

Certificate of Analysis

PI3K-C2 β , active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-907, 14-907-K, 14-907M

Parent Lot # D9MN050N

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 PI3K-C2 β full length, expressed by baculovirus in Sf21 insect cells. Purified using Ni²⁺/NTA agarose.

Purity 40% by SDS-PAGE and Coomassie blue staining. MW = 189kDa.

Specific Activity (Parent lot# D9MN050N): 153U/mg, where one unit of PI3K-C2 β , active activity is defined as 1nmol phosphatidylinositol 3-phosphate (PI(3)P) formed per minute at room temperature with a final ATP concentration of 200 μ M.

Formulation: 0.323mg/ml of enzyme in 50mM Tris/HCl pH8.0, 300mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 0.2mM PMSF, 1mM benzamidine, 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.

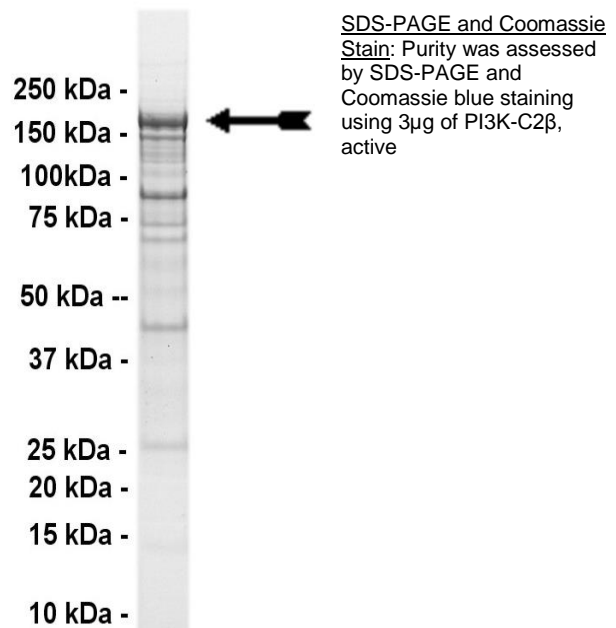
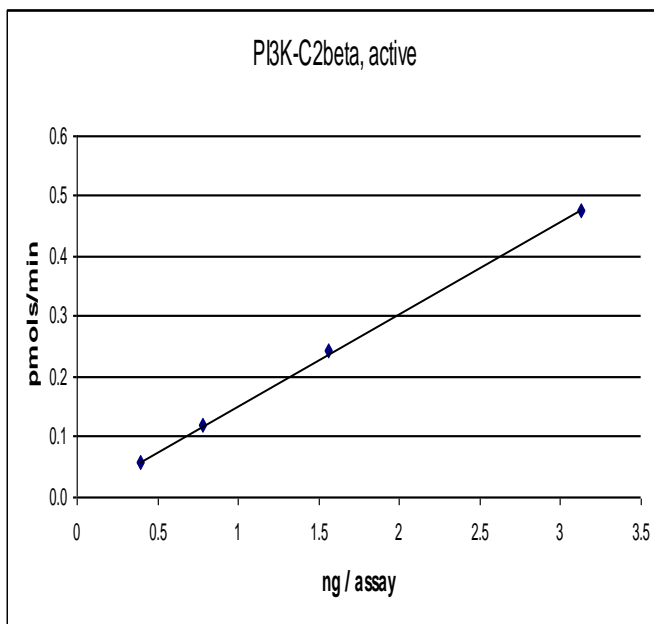
Handling Recommendations: Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled micro-centrifuge tubes and immediately snap-freeze the vials in liquid nitrogen prior to re-storage at -70°C.

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

Quality Control Testing

Kinase Assay: 0.4–3.2ng of this enzyme phosphorylated 25 μ M phosphatidylinositol in the assay referenced on page two.

MS Tryptic Fingerprint: Confirmed product identity as PI3K-C2 β , active with the translated sequence listed on page three



Certificate of Analysis

Eurofins PI3 Kinase Homogeneous Time-resolved Fluorescence (HTRF) Class II Reagent Kits

The following Eurofins kits are suitable for use with this enzyme. Assay protocol can be found in the corresponding Certificate of Analysis.

Cat. No	Kit Description
33-038	PI 3-Kinase (Class II) HTRF Assay 1-plate kit
33-039	PI 3-Kinase (Class II) HTRF Assay 5-plate kit

Certificate of Analysis

PI3K-C2 β Sequence Information

Protein	human PI3K-C2 β
Tags	N-terminal 6His
Native sequence	M31 of the recombinant protein is equivalent to M1 of human PI3K-C2 β
Accession number	GenBank NM_002646

Recombinant PI3K-C2 β amino acid sequence:

```

1 MSYYHHHHH DYDIPTTENL YFQGAMDPEF MSSTQDNGEH WKSLESVGIS RKELAMAEAL
61 QMEYDALSR L RHDKEENRAK QNADPSLISW DEPGVDFYSK PAGRRTDLKL LRGLSGSDPT
121 LNYNSLSPQE GPPNHSTSQG PQPGSDPWPK GSLSGDYLYI FDGSDGGVSS SPGPGDIEGS
181 CKKLSPPPLP PRASIWDTTP LPPRKGSPSS SKISQPSDIN TFSLVEQLPG KLEHRILEE
241 EEVLGGGGQG RLLGSVDYDG INDAITRLNL KSTYDAEMLR DATRGWKEGR GPLDFSKDTS
301 GKPVARSKTM PPQVPPRTYA SRYGNRK NAT PGKNRRISAA PVGSRPHTVA NGHELFEVSE
361 ERDEEVA AFC HMLDILRSGS DIQDYFLTGY VSAVTPSPE HLGDEVNLKV TVLCDRLQEA
421 LTFTCNCSST VDLLIYQ TLC YTHDDL RNVD VGDFVLKPCG LEEFLQNKHA LGSHEYIQYC
481 RKFDIDIRLQ LMEQKVRS D LARTVND DQS PSTLNYL VHL QERPVKQTIS RQALSLLFDT
541 YHNEVDA FLL ADGDFPLKAD RVVQSVK AIC NALAAVETPE ITSALNQLPP CPSRMQPKIQ
601 KDPSVLAVRE NREKVVEAL T AAIDLVELY CNTFNADFQT AVPGSRKHDL VQEACHFARS
661 LAFTVYATHR IPIIWATSYE DFYLSCSLSH GGKELCSPLQ TRRAHFSKYL FHLIVWDQOI
721 CFPVQVNR LP RETLLCATLY ALPIPPGSS SEANKQRRVP EALGWVTPL FNRQVLT CG
781 RKL LGLWPAT QENPSARWSA PNFHQPSVI LQIDFPTS AF DIKFTSPPGD KFSPRYEF GS
841 LREEDQRK LK DIMQKESLYW LTDADKRLW EKRYYCHSEV SSLPLVLASA PSWEWACL PD
901 IYVLLKQW TH MNHQDALGLL HATFPDQEV R RMAVQWIGSL SDAELLDYLP QLVQALKYEC
961 YLDSPLVR FL LKRAVSDLRV THYFFWLLKD GLKDSQFSIR YQYLLAALLC CCGKGLREEF
1021 NRQCWL VNAL AKLAQQVREA APSARQGILR TGLEEVKQFF ALNGSCLPL SPSLLVKGIV
1081 PRDCSYFN SN AVPLKLSFQN VDPLGENIRV IFKCGDDL RQ DMLTLQMIRI MSKIWVQEGL
1141 DMRMVI FRCF STGRGRGMVE MIPNAETLRK IQVEHGV TGS FKDRPLADWL QKHNPGEDEY
1201 EKAVENFI YS CAGCCVATYV LGICDRHNDN IMLKTTGHMF HIDFGRFLGH AQMFGNIKRD
1261 RAPFVFTSD M AYVINGGDKP SSRFHDFVDL CCQAYNLIRK HTHLFLNLLG LMLSCGIPEL
1321 SDLEDLKY VY DALRPQDTEA NATTYFTRLI ESSLGSVATK LNFFIHNL AQ MKFTGSDDRL
1381 TLSFASR TH T LKSSGRISDV FLCRHEKIFH PNKGYYVVK VMRENTHEAT YIQRTFEEFQ
1441 ELHNKLR L LF PSSHLPSFPS RFVIGRSRGE AVAERRREEL NGYIWHLIHA PPEVAECDLV
1501 YTFHFPL PRD EKAMGTSPAP KSSDGTWARP VGKVGGEVKL SISYKNNKLF IMVMHIRGLQ
1561 LLQDGNPD P YVKIYLLPDP QKTKRRTKV ARKTCNPTYN EMLVYDGIPK GDLQQRELQL
1621 SVLSEQGF WE NVLLGEVNIR LRELDLAEQEK TGWFALGSR S HGTL

```

Recombinant PI3K-C2 β nucleotide sequence:

```

1 atgtcgtact accatcacca tcaccatcac gattacgata tcccaacgac cgaaaacctg
61 tattttcagg gcgccatgga tccggaattc atgtcttcga ctcaggacaa tggggaacac
121 tggaagtccc tggagtcagt gggcatcagc cgcaaagagc tagcgatggc cgaagccctg
181 cagatggagt atgatgccct gtccccgctc cggcatgaca aggaggagaa cagagccaag
241 cagaacgcag acccctctct catcagctgg gatgagcctg gggtagactt ttacagcaag
301 ccagcaggaa ggcggaccga cctcaagctg ttacgcggtc tctctggctc tgatcctacc
361 ctaactaca actcactatc cccacaggaa gggccgcca accactctac ctcccaaggg
421 ccacagcctg gctcagatcc ctggcccaaa ggctcccctgt ctggagacta tctctacatt
481 tttgatggtt cagatggggg agtctcttcg tccccaggac caggggacat agagggtctc
541 tgcaagaaac tatccccacc tcctctgcct ccccgagctt ctatctggga taccctccc
601 ctgcctccca gaaaggggtc cccctcatcc tccaagatct cccagcccag tgacatcaac

```

Certificate of Analysis

```

661 actttctctt tggtcgaaca attgccaggc aaactgctag agcatcggat cctagaagag
721 gaagagggtgc tgggagggtgg gggtcagggg cgcctactgg ggtctgtgga ctatgatggt
781 atcaatgatg caattactag gctcaacttg aatcgacct atgatgcgga gatgttgagg
841 gatgccacca ggggctggaa ggagggccga gggccgctgg acttcagcaa agacacctct
901 ggaaaacccg tggccaggag caagactatg cccctcagg tgccccccg cacctatgcc
961 tcccgtctatg gcaaccgaaa gaatgcgacg cctggcaaga accgccggat ttctgcagcc
1021 ccgggtgggct cccggcccca cactgttgcc aatggccatg agttgtttga ggtctcagaa
1081 gagagagatg aggaggttgc tgcattttgc cacatgctgg atatccttcg atctggctct
1141 gacatccaag actacttctt cactggctat gtctggagtg ctgtcaccct tagcccagag
1201 cacctcgggg atgaggtcaa cctgaagggtg actgtgttgt gtgacaggct tcaagaggca
1261 ctacttttca cctgcaactg ttcctccact gtagacttgc ttatctacca gaccctgtgc
1321 tacacccatg atgacctgag gaatgtggac gtgggtgact ttgtgctaaa gccctgcggg
1381 ctggaggagt tcctgcagaa caagcatgcc ttgggcagtc atgagtacat ccaatactgc
1441 cgcaagtttg acattgacat tcggctacag ctgatggagc agaaggttgt gcgcagtgac
1501 ctggcccgga cggtgaatga tgaccagagc cctccacct tgaactacct cgtccatctc
1561 caagagaggc ctgtcaagca gaccatcagc aggcaggccc tgagtcttct gttcagacct
1621 taccacaatg aggtggatgc cttctgtctg gctgatggag acttccact gaaggctgac
1681 aggggtggtcc agtccgtcaa ggccatctgc aaccgctgg ccgccgtgga aaccctgag
1741 atcaccagtg ctctcaacca gctgccccc tgccccctcc gcatgcagcc taaaattcag
1801 aaggatccca gtgtcttggc tgtgagggaa aaccgagaga aggtcgtgga agccctgacc
1861 gctgccatct tggacctggt ggagctgtac tgcaacacat tcaacgcaga cttccagacg
1921 gcagtgcctg ggagccgcaa gcatgacctg gtccaggagg cctgccattt cgccaggtcc
1981 ctggccttca ctgtctatgc caccaccgc atccccatca tctgggctac cagctatgaa
2041 gatttctacc tctctgctc cctcagccat ggccggcaagg agctgtgcag ccccctgcag
2101 acccgaagag ctacttctc caagtacctc ttccacctca tcgtctggga ccagcagatc
2161 tgcttcccag tgcaggtgaa ccggctgcct cgggagacac tgctgtgtgc cactctctat
2221 gctctgccc tccccacc ggggagctcc tcagaggcca ataagcagcg gcgggtgcct
2281 gaagccctgg gctgggtcac taccctctc ttcaacttca ggcaggtcct gacctgtggc
2341 cggaaagctt tgggtttgtg gccagcaaca caggaaaatc ccagcgcctg ttggagtgca
2401 cctaatttcc accagccaga cagtgtcatc ctgcagattg acttccccac ctgcgctttt
2461 gacatcaagt tcaccagccc ccctggagac aagttcagcc cccgctatga gtttggcagc
2521 ctccgggaag aagaccagcg caagctaaa gacatcatgc agaaagagtc cttgtactgg
2581 ctactgatg ctgacaagaa ggcctgtgg gagaagcagc attactgcca ctgcggaggtg
2641 agctgcctcc ccctgggtgct cgccagcggc cccagctggg agtgggcttg cctgcctgac
2701 atctatgttc tcctgaagca gtggaccac atgaaccacc aggatgccct ggggctcctg
2761 catgccacct tcccggacca ggaggtgcgt cgtatggctg tgcatgggat tggctcactc
2821 tcagatgctg agctgctaga ctacctgcc cagctggtag aggcctgaa gtatgaatgc
2881 tacctggaca gcccgttggg gcgcttctc ctgaaacgag ctgtgtctga cttgagagtg
2941 actactact tcttctggtt actgaaggac ggcctcaagg actctcagtt cagcatccgc
3001 taccagtatc tgctggcagc cttactgtgc tgctgtggca aggggctgag agaagagttt
3061 aaccgccagt gctggcttgt caatgcctg gccaaactgg cccagcaggt ccgggaggca
3121 gccccatctg caaggcaggg aatcctccgc acgggcctgg aggaggtgaa gcagttcttt
3181 gccctcaatg gctcgtgccg cttgccactc agccccagtc tgctgggtaa ggggaattgtg
3241 cccagggact gttctactt caactcaat gctgtcccc tcaaactctc cttccaaaat
3301 gtggatcccc tgggtgagaa catccgtgtc atcttcaagt gtggggacga cttcggccag
3361 gacatgctaa cgctgcagat gattcgcac atgagcaaga tctgggtcca ggaggggctg
3421 gacatgcgca tggatcatct ccgctgctt tccaccggcc ggggcagagg gatggtggag
3481 atgatcccta atgctgagac cctgcgtaag atccaggtgg agcatggggg gaccggctcg
3541 ttcaaggacc ggcccctggc agactggctg cagaaacaca accctgggga ggacgagtat
3601 gagaaggctg tggagaactt tatctactc tgcgctggct gctgcgtggc cacgtacgtc
3661 ttgggcatct gtgaccgaca taatgacaac atcatgctga agaccactgg tcacatgttc
3721 cacattgatt ttggccgctt cctgggcat gcccagatgt ttggcaacat caagcgggac
3781 cgtgccccct ttgtcttcac ctccgacat gcgtatgtca tcaacggggg tgacaagcct
3841 tccagccgct tccatgattt tgttgacctt tgctgccaag cctacaacct cattcgcaag
3901 cacaccacc tcttctcaa cttctgggc ctgatgttgt cctgtgggat ccctgaactc
3961 tcagacctgg aggacctcaa gtatgtgtac gatgccctga ggcctcagga tacagaggcc
4021 aatgccacta cctacttcac taggttgatt gattccagcc tgggcagtg agccacaaag
    
```

Certificate of Analysis

```
4081 ctcaatTTTT tcatcataa tctggctcag atgaagttca cgggctcaga tgaccggctg
4141 accctctcct ttgcctccc aacacacact ctcaagagct ctggccgaat cagtgatggt
4201 ttctctgccc gccatgagaa gatcttccac cccaacaaag gctatatata tgtggtaaag
4261 gtgatgagag agaacactca cgaggccacc tacatccagc ggacctttga ggagttccag
4321 gaattacaca ataagttgag gctgctcttc ctttcttccc acttgcccag cttccctagt
4381 cgcttcgtga tcggccgctc ccggggagag gcggtggccg agcggcggag ggaggagcta
4441 aacggttaca tctggcactt gatccacgca ccccctgagg tggccgagtg tgatttgggt
4501 tacaccttct tccaccactt gccccgggat gagaaggcta tgggtaccag cccagctcct
4561 aagtcctcag atggcacatg ggccccggcc gtcggaaaagg tgggagggga ggtgaagctg
4621 tccatctcct acaaaaacaa taaactctt atcatggtga tgcatttcg gggcttgcaa
4681 ctgctccagg atggaatga ccctgacccc tatgtgaaaa tttacctcct tcctgaccct
4741 cagaaaacca ctaagaggaa aaccaaagt gcccggaaaa cctgcaatcc tacctacaat
4801 gagatggttg tatatgatgg gatccccaag ggtgacctgc agcagcggga gctccagctg
4861 agcgtgctga gtgagcaggg attctgggag aacgtcctc tcggtgaggt gaacatccgc
4921 ctgcgagagc tggacctggc tcaggagaag accggctggt tcgccctggg atctcgaagt
4981 catggcacct tgtga
```

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