

Certificate of Analysis

RSK1, active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-509, 14-509-K, 14-509M

Parent Lot # 1765772

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 full-length human RSK1 expressed in Sf21 insect cells. Purified using Ni²⁺/NTA agarose. Activated with PDK1 (cat# 14-280) and MAPK2 (cat# 14-173) and repurified using glutathione-Sepharose, heparin-Sepharose and Ni²⁺/NTA agarose. Purity 94.6% by SDS-PAGE and Coomassie blue staining. MW = 84kDa.

Specific Activity (Parent lot# 1765772): 3145U/mg, where one unit of RSK1 activity is defined as 1nmol phosphate incorporated into 30µM (KKNRRLSVA) per minute at 30°C with a final ATP concentration of 100µM.

Formulation: 1.632mg/ml of enzyme in 50mM Tris/HCl pH7.5, 150mM NaCl pH7.5, 0.03% Brij-35, 0.1mM EGTA, 270mM sucrose, 0.2mM PMSF, 1mM benzamidine, 0.1% 2-mercaptoethanol. Frozen solution.

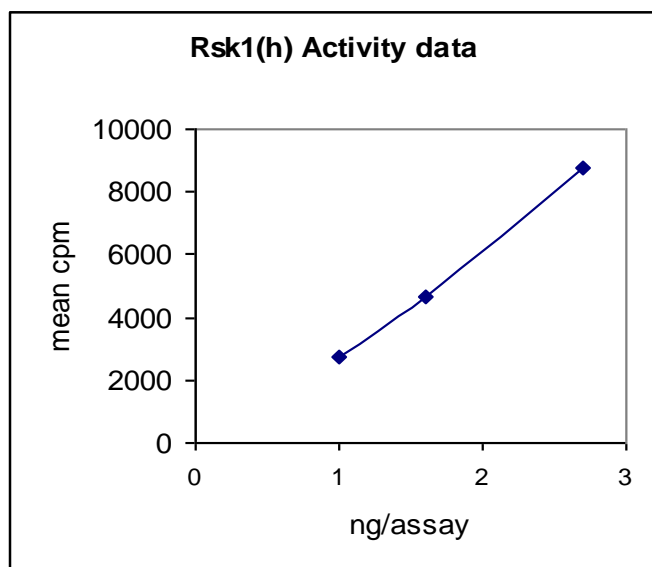
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 6 months from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

Handling Recommendations: Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled microcentrifuge tubes and immediately snap-freeze the vials in liquid nitrogen prior to re-storage at -70°C.

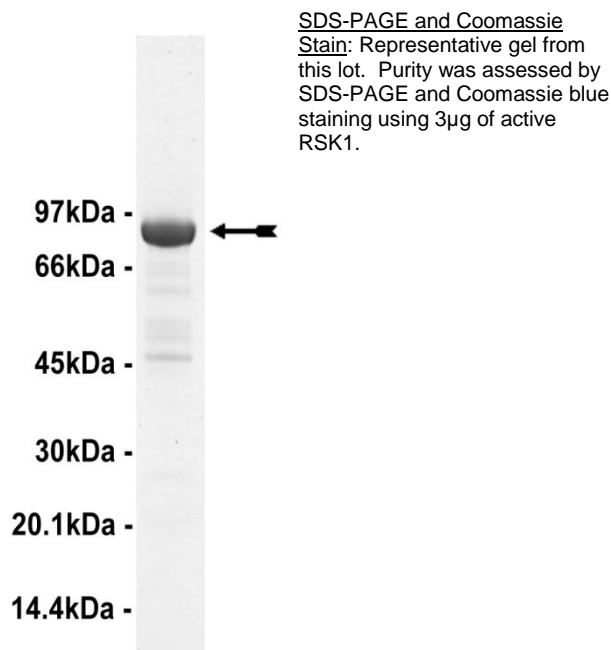
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 1.0–2.7ng of this lot of enzyme phosphorylated 30µM (KKNRRLSVA) 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 RSK1 with the translated sequence listed on page three.



Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

1. **5x Reaction Buffer:** 40mM MOPS/NaOH pH7.0, 1mM EDTA.
2. **(KKNRSLVA):** Use at a final assay concentration of 30µM. Prepare a 300µM stock. Add 2.5µl of stock per assay point.
3. **RSK1, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 1.0–2.7ng 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 **(KKNRSLVA)**.
3. Add **2.5µl (1.0–2.7ng) RSK1, active**.
4. Add 5µl of 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

RSK1 Sequence Information

<u>Protein</u>	human RSK1
<u>Tags</u>	N-terminal 6His
<u>Native sequence</u>	M10 of the recombinant protein is equivalent to M1 of human RSK1
<u>Accession number</u>	GenBank NM_002953

Recombinant RSK1 amino acid sequence:

```

1 MHHHHHHEFM PLAQLKEPWP LMELVPLDPE NGQTSGEEAG LQPSKDEGVL KEISITHHVK
61 AGSEKADPSH FELLKVLGQG SFGKVFLVRK VTRPDSGHLV AMKVLKKATL KVRDRVRTKM
121 ERDILADVNH PFVVKLHYAF QTEGKLYLIL DFLRGGDLFT RLSKEVMFTE EDVKFYLAEL
181 ALGLDHLHSL GIIYRDLKPE NILLEEGHI KLDFGLSKE AIDHEKKAYS FCGTVEYMAP
241 EVVNRQGHSH SADWWSYGVL MFEMLTGSLP FQGKDRKEM TLILKAKLGM PQFLSTEAQS
301 LLRALFKRNP ANRLGSGPDG AEEIKRHVYF STIDWNKLYR REIKPPFKPA VAQPDDTFYF
361 DTEFTSRTPK DSPGIPPSAG AHQLFRGFSF VATGLMEDDG KPRAPQAPLH SVVQQLHGKN
421 LVFSDGYVVK ETIGVGSYSE CKRCVHKATN MEYAVKVIDK SKRDPSEEIE ILLRYGQHPN
481 IITLKDVIYDD GKHVYLVTEL MRGSELLDKI LRQKFFSERE ASFVLHTIGK TVEYLHSQGV
541 VHRDLKPSNI LYVDESGNPE CLRICDFGFA KQLRAENGLL MTPCYTANFV APEVLKRQGY
601 DEGCDIWSLQ ILLYTMLAGY TPFANGPSDT PEEILTRIGS GKFTLSGGNW NTVSETAKDL
661 VSKMLHVDPH QRLTAKQVLQ HPWVTQKDKL PQSQLSHQDL QLVKGAMAAT YSALNSSKPT
721 PQLKPIESSI LAQRRVRKLP STTL

```

Recombinant RSK1 nucleotide sequence:

```

1 atgcatcatc accatcacca tgaattcatg ccgctcgccc agctcaagga gccctggccg
61 ctcatggagc tagtgccgct ggaccgagag aatggacaga cctcagggga agaagctgga
121 cttcagccgt ccaaggatga gggcgctcctc aaggagatct ccatcacgca ccacgtcaag
181 gctggctctg agaaggctga tccatcccat ttcgagctcc tcaaggttct gggccagggga
241 tccttgggca aagtcttcct ggtgcggaaa gtcaccgagc ctgacagtgg gcacctgat
301 gctatgaagg tgctgaagaa ggcaacgctg aaagtacgtg accgctccg gaccaagatg
361 gagagagaca tcctggctga tgtaaatcac ccattcgtgg tgaagctgca ctatgccttc
421 cagaccgagg gcaagctcta tctcattctg gacttctctg gtgggtggga cctcttcacc
481 cggctctcaa aagaggtgat gttcacggag gaggatgtga agttttacct ggccgagctg
541 gctctgggcc tggatcacct gcacagcctg ggtatcattt acagagacct caagcctgag
601 aacatccttc tggatgagga gggccacatc aaactcactg actttggcct gagcaaagag
661 gccattgacc acgagaagaa ggctattctt ttctgcgagg cagtggagta catggccctt
721 gaggtcgtca accgccaggg ccactcccat agtgcgagc ggtggtccta tggggtgttg
781 atgtttgaga tgctgacggg ctccctgccc ttccagggga aggaccgga ggagaccatg
841 aactgattc tgaaggcga gctaggcatg ccccagtttc tgagcactga agcccagagc
901 ctcttgccgg cctgttcaa gcggaatcct gcccaaccggc tcggtccgg ccctgatggg
961 gcagaggaaa tcaagcggca tgtcttctac tccaccattg actggaataa gctataccgt
1021 cgtgagatca agccaccctt caagccagca gtggctcagc ctgatgacac cttctacttt
1081 gacaccgagt tcacgtcccg cacaccaag gattccccag gcatcccccc cagcgtggg
1141 gcccatcagc tgttccgggg cttcagcttc gtggccaccg gcctgatgga agacgacggc
1201 aagcctcgtg ccccgaggc acccctgcac tcggtggtac agcaactcca tgggaagaac
1261 ctggttttta gtgacggcta cgtggtaaag gagacaattg gtgtgggctc ctactctgag
1321 tgcaagcgtg gtgtccacaa ggccaccaac atggagtatg ctgtcaaggt cattgataag
1381 agcaagcggg atccttcaga agagattgag attcttctgc ggtatggcca gcacccaac
1441 atcatcactc tgaagatgt gtatgatgat ggcaaacacg tgtacctggt gacagagctg
1501 atgcgggggt gggagctgct ggacaagatc ctgcggcaga agttcttctc agagcgggag
1561 gccagctttg tcctgcacac cattggcaaa actgtggagt atctgcactc acagggggtt

```

Certificate of Analysis

```
1621 gtgcacaggg acctgaagcc cagcaacatc ctgtatgtgg acgagtccgg gaatcccgag
1681 tgcctgcgca tctgtgactt tggttttgcc aaacagctgc gggctgagaa tgggctcctc
1741 atgacacctt gctacacagc caactttgtg gcgcctgagg tgctgaagcg ccagggctac
1801 gatgaaggct gcgacatctg gagcctgggc attctgctgt acaccatgct ggcaggatat
1861 actccatttg ccaacgggcc cagtgcacac ccagaggaaa tcctaaccg gatcggcagt
1921 gggaaagtta ccctcagtgg gggaaattgg aacacagttt cagagacagc caaggacctg
1981 gtgtccaaga tgctacacgt ggatccccac cagcgcctca cagctaagca ggttctgcag
2041 catccatggg tcaccagaa agacaagctt ccccaaagcc agctgtccca ccaggaccta
2101 cagcttgtga agggagccat ggctgccacg tactccgcac tcaacagctc caagcccacc
2161 cccagctga agccatcga gtcatccatc ctggcccagc ggcgagtgag gaagttgcca
2221 tccaccacc tgtga
```

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.