

Certificate of Analysis

PhKy2, active

(Recombinant enzyme expressed in *E. coli* cells)

Item # 14-698, 14-698-K, 14-698M

Parent Lot # D8CN057U

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, PhKy2, amino acids (1–301) containing the mutation F290L, expressed by *E. coli* cells. Purified using glutathione-agarose. Purity 89.1% by SDS-PAGE and Coomassie blue staining. MW = 61.37kDa.

Specific Activity (Parent lot# D8CN057U): 134U/mg, where one unit of PhKy2, active activity is defined as 1nmol phosphate incorporated into 250 μ M (KKLNRTLSFAEPG) per minute at 30°C with a final ATP concentration of 100 μ M.

Formulation: 1.138mg/ml of enzyme in 50mM Tris/HCl pH7.5, 300mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 1mM benzamide, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution.

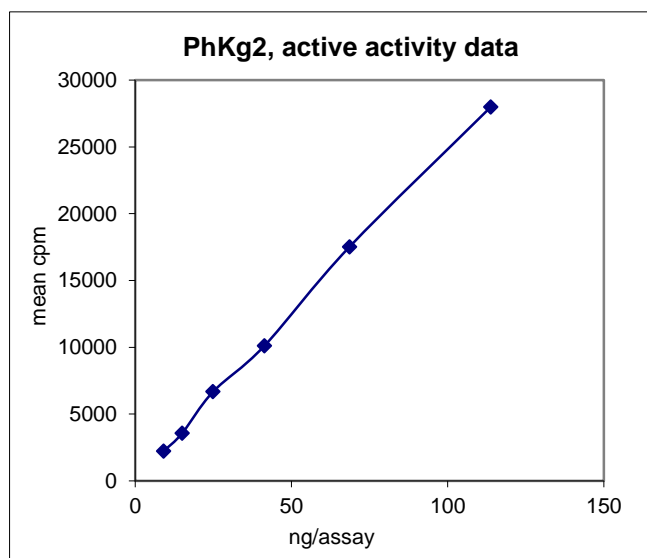
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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.

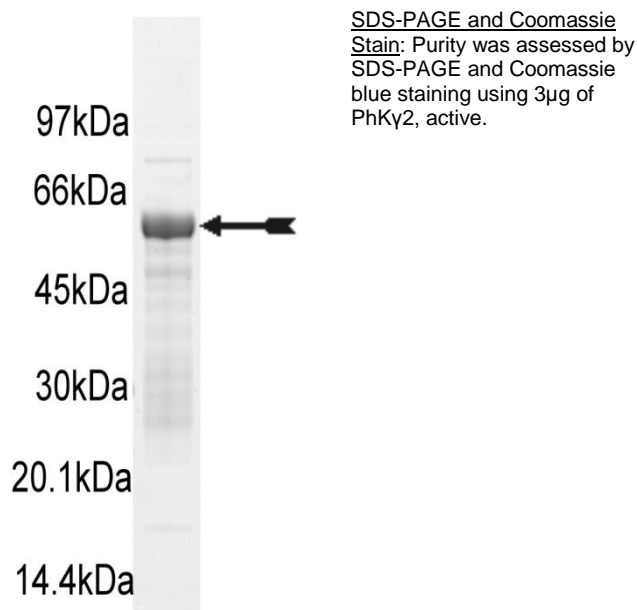
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 9–114ng of this lot of enzyme phosphorylated 250 μ M (KKLNRTLSFAEPG) 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 PhKy2 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. **(KKLNRTL^SFAEPG):** Use at a final concentration of 250µM. Make up a 2.5mM stock. Use 2.5µl of stock per assay point.
3. **PhKγ2, active:** Dilute with 20mM MOPS-NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 9–114ng 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 **(KKLNRTL^SFAEPG)**.
3. Add **2.5µl (9–114ng), PhKγ2 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

PhKy2 Sequence Information

Protein	Human PhKy2
Tags	N-terminal GST
Native sequence	M230 of recombinant sequence is equivalent to M1 of human PhKy2

Accession number GenBank NM_000294. The recombinant protein contains the amino acid substitution F290L with respect to this accession number. Inhibition studies on recombinant PhKy2 F290L exhibit a trend similar to that observed with the wild-type PhK holoenzyme using the inhibitors K252a, Bis-5 and Ro 318220 (Davies *et al.*, (2000), *Biochem. J.* **351**, 95-105; Elliott, *et al.*, (1990), *Biochem. Biophys. Res. Commun.* **171**,148-154)

Recombinant PhKy2 amino acid sequence:

```

1 MSPILGYWKI KGLVQPTRL L LEYLEEKYEE HLYERDEGDK WRNKKFELGL EFPNLPYYID
61 GDVKLTQ SMA IIRYIADKHN MLGGCPKERA EISMLEGAVL DIRYGVSR IA YSKDFETLKV
121 DFLSKLP EML KMFEDRLCHK TYLNGDHVTH PDFMLYDALD VVLYMDPMCL DAFPKLVCFK
181 KRIEAI PQID KYLKSSKYIA WPLQGWQATF GGGDHPPKSD LVPRGSPEFM TLDVGPEDL
241 PDWAAA KEFY QKYDPKDVIG RGVS SVVRR C VHRATGHEFA VKIMEVTAER LSPEQLEEV R
301 EATRRE THIL RQVAGHPHII TLIDSYESSS FMFLVFDLMR KGELFDYLTE KVALSEKET R
361 SIMRSL LEAV SFLHANNIVH RDLKPENILL DDNMQIRLSD FGFSCHLEPG EKLRELCGTP
421 GYLAPE ILC K SMDETHPGYG KEVDLWACGV ILFTLLAGSP PFWHRRQILM LRMIMEGQYQ
481 FSSPEW DDRS STVKDLISRL LQVDPEARLT AEQALQHPLF ERCEGSQPWN
  
```

Recombinant PhKy2 nucleotide sequence:

```

1 atgtccccta tactaggtta ttggaaaatt aagggccttg tgcaaccac tcgacttctt
61 ttggaatata ttgaagaaaa atatgaagag catttgtatg agcgcgatga aggtgataaa
121 tggcga aaca aaaagtttga attgggtttg gagtttccca atcttcctta ttatattgat
181 ggtgatg tta aattaacaca gtctatggcc atcatacgtt atatagctga caagcacaac
241 atgttgg gttg tttccaaa agagcgtgca gagatttcaa tgcttgaagg agcggttttg
301 gatattag at acggtgtttc gagaattgca tatagtaaag actttgaaac tctcaaagtt
361 gattttct ta gcaagctacc tgaatgctg aaaatgttcg aagatcgttt atgtcataaa
421 acatattt aa atggtgatca tgaacctat cctgacttca tgttgtatga cgctcttgat
481 gttgtttt at acatggacc aatgtgcctg gatgcgttcc caaaattagt ttgttttaa
541 aaacgtatt g aagctatccc acaaattgat aagtacttga aatccagcaa gtatatagca
601 tggccttt gc agggctggca agccacgttt ggtggtggcg accatcctcc aaaatcggat
661 ctggttcc gc gtggatcccc ggaattcatg acgctggacg tggggccgga ggatgagctg
721 cccgactgg g cccgcccaa agagttttac cagaagtacg accctaagga cgtcatcggc
781 agaggagt ga gctctgtggt cccgcttgt gttcatcgag ctactggcca cgagtttgcg
841 gtgaagatt a ttggaagtga agctgagcgg ctgagtcctg agcagctgga ggaggtgagg
901 gaagccac ac ggcgagagac acacatcctt cgccaggtcg ccggccacc ccacatcatc
961 accctcat cg attcctacga gtcttctagc ttcattgttcc tgggtgttga cctgatgagg
1021 aagggag agc tgtttgacta tctcacagag aagggtggccc tctctgaaa ggaaaccagg
1081 tccatcat gc ggtctctgct ggaagcagtg agctttctcc atgccaaca cattgtgcat
1141 cgagatct ga agcccagaaa tattctccta gatgacaata tgcagatccg actttcagat
1201 ttcgggtt ct cctgccactt ggaacctggc gagaagcttc gagagttgtg tgggacccca
1261 gggatatc tag cgccagagat ccttaaatgc tccatggatg aaaccacc ccaggctatggc
1321 aaggagg tgc acctctgggc ctgtgggggtg atcttgttca cactcctggc tggctcgcca
1381 cccttctg gc accggcggca gatcctgatg ttacgcatga tcatggaggg ccagtaccag
  
```

Certificate of Analysis

```
1441 ttcagttccc cggagtggga tgaccgttcc agcactgtca aagacctgat ctccaggctg  
1501 ctgcaggtgg atcctgaggc acgcctgaca gctgagcagg ccctacagca cccctctttt  
1561 gagcgttggt aaggcagcca accctggaac taa
```

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