

PPAR delta Transcription Factor Assay Kit ab133106

[3 References](#) [1 Image](#)

Overview

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

Predicted to work with: Mammals 

Product overview

Abcam's PPAR delta Transcription Factor Assay Kit (ab133106) 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 delta is detected by addition of specific primary antibody directed against PPAR delta. A secondary antibody conjugated to HRP is added to provide a sensitive colorimetric readout at 450 nm. ab133106 detects human PPAR delta.

This assay does not cross-react with PPAR gamma 1 and PPAR gamma 2 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 synthetic ligands.

PPAR delta is the least understood; however, identification of a high-affinity PPAR delta synthetic ligand suggests a potential role for this receptor in lipid metabolism and obesity. PPAR delta is ubiquitously expressed, but it is particularly abundant in tissues such as liver, intestine, kidney, abdominal adipose, and skeletal muscle.

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

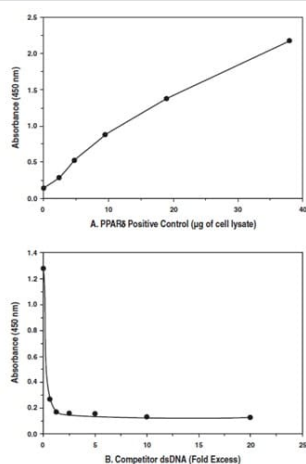
Properties

| | |
|-----------------------------|----------------------------|
| 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 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 PPARδ Positive Control | 1 vial |
| Transcription Factor PPARδ 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 | Ligand-activated transcription factor. Receptor that binds peroxisome proliferators such as hypolipidemic drugs and fatty acids. Has a preference for poly-unsaturated fatty acids, such as gamma-linoleic acid and eicosapentanoic acid. Once activated by a ligand, the receptor binds to promoter elements of target genes. Regulates the peroxisomal beta-oxidation pathway of fatty acids. Functions as transcription activator for the acyl-CoA oxidase gene. Decreases expression of NPC1L1 once activated by a ligand. |
| Tissue specificity | Ubiquitous with maximal levels in placenta and skeletal muscle. |
| Sequence similarities | Belongs to the nuclear hormone receptor family. NR1 subfamily. Contains 1 nuclear receptor DNA-binding domain. |
| Cellular localization | Nucleus. |

Images



Panel A: Increasing amounts of positive control (total lysate) are assayed for PPAR delta DNA-binding activity using ab133106.

Panel B: PPAR delta 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.

Functional Studies - PPAR delta Transcription
Factor Assay Kit (ab133106)

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