



## Cancer MicroRNA Array

Catalog Number AP-0003

(For Research Use Only)

### Introduction

Newly discovered microRNAs (miRNAs) are important to the regulation of gene expression (1). Up to 30% of mammalian genes might be regulated by miRNAs. So far, more than 400 miRNAs have been identified in the human genome and many of them are different only in one or a few nucleotides. Expression of mature miRNAs is tissue-specific (2) and the abundance of miRNAs varies in several orders of magnitude (3). More importantly, mis-regulation of miRNA expression might contribute to human cancers (4). Systematic profiling of miRNA expression displays unique signatures in a number of cancers (5).

Based on its proprietary technology, Signosis has developed a highly sensitive and discriminative array technology for monitoring miRNA. It combines oligo-ligation assay-based detection and T7 transcription-based linear amplification, providing a highly sensitive and specific array assay. Via ligation, selection, and transcription of a pair of oligos that are hybridized to each specific miRNA, the targets are amplified and array monitored. By assigning unique tag sequences to individual isoforms, they can be easily differentiated. The whole procedure is simple and straightforward. We are currently offering an array that targets the 60 most well studied miRNAs, each with cited literature, which is able to facilitate the comparison and discovery of differentiated miRNA expression in different samples.

### Principles of the technology

miRNAs are different from large messenger RNAs in three aspects; (1) miRNAs are small size molecules with quite a big difference in abundance, (2) mature miRNAs co-exist with their precursor pre-miRNA and pri-miRNA, differing only in length, and (3) many miRNAs are very closely related in sequences, such as isoforms, differing by only one or a few nucleotides. Therefore, the conventional microarray technologies cannot directly be applied to analyzing these molecules. A number of miRNA microarray products are commercially available, but they are either tedious in requiring pre-isolation of microRNA, lack discriminative power to differentiate isoforms, or are not sensitive enough to monitor low abundant miRNAs.

In our array assay, each miRNA molecule is targeted by two oligos, each that hybridizes a half molecule of the

target miRNA to form a RNA/DNA duplex. When the sequences are perfectly matched, they are aligned with the miRNA and the joint can be ligated by DNA ligase (figure 1). A single nucleotide difference among miRNAs will block either the hybridization or the ligation, so that miRNA isoforms can be differentiated. Due to the small size of miRNA, the hybrid might not be stable; therefore we introduce the stacking sequences. By extending these two oligos along with their complementary oligos the stability is increased. Once the pair of oligos is ligated, the ligated molecules are subjected to linear amplification via T7 transcription into RNA in the presence of biotin-UTP, which are used as probes for array hybridization. To differentiate each isoform, we assigned unique tag sequences to the ligation oligos, so that single nucleotide differences are converted into unique tag sequences. Therefore, each isoform can be easily distinguished by array hybridization.

We offer the miRNA profiling assay kit to profile the expression of the 60 most popular miRNAs and their isoforms. The procedure is simple and straight forward, including three steps: (1) mix the total RNA with provided oligos to form miRNA/oligo hybrids; (2) select the hybrids and remove free oligos, and ligate miRNA-directed pairing of oligos to become a single DNA; and (3) amplify the ligated DNA with T7 transcription.

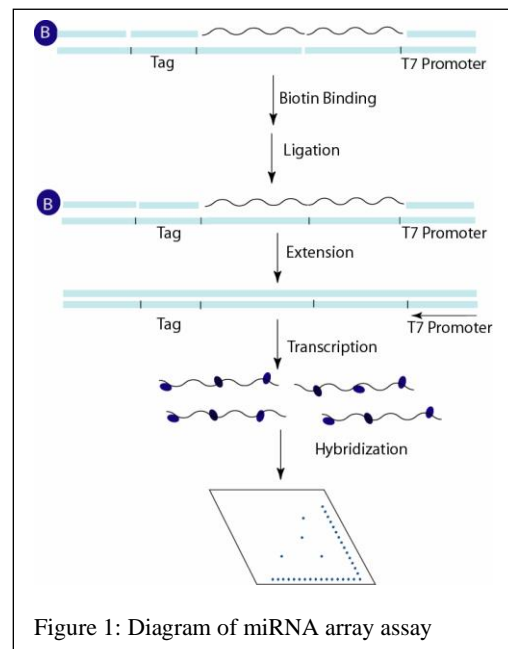


Figure 1: Diagram of miRNA array assay

## Materials provided with the kit

15 µl Cancer oligo mix (-20°C)  
5 µl Array Detection Oligo (-20°C)  
400 µl Annealing buffer (RT)  
15 µl Magnetic streptavidin beads (4 °C)  
120 µl Beads binding buffer (RT)  
1.0 ml Bead wash buffer (RT)  
6 µl Ligase (-20°C)  
250 µl Ligation buffer (-20°C)  
60 µl Extension mix (-20°C)  
60 µl Labeling mix (-20°C)  
6 µl T7 RNA polymerase (-20°C)  
30ml 1x Hybridization buffer (RT)  
30ml 5x Hybridization wash buffer (RT)  
60ml Blocking buffer (4 °C)  
50 µl Streptavidin-HRP conjugate (4 °C)  
40ml 5x Detection wash buffer (RT)  
1.8ml Substrate A (4 °C)  
1.8ml Substrate B (4 °C)  
3 Array membranes (RT)  
3 Detection sheets (RT)

## Material required but not provided

Magnetic stand (96 well plate)  
RNase free water  
PCR machine  
Hybridization oven  
Washing tray  
50ml Centrifuge tubes (Corning tubes are recommended, cat#430290) or hybridization bottles  
0.2ml PCR tube  
Alpha Innotech image or equivalent image system or X-ray film

## Reagent preparation before starting experiment

- Dilute the 5x Hybridization washing buffer and 5x Detection washing buffer to 1x buffer  
1x Hybridization washing buffer:  
30ml 5x Hybridization washing buffer  
120ml ddH<sub>2</sub>O  
1x Detection washing buffer:  
40ml 5x Detection washing buffer  
160 ml ddH<sub>2</sub>O
- Prewarm 1x Hybridization buffer, Blocking Buffer, 1x Hybridization wash buffer at 42°C for 1 hour or until the buffers are clear without visible precipitation before using.
- Pre-hybridization can be done during T7 RNA transcription at the Step 4.

11. Remove the sealing film from the plate.

## 1. Annealing of miRNA with Oligo mix

(1) Sample preparation  
X µl 5ug total RNA or 10ng isolated miRNA  
5 µl Cancer oligo mix  
1 µl Array Detection Oligo  
20 µl Annealing buffer  
X µl ddH<sub>2</sub>O  
-----  
40ul

(2) Incubate on PCR machine at 72°C for 5 minutes and 53°C for 90 minutes.

## 2. Selection of miRNA/oligo hybrids

- (1) Beads washing
- Resuspend the beads by gently tapping the tube to obtain a homogeneous suspension.
  - Transfer 5 µl of the beads to a 0.2ml PCR tube (the size of the tube that should fit into the magnetic stand.
  - Add 100 µl of annealing buffer to the tube and then place onto the magnetic stand for 30 seconds.
  - Aspirate out the liquid.
  - Remove the tube from the magnetic stand.

- (2) Beads selection
- Add 40 µl of the Bead binding buffer to 40 µl annealed miRNA/oligo hybrid mix from Step 1, transfer to the tube containing the washed beads from Step 2(1) and resuspend the beads in the solution.
  - Incubate at 37°C for 30 minutes.
  - Place the bead mixture on the magnetic stand for 30 seconds, and aspirate out the buffer. The beads will remain on the side of the tube.
  - Remove the tube from the magnetic stand and add 100 µl of the Bead wash buffer to resuspend the beads, pipette gently up and down, and place the tube on the magnetic stand for 30 seconds, aspirate the buffer.
  - Repeat the washing step once.

## 3. Ligation of miRNA-directed oligos to form a single molecule

(1) Add 50 µl of the Ligation buffer to resuspend the beads, pipette gently up and down, then place the tube on the magnetic stand for 30 seconds and aspirate the buffer.

(2) Remove the tube from the magnetic stand. Add 20 µl of ligation buffer to resuspend the beads, then add 2µl of Ligase to the resuspended beads and incubate at 37° for 90 minutes.

#### 4. T7 RNA transcription of ligated molecule

- (1) Add 100  $\mu$ l Bead washing buffer directly to 20  $\mu$ l ligat reaction mix from Step 3, place the tube on the magnetic stand 30 seconds, and aspirate the buffer.
- (2) Remove the tube from the magnetic stand and add 20  $\mu$ l of Extension mix to resuspend the beads.
- (3) Incubate the mixture on PCR machine at 94°C for 2 minutes, 54°C for 1 minute, 72°C for 1.5 minute, and 94°C for 30 seconds.
- (4) Place the reaction tube on the magnetic stand for 30 seconds. Immediately transfer the 20  $\mu$ l of the extension mix to a fresh tube (keep the solution and toss the beads).
- (5) Add 20  $\mu$ l of Labeling mix and 1  $\mu$ l of T7 RNA polymerase to the tube.
- (6) Incubate the mixture at 37°C for 1 hour.
- (7) The transcribed RNA is ready for hybridization.

#### 5. Pre-hybridization and hybridization

- (1) Place each array membrane into a 50 ml tube. Wet the membrane by filling the tube with dH<sub>2</sub>O, then carefully decant the water. The side of the membrane with the spotted oligos should face into the middle of the tube.
- (2) Add 4 ml of prewarmed 1x Hybridization buffer to each tube. Incubate the tubes in a hybridization oven at 42°C for at least 30-60 minutes.
- (3) Decant the hybridization buffer and replace with 4ml of prewarmed 1x Hybridization buffer. Add 40  $\mu$ l of transcribed RNA to prehybridized membrane and incubate overnight in a hybridization oven at 42°C.
- (4) Decant the hybridization mixture from each tube and wash each membrane as follows:
  - Rinse the membrane with 20 ml Hybridization washing buffer, and decant liquid.
  - Incubate the membrane with 20 ml Hybridization wash buffer at 42°C for 20 minutes. Decant liquid.

#### 6. Detection

- (1) Using forceps, carefully transfer the membrane from the hybridization tube to a container (an empty 200  $\mu$ l pipette tips box). Each box could have two membranes, one at each side of the box.
- (2) Rinse the membrane with 10 ml of 1X Detection wash buffer.
- (3) Block the membrane with 15 ml of Blocking buffer for 30 minutes at room temperature with moderate shaking.
- (4) Dilute 15  $\mu$ l of Streptavidin-HRP conjugate with 1 ml of the 1X Blocking buffer and transfer to the container. **Do not** add HRP diluted solution directly onto the membrane.
- (5) Continue shaking the membrane for 45 min at room temperature.
- (6) Decant the Blocking buffer and wash three times at room temperature with 15 ml of 1x Detection washing buffer, 10 minutes each wash.

- (7) Mix equal amounts of Substrate A and B. Place the membrane on the bottom side of detection sheet on a flat surface and overlay the membrane with 1 ml of substrate solution. To ensure that the solution remains evenly distributed over the membrane when enveloped by the detection sheet: gently lower the top side of the detection sheet halfway over the membrane then pull back up slightly to allow the solution to flow back over the membrane. Then slowly lay the top sheet down completely without trapping air bubbles. Incubate at room temperature for 5 minutes.
- (8) Remove excess substrate by gently applying pressure over the top sheet using a paper towel. Expose the membranes using either Hyperfilm ECL (2-10 min) or a chemiluminescence imaging system (i.e., FluorChem imager from Alpha Innotech). With either method, experiment with different exposure times.
- (9) Use the schematic diagram of human miRNA array I to identify the spots on the array.

#### Example of miRNA array analysis

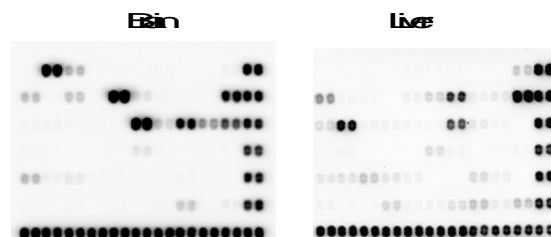
#### Trouble Shooting

##### Signals are too weak

- Total RNA may not contain small RNA
- RNA may be degraded
- If signals of the alignment spots are weak as well, the incubation of the membrane with Streptavidin HRP conjugate may be too short or the exposure time may be too short.

##### Uneven background

- Substrate was not evenly overlaid on the membrane



5 $\mu$ g total RNA was used for miRNA array assay and hybridization was detected with a chemiluminescence imaging system.

## References

1. Ambros, V. (2003), MicroRNA pathways in flies and worms: growth, death, fat, stress, and timing. *Cell*. 113:673–676.
2. Poy, et. al. (2004) A pancreatic islet-specific microRNA regulates insulin secretion. *Nature*. 432:226–230.
3. Berezikov, et. al. (2006) Many novel mammalian microRNA candidates identified by extensive cloning and RAKE analysis. *Genome Res*. 16:1289–1298.
4. Calin, et. al. (2002) Frequent deletions and down-regulation of micro-RNA genes *miR15* and *miR16* at 13q14 in chronic lymphocytic leukemia. *Proc. Natl. Acad. Sci. USA*. 99:15524–15529.
5. Calin, et. al. (2004) MicroRNA profiling reveals distinct signatures in B cell chronic lymphocytic leukemias. *Proc. Natl Acad. Sci. USA*. 101:11755–11760.

Schematic diagram of cancer miRNA array

let7a	let7b	let7c	let7d	let7e	let7f	let7g	let7i	miR1	miR7
miR9	miR9a	miR9a	miR9b	miR9c	miR9d	miR9e	miR9f	miR9g	miR9h
miR10	miR10a	miR10b	miR10c	miR10d	miR10e	miR10f	miR10g	miR10h	miR10i
miR125	miR125a	miR125b	miR125c	miR125d	miR125e	miR125f	miR125g	miR125h	miR125i
miR133a	miR133b	miR133c	miR133d	miR133e	miR133f	miR133g	miR133h	miR133i	miR133j
miR143a	miR143b	miR143c	miR143d	miR143e	miR143f	miR143g	miR143h	miR143i	miR143j
miR145	miR145a	miR145b	miR145c	miR145d	miR145e	miR145f	miR145g	miR145h	miR145i
miR147a	miR147b	miR147c	miR147d	miR147e	miR147f	miR147g	miR147h	miR147i	miR147j
miR150	miR150a	miR150b	miR150c	miR150d	miR150e	miR150f	miR150g	miR150h	miR150i
miR154	miR154a	miR154b	miR154c	miR154d	miR154e	miR154f	miR154g	miR154h	miR154i
miR155	miR155a	miR155b	miR155c	miR155d	miR155e	miR155f	miR155g	miR155h	miR155i
miR157	miR157a	miR157b	miR157c	miR157d	miR157e	miR157f	miR157g	miR157h	miR157i
miR159	miR159a	miR159b	miR159c	miR159d	miR159e	miR159f	miR159g	miR159h	miR159i
miR160	miR160a	miR160b	miR160c	miR160d	miR160e	miR160f	miR160g	miR160h	miR160i
miR161	miR161a	miR161b	miR161c	miR161d	miR161e	miR161f	miR161g	miR161h	miR161i
miR162	miR162a	miR162b	miR162c	miR162d	miR162e	miR162f	miR162g	miR162h	miR162i
miR163	miR163a	miR163b	miR163c	miR163d	miR163e	miR163f	miR163g	miR163h	miR163i
miR164	miR164a	miR164b	miR164c	miR164d	miR164e	miR164f	miR164g	miR164h	miR164i
miR165	miR165a	miR165b	miR165c	miR165d	miR165e	miR165f	miR165g	miR165h	miR165i
miR167	miR167a	miR167b	miR167c	miR167d	miR167e	miR167f	miR167g	miR167h	miR167i
miR168	miR168a	miR168b	miR168c	miR168d	miR168e	miR168f	miR168g	miR168h	miR168i
miR169	miR169a	miR169b	miR169c	miR169d	miR169e	miR169f	miR169g	miR169h	miR169i
miR17	miR17a	miR17b	miR17c	miR17d	miR17e	miR17f	miR17g	miR17h	miR17i