

## PI3K (p85β) Protein

Full length recombinant protein expressed in Sf9 cells

**Catalog # P31-30BH**

Lot # N104-3

### Product Description

Recombinant full length human PI3K (p85β) was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag. The gene accession number is [NM\\_005027](#).

### Gene Aliases

PIK3R2, P85B, p85-BETA

### 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.

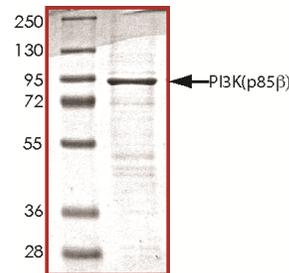
### Scientific Background

Phosphatidylinositol 3-kinase (PI3K) is a lipid kinase that phosphorylates the inositol ring of phosphatidylinositol and related lipid products at the 3-prime position. PI3K p85 subunit which is the regulatory subunit of PI3K and consists of 2 closely related proteins, p85-alpha and p85-beta (1). The activation of PI3K signaling in influenza A virus-infected cells is important for efficient virus replication (2). PI3K signaling in T cells can lead to organ-specific autoimmunity and the class IA PI3K-deficient mice manifest the cardinal features of human primary Sjogren's syndrome-like disease (3).

### References

1. Otsu, M., et.al: Characterization of two 85 kd proteins that associate with receptor tyrosine kinases, middle-T/pp60(c-src) complexes, and PI3-kinase. Cell 65: 91-104, 1991.
2. Hale, B. G. et.al: Influenza A virus NS1 protein binds p85-beta and activates phosphatidylinositol-3-kinase signaling. Proc. Nat. Acad. Sci. 103: 14194-14199, 2006.
3. Oak, J. S., et.al: Sjogren's syndrome-like disease in mice with T cells lacking class 1A phosphoinositide-3-kinase. Proc. Nat. Acad. Sci. 103: 16882-16887, 2006.

### Purity



The purity of PI3K (p85β) was determined to be **>85%** by densitometry. Approx. MW **88 kDa**.

## PI3K (p85β) Protein

Full length recombinant protein expressed in Sf9 cells

Catalog #	P31-30BH
Lot #	N104-3
Purity	>85%
Concentration	0.1µg/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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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