

# PRODUCT INFORMATION



## PPAR $\gamma$ FL (human recombinant from *E. coli*)

Item No. 61700 • Batch No. XXXXX

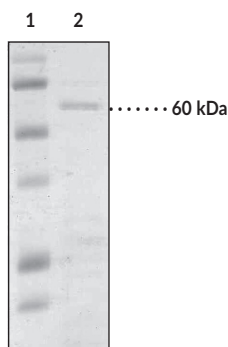
### Overview and Properties

**Synonym:** Peroxisome Proliferator-activated Receptor  $\gamma$  Full Length  
**Source:** Human recombinant N-terminal hexahistidine-tagged protein expressed in *E. coli*  
**Uniprot No.:** P37231

Batch specific information can be found on the Certificate of Analysis or by contacting Technical Support

**Molecular Weight:** ~60 kDa  
**Storage:** -80°C (as supplied)  
**Stability:** As supplied, 6 months from the QC date provided on the Certificate of Analysis, when stored properly  
**Purity:** *batch specific* (estimated by SDS-PAGE)  
**Supplied in:** 20 mM Tris HCl, pH 8.0, containing 250 mM KCl, 20% glycerol, 5 mM DTT, and 0.5 mM EDTA  
**Protein Concentration:** *batch specific* mg/ml

### Image



Lane 1: MW Standards  
Lane 2: PPAR $\gamma$  FL (2  $\mu$ 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, 07/11/2017

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 to regulate gene expression and differentiation.<sup>1</sup> The PPAR family of nuclear hormone receptors consists of three subtypes encoded by separate genes: PPAR $\alpha$ , PPAR $\delta$  (also referred to as hNUC1, PPAR $\beta$ , or FAAR), and PPAR $\gamma$ . Among them PPAR $\gamma$  is the most widely studied and has been implicated in the pathology of numerous diseases including obesity, diabetes, atherosclerosis, and cancer.<sup>2</sup> PPAR $\gamma$  is primarily expressed in adipose tissue and to a lesser extent in the colon, the immune system, and the retina.<sup>3</sup> PPAR $\gamma$  FL (human recombinant from *E. coli*) 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**, 421-424 (2000).
2. Vidal-Puig, A., Jimenez-Linan, M., Lowell, B.B., *et al.* Regulation of PPAR  $\gamma$  gene expression by nutrition and obesity in rodents. *J. Clin. Invest.* **97**, 2553-2561 (1996).
3. Clark, R.B. The role of PPARs in inflammation and immunity. *J. Leukoc. Biol.* **71**, 388-400 (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