

Certificate of Analysis

MSK2, unactive

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-625, 14-625-K, 14-625M

Parent Lot # D7NN001U

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 MSK2, amino acids 2–end, expressed by baculovirus in Sf21 insect cells. Purified using Ni²⁺/NTA-agarose. Purity 59.7% by SDS-PAGE and Coomassie blue staining. MW = 89.9kDa.

Formulation: 1.034mg/ml of enzyme in 50mM Tris/HCl pH7.5, 300mM NaCl, 0.1mM EGTA, 1mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol, 0.03% Brij-35, 270mM sucrose. 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# D7NN001U): as provided, this lot demonstrated <1% of maximum activity. Activated by phosphorylation with MAPK2.

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.

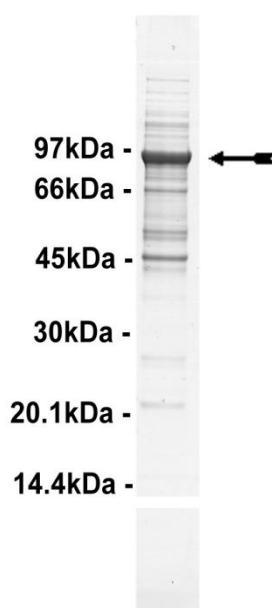
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: etc: Unactive MSK2 was activated using MAPK2 and the increased activity against Crosstide determined. The activation and subsequent assay are described on page two. Results of this assay are shown below.

MAPK2	Unactive MSK2	Mean cpm	Comments
85ng	180ng	41679	Kinase activity
None	360ng	482	Background

MS Tryptic Fingerprint: Confirmed identity as MSK2 with the translated native sequence listed on page three.



SDS-PAGE and Coomassie Stain: Purity was assessed by SDS-PAGE and Coomassie blue staining using 3µg of MSK2, unactive

Certificate of Analysis

Kinase Cascade Assay Protocol

Stock Solutions:

1. **10 x MSK2 activation buffer:** 500mM Tris/HCl pH 7.5, 1mM EGTA, 1% 2-mercaptoethanol.
2. **5 x MSK2 assay buffer:** 40mM MOPS/NaOH pH7.0, 1mM EDTA.
3. **Enzyme Dilution buffer:** Dilute in 20mM MOPS/NaOH pH7.0, 1mM EDTA, 5% glycerol, 0.01% Brij-35, 0.1% 2-mercaptoethanol, 1mg/ml BSA.
4. **Magnesium/ATP Cocktail (5 x stock):** 500 μ M cold ATP, and 50mM magnesium acetate.
5. **[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.)
6. **MAPK2, active (Catalogue# 14-173):** Use at a final assay concentration of 0.5 μ M (0.0339mg/ml). Prepare a 0.339mg/ml stock and add 2.5 μ l of stock per assay point.
7. **MSK2, inactive:** Use at a final concentration of 2 μ M (0.1798mg/ml). Prepare a 0.899mg/ml stock and add 5 μ l of stock per assay point.
8. **Modified Crosstide (GRPRTSSFAEGKK):** Use a final assay concentration of 30 μ M. Make a 300 μ M stock. Add 2.5 μ l of stock per assay point.

Assay Procedure:

Stage One: *Activation of MSK2 by MAPK2*

1. Add 2.5 μ l of MSK2 activation buffer to a microcentrifuge tube.
2. Add 2.5 μ l of **MAPK2, active**.
3. Add **5 μ l of MSK2, inactive**.
4. Add 10 μ l of dH₂O.
5. Add 5 μ l of Magnesium/ATP Cocktail.
6. Incubate for 45 minutes at 30°C.
7. Stop reaction by diluting 2.5–50 fold and placing on ice.

Stage Two: *Phosphorylation of Crosstide by activated MSK2*

1. Add 5 μ l of MSK2 assay buffer into a microcentrifuge tube.
2. Add 2.5 μ l of modified Crosstide (**GRPRTSSFAEGKK**).
3. Add 5 μ l of **Stage One** reaction product diluted 1–50 fold.
4. Add 2.5 μ l of dH₂O.
5. Add 10 μ l of the diluted [γ -³³P] ATP.
6. Incubate for 10 minutes at 30°C.
7. Slowly transfer 20 μ l onto the centre of a 2cm x 2cm **P81** paper.
8. Wash assay squares twice for 5 minutes with 75mM phosphoric acid
9. Wash assay squares once with acetone for 2 minutes.
10. Transfer assay squares to scintillation vials and add 1ml scintillation cocktail.
11. Read in scintillation counter. Compare cpm of enzyme samples with cpm of control samples that contain all appropriate assay components plus 1 μ l of 30 % phosphoric acid.

Certificate of Analysis

MSK2 Sequence Information

<u>Protein</u>	Human MSK2
<u>Tags</u>	N-terminal 6His
<u>Native sequence</u>	G37 of the recombinant protein is equivalent to G2 of human MSK2
<u>Accession number</u>	GenBank AJ010119

Recombinant MSSK2 amino acid sequence:

```

1 MSYYHHHHHH DYDIPTTENL YFQGAMGIRN SKAYVDGDED DDESCAVELR ITEANLTGHE
61 EKVSVENFEL LKVLGTGAYG KVFLVRKAGG HDAGKLYAMK VLRKAALVQR AKTQEHTRTE
121 RSVLELVRQA PFLVTLHYAF QTDAKLHLIL DYVSGGEMFT HLYQRQYFKE AEVRVYGGEI
181 VLALEHLHLK GIIYRDLKLE NVLLDSEGHI VLTFDGLSKE FLTEEKERTF SFCGTIEYMA
241 PEIIRSKTGH GKAVDWWSLG ILLFELLTGA SPFTLEGERN TQAEVSRRL KCSPPFPRI
301 GPVAQDLLQR LLCKDPKKRL GAGPQGAQEV RNHPFFQGLD WVALAARKIP APFRPQIRSE
361 LDVGNFAEEF TRLEPVYSPP GSPPPDPRI FQGYSFVAPS ILFDHNNAVM TDGLEAPGAG
421 DRPGRAAVAR SAMMQDSPFF QQYELDLREP ALGQGSFVVC RRCRQRQSGQ EFAVKILSRR
481 LEANTQREVA ALRLCQSHPN VVNLHEVHHD QLHTYLVLEL LRGGELLEHI RKKRHFSESE
541 ASQILRSLVS AVSFMHEEAG VVHRDLKPEN ILYADDTPGA PVKIIDFGFA RLRPQSPGVP
601 MQTPCFTLQY AAPELLAQQG YDESCDLWSL GVILYMMLSG QVPFQGASGQ GGQSQAAEIM
661 CKIREGRFSL DGEAWQGVSE EAKELVRGLL TVDPAKRLKL EGLRGSSWLQ DGSARSSPPL
721 RTPDVLESSG PAVRSGLNAT FMAFNRGKRE GFFLKSVENA PLAKRRKQKL RSATASRRGS
781 PAPANPGRAP VASKGAPRA NGPLPPS

```

Recombinant MSSK2 nucleotide sequence:

```

1 atgtcgtact accatcacca tcaccatcac gattacgata tcccaacgac cgaaaacctg
61 tattttcagg gcgccatggg gatccggaat tcaaaggcct acgtcgacgg ggacgaggac
121 gacgatgaga gctgcgccgt ggagctgcgg atcaccgaag ccaacctgac cgggcacgag
181 gagaaggtga gcgtggagaa cttcgagctg ctcaaggtgc tgggcacggg agcctacggc
241 aaggtgttcc tgggtgaggaa ggcgggcggg cacgacgagg ggaagctgta cgccatgaag
301 gtgctgcgca aggcagcgcct ggtgcagcgc gccaaagacg aagagcacac ggcgaccgag
361 cgctcggtgc tggagctggt gcgccaggcg cccttctctg tcacgctgca ctacgcttcc
421 cagacggatg ccaagctgca cctcatctg gactatgtga gcggcgggga gatgttcacc
481 cacctctacc agcgccagta cttcaaggag gctgaggtgc gcgtgtatgg ggtgtgatgc
541 gtgctgcccc tggaaacct gcacaagctc ggcatacatt accgagacct gaaactggag
601 aatgtgctgc tggactccga gggccacatt gtctctacgg acttcgggct gagcaaggag
661 ttctgacgag aggagaaaga gcggaccttc tccttctgtg gcaccatcga gtacatggcc
721 cccgaaatca tccgtagcaa gacggggcat ggcaaggctg tggactggtg gagcctgggc
781 atcttgcctc tcgagctgct gacgggggccc tcgcccttca ccctggaggg cgagaggaac
841 acgcaggctg aggtgtctcg acggatcctg aagtgtctcc ctcccttccc ccctcggatc
901 gggcccgtgg cgcaggacct gctgcagcgg ctgctttgta aggatcctaa gaagcgattg
961 ggcgcggggc cccagggggc acaagaagtc cggaaccatc ctttcttcca gggcctcgat
1021 tgggtggctc tggctgccag gaagattcca gccccattcc ggccccaaat ccgctcagag
1081 ctggatgtgg gcaactttgc ggaggaattc actcggctgg agcctgtcta ctacccccct
1141 ggcagcccc cacctgggga cccccgaatc tttcagggat actcctttgt ggcaccctcc
1201 attctctttg accacaacaa cgcggtgatg accgatgggc tggaaagcgc tgggtgctgga
1261 gaccggccag gtcgggcagc ggtggccagg agcgctatga tgcaggactc gcccttcttc
1321 cagcagtagc agctggacct gcgggagcct gcgctgggccc agggcagctt ttctgtgtgt
1381 cgccgctgcc gccagcgcca gacgggccag gagttcgcag tcaagatcct cagtcgcagg
1441 ctggaggcga acacgcagcg cgaagtggct gccctgcgcc tgtgccagtc acacccaac
1501 gtggtgaatc tgcacgaggt gcatcacgac cagctgcaca cgtacctggt cctggagctg
1561 ctgcggggcg gggagctgct ggagcacatc cgcaagaagc ggcacttcag cgagtcggaa
1621 gcaagccaga tcctgcgcag cctcgtgtcg gccgtgagct tcatgcacga ggaggcgggc

```

Certificate of Analysis

```
1681 gtggtgcacc gcgacctcaa gccggagaac atcctgtacg ccgacgacac gcccggggcc
1741 ccggtgaaaa tcatcgactt cgggttcgcg cggttgcggc cgcagagtcc cggggtgcc
1801 atgcagacgc cctgcttcac gctgcagtac gctgccccg agctgctggc gcagcagggc
1861 tacgacgagt cctgcgacct ctggagcctg ggcgtcattc tgtacatgat gctgtcgggg
1921 caggtccctt tccagggggc ctctggccag ggcgggcaga gccaggcggc cgagatcatg
1981 tgcaaaatcc gcgagggggc cttctccctt gacggggagg cctggcaggg tgtatccgag
2041 gaagccaagg agctggtccg agggctcctg accgtggacc ccgccaagcg gctgaagctc
2101 gagggactgc ggggcagctc gtggctgcag gacggcagcg cgcgctcctc gcccccgctc
2161 cggacgcccg acgtgctcga gtctctggg cccgcagtgc gctcgggtct caacgccacc
2221 ttcattggcat tcaaccgggg caagcgggag ggcttcttcc tgaagagcgt ggagaatgca
2281 cccctggcca agcggcggaa gcagaagctg cggagcgcca ccgcctccc cgggggctcc
2341 cctgcaccag ccaaccggg ccgagcccc gtcgcctcca aaggggcccc ccgccgagcc
2401 aacggcccc tgccccctc ctaa
```

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.