

Certificate of Analysis

ROCK-I, active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-601, 14-601-K, 14-601M

Parent Lot # D8NN032U

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 ROCK-I amino acids 17–535, expressed by baculovirus in Sf21 insect cells. Purified using Ni²⁺/NTA-agarose. Autoactivated by incubating with Mg/ATP and redialysed to remove excess ATP. Purity 71.1% by SDS-PAGE and Coomassie blue staining. MW = 61.4kDa.

Specific Activity (Parent lot# D8NN032U): 63U/mg, where one unit of ROCK-I activity is defined as 1nmol phosphate incorporated into 30µM Long S6 peptide per minute at 30°C with a final ATP concentration of 100µM.

Formulation: 0.367mg/ml of enzyme in 50mM Tris/HCl pH7.5, 300mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 0.2mM PMSF, 1mM benzamidine, 0.1% 2-mercaptoethanol. Frozen solution.

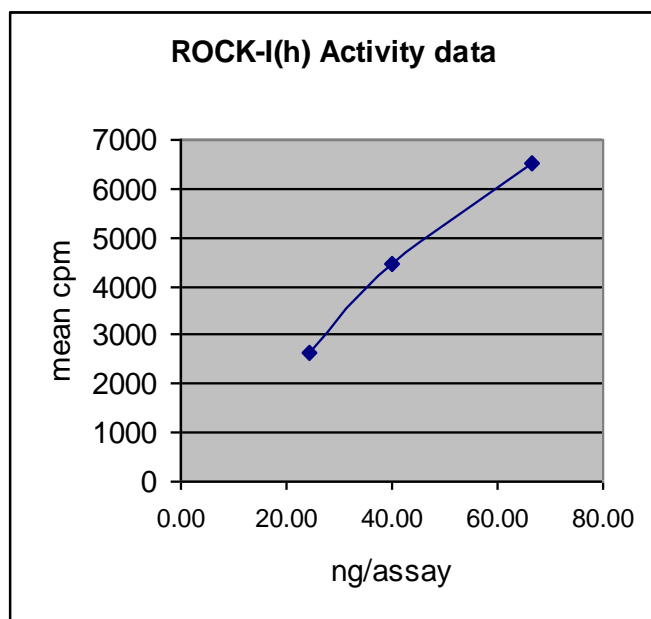
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 6 months from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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.

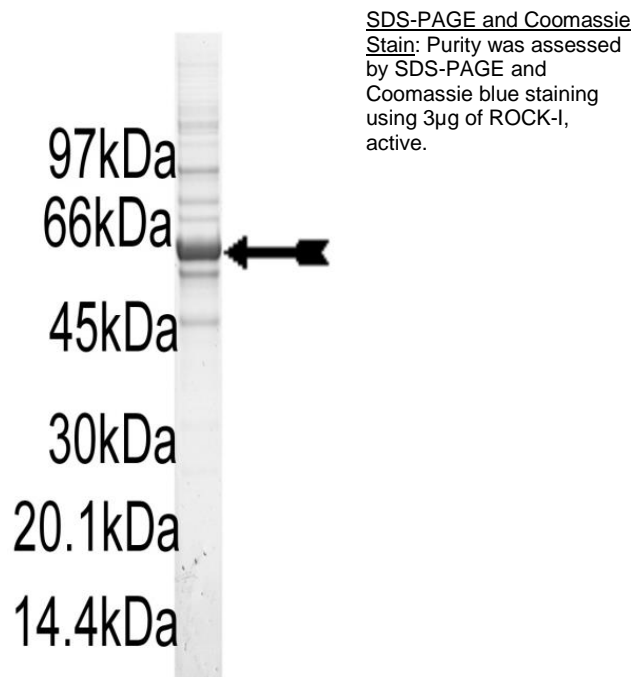
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 24.2–66.6ng of this lot of enzyme phosphorylated 30µM Long S6 peptide 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 ROCK-I with the translated native sequence listed on page three.



Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

1. **5 x Reaction Buffer:** 40mM MOPS/NaOH pH7.0, 1mM EDTA.
2. **Long S6 substrate peptide:** Use at a final assay concentration of 30 μ M. Make up a 300 μ M stock. Add 2.5 μ l of stock per assay point.
3. **ROCK-I, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 24.2–66.6ng 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 5 μ l of 5 x reaction buffer per assay to wells.
2. Add 2.5 μ l of **Long S6 substrate peptide**.
3. Add **2.5 μ l (24.2–66.6ng) ROCK-I, active**.
4. Add 5 μ l of dH₂O.
5. Add 10 μ l of diluted [γ -³³P]ATP mixture
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5 μ l of 3% phosphoric acid.
8. Transfer a 10 μ l 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 4 ml 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

ROCK-I Sequence Information

<u>Protein</u>	human ROCK-I
<u>Tags</u>	N-terminal 6His
<u>Native sequence</u>	L8 of the recombinant protein is equivalent to L17 of human ROCK-I
<u>Accession number</u>	GenBank NM_005406

Recombinant ROCK-I amino acid sequence:

```

1 MHHHHHLLR DPKSEVNSDC LLDGLDALVY DLDFPALRKN KNIDNFLSRY KDTINKIRDL
61 RMKAEDYEVV KVIGRGAFGE VQLVRHKSTR KVIYAMKLLSK FEMIKRSDSA FFWEERDIMA
121 FANSPWVVL FYAFQDDRYL YMVMEYMPGG DLVNLMSNYD VPEKWARFYT AEVVLALDAI
181 HSMGFIHRDV KPDNMLLDKS GHLKLADFGT CMKMNKEGMV RCDTAVGTPD YISPEVLKSQ
241 GGDGYYGREC DWWSVGVFLY EMLVGDTPFY ADSLVGTYSK IMNHKNSLTF PDDNDISKEA
301 KNLICAFSLD REVRLGRNGV EEIKRHLFFK NDQWAWETLR DTVAPVVPDL SSDIDTSNFD
361 DLEEDKGEEE TFPIPKAFVG NQLPFVGFY YSNRRYLSSA NPNDNRTSSN ADKSLQESLQ
421 KTIYKLEEQL HNEMQLKDEM EQKCRTSNIK LDKIMKELDE EGNQRRNLES TVSQIEKEKM
481 LLQHRINEYQ RKAEQENEKR RNVENEVSTL KDQLEDLKKV SQNSQL
  
```

Recombinant ROCK-I nucleotide sequence:

```

1 atgcatcacc atcaccatca tctgctgctg gatcccaaat cggaagtgaa ttcggattgt
61 ttgctggatg gattggatgc tttggatatat gatttggatt ttcctgcctt aagaaaaaac
121 aaaaatattg acaacttttt aagcagatat aaagacacaa taaataaaat cagagattta
181 cgaatgaaag ctgaagatta tgaagtagtg aaggtgattg gtagagggtgc atttgagaa
241 gttcaattgg taaggcataa atccaccagg aaggtatatg ctatgaagct tctcagcaaa
301 tttgaaatga taaagagatc tgattctgct tttttctggg aagaaagga catcatggct
361 tttgccaaca gtccttgggt tgttcagctt ttttatgcat tccaagatga tcggtatctc
421 tacatggtga tggatacat gcctgggtga gatcttgtaa acttaatgag caactatgat
481 gtgcctgaaa aatgggcacg attctatact gcagaagtag ttcttgcatt ggatgcaatc
541 cattccatgg gttttattca cagagatgtg aagcctgata acatgctgct ggataaatct
601 ggacatttga agttagcaga ttttgggtact tgtatgaaga tgaataagga aggcattgta
661 cgatgtgata cagcggttgg aacacctgat tatatttccc ctgaagtatt aaaatcccaa
721 ggtggtgatg gttattatgg aagagaatgt gactggtggg cggttggggg atttttatac
781 gaaatgcttg taggtgatac acctttttat gcagattctt tggttggaac ttacagtaaa
841 attatgaacc ataaaaattc acttaccttt cctgatgata atgacatata aaaagaagca
901 aaaaacctta tttgtgcctt cttactgac agggaagtga ggttagggcg aaatggtgta
961 gaagaaatca aacgacatct cttcttcaaa aatgaccagt gggcttggga aacgctccga
1021 gacactgtag caccagtgtg acccgattta agtagtgaca ttgatactag taattttgat
1081 gacttggaa gataaaagg agaggaagaa acattcccta ttcctaaagc tttcgttggc
1141 aatcaactac cttttgtagg atttacatat tatagcaatc gtagatactt atcttcagca
1201 aatcctaata ataacagaac tagctccaat gcagataaaa gcttgcagga aagtttgcaa
1261 aaaacaatct ataagctgga agaacagctg cataatgaaa tgcagttaaa agatgaaatg
1321 gagcagaagt gcagaacctc aaacataaaa ctagacaaga taatgaaaga attggatgaa
1381 gagggaaatc aaagaagaaa tctagaatct acagtgtctc agattgagaa ggagaaaatg
1441 ttgctacagc atagaattaa tgagtaccaa agaaaagctg aacaggaaaa tgagaagaga
1501 agaaatgtag aaaatgaagt ttctacatta aaggatcagt tggaagactt aaagaaagt
1561 agtcagaatt cacagctttaa
  
```

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.