

# PRODUCT INFORMATION



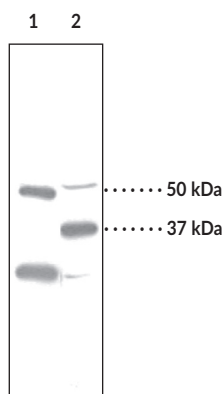
## PPAR $\delta$ Polyclonal Antibody

Item No. 101720

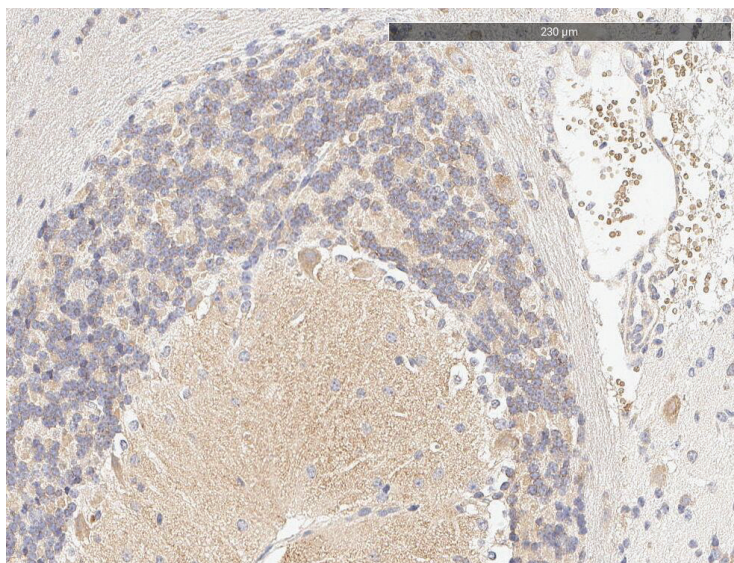
### Overview and Properties

<b>Contents:</b>	This vial contains 500 $\mu$ l peptide affinity-purified polyclonal antibody.
<b>Synonyms:</b>	FAAR, NUC1, Nuclear Hormone Receptor 1, PPAR $\beta$
<b>Immunogen:</b>	Synthetic peptide from the N-terminal region of human PPAR $\delta$
<b>Species Reactivity:</b>	(+) Human and mouse; other species not tested
<b>Uniprot No.:</b>	Q03181
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	$\geq 3$ years
<b>Storage Buffer:</b>	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
<b>Host:</b>	Rabbit
<b>Applications:</b>	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:200 for IHC and WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



Lane 1: Human cerebral cortex (30  $\mu$ g)  
Lane 2: Mouse liver (30  $\mu$ g)



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) mouse brain tissue after heat-induced antigen retrieval in pH 6.0 citrate buffer. After incubation with PPAR $\delta$  polyclonal antibody (Item No. 101720) at a 1:200 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.

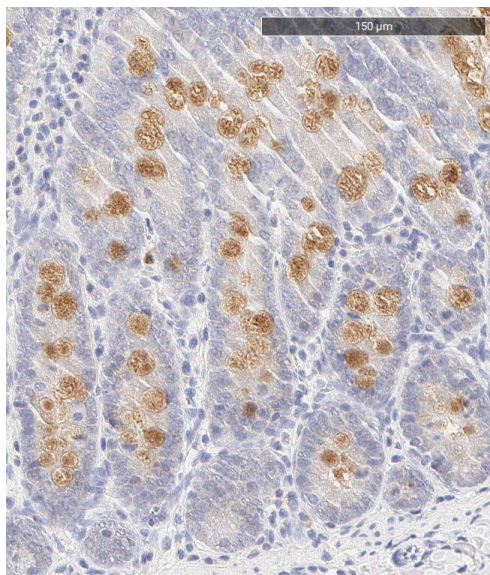
**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email 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, 11/02/2023

**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



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) rat colon tissue after heat-induced antigen retrieval in pH 6.0 citrate buffer. After incubation with PPAR $\delta$  polyclonal antibody (Item No. 101720) at a 1:200 dilution, slides were incubated with biotinylated secondary antibody, followed by alkaline phosphatase-streptavidin and chromogen (DAB).

## Description

PPAR $\delta$  is one of three peroxisome proliferator-activated receptor (PPAR) subtypes that possess a domain structure common to other members of the nuclear receptor gene family. It was first cloned from *Xenopus laevis* and named PPAR $\beta$ .<sup>1</sup> PPAR $\delta$  is ubiquitously expressed but is particularly abundant in tissues such as liver, intestine, kidney, abdominal adipose, and skeletal muscle, all of which are involved in lipid metabolism.<sup>2</sup> PPAR $\delta$  is a mediator of diverse physiological functions including lipid and cholesterol homeostasis, embryo implantation, and cancer development.<sup>3-6</sup> Most recently, attention has been focused on the role of PPAR $\delta$  in obesity.<sup>7</sup>

Cayman's PPAR $\delta$  Polyclonal Antibody can be used for WB and IHC to study the expression and functions of this protein. The antibody recognizes PPAR $\delta$  at 50 kDa from human samples. An additional smaller size of PPAR $\delta$  (~40 kDa)<sup>8</sup> is also detected in certain mouse tissues.

## References

1. Dreyer, C., Krey, G., Keller, H., *et al.* *Cell* **68**, 879-887 (1992).
2. Willson, T.M., Brown, P.J., Sternbach, D.D., *et al.* *J. Med. Chem.* **43**(4), 528-550 (2000).
3. Amri, E.-Z., Bonino, F., Ailhaud, G., *et al.* *J. Biol. Chem.* **270**, 2367-2371 (1995).
4. Berger, J., Leibowitz, M.D., Doebber, T.W., *et al.* *J. Biol. Chem.* **274**, 6718-6725 (1999).
5. Lim, H. and Dey, S.K. *TEM* **11**(4), 137-142 (2000).
6. He, T.-C., Chan, T.A., Vogelstein, B., *et al.* *Cell* **99**, 335-345 (1999).
7. Wang, Y.-X., Lee, C.-H., Tiep, S., *et al.* *Cell* **113**, 159-170 (2003).
8. Mammalian Gene Collection (MGC) Program Team *Proc. Natl. Acad. Sci. USA* **99**(26), 16899-16903 (2002).

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