
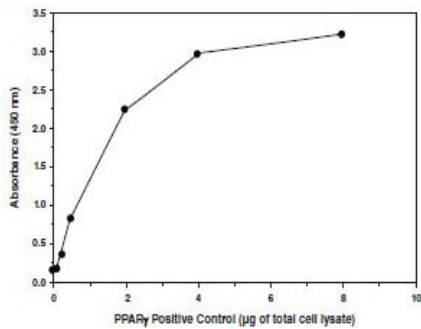


PPAR (alpha, delta, gamma) Transcription Factor Assay Kit ab133113

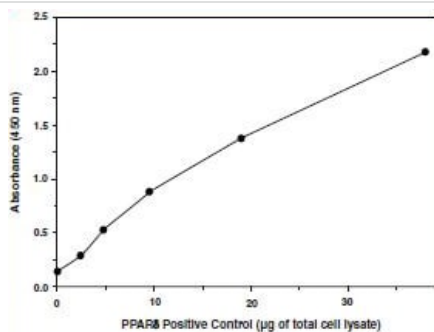
4 References 3 图像

概述	
产品名称	PPAR (alpha, delta, gamma) 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: Human 预测可用于: Mammals 
产品概述	<p>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.</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 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.</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. PPARs play pivotal roles in the regulation of lipid metabolism and homeostasis and are important indirect and direct regulators of cellular insulin sensitivity.</p>
平台	Microplate reader
性能	
存放说明	Please refer to protocols.



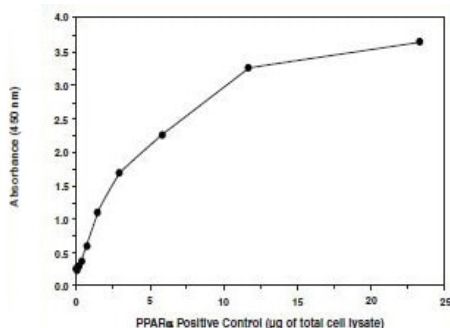
Transcription Factor Binding Assay for PPAR
gamma

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



Transcription Factor Binding Assay for PPAR delta

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



Transcription Factor Binding Assay for PPAR alpha

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

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