



Isolation

Quantification

Functionalization



## The miRNA Revolution

Fall miRNA webinar series

miRNA quantification:  
From experimental design through data analysis

Subu Yerramilli, Ph.D.  
Associate Director R&D  
[subu.yerramilli@qiagen.com](mailto:subu.yerramilli@qiagen.com)

*The second in a series of QIAGEN webinars on miRNA 2011*



# Welcome to the three-part webinar series on The miRNA Revolution



**Webinar 1 :** Meeting the challenges of miRNA research - An introduction to miRNA function & analysis  
**Date:** October 5<sup>th</sup>, 2011  
**Speaker:** Jonathan Shaffer, Ph.D.

**Webinar 2 :** miRNA Quantification  
From experimental design through data analysis  
**Date:** October 12<sup>th</sup>, 2011  
**Speaker:** Subu Yerramilli, Ph.D.

**Webinar 3 :** Profiling miRNA expression in Cells, FFPE, and Serum: on the road to biomarker development  
**Date:** October 19<sup>th</sup>, 2011  
**Speaker:** Eric Lader, Ph.D.



## miRNA quantification:

From experimental design through data analysis

### Topics to be covered

Challenges in miRNA quantification

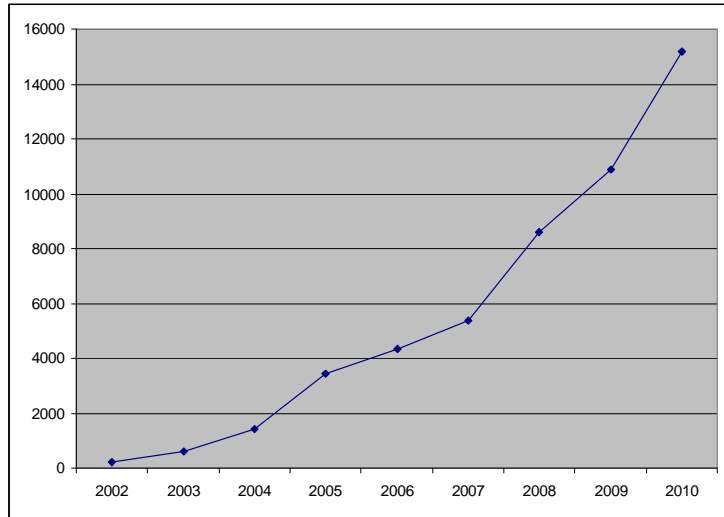
The miScript System

Profiling the miRNome

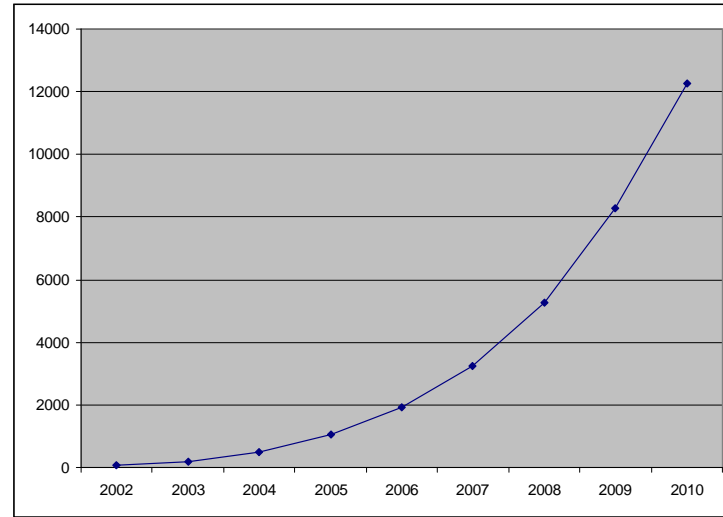
Data Analysis



# Why Quantify miRNA?



Known miRNAs in Sanger DB



miRNA publications in Medline

Virtually every publication included quantification experiments

Changes in miRNA correlated with gene expression changes in development, differentiation, signal transduction, infection, aging, and disease



# miRNA Quantification Challenges

- Short sequences: ~16-26 nt
- Presence of miRNA isoforms
  - Near identical with ~1-3 nt differences
  - Can be anywhere along the sequence
- miRNAs with less than ideal GC% (very low or very high)
- Presence of repetitive sequences
  - hsa-miR-296-5p: `AGGGCCCCCCUCAAUCCUGU`
  - hsa-miR-638: `AGGGAUCGCGGGCGGGUGGCGGCCU`
  - hsa-miR-940: `AAGGCAGGGCCCCCGCUCCCC`
  - mmu-miR-703: `AAAACCUUCAGAAGGAAAGAA`
  - mmu-miR-466b-5p: `UAUGUGUGUGUACAUGUACAUA`
  - mmu-miR-574-5p: `UGAGUGUGUGUGUGUGAGUGUGU`
- 3'-and 5'-end variations (Next slide)




# 3'- End & 5'-End Variations in miRNA

Home Search Browse Help Download Blog Submit

## Deep sequencing reads for stem-loop sequence MI0000060

Stem-loop ID [hsa-let-7a-1](#)

Reads	<a href="#">hsa-let-7a</a>	<a href="#">hsa-let-7a*</a>	Count
	..... <a href="#">AUGAGGUAGUAGGUUGUAUAGUU</a> .....		39
	..... <a href="#">AUGAGGUAGUAGGUUGUAUAGUUU</a> .....		2
↗	..... <a href="#">UGAGGUAGUAGGUUGUAUAGU</a> .....		76197
	..... <a href="#">UGAGGUAGUAGGUUGUAUA</a> .....		1740
↗	..... <a href="#">UGAGGUAGUAGGUUGUAUAGUUUU</a> .....		795
	..... <a href="#">UGAGGUAGUAGGUUGUA</a> .....		615
	..... <a href="#">UGAGGUAGUAGGUUG</a> .....		46
	..... <a href="#">UGAGGUAGUAGGUUGUAUAGUUUA</a> .....		19
	..... <a href="#">UGAGGUAGUAGGUUGUAUAGUUUUAGG</a> .....		8
	..... <a href="#">UGAGGUAGUAGGUUGUAUAGUUUUAG</a> .....		7
↗	..... <a href="#">GAGGUAGUAGGUUGUAUAG</a> .....		12
	..... <a href="#">GAGGUAGUAGGUUGUAUAGUUUU</a> .....		6
	..... <a href="#">AGGUAGUAGGUUGUAUAGUUUU</a> .....		4
↗	..... <a href="#">GGUAGUAGGUUGUAUAGUU</a> .....		6
↗	..... <a href="#">GUAGUAGGUUGUAUAGUU</a> .....		5
	.....	..... <a href="#">CUAUACAAUCUACUGUCUUU</a> .....	16
	.....	..... <a href="#">UAUACAAUCUACUGUCUUUCCU</a> .....	16
	.....	..... <a href="#">UAUACAAUCUACUGUCUUUCC</a> .....	3
	.....	..... <a href="#">AUACAAUCUACUGUCUUUCCU</a> .....	6
	UGGGAUGAGGUAGUAGGUUGUAUAGUUUUAGGGUCACACCCACCACUGGGAGAUAA	CUAUACAAUCUACUGUCUUUCUA	



# miRNA Quantification Strategies



- Micro array based approaches
  - Low specificity, needs further validation
- Expression analysis by deep sequencing
  - Expensive, and low throughput
  - Not practical for large number of samples
- Real time PCR based approaches
  - Fast, flexible, scalable, accurate, sensitive



# Real-Time PCR Based Approach miRNA 3'-End Dependent Reverse Transcription

e.g. hsa-Let-7a

5' -UGAGGUAGUAGGUUGUAUAGUU- 3'



→ miRNA specific cDNA

```
hsa-let-7a
..AUGAGGUAGUAGGUUGUAUAGUU.....
..AUGAGGUAGUAGGUUGUAUAGUU.....
...UGAGGUAGUAGGUUGUAUAGU.....
...UGAGGUAGUAGGUUGUAUA.....
...UGAGGUAGUAGGUUGUAUAGUUUU.....
...UGAGGUAGUAGGUUGUA.....
...UGAGGUAGUAGGUUG.....
...UGAGGUAGUAGGUUGUAUAGUUUUUA.....
...UGAGGUAGUAGGUUGUAUAGUUUUAGG.....
...UGAGGUAGUAGGUUGUAUAGUUUUAG.....
...GAGGUAGUAGGUUGUAUAG.....
...GAGGUAGUAGGUUGUAUAGUUUU.....
...AGGUAGUAGGUUGUAUAGUUUU.....
...GGUAGUAGGUUGUAUAGUU.....
...GUAGUAGGUUGUAUAGUU.....
```

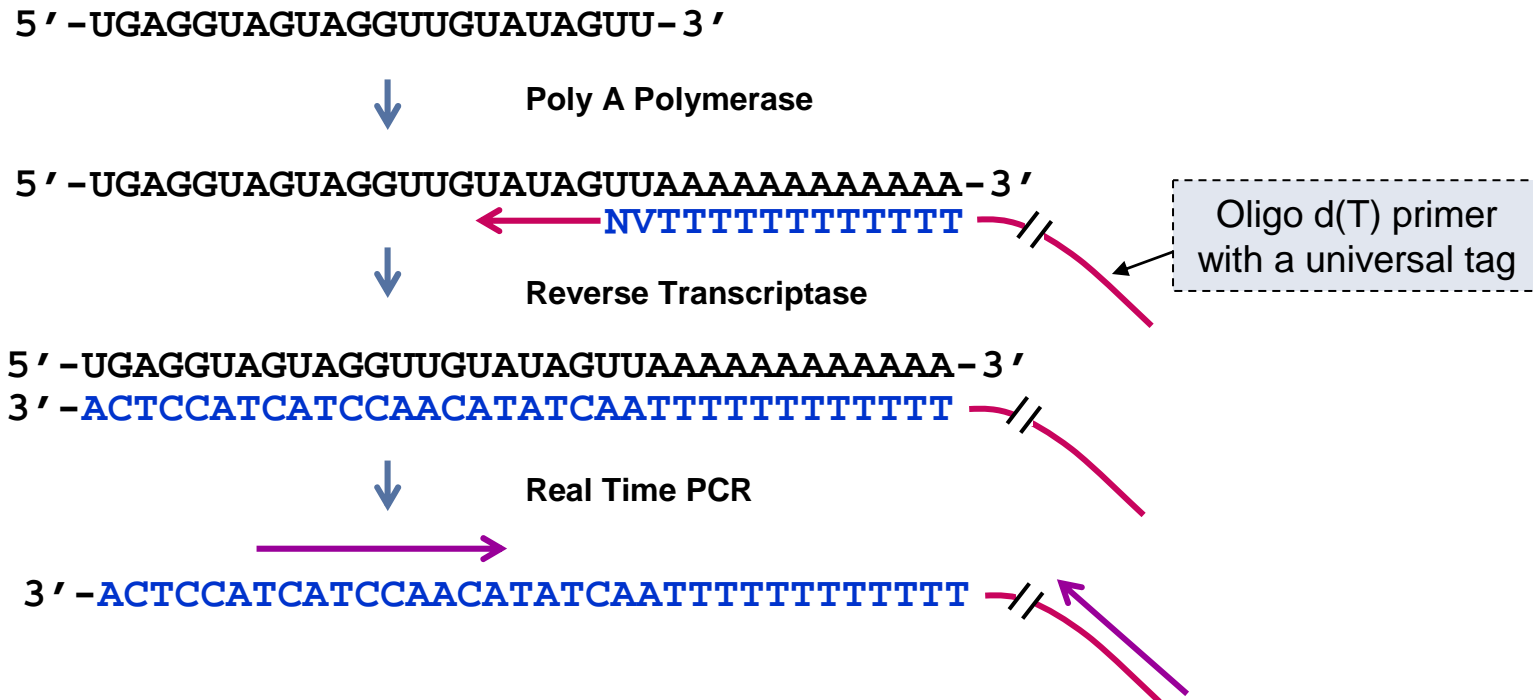
- Each miRNA detection requires a separate cDNA synthesis reaction
- A given RT primer could prime multiple miRNAs if these share the 3'-sequences
- Doesn't address 3'-end Polymorphism

[www.mirbase.org](http://www.mirbase.org)



# Real-Time PCR-Based Approaches

## Universal Reverse Transcription



qPCR with  
miRNA specific FWD primer  
Universal Reverse Primer

- A single cDNA prepared per sample is sufficient to interrogate many miRNAs
- 3'-end polymorphism tolerated



## 1. miScript II RT kit

- miScript Reverse Transcriptase Mix
- 5x HiSpec Buffer
- 5x HiFlex Buffer
- 10x Nucleics Mix
- Water

## 2. miScript SYBR Green PCR Kit

- QuantiTect SYBR Green PCR MM
- Universal Primer

## 3. Assays

- miScript Primer Assays
- miScript Precursor Assays
- QuantiTect Primer Assays
- miScript Controls (NEW)

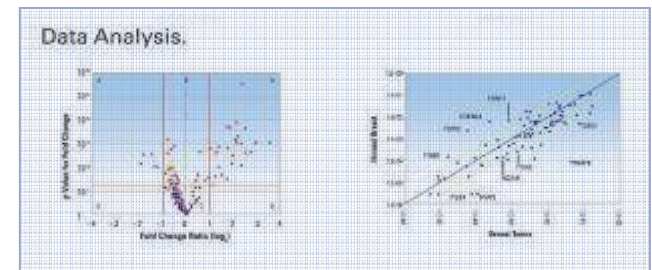
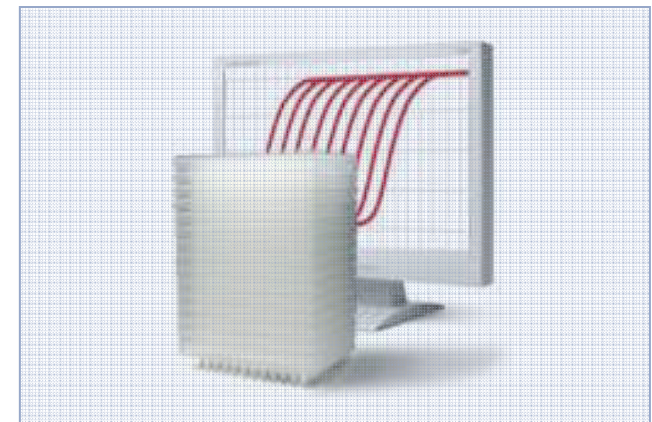
## 4. miScript PCR Arrays

- Whole miRNome PCR Arrays
- miScript miRNA PCR (Focused) Arrays



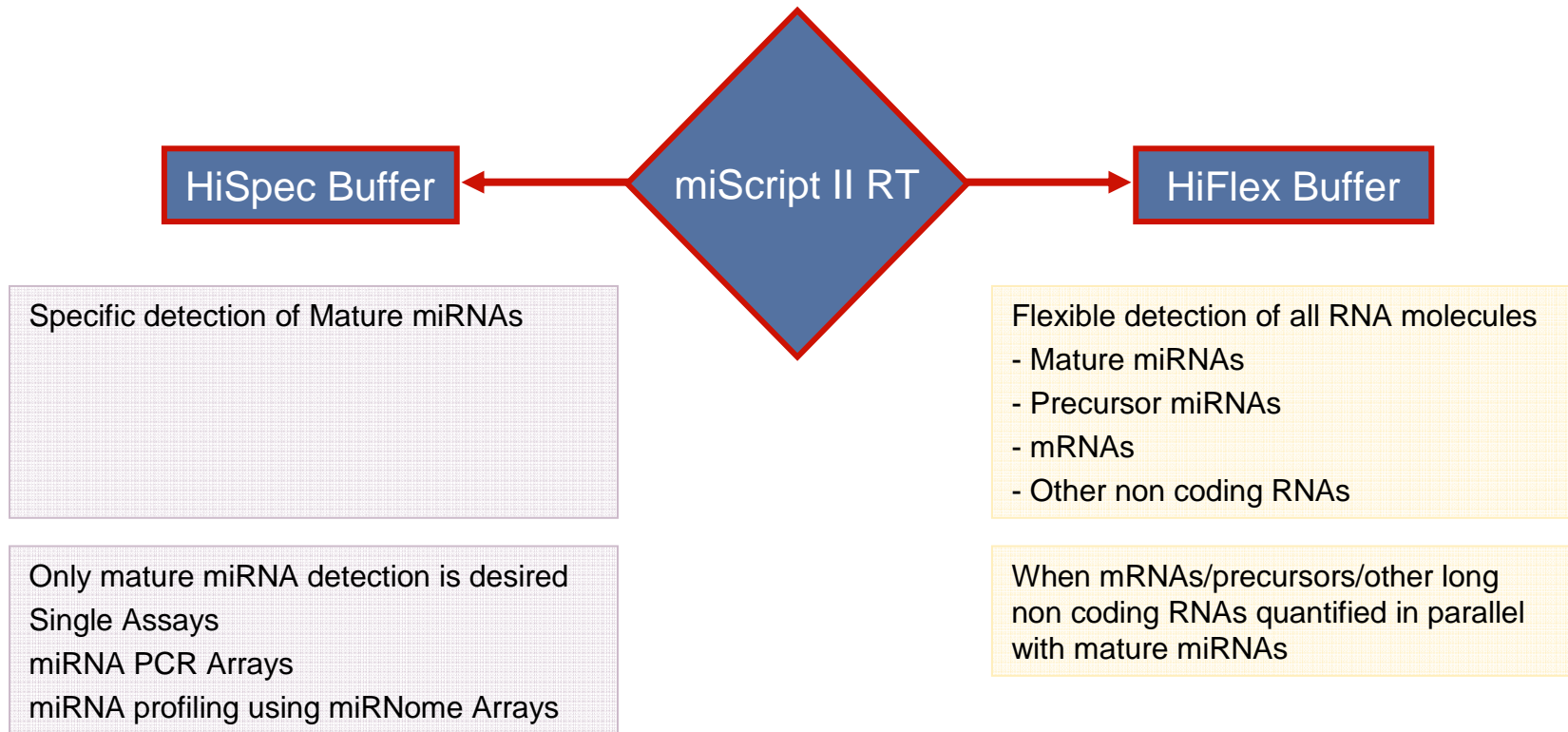
## 5. Online Array Data analysis software

- ACCESS AT NO COST**
- **Data analysis made easy**





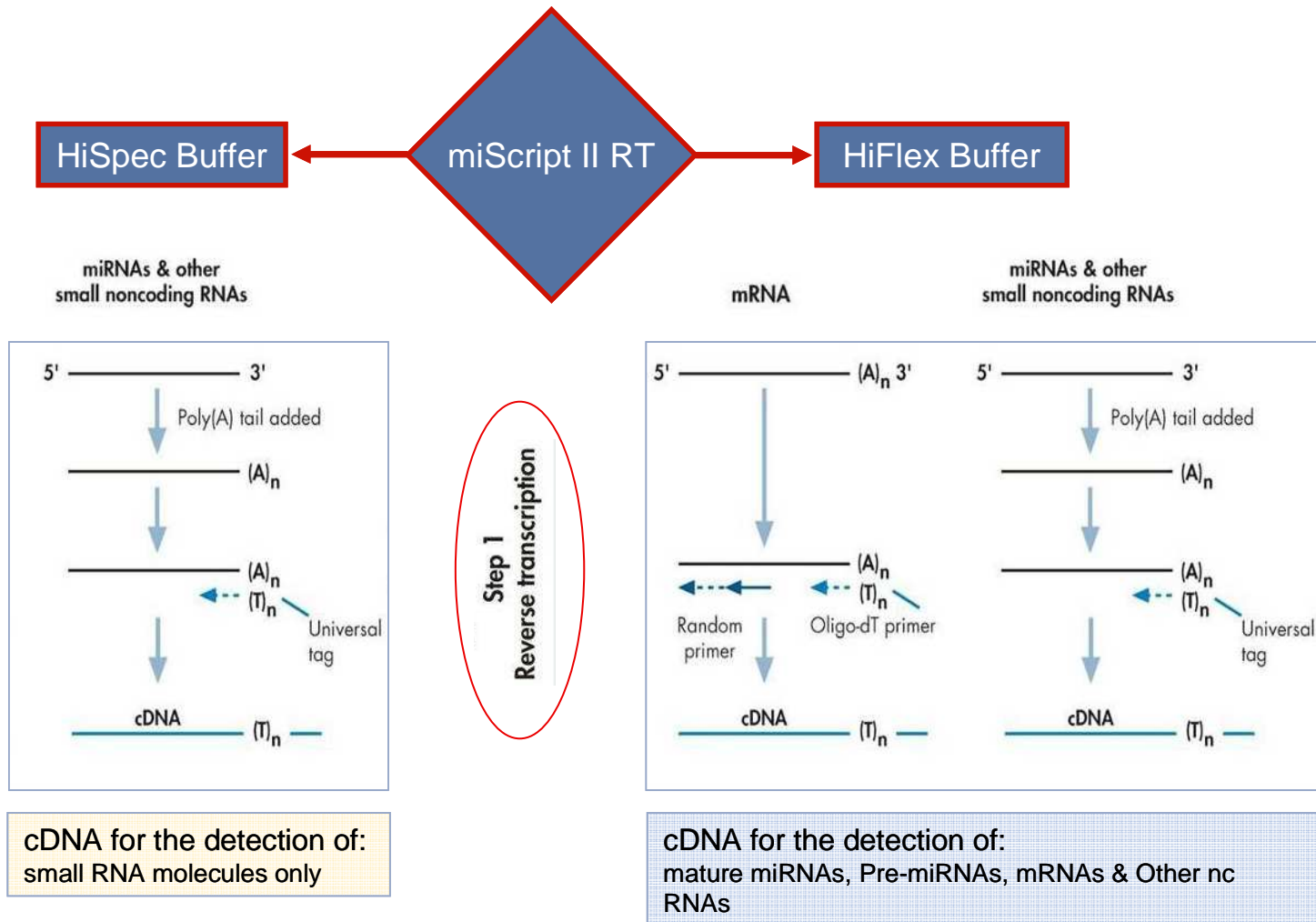
# miScript II RT: Dual Buffer System Offers Flexibility & Specificity



**New** miScript Controls work equally well in both buffer systems

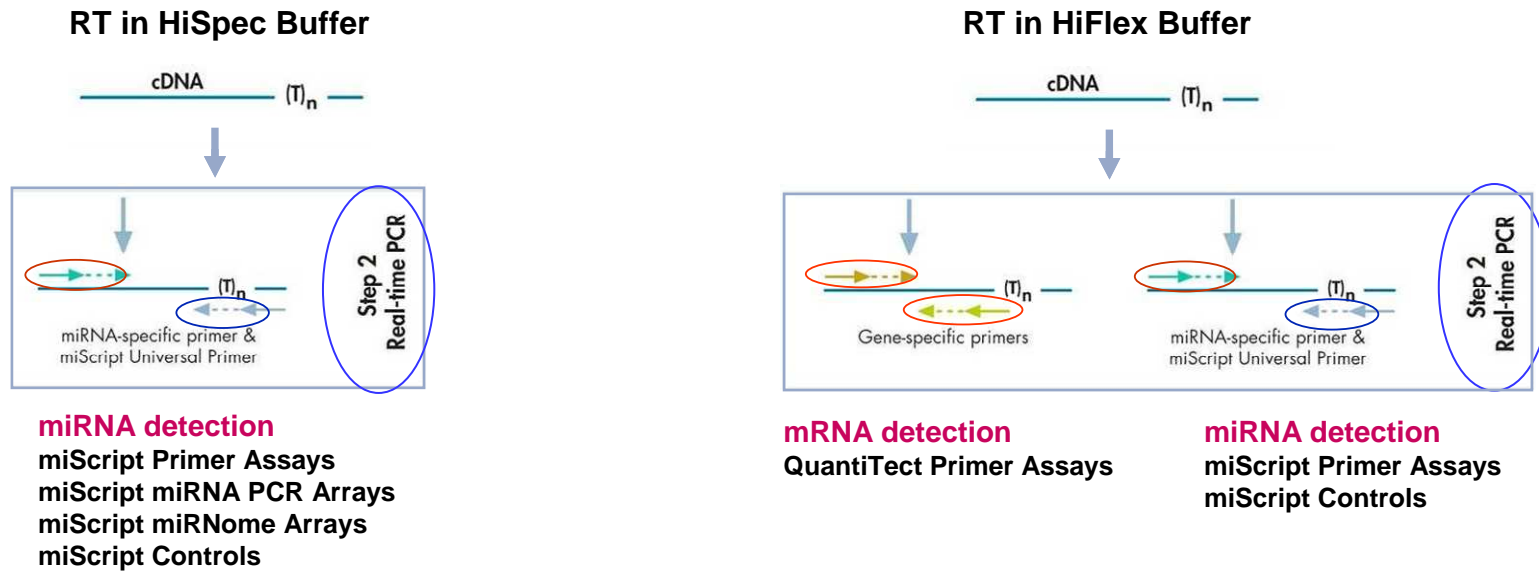


# Unique RT Buffer Formulation: Offers Flexibility, Sensitivity & Specificity

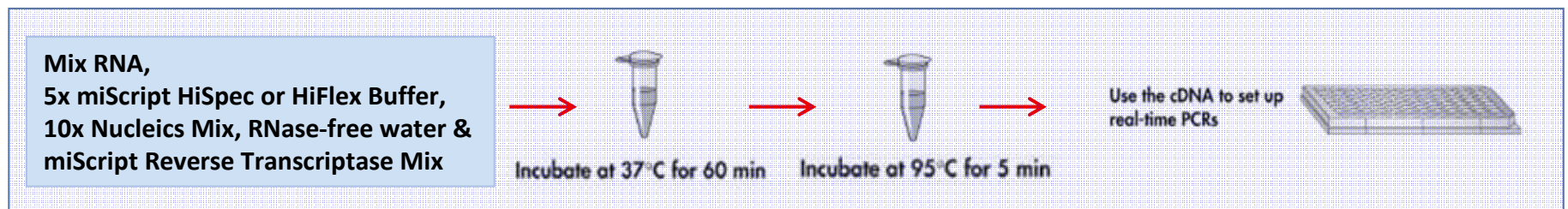




# miScript PCR System: Real time PCR Step



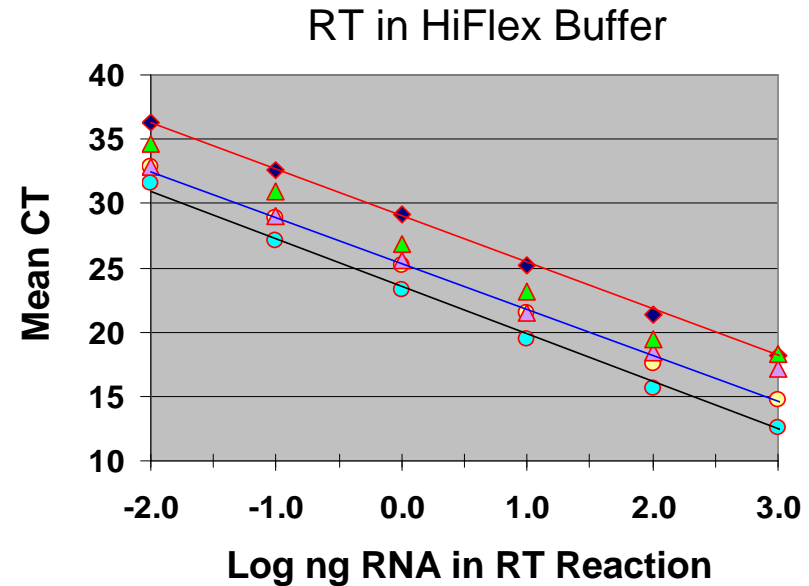
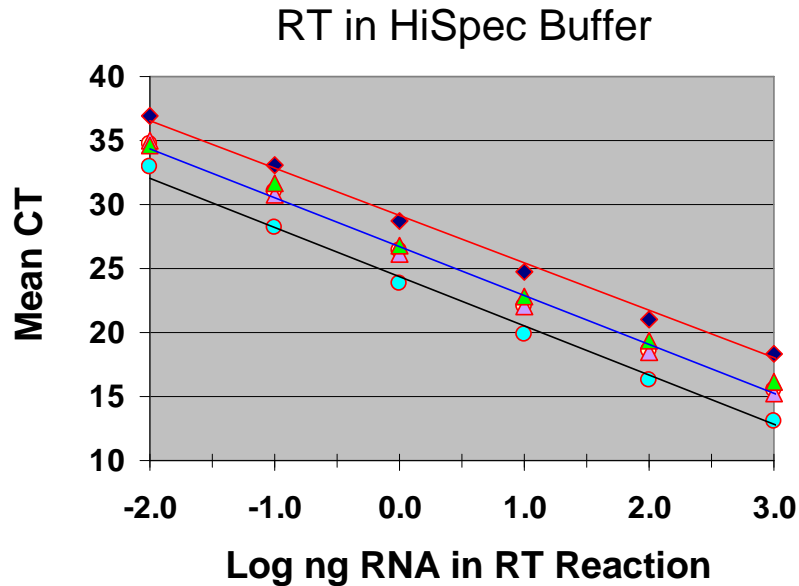
## miScript PCR System workflow: Simple & Easy steps





# Linear cDNA Synthesis Over Five Logs Of Input RNA: With miScript II RT

Linear in the range of 10 pg to 1 µg



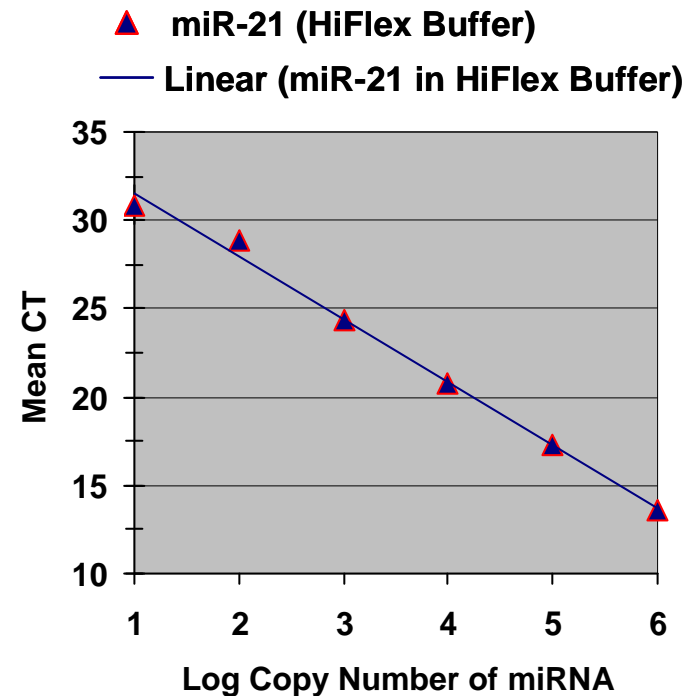
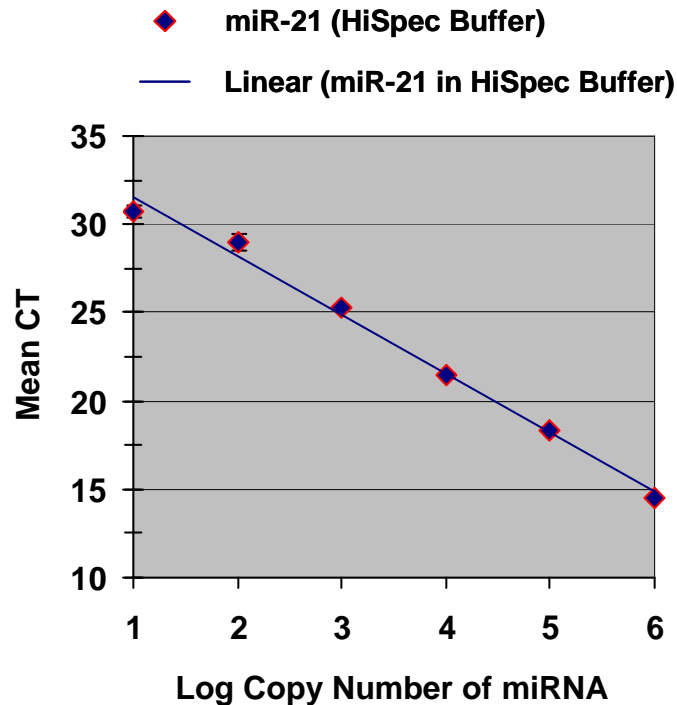
RNA input range (Recommended)

- 5x HiSpec Buffer → 10 pg to 2 µg
- 5x HiFlex Buffer
  - For pre-miRNA detections → max of 500 ng
  - For all the other RNAs → 10 pg to 1 µg

- ◆ miR-100
- miR-16
- miR-21
- △ miR-27a
- ▲ miR-20a
- Linear (miR-100)
- Linear (miR-16)
- Linear (miR-21)



# High Detection Sensitivity with miScript PCR System: From 10 to 10<sup>6</sup> Copies.



cDNA prepared either in HiSpec Buffer or in HiFlex Buffer using synthetic miR-21 in the presence of a carrier RNA was used as the template.

Real time PCR performed using log dilutions of the cDNA template representing various copy numbers.



## When & Why To Use HiSpec Buffer?

- When to use HiSpec Buffer for RT reaction ?
  - cDNA is used exclusively for mature miRNA detection
    - miRNA profiling using miScript miRNome Arrays
    - miScript miRNA PCR Arrays
    - Individual miScript Primer Assays for mature miRNA detection
  
- Why Use HiSpec Buffer ?
  - Increases sensitivity
  - Eliminates the need to verify the dissociation curves
    - A tedious step in miRNA profiling
  - Higher input amount of RNA tolerated (up to 2 µg of total RNA/RT reaction)
  - Only mature miRNAs & miScript controls are converted into cDNA
  - Any other longer RNA species are not converted into cDNA
    - Hence reduces background noise
    - Higher input amount of RNA tolerated (up to 2 µg of total RNA)
  - All miScript Primer Assays & Arrays are bench validated with cDNA made using HiSpec Buffer



## Why To Use HiFlex Buffer?

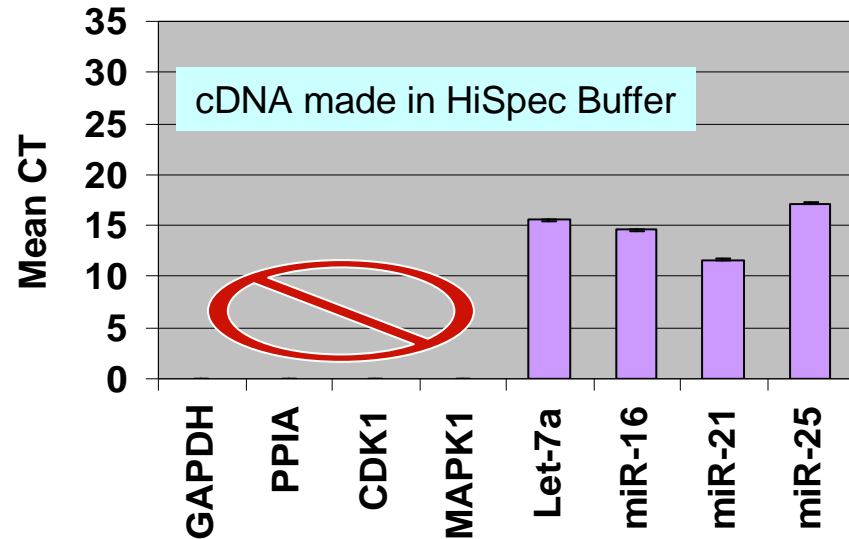
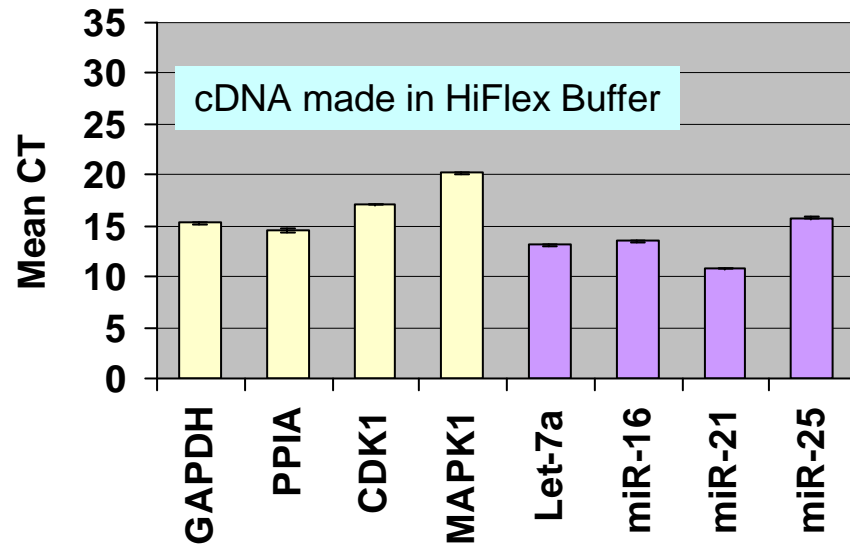
- Why use HiFlex Buffer in RT reaction ?
  - Flexible detection of all RNA molecules is desired
  - Same cDNA can be used for mature miRNAs, Pre-miRNAs, mRNAs & other non coding RNAs
    - Simultaneous detection of mature miRNAs along with its precursors or
    - Simultaneous detection of mature miRNAs along with their mRNA targets
  - Most helpful when starting RNA is a limiting factor when multiple RNA species are to be quantified



# HiFlex Buffer: Detection Of microRNAs & mRNAs From The Same cDNA

*HiFlex Buffer* should be used to prepare cDNA for the quantification of both mature microRNAs & mRNAs using appropriate Primer Assays.

Long RNAs such as mRNAs are not converted into cDNA during RT reaction in *HiSpec Buffer*



Primer Assays used

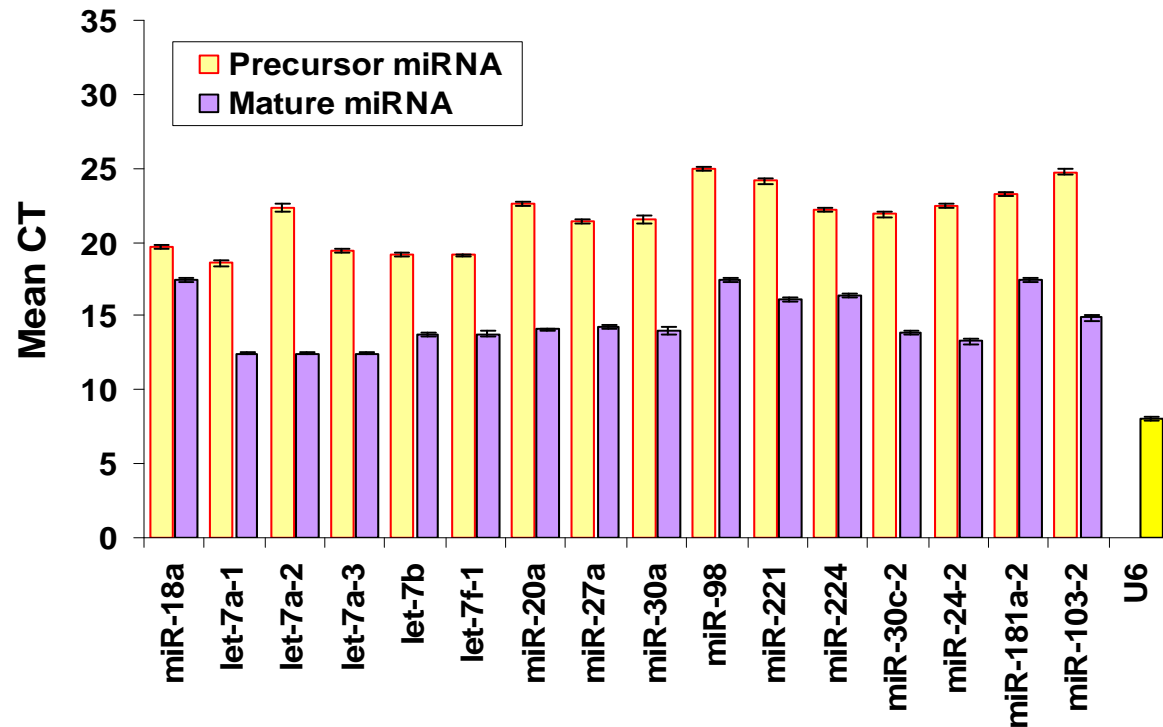
- mRNA quantification
- microRNA quantification

- QuantiTect Primer Assays
- miScript Primer Assays



# HiFlex Buffer Allows Parallel Detection of Precursor & Mature miRNAs From The Same cDNA

HiFlex Buffer should be used to prepare cDNA for the quantification of both mature miRNAs and their precursors using appropriate Primer Assays.



Relative levels of various precursor & mature miRNAs in HeLa S3 cells



## miScript miRNome PCR Array Content: miRBase v16

Species	Number of Assays (miRBase v16)	Total Number of Plates	
		96 well format	384 well format
Human	<b>1068</b>	<b>12</b>	<b>3</b>
Mouse	<b>947</b>	<b>12</b>	<b>3</b>
Rat	<b>653</b>	<b>8</b>	<b>2</b>
Dog	<b>277</b>	<b>4</b>	<b>1</b>

### miScript miRNA Assays & Arrays:

- Each Assay has been bench validated using synthetic miRNA pool (v16)  
miRNA specific FWD Primers only & Universal REV Primer is a part of the miScript SYBR Green PCR Kit



# miScript Primer Assays: Highly Specific. Detection of Let-7 Family of Isoforms With High Specificity

CLUSTAL 2.1 multiple sequence alignment

```

hsa-let-7b  UGAGGUAGUAGGUUGUGUGUU 22
hsa-let-7c  -----A----- 22
hsa-miR-98  -----A---A-U--- 22
hsa-let-7d  A-----CA-A--- 22
hsa-let-7e  -----G-----A-A--- 22
hsa-let-7a  -----A-A--- 22
hsa-let-7f  -----A---A-A--- 22
hsa-let-7g  -----U---ACA--- 22
hsa-let-7i  -----U---CU--- 22
  
```

\*\*\*\*\* \* \*\*\*

% Activity relative to perfect match primer assay as 100%\*

cDNA Used in qPCR	Primer Assay Used								
	Let-7b	Let-7c	miR-98	Let-7d	Let-7e	Let-7a	Let-7f	Let-7g	Let-7i
<b>Let-7b</b>	100.0	1.8	0.0	0.0	0.0	0.0	0.0	0.0	0.0
<b>Let-7c</b>	0.5	100.0	0.0	0.0	2.4	0.1	0.0	0.0	0.0
<b>miR-98</b>	0.0	0.2	100.0	0.1	0.0	0.1	0.0	0.0	0.1
<b>Let-7d</b>	0.1	0.0	0.0	100.0	0.0	0.4	0.0	0.0	0.0
<b>Let-7e</b>	0.1	0.0	0.0	0.0	100.0	0.2	0.0	0.0	0.0
<b>Let-7a</b>	0.1	0.6	0.0	0.5	3.1	100.0	0.1	0.0	0.0
<b>Let-7f</b>	0.6	0.1	0.0	0.1	0.0	1.0	100.0	0.1	0.1
<b>Let-7g</b>	0.6	0.2	0.0	0.1	0.0	0.0	0.0	100.0	0.2
<b>Let-7i</b>	0.1	0.0	0.0	0.0	0.0	0.0	0.0	0.1	100.0

cDNA prepared using HiFlex Buffer

\*Similar results obtained with cDNA made using HiSpec Buffer



# New miScript Controls

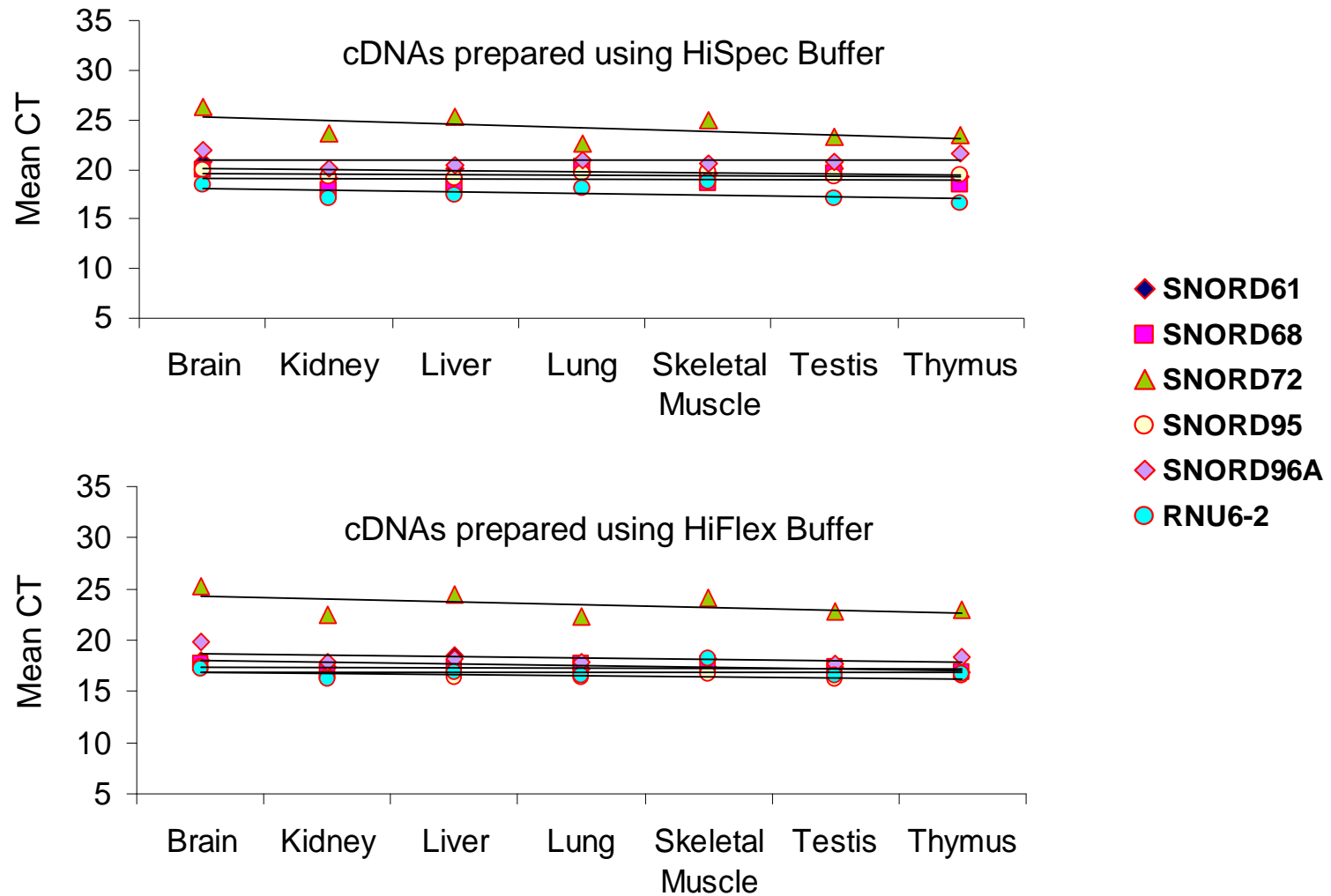
- Highly conserved across multiple species
  - Human, Mouse, Rat, dog (and possibly Rhesus, and pig)
  - The same controls work in all these species
- Relatively stable expression in many tissues
- Amplification efficiencies of these assays are 100%
- ➔ ■ Consistent performance in both HiSpec and HiFlex Buffers
- Ideal normalizers for  $\Delta\Delta\text{CT}$  method of data analysis

Official Symbol	Gene ID	Accession	Alias	PCR Product Size (bp)
SNORD61	26787	NR_002735	U61; RNU61; HBII-342	83
SNORD68	606500	NR_002450	HBII-202	84
SNORD72	619564	NR_002583	HBII-240	89
SNORD95	619570	NR_002591	U95	81
SNORD96A	619571	NR_002592	U96A	84
RNU6-2	26826	NR_002752	U6; RNU6B	85

MIQE Guidelines recommend the use of multiple Normalization controls

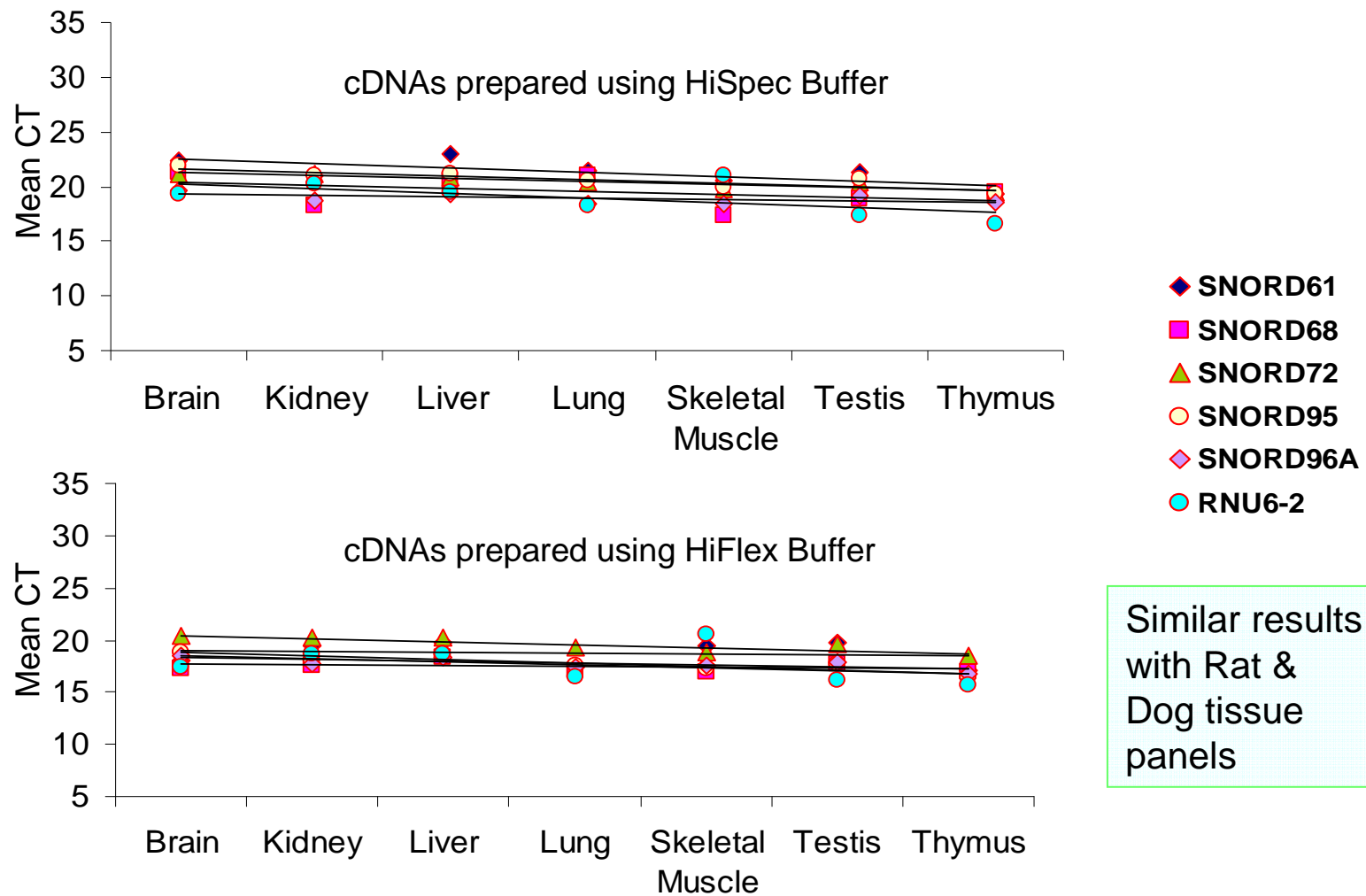


# miScript Controls: Constant Level Of Expression Across A Panel of Human Tissues





# miScript Controls: Constant Level Of Expression Across A Panel of Mouse Tissues





# miScript Controls: Amplification Efficiencies Close To 100%



Similar results with Rat & Dog Universal cDNA

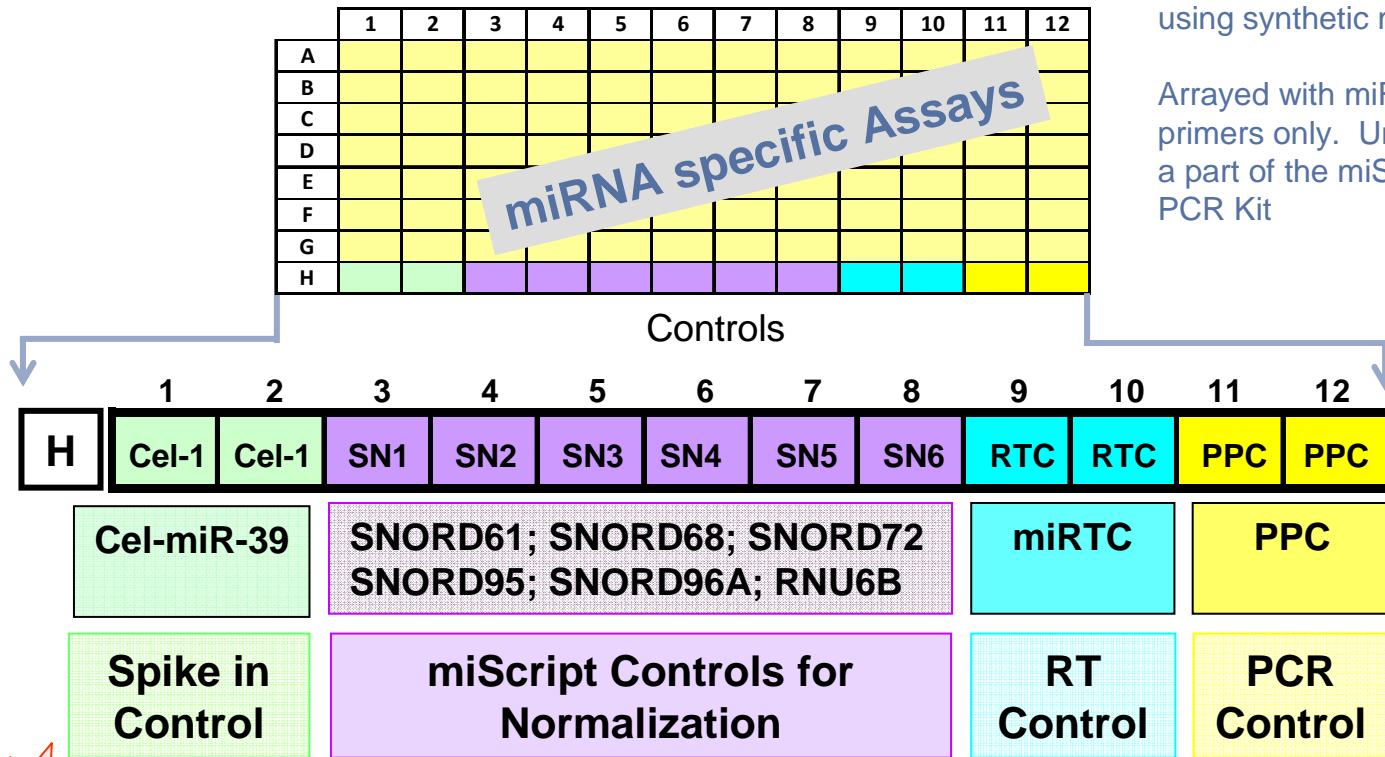


# miScript miRNA array format: Standardized controls (96 well plate)

## miScript miRNA Assays & Arrays

Each assay has been bench validated using synthetic miRNA pool (v16)

Arrayed with miRNA specific FWD primers only. Universal REV primer is a part of the miScript SYBR Green PCR Kit



miScript Controls: Common for Human, Mouse, Rat and Dog Arrays



# miScript miRNA array format: Standardized controls (4X96 well plate)

	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	
A	1	2																							
B	3	4																							
C																									
D																									
E																									
F																									
G																									
H																									
I																									
J																									
K																									
L																									
M																									
N																									
O	1	2	1	2	1	2	1	2	1	2	1	2	1	2	1	2	1	2	1	2	1	2	1	2	
P	3	4	3	4	3	4	3	4	3	4	3	4	3	4	3	4	3	4	3	4	3	4	3	4	

Arrayed with miRNA specific FWD primer only

Universal REV primer is a part of the miScript SYBR Green PCR MM

O	1	2	1	2	1	2	1	2	1	2	1	2	1	2	1	2	1	2	1	2	1	2	1	2
P	3	4	3	4	3	4	3	4	3	4	3	4	3	4	3	4	3	4	3	4	3	4	3	4

Cel-miR-39

SNORD61; SNORD68; SNORD72  
SNORD95; SNORD96A; RNU6B

miRTC

PPC

Spike in Control

miScript Controls for Normalization

RT Control

PCR Control



miScript Controls: Common for Human, Mouse, Rat and Dog Arrays

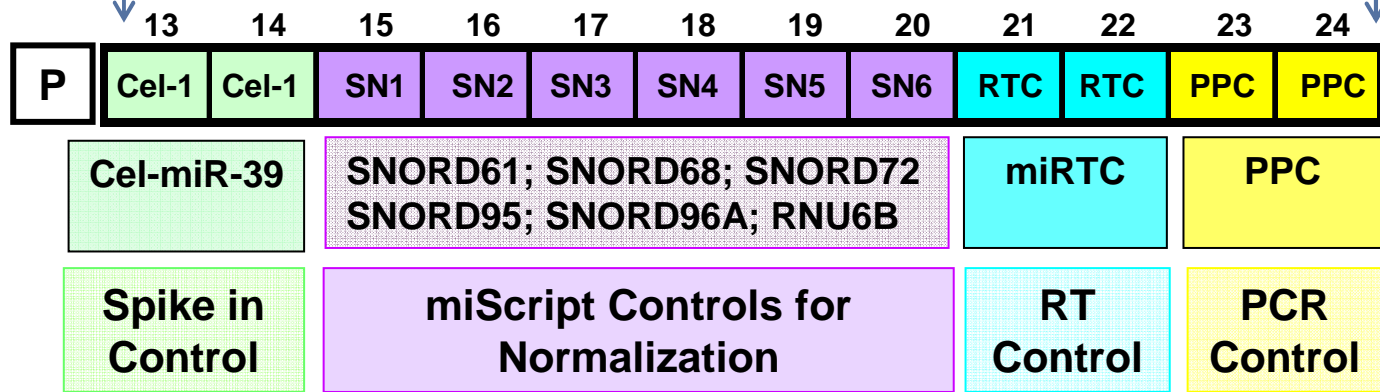


# miScript miRNA array format: Standardized controls (384 well plate)

	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24
A																								
B																								
C																								
D																								
E																								
F																								
G																								
H																								
I																								
J																								
K																								
L																								
M																								
N																								
O																								
P																								

**miRNA specific Assays**

Arrayed with miRNA specific FWD primer only  
Universal REV primer is a part of the miScript SYBR Green PCR MM



miScript Controls: Common for Human, Mouse, Rat and Dog Arrays



# miScript miRNA PCR Array portfolio

## miRNome, Disease, & Pathway-Focused Analysis



■ miRNome miScript miRNA PCR Arrays (v16)	✓	✓	✓	✓
■ Pathway Focused miScript miRNA PCR Arrays:				
<input type="checkbox"/> miFinder Array	✓	✓	✓	✓
<input type="checkbox"/> Serum & Plasma	✓	✓	✓	
<input type="checkbox"/> Cancer PathwayFinder Array	✓	✓	✓	
<input type="checkbox"/> Brain Cancer Array	✓	✓	✓	
<input type="checkbox"/> Breast Cancer Array	✓	✓	✓	
<input type="checkbox"/> Ovarian Cancer Array	✓	✓	✓	
<input type="checkbox"/> Cell Development & Differentiation Array	✓	✓	✓	
<input type="checkbox"/> Immunopathology	✓	✓	✓	
<input type="checkbox"/> Inflammatory Response & Autoimmunity Array	✓	✓	✓	
<input type="checkbox"/> Neurological Development & Disease	✓	✓	✓	



## QIAGEN Solutions for miRNA Research



### Characterizing the miRNome Using the miScript System



# miRNA quantification: Experimental Design & Key Considerations

## Scientific question

- A well defined biological question

## Experimental sample set

- Sample set containing experimental samples along with appropriate controls
- Statistically meaningful number of replicates (biological replicates)
  - A minimum of 3 replicates recommended
  - Replicates enhance the statistical power
- Inclusion of proper normalization controls & spike in controls in the analysis

## Hallmarks of a good experimental system

- Simple & easy to carry out
- High sensitivity & specificity
- Availability of companion research tools

## Data Analysis & statistical considerations

- Easy and simple data analysis tools → online data analysis software

## Confirmation of the results

- Customizable solutions & single assays



## How much RNA is Needed for miRNA Quantification? Depends on the Detection Limit of the System and miRNA Abundance

- System sensitivity:
  - miScript RT-PCR can detect 10 copies per PCR
- RNA abundance:
  - A 'typical' cell contains ~15-30 pg of total RNA
  - miRNA expression levels vary from as low as 10 copies/cell to as high as 35,000 copies/cell
- How much template is needed to detect an abundant miRNA?
  - An abundant miRNA should be detectable with 10-15 pg of template



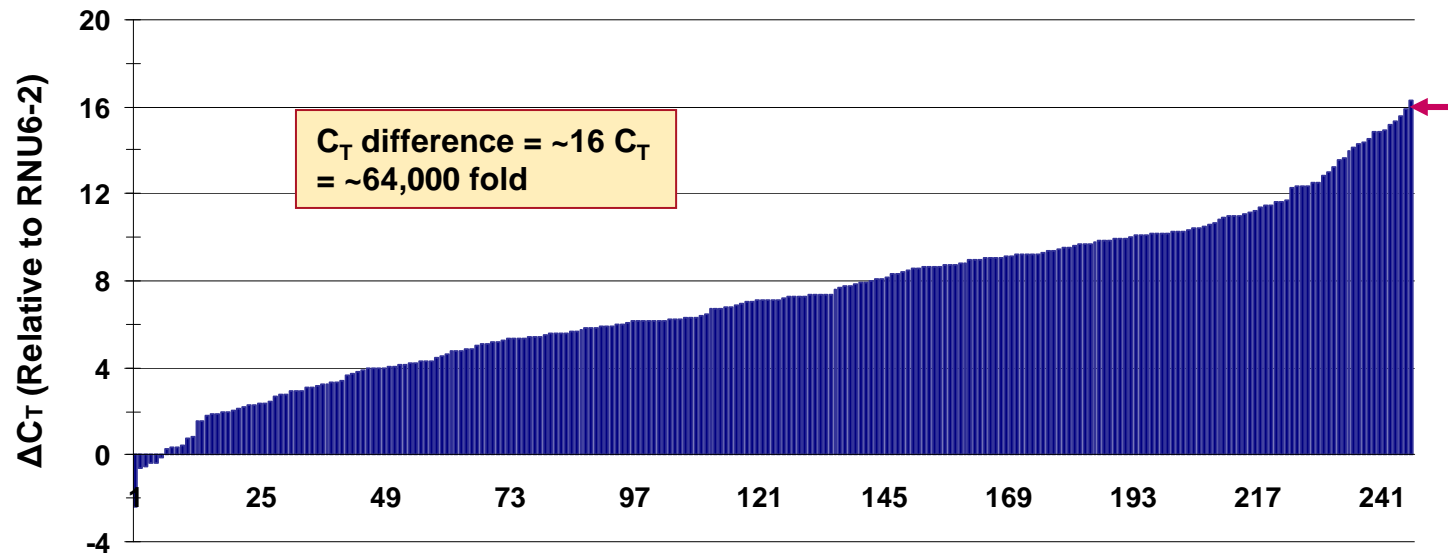
# How much RNA is Needed for miRNA Quantification?

Depends on the Detection Limit of the System and miRNA Abundance

- What about miRNAs expressed at ~10-100 copies/cell?
  - Considering that 10x more template result in a 3.3 CT shift
  - Use 10-20 times more template (0.1 ng-0.2 ng) to detect rare miRNAs
- To be able to 'call with certainty' a miRNA is 'not expressed'
  - use >20x more RNA (0.5 to 1.5 ng)
- As a guideline, consider using ~ 50 pg to 500 pg template/PCR
  - Up to 20,000 PCR reactions per 1  $\mu$ g miScript RT reaction
  - Use an RT system that can work in a broad input RNA range.
    - miScript allows 10 pg - 1 $\mu$ g in put RNA into RT reaction



# miRNA Expression In HeLa Cells

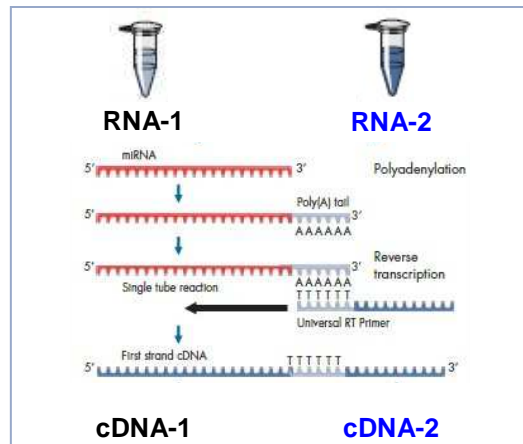


Deep sequencing results show that miRNA reads range from 20 to 292,000 (in *HeLa* cells)

*Shin, C., et. al. 2010; Mol. Cell, 38, 789*

# miScript miRNA PCR Array workflow: Genome-wide, disease, & pathway-focused analysis

## 1. Convert miRNA into cDNA



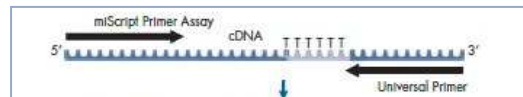
- cDNA Synthesis in HiSpec Buffer
  - 1 hour

## 2. Assemble Real time PCR



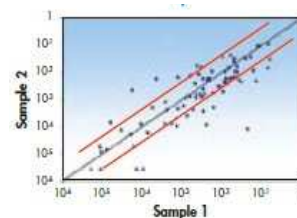
- Load Plates
  - 2 minutes

## 3. Run in real-time PCR cycler

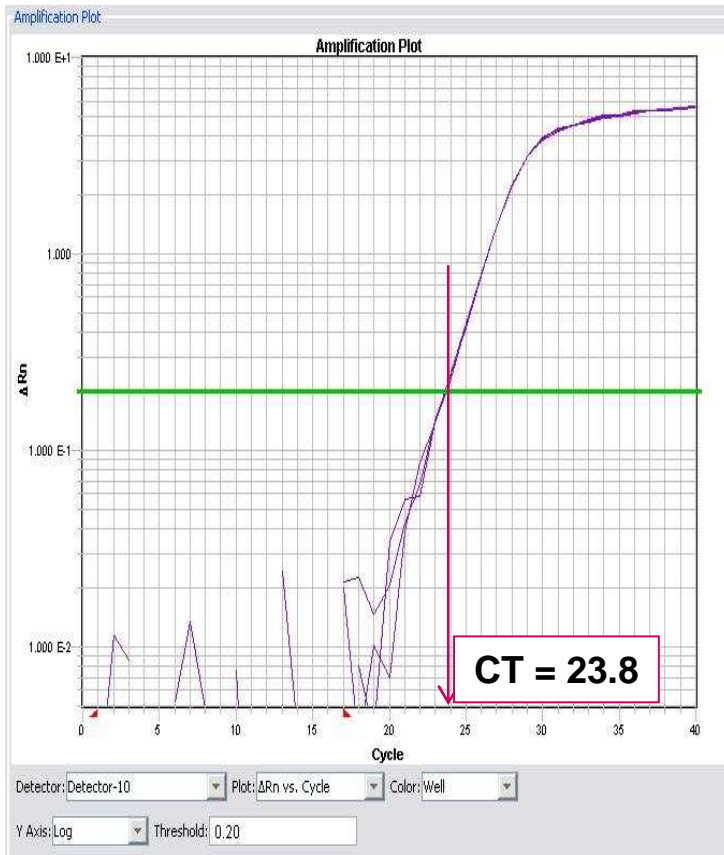


- Run 40 cycle qPCR Program
  - 2 hours

## 4. Data Analysis



- Upload and analyze Data
  - 15 minutes



- Absolute quantification
  - Absolute input copies, based on a standard curve
- Relative quantification
  - Comparative CT method (also known as the  $2^{-\Delta\Delta CT}$  method)
  - Selection of internal control (e.g. untreated control or normal sample)
  - Selection of calibrator
  - Assumes that the PCR efficiency of the target gene is similar to the internal control gene (and that the efficiency of the PCR is close to 1)
  - Fold change =  $2^{-\Delta\Delta CT}$

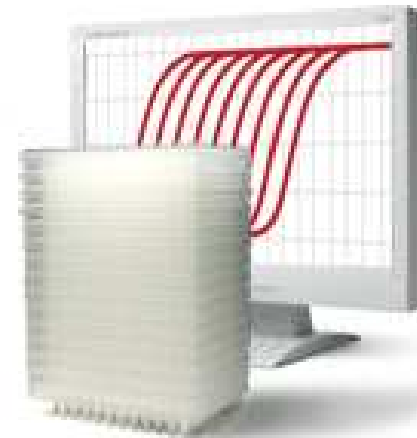
$$= \left[ \{ \text{gene of interest (CT)} - \text{internal control (CT) sample A} \} - \{ \text{gene of interest (CT)} - \text{internal control (CT) sample B} \} \right]$$

- (1) Schmittgen TD, Livak KJ.(2008):Analyzing real-time PCR data by the comparative C(T) method. Nat Protoc.;3(6):1101-8
- (2) Livak, KJ, and Schmittgen, TD.(2001): Analysis of Relative Gene Expression Data Using Real-Time Quantitative PCR and the  $2^{-\Delta\Delta CT}$  Method METHODS 25, 402–408
- (3) [www.Gene-Quantification.info](http://www.Gene-Quantification.info)



# miScript miRNA PCR Array Data Analysis tool

- Web-Based Software
  - No installation needed
- From Raw Ct Values to Fold Change Results
  - Using  $\Delta\Delta C(t)$  Method
- Multiple Analysis Formats
  - Scatter Plot
  - Volcano Plot
  - Multi-Group Plot
  - Clustergram

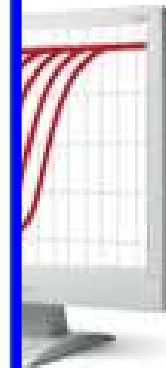
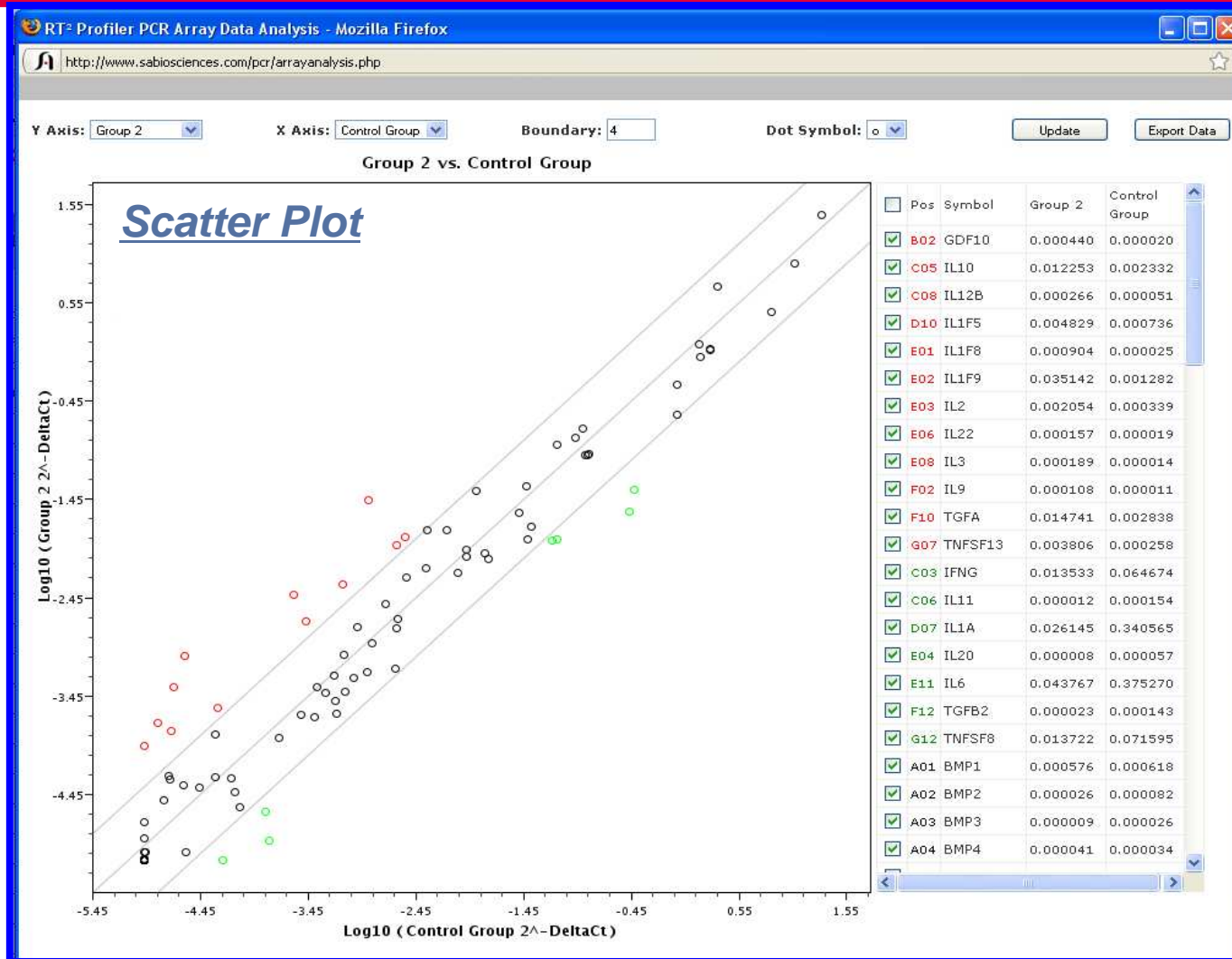


[Also available as downloadable Excel templates](#)



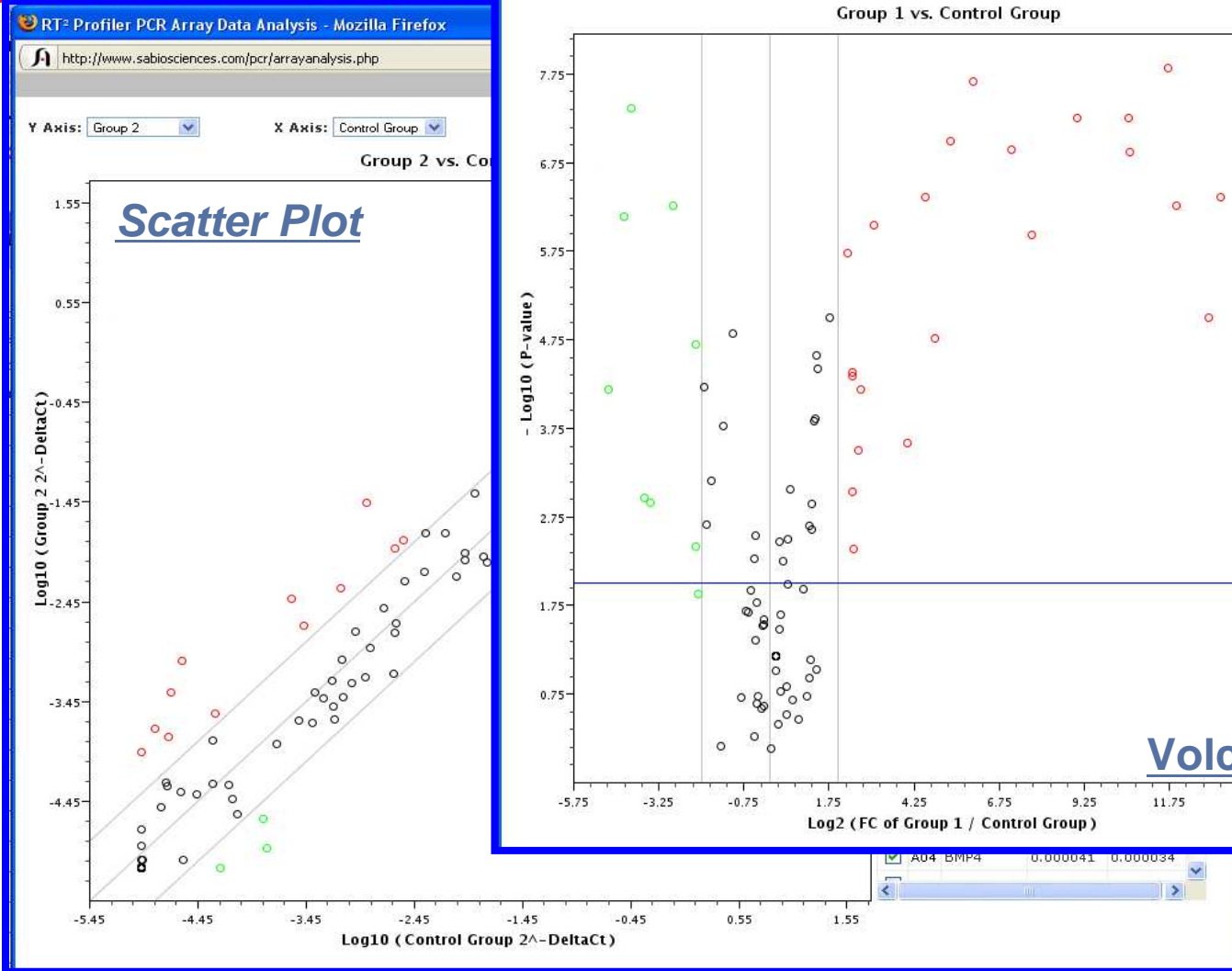
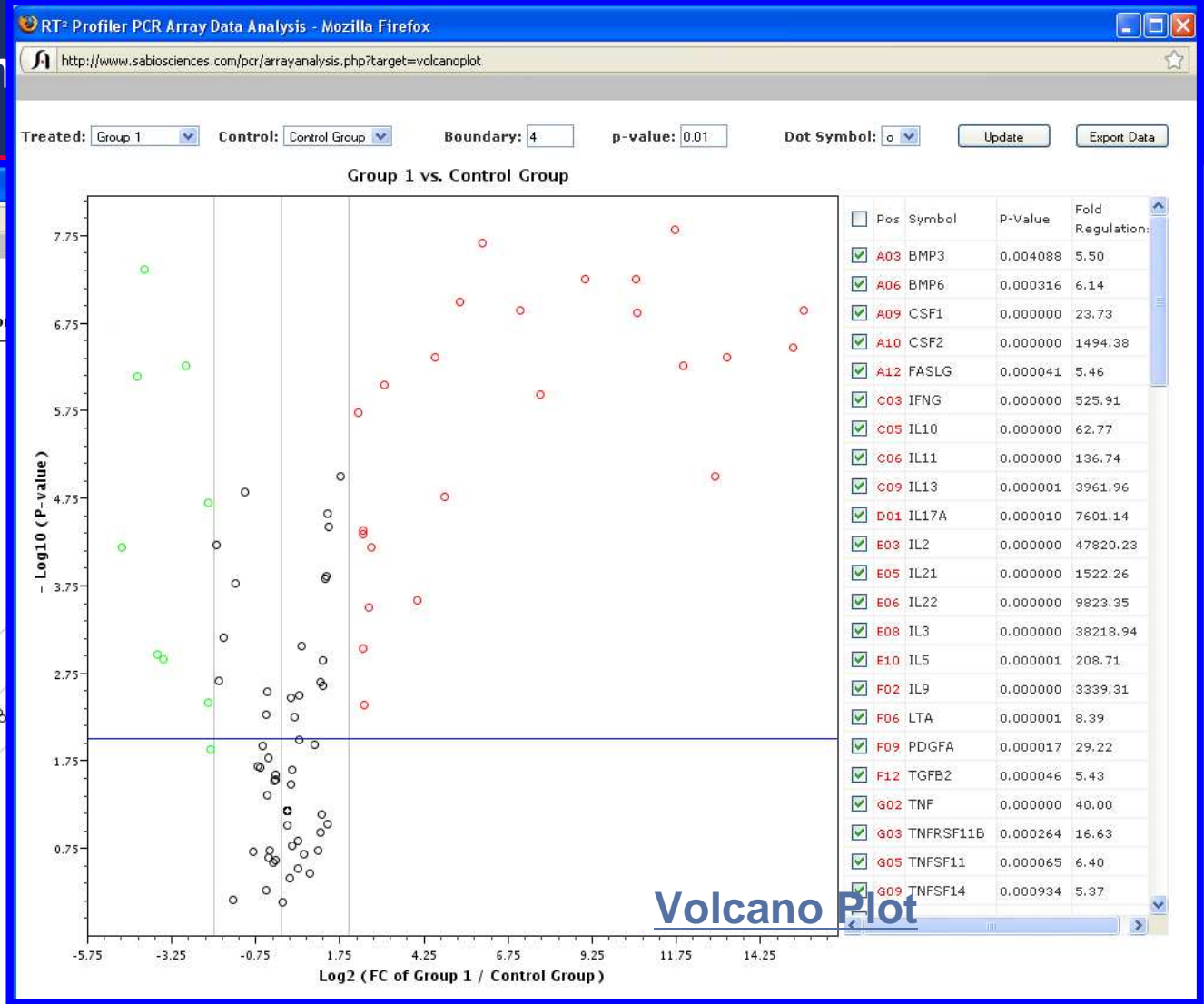


# miScript miRNA PCR Array Data Analysis





miScript m







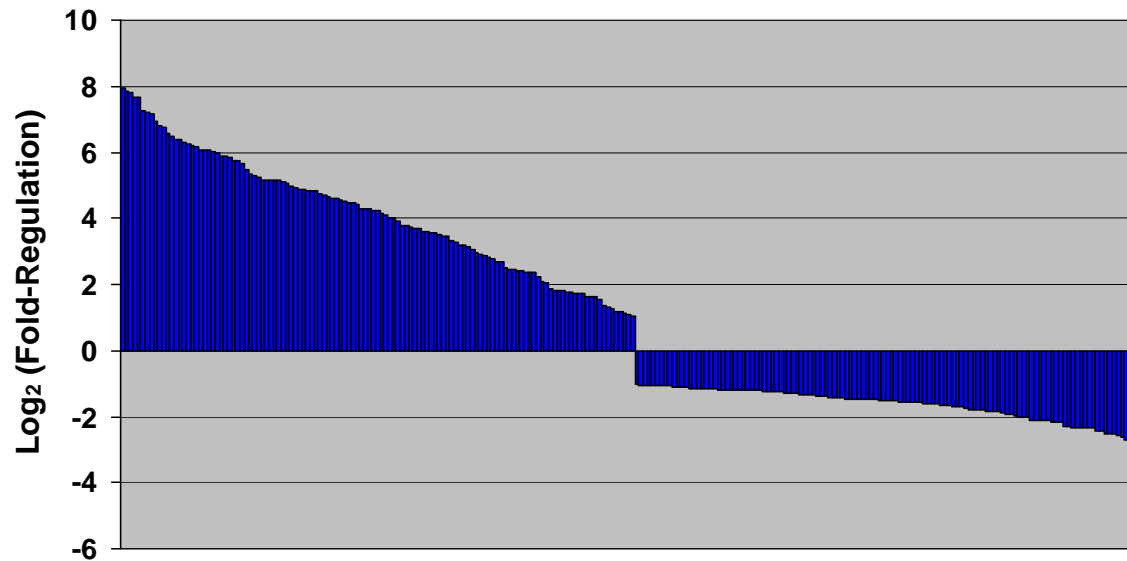
## miRNA profiling of HCT 116 Colorectal Cancer Cells: Following treatment with 5-aza-2'-deoxycytidine (AZDC)

- 5-aza-2'-deoxycytidine (AZDC) is an irreversible inhibitor of DNA methylation.
- AZDC incorporates into DNA & covalently bind to the active site of the DNMT
- HCT116 cells were treated with AZDC, and Untreated cells were used as controls
- Total RNA from each was used to make cDNA with miScript II RT in HiSpec Buffer
- Human whole-miRNome miScript miRNA PCR Array set was used to profile of mature miRNA expression
- Data analysis was performed using the online miScript miRNA PCR Array data analysis tool.

DNMT: DNA Methyl Transferase enzyme



# miRNA profiling of HCT 116 Colorectal Cancer Cells: Following treatment with 5-aza-2'-deoxycytidine (AZDC)

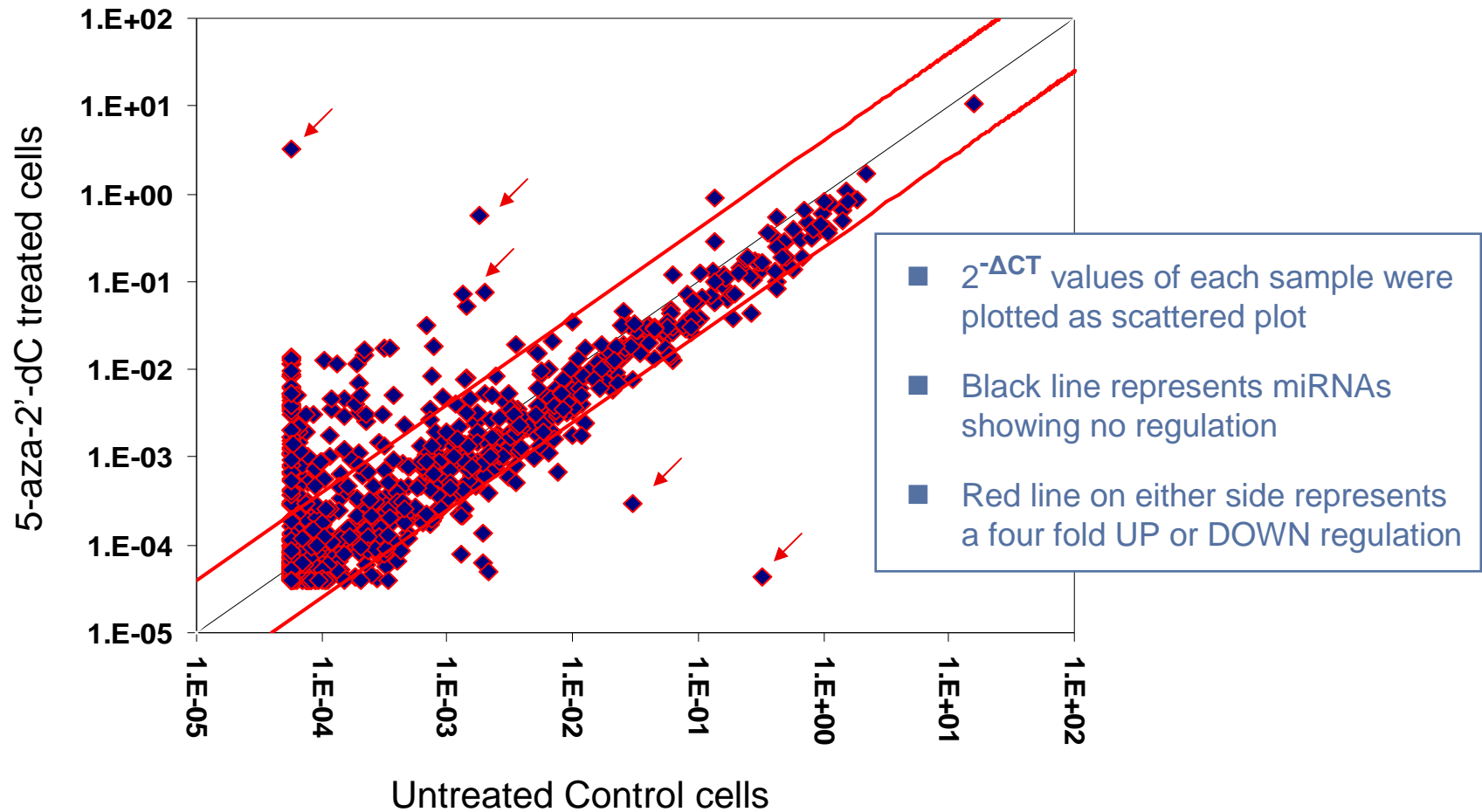


>2 fold UP:	250 miRNAs
>10 fold UP:	164 miRNAs
>2 fold DOWN:	369 miRNAs
>10 fold DOWN:	51 miRNAs

miRNA up regulation likely due to demethylation associated promoter activation

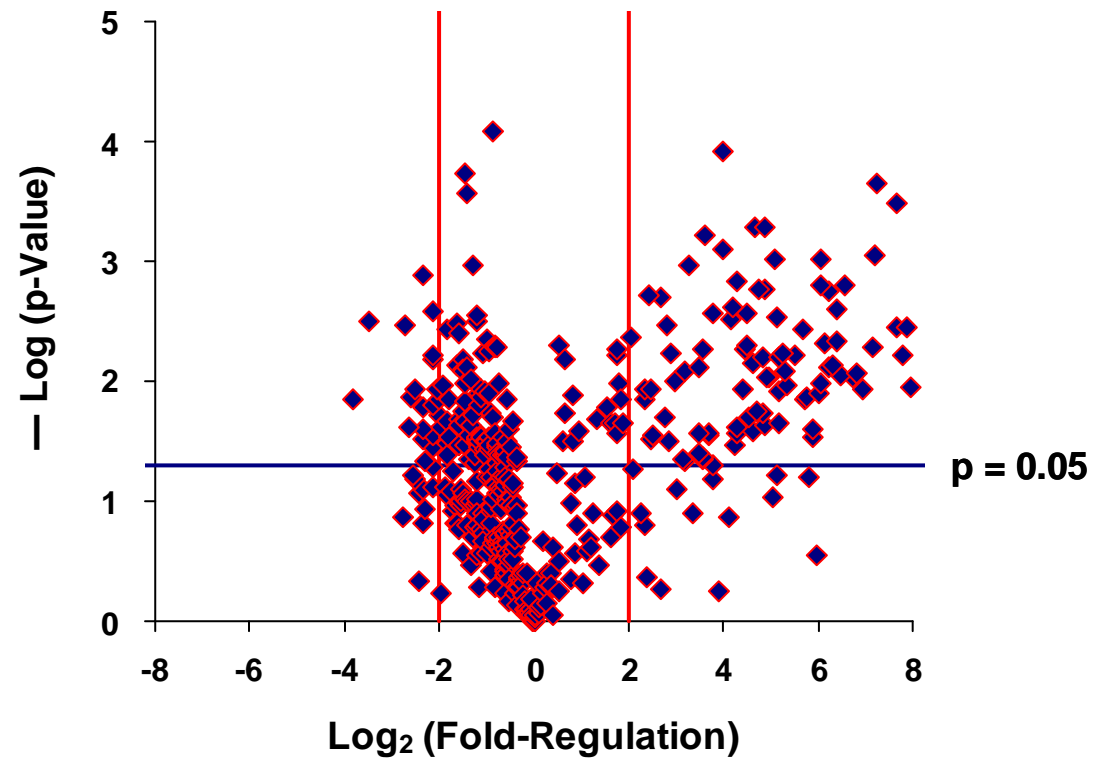


# Scatter Plot: AZDC Treated HCT 116 Colorectal Cancer Cells



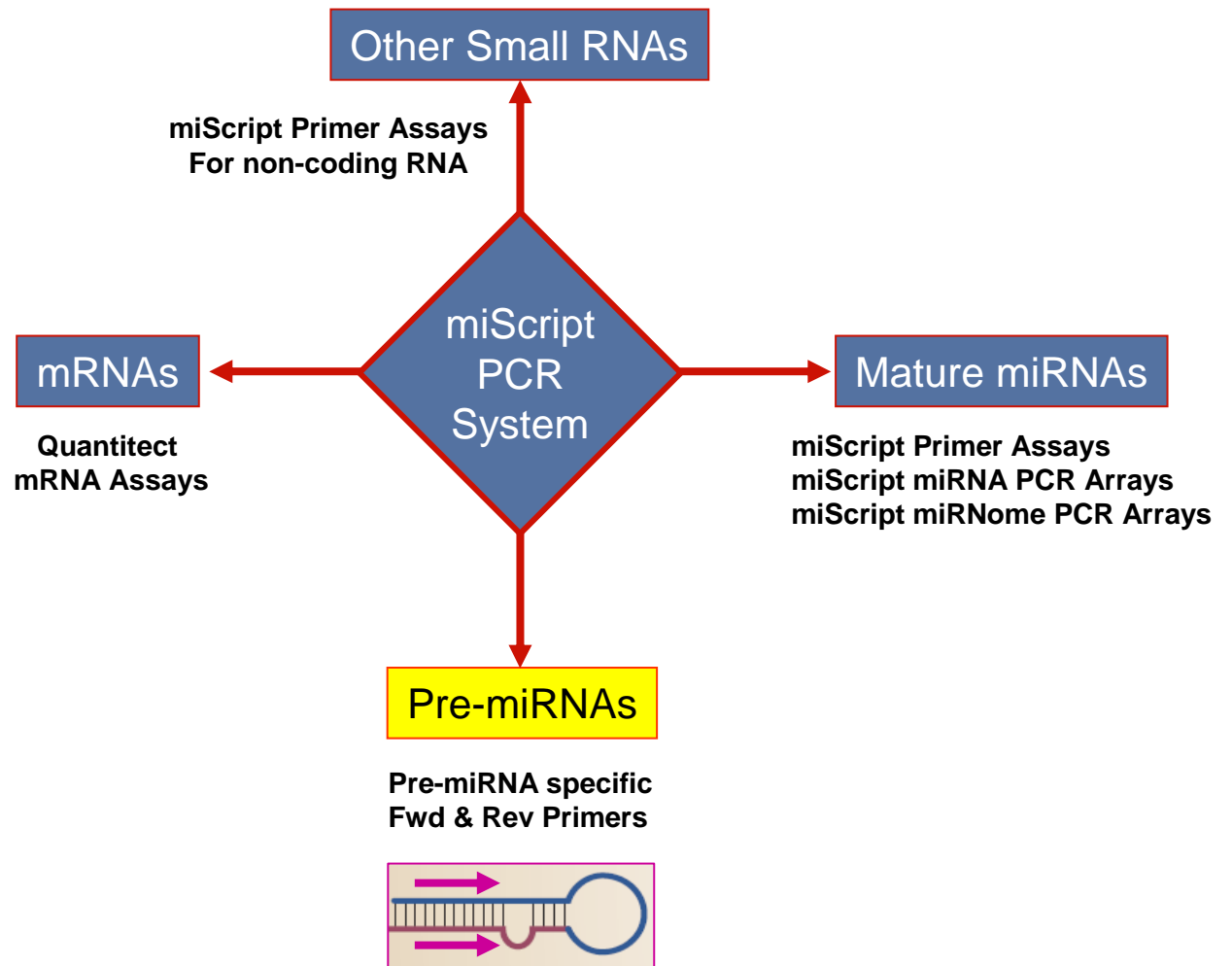


# Volcano Plot: AZDC Treated HCT 116 Colorectal Cancer Cells





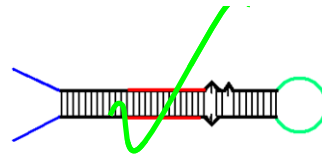
# miScript PCR System: Pre-miRNA Quantification



# Pre-miRNA Stem-Loop Quantification

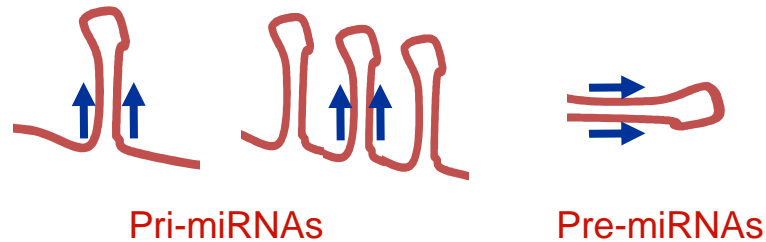
- Study transcriptional and/or post-transcriptional regulation in miRNA biogenesis
- Definitively identify which closely related genomic loci are expressed

## miScript Precursor Assays



- Precursor specific FWD & REV primer
- Specific for precursor miRNAs only

- Precursor assays detect all RNA species that contain the specific stem loop

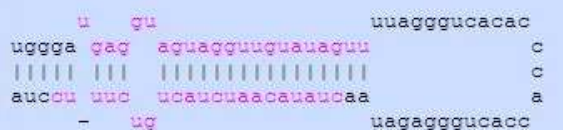




# Expression of Let-7a precursors in various tissues

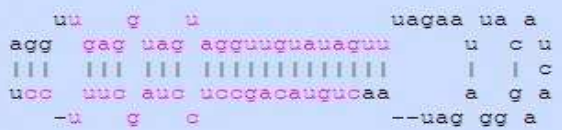
**CHR-9, Intergenic**

Homo sapiens let-7a-1 stem-loop



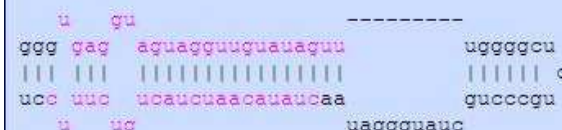
**CHR-11, NR\_024430.1: mir-100-let-7a-2 cluster host gene (non-protein coding)**

Homo sapiens let-7a-2 stem-loop

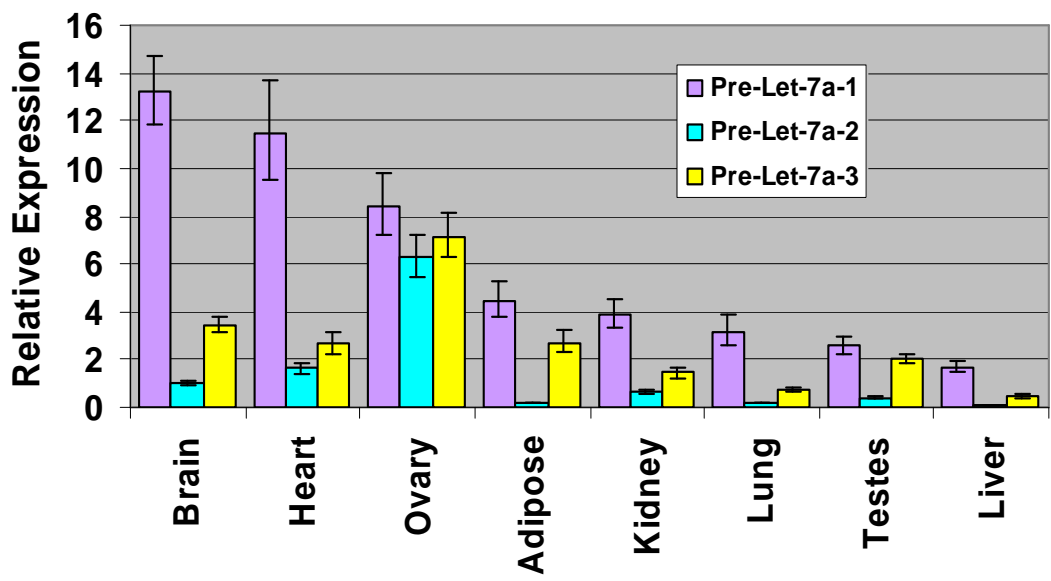


**CHR-22, NR\_027033.1: MIRLET7B host gene (non-protein coding)**

Homo sapiens let-7a-3 stem-loop



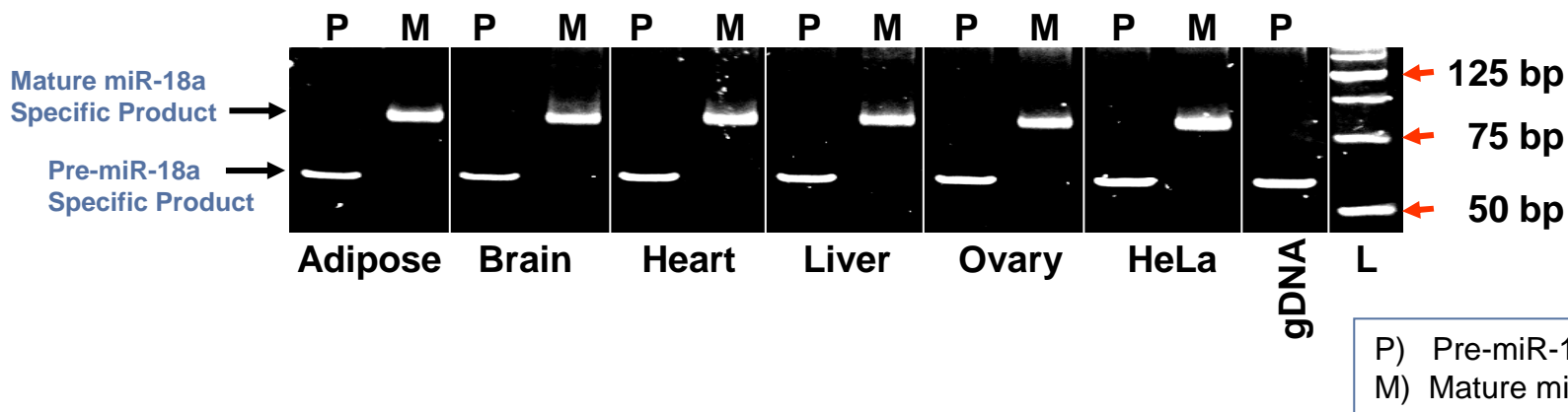
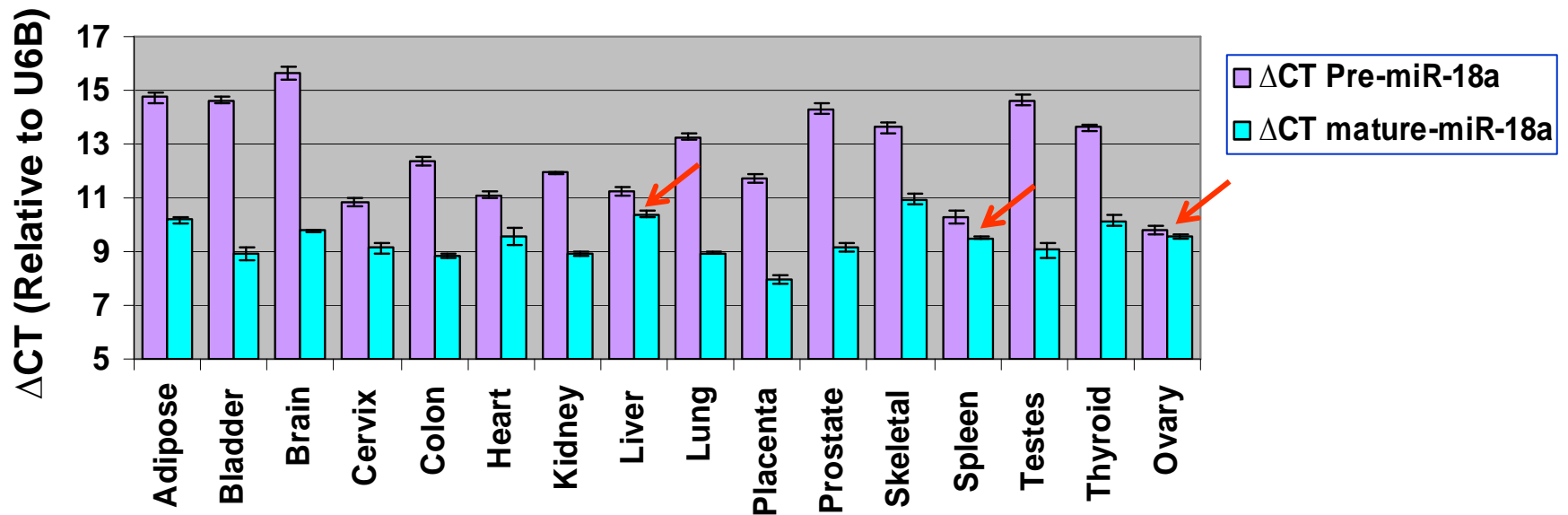
Mature hsa-Let-7a: 5'-UGAGGUAGUAGGUUGUAUAGUU-3'



Relative Expression =  $2^{-\Delta CT} \times 10^5$   
 Normalizer Used: U6B  
 10 ng cDNA /PCR

# Expression of miR-18a and its Precursor

## A member of the miR-17-92 Cluster (13q31.3)





## Summary

- Challenges in miRNA quantification
- miRNA quantification by real time PCR
- Available tools for miRNA profiling experiments
- Critical aspects about template amount requirements
- Examples of miRNA profiling studies
- Precursor detection
  - Measuring the relative levels of mature miRNA and its precursor
  - Identifying the genetic locus responsible for a given miRNA expression in a cell
- miScript miRNA arrays for simple and easy miRNA Profiling
- Data analysis



## GeneGlobe Search Center

Explore and order gene- and pathway-specific products for the following applications:

- ▶ Gene silencing
- ▶ **New!** Gene expression analysis
- ▶ microRNA research
- ▶ Protein research
- ▶ **Updated!** DNA methylation analysis
- ▶ Next-generation sequencing
- ▶ Pathway analysis
- ▶ **New!** PCR arrays for disease research

Search Custom Products New [Guided Tour](#)

Enter search term(s)

- Entrez Gene IDs (e.g., 835)
- RefSeq IDs (e.g., NM\_032983, NP\_116765)
- Gene symbols (e.g., CASP2)
- Cat. No. (e.g., SI00299551, QT01342509)
- Sanger ID or Accession (e.g., hsa-let-7b, MI0000063)

▶ or upload a list of search terms in a file

Species

- Human (Homo sapiens)
- Mouse (Mus musculus)
- Rat (Rattus norvegicus)
- Arabidopsis (Arabidopsis thaliana)
- Chicken (Gallus gallus)
- Dog (Canis familiaris)
- Fruit fly (Drosophila melanogaster)
- Nematode (Caenorhabditis elegans)
- Zebrafish (Danio rerio)
- Frog (clawed frog) (Xenopus laevis)
- Cow (Bos taurus)
- Epstein Barr virus
- Human cytomegalovirus
- Human immunodeficiency virus 1
- Kaposi sarcoma-associated herpesvirus
- Mouse cytomegalovirus
- Mareks disease virus
- Mareks disease virus type 2
- Mouse gammaherpesvirus 68
- Rhesus lymphocryptovirus
- Rhesus monkey rhadinovirus



# Try free miScript PCR Array Starter Pack



Experience PCR Array Performance → Try them today!

## miScript miRNA PCR Array Starter Pack

Promo Code: MS 11

FREE PCR Arrays of any Pathway

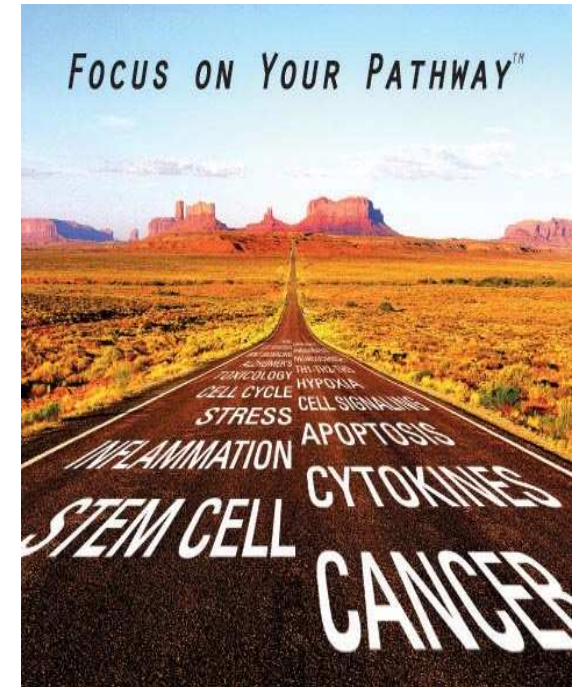
- 2 96-well/100-well (2 samples) OR
- 1 384-well (4 samples)
- With Purchase of:
  - miScript II Reverse Transcription Kit
  - miScript SYBR Green PCR Mastermix (2-Pack)

To try the free Starter Pack, click the link below

<http://www.sabiosciences.com/promotion/miscriptdemo.php>

## Receive your Free iPod Shuffle by

- Sharing your data with us
- Contact: [sab.us@qiagen.com](mailto:sab.us@qiagen.com) to claim your gift



Thousands of scientists have discovered the power of PCR Arrays. Join them on the road to success!



## Questions

**Webinar 3:** Profiling miRNA Expression:  
On the Road To Biomarker Development

**Date:** October 19, 2011, 1:00 pm EST

**Speaker:** Eric Lader Ph. D.  
Senior Director, QIAGEN R&D

**Thank you!**

**Subu Yerramilli**  
**Subu.yerramilli@qiagen.com**

For up-to-date licensing information and product-specific disclaimers, see the respective QIAGEN kit handbook or user manual. QIAGEN kit handbooks and user manuals are available at [www.qiagen.com](http://www.qiagen.com) or can be requested from QIAGEN Technical Services or your local distributor