

Certificate of Analysis

PKD2, active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-506, 14-506-K, 14-506M

Parent Lot # 33189U

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, full length, human PKD2, expressed by baculovirus in Sf21 insect cells. Purified using Ni²⁺-NTA agarose. Purity 52.3% by SDS-PAGE and Coomassie blue staining. MW = 100kDa.

Specific Activity (Parent lot# 33189U): 215U/mg, where one unit of PKD2 activity is defined as 1nmol phosphate incorporated into 30µM (KKLNRTLSSVA) per minute at 30°C with a final ATP concentration of 100µM.

Formulation: 1.004mg/ml of enzyme in 50mM Tris/HCl pH7.5, 150mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 1mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution.

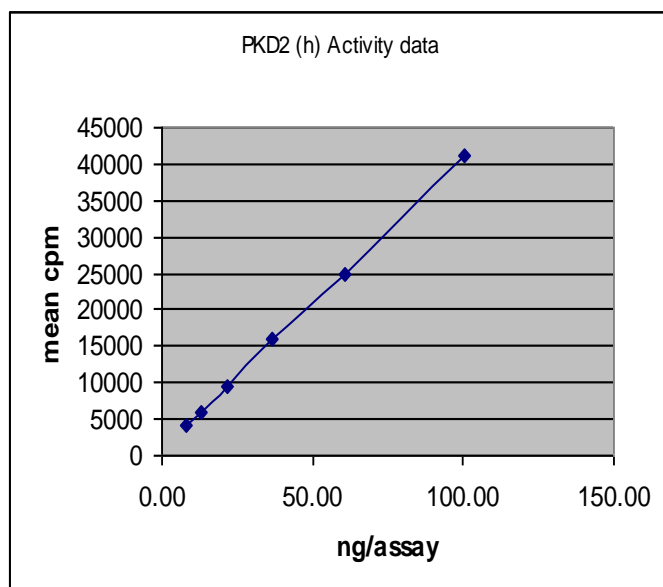
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 6 months from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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.

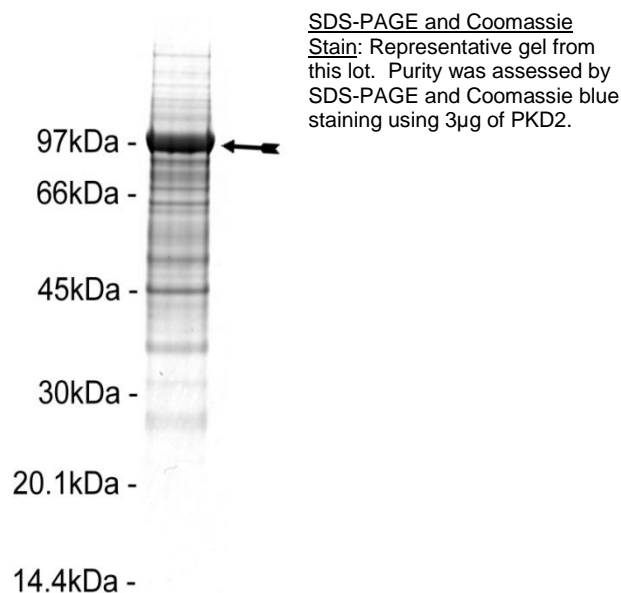
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 8–36ng of this lot of enzyme phosphorylated 30µM (KKLNRTLSSVA) 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 PKD2 with the translated sequence listed on page three.



Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

1. **10 x Reaction Buffer:** 200mM HEPES/NaOH pH7.4, 0.3% Triton X-100.
2. **(KKLNRTLVA):** Use at a final assay concentration of 30 μ M. Prepare a 300 μ M stock. Add 2.5 μ l of stock per assay point.
3. **PKD2, active:** Dilute with 20mM HEPES/NaOH pH7.4, 0.03% Triton X-100. Use 8–36ng 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 2.5 μ l of 10 x reaction buffer per assay to wells.
2. Add 2.5 μ l of **(KKLNRTLVA)**.
3. Add **2.5 μ l (8–36ng) PKD2, active**.
4. Add 7.5 μ l 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

PKD2 Sequence Information

<u>Protein</u>	Human PKD2
<u>Tags</u>	N-terminal 6His
<u>Native sequence</u>	M29 of the recombinant protein is equivalent to M1 of human PKD2
<u>Accession number</u>	GenBank NM_016457

Recombinant PKD2 amino acid sequence:

```

1 MSYYHHHHH DYDIPTTENL YFQGAMGMSMA TAPSYAGLP GSPGPGSPPP PGGLELQSP
61 PLLPQIPAPG SGVSFHIQIG LTREFVLLPA ASELAHVKQL ACSIVDQKFP ECGFYGLYDK
121 ILLFKHDPTS ANLLQLVRSS GDIQEGDLVE VVLSASATFE DFQIRPHALT VHSYRAPAFC
181 DHCGEMLFGL VRQGLKCDGC GLNYHKRCFA SIPNNCSGAR KRRLSSTSLA SGHSVRLGTS
241 ESLPCTAEEL SRSTTELLPR RPPSSSSSSS ASSYTGRPIE LDKMLLSKVK VPHTFLIHSY
301 TRPTVCQACK KLLKGLFRQG LQCKDCKFNC HKRCATRVPN DCLGEALING DVPMEEATDF
361 SEADKSALMD ESEDSGVIPG SHSENALHAS EEEEEGGKA QSSLGYIPLM RVVQSVRHTT
421 RKSSTTLREG WVVHYSNKDT LRKRHYWRLD CKCITLFQNN TTNRYYKEIP LSEILTVESA
481 QNFSLVPPGT NPHCFEIVTA NATYFVGEMP GGTPGGPSGQ GAEAARGWET AIRQALMPVI
541 LQDAPSAPGH APHRQASLSI SVSNSQIQEN VDIATVYQIF PDEVLGSGQF GVVYGGKHRK
601 TGRDVAVKVI DKLRFPKQE SQLRNEVAIL QSLRHPGIVN LECMFETPEK VFVMEKHLG
661 DMLEMILSSE KGRLPERLTK FLITQILVAL RHLHFKNIVH CDLKPENVLL ASADPFQVK
721 LCDFGFARII GEKSFRRSVV GTPAYLAPEV LLNQGYNRS LDMWSVGVIMY VLSLGTFFFN
781 EDEDINDQIQ NAAFMYASP WSHISAGAI D LINNLLQVKM RKRYSDKSL SHPWLQEYQT
841 WLDLRELEGK MGERYITHES DDARWEQFAA EHPLPGSGLP TDRDLGGACP PQDHDMQGLA
901 ERISVL

```

Recombinant PKD2 nucleotide sequence:

```

1 atgtcgtact accatcacca tcaccatcac gattacgata tcccaacgac cgaaaacctg
61 tattttcagg ggccatggg atccatggc accgcccct cttatcccgc cgggctccct
121 ggctctccc ggccggggtc tcctccgcc cccggcggcc tagagctgca gtcgccgcca
181 ccgctactgc ccagatccc ggccccgggt tccggggtct ctttcacat ccagatcggg
241 ctgaccgcg agttcgtgct gttgccgcc gcctccgagc tggctcatgt gaagcagctg
301 gcctgttcca tcgtggacca gaagttccct gagtgtggct tctacggcct ttacgacaag
361 atcctgcttt tcaaacatga ccccacgtcg gccaacctcc tgcagctggt gcgctcgtcc
421 ggagacatcc aggagggcga cctggaggag gtgggtgctgt cggcctcggc cacttcgag
481 gacttcaga tccgccgca cgccctcacg gtgcactcct atcgggcgcc tgccttctgt
541 gatcactgcg gggagatgct cttcggccta gtgcgccagg gcctcaagtg cgatggctgc
601 gggctgaact accacaagcg ctgtgcctt agcatccca acaactgtag tggggcccgc
661 aaacggcgcc tgtcatccac gtctctggcc agtgccact cggctgcct cggcacctcc
721 gaggccctgc cctgcacggc tgaagagctg agccgtagca ccaccgaact cctgcctcgc
781 cgtccccgt catcctctt ctcctcttct gcctcatcgt atacgggccc cccattgag
841 ctggacaaga tgctgctctc caaggtcaag gtgccgcaca cttcctcat ccacagctat
901 acacggccca ccggttgcca ggcttgcaag aaactcctca agggcctctt ccggcagggc
961 ctgcaatgca aagactgcaa gtttaactgt cacaacgct gcgccaccg cgtccctaat
1021 gactgcctgg gggaggccct tatcaatgga gatgtgccga tggaggaggc caccgatttc
1081 agcaggctg acaagagcgc cctcatggat gactcagagg actccggtgt catccctggc
1141 tcccactcag agaatgcgct ccacgccagt gaggaggagg aaggcgaggg aggcaaggcc
1201 cagagctccc tgggtacat ccccctaat aggggtggtc aatcgggtcg acacacgacg
1261 cggaatcca gcaccagct gcgggagggt tgggtggttc attacagcaa caaggacacg
1321 ctgagaaagc ggcactattg gcgcctggac tgcaagtgta tcacgctctt ccagaacaac
1381 acgaccaaca gatactataa ggaaattccg ctgtcagaaa tcctcacggt ggagtccgcc

```

Certificate of Analysis

```
1441 cagaacttca gccttgtgcc gccgggcacc aaccacact gctttgagat cgtcactgcc
1501 aatgccacct acttcgtggg cgagatgcct ggccgggactc cgggtgggccc aagtgggagc
1561 ggggctgagg cgcgccgggg ctgggagaca gccatccgcc aggccctgat gcccgatcgc
1621 cttcaggacg caccagcgc cccaggccac gcgccccaca gacaagcttc tctgagcatc
1681 tctgtgtcca acagtcagat ccaagagaat gtggacattg cactgtctta ccagatcttc
1741 cctgacgaag tgcctgggctc agggcagttt ggagtggctc atggaggaaa acaccggaag
1801 acaggccggg acgtggcagt taagtcattt gacaaactgc gcttccctac caagcaggag
1861 agccagctcc ggaatgaagt ggccattctg cagagcctgc ggcatcccgg gatcgtgaac
1921 ctggagtgca tggcagagac gcctgagaaa gtgtttgtgg tgatggagaa gctgcatggg
1981 gacatgttgg agatgatcct gtccagttag aaggccggc tgcctgagcg cctcaccaag
2041 ttcctcatca ccagatcctt ggtggctttg agacacctc acttcaagaa cattgtccac
2101 tgtgacttga aaccagaaaa cgtgtgtctg gcatcagcag acccatttcc tcagggtgaag
2161 ctgtgtgact ttggctttgc tcgcatcatc ggcgagaagt cgttcccgcc ctgagtggg
2221 ggcacgccgg cctactggc accgaggtg ctgctcaacc agggctacaa ccgctcgtg
2281 gacatgtggt cagtggcggt gatcatgtac gtcagcctca gcggcacctt cccttcaac
2341 gaggatgagg acatcaatga ccagatccag aacgccgcct tcatgtaccg gccagcccc
2401 tggagccaca tctcagctgg agccattgac ctcatcaaca acctgctgca ggtgaagatg
2461 cgcaaacgct acagcgtgga caaatctctc agccaccctt ggttacagga gtaccagacg
2521 tggctggacc tccgagagct ggaggggaaag atgggagagc gatacatcac gcatgagagt
2581 gacgacgcgc gctgggagca gtttgagca gagcatccgc tgcctggggtc tgggctgccc
2641 acggacaggg atctcgggtg ggcctgtcca ccacaggacc acgacatgca ggggctggcg
2701 gagcgcacga gtgttctctg a
```

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.