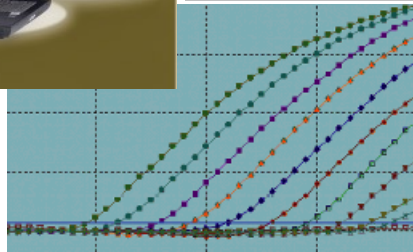
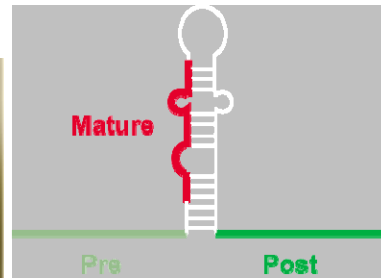


Microarray and QPCR applications for miRNAs



Cathy Cutler
Field Application Scientist
Stratagene Products Division

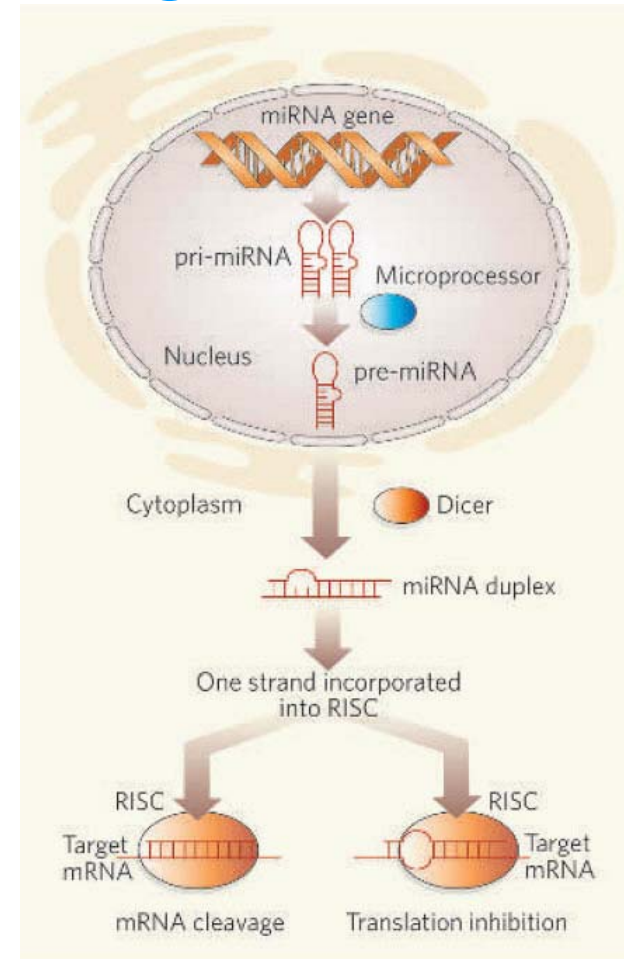


Introduction to miRNA

What are microRNAs

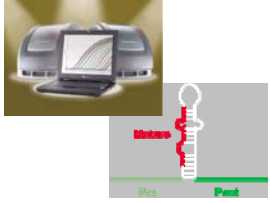
- **Non-coding small RNAs (19-30 nt)**
- **3 different forms:**
 - Pri- and pre-miRNA: hairpin shaped**
 - mature miRNA: bound to RISC complex**
- **Act in repression of translation or direct mRNA degradation**
- **Implicated heavily in disease (eg. cancer), development, apoptosis, proliferation and differentiation**
- **Involved in viral infections**
- **More than 700 human miRNAs identified (Sanger miRBase 10.0)**
- **Implicated in regulation of up to 30% of all human genes**
- **Projected \$100M to be awarded by the NIH for miRNA-related research in 2008***

*Source: NIH CRISP database at <http://crisp.cit.nih.gov/>



Cancer Genomics: Small RNAs with big impacts
from Nature 435: 745-746 (9 June 2005)



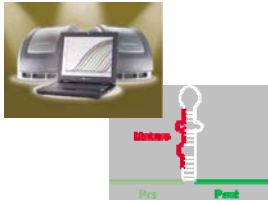


Introduction to miRNA

Why Are miRNAs Important?

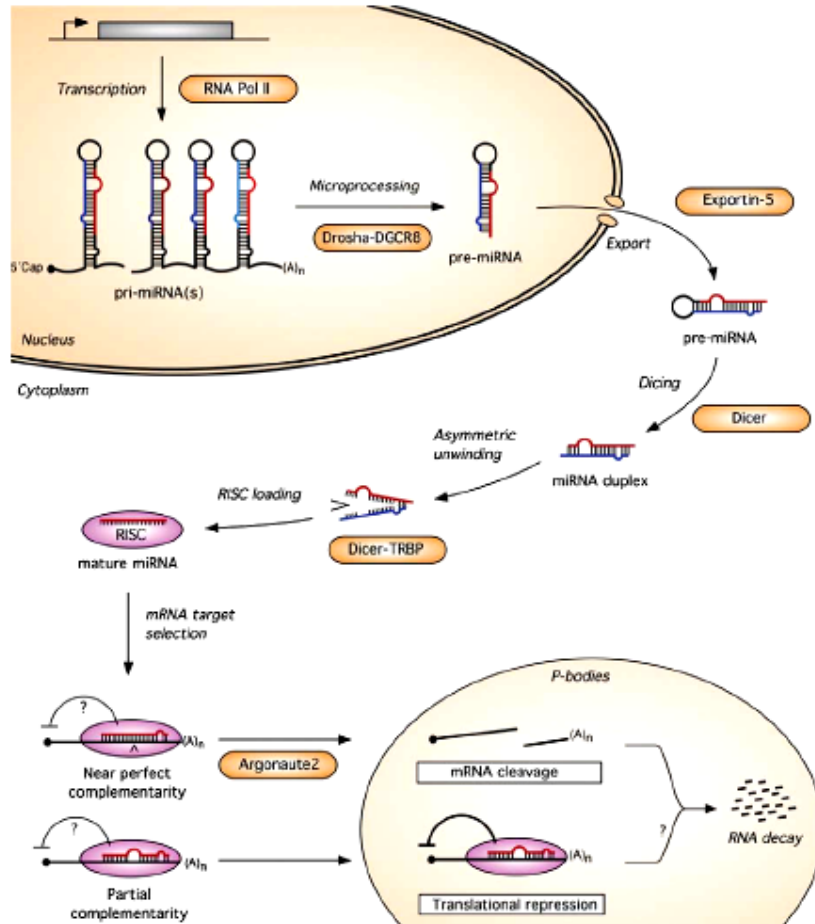
- Predicted to regulate expression of 30% of human genes
- Act as key regulators of many cellular processes, such as:
 - **Early development**
 - **Cell proliferation and apoptosis**
 - **Fat metabolism**
 - **Cell differentiation**
- Involved in viral infection processes
- Differentially expressed miRNAs are reported in many human cancers
- > 50% of miRNA genes are located within regions associated with amplification, deletion, and translocation in cancer (Calin, G.A. et al PNAS 2004 101:2999-3004)





Introduction to miRNA Biogenesis in mammals

E. Wienholds, R.H.A. Plasterk / FEBS Letters 579 (2005) 5911-5922



- Expression may be regulated by transcription factors
- monocistronic and polycistronic (40%) miRNA genes expressed by pol II promoter
- Intronic miRNA genes (~50%) regulated with mRNA that flanks miRNA
- miRNA transcripts can contain more than one miRNA



Introduction to miRNA

Mode of Action – Repression of Translation



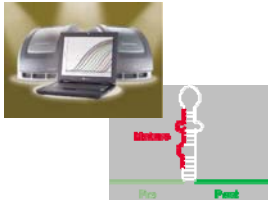
miRNA

miRNA bound by RISC complex

RISC bound miRNA binds target mRNA

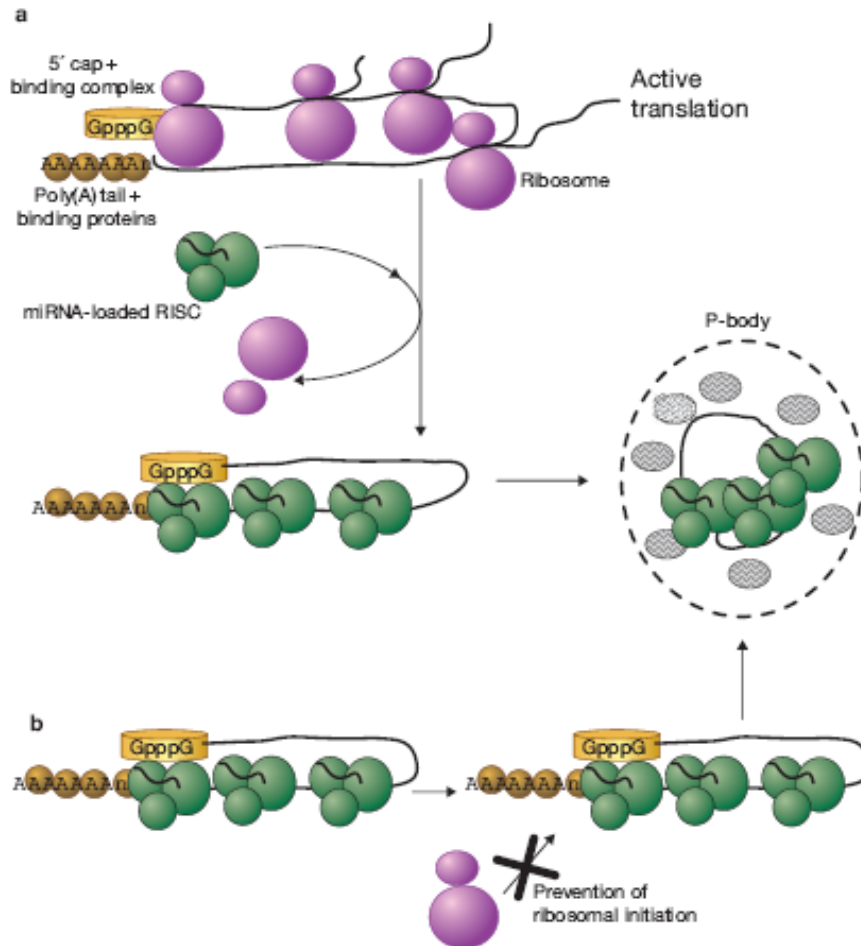
Bound miRNA-RISC blocks translation of mRNA by ribosome





Introduction to miRNA

Mode of Action – RNA Degradation



P-body
contains RNA that can no longer be translated and enzymes that are required for RNA decay and degradation

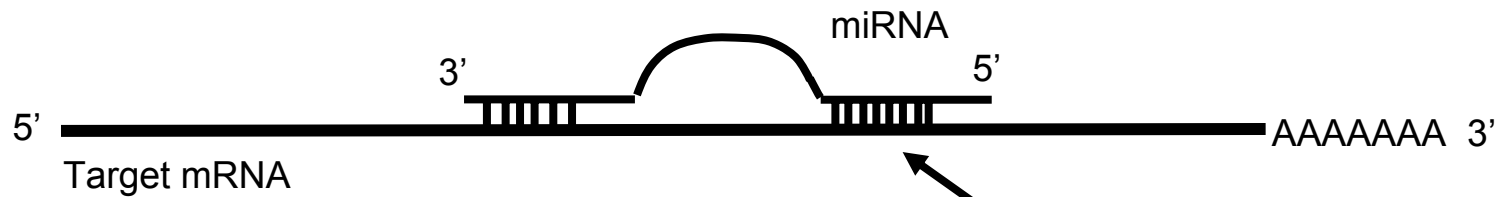
mRNA degradation allows for detecting effects of miRNA using microarrays

Rossi, J J 2005 Nature Cell Biol 7(7):643-644



Interaction of miRNA with mRNA

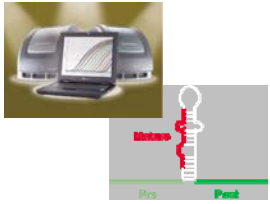
Family Grouping by Seed Region



Perfect homology, 5'
6-8nt of miRNA (seed)

- one miRNA may regulate hundreds of genes
- a gene may be regulated by more than one miRNA

miRNA	nucleotide sequence seed (5'-3')
hsa-let-7a	UGAGGUAGUAGGUUGUAUAGUU
hsa-let-7b	UGAGGUAGUAGGUUGUGUGGUU
hsa-let-7c	UGAGGUAGUAGGUUGUAUGGUU
hsa-let-7d	AGAGGUAGUAGGUUGCAUAGU
hsa-let-7e	UGAGGUAGGAGGUUGUAUAGU
hsa-let-7f	UGAGGUAGUAGAUUGUAUAGUU
hsa-miR-98	UGAGGUAGUAAGUUGUAUUGUU
hsa-let-7g	UGAGGUAGUAGUUUGUACAGU
hsa-let-7i	UGAGGUAGUAGUUUGUGCUU



Introduction to miRNA

miRNA Registry

- <http://www.sanger.ac.uk>
- Searchable database of all published miRNA sequences and annotation
- Mouse and human are highly conserved
- Human is not conserved with plants
- Data can be downloaded from ftp site:
 - <ftp://ftpsanger.ac.uk/pub/mirbase>

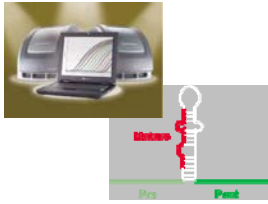




miRNA qRT-PCR Methods and Microarray Methods

- QRT-PCR miRNA detection methods are more specific than microarray methods.
- Microarray methods tend to show higher cross-hybridization which can lead to false positives. Data should be confirmed by QRT-PCR
- Microarray methods are ideal for profiling a small number of samples for many miRNAs
 - QRT-PCR methods are ideal for:
 - Profiling a large number of samples against a small number of miRNAs
 - Profiling a moderate number of samples against a moderate number of miRNAs

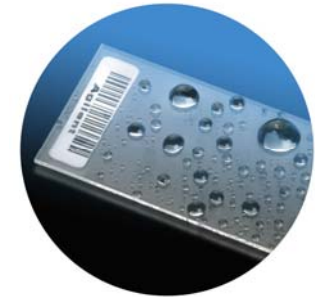




Choose from a comprehensive set of microarray applications

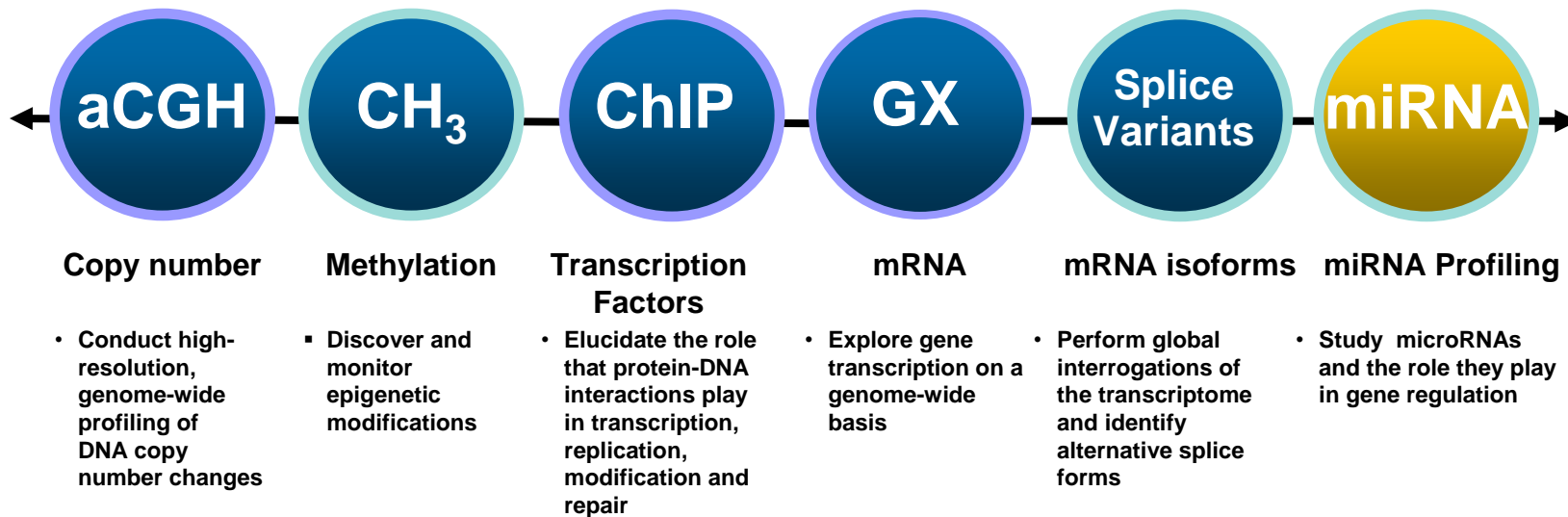
From one-dimensional to multi-dimensional ...

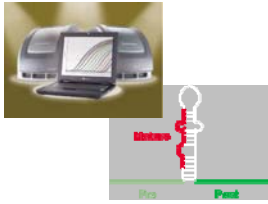
with multiple applications ...



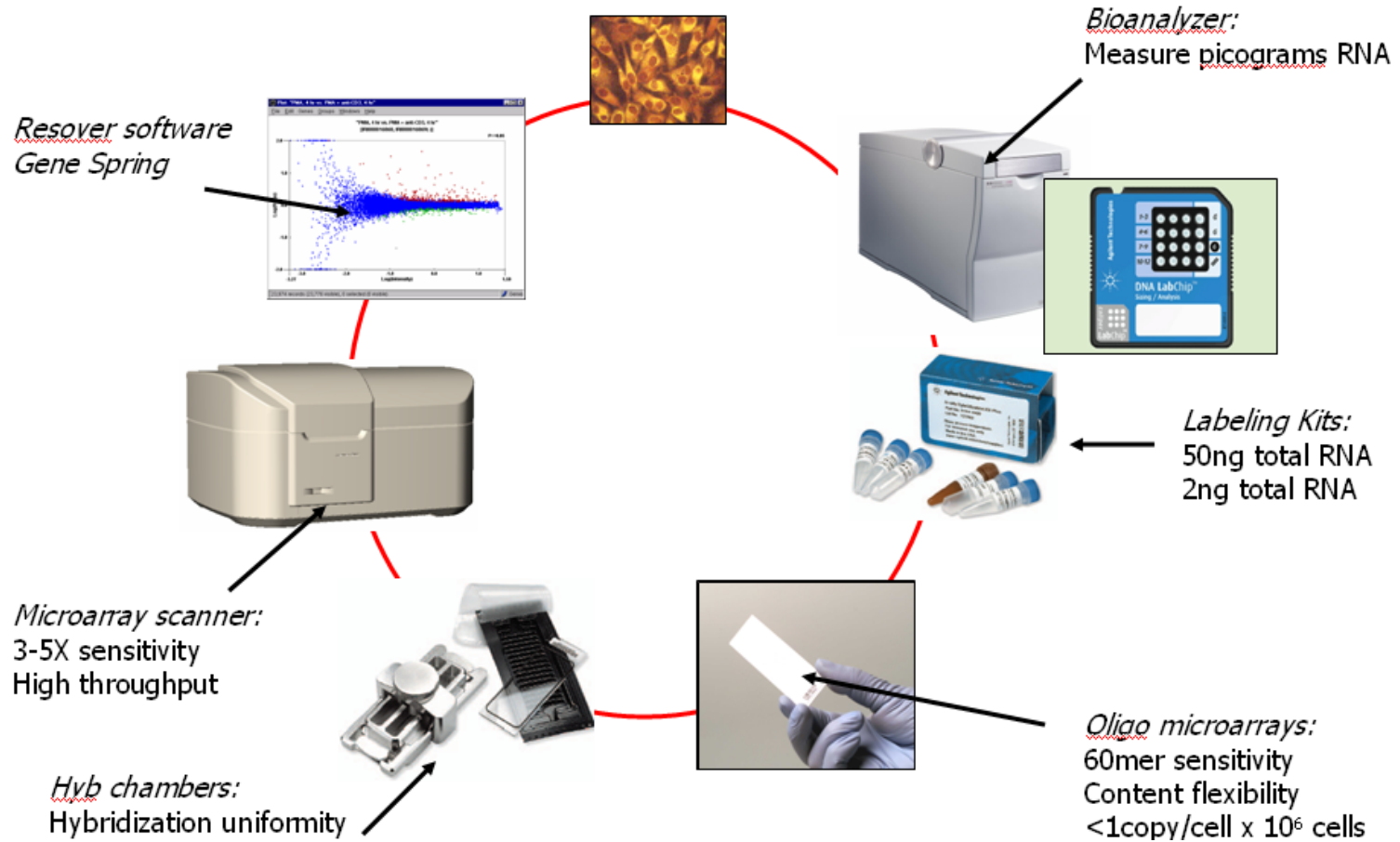
DNA

RNA

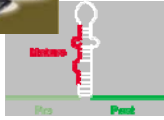




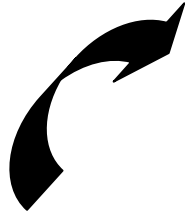
Microarray work flow



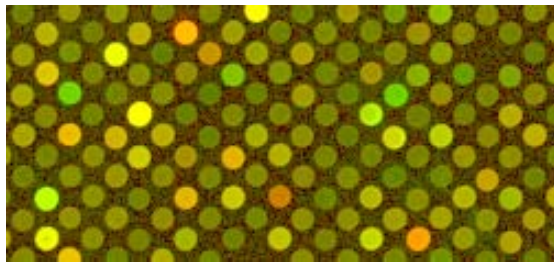
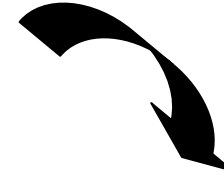
Data Workflow



Grid Template
Grid File
FE Protocol



Grid and Measure Spots
Reject Outlier Pixels
Subtract Background
Correct Dye Biases



16-bit Tiff Image
(uncompressed)



Feature Extraction Software



CGH analytics



GeneSpring



Chip analytics



QC Report
Grid
Shape
Text
JPEG
MAGE-ML
GEML

Feature Extraction
Result Files

Import

FTP
Export

Agilent and non-Agilent microarrays
scanned on [Agilent Scanner](#)

Agilent microarrays scanned on
[GenePix Scanner](#)





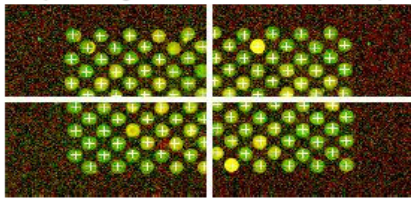
QC Metrics

Feature Extraction QC metrics

QC Report - Agilent Technologies : 2 Color CGH

Date:	Thursday, February 23, 2006 - 11:12	BG Method:	No Background
Image:	US45102826_251275011095_S01	Spatial Detrend:	On
Protocol:	CGH_44k_1005 (Read Only)	Global Adjust:	Off
User Name:	ikishawi	Dye Norm:	Linear
Grid:	012750_D_20050614	Linear DyeNorm Factor:	18.2 (Red) 28.4(Green)
FE Version:	8.5.1.1	Additive Error:	33(Red)104(Green)
		Derivative of Log Ratio Spread:	0.42

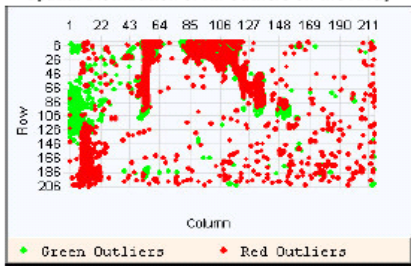
Spot Finding of the Four Corners of the Array



Feature	Local Background	
	Red	Green
Non Uniform	102	53
Population	39	82

Feature	Local Background	
	Red	Green
Non Uniform	4	4
Population	1877	2566

Spatial Distribution of All Outliers on the Array



Local Bkg (inliers)

Net Signal statistics

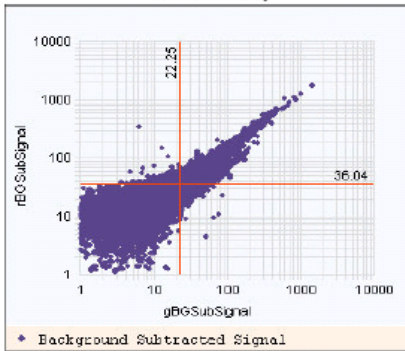
Non-Control probes:

	Red	Green
NumSat	0	1
99%	271	220
50%	68	68
1%	34	37

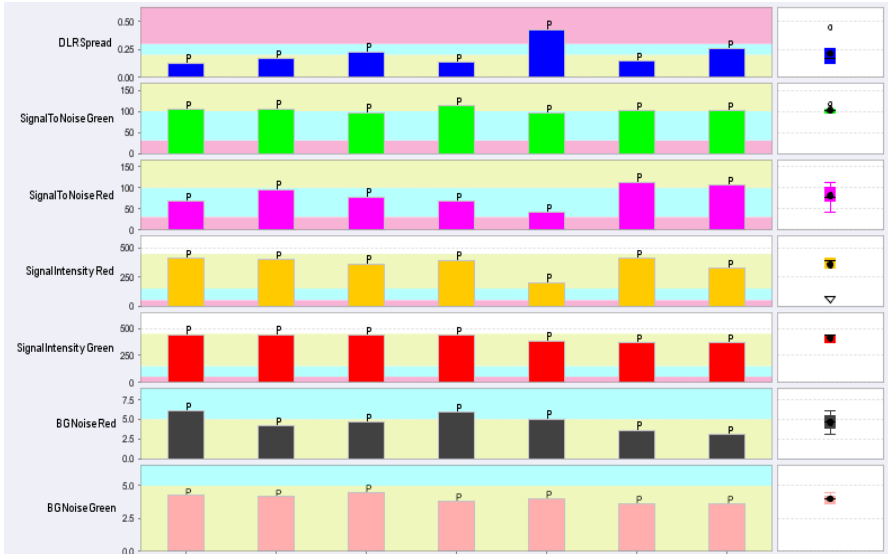
Negative Control Stats

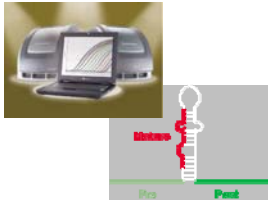
	Red	Green
Average Net Signals	32.36	45.92
StdDev Net Signals	1.93	4.45
Average BG Sub Signal	-0.03	-0.08
StdDev BG Sub Signal (BG Noise)	1.83	3.67

Red and Green Background Corrected Signals (Non-Control Inliers)



DNA Analytics QC metrics





Data quality: feature sizes and density

- Features should be far enough apart to prevent light leakage.
- Features should be big enough for perfect registration of each layer (no blurry edges), for individual feature statistics and higher confidence
- An ability to increase density while maintaining large enough features without compromising statistical significance

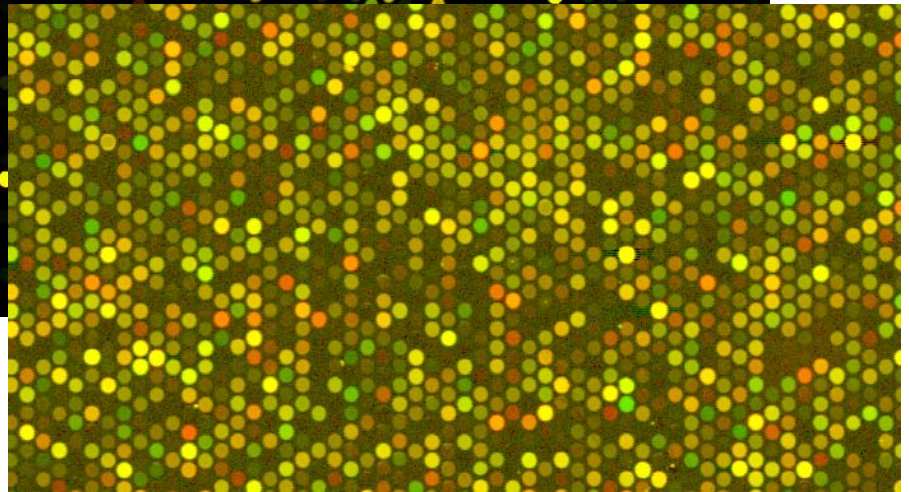
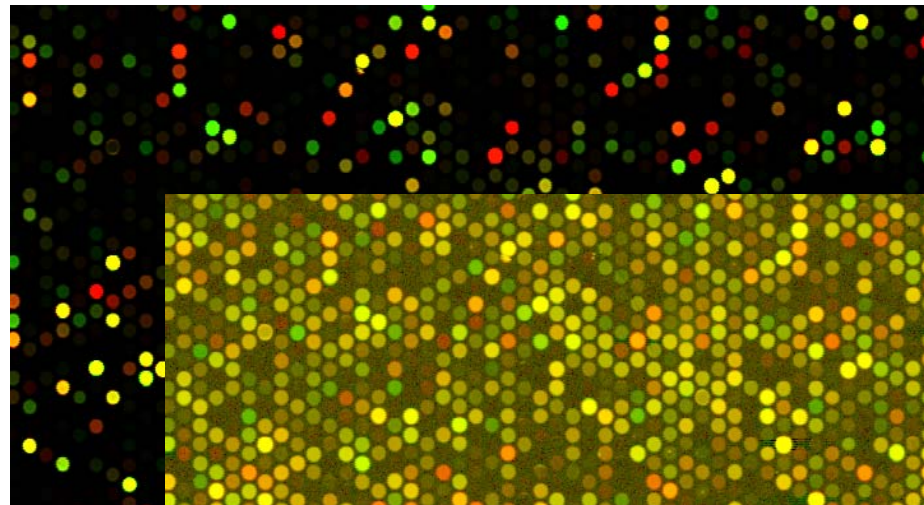
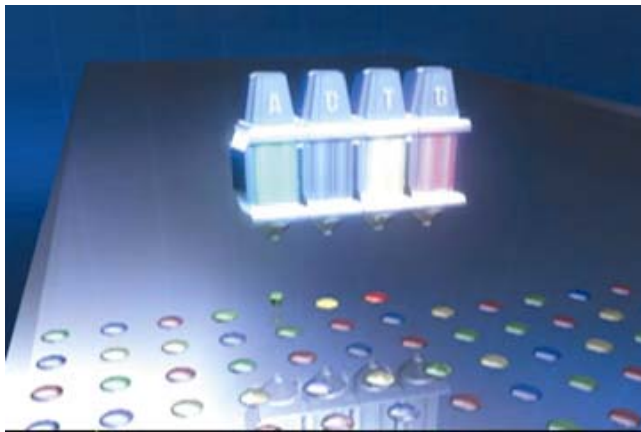
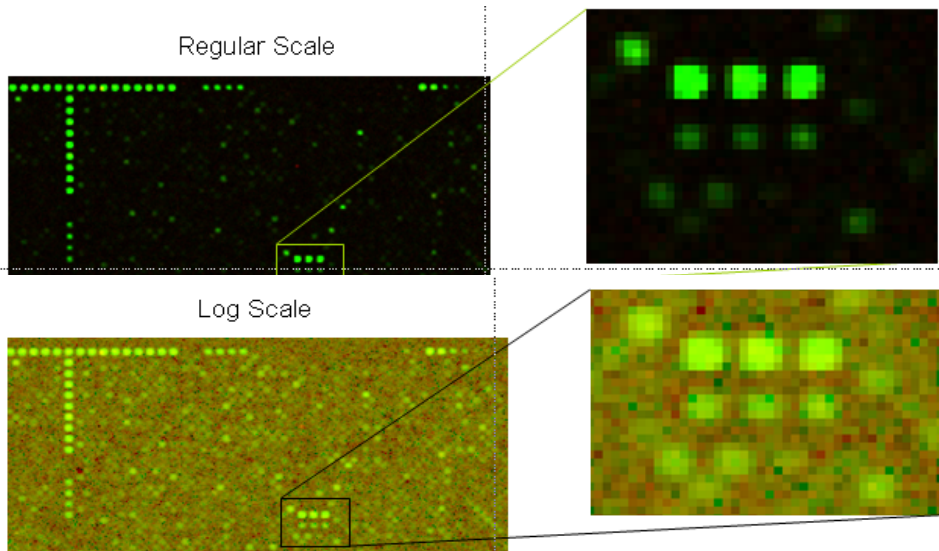


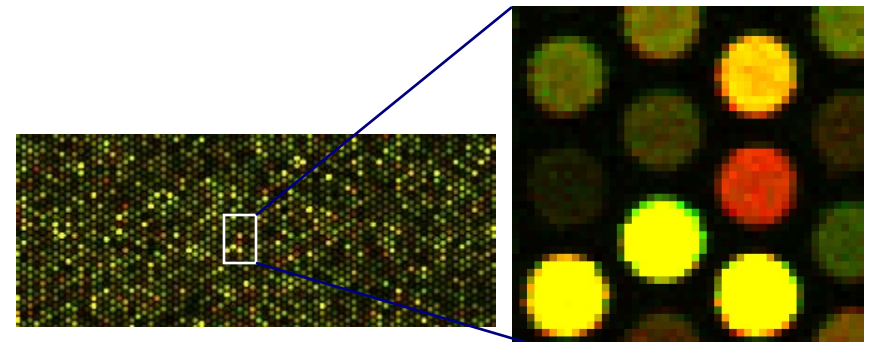


Image Value in Data quality

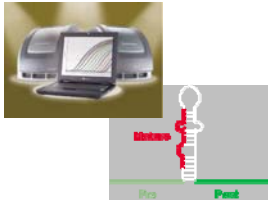
Agilent scanner at 5 micron resolution



Other array image

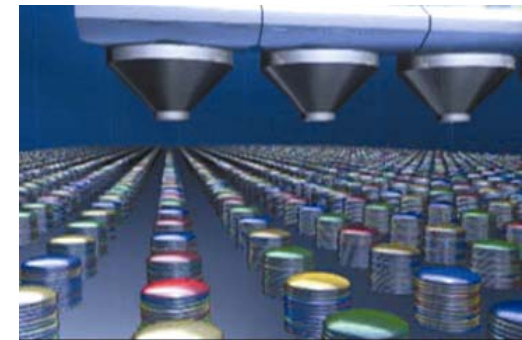
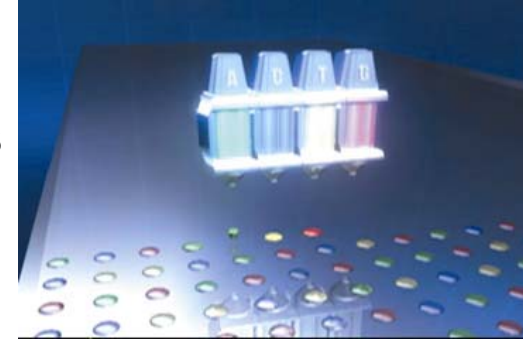


Agilent 224K image



Agilent's Microarray Platform

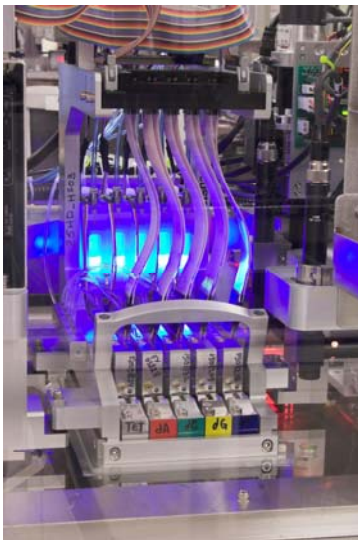
- **Reliable inkjet printing of phosphoramidite bases**
- **Sensitivity:** The chemistry reliability allows synthesis of long oligos that are highly sensitive and specific
- **Flexibility of microarrays, array is defined by electronic file.**
- **Ease of implementation**
- **Probe Design:** Narrow TM, Unique Probes, Self Structure, and GC content, All probes are empirically tested.
- **Probe Fidelity** can synthesize up to 200mer





Santa Clara Manufacturing Facility

- Manufacturing Process Development
- Bioinformatics



- Industrial manufacturing – Class 10,000 clean-room
- Wired directly into eArray, allowing direct customer access to fully customizable products
- High-performance inkjet printing enables long oligo manufacturing

Our measure is your success.



Agilent Technologies



Challenges in miRNA Profiling

- Small size – difficult to label with high efficiency
- High sequence homology – difficult to design probes with high specificity
- Presence of larger RNAs with highly homologous sequences
- Expressed with large dynamic range – difficult to avoid compression
- Growing & changing database

miRNA	Sequence	#NT
hsa-let-7a	ugagguaguagguuguauagu	22
hsa-let-7b	ugagguaguagguugu <u>g</u> guu	22
hsa-let-7c	ugagguaguagguuguau <u>g</u> guu	22
hsa-let-7d	<u>a</u> gagguaguagguug <u>c</u> auagu	22
hsa-let-7e	ugagguag <u>g</u> agguuguauagu	22
hsa-let-7f	ugagguaguag <u>a</u> uuguauagu	22
hsa-let-7g	ugagguaguag <u>u</u> uugua <u>c</u> agu	22
hsa-let-7i	ugagguaguag <u>u</u> uugu <u>gcu</u> guu	22

Sanger
miRBase 12.0



miRNA Arrays:

- Human miRNA Microarray, v2.0:
 - - 799 distinct probe sets (723 human and 76 human viral)
- Human miRNA Microarray, v1.0:
 - - 450 human distinct probe sets





Direct and Sensitive miRNA Profiling

www.rnajournal.org

RNA

VOL. 13, NO. 1 MAY 2007

In this issue:

- tRNA competition and translational misreading
- Conversion of pre-RISC to holo-RISC by Ago2
- Protein composition of human miRNAs spliced in vitro
- Role of metal ions in the tetraloop-receptor complex
- Direct and sensitive miRNA profiling

CSH PRESS COLD SPRING HARBOR, NY




Hui Wang, PhD
Senior Scientist
Agilent Technologies

METHOD

Direct and sensitive miRNA profiling from low-input total RNA

HUI WANG, ROBERT A. ACH, and BO CURRY
Agilent Technologies, Inc., Agilent Laboratories, Santa Clara, California 95051, USA

ABSTRACT

We have developed a sensitive, accurate, and multiplexed microRNA (miRNA) profiling assay that is based on a highly efficient labeling method and novel microarray probe design. The probes provide both sequence and size discrimination, yielding in most cases highly specific detection of closely related mature miRNAs. Using a simple, single-vial experimental protocol, 120 ng of total RNA is directly labeled using Cy3 or Cy5, without fractionation or amplification, to produce precise and accurate measurements that span a linear dynamic range from 0.2 amol to 2 fmol of input miRNA. The results can provide quantitative estimates of the miRNA content for the tissues studied. The assay is also suitable for use with formalin-fixed paraffin-embedded clinical samples. Our method allows rapid design and validation of probes for simultaneous quantitative measurements of all human miRNA sequences in the public databases and to new miRNA sequences as they are reported.

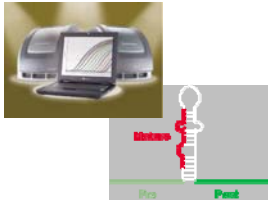
Keywords: miRNA profiling; microarray; RNA labeling; probe design; microRNA

INTRODUCTION

miRNAs (miRNAs) are a class of small single-stranded coding RNAs [19–30 nucleotides (nt)] that have been identified in animals, plants, and viruses, with over 400 identified in humans (Bartel 2004; Griffiths-Jones 2004; Kore and Haley 2005; Griffiths-Jones et al. 2006; Kim Nam 2006). MicroRNA genes are transcribed as pri-miRNAs, which are then processed to the shorter hairpin and pre-miRNAs (~70–90 nt) before they are cleaved to the mature single-stranded miRNAs (~22 nt) (Bartel; Zamore and Haley 2005; Kim and Nam 2006). As a multiprotein RNA-induced silencing complex, the miRNAs repress translation by forming imperfect base pairing with the 3' untranslated region of target messenger RNAs (mRNAs). Bioinformatics and cloning studies have estimated that miRNAs may regulate 30% of all human genes (Lencioni et al. 2005; Lim et al. 2005). Recent studies have shown that distinct miRNA expression patterns are associated with various cancer types (Cahn et al. 2004; Liu et al. 2005; Cassanese et al. 2006; Esquela-Kerscher and...

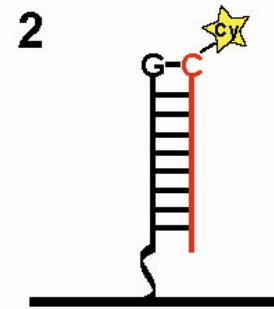
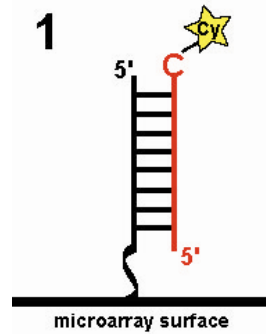
The ardent interest in profiling miRNA expression has led to developments in Northern blot (Cummins et al. 2006), cloning (Lau et al. 2001), PCR (Chen et al. 2005), bead-based (Lu et al. 2005), SAGE-based (Cummins et al. 2006), and microarray-based methods (Babak et al. 2004; Band et al. 2004; Galin et al. 2004; Liu et al. 2004; Nelson et al. 2004; Baskerville and Bartel 2005; Liang et al. 2005; Lim et al. 2005; Shingara et al. 2005; Castoldi et al. 2006). The ideal method would require submicrogram quantities of total RNA, have minimal sequence bias, be easy to multiplex, and involve simple experimental protocols. Microarray-based assays offer an efficient method for measuring the expression profile of large numbers of miRNAs simultaneously. However, the small size and high sequence homology of miRNAs present major challenges to sample labeling and microarray probe design. Here we present a novel miRNA assay employing simple high-efficiency direct labeling of submicrogram quantities of total RNA, without amplification or size fractionation. Our labeling protocol has little sequence bias, and our in situ synthesized DNA microarray probes (Hughes et al. 2001) are both sequence and size selective for mature miRNAs. The enzymatic labeling attaches a single fluorophore-labeled nucleotide to the 3' end of each miRNA with high yield and minimal sample manipulation. Hybridization to the microarray is carried out under conditions that result in near-equilibrium binding and high (>25%) hybridization yields for most miRNAs. The assay is easy to perform, has





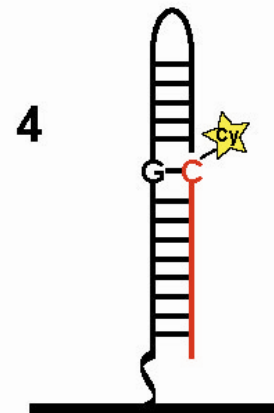
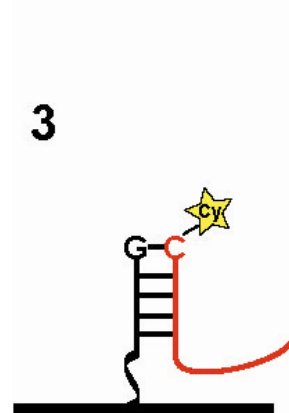
miRNA Probe Design Strategy

1. Start design with full-length miRNA-probe sequence, attached to a stilt sequence.



2. Utilize the **C** incorporated during labeling for additional **G-C** base pair on 3' end of miRNA to increase stability

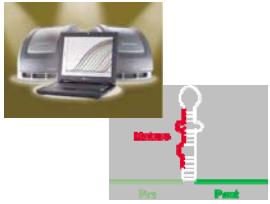
3. Sequentially shorten target-probe base pairing from 5' end of miRNA during preliminary T_m balancing by calculation.



4. Incorporate hairpin structure on probes to increase size specificity and probe:target stability.

FINAL STEP:

Select T_m -balanced probes for each miRNA empirically using microarray data.



Work Flow

Total RNA
(100 ng)

Direct Label with Cy-dye

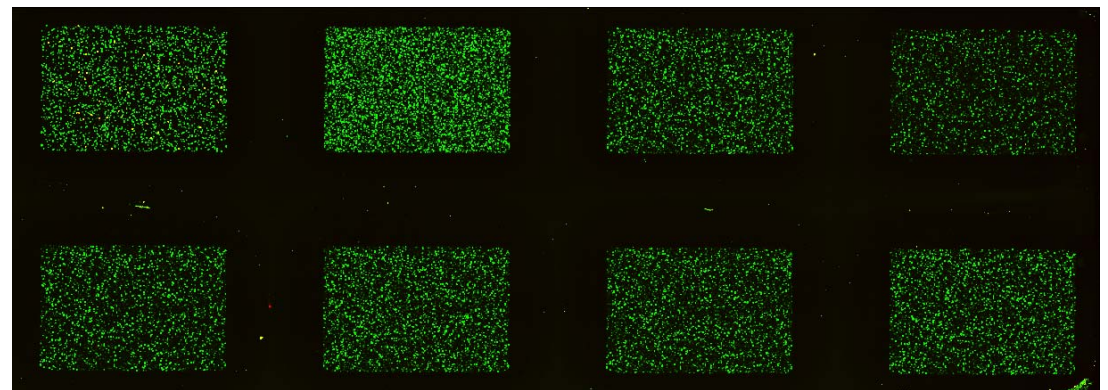
Labeled RNA

Hybridize

miRNA Profile

8-pack

100ng sample input
Direct probe labeling
High specificity & sensitivity





Specificity to Distinguish Homologous

miRNAs

	7a	7b	7c	7d	7e	7f	7g	7i
7a	100	10	51	3	1	5	2	0
7b	0	100	1	0	0	0	7	0
7c	6	70	100	1	0	0	3	0
7d	1	2	1	100	39	0	3	0
7e	4	1	1	2	100	0	1	0
7f	62	3	5	1	0	100	1	0
7g	1	0	0	0	0	0	100	1
7i	0	1	0	0	0	0	4	100

85 - 100%

70 - 84%

55 - 69%

40 - 54%

25 - 39%

10 - 24%

5 - 9%

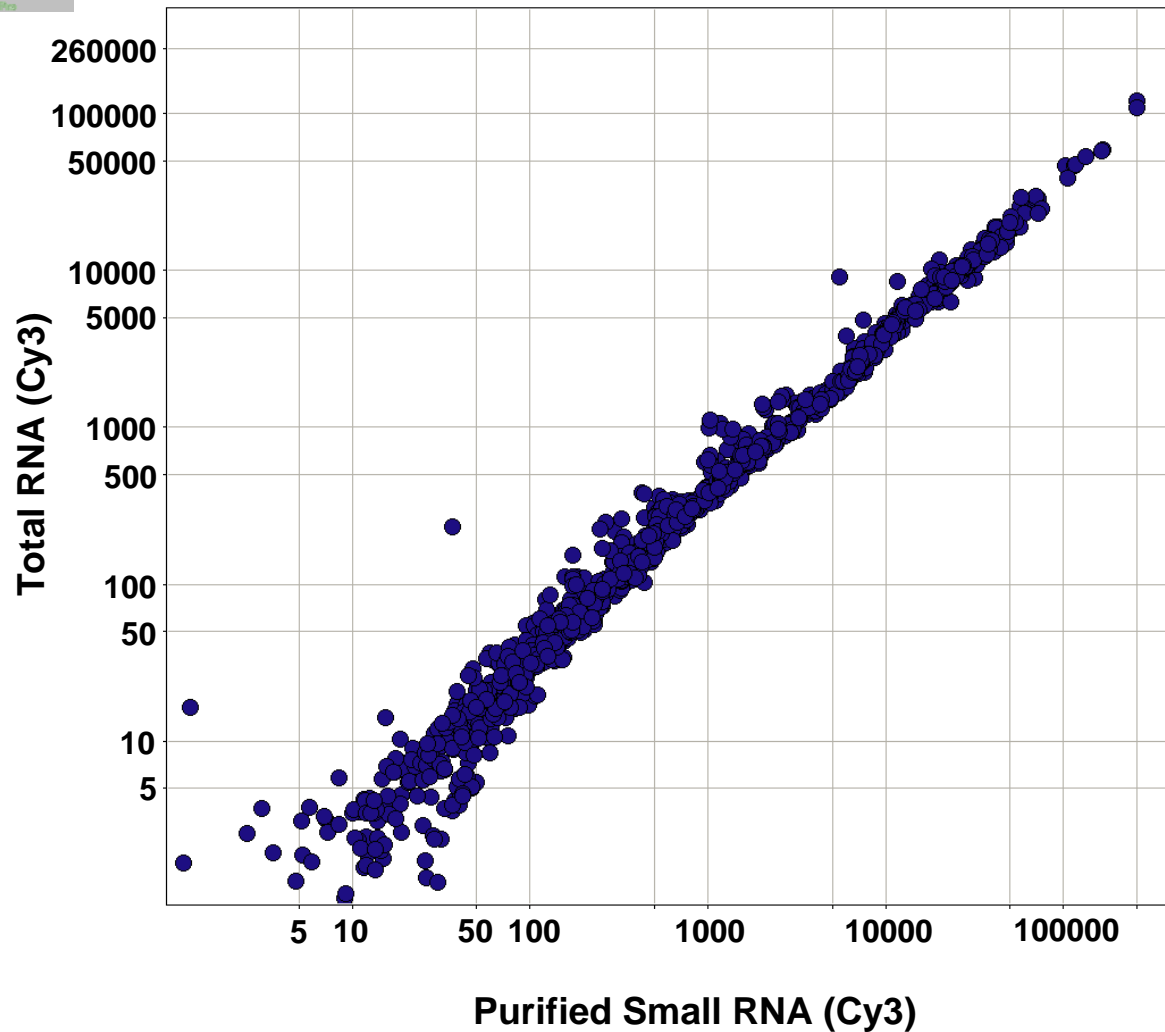
0 - 4% Grey Number

hsa-let-7 family

miRNA	Sequence	Length (nts)
hsa-let-7a	UGAGGUAGUAGGUUGUAUAGUU	22
hsa-let-7b	UGAGGUAGUAGGUUGUGUGGUU	22
hsa-let-7c	UGAGGUAGUAGGUUGUAUGGUU	22
hsa-let-7d	AGAGGUAGUAGGUUGCAUAGU	21
hsa-let-7e	UGAGGUAGGAGGUUGUAUAGU	21
hsa-let-7f	UGAGGUAGUAGAUUGUAUAGUU	22
hsa-let-7g	UGAGGUAGUAGUUUGUACAGU	21
hsa-let-7i	UGAGGUAGUAGUUUGUGCUGU	21



Direct miRNA Measurement From Total RNA



Data shown are background-subtracted signals with no filtering or normalization.

Each miRNA has signals from multiple probes.

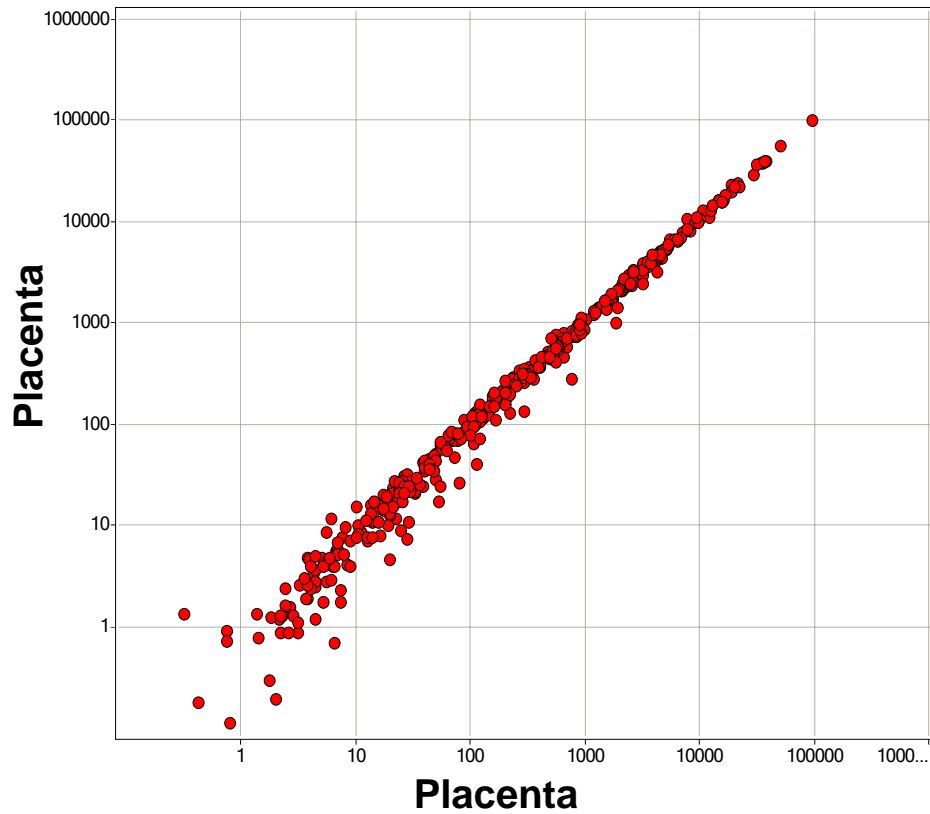




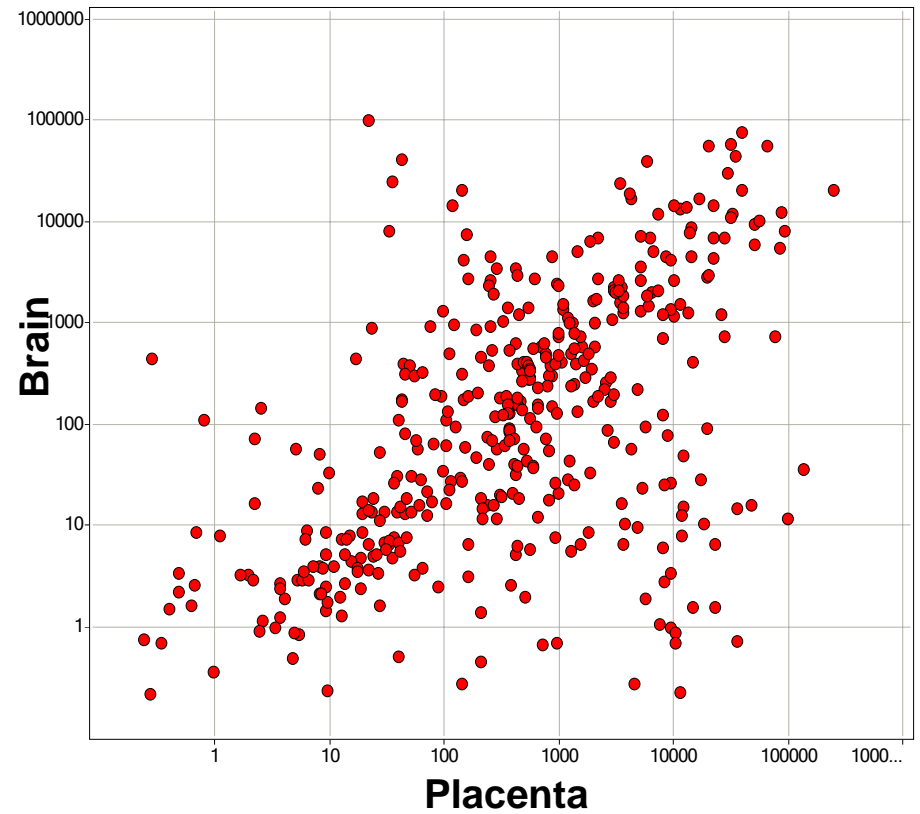
miRNA Profiles Using 100ng Total RNA



Reproducibility

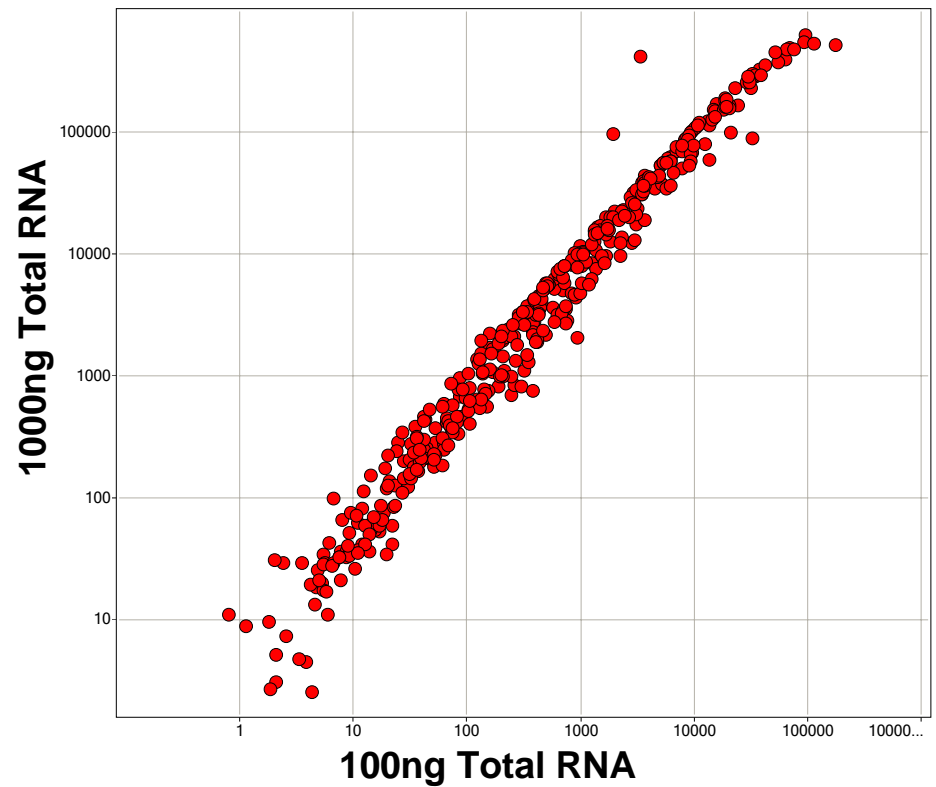
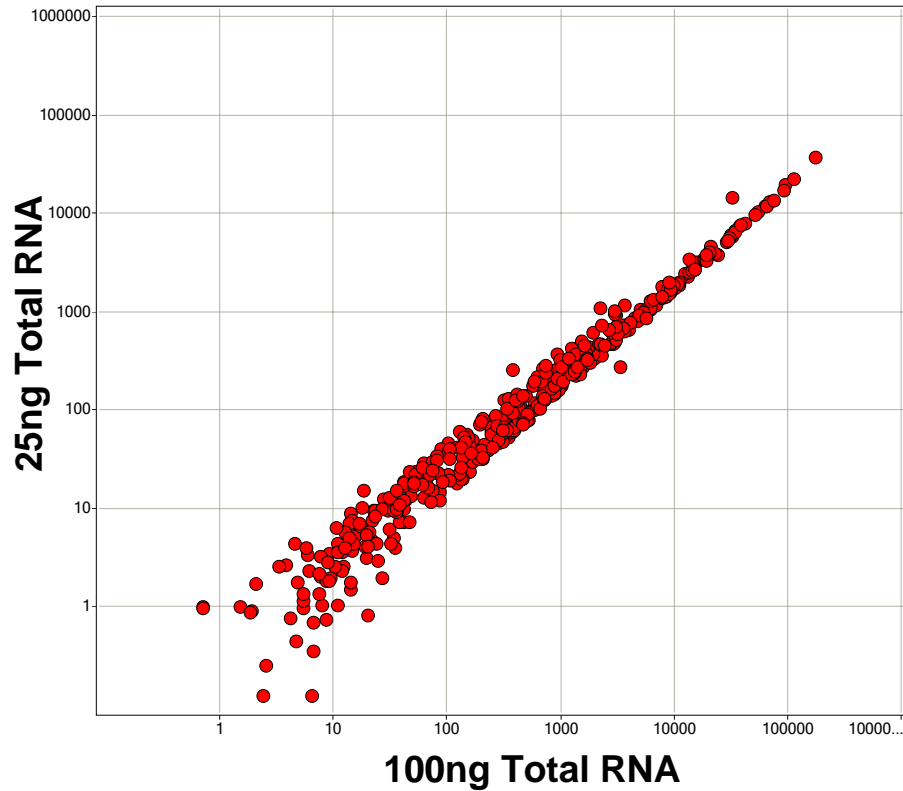


Differential Expression





miRNA Profiles Independent of Sample Input



Consistent miRNA profile for low (25ng) vs high (1000ng) total RNA input

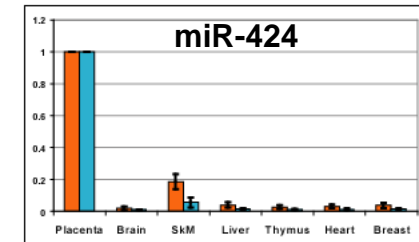
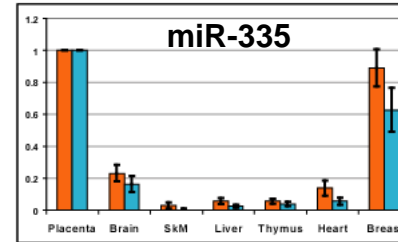
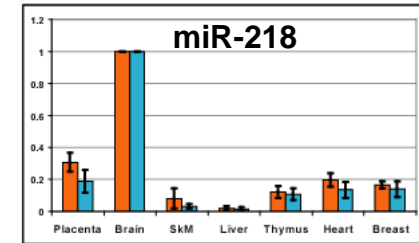
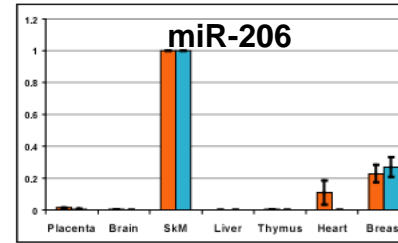
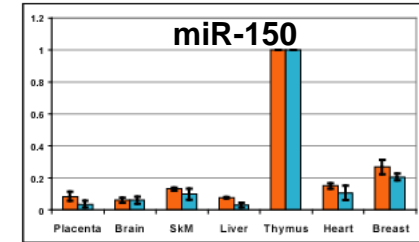
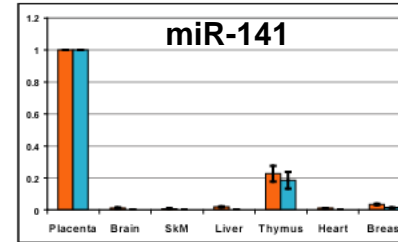
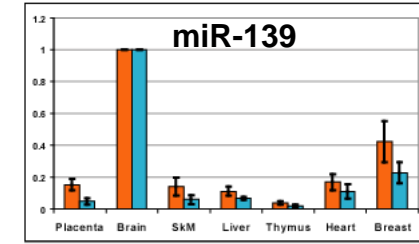
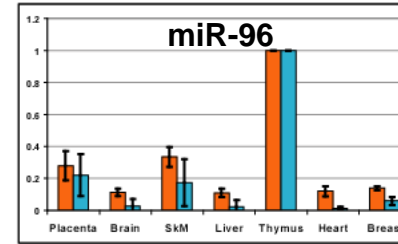
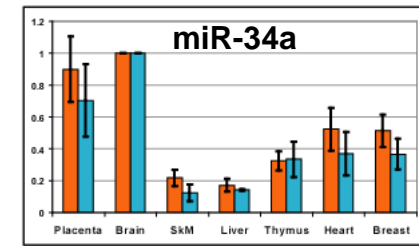
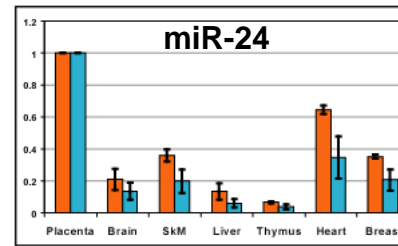




Correlation between microarray and quantitative RT-PCR results

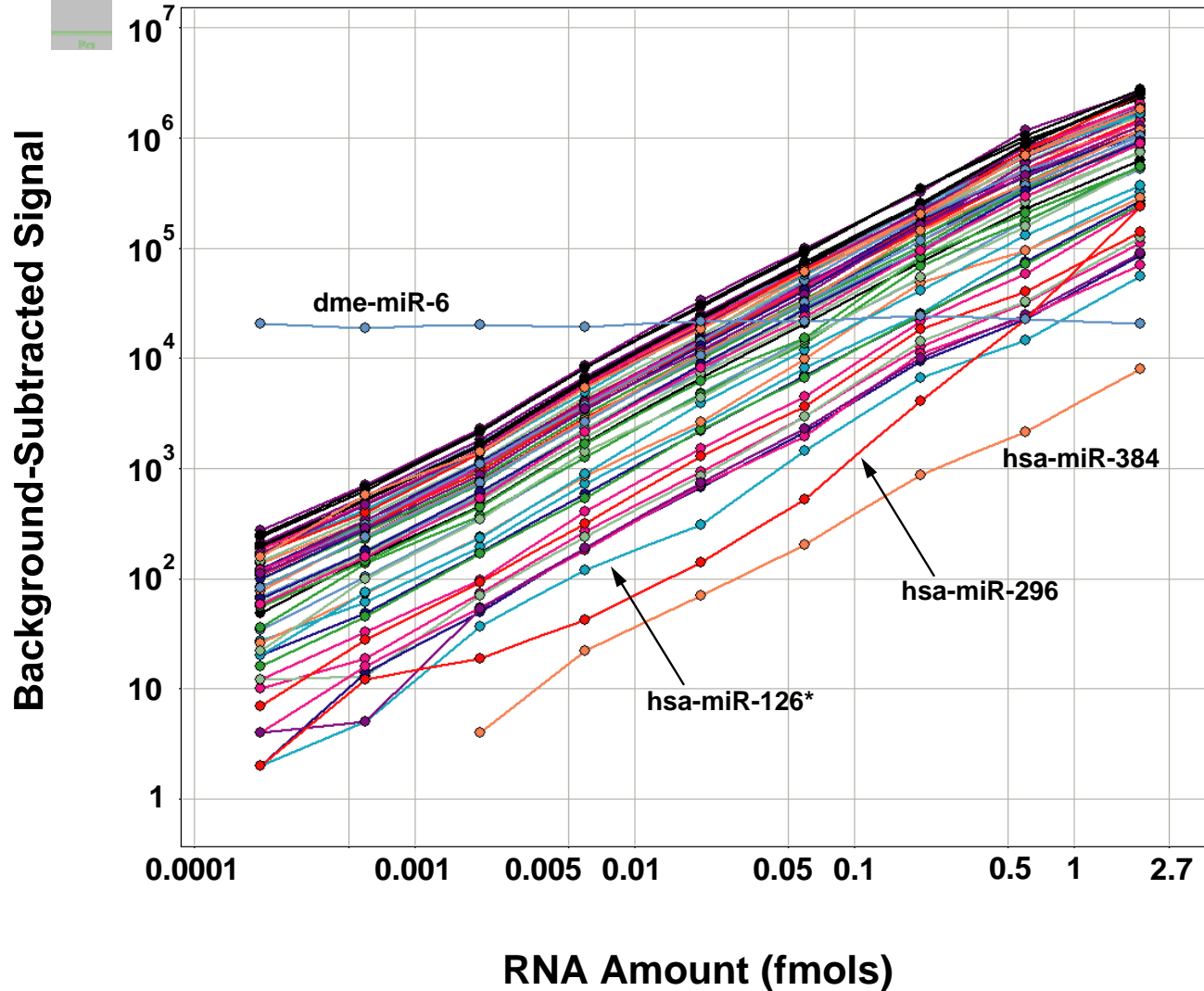
Quantitative RT-PCR (qPCR) reactions (orange bars) and microarray hybridizations (blue bars) were run in quadruplicate for each miRNA species.

 qPCR Data  Microarray Data





10⁴ Linear Dynamic Range



Synthetic miRNAs were selected for typical and atypical sequence representations.

Lowest and highest calculated Tms are represented.

Most miRNAs are detectable at the 0.2 α mol-2fmol range.





Platform Features, Benefits & Advantages

Feature	Benefit	Advantage to Investigator
8 x 15K format	Process multiple samples on the same slide	Reduce costs Run parallel experiments
Optimized probe design and direct labeling	Greater sensitivity and specificity	Detect low abundance miRNAs Detect highly homologous miRNAs
Small sample input	Ability to work with precious samples	Analyze new sample types Conserve precious samples
Use of total RNA	Does not require size separation	Does not introduce any miRNA isolation bias Easy to use Reduce hands-on time
Broad dynamic range	Detect all possible miRNAs in sample	Comprehensive miRNA profiling
Compatible with standard microarray platform	Correlate miRNA profiles with GE and CGH data	Requires only single capital investment Enables integrated genomics



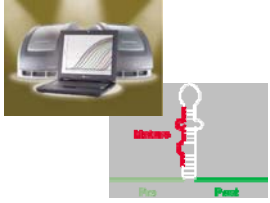


Response from Early Access Customer

We have shown for the first time that let-7 expression is frequently reduced in lung cancers and that alterations in the miRNA expression may have a prognostic impact on the survival of surgically treated lung cancer patients. Agilent miRNA arrays give us the comprehensive miRNA expression profile with excellent performance on sensitivity and accuracy. I expect that the studies of Agilent miRNA array may ultimately provide a foundation for a new paradigm of the involvement of miRNA in human oncogenesis.

Dr. Takashi Takahashi
Professor of Oncology
Molecular Carcinogenesis
Nagoya University





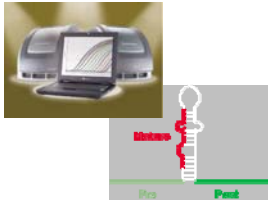
Stratagene QPCR MicroRNA Studies

Featuring:
High-Specificity miRNA QRT-PCR
Detection Kit

Our measure is your success.



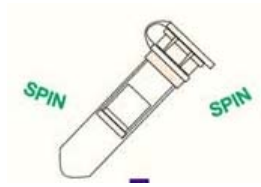
Agilent Technologies



Purification of miRNA

Challenges in small RNA Purification

Spin columns ➤ Retention of small RNAs on standard columns not sufficient: → Low yield of small RNAs



➤ Specialized solutions available:
miRACLE™ miRNA Isolation kit

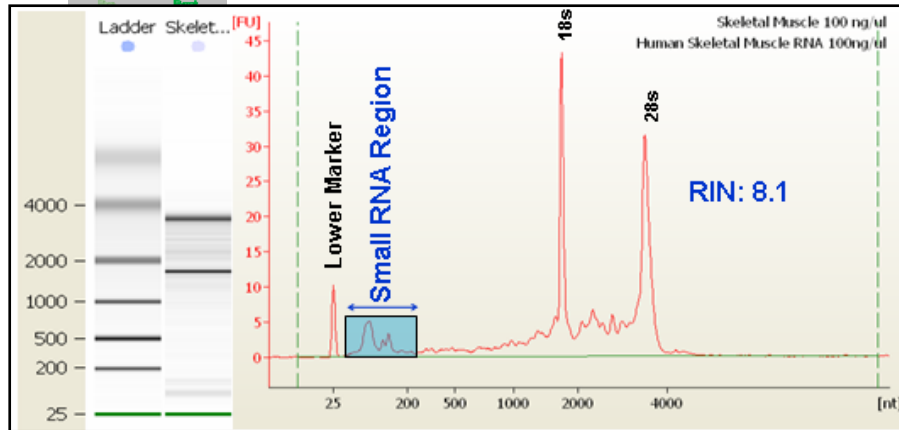
- Enrichment of miRNA species for increased sensitivity using specialized solutions like miRACLE™
- Usually Phenol-Chloroform based clean up to avoid loss of small RNAs
- Alternative methods like SideStep™ (Cells-to-PCR) available





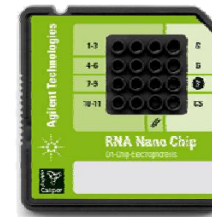
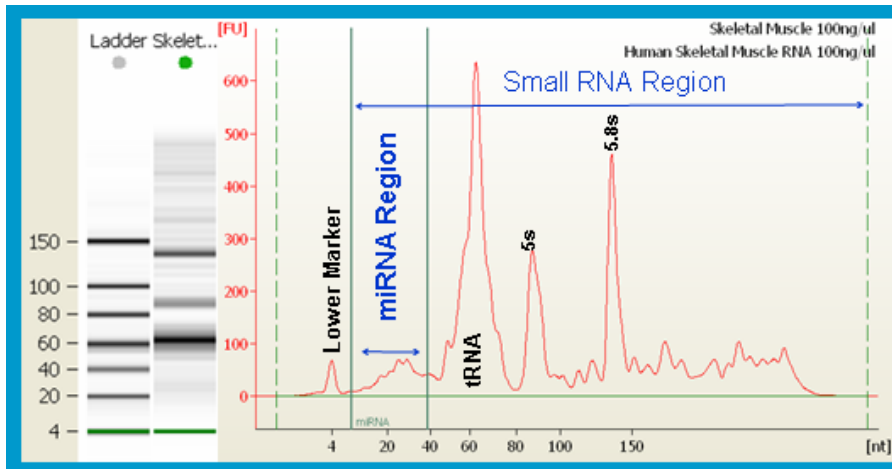
Purification of miRNA

Sample Verification – Bioanalyzer Small RNA kit



RNA 6000Nano kit

Small RNA Kit



RNA 6000 Nano Kit

Size range: 25-6000nt

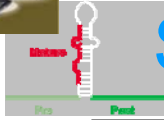
Results: Integrity, Total RNA amount, gDNA contamination



NEW! Small RNA Kit

Size range: 6-150nt

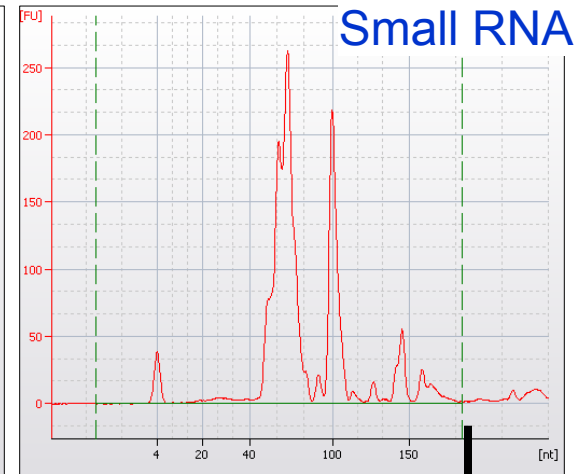
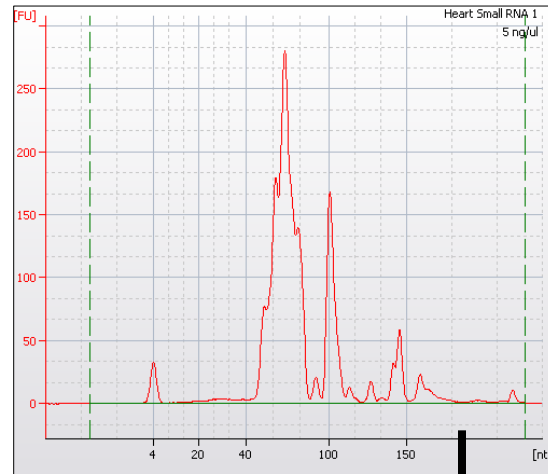
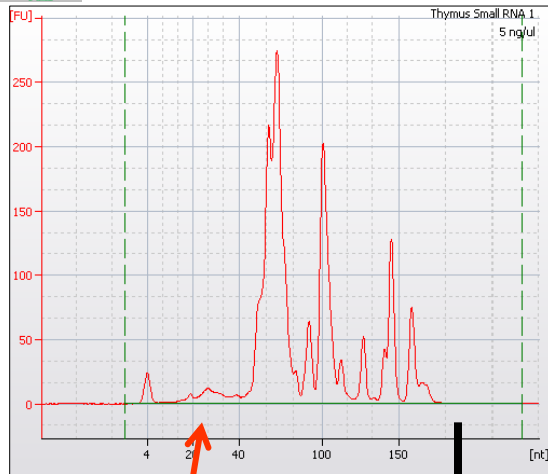
Results: miRNA amount, Ratio and amount of other Small RNA



Purification of miRNA

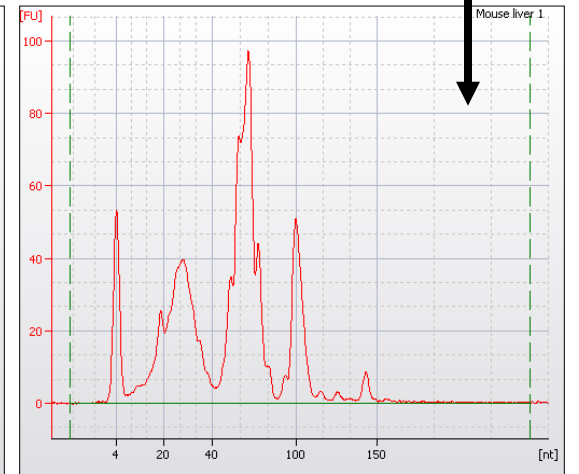
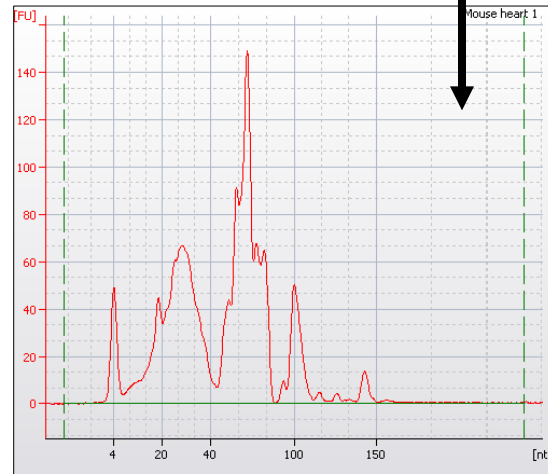
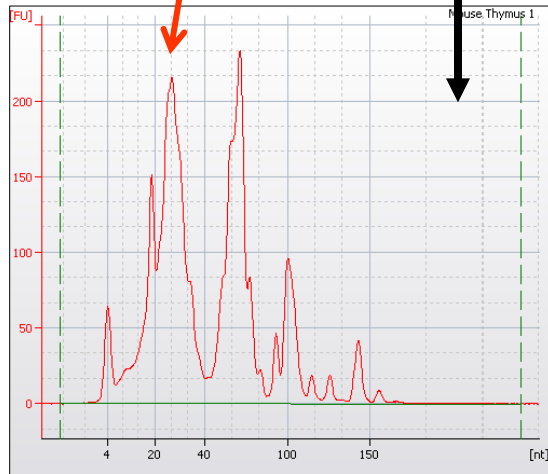
Sample Verification – Bioanalyzer Small RNA kit

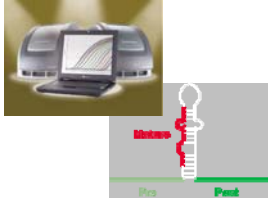
Pre-purified
small RNA
(0-150 nt)



miRNA Peak

enriched
miRNA
fraction



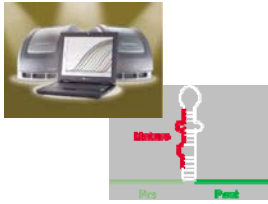


miRNA Detection Challenges in QPCR

Design challenges:

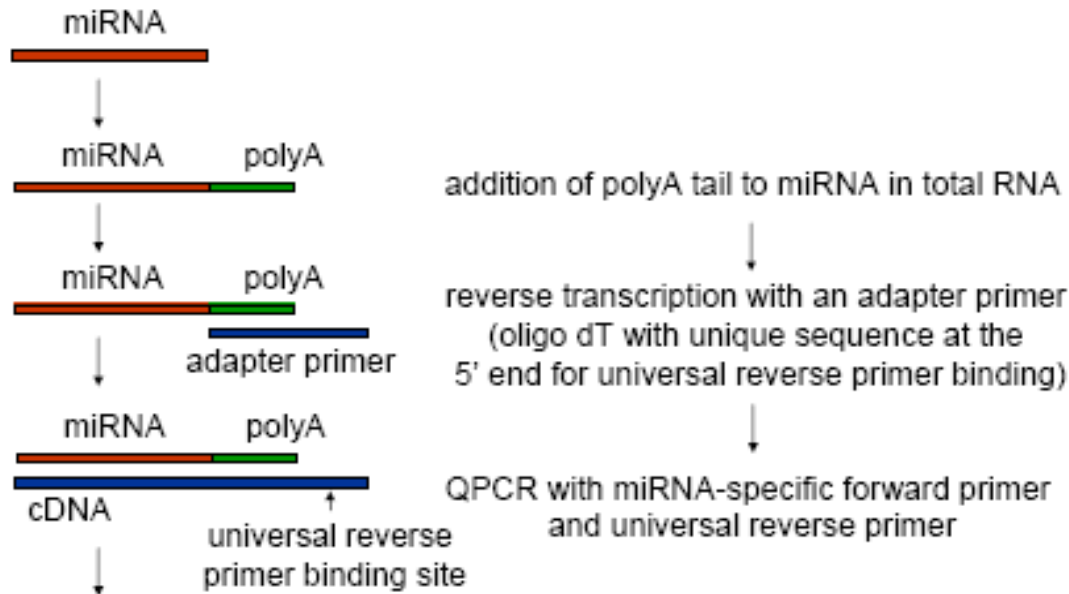
- Sequence length of miRNAs is very short
- Lack of common nucleotide sequence to be used as priming site
- High sequence homology between different miRNAs
- Presence of different forms of miRNA in a cell



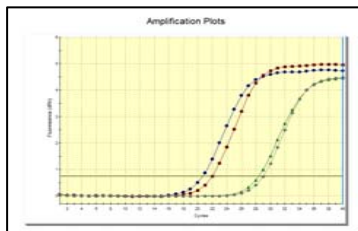


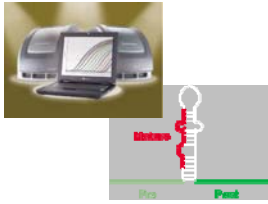
QRT-PCR Detection

High-Specificity miRNA QRT-PCR Detection Kit



- High-Specificity miRNA QRT-PCR Detection Kit
 - miRNA 1st Strand cDNA Synthesis Kit
 - High-Specificity miRNA QPCR Core Kit





Specificity of High-Specificity miRNA QRT-PCR Detection Kit

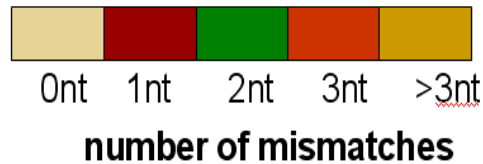
Discrimination between all let-7 family members using miRNA-specific PCRprimers.

miRNA-specific assay

miRNA synthetic templates

	let-7a	let-7b	let-7c	let-7d	let-7e	let-7f	miR-98	let-7g	let-7i
let-7a	100	0	0	0	0.7	0.2	0	0	0
let-7b	0.0	100	0.6	0	0	0	0	0	0
let-7c	0.8	0.1	100	0	0	0	0	0	0
let-7d	0.1	0	0	100	0	0	0	0	0
let-7e	0.6	0	0	0	100	0	0	0	0
let-7f	0.2	0	0	0	0.1	100	0	0	0
miR-98	0	0	0	0	0	0	100	0	0
let-7g	0	0	0	0	0	0	0	100	0
let-7i	0	0	0	0.1	0	0.1	0	0.1	100

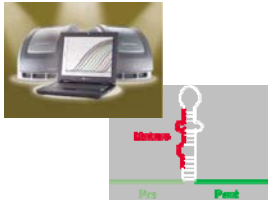
% relative detection



➤ The same number of let-7 miRNA (10^{10} miRNA synthetic templates) were converted to cDNA and detected with each of the let-7 miRNA-specific primers (miRNA-specific assay).

- Detection with each of the let-7 miRNA-specific primers (miRNA-specific assay)
- Copy number of the matching primer and template was 100%.
- Copy number of a non-matching primer and template determined as a percentage of the matching primer and template (% relative detection)
- The number of mismatched nucleotides is indicated in color as shown in the key (number of mismatches).





High-Specificity miRNA QRT-PCR Detection Kit Features

- Detects mature miRNA in 3 hours
- As little as 15 ng of total RNA input
- Variety of sample types
 - Lysed cells
 - Total RNA isolated from tissues
 - Cell cultures
 - FFPE tissues
- Sensitive detection down to 10 copies
- Single-nucleotide discrimination
- Detection of up to 6,000 different miRNA in a single sample preparation





High-Specificity miRNA QRT-PCR Detection Kit Features

Feature	High-Specificity miRNA QRT-PCR Detection Kit
Sensitivity	33 copies in 3.3ng total RNA (10 copies/ng)
Template (lowest)	30pg total RNA/600-6,000 miRNA
Template (recommended)	250-1000ng/600-6,000 miRNA
Linearity	7 logs
# human assays	50 human





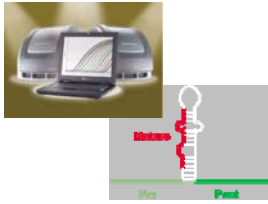
High-Specificity miRNA Detection Kit

- 50 miRNA-specific primers based on human miRNA
- selection was based on differential expression in cancer and during development most will also detect mouse and rat miRNA

human	mouse	rat
50	41 (miR-106a)	40 (miR-15a, miR-106a)

- will be adding another 8 miRNA-specific primers to the list
 - human and mouse U6 primers for normalization
- primer design rules will be available to customers on our website
lot number from RT adapter primer has to be input to access primer design rules

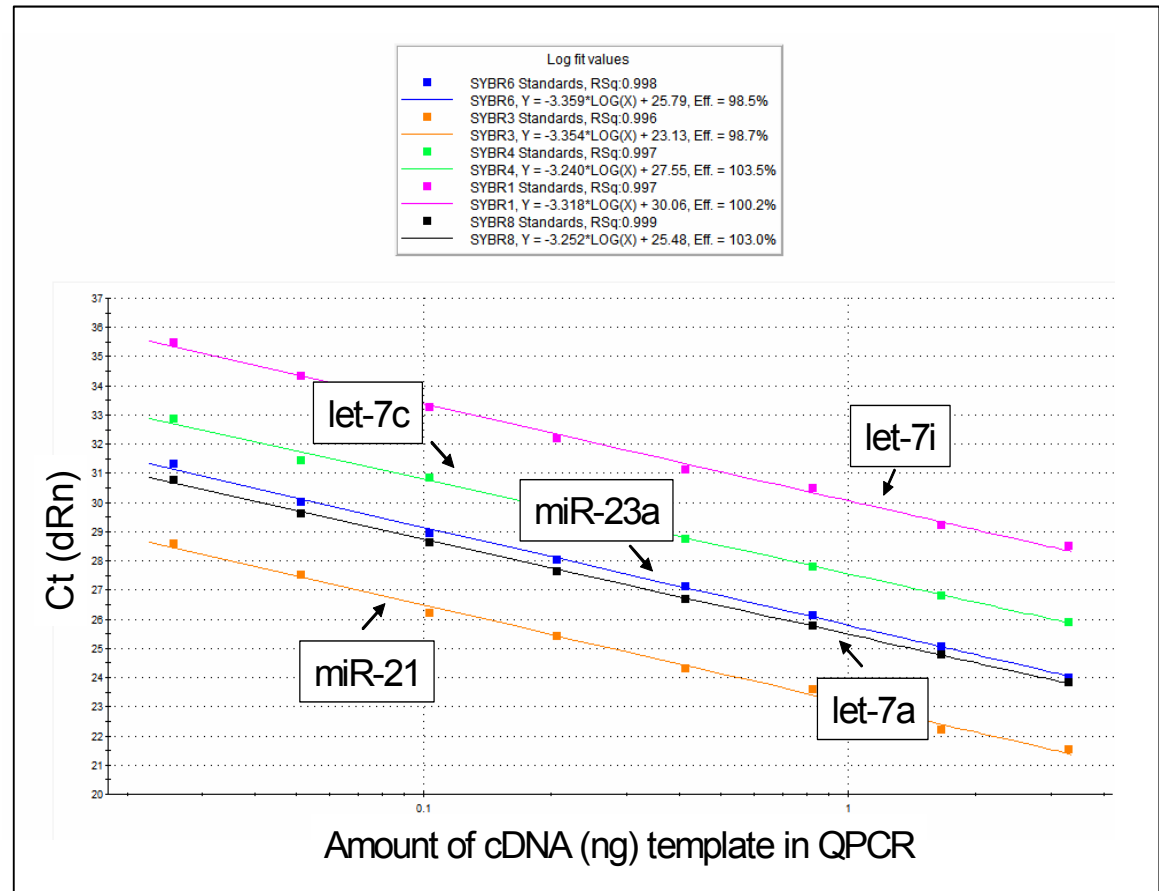




Kit Performance: Assay Linearity

➤ cDNA prepared from 1ug HeLa total RNA and using varying amounts as template in QPCR.

➤ Five different miRNA of varying abundance, let-7a, let-7c, let-7i, miR-21 and miR-23a, were detected.



- 25 pg – 3.3 ng cDNA input. Detection of miRNAs of varying abundance
- Linearity over 7 logs (after Poly A tailing and cDNA synthesis)



Detection of miRNA Expression

Breast Cancer Biomarker study

Introduction

- Preserved breast tumor tissue samples were obtained from ER positive and PR positive pre-menopausal women
- Women were treated post-surgery with tamoxifen. Clinical data indicated a wide range of mortalities (full responder, partial responder, non responder).
- All samples were from Caucasian women diagnosed with infiltrating duct carcinoma and had been preserved 8 years prior to RNA extraction (FFPE tissues)

Purpose

Retrospective study to demonstrate our High-Specificity miRNA QRT-PCR Detection kit successfully detects miRNA isolated from FFPE samples.





Detection of miRNA Expression in Breast Cancer

- Scott Basehore
- Natalia Novoradovskaya

Sample Preparation

Deparaffinize and digest FFPE tissues

Pulverize frozen tissues in liquid nitrogen

homogenize in Absolutely RNA lysis buffer: Lysate can be stored at -80°C

separate RNA and gDNA with acidified phenoll

Isolate total RNA using Absolutely RNA® kit

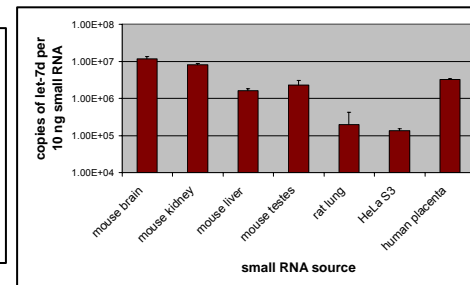
Isolate gDNA from acidified phenol phase

Isolate miRNA from aqueous phase

Nucleic Acid Isolation and QC

Results:

RNA sample, 50 ng	Ct (dRn)					
	B2M	GAPDH	HER2	PGR	Ki67	BIRC5
Frozen breast, normal	18.89	20.01	30.57	30.47	33.83	48.7
FFPE breast cancer 1	20.63	23.41	29.6	28.85	34.86	No Ct
FFPE breast cancer 2	20.22	23.23	31.48	26.02	32.33	36.12
FFPE breast cancer 3	20.43	24.82	30.19	36.13	37.3	No Ct
UHRR	19.44	16.35	32.96	28.01	25.75	26.24

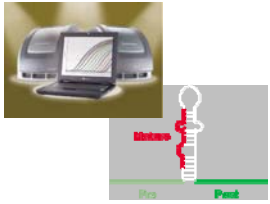




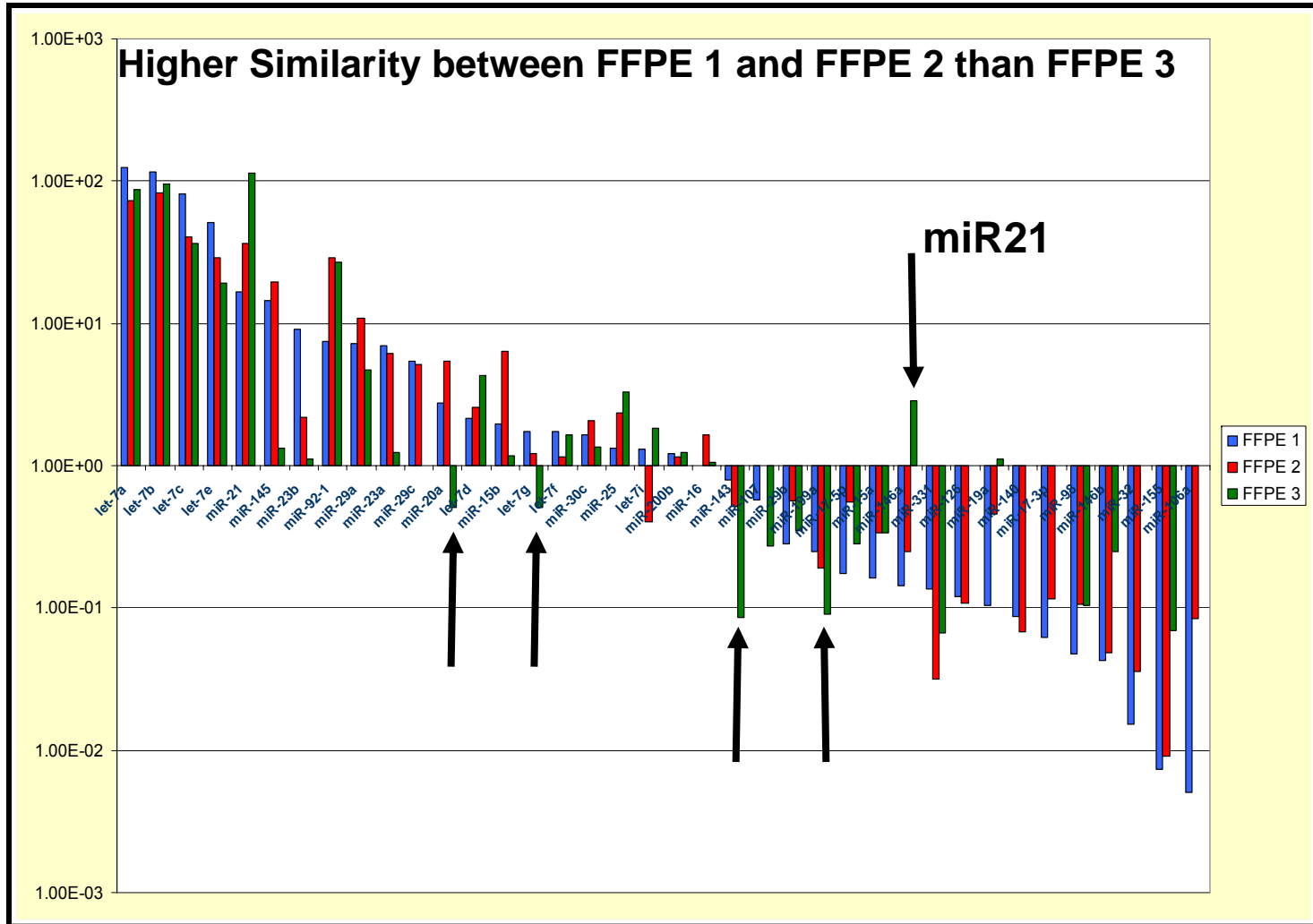
Detection of miRNA Expression in Breast Cancer

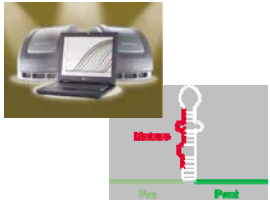
FFPE	Age	Differentiation Grade	TNM	Stage	No. Pos. lymph nodes	No of analyzed lymph nodes	Tumor size
1	47	grade 2	T2N0M0	IIA	0	6	5 cm
2	42	grade 2	T1N0M0	IIA	0	5	1 cm
3	51	grade 2-3	T2N1M0	IIB	2	7	3 cm

FFPE	Location of distant metastasis	Relapse (Y/N) w/in observation time	Disease-free survival (mos)	Death?	Death from main disease
1	no	no	84 mos	no	no
2	no	no	65 mos	no	no
3	lung, bone	yes	54 mos	yes	yes



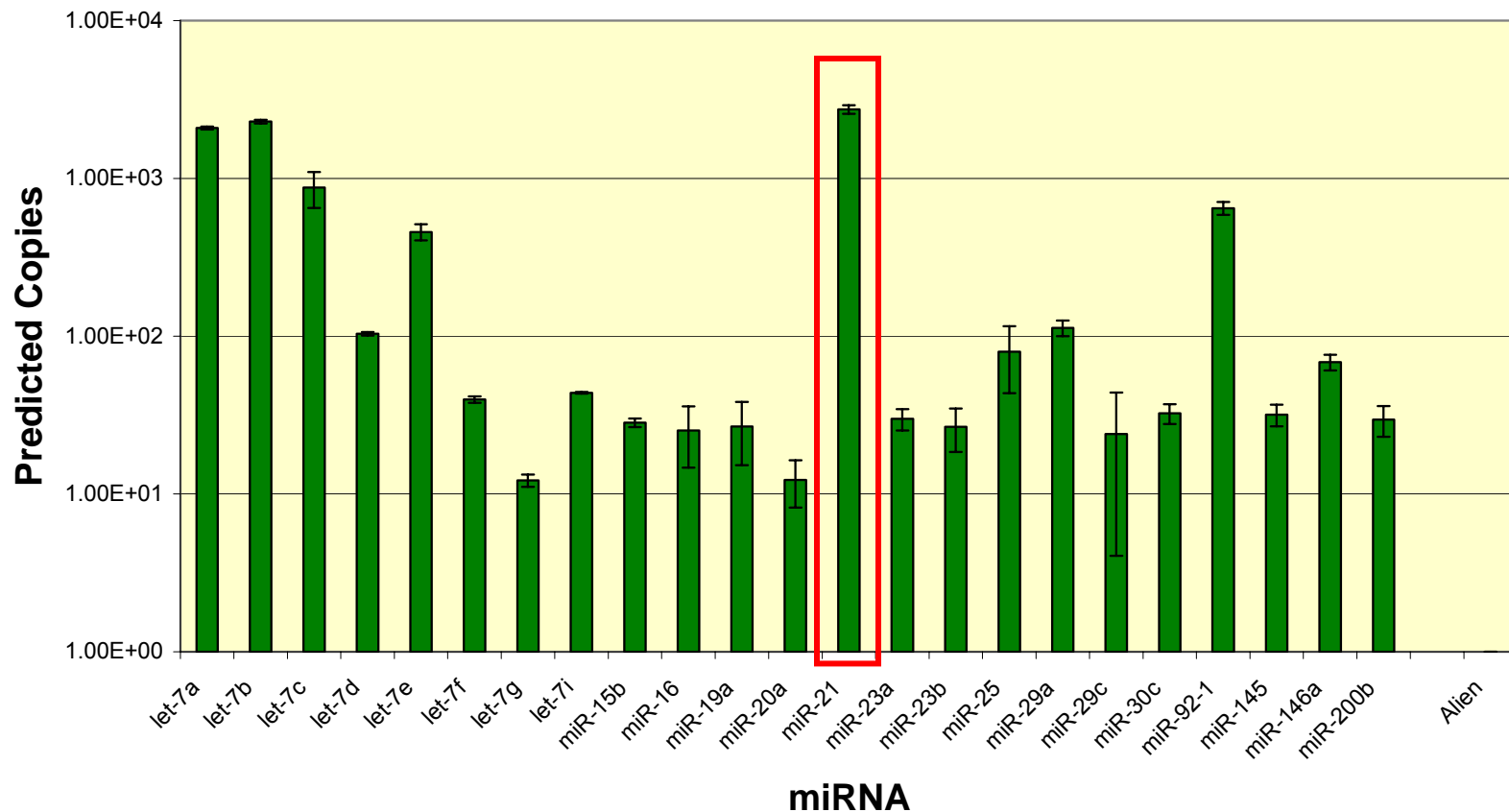
Detection of miRNA Expression in Breast Cancer





Detection of miRNA Expression in Breast Cancer

miRNA Profiling in Breast Cancer FFPE Tissue RNA





Detection of miRNA Expression in Breast Cancer

Conclusion:

- mRNA and miRNA RNA isolated using the Absolutely RNA FFPE kit from FFPE tissues that had been preserved 8 years were of high quality and suitable for QPCR analysis
- Fold differences of miRNA levels during cancer progression from stage I to IIA and to IIB show significant up-regulation of miR-21. This is in agreement with various published results on solid tumors.*

*Volinia, S., et al (2006) Proc Natl Acad Sci 103:2257-2261



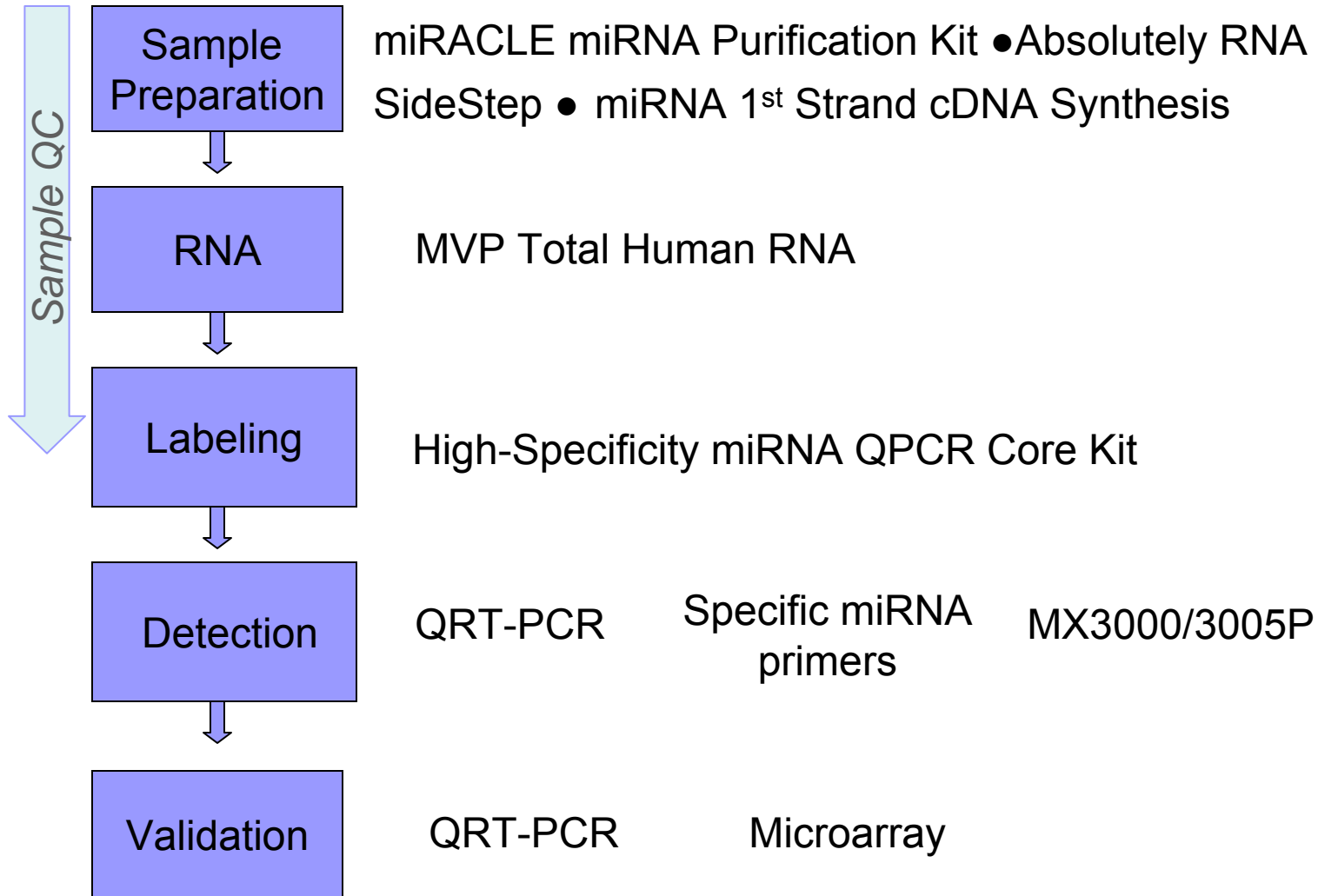


miRNA Forward Primers Relative to Disease

Types of Cancer	
Brain	miR-21
Bladder	miR-21, miR-126, miR-143, miR-145, miR-188, miR-200b, miR-219, miR-331
Breast	miR-17-5p, miR-21, miR-29b-2, miR-126, miR-143, miR-145, miR-146, miR-155, miR-181b-1, miR-188, miR-200b, miR-210, miR-213, miR-219 miR-331
Colon	miR-17-3p, miR-17-5p, miR-20a, miR-21, miR-24-1, miR-24-2, miR-29b-2, miR-30c, miR-32, miR-106a, miR-107, miR-126, miR-128b, miR-130a, miR-143, miR-145, miR-155, miR-188miR-191, miR-195, miR-200b, miR-218-2, miR-219, miR-221, miR-331, miR-223
Lung	miR-17-5p, miR-21, miR-30a, miR-126, miR-128b, miR-143, miR-145, miR-155, miR-188, miR-189, miR-191, miR-199a-1, miR-200b, miR-210,miR-213, miR-219, miR-223, miR-331
Pancreas	miR-17-5p, miR-20a, miR-21, miR-24-1, miR-24-2, miR-25, miR-29b-2, miR-30c, miR-32, miR-92-2, miR-106a, miR-107, miR-126, miR-128b, miR-143, miR-145, miR-146, miR-181b-1, miR-188, miR-191, miR-199a-1, miR-200b, miR-214, miR-218-2, miR-219, miR-221, miR-331
Prostate	miR-17-5p, miR-20a, miR-21, miR-25, miR-29b-2, miR-30c, miR-32, miR-92-2, miR-106a, miR-126, miR-143, miR-145, miR-146, miR-181b-1, miR-188, miR-191, miR-200b, miR-214, miR-218-2, miR-219, miR-223, miR-331
Stomach	miR-21, miR-24-1, miR-24-2, miR-25, miR-92-2, miR-107, miR-191, miR-199a-1, miR-214, miR-218-2, miR-221, miR-223
Thymus	miR-21, miR-126, miR-143, miR-145, miR-188, miR-200b, miR-219, miR-331

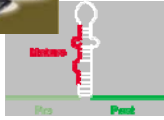


Stratagene's miRNA Workflow Overview





miRNA products from Agilent Technologies



- Microarray Platform:
 - Human miRNA microarray kit (version 1.0)
 - miRNA labeling reagent and hybridization kit
 - March 1: Mouse, Rat and Human (version 2.0) microarray kits
- 2100 Bioanalyzer:
 - Total RNA Assays (for RNA integrity)
 - RNA 6000 Nano Kit
 - RNA 6000 Pico Kit
 - Small RNA Kit (for analysis of small RNAs)
- Stratagene's qPCR:
 - High-Specificity miRNA QRT-PCR Detection Kit
 - miRNA Specific Forward Primers

