

## Certificate of Analysis

### MSK1 (2–end), active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-438, 14-438-K, 14-438M

Parent Lot # D8AN041U

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 6His-tagged recombinant human MSK1 residues 2–end. Expressed by baculovirus in Sf21 insect cells. Purified using Ni<sup>2+</sup>/NTA agarose, activated using human MAPK2, and then re-purified using Ni<sup>2+</sup>/NTA agarose. Purity 93.5% by SDS-PAGE and Coomassie blue staining. MW = 94.4kDa.

**Specific Activity (Parent lot# D8AN041U):** 734U/mg, where one unit of MSK1, active activity is defined as 1nmol phosphate incorporated into 30µM modified crosstide (GRPRTSSFAEGKK) per minute at 30°C with a final ATP concentration of 100µM.

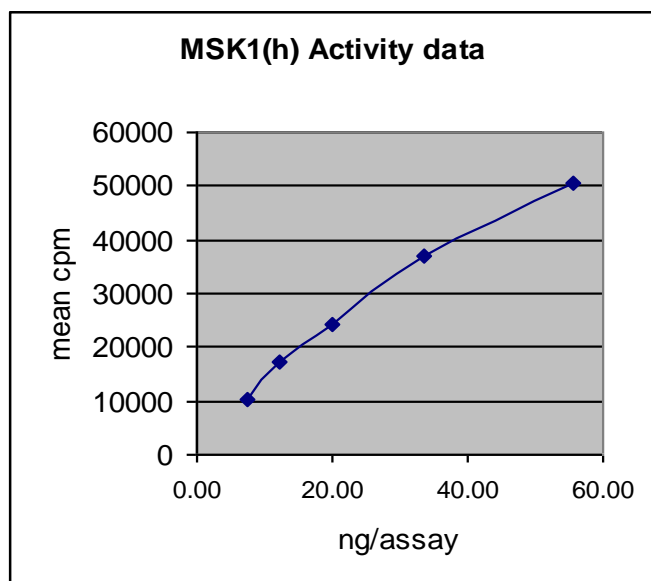
**Formulation:** 2.222mg/ml of enzyme in 50mM Tris/HCl pH7.5, 50% glycerol, 150mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 1mM benzamidine, 0.2mM PMSF, 0.1 % 2-mercaptoethanol. Frozen solution.

**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**

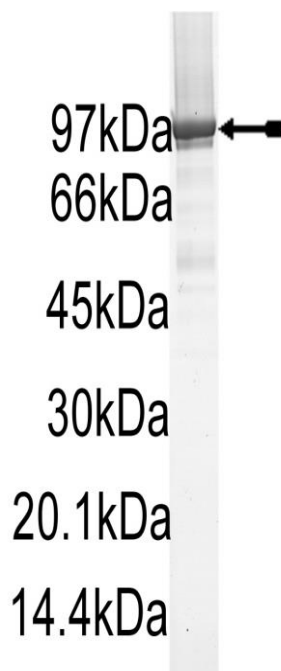
### Quality Control Testing

**Kinase Assay:** 7–55ng of this lot of enzyme phosphorylated 30µM modified crosstide 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 identity as MSK1 with the translated sequence listed on page three.

**SDS-PAGE and Coomassie Stain:** Purity was assessed by SDS-PAGE and Coomassie blue staining using 3µg of MSK1, active.



## Certificate of Analysis

### Kinase Assay Protocol

#### Stock Solutions:

1. **5 x Reaction Buffer:** 40mM MOPS/NaOH pH7.0, 1mM EDTA.
2. **Modified Crosstide (GRPRTSSFAEGKK):** Use at a final assay concentration of 30 $\mu$ M. Prepare a 300 $\mu$ M stock and add 2.5 $\mu$ l of stock per assay point.
3. **MSK1, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 5% glycerol, 0.01% Brij-35, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 7–55ng per assay point.
4. **[ $\gamma$ -<sup>33</sup>P]ATP:** 2.5 x magnesium acetate/[ $\gamma$ -<sup>33</sup>P]ATP cocktail: 25mM MgAc and 0.25mM ATP to which is added [ $\gamma$ -<sup>33</sup>P]ATP (specific activity approximately 500 - 800cpm/pmol as required.)

#### Assay Procedure (96 well plate format):

1. Add 5 $\mu$ l of 5 x reaction buffer per assay to wells.
2. Add 2.5 $\mu$ l of **modified crosstide**.
3. Add **2.5 $\mu$ l (7–55ng) MSK1, active**.
4. Add 5 $\mu$ l of dH<sub>2</sub>O.
5. Add 10 $\mu$ l of diluted [ $\gamma$ -<sup>33</sup>P] ATP mixture.
6. Incubate for 10 minutes at 30°C.
7. Stop reaction by adding 5 $\mu$ l of 3% phosphoric acid.
8. Transfer a 10 $\mu$ l aliquot onto the appropriate area of a **P30 Filtermat**.
9. Wash the filtermat three times for 5 minutes with 50mM 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 1 $\mu$ l of 30% phosphoric acid.

## Certificate of Analysis

### MSK1 Sequence Information

<b>Protein</b>	human MSK1 (2–end)
<b>Tags</b>	N-terminal 6His
<b>Native sequence</b>	E39 of the recombinant protein is equivalent to E2 of human MSK1
<b>Accession number</b>	GenBank AF074393

#### Recombinant MSK1 amino acid sequence:

```

1 MSYYHHHHHH DYDIPTTENL YFQGAMGSAT MDYKDDDDKE EEGSSGGAA GTSADGGDGG
61 EQLLTVKHEL RTANLTGHAE KVGIENFELL KVLGTGAYGK VFLVRKISGH DTGKLYAMKV
121 LKKATIVQKA KTTEHTRTER QVLEHIRQSP FLVTLHYAFQ TETKLHLILD YINGGELFTH
181 LSQRERFTEH EVQIYVGEIV LALEHLHKLK IYRDIKLEN ILLDSNGHVV LTDFGLSKEF
241 VADETERAYS FCGTIEYMAP DIVRGGDSGH DKAVDWWSLG VLMYELLTGA SPFTVDGEKN
301 SQAEISRRIL KSEPPYPQEM SALAKDLIQR LLMKDPKKRL GCGPRDAEI KEHLFFQKIN
361 WDDLAAKKVP APFKPVIRDE LDVSNFAEEF TEMDPTYSPA ALPQSSEKLF QGYSFVAPSI
421 LFKRNAVID PLQFHMVER PGVTNVARSA MMKDSPFYQH YDLDLKDKPL GEGSFSICRK
481 CVHKSNQAF AVKIIISKRME ANTQKEITAL KLCEGHPNIV KLHEVFHDQL HTFLVMELLN
541 GGELFERIKK KKHFSETEAS YIMRKLVS AVSHMHVGVVH RDLKPENLLF TDENDNLEIK
601 IIDFGFARLK PPDNQPLKTP CFTLHYAAPE LLNQNGYDES CDLWSLGVIL YTMLSGQVVP
661 QSHDRSLTCT SAVEIMKKIK KGDFSFEAEA WKNVSQEAKD LIQGLLTVDP NKRLKMSGLR
721 YNEWLQDGSQ LSSNPLMTPD ILGSSGA AVHTCVKATFHAF NKYKREGFCL QNVDKAPLAK
781 RRMKKTSTS TETRSSSES SHSSSSHSHG KTTPTKTLQP SNPADSNNPE TLFQFSDSVA
  
```

#### Recombinant MSK1 nucleotide sequence:

```

1 atgtcgtact accatcacca tcaccatcac gattacgata tcccaacgac cgaaaacctg
61 tattttcagg gcgccatggg atctgccacc atggactaca aggacgacga tgacaaggag
121 gaggagggtg gcagcagcgg cggcgccgcg gggaccagcg cggacggcgg cgacggagga
181 gagcagctcc tactgtcaa gcacgagctg cggactgcta atttgacagg acatgctgag
241 aagtgaggaa tagaaaatth tgagctcctg aaggtcctag gaactggagc ttatggaaaa
301 gtatttctag ttcgtaaaat aagtggccat gatactggaa agctgtatgc catgaaagtt
361 ttgaaaaaag caacaatcgt tcaaaaggcc aaaaccacag agcatacaag gacagaacga
421 caagtctcgg aacacattag gcagtcgcca tttttggtaa cattacatta tgctttccag
481 acagaaacca aacttcatct catttttagat tatataaatg gtggtgaact ttttactcat
541 ctttctcaaa gagagcgttt cacagagcat gaggtgcaga tttatggttg agagattgtg
601 cttgccctcg aacatctcca caagttgggg attatataatc gtgatattaa gcttgagaat
661 atttactctg attctaattg ccatgtggtg ctgacagatt ttggtctgag taaggagttt
721 gtggctgatg aaactgaaag agcatattcc ttttgggaa ctattgaata ctggcacca
781 gatattgtca gagggggaga ttcaggacat gacaaggcag ttgactggtg gatgtgggt
841 gttctaattg atgaattact aactggagca tctcctttca ctggtgatgg agaaaaaat
901 tcccaagctg agatatctag gagaatatta aaaagtgagc ctccatatcc ccaagaaatg
961 agtgctttag cgaaagacct aattcagcgt cttttgatga aagatcccaa gaagagattg
1021 ggatgtggtc cacgtgatgc agatgaaatc aaagaacatc tcttctttca gaaaataaat
1081 tgggatgatt tagccgcaa aaaagtgcct gcaccattta agccagtcac tcgagatgaa
1141 ttagatgtga gtaactttgc agaagagttc acagaaatgg atcccactta ttctcccgca
1201 gccctgcccc agagttctga gaagctggtt cagggtatt ctttggttgc tccttccatc
1261 ctattcaagc gtaatgcagc tgtcatagac cctcttcagt ttcacatggg agttgaacgt
1321 cctggagtga caaatggtgc caggagtgca atgatgaagg actctccatt ctatcaacac
1381 tatgacctag atttgaagga caaacccctg ggagaaggta gtttttcaat ttgtcgaaag
1441 tgtgtgcata aaaaaagtaa ccaagctttt gcagtcaaaa taatcagcaa aaggatggaa
1501 gccaatactc aaaaggaaat aacagctctg aaactctgtg aaggacaccc caatattgtg
1561 aagttgcatg aagtttttca tgatcagctt cacacgtttc tagtgatgga acttctgaat
1621 ggaggagaac tgtttgagcg cattaagaaa aagaagcact tcagtgagac ggaagccagc
  
```

## Certificate of Analysis

```
1681 tacatcatga ggaagcttgt ttcagctgta agccacatgc atgatgttgg agtgggtgac
1741 agggatctga aacctgagaa tttattgttc accgatgaaa atgacaattt ggaaattaaa
1801 ataattgatt ttggatttgc acggctaaag ccaccggata atcagcccct gaagactcca
1861 tgcttcaccc ttcattatgc cgccccagag ctcttgaatc agaacggcta cgatgagtcc
1921 tgtgacctgt ggagcttggg cgtcattttg tacacaatgt tgtcaggaca ggttcccttc
1981 caatctcatg accgaagttt gacgtgtacc agcgcggtgg aaatcatgaa gaaaattaaa
2041 aaggagatt tctcctttga aggagaagcc tggagaatg tatcccaaga ggctaaagat
2101 ttgatccaag gacttctcac agtagatcca aacaaaaggc ttaaaatgtc tggcttgagg
2161 tacaatgaat ggctacaaga tggagtcag ctgtcctcca atcctctgat gactccggat
2221 attctaggat cttccggagc tgccgtgcat acctgtgtga aagcaacctt ccacgccttt
2281 aacaaataca agagagaggg gttttgcctt cagaatgttg ataaggcccc tttggctaag
2341 agaagaaaaa tgaaaaagac tagcaccagt accgagacgc gcagcagttc cagtgagagt
2401 tcccattctt cttcctctca ttctcacggt aaaactacac ccaccaagac actgcagccc
2461 agcaatcctg ccgacagcaa taaccggag accctcttcc agttctcggg ctcagtagct
2521 tag
```

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.