

Certificate of Analysis

PEK, active

(Recombinant enzyme expressed in Sf21 cells)

Item # 14-916, 14-916-K, 14-916M

Parent Lot # 2042626

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 PEK amino acids 536 – end, expressed by baculovirus in Sf21 insect cells. Purified using immobilized metal affinity chromatography. Purity 81.7% by SDS-PAGE and Coomassie blue staining. MW = 71kDa.

Specific Activity (Parent lot# 2042626): 383U/mg, where one unit of PEK, active activity is defined as 1nmol phosphate incorporated into 300µM RSRSRSRSRSRSRSR per minute at 30°C with a final ATP concentration of 100µM.

Formulation: 0.82mg/ml of enzyme in of 50mM Tris/HCl pH7.5, 300mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 1mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution

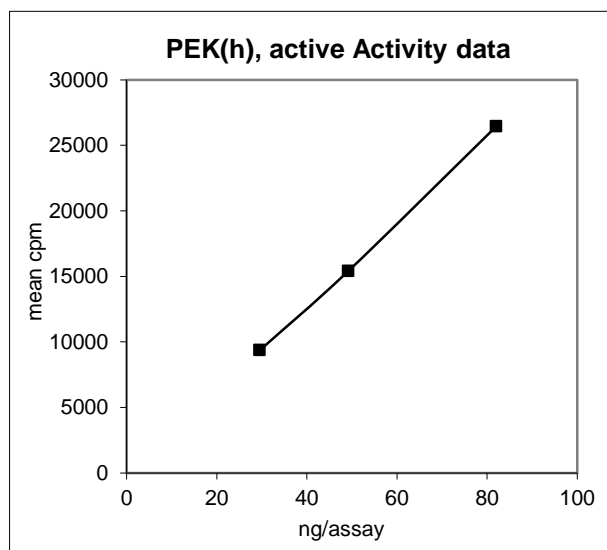
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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.

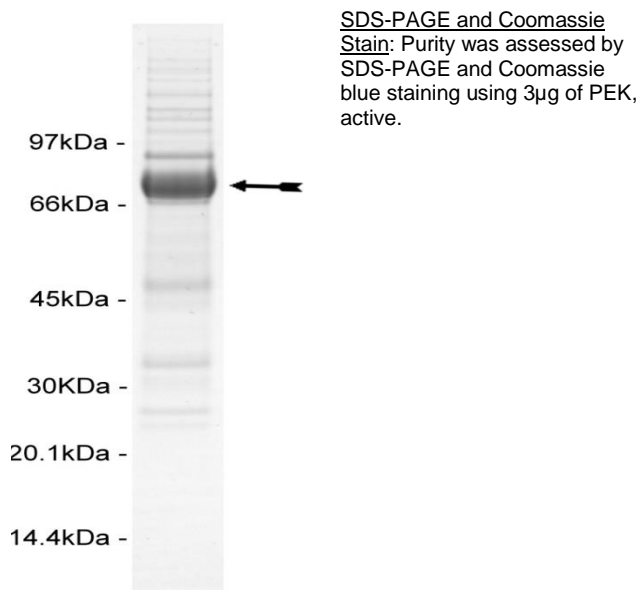
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 30–82ng of this lot of enzyme phosphorylated 300µM RSRSRSRSRSRSRSR 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 PEK 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, other components.
2. **RSRSRSRSRSRSRSR:** Use at a final assay concentration of 300 μ M. Prepare a 3mM stock and add 2.5 μ l of stock per assay point.
3. **PEK(h), active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 30–82ng 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 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 **RSRSRSRSRSRSRSR**.
3. **Add 2.5 μ l (30–82ng) PEK(h), 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

PEK Sequence Information

Protein	Human PEK
Tags	N-terminal 6His
Native sequence	T37 of the recombinant protein is equivalent to T536 of human PEK
Accession number	GenBank NM_004836

Recombinant PEK amino acid sequence:

```

1 MSYYHHHHHH DYDIPTTENL YFQGAMDPEF KGLRRQTTFI VRRLFHPPHP RQRKESETQC
61 QTENKYDSVS GEANDSSWND IKNSGYISRY LTDFEPIQCL GRGGFGVVFE AKNKVDDCNY
121 AIKRIRLPNR ELAREKVMRE VKALAKLEHP GIVRYFNAWL EAPPEKWQEK MDEIWLKDES
181 TDWPLSSPSP MDAPSVKIRR MDPFATKEHI EIIAPSPQRS RSFVSGISCD QTSSSESQFS
241 PLEFSGMDHE DISESVDAAY NLQDSCLTDC DVEDGTMDGN DEGHSFELCP SEASPYVRSR
301 ERTSSSIVFE DSGCDNASSK EEPKTNRLHI GNHCANKLTA FKPTSSKSSS EATLSISPPR
361 PTTLSLDLTK NTTEKLQPSK PKVYLYIQMQ LCRKENLKDW MNGRCTIEER ERSVCLHIFL
421 QIAEAVEFLH SKGLMHRDLK PSNIFFTMDD VVKVGDFGLV TAMDQDEEEQ TVLTPMPAYA
481 RHTGQVGTKL YMSPEQIHGN SYSHKVDIFS LGLILFELLY PFSTQMERVER TLT DVRNLKF
541 PPLFTQKYPC EYVMVQDMLS PSPMERPEAI NIIENAVFED LDFPGKTVLR QRSRSLSSSG
601 TKHSRQSNNS HSPLPSN
  
```

Recombinant PEK nucleotide sequence:

```

1 atgtcgtact accatcacca tcaccatcac gattacgata tcccaacgac cgaaaacctg
61 tattttcagg ggcgatgga tccggaattc aaaggcctac gtcgacaaac aacgtttatt
121 gtgcgagcgc ttttccatcc tcacctcac aggcaaagga aggagtctga aactcagtgt
181 caaactgaaa ataaatatga ttctgtaagt ggtgaagcca atgacagtag ctggaatgac
241 ataaaaaact ctggatata atcacgatata ctaactgatt ttgagccaat tcaatgcctg
301 ggacgtgggtg gctttggagt tgtttttgaa gctaaaaaca aagtagatga ctgcaattat
361 gctatcaaga ggatccgtct ccccaatagg gaattggctc gggaaaagg aatgcgagaa
421 gttaaagcct tagccaagct tgaacacccg ggcattgta gatatttcaa tgctggctc
481 gaagcaccac cagagaagtg gcaagaaaag atggatgaaa ttggctgaa agatgaaagc
541 acagactggc cactcagctc tcctagccca atggatgcac catcagttaa aatacgcaga
601 atggatcctt tcgctacaaa agaactatatt gaaatcatag ctccttcacc acaagaagc
661 aggtcttttt cagtaggat ttctgtgac cagacaagtt catctgagag ccagttctca
721 ccaactggaat tctcaggaat ggaccatgag gacatcagtg agtcagtgga tgcagcatac
781 aacctccagg acagttgcct tacagactgt gatgtggaag atgggactat ggatggcaat
841 gatgaggggc actcctttga actttgtcct tctgaagctt ctccttatgt aaggtcaagg
901 gagagaacct cctcttcaat agtatttgaa gattctggct gtgataatgc ttccagtaaa
961 gaagagccga aaactaatcg attgcatatt ggcaaccatt gtgctaataa actaactgct
1021 ttcaagccca ccagtagcaa atcttcttct gaagctacat tgtctatttc tcctccaaga
1081 ccaaccactt taagtttaga tctcactaaa aacaccacag aaaaactcca gccagttca
1141 ccaagggtgt atctttacat tcaaatgcag ctgtgcagaa aagaaaacct caaagactgg
1201 atgaatggac gatgtaccat agaggagaga gagaggagcg tgtgtctgca catcttcctg
1261 cagatcgagc aggcagtgga gtttcttcac agtaaaggac tgatgcacag ggacctcaag
1321 ccatccaaca tattctttac aatggatgat gtggtcaagg ttggagactt tgggttagtg
1381 actgcaatgg accaggatga ggaagagcag acggttctga ccccaatgcc agcttatgcc
1441 agacacacag gacaagtagg gaccaaactg tatatgagcc cagagcagat tcatggaaac
1501 agctattctc ataaagtgga catcttttct ttaggcctga ttctatttga attgctgtat
1561 ccattcagca ctcagatgga gagagtcagg accttaactg atgtaagaaa tctcaaattt
1621 ccaccattat ttactcagaa atatccttgt gagtacgtga tggttcaaga catgctctct
1681 ccatccccc a tggaaacgacc tgaagctata aacatcattg aaaatgctgt atttgaggac
  
```

Certificate of Analysis

1741 ttggactttc caggaaaaac agtgctcaga cagaggtctc gctccttgag ttcacgga
1801 aaaaaacatt caagacagtc caacaactcc catagccctt tgccaagcaa ttag

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.