


PPAR gamma Transcription Factor Assay Kit ab133101

★★★★★ 3 Abreviews 21 References 4 Images

Overview

Product name	PPAR gamma Transcription Factor Assay Kit
Detection method	Colorimetric
Sample type	Adherent cells, Suspension cells, Nuclear Extracts
Assay type	Semi-quantitative
Species reactivity	Reacts with: Mouse, Rat, Human

Product overview

Predicted to work with: Mammals 

PPAR gamma Transcription Factor Assay kit ab133101 is a non-radioactive, sensitive method for detecting specific transcription factor DNA binding activity in nuclear extracts.

A 96 well enzyme-linked immunosorbent assay (ELISA) replaces the cumbersome radioactive electrophoretic mobility shift assay (EMSA). A specific double stranded DNA (dsDNA) sequence containing the peroxisome proliferator response element (PPRE) is immobilized onto the bottom of wells of a 96 well plate. PPARs contained in a nuclear extract bind specifically to the PPRE. PPAR gamma is detected by addition of specific primary antibody directed against PPAR gamma. A secondary antibody conjugated to HRP is added to provide a sensitive colorimetric readout at 450 nm. PPAR gamma will not crossreact with PPAR delta or PPAR alpha.

Notes

Peroxisome proliferator-activated receptors (PPARs) are ligand activated nuclear receptors. Three PPAR subtypes have been identified: alpha, delta and gamma. PPARs can be activated by polyunsaturated fatty acids, eicosanoids and various synthtic ligands.

PPAR gamma is primarily expressed in adipose tissue and to a lesser extent in the colon, immune system and the retina. PPAR gamma was first identified as regulator of adipogenesis, but also plays an important role in cellular differentiation, insulin sensitization , atherosclerosis and cancer.

Platform	Microplate reader
----------	-------------------

Properties

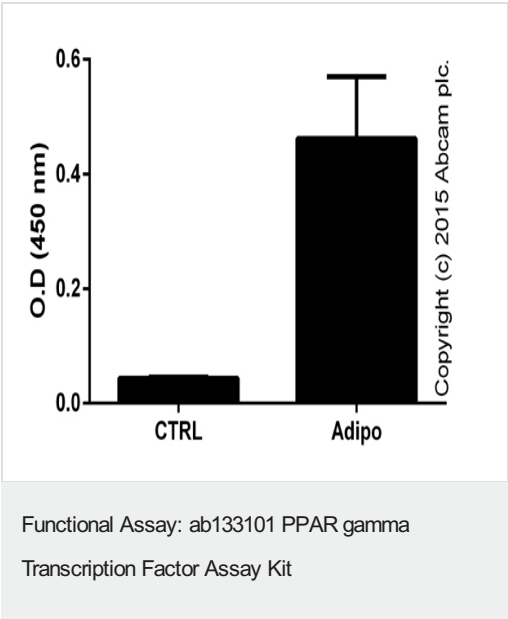
Storage instructions	Please refer to protocols.
----------------------	----------------------------

Components	96 tests
96-Well Plate Cover	1 unit

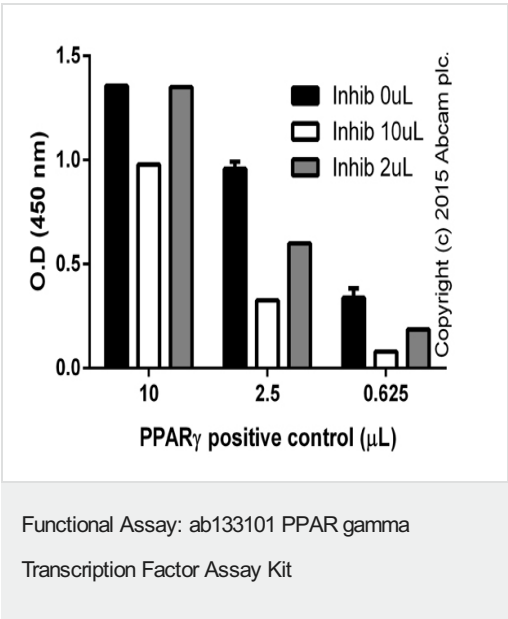
Components	96 tests
Polysorbate 20	1 vial
Transcription Factor Antibody Binding Buffer (10X)	1 x 3ml
Transcription Factor Binding Assay Buffer (4X)	1 x 3ml
Transcription Factor Developing Solution	1 x 12ml
Transcription Factor Goat Anti-Rabbit HRP Conjugate	1 x 100µl
Transcription Factor PPAR 96-Well Strip Plate	1 unit
Transcription Factor PPAR Competitor dsDNA	1 vial
Transcription Factor PPAR gamma Positive Control	1 vial
Transcription Factor PPAR gamma Primary Antibody	1 vial
Transcription Factor Reagent A	1 x 120µl
Transcription Factor Stop Solution	1 x 12ml
Wash Buffer Concentrate (400X)	1 x 5ml

Function	Receptor that binds peroxisome proliferators such as hypolipidemic drugs and fatty acids. Once activated by a ligand, the receptor binds to a promoter element in the gene for acyl-CoA oxidase and activates its transcription. It therefore controls the peroxisomal beta-oxidation pathway of fatty acids. Key regulator of adipocyte differentiation and glucose homeostasis.
Tissue specificity	Highest expression in adipose tissue. Lower in skeletal muscle, spleen, heart and liver. Also detectable in placenta, lung and ovary.
Involvement in disease	<p>Note=Defects in PPARG can lead to type 2 insulin-resistant diabetes and hypertension. PPARG mutations may be associated with colon cancer.</p> <p>Defects in PPARG may be associated with susceptibility to obesity (OBESITY) [MIM:601665]. It is a condition characterized by an increase of body weight beyond the limitation of skeletal and physical requirements, as the result of excessive accumulation of body fat.</p> <p>Defects in PPARG are the cause of familial partial lipodystrophy type 3 (FPLD3) [MIM:604367]. Familial partial lipodystrophies (FPLD) are a heterogeneous group of genetic disorders characterized by marked loss of subcutaneous (sc) fat from the extremities. Affected individuals show an increased preponderance of insulin resistance, diabetes mellitus and dyslipidemia.</p> <p>Genetic variations in PPARG can be associated with susceptibility to glioma type 1 (GLM1) [MIM:137800]. Gliomas are central nervous system neoplasms derived from glial cells and comprise astrocytomas, glioblastoma multiforme, oligodendrogliomas, and ependymomas.</p> <p>Note=Polymorphic PPARG alleles have been found to be significantly over-represented among a cohort of American patients with sporadic glioblastoma multiforme suggesting a possible contribution to disease susceptibility.</p>
Sequence similarities	<p>Belongs to the nuclear hormone receptor family. NR1 subfamily.</p> <p>Contains 1 nuclear receptor DNA-binding domain.</p>

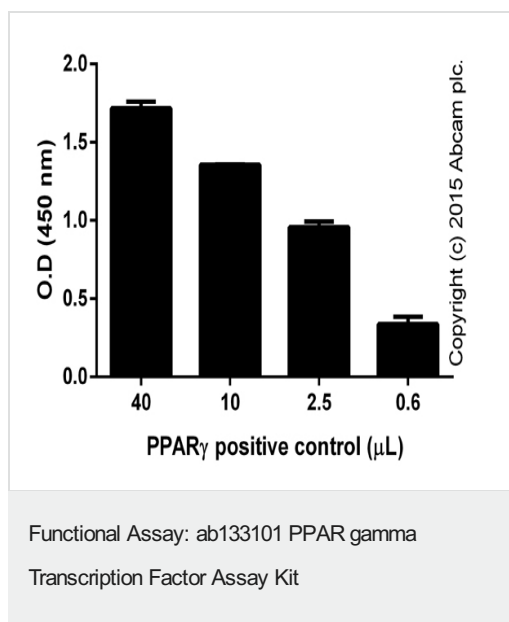
Images



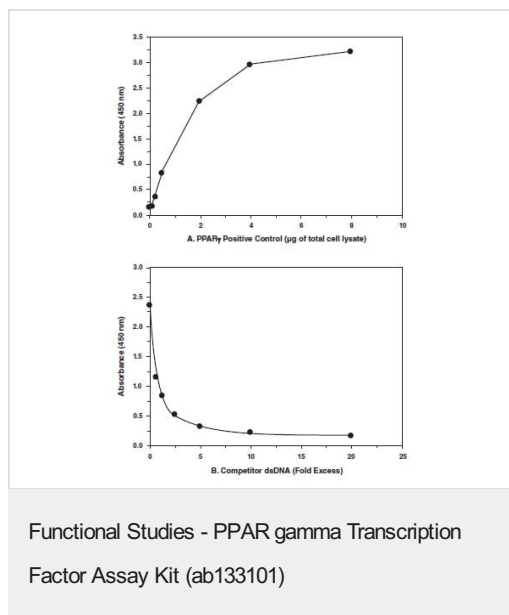
3T3-L1 cells were differentiated to adipocytes with 1 uM dexamethasone ([ab120743](#)), 1 ug x mL⁻¹ insulin ([ab123768](#)) and 0.5 mM IBMX ([ab120840](#)). From the third day, the cells were grown in normal medium with the addition of only 1 ug x mL⁻¹ insulin. From day six, the cells were cultured in normal medium for an additional two days. 40 uL of nuclear extracts were tested in duplicates (+/- SD).



Different volumes of positive control with inhibitor (duplicates, +/- SD).



Different volumes of positive control (duplicates, +/-SD).



Panel A: Increasing amounts of positive control (total lysate) are assayed for PPAR gamma DNA-binding activity using ab133101.

Panel B: PPAR gamma DNA-binding assays are performed in the presence of competitive dsDNA. The decrease in signal caused by addition of competitive dsDNA confirms the assay specificity.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors