

Certificate of Analysis

CRIK, active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 16-040, 16-040-K, 16-040M

Parent Lot # D17HP011N

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 CRIK amino acids 1-449 expressed by baculovirus in Sf21 insect cells. Purified using glutathione agarose.

Purity 88% by SDS-PAGE and Coomassie blue staining. MW = 77kDa.

Specific Activity (Parent lot# D17HP011N): 96U/mg, where one unit of CRIK activity is defined as 1nmol phosphate incorporated into 150 μ M KKLRRRLSVA per minute at 30°C with a final ATP concentration of 100 μ M.

Formulation: 1.24mg/ml of enzyme in 50mM Tris/HCl pH7.5, 300mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 1mM benzimidazole, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution.

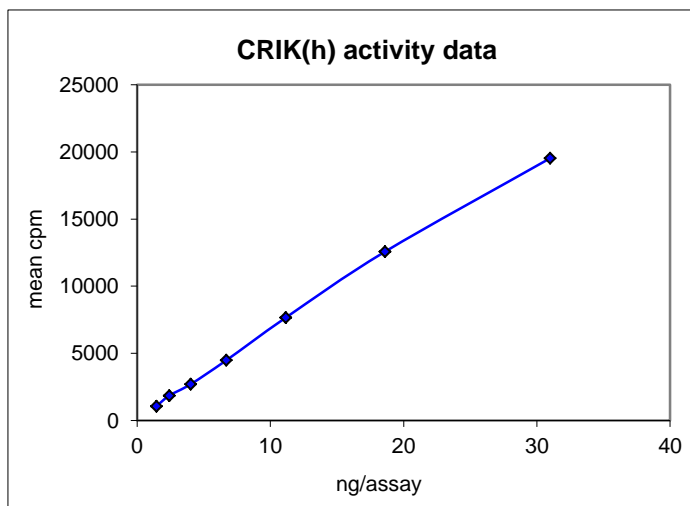
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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.

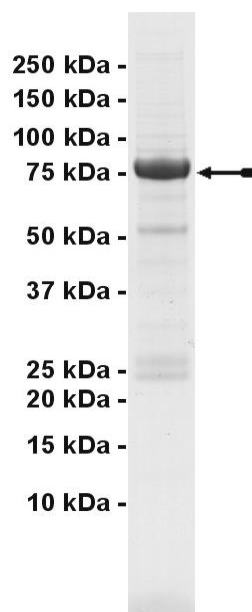
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 1–31ng of this lot of enzyme phosphorylated 150 μ M KKLRRRLSVA 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 CRIK with the translated sequence listed on page three.



SDS-PAGE and Coomassie Stain: Purity was assessed by SDS-PAGE and Coomassie blue staining using 3 μ g of CRIK, active..

Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

- 1. 5 x Reaction Buffer:** 40mM MOPS/NaOH pH7.0, 1mM EDTA.
- 2. KCLRRTLSVA:** Use at a final assay concentration of 150 μ M. Prepare a 1.5mM stock and add 2.5 μ l of stock per assay point.
- 3. CRIK, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 1–31ng per assay point.
- 4. [γ -³³P]ATP:** 2.5 x MgAc/[γ -³³P]ATP cocktail: 25mM MgAc and 0.25mM ATP to which is added [γ -³³P]ATP (specific activity approximately 1500 - 2400cpm/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 KCLRRTLSVA.
3. Add **2.5 μ l (1–31ng) CRIK, active.**
4. Add 5 μ l of dH₂O.
5. Add 10 μ l of diluted [γ -³³P]ATP mixture.
6. Incubate for 30 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

CRIK, active Sequence Information

<u>Protein</u>	Human CRIK
<u>Tags</u>	N-terminal GST
<u>Native sequence</u>	M230 of the recombinant protein is equivalent to M1 of human CRIK
<u>Accession number</u>	GenBank NM_007174.2

Recombinant CRIK amino acid sequence:

```

1 MSPILGYWKI KGLVQPTRLL LEYLEEKYEE HLYERDEGDK WRNKKFELGL EFPNLPYYID
61 GDVKLTQSMA IIRYIADKHN MLGGCPKERA EISMLEGAVL DIRYGVSRIA YSKDFETLKV
121 DFLSKLPPEML KMFEDRLCHK TYLNGDHVTH PDFMLYDALD VVLYMDPMCL DAFPKLVCFK
181 KRIEAIPOID KYLKSSKYIA WPLQGWQATF GGGDHPPKSD LVPRGSKELM LKFKYGARNP
241 LDAGAAEPIA SRASRLNLFQ QGKPPFMTQQ QMSPLSREGI LDALFVLFEE CSQPALMKIK
301 HVSNFVRKYS DTIAELQELQ PSAKDFEVRS LVGCGHFAEV QVVREKATGD IYAMKVMKKK
361 ALLAQEQVSF FEEERNILSR STSPWIPQLQ YAFQDKNHLY LVMEYQPGGD LLSLLNRYED
421 QLDENLIQFY LAELILAVHS VHLMGYVHRD IKPENILVDR TGHIKLVDFG SAAKMNSNKM
481 VNAKLPIGTP DYMAPEVLTV MNGDGKGTYG LDCDWWSVGV IAYEMIYGRS PFAEGTSART
541 FNNIMNFQRF LKFPDDPKVS SDFLDLIQSL LCGQKERLKF EGLCCHPFFS KIDWNNIRNS
601 PPPFVPTLKS DDDTSNFDEP EKNSWVSSSP CQLSPSGFSG EELPFVGFYSY SKALGILGRS
661 ESVVSGLDSP AKTSSMEK
    
```

Recombinant CRIK nucleotide sequence:

```

1 atgtccccta tactaggtta ttggaaaatt aagggccttg tgcaaccacac tcgaacttctt
61 ttggaatattc ttgaagaaaa atatgaagag catttgatag agcgcgatga aggtgataaaa
121 tggcgaaaca aaaagtttga attgggtttg gagtttccca atcttccctta ttatattgat
181 ggtgatgta aattaacaca gtctatggcc atcatacgtt atatagctga caagcacaac
241 atgttgggtg gttgtccaaa agagcgtgca gagatttcaa tgcttgaagg agcgggttttg
301 gatattagat acggtgtttc gagaattgca tatagtaaag actttgaaac tctcaaagtt
361 gattttctta gcaagctacc tgaaatgctg aaaatgttcg aagatcgttt atgtcataaaa
421 acatatttaa atggtgatca tgtaaccat cctgacttca tgttgatga cgctcttgat
481 gttgttttat acatggacc aatgtgcctg gatgcttcc caaaattagt ttgttttaaa
541 aaacgtattg aagctatccc acaattgat aagtacttga aatccagcaa gtatatagca
601 tggcctttgc agggctggca agccacgttt ggtggtggcg accatcctcc aaaatcggat
661 ctggttccgc gtggatcaa ggaattgat ttgaagttca aatatggagc gcggaatcct
721 ttggatgctg gtgctgctga acccattgcc agccgggcct ccaggctgaa tctgttcttc
781 caggggaaac caccctttat gactcaacag cagatgtctc ctcttcccag agaagggata
841 ttagatgccc tctttgttct ctttgaagaa tgcagtcagc ctgctctgat gaagattaag
901 cacgtgagca actttgtccg gaagtattcc gacaccatag ctgagttaca ggagctccag
961 cttcggcaa aggacttcca agtcagaagt cttgtagggt gtggtcactt tgctgaagtg
1021 caggtggtaa gagagaaagc aaccggggac atctatgcta tgaaagtgat gaagaagaag
1081 gctttatttg cccaggagca ggtttcattt tttgaggaag agcgggaacat attatctcga
1141 agcacaagcc cgtggatccc ccaattacag tatgcctttc aggacaaaaa tcacctttat
1201 ctggtcatgg aatatcagcc tggaggggac ttgctgtcac ttttgaatag atatgaggac
1261 cagttatagat aaaacctgat acagttttac atagctgagc tgattttggc tgttcacagc
1321 gttcatctga ttggatacgt gcatcgagac atcaagcctg agaacattct cgttgaccgc
1381 acaggacaca tcaagctggt ggattttggga tctgccgcga aatgaattc aaacaagatg
1441 gtgaatgcc aactcccgat tgggacccca gattacatgg ctctgaagt gctgactgtg
1501 atgaacgggg atggaaaagg cacctacggc ctggactgtg actggtggtc agtggggctg
1561 attgcctatg agatgattta tgggagatcc cccttcgcag agggaaacctc tgccagaacc
    
```

Certificate of Analysis

```
1621 ttcaataaca ttatgaatth ccagcggtht ttgaaatthc cagatgacc caaagtgagc
1681 agtgactthc ttgatctgat tcaaagcttg ttgtgcggcc agaaagagag actgaagtht
1741 gaaggtctth gctgccatcc thtctthctc aaaattgact ggaacaacat tcgtaactct
1801 cctccccctc tcgttcccac cctcaagtct gacgatgaca cctccaatth tgatgaacca
1861 gagaagaatt cgtgggthtc atcctctccg tgccagctga gccctcagg cttctcgggt
1921 gaagaactgc cgtthtgthg gththctgtac agcaaggcac tggggattct tggtagatct
1981 gagtctgthg thtccggthc ggactcccct gccaaagacta gctccatgga aaagtaa
```

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.