

PRODUCT INFORMATION



PPAR α Ligand-binding Domain (human, recombinant)

Item No. 10009088

Overview and Properties

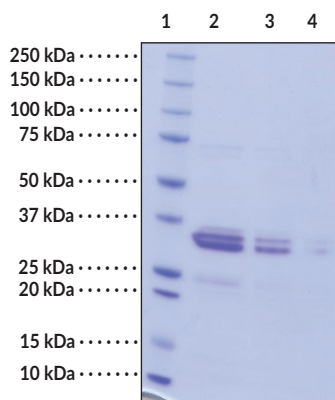
Synonyms: Peroxisome Proliferator-activated Receptor α Ligand Binding Domain, PPAR α LBD
Source: Recombinant human His-tagged PPAR α ligand-binding domain expressed in *E. coli*
Amino acids: 170-430
Uniprot No.: Q07869
Molecular Weight: ~34 kDa/subunit
Storage: -80°C (as supplied)
Stability: ≥ 1 year
Purity: *batch specific* ($\geq 55\%$ estimated by SDS-PAGE)
Supplied in: 50 mM sodium phosphate, pH 7.2, containing 20% glycerol and 100 mM sodium chloride

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 α LBD (human recombinant) (4 μ g)

Lane 3: PPAR α LBD (human recombinant) (2 μ g)

Lane 4: PPAR α LBD (human recombinant) (1 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, 02/16/2023

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



Description

Peroxisome proliferator-activated receptor α (PPAR α) is a member of the nuclear receptor family of ligand-activated transcription factors that regulates a variety of metabolic functions and inflammation.¹ It contains an N-terminal domain that is subject to phosphorylation, a DNA-binding domain, and a C-terminal ligand-binding domain.² PPAR α is highly expressed in tissues with high fatty acid oxidation rates, including the liver, heart, skeletal muscle, brown adipose tissue, and kidney, as well as in macrophages and T cells.^{2,3} It is activated by a variety of endogenous ligands such as fatty acids, eicosanoids, and endocannabinoids, as well as synthetic agents, including fenofibrate (Item No. 10005368) and gemfibrozil (Item No. 14835).⁴ Upon activation, PPAR α heterodimerizes with the retinoid X receptor (RXR) and binds to PPAR response elements in PPAR α target genes, recruiting RNA polymerase II and initiating gene transcription.¹ PPAR α transcriptionally regulates a variety of genes involved in several cellular processes, including lipid and hormone transport, peroxisomal and mitochondrial β -oxidation, amino acid metabolism, and inflammation.^{1,2} Genome-wide deletion of *Ppara* protects mice from high-fat diet-induced hyperinsulinemia and insulin resistance.⁵ PPARA SNPs have been found in individuals with a variety of cardiovascular conditions, including hypertension, atherosclerosis, coronary artery disease, left ventricular hypertrophy, or myocardial infarction.¹ Formulations containing PPAR α agonists have been used in the treatment of hyperlipidemia. Cayman's PPAR α Ligand-binding Domain protein can be used for Western blot (WB) applications.

References

1. Li, S., Yang, B., Du, Y., *et al.* Targeting PPAR α for the treatment and understanding of cardiovascular diseases. *Cell Physiol. Biochem.* **51(6)**, 2760-2775 (2018).
2. Pawlak, M., Lefebvre, P., and Staels, B. Molecular mechanism of PPAR α action and its impact on lipid metabolism, inflammation and fibrosis in non-alcoholic fatty liver disease. *J. Hepatol.* **62**, 720-733 (2015).
3. Rakhshandehroo, M., Knoch, B., Müller, M., *et al.* Peroxisome proliferator-activated receptor alpha target genes. *PPAR Res.* (2010).
4. Ruscica, M., Busnelli, M., Runfola, E., *et al.* Impact of PPAR-Alpha polymorphisms-the case of metabolic disorders and atherosclerosis. *Int. J. Mol. Sci.* **20(18)**, 4378 (2019).
5. Guerre-Millo, M., Rouault, C., Poulain, P., *et al.* PPAR- α -null mice are protected from high-fat diet-induced insulin resistance. *Diabetes* **50(12)**, 2809-2814 (2001).

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