

Certificate of Analysis

PI3K-C2 γ , active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-910, 14-910-K, 14-910M

Parent Lot # 2023057

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 human PI3K-C2 γ amino acids 468–1203, expressed by baculovirus in Sf21 insect cells. Purified using glutathione agarose.

Purity 70.5% by SDS-PAGE and Coomassie blue staining. MW = 112kDa.

Specific Activity (Parent lot# 2023057):

953U/mg, where one unit of PI3K-C2 γ , 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: 0.406mg/ml of enzyme in 50mM Tris/HCl pH7.5, 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.

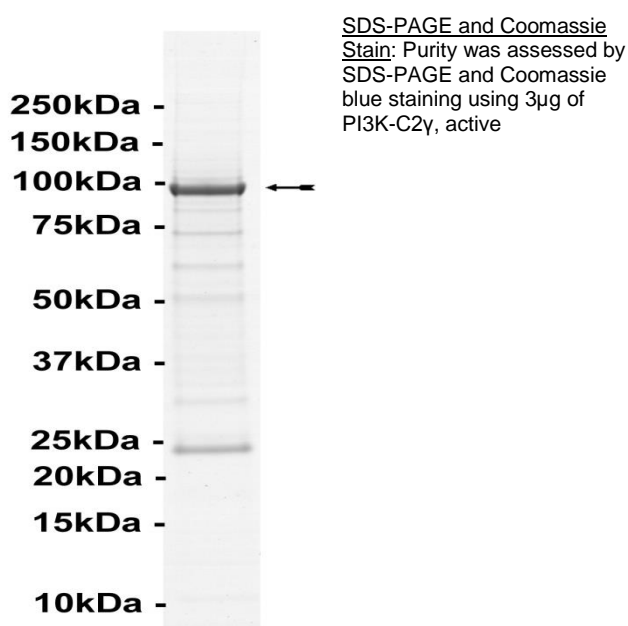
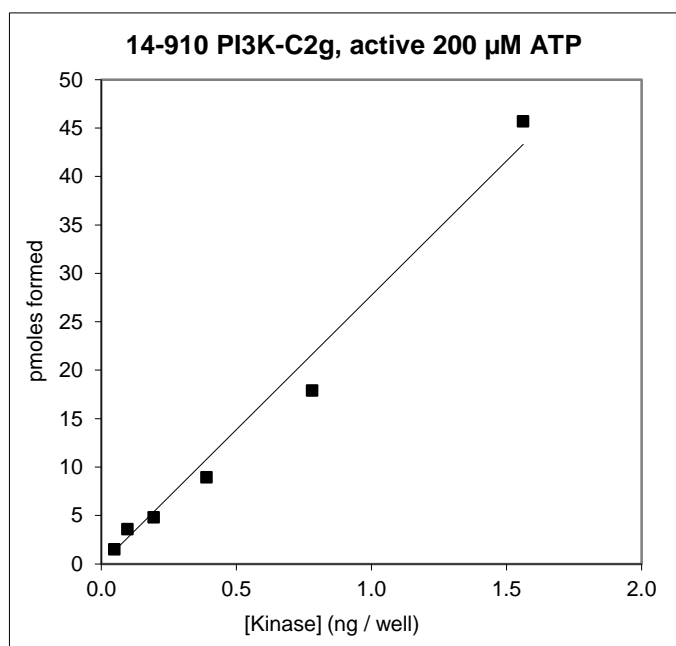
Handling Recommendations: Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled micro-centrifuge tubes and immediately snap-freeze the vials in liquid nitrogen prior to re-storage at -70°C.

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

Quality Control Testing

Kinase Assay: 0.05–1.56ng of this enzyme phosphorylated 25 μ M phosphatidylinositol in the assay referenced on page two.

MS Tryptic Fingerprint: Confirmed identity as PI3K-C2 γ 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 γ Sequence Information

Protein	human PI3K-C2 γ
Tags	N-terminal GST
Native sequence	Q237 of the recombinant protein is equivalent to Q468 of human PI3K-C2 γ
Accession number	GenBank BC130277

Recombinant PI3K-C2 γ amino acid sequence:

```

1 MSPILGYWKI KGLVQPTRL L LEYLEEKYEE HLYERDEGDK WRNKKFELGL EFPNLPYYID
61 GDVKLTQSMA IIRYIADKHN MLGGCPKERA EISMLEGAVL DIRYGVSRIA YSKDFETLKV
121 DFLSKLPEML KMFEDRLCHK TYLNGDHVTH PDFMLYDALD VVLYMDPMCL DAFPKLVCFK
181 KRIEAIQID KYLKSSKYIA WPLQGWQATF GGDHPPKSD LEVLFQGPEF KGLRRHQSSSE
241 TSAKGLIEKV TTELSTSIYQ LINVYCNSFY ADFQPVNVPR CTSYLNPLGLP SHLSFTVYAA
301 HNIPETWVHR INFPLEIKSL PRESMLTVKL FGIACATNNA NLLAWTCLPL FPKEKSILGS
361 MLFSMTLQSE PPVEMITPGV WDVSQSPSPT LQIDFPATGW EYMKPDSEEN RSNLEEPLKE
421 CIKHIARLSQ KQTPLLLSEE KKRYLWFYRF YCENNENCSLP LVLGSAPGWD ERTVSEMHTI
481 LRRWTFSQL EALGLLTSSF PDQEIRKQVAV QQLDNLLNDE LLEYLPQLVQ AVKFEWNLES
541 PLVQLLLHRS LQSIQVAHRL YWLLKNAENE AYFKSWYQKL LAALQFCAGK ALNDEFSKEQ
601 KLIKILGDIG ERVKSASDHQ RQEVLKKEIG RLEEFFQDVN TCHLPLNPAL CIKGIDHDAC
661 SYFTSNALPL KITFINANLM GKNISIFKA GDDLQDMLV LQLIQVMDNI WLQEGLDMMQ
721 IYRCLSTGK DQGLVQMPVD AVTLAKIHRH SGLIGPLKEN TIKKWFSQHN HLKADYEKAL
781 RNFYSCAGW CVVTFILGVC DRHNDNIMLT KSGHMFHIDF GKFLGHAQTF GGIKRDRAPF
841 IFTSEMEYFI TEGGKNPQHF QDFVELCCRA YNIIRKHSQL LLNLLEMMLY AGLPELSGIQ
901 DLKYVYNNLR PQDTDLEATS HFTKKIKESL ECFPVKLNLL IHTLAQMSAI SPAKSTSQTF
961 PQESCLLSTT RS

```

Recombinant PI3K-C2 γ nucleotide sequence:

```

1 atgtccccta tactaggtta ttggaaaatt aagggccttg tgcaaccac tcgacttctt
61 ttggaatata ttgaagaaaa atatgaagag catttgtatg agcgcgatga aggtgataaa
121 tggcgaaca aaaagtttga attgggtttg gagtttccca atcttcctta ttatattgat
181 ggtgatgta aattaacaca gtctatggcc atcatacgtt atatagctga caagcacaac
241 atgttgggtg gttgtccaaa agagcgtgca gagatttcaa tgcttgaagg agcggttttg
301 gatattagat acggtgtttc gagaattgca tatagtaaag actttgaaac tctcaaagtt
361 gattttctta gcaagctacc tgaatgctg aaaatgttcg aagatcgttt atgtcataaa
421 acatatttaa atggtgatca tgaaccat cctgacttca tgttgatga cgctcttgat
481 gttgttttat acatggacc aatgtgcctg gatgcgttcc caaaattagt ttgttttaa
541 aaacgtattg aagctatccc acaaattgat aagtacttga aatccagcaa gtatatagca
601 tggcctttgc agggctggca agccacgttt ggtggtggcg accatcctcc aaaatcggat
661 ctggaagttc tgttcaggg gcccgaattc aaaggcctac gtcgacatca aagttcagag
721 acttcagcaa aaggcttgat agagaaggta acaactgaac tatccacatc catctaccag
781 ctaatcaatg tctactgtaa cagcttttat gcagattttc agcctgtaaa tgtacctaga
841 tgacttctct atctaaatcc cgggcttctt tcccacctca gcttcacagt gtatgcagca
901 cacaacattc cagaaacctg ggtgcacagg atcaattttc cccttgaaat aaagtcactt
961 ccaagggat ccatgctcac tgtaaaactg tttgggattg cctgtgcaac caacaatgca
1021 aatttactgg cgtggacttg tcttccactg tttccaaaag aaaaatccat tctcgggtct
1081 atgctgttca gcatgacatt acagagttag cctcccgtag aatgataac tccaggagtg
1141 tgggatgtaa gtcagccatc cccggtgacc ctgcagattg attttccagc tactgggtgg
1201 gagtatatga aacctgattc tgaagagaat agaagtaatc ttgaagagcc actaaaggag
1261 tgtataaaac atattgccag actttcacag aacagactc ccctactact ctctgaagaa
1321 aagaaaagat atttatggtt ttatcgcttc tactgcaata atgaaaactg ctcccttctt

```

Certificate of Analysis

```

1381 ttagtcctgg gtagtcccc tggatgggat gaaaggactg tttcagaaat gcataccatt
1441 ttgagaagat ggacattttc tcaaccttta gaggctcttg ggcttttgac ttccagtttt
1501 ccagatcaag aaattcgtaa agtggcagtt caacaattag acaacctctt gaatgatgaa
1561 ctactggaat atctcccaca gctagtctag gctgtcaagt ttgaatggaa ccttgagagt
1621 ctttagtgct aacttctact ccaccgctcc ttgcagagca tccaggttgc ccatcgctct
1681 tactggctgc taaaaaatgc agaaaatgaa gcttatttta aaagctggta tcagaagcta
1741 ctactgctc tccaattctg tgcaggtaaa gccttgaatg atgagttttc caaggagcag
1801 aaacttatca aaattctggg agatattggg gaaagagtca agtctgccag tgaccatcaa
1861 agacaggagg tactgaagaa agaaattggc agactagaag agttctttca agatgtaaat
1921 acttgtcatc ttcctctgaa ccctgcccta tgtataaaaag ggattgatca cgatgcatgt
1981 tcatatttta catctaatac tttgccattg aagattactt tcatcaatgc taatctgatg
2041 ggcaaaaaca tcagcattat ttttaaggct ggagatgatc ttcgtcagga tatgcttggt
2101 ctgcagctta ttcaagtgat ggacaatatt tggctgcagg aaggcttgga tatgcaaagt
2161 atcatttata gatgtctatc cacagaaaaa gaccaaggat tggatgcagat ggtacctgat
2221 gctgtgacct tagcaaagat tcatcgccat tctggactga taggaccatt gaaagaaaaa
2281 acaattaaaa agtggttcag tcagcacaac cacttaaagg cagattatga aaaggccttg
2341 aggaactttt tctactcctg tgctggctgg tgtgtggtaa cattcatcct gggagtatgt
2401 gaccgtcaca atgataatat catgctgaca aagtcgggccc acatgtttca tattgacttt
2461 ggaaaattct taggtcatgc acaaacattt ggagggataa aaagggaccg agtcctttc
2521 atttttactt cagagatgga atactttatt acagagggtg ggaaaaacc acagcatttt
2581 caagattttg tggaaacttg ctgtcgtgct tataatatta tcagaaagca cagccaactg
2641 ctcttgaacc tgctggaat gatgctgtat gcaggactgc ctgagctaag tggaattcaa
2701 gacctgaaat atgtgtataa taatcttctg ccacaagaca cagacctgga agcaacaagt
2761 cattttacca agaaaataaa ggaaagtctg gagtgtttcc ctgttaaatt gaataacttg
2821 atccacacac ttgcacaaat gtcagccata agccctgcca aatctacttc acagactttt
2881 cctcaggaat cctgtttgct gagtacaact aggtcgtaa
  
```

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