

Certificate of Analysis

MKK6/SKK3, unactive

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

Item # 14-304, 14-304-K, 14-304M

Parent Lot # 1604418

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 maltose binding protein tagged, recombinant MKK6 amino acids 4–end, expressed in *E.coli* cells. Purified using amylose agarose. Purity 99% by SDS-PAGE and Coomassie blue staining. MW = 80kDa.

Formulation: 3.19mg/ml of enzyme in 50mM Tris/HCl pH7.5, 150mM NaCl, 0.1mM EGTA, 0.1% 2-mercaptoethanol, 0.03% Brij-35, 20% glycerol. Frozen solution.

Specific Activity (Parent lot# 1604418): As provided, this lot demonstrated 3% of maximum activity. Activated by phosphorylation with MEKK (cat# 14-196).

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.

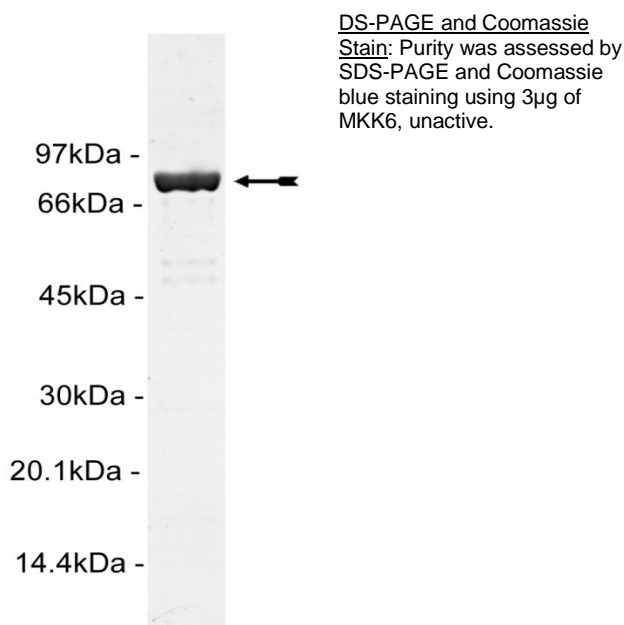
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Activation Assay: 4µM MKK6, unactive was activated using 200µg/ml MEKK (Catalogue# 14-196) which in turn was used to activate 2µM SAPK2a, and the increased activity against MBP determined. The activation and assay are described on pages two and three. Results of this assay are shown below.

MS Tryptic Fingerprint: Confirmed product identity as MKK6 with the translated sequence listed on page four

Active MEKK	Unactive MKK6	Mean cpm	Comments
5µg	8.06µg	20658	Kinase activity
None	8.06µg	914	Background



Certificate of Analysis

Kinase Assay Protocol

Stock Solutions

1. **10 x Activation Buffer:** 500mM Tris/HCl pH7.5, 1mM EGTA, 1% 2-mercaptoethanol, 1mM Na₃VO₄, 10mg/ml BSA.
2. **10 x Reaction Buffer:** 250mM Tris/HCl pH7.5, 1mM EGTA.
3. **Enzyme Dilution Buffer:** 50mM Tris/HCl pH7.5, 0.1mM EGTA, 0.1% 2-mercaptoethanol, 0.1mM Na₃VO₄, 1mg/ml BSA.
4. **5 x Mg/ATP (Stages One and Two):** 50mM MgAc, 0.5mM ATP.
5. **MEKK:** Use at a final assay concentration of 200µg/ml. Prepare a 1mg/ml stock. Use 5µl of stock per assay point.
6. **MKK6 unactive:** Use at a final assay concentration of 4µM (0.322mg/ml). Prepare a 1mg/ml stock. Use 8µl of stock per assay point.
7. **SAPK2a unactive:** Use at a final assay concentration of 2µM (0.135mg/ml). Prepare a 0.675mg/ml stock. Use 5µl of stock per assay point.
8. **MBP:** Use at a final assay concentration of 0.33mg/ml. Prepare a 3.33mg/ml stock. Use 2.5µl of stock per assay point.
9. **[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:

Stage One: *Activation of MKK6 by MEKK*

1. Add 2.5µl of 10 x activation buffer to a microcentrifuge tube.
2. Add 5µl of **MEKK**.
3. Add **8µl MKK6, unactive**.
4. Add 4.5µl of dH₂O.
5. Add 5µl of stage one Mg/ATP mixture. (In appropriate controls add dilution buffer to a final volume of 25µl).
6. Incubate for 1 hour at 30°C.
7. Stop the reaction by diluting the MKK6 250-1000 fold and storing on ice.

Stage Two: *Activation of SAPK2a by MKK6*

1. Add 2.5µl of 10 x activation buffer to a microcentrifuge tube.
2. Add 5µl of **SAPK2a unactive**.
3. Add **2.5µl diluted MKK6 (activated) from Stage One**.
4. Add 10µl of dH₂O.
5. Add 5µl of stage two Mg/ATP mixture.
6. Incubate at 30°C for 15 minutes
7. Stop reaction by immediately transferring 5µl into **Stage Three** reaction mixture.

Certificate of Analysis

Stage Three: *Phosphorylation of MBP by SAPK2a*

1. Add 2.5µl of reaction buffer to a microcentrifuge tube.
2. Add 2.5µl of **MBP**.
3. Add 5µl of **activated SAPK2a from Stage Two**.
4. Add 5µl of dH₂O.
5. Add 10µl of the diluted [γ -³³P]ATP mixture.
6. Incubate for 15 minutes at 30°C.
7. Transfer a 20µl aliquot onto the centre of a 2cm x 2cm **P81** paper square.
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 vial 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 plus 1µl of 30% phosphoric acid.

Certificate of Analysis

MKK6 (4-end) Sequence Information

Protein	human MKK6 (4-end)
Tags	N-terminal maltose binding protein
Native sequence	S396 of the recombinant protein is equivalent to S4 of human MKK6
Accession number	EMBL U39657

Recombinant MKK6 amino acid sequence:

```

1 MKTEEGKLVI WINGDKGYNG LAEVGKKFEK DTGIKVTVEH PDKLEEKFPQ VAATGDGPDI
61 IFWAHDRFGG YAQSGLLAEI TPKAFQDKL YPFTWDAVRY NGKLIAYPIA VEALSIIYNK
121 DLLPNPPKTW EEIPALDKEL KAKGKSALMF NLQEPYFTWP LIAADGGYAF KYENKDYDIK
181 DVGVDNAGAK AGLTFLVDLI KNKHMNADTD YSIAEAAFNK GETAMTINGP WAWSNIDTSK
241 VNYGVTVLPT FKGQPSKPFV GVLSAGINAA SPNKELAKEF LENYLLTDEG LEAVNKDKPL
301 GAVALKSYEE ELAKDPRIAA TMENAQGEI MPNIPQMSAF WYAVRTAVIN AASGRQTVDE
361 ALKDAQTNSS SNNNNNNNNN NLGIEGRISE FGSSRSKGKK RNPGLKIPKE AFEQPQTSST
421 PPRDLDSKAC ISIGNQNFV KADDLEPIME LGRGAYGVVE KMRHVPSGQI MAVKRIRATV
481 NSQEQKRLLM DLDISMRTVD CPFTVTFYGA LFREGDVVIC MELMDTSLDK FYKQVIDKGQ
541 TIPEDILGKI AVSIVKALEH LHSKLSVIHR DVKPSNVLIN ALGQVKMCDF GISGYLVDSV
601 AKTIDAGCKP YMAPERINPE LNQKGYSVKS DIWSLGITMI ELAILRFPYD SWGTPFQQLK
661 QVVEEPSPQL PADKFSAEFV DFTSQCLKKN SKERPTYPEL MQHPFFTLHE SKGTDVASFV
721 KLILGD
  
```

Recombinant MKK6 nucleotide sequence:

```

1 atgaaaactg aagaaggtaa actgtaatc tggattaacg gcgataaagg ctataacggt
61 ctcgctgaag tcggtaaгаа attcgagaaa gataccggaa ttaaagtcac cgttgagcat
121 ccggataaac tggaaгagaa attcccacag gttgcggcaa ctggcgatgg ccctgacatt
181 atcttctggg cacacgaccg ctttggtggc tacgctcaat ctggcctggt ggctgaaatc
241 acccggaca aagcgttcca ggacaagctg tatccgttta cctgggatgc cgtacgttac
301 aacggcaagc tgattgctta cccgatcgct gttgaagcgt tatcgctgat ttataacaaa
361 gatctgctgc cgaaccgccc aaaaacctgg gaagagatcc cggcgctgga taaagaactg
421 aaagcгааag gtaagagcgc gctgatgttc aacctgcaag aaccgtactt cacctggccg
481 ctgattgctg ctgacggggg ttatgcgttc aagtatgaaa acggcaagta cgacattaaa
541 gacgtgggcg tggataacgc tggcgcgaaa gcgggctctga cttcctggtg tgacctgatt
601 aaaaacaaac acatgaatgc agacaccgat tactccatcg cagaagctgc ctttaataaa
661 ggcгааacag cgatgaccat caacggcccг tgggcatggt ccaacatcga caccagcaaa
721 gtgaattatg gtgtaacggт actgccgacc ttcaagggtc aaccatccaa accgttcggt
781 ggcgtgctga gcgcaggtat taacgccgcc agtccgaaca aagagctggc aaaagagttc
841 ctcгaaaact atctgctgac tgatgaaggt ctggaagcgg ttaataaaga caaaccgctg
901 ggtgccgtag cgtgaagtc ttacgaggaa gagttggcga aagatccacg tattgccgcc
961 accatggaaa acgccagaa aggtgaaatc atgccgaaca tcccgcagat gtccgctttc
1021 tggtatgccg tgcgtactgc ggtgatcaac gccgccagcg gtcgtcagac tgtcgatgaa
1081 gccctгааag acgгcгagac taattcgagc tcгaaacaaca acaacaataa caataacaac
1141 aacctгggga tcгaggгaaг gatttcгaa тtcгgatcct ctagatcгaa agгcaagaag
1201 cгaaaccctг gccttaaaat tccaaaгaa гcatttгaaс aacctгagac cгttccaca
1261 ccacctгgag atttagactc caaggctгc atttctattg gaaatcгaa ctttgagggtg
1321 aaggcagatg acctгgagcc tataatггaa ctgggacгag gtгcгtacгg ggtгgtгgгag
1381 aagatгcггc acгtгcccag cgggгagatc atггcгagtг agcгgгatccг agccacagta
1441 aatagccagг aacгaaacг gctactгatг gatttгgгata tttccatгag gacгgtгgгac
1501 gtгccattca ctгtcacctt ttatггcгca ctгtttcггg agгgtгatгt gtгgгatctгc
1561 atгgгagctca tгgгatacatc actagataaa ttctacaaac aagttattgг taaagгccag
  
```

Certificate of Analysis

```
1621 acaattccag aggacatctt agggaaaata gcagtttcta ttgtaaaagc attagaacat
1681 ttacatagta agctgtctgt cattcacaga gacgtcaagc cttctaattgt actcatcaat
1741 gctctcggtc aagtgaagat gtgcgatfff ggaatcagtg gctacttggg ggactctggt
1801 gctaaaacia ttgatgcagg ttgcaaacca tacatggccc ctgaaagaat aaaccagag
1861 ctcaaccaga agggatacag tgtgaagtct gacatttggg gtctgggcat cacgatgatt
```

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