

PPAR delta Transcription Factor Assay Kit ab133106

[3 References](#) [1 图像](#)

概述

产品名称	PPAR delta Transcription Factor Assay试剂盒
检测方法	Colorimetric
样品类型	Adherent cells, Suspension cells, Nuclear Extracts
检测类型	Semi-quantitative
种属反应性	与反应: Mouse, Human

预测可用于: Mammals 

产品概述

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.

说明

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.

平台

Microplate reader

性能

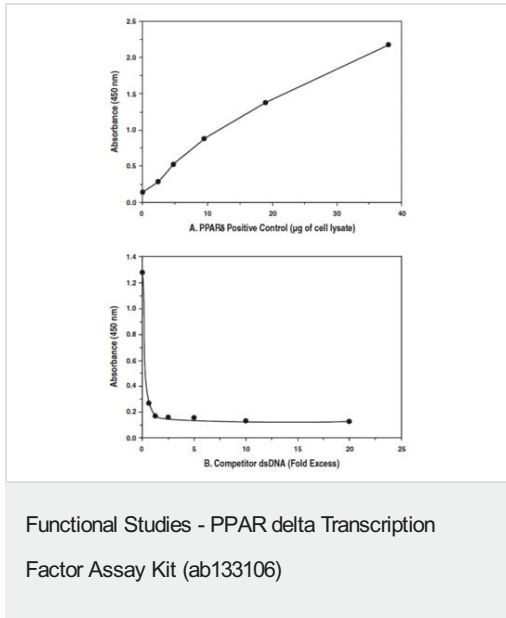
存放说明

Please refer to protocols.

组件	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 PPAR δ Positive Control	1 vial
Transcription Factor PPAR δ 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

功能	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.
组织特异性	Ubiquitous with maximal levels in placenta and skeletal muscle.
序列相似性	Belongs to the nuclear hormone receptor family. NR1 subfamily. Contains 1 nuclear receptor DNA-binding domain.
细胞定位	Nucleus.

图片



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.

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.cn/abpromise> or contact our technical team.

Terms and conditions

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