

## PPAR- $\gamma$ (Ab-112) Antibody [#B0557]

<b>Catalog Number:</b>	B0557
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P37231
<b>Other Names:</b>	NR1C3, PPAR-gamma-1, Peroxisome proliferator activated receptor gamma
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized non-phosphopeptide derived from human PPAR- $\gamma$ around the phosphorylation site of serine 112 (P-A-S <sup>P</sup> -P-P).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	PPAR- $\gamma$ (Ab-112) antibody detects endogenous levels of total PPAR- $\gamma$ protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
<b>References:</b>	Filiz Akbiyik, Blood, Sep 2004; 104: 1361 - 1368. Robert B. Clark, J. Immunol., Feb 2000; 164: 1364. JOSUÉ PADILLA, Ann. N.Y. Acad. Sci., Apr 2000; 905: 97. Chinyu G. Su, J. Clin. Invest., Aug 1999; 104: 383.