

Recombinant PPAR γ protein

Catalog No: 81214, 81914

Expressed In: Baculovirus

Quantity: 20, 1000 μ g

Concentration: 0.4 μ g/ μ l

Source: Human

Buffer Contents: Recombinant PPAR γ protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, 0.5 mM TCEP.

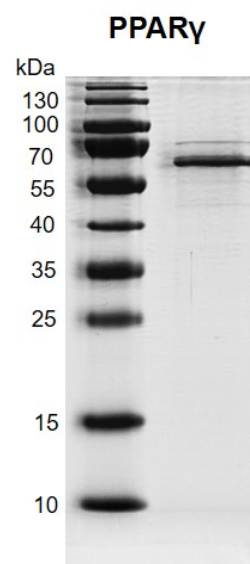
Background: PPAR γ (Peroxisome Proliferator Activated Receptor Gamma), also called as PPARG or NR1C3, is a member of the peroxisome proliferator-activated receptor (PPAR) subfamily of nuclear receptors. PPARs form heterodimers with retinoid X receptors (RXRs) and these heterodimers regulate transcription of various genes. Three subtypes of PPARs are known: PPAR-alpha, PPAR-delta, and PPAR-gamma.

PPAR γ is a nuclear receptor that binds peroxisome proliferators such as hypolipidemic drugs and fatty acids. Once activated by a ligand, it can bind to DNA specific PPAR response elements (PPRE) and modulate the transcription of its target genes, such as acyl-CoA oxidase. It therefore controls the peroxisomal beta-oxidation pathway of fatty acids. It's also a key regulator of adipocyte differentiation and glucose homeostasis. PPAR γ acts as a critical regulator of gut homeostasis by suppressing NF-kappa-B-mediated proinflammatory responses. It plays a role in the regulation of cardiovascular circadian rhythms by regulating the transcription of ARNTL/BMAL1 in the blood vessels (By similarity).

Protein Details: Recombinant PPAR γ protein was expressed in baculovirus expression system as the full length protein (accession number NP_056953.2) with an N-terminal FLAG tag. The molecular weight of the protein is 58.7 kDa.

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data this product is shown.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.



Recombinant PPAR γ protein gel
10% SDS-PAGE Coomassie staining
MW: 58.7 kDa
Purity: >85%