

PRODUCT INFORMATION



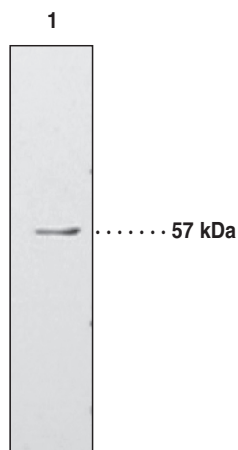
PPAR γ Polyclonal Antibody

Item No. 101700

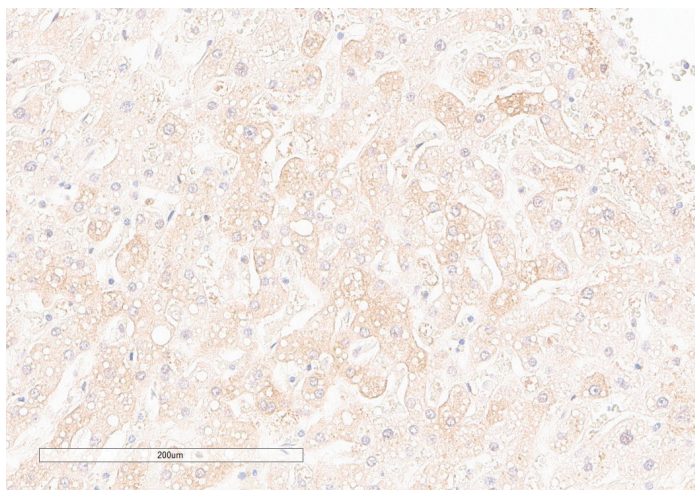
Overview and Properties

| | |
|----------------------------|---|
| Contents: | This vial contains 500 μ l of peptide affinity-purified polyclonal antibody. |
| Synonym: | Peroxisome Proliferator-activated Receptor γ |
| Immunogen: | Synthetic peptide from the internal region of PPAR γ 1 |
| Cross Reactivity: | (+) PPAR γ 1 and PPAR γ 2 |
| Species Reactivity: | (+) Human, mouse, rat; other species not tested |
| Uniprot No.: | P37231 |
| Form: | Liquid |
| Storage: | -20°C (as supplied) |
| Stability: | \geq 3 years |
| Storage Buffer: | PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide |
| Host: | Rabbit |
| Applications: | Immunohistochemistry (IHC), immunoprecipitation (IP), and Western blot (WB); the recommended starting dilutions are 1:80, 12 μ l of IgG per reaction, and 1:200, respectively. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically. |

Images



Lane 1: Rat adipose homogenate (30 μ g)



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) human liver tissue after heat induced antigen retrieval in pH 6.0 citrate buffer. After incubation with PPAR γ Polyclonal Antibody, (Item No. 101700), at a 1:80 dilution, slides were incubated with biotinylated secondary antibody, followed by alkaline phosphatase-streptavidin and chromogen (DAB).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/01/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

PPAR γ is a ligand-activated transcription factor involved in the regulation of lipid homeostasis and may function as a master regulator of adipogenesis.¹⁻⁴ PPAR γ mRNA is expressed prominently in adipose tissue, but is also found in large intestine, kidney, liver, and small intestine.⁵ Alternative splicing of the PPAR γ gene results in at least two mRNA species that differ at their 5' ends.^{1,6} Human PPAR γ 1 and PPAR γ 2 proteins are 53 and 57 kDa, respectively, based on the deduced amino acid sequences.⁶ PPAR γ 2 is the major PPAR γ isoform found in both the cytosolic and nuclear fractions of undifferentiated 3T3-L1 cells.⁷ Total cellular PPAR γ 2 protein increases approximately 2-fold following differentiation of 3T3-L1 cells, whereas only small quantities of PPAR γ 1 are detected in the nuclear fraction following differentiation.⁷

References

1. Greene, M.E., Blumberg, B., McBride, O.W., et al. *Gene Expression* **4**, 281-299 (1995).
2. Elbrecht, A., Chen, Y., Cullinan, C.A., et al. *Biochem. Biophys. Res. Commun.* **224**, 431-437 (1996).
3. Spiegelman, B.M. and Flier, J.S. *Cell* **87**, 377-389 (1995).
4. Lemberger, T., Desvergne, B., and Wahli, W. *Annu. Rev. Cell Dev. Biol.* **12**, 335-363 (1996).
5. Fajas, L., Auboeu, D., Raspé, E., et al. *J. Biol. Chem.* **272**, 18779-18789 (1997).
6. Mukherjee, R., Jow, L., Croston, G.E., et al. *J. Biol. Chem.* **272**, 8071-8076 (1997).
7. Thuillier, P., Baillie, R., Sha, X., et al. *J. Lipid Res.* **39**, 2329-2338 (1998).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM