

Data Sheet

p110 γ /PIK3R5

Human, recombinant

Catalog #: 40626

Lot #: 120605 **Conc.** 0.14 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 16 mM
Glutathione, 3 mM DTT, 20% Glycerol.

Stability: >6 months at -80°C

References:

1. Shymanets, A. *et al.*, *Biochem. J.*
2012 Feb 1;**441(3)**: 851-8.
2. Brazzatti, J.A., *et al.*, *Oncogene.*
2012 May 3;**31(18)**:2350-61.

Description:

Complex of N-terminal His-tag recombinant human p110 γ , or PIK3CG, (GenBank Accession No. NM_002649) a.a. 2-1102 (end), and N-terminal GST-tag recombinant human PIK3R5 (GenBank Accession No. NM_001142633) a.a. 2 - 880 (end). Co-expressed in a Baculovirus infected Sf9 cell expression system. p110 γ MW=127 kDa, PIK3R5 MW=124 kDa.

Application:

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

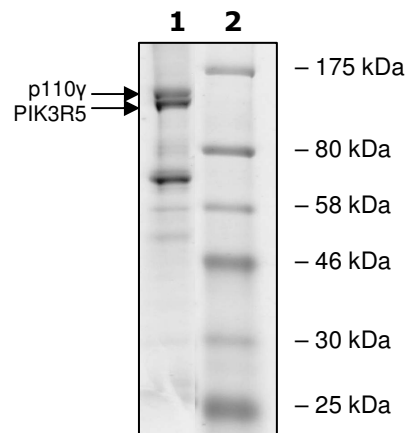
10% SDS-PAGE
Coomassie staining

Lane 1:
1.4 μ g p110 γ /PIK3R5

Lane 2:
Protein Marker
BioLabs (#P7708L)

MW:
p110 γ : 127 kDa
PIK3R5: 124 kDa

Purity: \geq 40%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Protein Sequence

p110y

MHHHHHHELENYKQPVVLRDNCRRRRRMKPRSAASLSSMELIPIEFVLPTSQRKC
KSPETALLHVAGHGNVEQMKAQVWLRALETSVAADFYHRLGPHHFLLLYQKKGQW
YEIYDKYQVVQTLDCRLRYWKATHRSPGQIHLVQRHPPSEESQAFQRQLTALIGYDVTD
VSNVHDELEFTRRGLVTPRMAEVASRDPKLYAMHPWVTSKPLPEYLWKKIANNCFI
VIHRSTTSQTIKVPDDTPGAILQSFFTKMAKKKSLMDIPESQSEQDFVLRVCGRDEYL
VGETPIKNFQWVRHCLKNGEEIHVVLDTPDPALDEVKKEEWPLVDDCTGVTGYHEQ
LTIHGKDHESVFTVSLWDCDRKFRVKIRGIDIPVLPNTDLTVFVEANIHQGQQVLCQR
RTSPKPFTEEVLWNVWLEFSIKIKDLPGALLNLQIYCGKAPALSSKASAESPSSSEK
GKVQLLYVNNLLIDHRFLLRRGEYVLHMWQISGKGEDQGSFNADKLTSAATNPDKEN
SMSISILLDNYCHPIALPKHQPTDPEGDRVRAEMPQLRKQLEAIIATDPLNPLTAED
KELLWHFRYESLKHKPKAYPKLFSSVKWGQQEIVAKTYQLLARREVWDQSALDVGLT
MQLLDCNFSDENVRAIAVQKLESLEDDDLVHLLQLVQAVKFEPYHDSALARFLLKR
GLRNKRIGHFLFWFLRSEIAQSRHYQQRFAVILEAYLRGCGTAMLHDFQQVQVIEML
QKVTLDIKLSAEKYDVSSQVISQLKQKLENLQNSQLPESFRVPYDPGLKAGALAEK
CKVMASKKKPLWLEFKCADPTALSNETIGIIFKHGDDLQDMLILQILRIMESIWETESL
DLCLLPYGCISTGDKIGMIEIVKDATTIAKIQQSTVGNTGAFKDEVLNHWLKEKSPTEE
KFQAAVERFVYSCAGYCVATFVLGIGDRHNDNIMITETGNLFHIDFGHILGNYKSFLGIN
KERVPFVLTDFLVMGTSGKKTSPHFQKFQDICVKAYLALRHHTNLLILFSMMLMTG
MPQLTSKEDIEYIRDALTVGKNEEDAKKYFLDQIEVCRDKGWTVQFNWFLHLVLGIKQ
GEKHS

PIK3R5

MSPILGYWKIKGLVQPTRLLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMAIRYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSRIAYSKDFETLK
VDFLSKLPEMLKMFKDRLLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIQIDKYLKSSKYIAWPLQGWQATFGGGDHPPKSDPLVPRGSPGEFQPGA
TTCTEDRIQHALLERCLHGLSLSRRSTSWAGLCLNCWSLQELVSRDPGHFLILLEQIL
QKTREVQEKGTDLLTPLALLFYSTVLCTPHFPPDSDLLLKAASYHRFLTWPVYCS
ICQELLTFIDAELKAPGISYQRLVRAEQGLPIRSHRSSTVTVLLNPNVEVQAEFLAVAN
KLSTPGHSPHSAYTTLLHAFQATFGAHCDDVPLHCRLLQAKTLAELEDIFTETAEAEQ
LASGIGDAAEARWLRTKLQAVGEKAGFPGLVDTAKPGKLHTIPIVARCYTYSWSQ
DSFDILQEILLKEQELLQPGILGDDEEEEEEEEEVEEDLETGDHCAERDSSLSTSSLAS
HDSTLSLASSQASGPALSRHLLTSFVSGLSGMDSGYVEDSESSSEWPWRRGSQE
RRGHRRPGQKFIRIYKLFKSTSQVLRRDRSLEGSSDTALPLRRAGSLCSPLDEPVS
PPSRAQRSRSLPQPKLGTQLPSWLLAPASRPQRRRPFLSGDEDPKASTLRVVVFGS
DRISGKVARAYSNLRRLENNRPLLTRFFKLQFFYVPVKRSHGTSPGACPPPRSQTSP
PPTDSPRHASPGELGTPWEESTNDISHYLGMLDPWYERNVGLMLHLPPEVLCQQSL
KAEAQALEGSPTQLPILADMLLYCRFAARPVLLQVYQTELTFTITGEKTTEIFIHSLELG
HSAATRAIKASGPGSKRLGIDGDREAVPLTLQIIYSKGAISGRSRWSNLEKVCTSVNLN
KACRKQEELDSSMEALTLNLTEVVKRQNSKSKKGFNQISTSIKVDKVIIGNSNCPF
AVCLDQDERKILQSVRCEVSPCYKPEKSDLSPPQTPDLPAQAAPDLCSLLCLPIM
TFSGALP

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone 1.858.829.3082 Fax 1.858.481.8694
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com