
PPAR gamma Polyclonal Antibody

(Catalog # A71563)

Background

This gene encodes a member of the peroxisome proliferator-activated receptor (PPAR) subfamily of nuclear receptors. PPARs form heterodimers with retinoid X receptors (RXRs) and these heterodimers regulate transcription of various genes. Three subtypes of PPARs are known: PPAR-alpha, PPAR-delta, and PPAR-gamma. The protein encoded by this gene is PPAR-gamma and is a regulator of adipocyte differentiation. Additionally, PPAR-gamma has been implicated in the pathology of numerous diseases including obesity, diabetes, atherosclerosis and cancer. Alternatively spliced transcript variants that encode different isoforms have been described.

Description

PPAR gamma Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.01% thimerosal, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P37231

Purification

Affinity Purification

Immunogen

A synthetic peptide corresponding to a sequence within amino acids 100-200 of human PPAR γ (NP_056953.2).

Storage

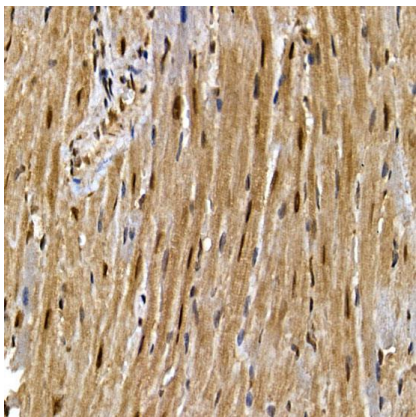
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

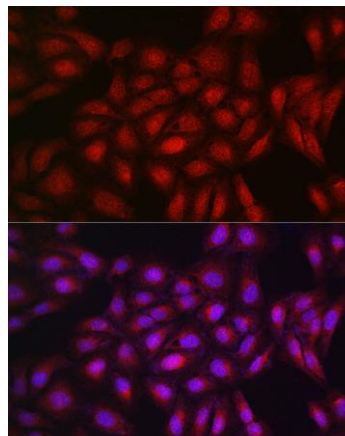
PPARG; CIMT1; GLM1; NR1C3; PPARG1; PPARG2; PPARgamma; peroxisome proliferator-activated receptor gamma

Application

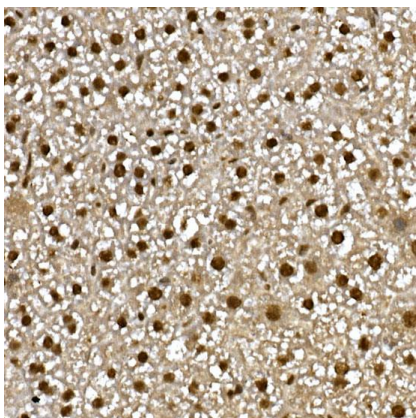
IHC, IF; Recommended dilution: IHC 1:50 - 1:200, IF 1:50 - 1:200



Immunohistochemistry analysis of PPAR γ in paraffin-embedded mouse heart using PPAR gamma Polyclonal Antibody at dilution of 1:200 (40x lens). Perform high-pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



Immunofluorescence analysis of U2OS cells using PPAR gamma Polyclonal Antibody at dilution of 1:200 (40x lens). Secondary antibody: Cy3 Goat Anti-Rabbit IgG (H+L) at 1:500 dilution. Blue: DAPI for nuclear staining.



Immunohistochemistry analysis of PPAR γ in paraffin-embedded mouse liver using PPAR gamma Polyclonal Antibody at dilution of 1:200 (40x lens). Perform high-pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.