

PPAR- γ antibody (mAb)

Catalog Nos: 61025, 61026

RRID: AB_2793477

Clone: 3PP-1H4

Isotype: IgG1

Application(s): WB

Reactivity: Human

Quantity: 100 μ g

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 μ g/ μ l

Molecular Weight: 55 kDa

Background: Peroxisome Proliferator-Activated Receptors (PPARs) are nuclear receptors involved in lipid transport and metabolism. As such, their roles in chronic diseases such as diabetes, obesity, atherosclerosis and cancer are heavily investigated. Transcriptional activity of PPARs is regulated by fatty acid binding. Three PPAR isotypes have been identified: α , δ and γ . PPAR- γ stimulates lipolysis of circulating triglycerides and the subsequent uptake of fatty acids into adipose cells. PPARs bind to peroxisome proliferator response elements (PPREs) as heterodimers with the retinoid X receptor (RXR).

Immunogen: This PPAR-gamma antibody was raised against a peptide within the N-terminal region of human PPAR gamma 1 and gamma 2.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Published Applications:

WB

See references for more information. Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Upon receipt, unconjugated antibodies may be stored at -20°C for up to 2 years. Fluorophore- & enzyme-conjugated antibodies should be stored at 4°C. Fluorophore-conjugated antibodies should be protected from light. Keep reagents on ice when not in storage; to avoid repeated freeze/thaw cycles, we recommend aliquoting items that will be stored frozen into single-use fractions prior to freezing. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.