



miProfile™ miRNA PCR Arrays (96-Well)

For high-throughput profiling of miRNA expression

User Manual

GeneCopoeia, Inc.
9620 Medical Center Drive, #101
Rockville, MD 20850
USA

301-762-0888
866-360-9531

support@genecopoeia.com

www.genecopoeia.com

USER MANUAL

miProfile™ miRNA PCR Array

- I. Introduction
- II. Kit Components and Array Format Options
- III. Preparation
- IV. Procedure
- V. Data Analysis
- VI. Appendix I
- VII. Appendix II
- VIII. Limited Use License and Warranty

I. Introduction

MicroRNAs (miRNAs) are small non-coding RNAs that regulate gene expression at the post-transcriptional level. Usually 21-23 nucleotides in length, miRNAs are important modulators in cellular pathways and are highly conserved in eukaryotic organisms. Irregularities in miRNA-regulated gene expression have been found to be associated with cancers, cardiovascular disorders and a variety of other diseases.

The miProfile miRNA PCR Arrays are designed for profiling the expressions of pre-defined or customized sets of miRNAs in various tissues or cells. The resulting differential expressions of profiled miRNAs help researchers to identify those miRNAs that are of biological significance and importance relevant to their research. Each 96-well plate contains up to 84 pairs of PCR primers (forward: miRNA-specific primer; reverse: universal primer), which have been pre-validated and deposited in designated wells. Each plate also has 12 wells that contain different types of controls for monitoring the efficiency of the entire experimental process: from reverse transcription to qPCR reaction.

The All-in-One™ miRNA First-Strand cDNA Synthesis Kits for miRNA qPCR array(QP017, QP018) and qPCR Mix Kits (QP001, QP002, QP004) are the recommended RT-PCR reagents for use with the miProfile miRNA qPCR arrays. These reagents have been optimized to produce high sensitivity, efficiency, and specificity. The All-in-One reverse transcriptase mix contains a novel and optimized blend of polyA polymerase and reverse transcriptase in a buffer that allows high activities and maximal performances of both enzymes. In such reactions, the polyA polymerase adds poly-A tails to mature miRNAs to generate polyA miRNAs. In the same reaction, m-MLV RTase and a unique oligo dT adaptor primer (compatible with the PCR universal reverse primer pre-deposited in the miRNA plates) reverse-transcribe the polyA miRNAs. The All-in-One qPCR Mix containing SYBR[®] Green is used to specifically detect the reverse-transcribed miRNA with the miRNA-specific forward primer and PCR universal reverse primer which are pre-deposited in the miRNA plates. Similar reagents from third-party vendors may be compatible for use. However, their uses are not supported.

Using a universal real-time PCR condition, one can easily profile and analyze the miRNA expression in a high-throughput fashion.

Small RNA is recommended as the input RNA to increase the specificity of detection, although use of total RNA can achieve similar results.

Key advantages

- **Genome-wide coverage, pre-arranged groups, or customized groups**
Largest genome-wide miRNA coverage
Cancer-related groups
Customized miRNA arrays for focused study
- **Robust performance**
Sensitive – Detect miRNAs from as little as 10 pg of input small RNA or 20 pg of total RNA
Specific – Capable of distinguishing miRNAs with single nucleotide mismatches. Each primer set has been experimentally validated for specific amplification
Broad linearity – Allow miRNAs at different expression levels to be detected simultaneously
Reproducible -High reproducibility ($R^2 > 0.99$) for inter-array and intra-array replicates

- Validated miRNA primers
Each miRNA primer is designed using a proprietary algorithm and has been experimentally validated

Protocol overview

A. Prepare cDNA from your RNA Samples



B. Add qPCR Mix and cDNA to the qPCR Array Plate



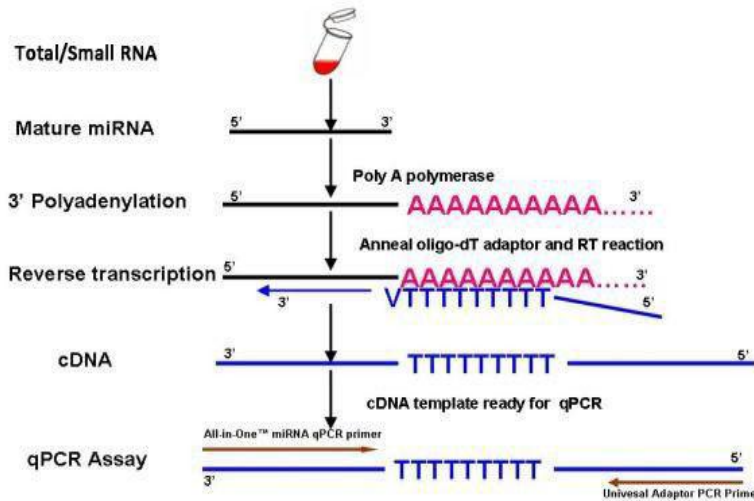
C. Perform real-time PCR



D. Analyze the qPCR Results with GeneCopoeia's Online Data Analysis System

qPCR Array Catalog#	HmiRWG-07	miRNA ID	Homo Brain Assay Data(Sample)			Homo Liver Assay Data(Control)			Fold Change	T Test
			Ct Value-1	Ct Value-2	Ct Value-3	Ct Value-1	Ct Value-2	Ct Value-3		
1	MIMAT0004992	hsa-miR-939	26.27	26.00	26.32	23.79	30.06	30.28	112.69	1.93E-06
2	MIMAT0004951	hsa-miR-807	27.16	27.15	27.30	30.84	30.79	30.95	98.51	3.97E-06
3	MIMAT0004804	hsa-miR-615-5p	32.75	31.99	32.55	33.93	34.02	33.87	22.32	9.52E-05
4	MIMAT0004636	hsa-miR-323-5p	20.32	20.35	20.82	27.69	27.81	27.59	1154.03	6.94E-06
5	MIMAT0003324	hsa-miR-661	27.08	27.25	27.12	31.49	31.89	31.05	156.93	6.38E-04
6	MIMAT0004983	hsa-miR-940	18.95	19.02	18.89	19.12	18.95	19.02	0.27	1.53E-04
7	MIMAT0003306	hsa-miR-636	20.40	20.39	20.45	24.97	24.78	25.11	182.60	1.49E-06
8	MIMAT0005967	hsa-miR-663b	27.41	27.39	27.42	22.80	22.48	22.59	0.28	1.81E-03
9	MIMAT0002178	hsa-miR-487a	23.93	23.94	24.01	31.83	31.56	31.78	1703.48	9.72E-07
10	MIMAT0003180	hsa-miR-487b	21.42	22.01	21.89	28.77	28.89	28.59	987.94	1.18E-05

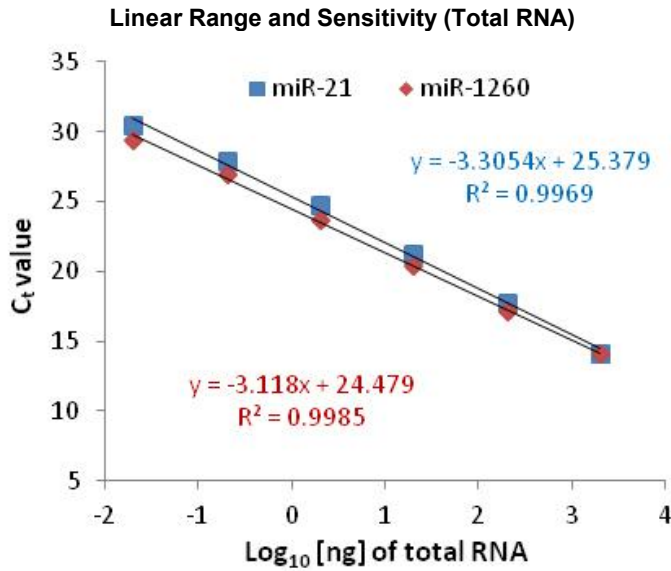
(A)



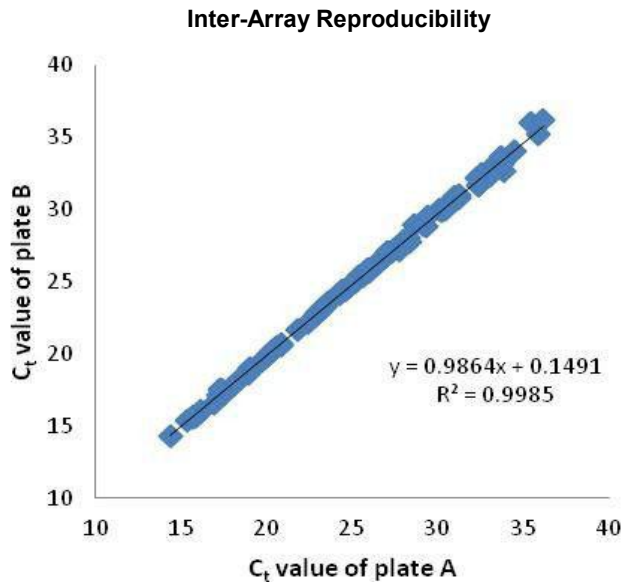
(B)

Figure 1. miRNA PCR array experiment work flow (A) and miRNA RT-PCR mechanism (B)

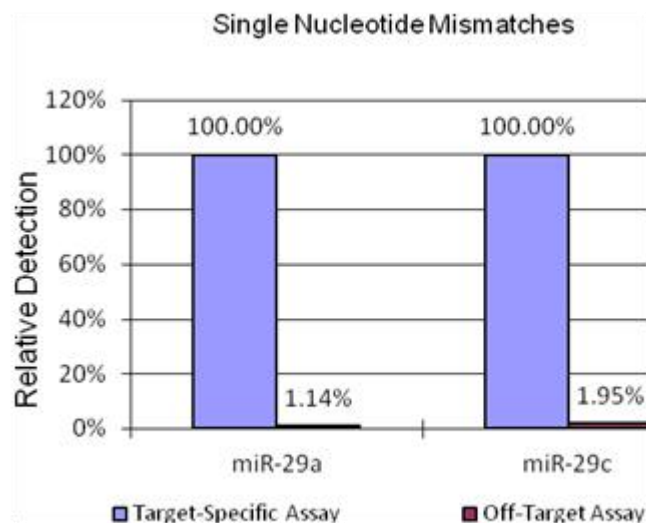
Performance data

**Figure 2. Broad linear range and high sensitivity**

Starting with serially diluted amounts of human colon cancer total RNA, miR-21 and miR-1260 were detected using All-in-One™ miRNA qRT-PCR Detection Kit. The resulting Ct values were plotted against the log₁₀ of the amounts of input total RNA. The data demonstrated a broad linear dynamic range from 20pg to 2 µg of input total RNA as well as high sensitivity. This allows the detection of miRNAs at varying expression levels, including low expressers.

**Figure 3. High inter-array reproducibility**

Two miProfile PCR array replicates (plate A and B) were analyzed using human total RNA (10-tissue mix) on the Bio-Rad iQ5. The Ct values of the replicate plates were plotted against each other. $R^2 > 0.99$ were observed for high inter-array reproducibility. $R^2 > 0.99$ is also observed for intra-array reproducibility (data not shown).



(A)

hsa-miR-29a	UAGCACCAUCUGAAAUCGGUUA
hsa-miR-29c	UAGCACCAU <u>U</u> UGAAAUCGGUUA

(B)

Figure 4. Specificity of miRNA detection

miRNA miR-29a and miR-29c with one single nucleotide mismatch (B) can be distinguished. Relative detection, defined as a percentage of the perfect match ($100\% \times 2^{-\Delta Ct}$), was calculated using the Ct values of on-target and off-target assays, which were performed to detect miRNA plasmid DNA templates using All-in-One miRNA qRT-PCR Detection Kits (A).

II. Product, Array Layout and Array Format Options

Catalog miProfile miRNA qPCR arrays

See the complete list in Appendix II or visit <http://www.genecopoeia.com/product/mirna-solutions/mirna-qpcr-arrays/>.

Array format options

GeneCopoeia provides five qPCR array formats (A, B, C, D, and E) suitable for use with the following real-time cyclers.

Important note: Upon receiving, please check to make sure that the correct array format was ordered to ensure the compatibility with your qPCR instrument.

Plate format	Instrument provider	qPCR instrument model
A (96-well)	Applied Biosystems	5700, 7000, 7300, 7500, 7700, 7900HT (Standard 96-well block), ViiA™7 (Standard 96-well block)
B (96-well)	Applied Biosystems	7500 (Fast block), 7900HT (Fast block), StepOnePlus™, ViiA™7 (Fast block)
C (96-well)	Bio-Rad Laboratories	iCycler iQ®, MyiQ™, iQ™5
D (96-well)	Bio-Rad Laboratories	CFX96™, DNA Engine Opticon™, DNA Engine Opticon 2™, Chromo4™
E (96-well)	Roche Applied Science	LightCycler® 480 (96-well block)

Array layout

	1	2	3	4	5	6	7	8	9	10	11	12
A	1	2	3	4	5	6	7	8	9	10	11	12
B	13	14	15	16	17	18	19	20	21	22	23	24
C	25	26	27	28	29	30	31	32	33	34	35	36
D	37	38	39	40	41	42	43	44	45	46	47	48
E	49	50	51	52	53	54	55	56	57	58	59	60
F	61	62	63	64	65	66	67	68	69	70	71	72
G	73	74	75	76	77	78	79	80	81	82	83	84
H	NC	NC	HK1	HK2	HK3	HK4	HK5	HK6	RT	RT	PCR	PCR

Figure 5. Illustration of miProfile miRNA qPCR array layout (96-well plate)

- **miRNA primer pairs:** Wells 1-84 are designated wells for pre-deposited miRNA primer pairs.
- **NC:** Negative controls, which only have the pre-deposited reverse universal primers
- **HK1-6:** Six pre-deposited housekeeping snRNAs (**HK1-6**) primer pairs, which can be used as endogenous positive controls as well as for array normalization.
- **RT:** Spike-in reverse transcription controls, which can be used to monitor the efficiency of the RT reactions. These pre-deposited primer pairs specifically amplify the cDNA template reversed transcribed from the spike-in exogenous RNA in the sample.
- **PCR:** Positive PCR controls, which are used to verify the PCR efficiency by amplifying the pre-deposited DNA template with its specific pre-deposited primer pairs.

RNA extraction and RT-PCR reagents required (sold separately)

Cat. No.	Products	Quantity/set	Shipping and storage condition
QP020	RNAzol® RT RNA isolation reagent	50 ml	Shipped at room temperature. Stable for at least two years when stored at room temperature.
QP017 QP018	All-in-One miRNA first-strand cDNA synthesis kit For miRNA qPCR array	20 reactions 60 reactions	Shipped with dry ice. Store at –20°C (Stable for at least 12 months). Alternatively, store at –80°C in aliquots. Avoid repeated freezing/ thawing.
QP001 QP002 QP004	All-in-One qPCR mix	200 reactions 1000 reactions 4000 reactions	Shipped with dry ice. Store at –20°C (Stable for at least 12 months). Alternatively, store at –80°C in aliquots. Avoid repeated freezing/ thawing.

Estimates of number of RT-PCR reactions required for EACH SAMPLE

Array format	Numbers of plates per sample	Numbers of RT reactions per sample	Numbers of PCR reactions per sample
96-well plate	1	1	110
	2	1	220
	5	3	550
	10	5	1100
	20	10	2200
	40	20	4400

Other materials required but not provided

Small/total RNA extraction kit (i.e. RNAzol® RT)
 DNase/RNase free tips, PCR reaction tubes, 1.5 ml microcentrifuge tubes
 5 ml and 10 ml graduated pipettes, beakers, flasks, and cylinders

10 µl to 1,000 µl adjustable single channel micropipettes with disposable tips
 5 µl to 20 µl adjustable multichannel micropipette, disposable tips, and reservoir
 qPCR instrument, compatible with miRNA qPCR arrays ordered

III. Preparation

Important notes

1. Before use, remove any condensation that has accumulated on the plate sealing surface and centrifuge plates briefly to collect the contents to the bottom of the plate wells.
2. Strictly follow the standard procedures for PCR to avoid nucleic acid contamination and non-specific amplifications.
3. Read the instructions thoroughly before attempting to perform the procedures.

Estimates of RNA and number of RT-PCR reactions required for EACH SAMPLE

Array format	Number of plates per sample	of Small RNA per recommended per sample	Total RNA recommended per sample	Number of RT reactions per sample	Number of qPCR reactions per sample
96-well plate	2	0.5-1 ug	1-2 ug	1	220
	5	1.5-3 ug	5-10 ug	3	550
	10	2.5-5 ug	5-10 ug	5	1,100
	20	5-10 ug	10-20 ug	10	2,200
	40	10-20 ug	20-40 ug	20	4,400

RNA quantification and quality control

1. Dilute the RNA sample with the RNase-free water and measure the absorbance at 260 nm and 280 nm. A260/280 should be greater than 1.8.
2. Use the formula $A_{260} \times \text{dilution} \times 40 = \mu\text{g RNA/mL}$ to determine the RNA concentration.
3. Check the RNA integrity by agarose electrophoresis.

IV. Procedure

First-strand cDNA synthesis

Note: High-quality cDNA is a prerequisite for accurate detection of miRNA expression. GeneCopoeia's All-in-One miRNA First-Strand cDNA Synthesis Kit is required for small cDNA synthesis.

1. Thaw the reagents in All-in-One miRNA First-Strand cDNA Synthesis Kit, mix by gently flicking the tube, briefly centrifuge to bring the contents to the bottom of the tubes and then place them on ice.
2. Prepare miRNA polyA polymerase (PAP) and reverse transcriptase (RT) reaction mix. Add the following reagents to the ice-chilled RNase-free reaction tubes to a final total volume of 25 µl.

Component	Volume ^a	Quantity
Small RNA		0.5~1.0 µg ^b
2.5U/µl PolyA Polymerase	1 µl	2.5 U
RTase mix	1 µl	
5×PAP/RT buffer	5 µl	1×
Spike-in RT control	1 µl	
ddH2O(RNase/DNase free)	to final 25µl	

- a. cDNA product from a standard miRNA reverse transcription reaction (25 µl) should be enough for 2 plates of 96-Well reactions. Prepare at least 10 standard miRNA reverse transcription reactions for the 19 plates of whole genome miRNA PCR Arrays.
- b. To increase the rate of positive detection, an input of 0.5~1.0µg of small RNA is recommended for

- the standard miRNA reverse transcription reaction (25 µl).
- Perform reverse transcription reaction:
Mix the prepared reaction mix gently by pipetting up and down. Incubate at 37°C for 60 minutes. Terminate the reaction by incubating at 85°C for 5 minutes. After the incubation, dilute the cDNA products 10 times by adding 225µl of sterile water to each RT reaction and use it for the subsequent qPCR reactions. The diluted cDNA can be stored at -20°C for several weeks.

qPCR reaction

Note: Be sure the miProfile miRNA PCR Array plate is compatible with your qPCR instrument before beginning this protocol.

- Thaw the reagents of All-in-One miRNA qPCR Mix Kit. Invert the tubes to mix gently but thoroughly. Briefly centrifuge to bring the contents to the bottom of the tubes and then place them on ice. Remove any condensation that has accumulated on the plate sealing surface and centrifuge briefly to collect the contents to the bottom of the plate wells. Carefully remove sealing film before use 96-Well-qPCR.
- Prepare qPCR solution on ice

Components	1 well	N well ^a
2×All-in-One qPCR Mix	10µl	11µl × N
miRNA cDNA (10 times dilution)	1µl	1.1µl × N
50 X Rox Reference Dye ^b	0.4µl	0.44µl× N
ddH2O		
■ Not using Rox Reference Dye	9 µl	9.9 µl× N
■ Using Rox Reference Dye	8.6 µl	9.5 µl× N
Final Volume	20µl	22µl× N

- miProfile miRNA PCR Array is used to detect multiple miRNAs simultaneously in the same sample. Ensure sufficient mix is available by preparing enough for the number of reactions to be used with a 10% additional volume for pipetting loss.
 - 50×Rox Reference Dye is added only for qPCR instruments that require ROX for calibration.
- Mix the qPCR solution thoroughly and centrifuge briefly. Accurately transfer exactly 20 µl reaction mix to each well. Change tips after each transfer to avoid cross-contamination.
 - Tightly seal the qPCR reaction plate with a new sealing film, Ensure that the film seals smoothly to prevent refraction of light. Centrifuge briefly to remove bubbles.
 - Run qPCR. The following three-step PCR program is recommended for running qPCR.

Cycles	Steps	Temperature	Duration	Detection
1 40	Initial denaturation	95°C	10min ^a	No
	Denaturation	95°C	10sec.	No
	Annealing	60°C ^b	20 sec.	No
	Extension	72°C ^c	10 sec.	Yes

- The DNA polymerase used in the 2X All-in-One qPCR Mix is a special chemically modified hot-start enzyme. The indicated initial denaturation is sufficient to activate the enzyme.
- The annealing temperatures of the cross-linked primers in All-in-One qPCR Primer Array are designed and optimized. For comparing the miRNAs with single nucleotide difference, a higher annealing temperature (65°C) might be necessary.
- The extension time indicated above is suitable for Bio-Rad's iQ5 real-time PCR instrument. Adjust the time duration according to the documentation provided with your instrument.

When using SYBR Green dye to monitor the qPCR reaction, a melting curve analysis should be performed immediately after qPCR cycling.

Temperature range	Heating rate	Constant temperature	Detection
66°C ~ 95°C	0.5°C/unit time	6sec./unit time	Yes

V. Data Analysis

1. Define the baseline

The baseline is the noise level in early cycles. Each real-time PCR instrument has algorithms to perform the baseline-setting. This may be a fixed number of cycles for all samples or adaptive for each sample, depending on the type of instrument that is being used. If the lowest Ct is less than the upper limit of the baseline setting, then the baseline should be manually adjusted. Use the “Linear View” of the amplification plot to determine the earliest visible amplification, and then set the baseline from cycle 2 to two cycles before the earliest visible amplification. Normally it is between 2 to 10 cycles. Do not use cycles greater than 15.

Ensure that baseline settings are the same across all PCR runs in the same analysis to allow comparison of results.

2. Set threshold

Correct placement of the threshold is the next crucial step in data analysis. To adjust the threshold properly, set the threshold value within the exponential phase of all amplification plots when viewed using the logarithmic scale for the y axis. Generally, the expression level of each reference gene should be higher than most other genes.

3. Obtain the Ct or Cp values

The Ct is defined as the cycle when sample fluorescence exceeds a chosen threshold above background fluorescence. This is also known as the Cp or crossing point.

4. Export the data. Most qPCR instruments provide a function for exporting Ct or Cp values to Excel.

5. Analyze the qPCR results using the $\Delta\Delta C_T$ method of relative quantification and interpretation of the control wells.

6. All Ct values reported as greater than 35 or as N/A (not detected) are considered as not detectable.

QC

1. Examined amplification and melting status of each gene using the qPCR instrument software. Each reference gene, RTC and PPC should exhibit only one melting peak per reaction.

2. Examined CT values of the positive PCR control wells (PCR). If the RNA sample is of high quality, the cycling program has been correctly run, and the thresholds have been correctly defined, the value of Ct of PCR should be ~~20-22~~ across all arrays or samples.

3. Examined CT values of the positive RT control wells (RT). If the RNA sample is of high quality, the cycling program has been correctly run, and the thresholds have been correctly defined, the value of Ct of RT should be ~~20-23~~ across all arrays or samples.

Data analysis

Analyze the qPCR result with GeneCopoeia’s online Data Analysis System (free), which is available at <http://www.genecopoeia.com/product/qpcr/analyse/>

This Data Analysis System uses the $\Delta\Delta C_T$ method to perform fold-change analysis or simple statistical analysis of the expression level (C_t or C_p values) for each gene.

1. Download and read the “Primer Array Data Analysis Operation Guide” before performing analysis.

2. Import the C_t or C_p values into the corresponding data analysis template form (*Sample Data.xls* and *Control Data.xls*). Upload the template form and choose the correct reference and analysis factors.

Note: The reference factor chosen for qPCR Primer Array for normalization with the $\Delta\Delta C_T$ method must not be influenced by the experimental design. Therefore use one or more factors that have been previously verified experimentally. A single value or an average of the C_t values for the reference factor can be used for normalization.

3. Perform the specified analysis. When a test is repeated at least three times, statistical results (p value) are provided. The analysis results allow genes of interest to be simply and rapidly selected for further study.

VI. Appendix I

ΔΔCt data analysis method

ΔΔCt data analysis, a relative quantitative analysis technique, is the most simple and direct method for gene expression analyses. The method requires stable expression from a reference gene to normalize the variation introduced by each step, including sample collection, RNA isolation, reverse transcription and amplification. Typically housekeeping genes are used as reference genes.

In qPCR, as in any amplification-based technique, the number of amplification products (N) is calculated as follows:

$$N = N_0 \times (1 + E)^{C_t}$$

N₀: number of template molecules

C_t: threshold cycle

E: amplification efficiency

When the amplification efficiency E is 100%, the number of template molecules in pre-amplification mix is calculated as follows:

$$N_0 = N \times 2^{-C_t}$$

To analyze the change in expression level for the gene of interest in multiple samples using the ΔΔCt method, the amount of the amplification template from different samples is normalized by dividing the expression level of the gene of interest (x) with the reference factor (r) as follows:

$$N_{rel} = N_0x/N_0r = N \times 2^{-C_{tx}} / N \times 2^{-C_{tr}} = 2^{-(C_{tx} - C_{tr})} = 2^{-\Delta C_t}$$

The change in normalized expression levels of the gene of interest (x) between experimental sample (sample 1) and the control sample (sample 2) is as follows:

$$N_{rel1}/N_{rel2} = 2^{-\Delta C_{t1}} / 2^{-\Delta C_{t2}} = 2^{-(\Delta C_{t1} - \Delta C_{t2})} = 2^{-\Delta\Delta C_t}$$

The value of $2^{-\Delta\Delta C_t}$ is the change in expression level of the gene of interest between different samples.

VII. Appendix II

Note: for most updated catalog array list, please visit

<http://www.genecopoeia.com/product/mirna-solutions/mirna-qpcr-arrays/>

Cat. No.	qPCR array products	Quantity/set	Shipping and storage condition
QM001	miProfile™ human miRNome miRNA qPCR arrays	1700 miRNAs 21 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM002	miProfile™ mouse miRNome miRNA qPCR arrays	834 miRNAs 10 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM003	miProfile™ human single-nucleotide mismatch miRNA qPCR arrays	61 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM004	miProfile™ human cancer miRNA qPCR arrays	420 miRNAs 5 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM005	miProfile™ human breast cancer miRNA qPCR arrays	168 miRNAs 2 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM006	miProfile™ human colorectal cancer miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM007	miProfile™ human endometrial cancer miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C

miProfile™ miRNA PCR Array User Manual

QM008	miProfile™ human gastric cancer miRNA qPCR arrays	80 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM009	miProfile™ human hepatocellular carcinoma miRNA qPCR arrays	168 miRNAs 2 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM010	miProfile™ human lung cancer miRNA qPCR arrays	168 miRNAs 2 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM011	miProfile™ human melanoma miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM012	miProfile™ human brain cancer miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM013	miProfile™ human ovarian cancer miRNA qPCR arrays	168 miRNAs 2 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM014	miProfile™ human pancreatic cancer miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM015	miProfile™ human prostate cancer miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM016	miProfile™ human head and neck cancer miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM017	miProfile™ human leukemia miRNA qPCR arrays	168 miRNAs 2 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM018	miProfile™ human bladder cancer miRNA qPCR arrays	79 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM019	miProfile™ human lymphoma miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM020	miProfile™ mouse breast cancer miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM021	miProfile™ mouse colorectal cancer miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM022	miProfile™ mouse hepatocellular carcinoma miRNA qPCR arrays	168 miRNAs 2 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM023	miProfile™ mouse lung cancer miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM024	miProfile™ mouse melanoma miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM025	miProfile™ mouse brain cancer miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C

miProfile™ miRNA PCR Array User Manual

QM026	miProfile™ mouse ovarian cancer miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM027	miProfile™ mouse pancreatic cancer miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM028	miProfile™ mouse prostate cancer miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM029	miProfile™ mouse head and neck cancer miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM030	miProfile™ mouse leukemia miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM031	miProfile™ human inflammatory miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM032	miProfile™ human heart disease miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM033	miProfile™ human immunopathology miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM034	miProfile™ human IPS (Stem cell) miRNA qPCR arrays	168 miRNAs 2 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM035	miProfile™ human muscle disease miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM036	miProfile™ human toxicology related miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM037	miProfile™ human serum and plasma miRNA qPCR arrays	168 miRNAs 2 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM038	miProfile™ mouse inflammatory miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM039	miProfile™ mouse heart disease miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM040	miProfile™ mouse immunopathology miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C
QM041	miProfile™ mouse serum & plasma miRNA qPCR arrays	84 miRNAs 1 x 96-well plate	Shipped at room temperature Stable for at least 6 months when stored at -20°C

VIII. Limited Use License and Warranty

Limited use license

Following terms and conditions apply to use of miProfile miRNA qPCR Arrays (the Products). If the terms and conditions are not acceptable, the Product in its entirety must be returned to GeneCopoeia within 5 calendar days. A limited End-User license is granted to the purchaser of the Product. The Product shall be used by the purchaser for internal research purposes only. The Product is expressly not designed, intended, or warranted for use in humans or for therapeutic or diagnostic use. The Product must not be resold, repackaged or modified for resale, or used to manufacture commercial products or deliver information obtained in service without prior written consent from GeneCopoeia. This Product should be used in accordance with the NIH guidelines developed for recombinant DNA and genetic research. Use of any part of the Product constitutes acceptance of the above terms.

Limited warranty

GeneCopoeia warrants that the Product meets the specifications described in the accompanying Product Datasheet. If it is proven to the satisfaction of GeneCopoeia that the Product fails to meet these specifications, GeneCopoeia will replace the Product. In the event a replacement cannot be provided, GeneCopoeia will provide the purchaser with a refund. This limited warranty shall not extend to anyone other than the original purchaser of the Product. Notice of nonconforming products must be made to GeneCopoeia within 30 days of receipt of the Product. GeneCopoeia's liability is expressly limited to replacement of Product or a refund limited to the actual purchase price. GeneCopoeia's liability does not extend to any damages arising from use or improper use of the Product, or losses associated with the use of additional materials or reagents. This limited warranty is the sole and exclusive warranty. GeneCopoeia does not provide any other warranties of any kind, expressed or implied, including the merchantability or fitness of the Product for a particular purpose.

GeneCopoeia is committed to providing our customers with high-quality products. If you should have any questions or concerns about any GeneCopoeia products, please contact us at 301-762-0888.

© 2018 GeneCopoeia, Inc.

GeneCopoeia, Inc.
9620 Medical Center Drive, Suite 101
Rockville, MD 20850
+1 (301) 762-0888
+1 (866) 360-9531
support@gencopoeia.com

GeneCopoeia Products are for Research Use Only

Copyright © 2018 GeneCopoeia, Inc.

Trademarks: GeneCopoeia™, miProfile™, All-in-One™, miProfile™ (GeneCopoeia Inc.); SYBR® (Molecular Probes); iCycler iQ®, MyiQ™, iQ™5, CFX96™, DNA Engine Opticon™, DNA Engine Opticon 2™, Chromo4™ (Bio-Rad); LightCycler® (Roche); Trizol™, ABi®, ROX®, ViiA™, StepOnePlus™ (Life Technologies); RNAzol® (Molecular Research Center, Inc.); NanoDrop™ (Thermo Scientific).

QM072318