggaaccag
1021 atcataagg g agctgcaggt tctgcacgag tgcaactccc cgtacatcgt gggcttctac
1081 ggggcatt ct acagcgatgg cgagatcagc atctgcatgg agcacatgga cgggggttcc
1141 ttggatca ag tcctgaagaa agctggacgg attcccagc aaattttggg gaaagtttagc
1201 attgctgt ga tcaagggcct gacgtatctg agggagaagc acaagatcat gcacagagac
1261 gtcaagcc ct ccaacatcct ggtcaactcc cgcggggaga tcaagctctg tgacttcggg
1321 gtcagtgg gc agctcatcga ctccatggcc aactccttcg tgggcaccag gtcttatatg
1381 tcgccgga ga gactccagg gacacactac tctgtgcagt cggacatctg gagcatgggg
1441 ctgtccct gg tggagatggc ggtggggcgg taccatcc cgcccccgga cgccaaggag
1501 ctggagct ga tgtttgggtg ccaggtggag ggcgatgcgg ccgagactcc gccaggccc
1561 aggaccctg ggcggcccct cagctcgtat ggaatggata gccggcctcc catggcgatt
  
```

Certificate of Analysis

```
1621 ttgagctgc tggattacat cgtcaatgag cctcctccga aactccccag cggagtcttc
1681 agcctggagt ttcaagattt tgtgaataaa tgcttaataa aaaaccccgc cgagagagca
1741 gacttgaagc agctcatggt tcatgctttt atcaagaggt ctgatgccga ggaggtggat
1801 ttgctggtt ggctgtgctc caccatcggc cttaccagc ccagcacgcc gacgcacgcg
1861 gccggtgtgc atcatcacca ccatcactaa
```

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.