

PI3K (p55 gamma) Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # P31-30CH

Lot # S298-1

Product Description

Recombinant full length human PI3K (p55 gamma) was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag. The gene accession number is [BC021622](#).

Gene Aliases

PIK3R3, p55, p55-GAMMA, FLJ41892, DKFZp686P05226

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

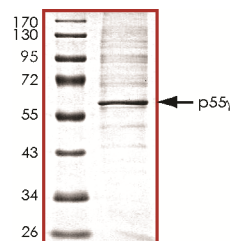
Scientific Background

PI3K (p55γ) also known as PI3K regulatory subunit 3 (PIK3R3) was first identified by 2-hybrid system as a protein that bind to the intracellular domain of IGF1R. PIK3R3 contains 2 Src homology-2 (SH2) domains, an inter-SH2/p110-binding site and a proline-rich motif. PIK3R3 transcript is detected in all fetal and adult tissues except peripheral blood leukocytes. Expression is low in adult liver and most abundant in adult testis. PIK3R3 can bind IGF1R, INSR, and INSR substrate-1 in vitro and the binding requires receptor activation (1).

References

1. Dey, B R. et al: Cloning of human p55-gamma, a regulatory subunit of phosphatidylinositol 3-kinase, by a yeast two-hybrid library screen with the insulin-like growth factor-1 receptor. Gene 209: 175-183, 1998.

Purity



The purity of PI3K (p55 gamma) was determined to be **>80%** by densitometry. Approx. MW **61 kDa**.

PI3K (p55 gamma) Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	P31-30CH
Specific Lot Number	S298-1
Purity	>80%
Concentration	0.2µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.