

## Anti PI 3-Kinase gamma (IP), cell culture supernatant anti-human phosphoinositide 3-kinase $\gamma$ , recognizes full length protein mouse, monoclonal

Cat. No.	Amount
ABD-027	1 ml

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

### Avoid freeze / thaw cycles

### Application

mAb p110 $\gamma$  can be used for IP and Immunocyto-chemistry

### Form

Liquid. Supplied as cell culture supernatant in Dulbecco's Modified Eagles Medium, 10% FCS, 0.02% sodium azide.

### Description

This Monoclonal Antibody (mAb) p110 $\gamma$  is a mouse IgG2a subtype antibody.

### Immunogen

Recombinant human full length protein PI3K $\gamma$  (p110 $\gamma$ ).

### Specificity

mAb p110 $\gamma$  recognizes human, rat and mouse PI3K $\gamma$  (p110 $\gamma$ ). The antibody does not cross-react with p110 $\alpha$ , p110 $\beta$  and p110 $\delta$ .

### 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.  
Lehmann *et al.* (2009) PI3K $\gamma$  controls oxidative bursts in neutrophils via interactions with PKC $\alpha$  and p47phox. *Biochem. J.* **419**:603-610.