

Product Information

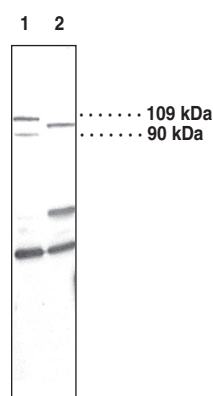


PGC-1 Polyclonal Antibody

Catalog No. 101707 • Lot No. XXXXXX

- Synonyms:** Peroxisome Proliferator-activated Receptor γ Coactivator 1, PPAR γ Coactivator 1
- Contents:** This vial contains XXX μ g peptide affinity-purified IgG in XXX μ l TBS, pH 7.4, containing 50% glycerol, 0.5 mg/ml, and 0.02% sodium azide
- Host:** Rabbit
- Antigen:** Human PGC-1 α amino acids 75-90; the antigen alignment with other known sequences is as follows:
Human PGC-1 α NIFEKIDEENEANLLA
Human PGC-1 β s e f q I D s E N E A 1 L a e
Murine and rat PGC-1 β s e f q I D s E N E A 1 L a A
- Cross-reactivity:** (+) Human, murine, rat PGC-1 α and PGC-1 β ; other species not tested
- Stability:** \geq 1 year at -20°C
- Applications:** The recommended starting dilution for western blot is 1:XXX and 2 μ g/ml for immunohistochemistry (paraffin-embedded tissue sections). Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Three peroxisome proliferator-activated receptor γ coactivator 1 (PGC-1) isoforms have been characterized - PGC-1 α , PGC-1 β , and PGC-1-related coactivator (PRC). PGC-1 α and PGC-1 β are products of two distinct genes with protein sequences sharing high homology at both the N- and C-terminus.¹ PGC-1 α is a tissue-specific and inducible transcriptional coactivator for several nuclear receptors and plays a key role in energy metabolism, hepatic gluconeogenesis, and cholesterol homeostasis.² PGC-1 β is also thought to activate oxidative metabolism in tissues, although it does so through a relatively restricted set of transcriptional partners compared to PGC-1 α . PGC-1 α is most abundant in kidney and heart whereas PGC-1 β is primarily expressed in heart, brown adipose, and skeletal muscle.^{3,4} Changes in PGC-1 α levels may play a role in metabolic disorders such as type II diabetes and obesity.⁵ Human PGC-1 α is 798 amino acids in length with a calculated molecular weight of 91 kDa. Cayman's PGC-1 polyclonal antibody detects the protein at around 96 kDa in murine kidney and liver samples. In human liver microsomes and HEK293 cells, both PGC-1 α and PGC-1 β are detected.



Lane 1: HEK293 cell lysate (40 μ g)
Lane 2: Murine kidney 100,000 x g pellet resuspension (40 μ g)

Laboratory Procedures

Immunoperoxidase staining procedure

A. Paraffin sections

1. Deparaffinize sections 3 times with xylene or xylene substitute, 5 minutes each.
2. Rehydrate sections 2 times with 100% ethanol, 5 minutes each, followed by 95%, 90%, 80%, 70%, and 50% ethanol, 5 minutes each.
3. Rinse sections in distilled water for 5 minutes.
4. Block endogenous peroxidase activity with 0.3% H₂O₂ in water (use methanol instead of water in case of strong endogenous peroxidase activity) for 15 minutes.
5. Wash sections 3 times in TBS containing 0.1% Tween 20 (TBST), pH 7.4, 5 minutes each.
6. Incubate sections with 5% normal serum (same species as the host of the secondary antibody) for 30 minutes.
7. Rinse sections 3 times with TBST, pH 7.4, 5 minutes each.
8. Incubate sections with a 1:300 dilution of PGC-1 polyclonal antibody (recommended starting dilution. Optimal dilution to be determined by end user) overnight at room temperature.
9. Wash sections 3 times in TBST, pH 7.4, 5 minutes each.

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, 03/06/2010

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

10. Incubate sections for 30 minutes with biotinylated secondary antibody, using a dilution as recommended by the provider.
 11. Wash sections 3 times in TBST, pH 7.4, 5 minutes each.
 12. Incubate sections for 30 minutes with ABC reagent, using a dilution as recommended by the provider.
 13. Wash sections 3 times in TBST, pH 7.4, 5 minutes each.
 14. Incubate sections in peroxidase substrate solution. Check staining under a microscope frequently. When desired staining intensity is achieved, rinse sections with distilled water thoroughly.
 15. Counter stain sections if desired. Rinse sections thoroughly after counter stain.
 16. Dehydrate sections through 50%, 70%, 80%, 90%, 95%, and 100% ethanol (2 times) for 5 minutes each.
 17. Clear sections 3 times with xylene or xylene substitute, 5 minutes.
 18. Mount sections with coverslips.
- B. Fresh Frozen Sections
1. After briefly fixing sections with an appropriate fixative (e.g., 10% formaldehyde for 2 minutes), wash sections 3 times with TBS, pH 7.4, 5 minutes each.
 2. Follow steps 4-18 of the procedure recommended for paraffin section.

References

1. Lin, J., Puigserver, P., Donovan, J., *et al.* Peroxisome proliferator-activated receptor γ coactivator 1 β (PGC-1 β), a novel PGC-1-related transcription coactivator associated with host cell factor. *J. Biol. Chem.* **277**(3), 1645-1648 (2002).
2. Lin, J., Handschin, C., and Spiegelman, B.M. Metabolic control through the PGC-1 family of transcription coactivators. *Cell Metabolism* **1**, 361-370 (2005).
3. Larrouy, D., Vidal, H., Andreelli, F., *et al.* Cloning and mRNA tissue distribution of human PPAR γ coactivator-1. *Int. J. Obes.* **23**, 1327-1332 (1999).
4. Meirhaeghe, A., Crowley, V., Lenaghan, C., *et al.* Characterization of the human, mouse, and rat PGC1 β (peroxisome-proliferator-activated receptor- γ co-activator 1 β) gene *in vitro* and *in vivo*. *J. Biochem.* **373**, 155-165 (2003).
5. Flier, J.S. Obesity wars: Molecular progress confronts an expanding epidemic. *Cell* **116**, 337-350 (2004).

Related Products

CD36 Polyclonal Antibody - Cat. No. 100011 • PPAR γ Polyclonal Antibody - Cat. No. 101700 • PPAR α Polyclonal Antibody - Cat. No. 101710 • PPAR δ Polyclonal Antibody - Cat. No. 101720 • MCAD Polyclonal Antiserum - Cat. No. 101730 • CD36 Monoclonal Antibody- Cat. No. 188150 • FABP4 Polyclonal Antibody - Cat. No. 10004944 • PPAR γ Transcription Factor Assay Kit - Cat. No. 10006855 • PPAR δ Transcription Assay Factor Kit - Cat. No. 10006914 • PPAR α Transcription Factor Assay Kit - Cat. No. 10006915

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