

Anti PI 3-Kinase gamma (IP), purified anti-human phosphoinositide 3-kinase γ , recognizes full length protein mouse, monoclonal

Cat. No.	Amount
ABD-027P	100 μ g

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Application

mAb p110g can be used for IP and Immunocyto-chemistry

Form

Liquid. Supplied in 50 mM Tris pH 8.0 and 50% glycerol.

Description

Monoclonal Antibody (mAb) p110 γ is a mouse IgG2a subtype purified from Hybridoma cell culture supernatant with Protein G-Sepharose.

Immunogen

Recombinant human full length protein PI3K γ (p110 γ).

Specificity

mAb p110 γ recognizes human, rat and mouse PI3K γ (p110 γ). The antibody does not cross-react with p110 α , p110 β and p110 δ .

Selected References:

Leopoldt *et al.* (1998) Gbetagamma stimulates phosphoinositide 3-kinase-gamma by direct interaction with two domains of the catalytic p110 subunit. *J. Biol. Chem.* **273**:7024.