

Certificate of Analysis

ACK1, active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-756, 14-756-K, 14-756M

Parent Lot # D7EN072U

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 GST-tagged recombinant, human ACK1, residues 1–389. Expressed by baculovirus in Sf21 insect cells. Purified using glutathione agarose. Purity 65% by SDS-PAGE and Coomassie blue staining. MW = 71.5kDa.

Specific Activity (Parent lot# D7EN072U): 249U/mg, where one unit of ACK1, active activity is defined as 1nmol phosphate incorporated into 400 μ M (EFPIYDFLPAKKK) per minute at 30°C with a final ATP concentration of 100 μ M.

Formulation: 0.822mg/ml of enzyme in 50mM Tris/HCl pH7.5, 150mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 0.1mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution.

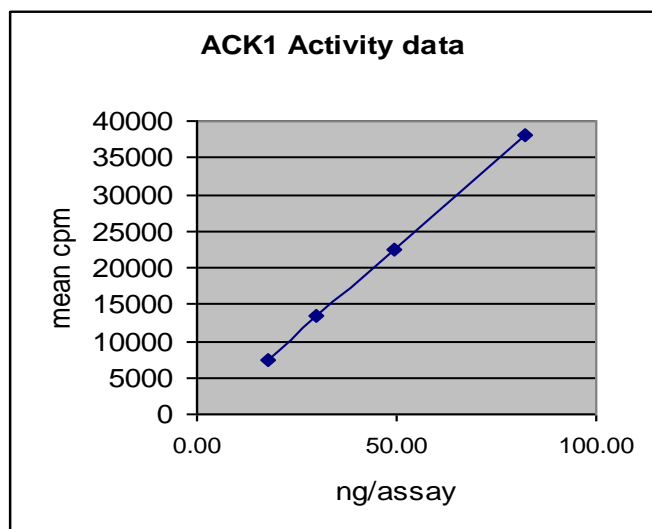
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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.

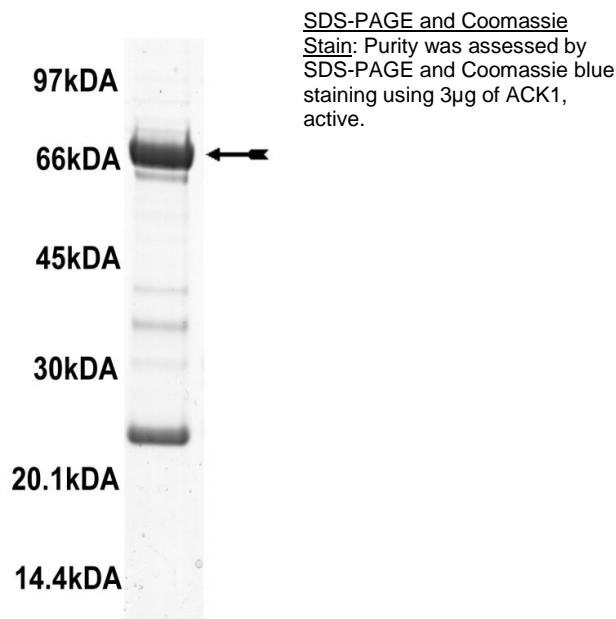
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 18–82ng of this lot of enzyme phosphorylated 400 μ M (EFPIYDFLPAKKK) 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 ACK1 with the translated 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. **(EFPIYDFLPAKKK):** Use at a final assay concentration of 400 μ M. Prepare a 4mM stock and add 2.5 μ l of stock per assay point.
3. **ACK1, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 5% glycerol, 0.01% Brij-35, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 18–82ng 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 **(EFPIYDFLPAKKK)**.
3. Add **2.5 μ l (18–82ng) ACK1, 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 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 μ l of 30% phosphoric acid.

Certificate of Analysis

ACK1 Sequence Information

<u>Protein</u>	human ACK1
<u>Tags</u>	N-terminal GST
<u>Native sequence</u>	M231 of recombinant sequence is equivalent to M1 of the native ACK1
<u>Accession number</u>	GenBank NM_005781

Recombinant ACK1 amino acid sequence:

```

1 MSPILGYWKI KGLVQPTRL L EYLEEKYEE HLYERDEGDK WRNKKFELGL EFPNLPYYID
61 GDVKLTQ SMA IIRYIADKHN MLGGCPKERA EISMLEGAVL DIRYGVSRIA YSKDFETLKV
121 DFLSKLPEML KMFEDRLCHK TYLNGDHVTH PDFMLYDALD VVLYMDPMCL DAFPKLVCFK
181 KRIEAIPQID KYLKSSKYIA WPLQGWQATF GGGDHPPKSD LEVLFQGP EF MQPEEGTGWL
241 LELLSEVQLQ QYFLRLRDDL NVTRLSHF EY VKNEDLEKIG MGRPGQRR LW EAVKRRKALC
301 KRKSWMSKVF SGKRLEAEFP PHHSQSTFRK TSPAPGGPAG EGPLQSLTCL IGEKDLR LLE
361 KLGDSFGV V RRGWDAPSG KTVSVAVKCL KPDVLSQPEA MDDFIREVNA MHSLDHRNLI
421 RLYGVVLT PP MKMVT ELAPL GSLLDRLRKH QGHFLLGTLS RYAVQVAEGM GYLESKRFIH
481 RDLAARNLL L ATRDLVKIGD FGLMRALPQN DDHYVMQEHR KVPFAWCAPE SLKTRTFSHA
541 SDTWMFGV TL WEMFTYGQEP WIGLNGSQIL HKIDKEGERL PRPEDCPQDI YNVMVQCWAH
601 KPEDRPTFVA LRDFLLEAQ
    
```

Recombinant ACK1 nucleotide sequence:

```

1 atgtccccta tactaggtta ttggaaaatt aagggccttg tgcaaccac tcgacttctt
61 ttggaatatac ttgaagaaaa atatgaagag catttgtatg agcgcgatga aggtgataaa
121 tggcgaaca aaaagtttga attgggtttg gagtttcca atcttcctta ttatattgat
181 ggtgatgta aattaacaca gtctatggcc atcatacgtt atatagctga caagcacaac
241 atgttgggtg gttgtccaaa agagcgtgca gagatttcaa tgcttgaagg agcggttttg
301 gatattagat acggtgtttc gagaattgca tatagtaaag actttgaaac tctcaaagtt
361 gattttctta gcaagctacc tgaatgctg aaaatgctc aagatcgtt atgtcataaa
421 acatatttaa atggtgatca tgaaccat cctgacttca tgttgtatga cgctcttgat
481 gttgttttat acatggacc aatgtgcctg gatgcgttcc caaaattagt ttgttttaa
541 aaacgtattg aagctatccc acaaattgat aagtacttga aatccagcaa gtatataaga
601 tggcctttgc agggctggca agccacgttt ggtgggtggcg accatcctcc aaaatcggat
661 ctggaagtctc tgttccagg gcccgaattc atgcagccag aggaggcac agcctggctg
721 ctggagctgc tgcctgaggt gcagctgcaa cagtacttcc tgcggctcc agacgacctc
781 aacgtcacc cctgtccca ctttgagtac gtcaagaatg aggacctgga gaagatcggc
841 atgggtcggc ctggccagcg gcggctgtgg gaggctgtga agaggaggaa ggccttgtgc
901 aaacgcaagt cgtggatgag taaggtgtc agtggaaagc gactggaggc tgagttccca
961 cctcatcact ctcaagcac cttccggaag acctcgccg cccctggggg cccagcaggg
1021 gaggggcccc tgcagagcct cacctgcctc attggggaga aggacctgcg cctcctggag
1081 aagctgggtg atggttcctt tggcgtgggt cgcaggggagc agtgggagc gccctcaggg
1141 aagacggtga gtgtggctgt gaagtgcctg aagcccgatg tcctgagcca gccagaagcc
1201 atggagcact tcatccggga ggtcaatgcc atgcactcgc tcgaccaccg aaacctcatc
1261 cgcctctacg ggggtgtgct cacgccccc atgaagatgg tgacagagct ggcacctctg
1321 ggatcgttgt tggaccggct acgtaagcac cagggccact tcctcctggg gactctgagc
1381 cgctacgctg tgcaggtggc tgagggcagtg ggctacctgg agtccaagcg ctttattcac
1441 cgtgacctgg ctgccgcaa tctgctgttg gctaccgcg acctggtcaa gatcggggac
1501 tttgggctga tgcgagcact acctcagaat gacgaccatt acgtcatgca ggaacatcgc
1561 aaggtgccct tgcctgggtg tcccccgag agcctgaaga cacgtacctt ctcccatgcc
1621 agcgacacct ggatgttcgg ggtgacctg tgggaaatgt tcacctacgg ccaggagccc
1681 tggatcggcc tcaacggcag tcagatcctg cataagatcg acaaggaggg ggagcggctg
    
```

Certificate of Analysis

1741 ccccgcccg aggactgtcc ccaggacatc tacaacgtca tgggccagtg ctgggctcac
1801 aagccagagg acagaccac gtttgtggcc ctgcgggact tcctgctgga ggcccagtaa

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