
PPAR gamma Polyclonal Antibody

(Catalog # A71564)

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

Recombinant fusion protein containing a sequence corresponding to amino acids 1-120 of human PPARγ (NP_056953.2).

Storage

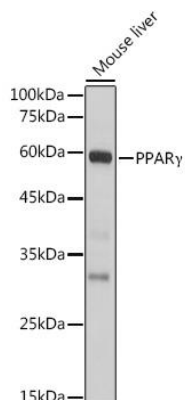
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

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

Application

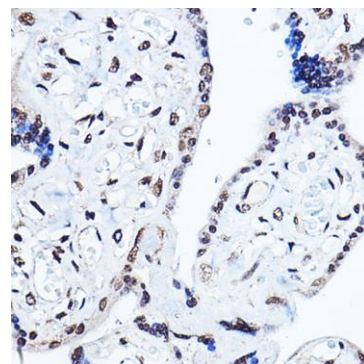
WB, IHC; Recommended dilution: WB 1:500 - 1:1000, IHC 1:100 - 1:500



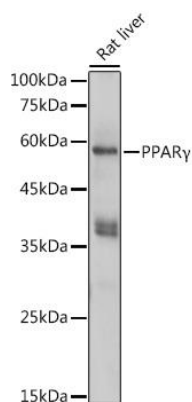
Western blot analysis of extracts of Mouse liver, using PPAR gamma Polyclonal Antibody at 1:1000 dilution. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25µg per lane.

Blocking buffer: 3% nonfat dry milk in TBST.



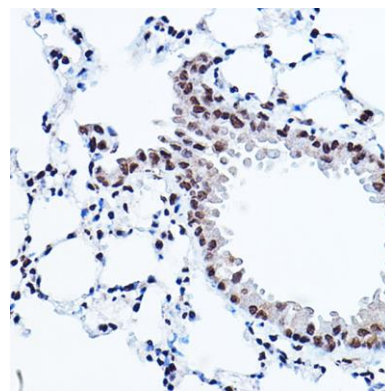
Immunohistochemistry analysis of paraffin-embedded human placenta using PPAR gamma Polyclonal Antibody at dilution of 1:500 (40x lens). Perform high-pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



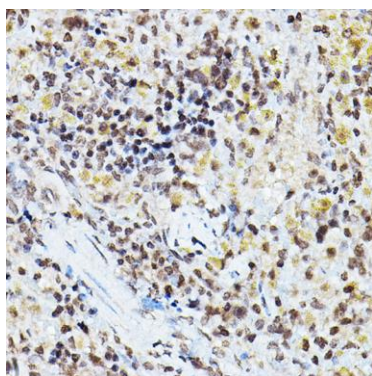
Western blot analysis of extracts of Rat liver, using PPAR gamma Polyclonal Antibody at 1:1000 dilution. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25µg per lane.

Blocking buffer: 3% nonfat dry milk in TBST.



Immunohistochemistry analysis of paraffin-embedded mouse lung using PPAR gamma Polyclonal Antibody at dilution of 1:500 (40x lens). Perform high-pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



Immunohistochemistry analysis of paraffin-embedded rat spleen using PPAR gamma Polyclonal Antibody at dilution of 1:500 (40x lens). Perform high-pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.
