

Certificate of Analysis

MAP kinase 1, active

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

Item # 14-439, 14-439-K, 14-439M

Parent Lot # 2492972

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: Recombinant, full-length, human MAP kinase 1 containing an *N*-terminal GST-tag. Expressed in *E.coli* cells and purified by glutathione-agarose. Activated with MEK1 and re-purified on Ni²⁺/NTA agarose. Purity = 88.7% by SDS-PAGE and Coomassie blue staining. MW = 69.9kDa.

Specific Activity (Parent lot# 2492972) 896U/mg, where one unit of MAP kinase 1 activity is defined as 1nmol phosphate incorporated into 0.33mg/ml myelin basic protein (MBP) per minute at 30°C with a final ATP concentration of 100µM.

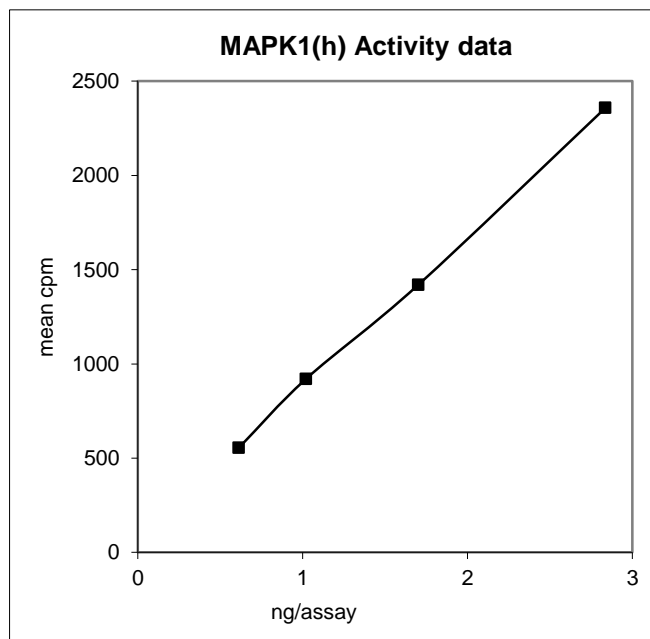
Formulation: 3.414mg/ml of enzyme in 20mM MOPS/NaOH pH7.2, 25mM sodium-β-glycerophosphate, 5mM EGTA, 1mM sodium orthovanadate, 1mM dithiothreitol, 50 % glycerol. Liquid at -20°C.

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: 0.6–2.8ng of this lot of enzyme phosphorylated 0.33mg/ml MBP 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 product identity as MAPK1 with the translated sequence listed on page three.



Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

- 1. 5 x Reaction Buffer:** 125mM Tris/HCl pH7.5, 0.1mM EGTA.
- 2. Myelin Basic Protein (MBP) substrate:** Use a final assay concentration of 0.33mg/ml. Make up a 3.3 mg/ml stock. Use 2.5µl of stock per assay point.
- 3. MAP kinase 1, active:** Dilute with 50mM Tris/HCl pH7.5, 0.1mM EGTA, 0.1mM Na₃VO₄, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 0.6–2.8ng per assay point.
- 4. [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.)

Assay Procedure (96 well plate format):

1. Add 5µl of 5 x reaction buffer per assay to wells.
2. Add 2.5µl of **MBP substrate**.
3. Add **2.5µl (0.6–2.8ng) MAP kinase 1, active**.
4. Make up to 15µl with dH₂O.
5. Add 10µl of 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. 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.
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 1µl of 30% phosphoric acid.

Certificate of Analysis

MAP kinase 1 Sequence Information

<u>Protein</u>	Human MAP kinase 1
<u>Tags</u>	N-terminal GST
<u>Native sequence</u>	M231 of the fusion protein is equivalent to M1 of human MAP kinase 1
<u>Accession number</u>	GenBank XM_055766

Recombinant MAP kinase 1 amino acid sequence:

```

1 MSPILGYWKI KGLVQPTRLL LEYLEEKYEE HLYERDEGDK WRNKKFELGL EFPNLPYYID
61 GDVKLTQSMA IIRYIADKHN MLGGCPKERA EISMLEGAVL DIRYGVSRIA YSKDFETLKV
121 DFLSKLPEML KMFEDRLCHK TYLNGDHVTH PDFMLYDALD VVLYMDPMCL DAFPKLVCFK
181 KRIEAIQID KYLKSSKYIA WPLQGWQATF GGGDHPPKSD LVPRGSPGIP MAAAAAQGGG
241 GGEPRRTEGV GPGVPGEVEM VKGQPFVGV RYTQLQYIGE GAYGMVSSAY DHVRKTRVAI
301 KKISPFEHQT YCQRTLREIQ ILLRFRHENV IGIRDILRAS TLEAMRDVYI VQDLMETDLY
361 KLLKSQQLSN DHICYFLYQI LRGLKYIHS NVLHRDLKPS NLLINTTCDL KICDFGLARI
421 ADPEHDHTGF LTEYVATRWY RAPEIMLNSK GYTKSIDIWS VGCILAEMLS NRPIFPKGHY
481 LDQLNHILGI LGSPSQEDLN CIINMKARNY LQSLPSKTKV AWAKLFPKSD SKALDLLDRM
541 LTFNPNKRIT VEEALAHPYL EQYYDPTDEP VAEEPFTFAM ELDDLPERL KELIFQETAR
601 FQPGVLEAP
  
```

Recombinant MAP kinase 1 nucleotide sequence:

```

1 atgtccccta tactaggtta ttggaaaatt aagggccttg tgcaaccac tcgacttctt
61 ttggaatatt ttgaagaaaa atatgaagag ctttgtatg agcgcgatga aggtgataaa
121 tggcgaaca aaaagtttga attgggtttg gagtttccca atcttcctta ttatattgat
181 ggtgatgta aattaacaca gtctatggcc atcacatgctt atatagctga caagcacaac
241 atgttgggtg gttgtccaaa agagcgtgca gagatttcaa tgcttgaagg agcggttttg
301 gatattagat acggtgtttc gagaattgca tatagtaaag actttgaaac tctcaaagtt
361 gattttctta gcaagctacc tgaatgctg aaaatgttcg aagatcgttt atgtcataaa
421 acatatttaa atggtgatca tgtaaccat cctgacttca tgttgtatga cgctcttgat
481 gttgttttat acatggacc aatgtgcctg gatgcgttcc caaaattagt ttgttttaaa
541 aaacgtattg aagctatccc acaaattgat aagtacttga aatccagcaa gtatatagca
601 tggcctttgc agggctggca agccacgttt ggtgggtggcg accatcctcc aaaatcggat
661 ctgcttccgc gtgcatcccc gggaaattccc atggcggcgg cggcggctca gggggcgggg
721 ggcgggggagc cccgtagaac cgagggggtc ggcccggggg tcccggggga ggtggagatg
781 gtgaaggggc agcgttcga cgtgggcccg cgctacacgc agttgcagta catcggcgag
841 ggcggtacg gcatggtcag ctcgccctat gaccacgtgc gcaagactcg cgtggccatc
901 aagaagatca gcccttcga acatcagacc tactgccagc gcacgctccg ggagatccag
961 atcctgctgc gcttccgcca tgagaatgtc atcggcatcc gagacattct gcgggcttcc
1021 accctggaag ccatgagaga tgtctacatt gtgcaggacc tgatggagac tgacctgtac
1081 aagttgctga aaagccagca gctgagcaat gaccatatct gctacttctt ctaccagatc
1141 ctgcggggcc tcaagtacat ccactccgcc aacgtgctcc accgagatct aaagccctcc
1201 aacctgctca tcaacaccac ctgcgacctt aagatttgtg atttcggcct ggcccggatt
1261 gccgatcctg agcatgacca caccggcttc ctgacggagt atgtggctac gcgctggtac
1321 cgggccccag agatcatgct gaactccaag ggctatacca agtccatcga catctggtct
1381 gtgggctgca ttctggctga gatgctctt aaccggcca tcttccctgg caagcactac
1441 ctggatcagc tcaaccacat tctgggcatc ctgggctccc catcccagga ggacctgaat
1501 tgtatcatca acatgaaggc ccgaaactac ctacagtctc tgccctcaa gaccaaggtg
1561 gcttgggcca agcttttccc caagtcagac tcaaagccc ttgacctgct ggaccggatg
1621 ttaaccttta accccaataa acggatcaca gtggaggaag cgctggctca cccctacctg
  
```

Certificate of Analysis

1681 gagcagtact atgacccgac ggatgagcca gtggccgagg agcccttcac cttcgccatg
1741 gagctggatg acctacctaa ggagcggctg aaggagctca tcttcagga gacagcacgc
1801 ttccagcccg gagtgctgga ggccccctag

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