

Certificate of Analysis

PKG1 α , active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-688, 14-688-K, 14-688M

Parent Lot # 2147495

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 PKG1 α , expressed by baculovirus in Sf21 insect cells. Purified using Ni²⁺/NTA agarose. Purity 86.7% by SDS-PAGE and Coomassie blue staining. MW = 80kDa.

Specific Activity (Parent lot# 2147495): 10980U/mg, where one unit of PKG1 α activity is defined as 1nmol phosphate incorporated into 200 μ M PAKtide (RRRLSFAEPG) per minute at 30°C with a final ATP concentration of 100 μ M.

Formulation: 1.163mg/ml of enzyme in 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.

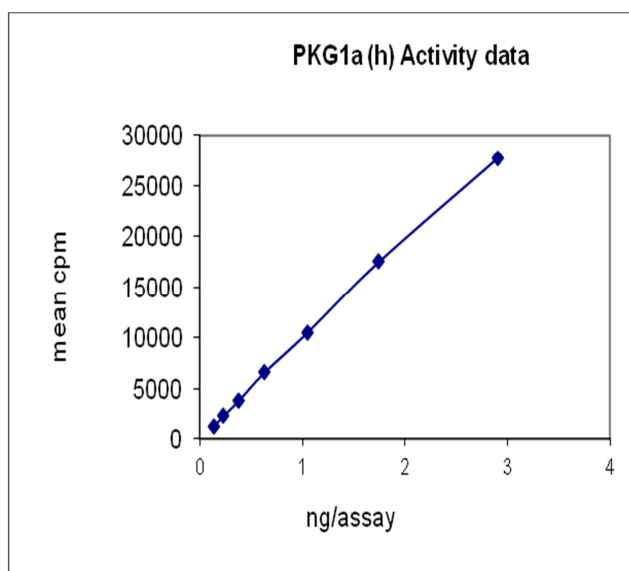
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

Handling Recommendations: Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled microcentrifuge tubes and immediately snap-freeze the vials in liquid nitrogen prior to re-storage at -70°C.

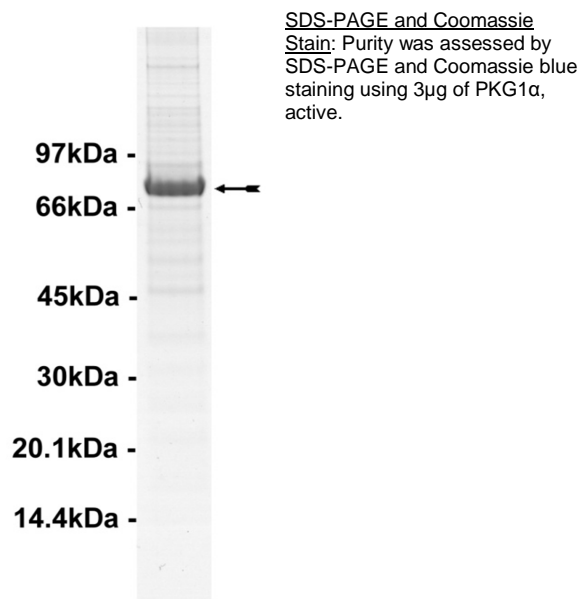
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 0.1–2.9ng of this lot of enzyme phosphorylated 200 μ M PAKtide (RRRLSFAEPG) 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 PKG1 α 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. **PAKtide (RRRLSFAEPG):** Use at a final assay concentration of 200 μ M. Make a 2mM stock. Add 2.5 μ l of stock per assay point.
3. **cGMP:** Use at a final assay concentration of 10 μ M. Make a 100 μ M stock. Add 2.5 μ l of stock per assay point.
4. **PKG1 α , active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 0.1–2.9ng per assay point.
5. **[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:

1. Add 5 μ l of 5 x reaction buffer per assay to wells.
2. Add 2.5 μ l of **PAKtide (RRRLSFAEPG)**.
3. Add 2.5 μ l of 100 μ M **cGMP**.
4. Add **2.5 μ l (0.1–2.9ng) PKG1 α , active**.
5. Add 2.5 μ l of dH₂O.
6. Add 10 μ l of diluted [γ -³³P]ATP mixture.
7. Incubate for 10 minutes at 30°C.
8. Stop the reaction by adding 5 μ l of 3% phosphoric acid.
9. Transfer a 10 μ l aliquot onto the appropriate area of a **P30 Filtermat**.
10. Wash the filtermat three times for 5 minutes with 75mM phosphoric acid.
11. Wash the filtermat once for 2 minutes with methanol.
12. Transfer the filtermat to a sealable plastic bag and add 4ml of scintillation cocktail.
13. 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

PKG1 Sequence Information

Protein	Human PKG1
Tags	N-terminal 6His
Native sequence	M31 of the recombinant protein is equivalent to M1 of native human PKG1 α
Accession number	GenBank D45864

Recombinant PKG1 α amino acid sequence:

```

1 MSYYHHHHHH DYDIPTTENL YFQGAMDPEF MSELEEDFAK ILMMLKEERIK ELEKRLSEKE
61 EEIQELKRKL HKCQSVLPVP STHIGPRTRR AQGISAEPQT YRSFHDLRQA FRKFTKSERS
121 KDLIKEAILD NDFMKNLELS QIQEIVDCMY PVEYGKDSCI IKEGDVGSVLV YVMEDGKVEV
181 TKEGVKLCMT GPGKVFGELE ILYNCTRAT VKTLVNVKLW AIDRQCFQTI MMRTGLIKHT
241 EYMEFLKSVP TFQSLPEEIL SKLADVLEET HYENGEYIIR QGARGDTFFI ISKGTVMVTR
301 EDSPSEDPVF LRTLKGDWF GEKALQGEDV RTANVIAAEA VTCLVIDRDS FKHLIGGLDD
361 VSNKAYEDAE AKAKYEAEEA FFANLKLSDF NIIDTLGVGG FGRVELVQLK SEESKTFAMK
421 ILKKRHIVDT RQQEHIRSEK QIMQGAHSDF IVRLYRTFKD SKYLYMLMEA CLGGELWTIL
481 RDRGSFEDST TRFYTACVVE AFAYLHSKGI IYRDLKPENL ILDHRGYAKL VDFGFACKIG
541 FGKKTWTFCG TPEYVAPEII LNKGHDISAD YWSLGILMYE LLTGSPPFSG PDPMKTYNII
601 LRGIDMIEFP KKIAKNAANL IKKLCRDNPS ERLGNLKNV KDIQKHKWE GFNWEGLRKG
661 TLTPPIIPSV ASPTDTSNFD SFPEDNDEPP PDDNSGWDID F
    
```

Recombinant PKG1 α nucleotide sequence:

```

1 atgtcgtact accatcacca tcaccatcac gattacgata tcccaacgac cgaaaacctg
61 tattttcagg ggcccatgga tccggaattc atgagcgagc tagaggaaga ctttgccaag
121 attctcatgc tcaaggagga gaggatcaaa gagctggaga agcggctgtc agagaaggag
181 gaagaaattc aggagctgaa gaggaaactc cacaaatgcc agtcggtgct cccagtgtccc
241 tcgaccacaca tcggcccccg gaccacccgg gcgcagggca tctcggccga gccgcagacg
301 tacaggtcct tccacgacct ccgacaggca ttccggaagt tcaccaagtc cgaaagggtcc
361 aaggatctta taaaggaagc tatccttgac aatgacttta tgaagaactt ggagctgtcg
421 cagatccagg agattgtgga ttgtatgtac ccggtggagt atggcaagga cagttgcatc
481 atcaaagaag gagacgtggg gtcactgggt tatgtcatgg aagatggtaa ggttgaagtt
541 acaaaagaag gtgtgaagtt gtgtaccatg ggtccaggaa aagtgtttgg ggaattggct
601 attctttaca actgtacccg gacagcgacc gtcaagactc ttgtaaagt aaaactctgg
661 gccattgatc gacaatgttt tcaacaata atgatgagga caggactcat caagcatacc
721 gagtatatgg aatttttaa aagcgttcca acattccaga gccttcctga agagatcctc
781 agcaagcttg ctgatgtcct tgaagagacc cactatgaaa atggagaata tattatcagg
841 caagggtgcaa gaggggacac cttctttatc atcagcaaag gaacggtaaa tgtcactcgt
901 gaagactcac cgagtgaaga cccagtcttt cttagaactt taggaaaagg agactggttt
961 ggagagaaaag cttgcaggg ggaagatgtg agaacagcaa acgtaattgc tgcagaagct
1021 gtaacctgcc ttgtgattga cagagactct tttaaacatt tgattggagg gctggatgat
1081 gtttctaata aagcatatga agatgcagaa gctaaagcaa aatatgaagc tgaagcggct
1141 ttcttcgcca acctgaagct gtctgatttc aacatcattg atacccttgg agttggagggt
1201 ttcggacgag tagaactggt ccagttgaaa agtgaagaat ccaaaacggt tgcaatgaag
1261 attctcaaga aacgtcacat tgtggacaca agacagcagg agcacatccg cttagagaag
1321 cagcatatgc agggggctca ttccgatttc atagttagac tgtacagaac atttaaggac
1381 agcaaatatt tgtatatggt gatggaagct tgtctagggt gagagctctg gaccattctc
1441 agggatagag gttcgtttga agattctaca accagatttt acacagcatg tgtgttagaa
1501 gcttttgctc atctgcattc caaaggaatc atttacaggg acctcaagcc agaaaatctc
1561 atcctagatc accgaggtta tgccaaactg gttgattttg gctttgcaaa gaaaatagga
1621 tttgaaaaga aaacatggac tttttgtggg actccagagt atgtagcccc agagatcatc
    
```

Certificate of Analysis

```
1681 ctgaacaaag gccatgacat ttcagccgac tactgggtcac tgggaatcct aatgtatgaa
1741 ctctgactg gcagcccacc tttctcaggc ccagatccta tgaaaaccta taacatcata
1801 ttgaggggga ttgacatgat agaatttcca aagaagattg ccaaaaatgc tgctaattta
1861 attaaaaaac tatgcaggga caatccatca gaaagattag ggaatttgaa aaatggagta
1921 aaagacattc aaaagcaca atggtttgag ggctttaact ggaaggctt aagaaaaggt
1981 accttgacac ctctataat accaagtgtt gcatcaccca cagacacaag taattttgac
2041 agtttccctg aggacaacga tgaaccacca cctgatgaca actcaggatg ggatatagac
2101 ttctaa
```

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