

PRODUCT INFORMATION



PPAR δ (human, recombinant)

Item No. 10007451

Overview and Properties

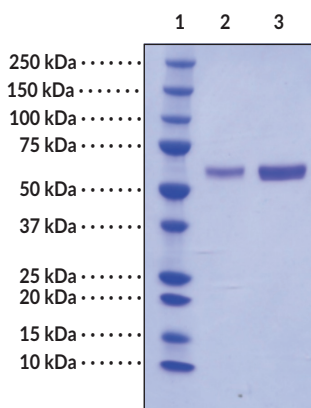
Synonyms: FAAR, NUC1, Nuclear Hormone Receptor 1, Peroxisome Proliferative Activated Receptor δ , PPAR β
Source: Recombinant protein isolated from a baculovirus overexpression system in Sf21 cells
Amino Acids: 2-441 (full length)
Molecular Weight: 54 kDa
Storage: -80°C (as supplied)
Stability: ≥ 1 year
Purity: *batch specific* ($\geq 95\%$ estimated by SDS-PAGE)
Supplied in: 50 mM of sodium phosphate, pH 7.2, containing 20% glycerol, 100 mM sodium chloride, and 1 mM DTT

Protein

Concentration: *batch specific* mg/ml

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



Lane 1: MW Markers
Lane 2: PPAR δ (2 μ g)
Lane 3: PPAR δ (4 μ g)

Representative gel image shown; actual purity may vary between each batch.

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, 12/02/2021

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

Peroxisome proliferator-activated receptors (PPARs) are members of the nuclear receptor family of ligand activated transcription factors that heterodimerize with retinoic acid like receptors, regulating gene expression and differentiation.¹ The PPAR family of nuclear hormone receptors consists of three subtypes encoded by separate genes: PPAR α , PPAR δ (also referred to as hNUC1, PPAR β , or FAAR), and PPAR γ . PPAR δ 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.² PPAR δ is a mediator of diverse physiological functions including lipid and cholesterol homeostasis, embryo implantation, and cancer development.³⁻⁶ Most recently, attention has been focused on the role of PPAR δ in obesity.⁷

Cayman's PPAR δ protein is purified from a baculovirus overexpression system. The purity was determined using gel electrophoresis followed by coomassie staining.

References

1. Kersten, S., Desvergne, B., and Wahli, W. Roles of PPARs in health and disease. *Nature* **405** (6785), 421-424 (2000).
2. Willson, T.M., Brown, P.J., Sternbach, D.D., *et al.* The PPARs: From orphan receptors to drug discovery. *J. Med. Chem.* **43**(4), 528-550 (2000).
3. Amri, E.-Z., Bonino, F., Ailhaud, G., *et al.* Cloning of a protein that mediates transcriptional effects of fatty acids in preadipocytes. Homology to peroxisome proliferator-activated receptors. *J. Biol. Chem.* **270**, 2367-2371 (1995).
4. Berger, J., Leibowitz, M.D., Doebber, T.W., *et al.* Novel peroxisome proliferator-activated receptor (PPAR) γ and PPAR δ ligands produce distinct biological effects. *J. Biol. Chem.* **274**, 6718-6725 (1999).
5. Lim, H. and Dey, S.K. PPAR δ functions as a prostacyclin receptor in blastocyst implantation. *Trends Endocrinol. Metab.* **11**(4), 137-142 (2000).
6. He, T.-C., Chan, T.A., Vogelstein, B., *et al.* PPAR δ is an APC-regulated target of nonsteroidal anti-inflammatory drugs. *Cell* **99**(3), 335-345 (1999).
7. Wang, Y.-X., Lee, C.-H., Tjep, S., *et al.* Peroxisome-proliferator-activated receptor δ activates fat metabolism to prevent obesity. *Cell* **113**(2), 159-170 (2003).

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