

PRODUCT INFORMATION



PPAR γ Blocking Peptide

Item No. 301700

Overview and Properties

Contents:	This vial contains 200 μ g of lyophilized peptide.
Synonym:	Peroxisome Proliferator-activated Receptor γ
Form:	Solid
Storage:	-20°C (as supplied)
Stability:	\geq 3 years

Procedures

This vial contains 200 μ g of lyophilized peptide derived from the human PPAR γ sequence. This peptide was used as an antigen for production of the PPAR γ Polyclonal Antibody (Item No. 101700) and can be used in conjunction with this antibody to block protein-antibody complex formation during immunochemical analysis for this protein

To block antibody/protein complex formation, the following procedure is recommended:

1. Mix the PPAR γ Polyclonal Antibody (Item No. 101700) and blocking peptide together in a 1:1 (v/v) ratio in a microfuge tube. For example, mix 20 μ l of antibody and 20 μ l of peptide.*
2. Incubate for one hour at room temperature with occasional mixing prior to further dilution and application of the mixture to the immunoblot.
3. Dilute the mixture to the final working antibody concentration and apply to the slide or membrane as usual.

*This is a recommended mixture. The minimum amount of peptide needed for complete blocking has not been precisely determined and may vary depending on the sample being analyzed. The amount of peptide required may need to be increased if sufficient blocking does not occur.

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, 12/07/2023

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