

For Research Use

---

# TAKARA

## PrimerArray™ Series

---

### Product Manual

Product Name	Cat. #
PrimerArray™ Cytokine-cytokine receptor interaction (Human)	PH001
PrimerArray™ Cell cycle (Human)	PH002
PrimerArray™ Cell adhesion molecules (Human)	PH003
PrimerArray™ Jak-STAT signaling pathway (Human)	PH004
PrimerArray™ Natural killer cell mediated cytotoxicity (Human)	PH005
PrimerArray™ Axon guidance (Human)	PH006
PrimerArray™ Focal adhesion (Human)	PH007
PrimerArray™ T cell receptor signaling pathway (Human)	PH008
PrimerArray™ TGF-beta signaling pathway (Human)	PH009
PrimerArray™ Wnt signaling pathway (Human)	PH010
PrimerArray™ Colorectal Cancer & Pancreatic Cancer (Human)	PH011
PrimerArray™ Prostate Cancer & Melanoma (Human)	PH012
PrimerArray™ Small Cell Lung Cancer & Non-small Cell Lung Cancer (Human)	PH013
PrimerArray™ Asthma & Rheumatoid arthritis (Human)	PH014
PrimerArray™ Diabetes mellitus, Typel & Typell (Human)	PH015
PrimerArray™ Cytokine-cytokine receptor interaction (Mouse)	PN001
PrimerArray™ Cell cycle (Mouse)	PN002
PrimerArray™ Cell adhesion molecules (Mouse)	PN003
PrimerArray™ Jak-STAT signaling pathway (Mouse)	PN004
PrimerArray™ Natural killer cell mediated cytotoxicity (Mouse)	PN005
PrimerArray™ Axon guidance (Mouse)	PN006
PrimerArray™ Focal adhesion (Mouse)	PN007
PrimerArray™ T cell receptor signaling pathway (Mouse)	PN008
PrimerArray™ TGF-beta signaling pathway (Mouse)	PN009
PrimerArray™ Wnt signaling pathway (Mouse)	PN010
PrimerArray™ Colorectal Cancer & Pancreatic Cancer (Mouse)	PN011
PrimerArray™ Prostate Cancer & Melanoma (Mouse)	PN012
PrimerArray™ Small Cell Lung Cancer & Non-small Cell Lung Cancer (Mouse)	PN013
PrimerArray™ Asthma & Rheumatoid arthritis (Mouse)	PN014
PrimerArray™ Diabetes mellitus, Typel & Typell (Mouse)	PN015

## Table of Contents

I. Description .....	3
II. Components .....	5
III. Materials Required but not Provided .....	5
IV. Storage .....	5
V. Protocol.....	6
VI. Appendix .....	13
VII. Related Products .....	15

## I. Description

The PrimerArray Series are sets of real-time RT-PCR primers for the analysis of the expression of genes associated with specific biological pathways. Each array contains 96 primer pairs representing 88 genes associated with a biological pathway and 8 housekeeping genes. When comparing an unknown sample to a control sample, gene expression differences can be expressed using the relative quantification method. In addition, expression levels of multiple genes can be screened simultaneously. The PrimerArray Analysis Tool Ver. 2.2\*<sup>1</sup> and/or Multiplate RQ\*<sup>2</sup> are useful for analyzing data obtained using these products.

- \* 1 : Please download the tool from the PrimerArray product page. Using this tool, it is possible to compare a control sample and one unknown sample.
- \* 2 : The Multiplate RQ software is used for relative quantification of gene expression data obtained with Thermal Cycler Dice Real Time Systems including the Thermal Cycler Dice Real Time System // MRQ (Cat. # TP960)\* and the Thermal Cycler Dice Real Time System Single MRQ (Cat. # TP870)\*. With Multiplate RQ, comparative analysis can be performed between numerous unknown samples and a control sample.

\*Not available in all geographic locations. Check for availability in your region.

### Procedural Overview :

Below is a flowchart, starting with RNA extraction and ending with data analysis. In parentheses is an estimate of time required for one real-time PCR experiment, and the entire process can be finished in approximately 2.5 to 3.5 hours. Multiple samples can be processed together through the reverse transcription step, and the synthesized cDNA should be stored at -20°C for subsequent experiments.

#### RNA extraction (approx. 30 min) :

Extract RNA from experimental materials (control sample and test sample using Nucleospin® RNA II (Cat. #740955.50) or RNAiso Plus (Cat. #9108)) and treat with DNase I. Use approximately 2.5 µg of total RNA for one experiment.



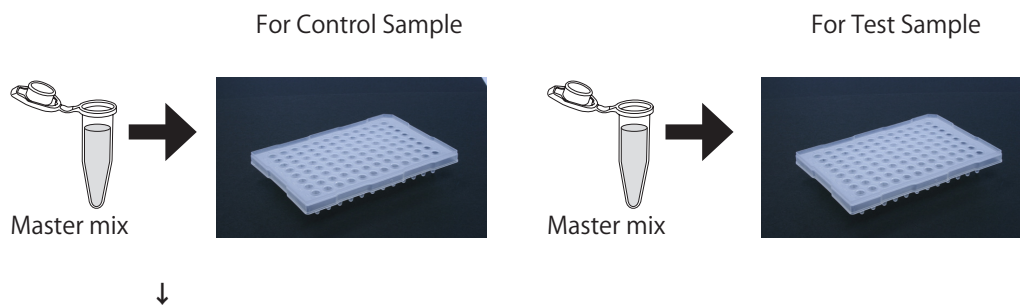
#### Reverse transcription (approx. 20 min) :

Synthesize cDNA from each of the total RNA samples. It is recommended that PrimeScript™ RT Master Mix (Perfect Real Time) (Cat. # RR036A) be used.



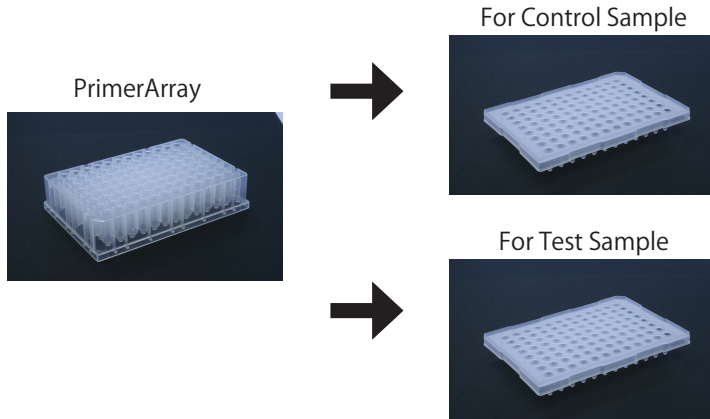
#### Dispense real-time PCR reaction solution (approx. 10 min) :

Combine the synthesized cDNA and SYBR® *Premix Ex Taq* II (Tli RNase Plus) (Cat. # RR820A) to prepare a master mix solution for control and test samples, then dispense into wells of a 96-well real-time PCR plate (see illustration below).



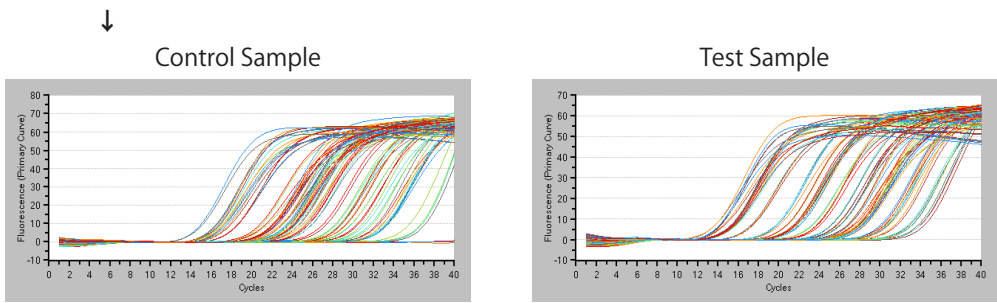
**Add primers (approx. 5 min) :**

Add PrimerArray primers to the real-time PCR plate using 8-multichannel pipette, etc. (see illustration below).



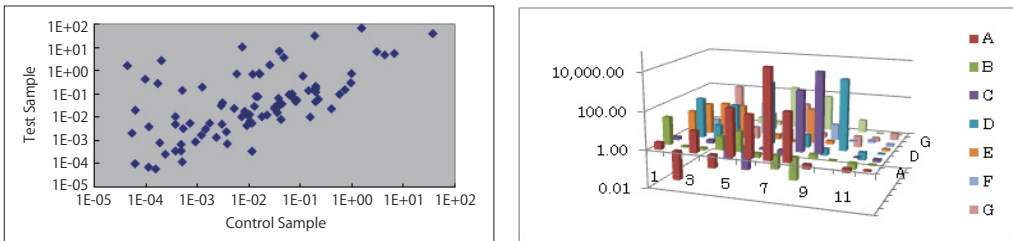
**Real-time PCR (1 - 2 hours) :**

The reaction is carried out with real-time PCR instrument.



**Data analysis (about 30 minutes) :**

Relative quantification analysis is achieved by the  $\Delta\Delta C_t$  method.



## II. Components

### PrimerArray

Primers for pathway-related genes 50  $\mu$ l x 88 wells

Primers for housekeeping genes 50  $\mu$ l x 8 wells

\* Forward & Reverse primer mix (2.5  $\mu$ M each) per well.



### NOTES:

The rubber mat (lid) is re-used repeatedly; therefore, do not discard.

Primer information can be downloaded in excel spreadsheet format from the PrimerArray product page.

### Primer 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	85	86	87	88	89	90	91	92	93	94	95	96

 Primers for pathway-related gene  
 Primers for housekeeping gene

## III. Materials Required but not Provided

Thermal cycler for real-time PCR

Thermal Cycler Dice Real Time System II (Cat. #TP900/TP960)\*

Reaction plate and seal for real-time PCR

8-multi-channel pipette, micropipette, and tips

Centrifuge for 96 well plate

RNA extraction reagent

NucleoSpin® RNA II (Clontech, Cat. # 740955.10); RNAiso Plus (Cat. #9108/9109)\*

Recombinant DNase I (RNase-free) (Cat. #2270A)

Reverse transcription reagents

PrimeScript RT Master Mix (Perfect Real Time) (Cat. #RR036A) is recommended

Real time PCR reagent

SYBR® Premix Ex Taq II (Tli RNaseH Plus) (Cat. #RR820A) is recommended

PrimerArray Analysis Tool Ver. 2.2 or Multiplate RQ

\*: Not available in all geographic locations. Check for availability in your region.

## IV. Storage -20°C

\* Expires 1 year from date of receipt when kept under proper storage conditions.

\* If this product is used over a short time period (~1 month), it should be stored at 4°C. This product does not contain a preservative, therefore caution should be used when handling the product to prevent contamination.

## V. Protocol

Please refer to the operation manual for instruments and instructions for specific reagents.

### 1. Preparation :

When the PrimerArray plate is stored at  $-20^{\circ}\text{C}$ , remove from the freezer and thaw at room temperature prior to use. After checking that the rubber mat (lid) is secure, gently shake the PrimerArray plate to uniformly mix the solution. Then briefly centrifuge the plate to collect the solution at the bottom of the well. When the PrimerArray plate is stored at  $4^{\circ}\text{C}$ , mix the plate gently and spin down briefly.

### 2. RNA extraction :

Please refer to "1. Preparation of RNA sample" in VI. Appendix for general considerations for RNA preparation.

#### 2-1. RNA Isolation

It is recommended that RNA isolation kits such as NucleoSpin® RNA II (Clontech, Cat. # 740955.10) or RNAiso Plus (Cat. #9108/9109)\* be used for isolating high-purity total RNA. Please refer to each instruction manual for detailed protocols. After removing the genomic DNA (see step 2-2), dissolve the RNA sample in sterile distilled water or TE buffer, and adjust the concentration to  $250\text{ ng}/\mu\text{l}$  in preparation for reverse transcription.

\*: Not available in all geographic locations. Check for availability in your region.

#### 2-2. Removal of genomic DNA

In some cases, total RNA samples may contain a small amount of genomic DNA, which could potentially be amplified by PCR and lead to inaccurate results. To avoid this situation, we recommend removing genomic DNA by DNase I treatment.

##### Removal of genomic DNA by DNase I treatment

After extracting total RNA, the RNA sample is treated with Recombinant DNase I (RNase-free) (Cat. #2270A). After the treatment, DNase I should be inactivated by either heat treatment or phenol/chloroform extraction.

(1) Prepare the following reaction mixture:

Total RNA	$x\ \mu\text{l}$	(Final) 20 - 50 $\mu\text{g}$
10X DNase I Buffer	5 $\mu\text{l}$	1X
RNase Inhibitor (40 U/ $\mu\text{l}$ )	0.5 $\mu\text{l}$	20 U
DNase I (RNase-free)	2 $\mu\text{l}$	10 U
DEPC-treated water	$(42.5-x)\ \mu\text{l}$	to 50 $\mu\text{l}$

(2) Incubate at  $37^{\circ}\text{C}$  for 20 min.

(3) Perform one of the following procedures to inactivate DNase I:

##### A. Heat treatment :

- i) Add 2.5  $\mu\text{l}$  0.5 M EDTA, and incubate at  $80^{\circ}\text{C}$  for 2 min.
- ii) Bring the reaction volume to 100  $\mu\text{l}$  with DEPC treated water.

##### B. Phenol/Chloroform extraction :

- i) Add 50  $\mu\text{l}$  of DEPC treated water and 100  $\mu\text{l}$  of phenol/chloroform/isoamyl alcohol (25 : 24 : 1), and mix well.
- ii) Centrifuge at 15,000 rpm for 5 min. at room temperature, and transfer the upper layer to a new tube.
- iii) Add equal amount of chloroform/isoamyl alcohol (24 : 1), and mix well.
- iv) Centrifuge at 15,000 rpm for 5 min. at room temperature, and transfer the upper layer to a new tube.

- (4) Add 10  $\mu$ l of 3 M sodium acetate, 250  $\mu$ l of cold ethanol, and mix well. Then incubate on ice for 10 min.
- (5) Centrifuge at 15,000 rpm for 15 min. at 4°C, and remove the supernatant.
- (6) Wash the pellet with 70% ethanol, centrifuge at 15,000 rpm for 5 min. at 4°C, and remove the supernatant.
- (7) Dry the pellet.
- (8) Dissolve the pellet in an appropriate amount of DEPC treated water.

### 3. Reverse Transcription :

Perform reverse transcription using the total RNA prepared above as a template. It is recommended that PrimeScript RT Master Mix (Perfect Real Time) (Cat. #RR036A) be used for reverse transcription.

- (1) Prepare the reverse transcription mixture on ice.

[For 1 reaction]

	Volume	Final.
5X PrimeScript RT Master Mix	10 $\mu$ l	1X
total RNA (250 ng/ $\mu$ l)	10 $\mu$ l	2.5 $\mu$ g
RNase Free dH <sub>2</sub> O	30 $\mu$ l	
Total	50 $\mu$ l	

\* It is possible to scale up the RT reaction as needed.

- (2) Incubate the reaction mixture under the following conditions
  - 37°C 15 min. (Reverse transcription reaction)
  - 85°C 5 sec. (Heat inactivation of reverse transcriptase)
  - 4°C

**NOTE :** The subsequent protocol depends on the specific real-time PCR instrument used. In this section, the protocol for the Thermal Cycler Dice Real Time System is described. For other instruments, please refer to the section, "2. How to use with various real-time PCR instruments" in VI. Appendix.

### 4. Preparation of Real-time PCR Reaction plates.

The preparation protocol shown in this section is for two real-time PCR reaction plates, one control sample and one unknown sample. Please scale up the number of reaction plates according to the number of experimental samples.

#### 4-1. Preparation and dispensing of Master Mix

The cDNA prepared in "3. Reverse Transcription" and SYBR® *Premix Ex Taq* II (Tli RNaseH Plus) (Cat. #RR820A) are combined to prepare a master mix, then the mixture is dispensed to real-time PCR plates.

- (1) Using cDNA of the control sample or of the unknown sample, prepare each master mix (PCR reaction) as outlined below.

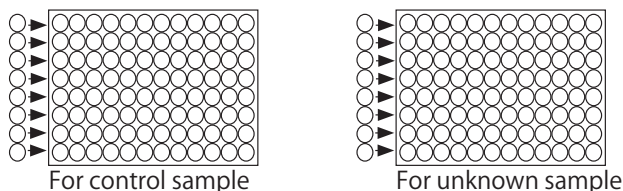
	For 1 well	For 110 wells
SYBR® Premix Ex Taq II (Tli RNaseH Plus) (2X)	12.5 $\mu$ l	1375 $\mu$ l
cDNA (50 ng/ $\mu$ l)*	0.4 $\mu$ l	44 $\mu$ l
sterile distilled dH <sub>2</sub> O	8.1 $\mu$ l	891 $\mu$ l
<b>Total</b>	<b>21 <math>\mu</math>l</b>	<b>2310 <math>\mu</math>l</b>

\* : total RNA equivalent

- (2) Dispense 21  $\mu$ l of the master mix into each well of the plate that will be used for the real-time reaction.

Example using an 8-multichannel pipette :

- 1) Dispense 273  $\mu$ l (sufficient for 13 wells) of the master mix into an appropriate 8-multichannel container.
- 2) Dispense 21  $\mu$ l of the mix to each well of the plate for real-time PCR with an 8-multichannel pipette.



#### 4-2. Addition of Primers:

The PrimerArray primers are added into the plate for real-time PCR that contains the master mix.

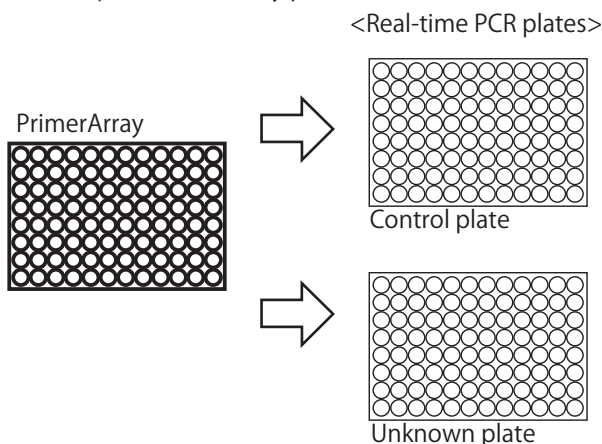
- (1) Remove the rubber mat (lid) of PrimerArray.

The rubber mat should be removed carefully, ensuring that primer solutions do not leak into adjacent wells. Please pay close attention to this as contamination will occur if primer solutions mix. (Do not throw away the rubber mat.)

- \* The primer solution should be mixed well, and spun down briefly prior to use.
- \* After using the PrimerArray, it should be covered with the rubber mat again and stored appropriately.

- (2) Dispense primers.

Add 4  $\mu$ l of PrimerArray primers to each well of the real-time PCR plates.





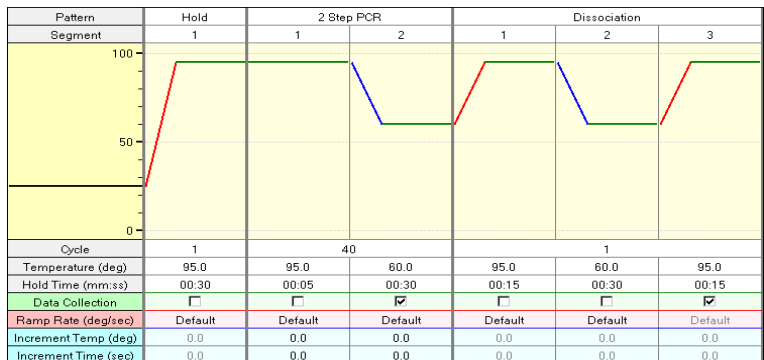
- (3) Cover the plate with Sealing Film for real-time PCR.  
 Firmly attach the Sealing Film to the plate using a special spatula (Plate Sealing Pad). If it is not attached firmly, the reaction mixtures will evaporate in the unsealed areas of the plate during real-time PCR reaction, making it possible to get inaccurate results .  
 \* Preparation of real-time PCR reaction plate is completed. If real-time PCR will not be started immediately, store the plate at 4°C protected from light. We highly recommend starting the reaction within 24 hours after plate preparation.
- (4) Cover the PrimerArray with the rubber mat.  
 The rubber mat should be firmly reattached to the plate. If it is not attached firmly, evaporation or contamination of the primer solution may occur during storage.

**5. Real-time PCR reaction :**

- (1) Centrifuge the real-time PCR plate briefly to collect the reaction mixture at the bottom of well.
- (2) Set the plate in the real-time PCR instrument, and start the reaction.
- (3) Perform PCR using the conditions below.

Initial denaturation  
 95°C 30 sec.  
 2 Step PCR  
 95°C 5 sec.  
 60°C 30 sec.  
 (40 cycles)

Melting curve analysis



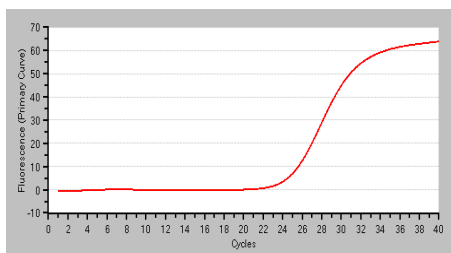
**6. Data Analysis :**

Ct values are calculated by the real-time PCR instrument's analysis software after the reaction. Relative quantification analysis is performed using PrimerArray Analysis Tool Ver. 2.2 and/or Multiplate RQ.

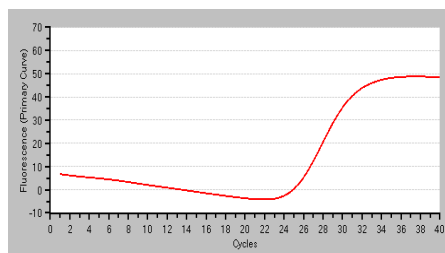
- (1) Setting analysis parameters : For most real-time PCR software, the analysis parameters are set automatically. Check that the setting parameters are correct, and reset them manually if necessary.

- Baseline region : Set the flat region of the curve on the left side of the graph as a baseline region. If the baseline is set incorrectly, the amplification curve will not be flat, but will tilt.

**Note :** The baseline correction will not work well if the region is narrow, and if the region is too wide, the curve will progressively slope right-down. (see below)

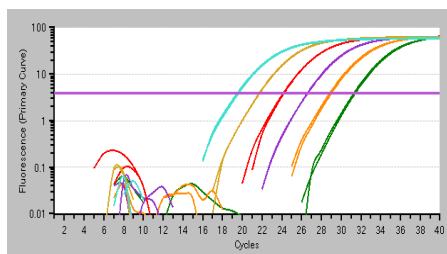


Baseline is set correctly



Baseline region is too wide

- Threshold : Set the threshold in the exponential region of the primary curve. The exponentially amplified region is the range where the primary curve becomes a straight line when the fluorescence intensity (vertical axis) is plotted is on a log scale.



Correctly set threshold

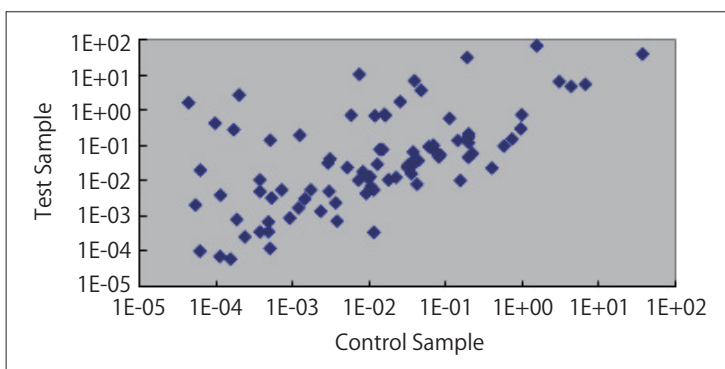
- (2) Calculation of Ct value and Tm value : Ct and Tm value are automatically calculated by the real-time PCR software.
  - (3) Output data : Export Ct values in either the Excel format or CSV format to perform relative quantification analysis using the PrimerArray Analysis Tool Ver. 2.2\*. When using the Thermal Cycler Dice Real Time System, you do not need to obtain Ct values for performing analysis with Multiplate RQ. For data obtained with other instruments, Multiplate RQ can not be used for analysis.
- \* : For some types of real-time PCR software, data generated from wells without sample information are omitted from analysis and may not be exported. This can lead to mistakes when the data is imported into PrimerArray Analysis Tool Ver. 2.2. Therefore, the data should be reviewed carefully to make sure that all of the well information is correct when imported into the Tool.

(4) Relative quantification analysis : Expression of each gene in the unknown sample is relative to the expression in control sample. Please refer to the Multiplate RQ manual and/or PrimerArray Analysis Tool Ver. 2.2 manual for details of the analysis procedure.

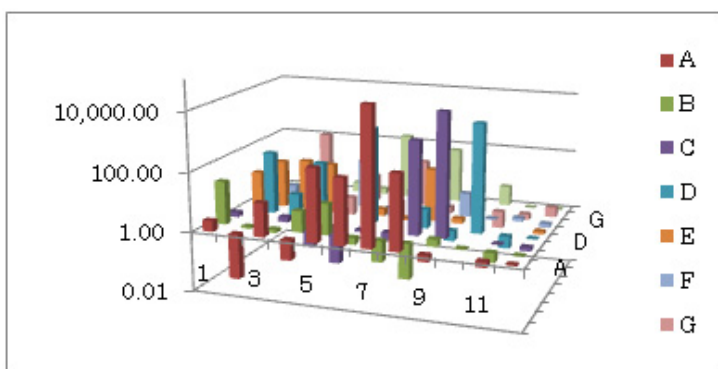
Analysis using PrimerArray Analysis Tool Ver. 2.2

The results comparing the control sample with one unknown sample are shown graphically as a Scatter Plot and a 3D Profile of Fold Difference.

- ScatterPlot : The expression level of the control sample is plotted on the horizontal axis, and the expression level of the unknown sample is plotted on the vertical axis.
- 3D Profile (Fold Difference) : The expression level (Fold Difference) of the unknown sample is normalized to the expression level of the control sample (i.e., control expression set to 1.0).



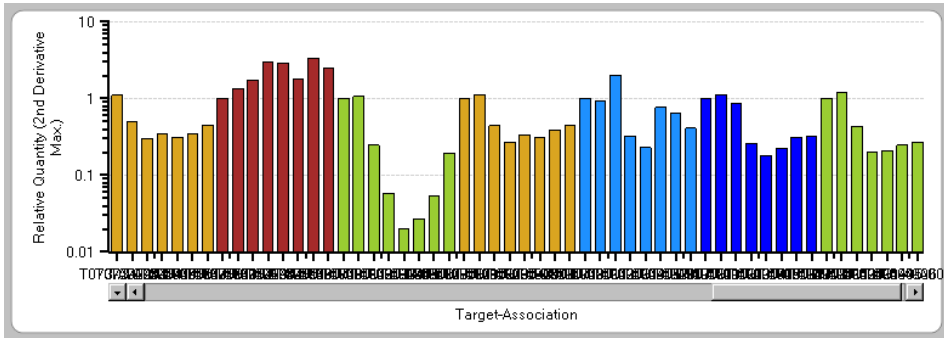
Scatter Plot



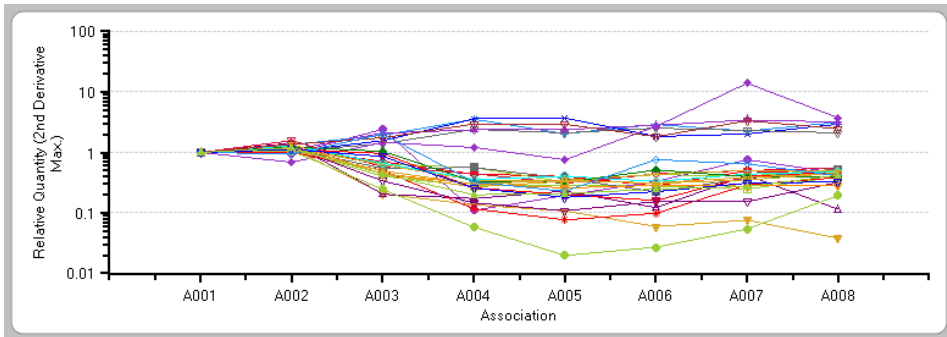
3D Profile

Analysis using Multiplate RQ

The results comparing numerous unknown samples and a control sample (relative quantification) are shown graphically as a bar chart (Relative Quantity Chart) or a line graph (Gene Profile).



Relative Quantity Chart



Gene Profile

## VI. Appendix

### 1. Preparation of RNA sample

These kits are designed to perform the reverse transcription of RNA to cDNA and subsequent amplification. It is important to use an RNA sample with high purity for better cDNA yields. In addition, it is essential to inhibit cellular RNase activity and to prevent contamination with RNase from the equipment and solutions used. Extra precautions should be taken during the sample preparation to prevent the RNase contamination from operators' sweat or saliva, including using disposable gloves and a workspace dedicated exclusively for RNA preparation, and avoiding unnecessary speaking during sample preparation.

#### Equipment to be used

- (1) Dry-heat sterilization (180°C, 60 min).
- (2) Treatment with 0.1% Diethylpyrocarbonate (DEPC) at 37°C for 1 hour, followed by autoclaving at 120°C for 30 min to remove DEPC. It is recommended that all the equipment be used as the exclusive use for RNA preparation.

Disposable plastic equipment should be used. Glass tools should be treated with either of the above protocols prior to use.

#### Reagents

All reagents to be used in this experiment must be prepared using tools that were treated as described in previous section (Dry-heat sterilization or DEPC treatment), and distilled water must be treated with 0.1% DEPC and autoclaved. All reagents and distilled water should be used exclusively for RNA experiments.

### 2. How to use with various real-time PCR instruments

Please operate according to the instruction manual of each instrument. Please refer to "V. Protocol" for details and precautions for the experimental procedure.

#### **[Applied Biosystems 7300/7500 Real Time PCR System (Life Technologies)]**

- (1) Prepare a master mix (PCR reaction) using each cDNA as shown below.

	For 1 well	For 110 wells
SYBR® <i>Premix Ex Taq</i> II (Tli RNaseH Plus) (2X)	25 $\mu$ l	2750 $\mu$ l
ROX Reference Dye or Dye II (50X)* <sup>1</sup>	1 $\mu$ l	110 $\mu$ l
cDNA (50 ng/ $\mu$ l)* <sup>2</sup>	0.8 $\mu$ l	88 $\mu$ l
dH <sub>2</sub> O (sterile distilled water)	15.2 $\mu$ l	1672 $\mu$ l
<b>Total</b>	<b>42 <math>\mu</math>l</b>	<b>4,620 <math>\mu</math>l</b>

\* 1 : ROX Reference Dye II is used for Applied Biosystems 7500 Real Time PCR System. ROX Reference Dye is used for 7300 Real Time PCR System.

\* 2 : total RNA equivalent

- (2) Dispense 42  $\mu$ l of the mixture into each well of a real-time PCR plate.

- (3) Add 8  $\mu$ l of each PrimerArray primer to wells of the real-time PCR plate.  
 (4) Perform real-time PCR reaction using the following conditions.

Initial denaturation

95°C                      30 sec.

2 Step PCR

95°C                      5 sec.  
 60°C                      31 or 34 sec.    \*3    ] 40 cycles

Melting curve analysis

\* 3 : For the Applied Biosystems 7300, set to 31 sec; for the Applied Biosystems 7500, set to 34 sec.

**[Applied Biosystems 7500 Fast Real Time PCR System (Life Technologies)]**

- (1) Prepare a master mix (PCR reaction) using each cDNA as shown below.

	For 1 well	For 110 wells
SYBR® <i>Premix Ex Taq</i> II (Tli RNaseH Plus) (2X)	10 $\mu$ l	1100 $\mu$ l
ROX Reference Dye II (50X)	0.4 $\mu$ l	44 $\mu$ l
cDNA (50 ng/ $\mu$ l) *	0.4 $\mu$ l	44 $\mu$ l
dH <sub>2</sub> O(sterile distilled water)	6 $\mu$ l	660 $\mu$ l
<b>Total</b>	<b>16.8 <math>\mu</math>l</b>	<b>1848 <math>\mu</math>l</b>

\* : total RNA equivalent

- (2) Dispense 16.8  $\mu$ l of the mixture into each well of a real-time PCR plate.  
 (3) Add 3.2  $\mu$ l of each PrimerArray primer to wells of the real-time PCR plate.  
 (4) Perform real-time PCR reaction using the following conditions.

Initial denaturation

95°C                      30 sec.

2 Step PCR

95°C                      3 sec.  
 60°C                      30 sec.    ] 40 cycles

Melting curve analysis

**VII. Related Products**

- PrimerArray™ Embryonic Stem Cells (Cat. #PH016, PN016)
  - PrimeScript™ RT Master Mix (Perfect Real Time) (Cat. #RR036A)
  - SYBR® *Premix Ex Taq*™ II (Tli RNaseH Plus) (Cat. #RR820A/B)
  - Real-time PCR systems:
    - Thermal Cycler Dice™ Real Time System // (Cat. #TP900)\*
    - Thermal Cycler Dice™ Real Time System // MRQ (Cat. #TP960)\*
    - Thermal Cycler Dice™ Real Time System Single (Cat. #TP850)\*
    - Thermal Cycler Dice™ Real Time System Single MRQ (Cat. #TP870)\*
  - 96well Hi-Plate for Real Time (Cat. #NJ400)\*
  - Sealing Film for Real Time (Cat. #NJ500)\*
  - NucleoSpin® RNA II (Clontech, Cat. #740955.10/.50/.250)
  - RNAiso Plus (Cat. #9108/9109)\*
  - Recombinant DNase I (RNase-free) (Cat. #2270A/B)
- \* Not available in all geographic locations. Check for availability in your region.

**NOTE:** This product is for research use only. It is not intended for use in therapeutic or diagnostic procedures for humans or animals. Also, do not use this product as food, cosmetic, or household item, etc.

Takara products may not be resold or transferred, modified for resale or transfer, or used to manufacture commercial products without written approval from TAKARA BIO INC.

If you require licenses for other use, please contact us by phone at +81 77 543 7247 or from our website at [www.takara-bio.com](http://www.takara-bio.com).

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product web page. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

All trademarks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions.