

PPARG polyclonal antibody

Other names: PPAR-gamma, NR1C3

Cat. No. C15310133

Type: Polyclonal

Source: Rabbit

Lot #: A576-001

Size: 100 µl

Concentration: not determined

Specificity: Human: positive

Other species: not tested

Purity: Whole antiserum from rabbit containing 0.05% azide.

Storage: Store at -20°C; for long storage, store at -80°C.

Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Description: Polyclonal antibody raised in rabbit against human PPARG (peroxisome proliferator-activated receptor gamma), using a KLH-conjugated synthetic peptide containing a sequence from the central part of the protein.

Applications

	Suggested dilution	Results
ELISA	1:500	Fig 1
Western blotting	1:500	Fig 2

Target description

PPARG (UniProtKB/Swiss-Prot entry P37231) is a nuclear hormone receptor which binds peroxisome proliferators such as hypolipidemic drugs and fatty acids. Like many other nuclear hormone receptors, PPARG forms a heterodimer with the retinoid X receptor (RXR) leading to transcriptional regulation of various genes including acyl-CoA oxidase and cytochrome P450 A6. PPARG has been implicated in adipocyte differentiation and glucose homeostasis and in various diseases such as obesity, diabetes, atherosclerosis and cancer.

Results

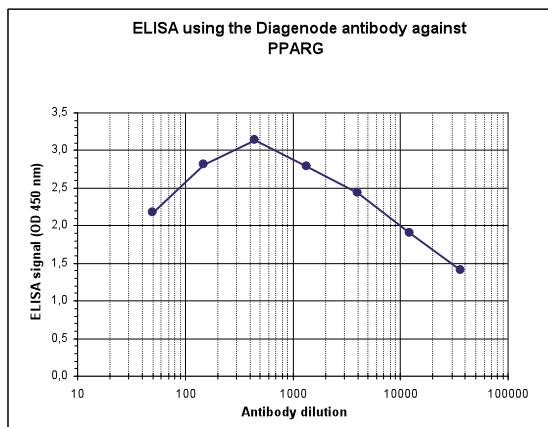


Figure 1. Determination of the titer

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against human PPARG (Cat. No. C15310133). The plates were coated with the peptide used for immunization of the rabbit. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:27,600.

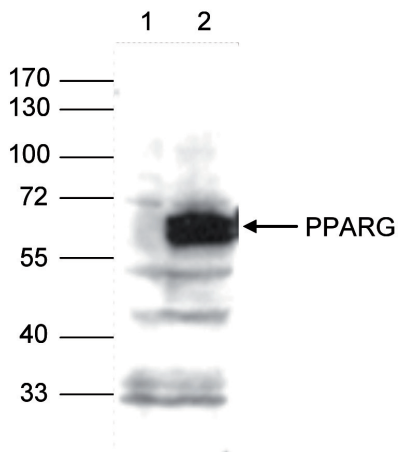


Figure 2. Western blot analysis using the Diagenode antibody directed against PPARG

293T cells were analysed by Western blot using the Diagenode antibody against PPARG (Cat. No. C15310133) diluted 1:500. Figure 2 shows the result of 293T cells transfected with pNTAP-PPARG (lane 2) or with an empty vector (lane 1). The position of the protein of interest is indicated on the right (expected size: 58 kDa); the marker (in kDa) is shown on the left.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue Bois Saint-Jean, 3
4102 Seraing (Ougrée) - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: March 3, 2015