

PPAR alpha Transcription Factor Assay Kit ab133107

5 References 1 图像

概述

产品名称	PPAR alpha 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, Rat, Human
产品概述	<p>Abcam's PPAR alpha Transcription Factor Assay Kit (ab133107) is a non-radioactive, sensitive method for detecting specific transcription factor DNA binding activity in nuclear extracts.</p> <p>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 alpha is detected by addition of specific primary antibody directed against PPAR alpha. A secondary antibody conjugated to HRP is added to provide a sensitive colorimetric readout at 450 nm. PPAR alpha Transcription Factor Assay will not cross-react with PPAR gamma or PPAR delta.</p>
说明	<p>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.</p> <p>PPAR alpha is predominantly expressed in liver, heart and kidneys. Upon activation by ligands such as free fatty acids, lipids and other small molecules, PPAR alpha forms a heterodimeric complex with RXR alpha, and this complex will bind to response elements allowing transcription of target genes involved in lipid metabolism.</p>
平台	Microplate reader

性能

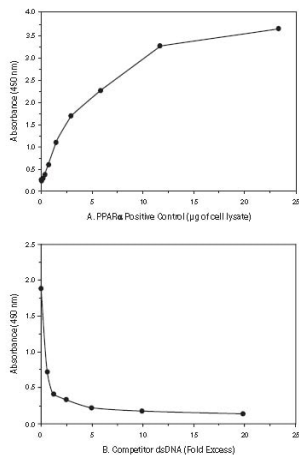
存放说明 Please refer to protocols.

组件	96 tests
96-Well Plate Cover	1 unit

组件	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 alpha Positive Control	1 vial
Transcription Factor PPAR alpha Primary Antibody	1 vial
Transcription Factor PPAR Competitor dsDNA	1 vial
Transcription Factor Reagent A	1 x 120µl
Transcription Factor Stop Solution	1 x 12ml
Wash Buffer Concentrate (400X)	1 x 5ml

<b>功能</b>	Ligand-activated transcription factor. Key regulator of lipid metabolism. Activated by the endogenous ligand 1-palmitoyl-2-oleoyl-sn-glycerol-3-phosphocholine (16:0/18:1-GPC). Activated by oleylethanolamide, a naturally occurring lipid that regulates satiety (By similarity). Receptor for peroxisome proliferators such as hypolipidemic drugs and fatty acids. Regulates the peroxisomal beta-oxidation pathway of fatty acids. Functions as transcription activator for the ACOX1 and P450 genes. Transactivation activity requires heterodimerization with RXRA and is antagonized by NR2C2.
<b>组织特异性</b>	Skeletal muscle, liver, heart and kidney.
<b>序列相似性</b>	Belongs to the nuclear hormone receptor family. NR1 subfamily. Contains 1 nuclear receptor DNA-binding domain.
<b>细胞定位</b>	Nucleus.

**图片**



Panel A: Increasing amounts of positive control (total lysate) are assayed for PPAR alpha DNA-binding activity using ab133107.

Panel B: PPAR alpha 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 alpha Transcription  
Factor Assay Kit (ab133107)

**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