



Mouse (monoclonal) Anti-PI3-Kinase p85 β Subunit

PRODUCT ANALYSIS SHEET

Catalog Number:	AHO0572
Lot Number:	See product label
Quantity:	2 mL
Clone Number:	T15
Isotype:	IgG1 (mouse)
Form of the Antibody:	Tissue culture supernatant.
Preservation:	0.02% sodium azide (Caution: sodium azide is a poisonous and hazardous substance. Handle with care and dispose of properly.)
Immunogen:	Recombinant bovine p85 β subunit of PI3-kinase, expressed in baculovirus.
Myeloma/Fusion Partners:	Immunized mouse splenocytes were fused with NS1 mouse myeloma cells.

Specificity: Phosphoinositide 3-kinase (PI3-kinase, PI3K) catalyzes the addition of a phosphate group to the 3' hydroxyl group of phosphatidylinositol, yielding PI 3,4 bisphosphate (PIP₂) and PI 3,4,5 trisphosphate (PIP₃). PIP₃ serves to localize proteins containing a Pleckstrin homology (PH) domain, a feature present in many classes of signaling proteins (serine/threonine kinases, tyrosine kinases, GTPase activating proteins, cytoskeletal proteins, etc.), to the plasma membrane. Once recruited to the plasma membrane by PIP₃, PH domain-containing proteins are co-localized with other membrane-associated signaling proteins, where they can influence signaling pathways.

PI3K is an integral part of several signaling cascades which culminate in cell proliferation, survival, cytoskeletal organization, membrane trafficking, membrane ruffling, differentiation, chemotaxis, and glucose homeostasis. In growth factor signaling, receptor ligand binding stimulates tyrosine phosphorylation of the receptor and downstream substrates. These regions of phosphorylated tyrosine residues serve as docking sites for PI3K via its SH2 domains. The association of PI3K with phosphorylated growth factor receptors localizes PI3K to the plasma membrane, bringing it in close proximity to its inositol lipid substrates. PI3K binding to downstream substrates, such as FAK and Pyk2, also allows these signaling proteins to direct PI3K.

Enzymes which are regulated by PI3K are numerous and include phosphoinositide dependent kinases (PDK's), integrin-linked kinase (ILK), MAPK, Bruton's tyrosine kinase, and PKC ξ . PI3K's regulation of PDK1 is of great interest because this kinase phosphorylates and thereby activates the serine/threonine kinase Akt/PKB. Once activated, Akt/PKB then phosphorylates and in turn inhibits the activities of many targets including GSK-3 α and β , Forkhead-related transcription factor 1 (FKHR-L1), and Bad.

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PI AHO0572

(Rev 10/08) DCC-08-1089

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.

Specificity (cont'd):

PIP₃ also localizes phospholipase C_γ to the plasma membrane through its PH domain. By recruiting PLC_γ to the membrane, PI3K increases the rate of hydrolysis of phosphatidylinositol 4,5-bisphosphate, generating the second messengers diacylglycerol (DAG), activating PKC, and inositol 1,4,5-trisphosphate (IP₃), releasing Ca²⁺ from intracellular stores which in turn activates Ca²⁺ sensitive protein such as the focal adhesion kinase family member Pyk2/RAFTK.

PI3K's comprise a family of homologous heterodimeric proteins composed of a catalytic subunit and a regulatory subunit. Three isoforms of the catalytic subunit have been described, p110 α , p110 β , and p110 δ , each of which is encoded by its own gene. Five isoforms of the regulatory subunit have been described, p85 α , p85 β , p55 α , p55 γ , and p50 α . The p85 α gene encodes three of these regulatory subunits, produced through alternative splicing of the transcripts. The three products of the p85 α gene are p85 α , p50 α , and p55 α regulatory subunits. The p85 β and the p55 γ regulatory subunits are each encoded by a single gene. All of the regulatory subunits possess two SH2 domains with an intervening catalytic subunit-binding domain. The regulatory subunits also possess variable numbers of Bcr, SH3, and PRM domains. The presence of these various domains allows the regulatory subunits to serve as adaptor proteins.

This antibody recognizes the p85 β subunit of PI3K.

Species Reactivity:

Human, monkey, mouse, rat, and bovine. Other species were not tested.

Applications:

This monoclonal antibody is suitable for use in immunohistochemistry with cryostat sections, immunoprecipitation, and Western blot analysis.

Suggested Working Dilutions:

The recommended dilution for Western blot analysis is 1:1000. The optimal concentration should be determined for each specific application.

Storage:

Store at 2-8°C. For long term storage, apportion into working aliquots and store at -20°C. Avoid repeated freeze-thaw cycles to prevent denaturing the antibody.

Expiration Date:

Expires one year from date of receipt when stored as instructed.

Related Products:

Rabbit (polyclonal) anti-AKT/PKB [pT308]	Cat. #	44-602G
Rabbit (monoclonal) anti-AKT/PKB [pS473]	Cat. #	44-621G
Rabbit (polyclonal) anti-GSK-3 β [pS9]	Cat. #	44-600G
Rabbit (polyclonal) anti-GSK-3 α [pY279]/GSK-3 β [pY216]	Cat. #	44-604G
Rabbit (polyclonal) anti-GSK-3 α / β	Cat. #	44-610
Rabbit (polyclonal) anti-PLC γ -1 [pY783]	Cat. #	44-696
Mouse (monoclonal) anti-PLC γ -1	Cat. #	AHO0792
Rabbit (polyclonal) anti-ERK1/2 [pTpY185/187]	Cat. #	44-680G
Rabbit (polyclonal) anti-BAD [pS112]	Cat. #	44-522
Rabbit (polyclonal) anti-BAD [pS136]	Cat. #	44-524

References:

Cantley, L.C. (2002) The phosphoinositide 3-kinase pathway. *Science* 296:1655-1657 (review).

Katso, R., K. Okkenhaug, K. Ahnadi, S. White, J. Timms, and M.D. Waterfield (2001) Cellular function of phosphoinositide 3-kinases: implications for development, homeostasis, and cancer. *Ann. Rev. Cell Dev. Biol.* 17:615-675 (review).

Wymann, M.P. and L. Pirola (1998) Structure and function of phosphoinositide 3-kinases. *Biochim. Biophys. Acta* 1436:127-150 (review).

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PI AHO0572

(Rev 10/08) DCC-08-1089

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.