


Product datasheet

PPAR (alpha, delta, gamma) Transcription Factor Assay Kit ab133113

[4 References](#) [3 Images](#)

Overview

Product name	PPAR (alpha, delta, gamma) Transcription Factor Assay Kit
Detection method	Colorimetric
Sample type	Adherent cells, Suspension cells, Nuclear Extracts
Assay type	Semi-quantitative
Species reactivity	Reacts with: Human Predicted to work with: Mammals 

Product overview Abcam's PPAR Transcription Factor Assay Kit (ab133113) is a non-radioactive, sensitive method for detecting specific transcription factor DNA binding activity in nuclear extracts and whole cell lysates.

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 PPAR response element is immobilized onto the bottom of wells of a 96-well plate. PPARs contained in a nuclear extract, bind specifically to the PPAR response element. PPAR alpha, delta, or gamma are detected by addition of specific primary antibodies directed against the individual PPARs. A secondary antibody conjugated to HRP is added to provide a sensitive colorimetric readout at 450 nm. ab133113 comes with a single plate that measures all three isoforms of PPAR alpha, delta, and gamma. There are enough reagents for one-third of a plate for each isoform.

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 synthetic ligands. PPARs play pivotal roles in the regulation of lipid metabolism and homeostasis and are important indirect and direct regulators of cellular insulin sensitivity.

Platform Microplate reader

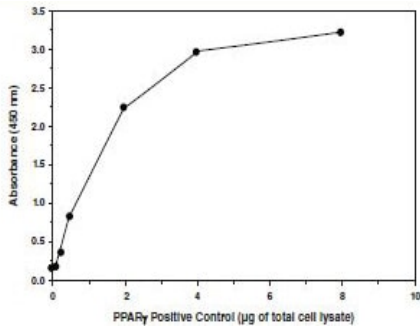
Properties

Storage instructions Please refer to protocols.

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

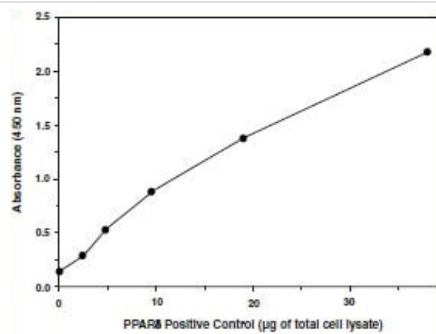
Cellular localization Cytoplasm, nucleus.

Images



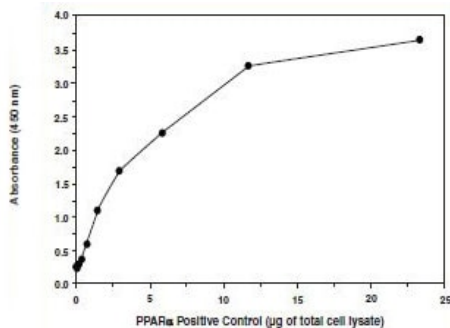
Increasing amounts of positive control (total lysate) are assayed for PPAR gamma DNA-binding activity using ab133113.

Transcription Factor Binding Assay for PPAR gamma



Increasing amounts of positive control (total lysate) are assayed for PPAR delta DNA-binding activity using ab133113.

Transcription Factor Binding Assay for PPAR delta



Increasing amounts of positive control (total lysate) are assayed for PPAR alpha DNA-binding activity using ab133113.

Transcription Factor Binding Assay for PPAR alpha

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