

Certificate of Analysis

MKK4/SKK1, unactive

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

Item # 14-378, 14-378-K, 14-378M

Parent Lot # 30901U

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, mouse MKK4, amino acids 34–end expressed in *E.coli* cells. Purified using glutathione-agarose. Purity 50%. MW = 67.8kDa.

Formulation: 3.945mg/ml of enzyme in 50mM Tris/HCl pH7.5, 150mM NaCl, 50% glycerol, 1mM benzamidine, 0.2mM PMSF, 0.1mM EGTA, 0.1% 2-mercapoethanol, 0.03% Brij-35. Liquid at -20°C.

Specific Activity (Parent lot# 30901U): As provided, this lot demonstrated <2.5% of maximum activity. Activated by phosphorylation with MEKK (cat# 14-196).

Storage and Stability: On receipt of material store at -20°C. Unopened reagent is stable for a minimum 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.

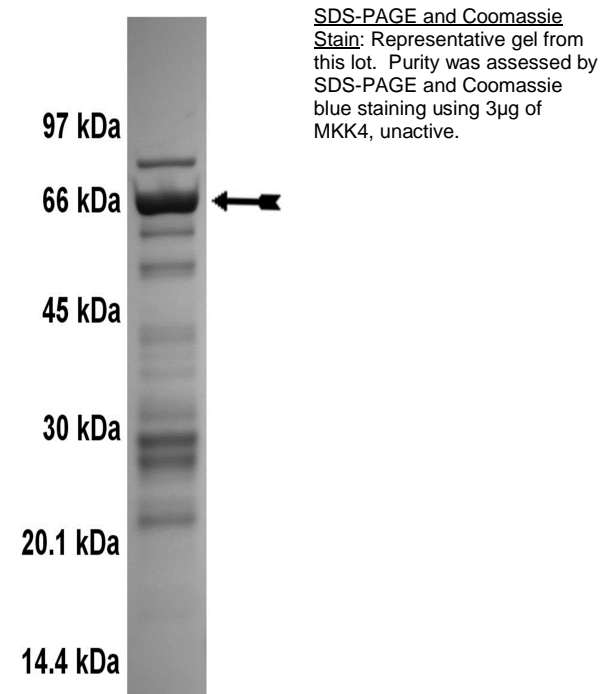
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Activation Assay: 4µM of MKK4, unactive was activated using 0.1mg/ml MEK kinase, then diluted 4-fold, used to activate JNK1α1, and the increased activity of the JNK 1α1 against ATF-2 determined. The activation of MKK4, subsequent activation of JNK 1α1, and assay are described on page two. Results of this assay are shown below.

Active MEKK	Unactive MKK4	Mean cpm	Comment
None	6.8µg	531	MKK4 unactive
2.5µg	6.8µg	22761	Activated MKK4

MS Tryptic Fingerprint: Confirmed identity as MKK4 with the translated native sequence listed on page four.



Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

1. **5 x Activation Buffer:** 250mM Tris/HCl pH7.5, 0.5mM EGTA, 0.5% 2-mercaptoethanol.
2. **10 x Assay Buffer:** 500mM Tris/HCl pH7.5, 1mM EGTA, 1mM Na₃VO₄, 1% 2-mercaptoethanol.
3. **Dilution Buffer:** 50mM Tris/HCl pH7.5, 0.1mM EGTA, 0.1mM Na₃VO₄, 0.1% 2-mercaptoethanol 1mg/ml BSA.
4. **Stages One and Two 5 x Mg/ATP:** 50mM magnesium acetate, 0.5mM ATP.
5. **MEKK, active (Catalogue# 14-196):** Use at a final assay concentration of 0.1mg/ml. Dilute with Dilution Buffer to 1.00mg/ml. Add 2.5µl of stock per assay point.
6. **MKK4, unactive:** Use at a final assay concentration of 4µM (0.271mg/ml). Dilute with Dilution Buffer to 2.71mg/ml. Use 2.5µl of stock per assay point.
7. **JNK1α1, unactive (Catalogue#14-328):** Use at a final assay concentration of 2µM (0.09mg/ml). Dilute with Dilution Buffer to 0.45mg/ml. Add 5µl of stock per assay point.
8. **ATF-2 (Catalogue# 12-367):** Use at a final assay concentration of 3µM (0.108mg/ml). Dilute with Dilution Buffer to 1.08mg/ml. Add 2.5µl of stock per assay point.
9. **[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 500 - 800cpm/pmol as required.)

Assay Procedure:

Stage One: *Activation of MKK4*

1. Add 5µl of 5 x activation buffer to a microcentrifuge tube.
2. Add **2.5µl (6.8µg) of MKK4 unactive.**
3. Add 2.5µl (2.5µg) **MEKK, active.**
4. Add 10µl of dH₂O.
5. Add 5µl stages one and two 5 x Mg/ATP.
6. In appropriate controls, add dilution buffer to a final volume of 25µl.
7. Incubate for 60 minutes at 30°C.
8. Stop the reaction by diluting 4-fold in dilution buffer. Store on ice.

Stage Two: *Activation of JNK 1α1.*

1. Add 5µl 5 x activation buffer to a microcentrifuge tube.
2. Add 5µl (2.25µg) of **JNK 1α1 unactive.**
3. Add 2.5µl of activated MKK4 from **Stage One.**
4. Add 7.5µl of dH₂O.
5. Add 5µl of stages one and two 5 x Mg/ATP.
6. Incubate for 15 minutes at 30°C.
7. Immediately transfer 2.5µl of the reaction product to the **Stage Three** component mixture.

Certificate of Analysis

Stage Three: *Phosphorylation of ATF-2 by Activated JNK 1 α 1.*

1. Add 2.5 μ l 10 x assay buffer to a microcentrifuge tube.
2. Add 2.5 μ l (2.7 μ g) **ATF-2**.
3. Add 7.5 μ l of dH₂O.
4. Add 10 μ l of the 2.5 x magnesium acetate/ [γ -³³P]ATP cocktail.
5. Add 2.5 μ l of JNK 1 α 1 from **Stage Two** to start the reaction.
6. Incubate for 15 minutes at 30°C.
7. Spot 20 μ l onto the centre of a 2cm x 2cm **P81** paper.
8. Wash the assay squares twice for 5 minutes with 75mM phosphoric acid.
9. Wash the assay squares once for 2 minutes with acetone.
10. Transfer the assay squares to a scintillation vial and add 1ml of scintillation cocktail.
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 1 μ l of 30% phosphoric acid.

Certificate of Analysis

MKK4 Sequence Information

Protein	Mouse MKK4
Tags	N-terminal GST
Native sequence	S226 of the recombinant protein corresponds to S34 of mouse MKK4
Accession number	EMBL U18310

Recombinant MKK4 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 KRIEAIPOID KYLKSSKYIA WPLQGWQATF GGDHPPKSD LVPRGSMQ GK RKALKLN FAN
241 PPVKSTARFT LNPNTTGVQN PHIERLRTHS IESSGKLIKIS PEQHWDF TAE DLKDLGEIGR
301 GAYGSVNKMV HKPSGQIMAV KRIRSTVDEK EQQLLMDLD VVMRSSDCPY IVQFYGALFR
361 EGDCWICMEL MSTSFDFKYK YVYSVLDDVI PEEILGKITL ATVKALNHLK ENLKI IHRDI
421 KPSNILLDRS GNIKLCDFGI SGQLVDSIAK TRDAGCRPYM APERIDPSAS RQGYDVRSDV
481 WSLGITLYEL ATGRFPYPKW NSVFDQLTQV VKGDPPQLSN SEEREFSPSF INFVNLCLTK
541 DESKRPKYKE LLKHPFILMY EERTVEVACY VCKILDQMPA TPSSPMYVD
  
```

Recombinant MKK4 nucleotide sequence:

```

1 atgtccccta tactaggtta ttggaaaatt aagggccttg tgcaaccac tcgacttctt
61 ttggaatata ttgaagaaaa atatgaagag catttgtatg agcgcgatga aggtgataaa
121 tggcgaaaaca aaaagtttga attgggtttg gagtttccca atcttcctta ttatattgat
181 ggtgatgtta aattaacaca gtctatggcc atcatacgtt atatagctga caagcacaac
241 atgttgggtg gttgtccaaa agagcgtgca gagatttcaa tgcttgaagg agcgggtttg
301 gatattagat acggtgttcc gagaattgca tatagtaaag actttgaaac tctcaaagtt
361 gattttctta gcaagctacc tgaaatgctg aaaatgttctg aagatcgttt atgtcataaa
421 acatatttaa atggtgatca tgtaacccat cctgacttca tgttgtatga cgctcttgat
481 gttgttttat acatggacc aatgtgcctg gatgcttcc caaaattagt ttgttttaaa
541 aaacgtattg aagctatccc acaaattgat aagtacttga aatccagcaa gtatatagca
601 tggcctttgc agggctggca agccacgttt ggtgggtggcg accatcctcc aaaatcggat
661 ctggttccgc gtggatccat gcagggttaag cgcaaagcac tgaagttgaa ttttgcaaat
721 ccacctgtca aatcgacagc acggtttacc ctgaatccta atactacagg agtccagaac
781 ccacacatag agagactgag aacacacagc attgagtcat caggaaaact gaagatctcc
841 cctgaacaac actgggattt cactgcagag gacttgaaag accttgagaa aattggacga
901 ggagcttatg gttctgtcaa caaaatggtc cacaaaacaa gtgggcagat aatggcagtt
961 aaaagaattc ggtcaactgt ggatgaaaaa gaacaaaaac aacttcttat ggatttggat
1021 gtagtaaatgc ggagtagtga ttgccatac attgttcagt tctatggtgc actcttcaga
1081 gagggcgact gttggatctg tatggagctc atgtctacct cgttcgataa gttttacaaa
1141 tatgtatata gtgtgttaga tgacgttatt ccggaagaga tcttaggcaa aatcacttta
1201 gcaactgtga aagcactaaa ccacttaaaa gaaaacttga aaattattca cagagacatc
1261 aaaccttcca atattcttct ggacagaagt ggaaatataa agctctgtga tttcggcatc
1321 agtggacagc ttgtggactc tattgccaag acaagagatg ctgggtgtag gccgtatatg
1381 gcacctgaaa gaatagacc aagtcatca agacaagggg atgatgtccg cctgatgtc
1441 tggagtttgg ggatcacatt gtacgagttg gccacaggcc gatttctta tccaaagtgg
1501 aatagtgtat ttgatcagct aacacaagtg gtgaaaggag accctccgca gctgagtaat
1561 tctgaagaaa gggagttctc cccagtttc atcaactttg tcaacttctg ccttacgaag
1621 gatgaatcca aaaggccaaa gtataaagag cttctgaaac atcccttat tttgatgat
  
```

Certificate of Analysis

1681 gaagaacgta ctgtagaggt cgcattgctat gtttgtaaaa tcttgatca gatgccagcc
1741 actcccagct cgccatgta tctcgactga

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.