



Mouse HIF-Regulated cDNA Plate Array

Catalog Number AP-1111

(For Research Use Only)

Introduction

Rapidly growing tumors result in hypoxic regions. Adaptive responses of most cells to hypoxia are (1) to produce VEGF and other hypoxia-induced angiogenic cytokines that promote increased tissue vascularization, thereby increasing tissue oxygenation, and (2) to switch metabolically from oxidative phosphorylation to anaerobic glycolysis. Hypoxia-inducible factor 1 (HIF-1) is an oxygen-regulated transcriptional activator that plays essential roles in the process. The HIF-1 α subunit is oxygen-dependent ubiquitination and proteasomal degradation. In addition, cytokines including interleukin-1 (IL-1) and tumor necrosis factor- α (TNF- α) stimulate HIF-1 dependent gene expression. HIF-1 increases the expression of several genes that promote blood flow and inflammation, such as vascular endothelial growth factor (VEGF). Signosis' HIF cDNA Plate Array allows profiling 20+ related 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-captured 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 23 different capture oligos (RT)
- Mouse HIF 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)
- Block buffer (RT)
- 5xDetection wash buffer (RT)
- Substrate A (4°C)
- Substrate B (4°C)
- Substrate dilution buffer (RT)

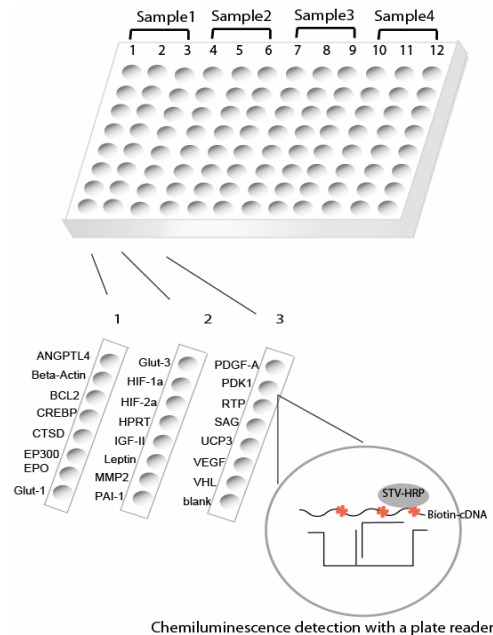


Diagram of Mouse HIF-regulated cDNA plate array

Material required but not provided

- PCR machine
- Incubator
- 0.2ml PCR tube
- Plate reader for chemiluminescence detection
- 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 before use.
- Dilute 1000 times of streptavidin-HRP with block buffer before use at Step 3(3).

Assay procedure

1. cDNA synthesis

Note: Briefly spin tubes before opening

(1) Sample preparation

- X µl 1-10µg total RNA
- 2 µl Mouse HIF 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, ready for hybridization on the plate.

2. Plate hybridization

- (1) Arrange the appropriate number of the wells of the plate based on your experiment by removing the top foil sealing film with a blade. Keep the unused well sealed. The whole plate is divided into 4 repeat sections, Column 1-3, 4-6, 7-9, 10-12 for 4 different samples.

Note: keep the removed sealing film for resealing during hybridization

- (2) 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.
- (3) Seal the wells with foil film 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.

3. Plate detection

- (1) Invert the plate over an appropriate container and expel the contents forcibly, and wash the plate by adding 300µl of warmed 1x Plate hybridization wash buffer. Repeat the washing process two times for a total of three washes. Complete removal of liquid at each wash by firmly tapping the plate against clean paper towels.
Optional: To ensure remove all traces of the buffer in last washing step, invert the plate on a clean, dry paper towel and centrifuge at 240xg for 1 minute at room temperature

- (2) Add 200µl of Block buffer incubate for 15 minutes at room temperature with gentle shaking.
- (3) Invert the plate over an appropriate container to remove block buffer. Add 100 µl of diluted streptavidin-HRP conjugate to each well and incubate for 45 min at room temperature with gentle shaking.
- (4) Wash the plate with 200ul 1X Detection wash buffer for 5 min at room temperature with gently shaking. Complete removal of liquid at each wash by firmly tapping the plate against clean paper towels.
- (5) Repeat Step (4) for additional 2 time washes.
- (6) Freshly prepare the substrate solution
For the whole plate:
1ml Substrate A
1ml Substrate B
8ml Substrate dilution buffer
- (7) Add 95µl substrate solution to each well and incubate for 1 minutes.
- (8) Determine the chemiluminescence of each well with a microplate reader within 5 minutes.

Example of Data Analysis

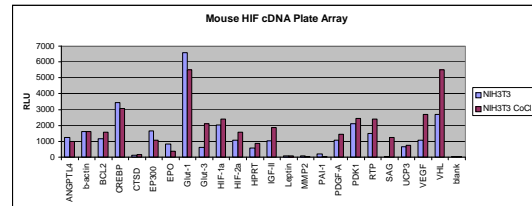


Figure 1: Mouse HIF-regulated cDNA plate assay for analysis of 21 genes.

NIH3T3 cells were treated with and without 200 uM cobalt chloride for 24 hours, RNAs prepared, cDNA synthesized with biotin label and subjected to cDNA plate array hybridization and detection.

Diagram of Mouse HIF-regulated cDNA plate array

	1	2	3	4	5	6	7	8	9	10	11	12
A	ANGPTL4	Glut-3	PDGF-A	ANGPTL4	Glut-3	PDGF-A	ANGPTL4	Glut-3	PDGF-A	ANGPTL4	Glut-3	PDGF-A
B	Beta-Actin	HIF-1a	PDK1	Beta-Actin	HIF-1a	PDK1	Beta-Actin	HIF-1a	PDK1	Beta-Actin	HIF-1a	PDK1
C	BCL2	HIF-2a	RTP	BCL2	HIF-2a	RTP	BCL2	HIF-2a	RTP	BCL2	HIF-2a	RTP
D	CREBP	HPRT	SAG	CREBP	HPRT	SAG	CREBP	HPRT	SAG	CREBP	HPRT	SAG
E	CTSD	IGF-II	UCP3	CTSD	IGF-II	UCP3	CTSD	IGF-II	UCP3	CTSD	IGF-II	UCP3
F	EP300	Leptin	VEGF	EP300	Leptin	VEGF	EP300	Leptin	VEGF	EP300	Leptin	VEGF
G	EPO	MMP2	VHL	EPO	MMP2	VHL	EPO	MMP2	VHL	EPO	MMP2	VHL
H	Glut-1	PAI-1	Blank	Glut-1	PAI-1	Blank	Glut-1	PAI-1	Blank	Glut-1	PAI-1	Blank