

Product Information

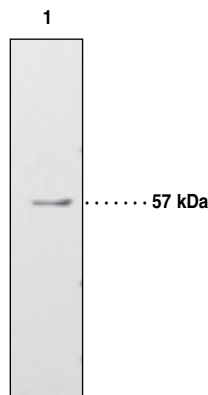


PPAR γ Polyclonal Antibody

Item No. 101700 • Lot No. XXXXX

- Contents:** This vial contains (100-500 μ g of peptide affinity-purified IgG, *lot specific*) in 500 μ l TBS, pH 8.0, containing 50% glycerol, 0.5 mg/ml BSA, and 0.02% sodium azide.
- Host:** Rabbit
- Antigen:** The PPAR γ polyclonal antibody was raised against a peptide corresponding to amino acids 82-101 of human PPAR γ 1 (PASPPYYSEKTQLYNKPHEE; amino acids 110-129 of PPAR γ 2).¹
- Cross-reactivity:** (+) Human and mouse PPAR γ 1 and PPAR γ 2; other species not tested but are expected to detect PPAR γ isoforms from:
- | | |
|----------------------------------|-----------------|
| canine, ape, and monkey | (100% identity) |
| porcine, guinea pig, and duck | (95% identity) |
| ovine, bovine, goat, and hamster | (90% identity) |
| chicken, goose, and crocodile | (85% identity) |
- Stability:** \geq 1 year at -20°C
- Application:** This polyclonal antibody can be used for detection of PPAR γ by western blot and immunohistochemistry at a dilution of 1:*lot specific*. Recommended for immunoprecipitation using 5 μ g IgG per reaction (PPAR γ Affinity Sorbent Item No. 401709, also available.)
- Concentration:** Varies by lot, from 0.2-1.0 mg/ml (100-500 μ g/vial). Always 100 ml final working volume for western blotting.

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



Lane 1: Rat adipose homogenate (30 μ g)

References

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

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/101700

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/12/2011

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com