

Human PPARr-Regulated cDNA Plate Array

Catalog Number AP-0201

(For Research Use Only)

Introduction

The peroxisome-proliferator activated receptor gamma (PPARr) is a member of the nuclear receptor superfamily of transcription factors. PPARr is important to a variety of biological processes, such as adipocyte differentiation, glucose homeostasis, lipid trafficking as well as vascular function and hypertension. PPARr exerts its effects by regulating target gene transcription in a ligand-dependent manner. Upon activation, the PPAR γ -RXR heterodimer binds to specific DNA sequences on the target gene promoter and leads to transcriptional regulation. Signosis developed a plate-based array for profiling 20+ PPARr-regulated genes.

Principle of the assay

Signosis' proprietary cDNA plate array is a plate-based hybridization profiling analysis for monitoring the expression of dozens of genes through reverse transcription of mRNA into cDNA. Like array analyses, total RNA is first reverse transcribed into cDNA in the presence of biotin-dUTP in the assay. Targeted genes are then specifically captured onto individual wells on a plate, instead of membranes, through a pre-coated gene-specific oligonucleotide. The captured cDNAs are further detected with streptavidin-HRP. Luminescence is reported as relative light units (RLUs) on a microplate luminometer. The expression level of genes is directly proportional to the luminescent intensity.

Materials provided with the kit

- A 96-well plate coated with 31+ different capture oligos (RT)
- Human RRARr Reg. Primer Mix (-20 °C)
- Reverse transcription buffer mix (-20 °C)
- Reverse transcriptase RT (-20 °C)
- Streptavidin-HRP conjugate (4°C)
- Plate hybridization buffer (RT)
- 5x Plate hybridization wash buffer (RT)
- Blocking buffer (RT)
- 5xDetection wash buffer (RT)
- Substrate A (4°C)
- Substrate B (4°C)
- Substrate dilution buffer (4°C)

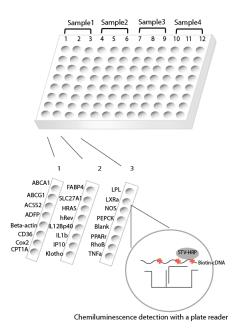


Diagram of human PPARr-regulated cDNA plate array

Material required but not provided

PCR machine Incubator 0.2ml PCR tube luminometer plate reader ddH2O (RNAase free)

Reagent preparation before starting experiment

- Dilute 30ml of 5x Plate hybridization wash buffer with 120 ml of dH₂O before use.
- Dilute 40ml of 5x Detection wash buffer with 160 ml of dH₂O before use.
- Warm up Plate hybridization buffer and Hybridization Wash buffer at 45 °C until no visible precipitate before use. Stir the solution with 10ml or 5ml pipette to facilitate the dissolving process.
- Dilute 500 times of streptavidin-HRP with blocking buffer before use.

Assay procedure

1. cDNA synthesis using PCR machine Note: Briefly spin tubes before opening

Sample preparation
X μl 1-10μg total RNA
μl Human PPARr Reg. Primer Mix
X μl ddH2O

11µl

- (2) Incubate for 5 minutes at 65 °C, and chill on ice.
- (3) Add 8 μl Reverse transcription buffer mix and 1μl RT to each reaction tube, and incubate for 1 hour at 45 °C.
- (4) Heat the reaction to 98 °C for 5 minutes, and chill on ice.
- (5) The 20ul cDNA is synthesized and labeled with biotin and ready for hybridization on the plate.

2. Plate hybridization

- (1) Remove the sealing film
- (2) Arrange the appropriate number of the wells of the plate based on your experiment. The whole plate is divided into 4 repeat sections, Column 1-3, 4-6, 7-9, 10-12 for 4 different samples.
- (3) Mix 20ul cDNA with 2.8ml pre-warmed Plate hybridization buffer, and dispense 95ul mixture to each well in a section **immediately**. A reagent reservoir can be used for dispensing cDNA mixture into the wells with a 8 multichannel pipette. Add 100ul Plate hybridization buffer without cDNA to the 'blank' well.
- (4) Seal the whole plate with foil film (provided) securely and incubate the plate at 45 °C for overnight. Ensure the numbers and letters on the plate are clearly visible from under foil seal by pressing the foil down on every single experimental well.

3. Plate detection

- (12) Remove the foil film from the experimental wells with a blade. Keep the unused well sealed.
- (13) Invert the Hybridization Plate over an appropriate container and expel the contents forcibly, and wash the plate 3 times by adding 200ul of pre-warmed 1x Plate hybridization wash buffer to each well. At each wash, incubate the plate for 5 minutes with gently shaking at room temperature.

- (3) Complete removal of liquid at each wash by firmly tapping the plate against clean paper towels.
- (4) Add 200μl of Blocking buffer to each well and incubate for 15 minutes at room temperature with gently shaking.
- (5) Invert the plate over an appropriate container to remove block buffer.
- (6) Add 20 μl of streptavidin-HRP conjugate in 10ml blocking buffer (1:500) dilution, enough for one plate. Add 95 μl of diluted streptavidin-HRP conjugate to each well and incubate for 45 min at room temperature with gently shaking.
- (7) Wash the plate 3 times by adding 200ul 1X Detection wash buffer to each well, At each wash, incubate the plate for 10 minutes with gently shaking at room temperature.
- (8) Complete removal of liquid at each wash by firmly tapping the plate against clean paper towels. At the last wash, invert plate on clean paper towels for 1-2 min to remove excessive liquid.
- (9) Freshly prepare the substrate solution:

For the whole plate:

1ml Substrate A

1ml Substrate B

8ml Substrate dilution buffer

- (10) Add 95 μ l substrate solution to each well and incubate for 1 min.
- (11) Place the plate in the luminometer. Allow plate to sit inside machine for 5 min before reading. Set integration time to 1 second with no filter position. For the best results, read the plate within 5-20 minutes.

Diagram of human PPAR-regulated cDNA plate array

	1	2	3	4	5	6	7	8	9	10	11	12
Α	ABCA1	FABP4	LPL	ABCA1	FABP4	LPL	ABCA1	FABP4	LPL	ABCA1	FABP4	LPL
В	ABCG1	FATP(SLC27A1)	LXRa	ABCG1	FATP(SLC27A1)	LXRa	ABCG1	FATP(SLC27A1)	LXRa	ABCG1	FATP(SLC27A1)	LXRa
С	ACSS2	HRASLS3	NOS2	ACSS2	HRAS	NOS	ACSS2	HRAS	NOS	ACSS2	HRAS	NOS
D	ADFP	hRev	PEPCK	ADFP	hRev	PEPCK	ADFP	hRev	PEPCK	ADFP	hRev	PEPCK
Е	Beta-actin	IL12B p40	Blank	B-actin	IL12B p40	Blank	B-actin	IL12B p40	Blank	B-actin	IL12B p40	Blank
F	CD36	IL-1b	PPARr	CD36	IL-1b	PPARr	CD36	IL-1b	PPARr	CD36	IL-1b	PPARr
G	Cox2	IP-10	RhoB	Cox2	IP-10	RhoB	Cox2	IP-10	RhoB	Cox2	IP-10	RhoB
Н	CPT1A	Klotho	TNFa	CPT1A	Klotho	TNFa	CPT1A	Klotho	TNFa	CPT1A	Klotho	TNFa