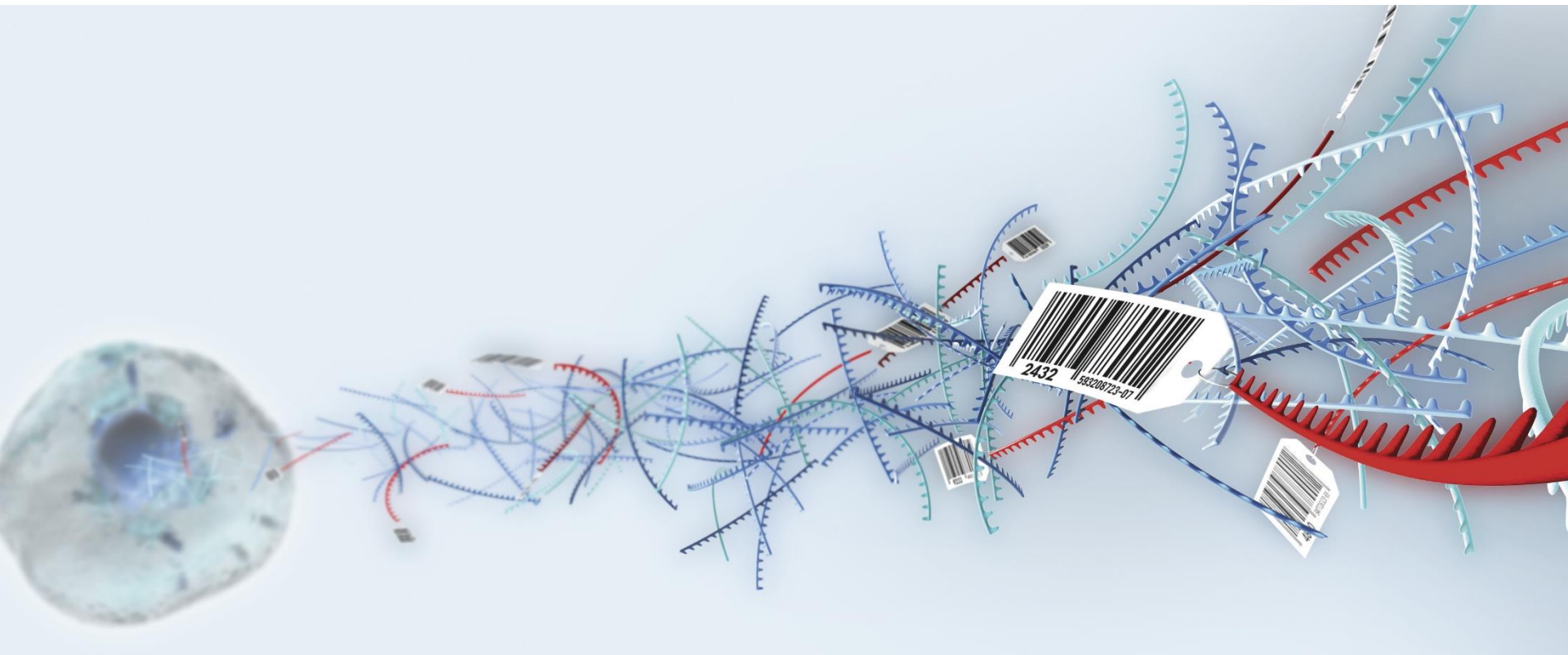


# QIAseq miRNA Library Kit

The newest solution for gel-free miRNA sequencing



Jonathan Shaffer, Ph.D.

[Jonathan.Shaffer@qiagen.com](mailto:Jonathan.Shaffer@qiagen.com)

Associate Director, Product Development



QIAGEN products shown here are intended for molecular biology applications. These products are not intended for the diagnosis, prevention or treatment of a disease.

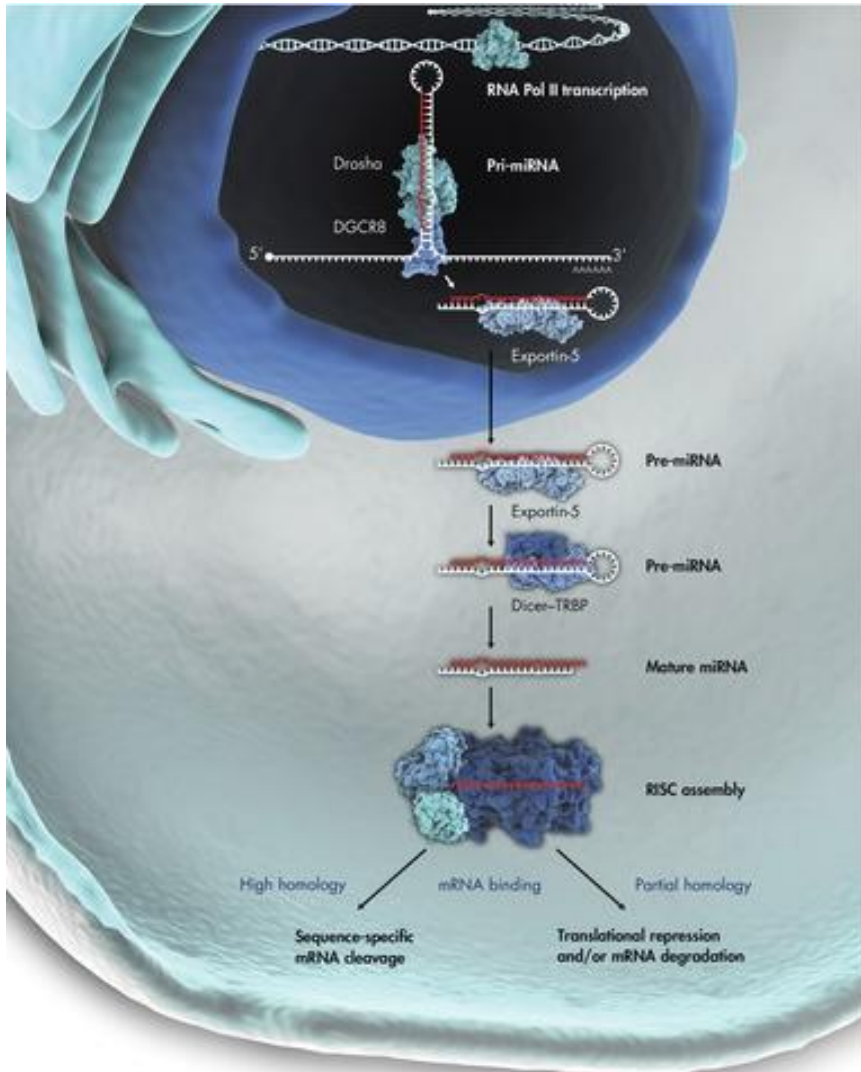
For up-to-date licensing information and product-specific disclaimers, see the respective QIAGEN kit handbook or user manual. QIAGEN kit handbooks and user manuals are available at [www.qiagen.com](http://www.qiagen.com) or can be requested from QIAGEN Technical Services or your local distributor.



- 1 miRNA background
- 2 QIAsiq miRNA sequencing: Product overview
- 3 Data analysis overview
- 4 Performance and application data
- 5 Summary and questions



- 1 miRNA background
- 2 QIAseq miRNA sequencing: Product overview
- 3 Data analysis overview
- 4 Performance and application data
- 5 Summary and questions



miRNAs are ~21 nucleotide, small non-coding RNAs that are expressed in most tissues

Changes in miRNA expression can be correlated with gene expression changes in:

- Development
- Differentiation
- Signal transduction
- Infection
- Aging
- Disease

Role in cancer: miRNAs are associated with cell proliferation, resistance to apoptosis, invasiveness and differentiation in cancer cells

# Why quantify miRNAs?

Using a signature of altered miRNA expression to differentiate cancer tissue from normal tissue

Using miRNA-based classifier to identify the tissue of origin for cancers of unknown primaries

Profiling circulating blood or tumor-derived exosomal miRNAs, surpassing the invasive procedures to aid in early detection of cancers



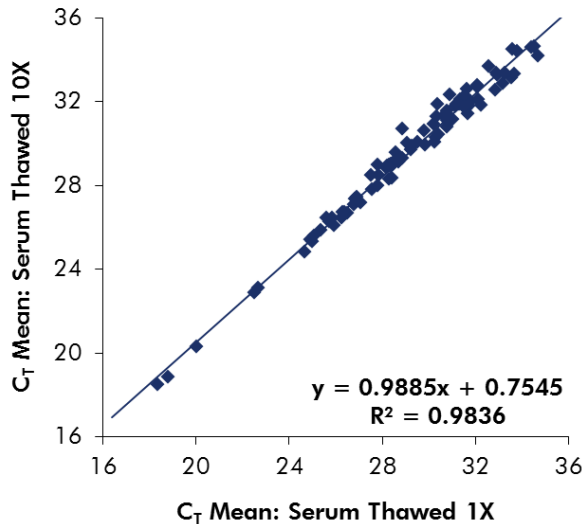
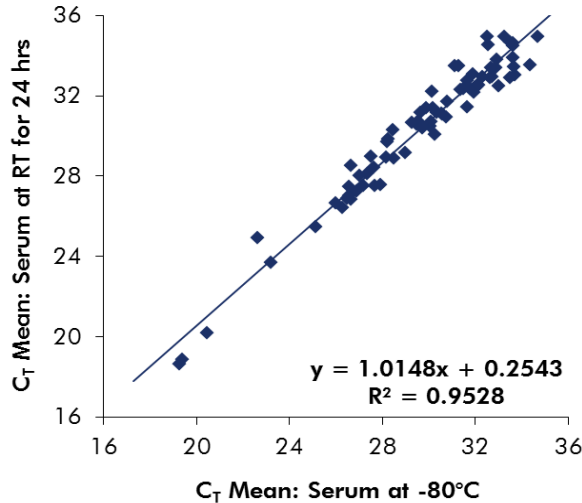
Distinguishing tumor subtypes using a panel of miRNAs that show differential expression within one cancer type

Studying SNPs in the miRNA genes, miRNA binding sites in the target mRNA genes or in the miRNA processing/machinery pathway genes to predict cancer predisposition

Source: Paranjape. et. al. (2009) Gut.

Quantifying miRNAs help identify novel expression patterns for miRNAs – both annotated (i.e. in miRBase) and non-annotated (i.e. identified using small RNA sequencing).

# Why are circulating miRNAs promising biomarker candidates?



## Expressed in most biofluids

- Serum, plasma, CSF, urine and saliva
- Robustly expressed in serum and plasma

## Minimally invasive and easily liberated

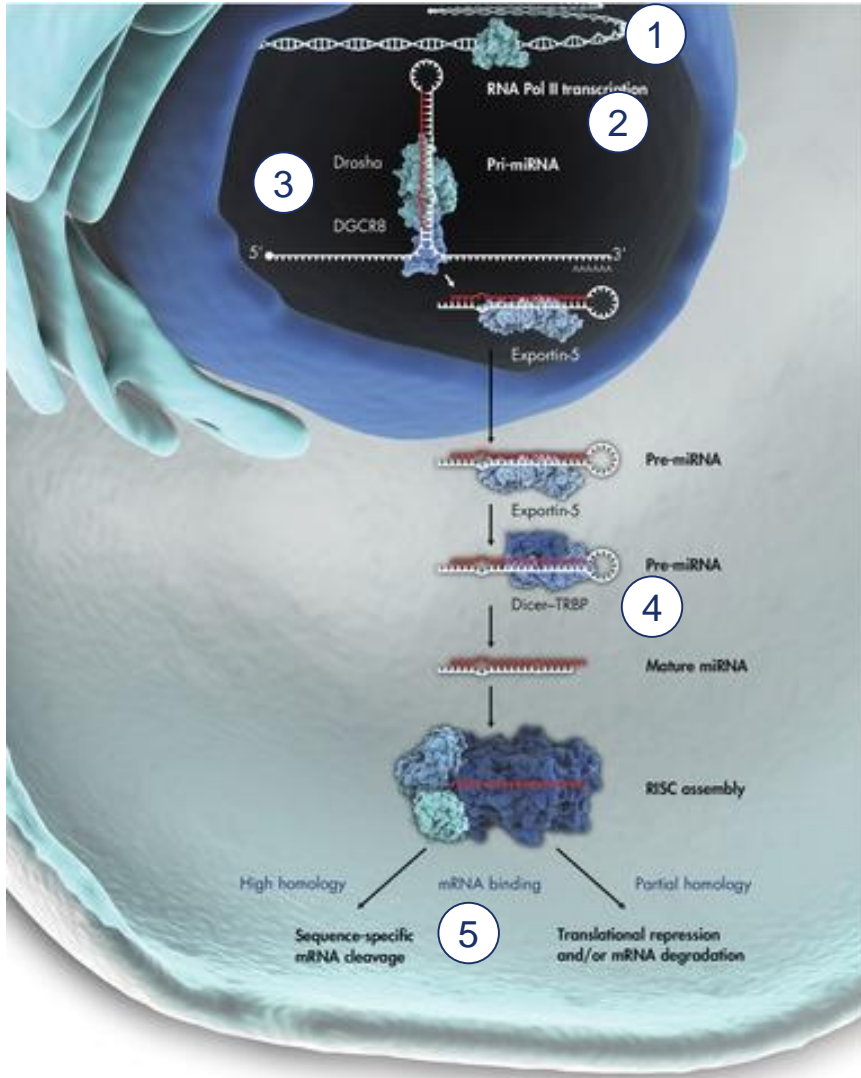
- Sensitive, specific and precise platforms exist for miRNA quantification

## Protected from degradation

- Resistant to freeze-thaw cycles
- Stable at room temperature for 24 hours

● miRNA expression is associated with normal and disease biology.

## Disruption of miRNA–mRNA interaction



- ① Genomic instability
  - Amplification or deletion
  - Translocation
  - Insertional mutagenesis
- ② Altered transcription
  - Methylation
  - Histone modification
  - Transcription factor
- ③ Drosha processing
- ④ Dicer processing
- ⑤ Loss of miRNA binding site in target mRNA
  - SNP or mutation
  - Alternative splicing
  - Loss or change of 3'-UTR



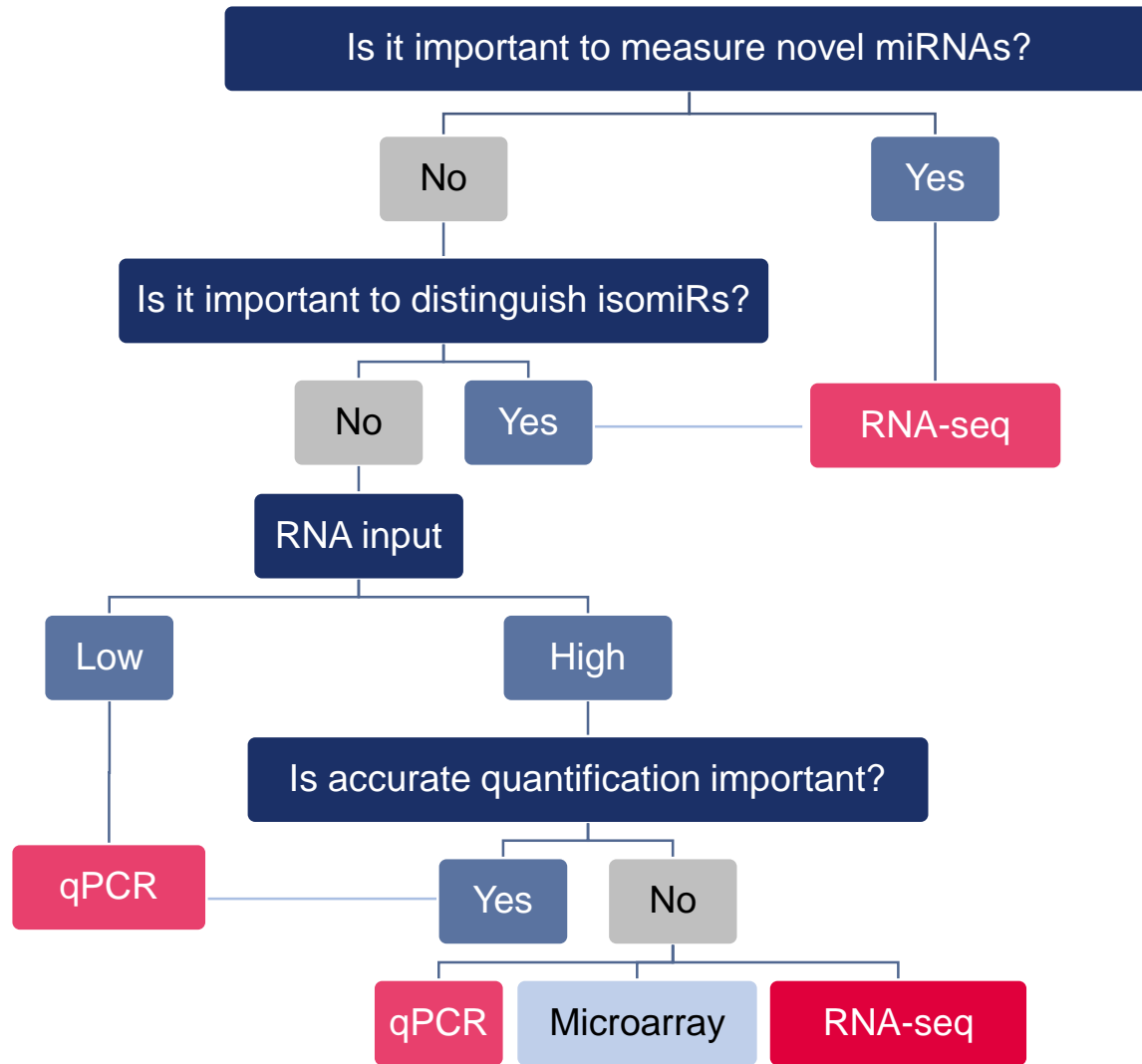
## miRNAs as biomarkers

- Tissue-based miRNA profiling for biomarker discovery
- miRNA profile-based classification of tissues of unknown origin
- **Circulating miRNA biomarkers**
- Forensics

## Understanding gene regulation

- Developmental biology
- Novel miRNA discovery
- Studying miRNA-mRNA and miRNA-protein interactions
- **Total RNA discovery** – integrative analyses of miRNAs in the context of gene regulatory networks

Source: Pritchard, C. C. et. al. (2012) Nature Rev. Genet. 13, 358-369



Source: Pritchard, C. C. et. al. (2012) Nature Rev. Genet. 13, 358-369

Inconvenient  
sample  
inputs

- Kits require up to 1  $\mu\text{g}$  of sample input and not ideal with lower inputs
- May not be compatible or optimized for biofluids

Poor data  
quality

- On target miRNA NGS reads account for only 20–30% of total reads
- Adapter dimers eat away your read budget
- Gel band excision is non-precise (not 100% miRNA)
- Biofluid samples may be contaminated with similar-sized RNA types

Tedious  
workflow

- Gel-based workflows take multiple days to get to sequencing
- Large sample numbers become tedious with gel-based workflows
- Optimization of sample isolation is not done

miRNA  
sequencing

Novel miRNA  
discovery and  
sample screening

Real-time  
PCR

NGS validation and  
expression analysis



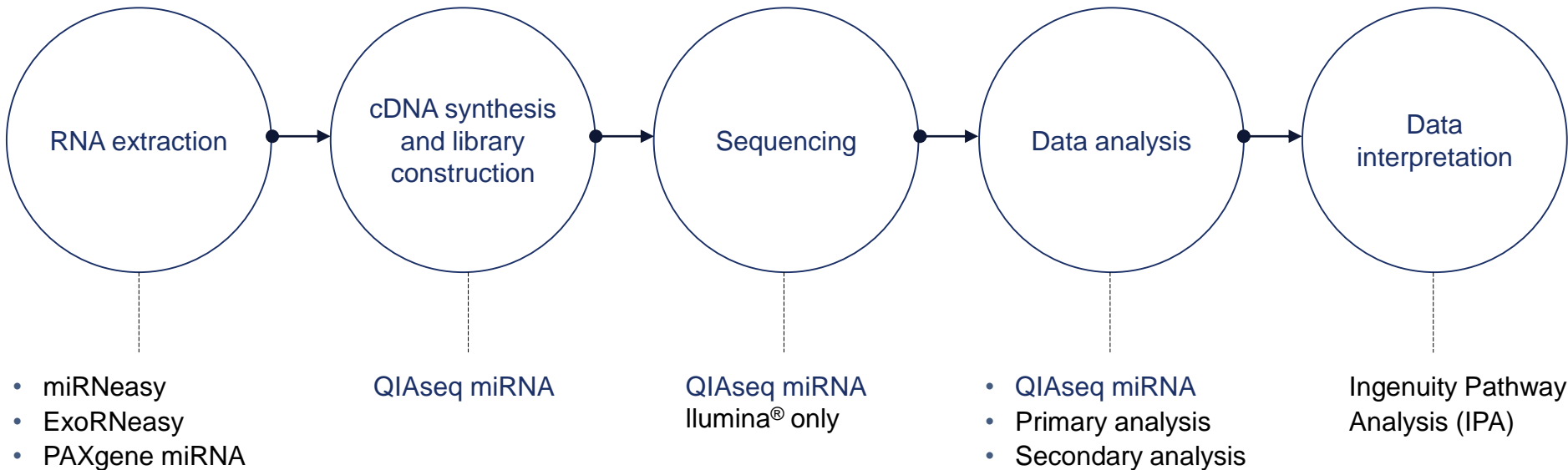
- Novel miRNA discovery
- Sample screening
- Accurate quantification
- Broad RNA input range
- No gels



- 1 miRNA background
- 2 QIAseq miRNA sequencing: Product overview**
- 3 Data analysis overview
- 4 Performance and application data
- 5 Summary and questions

Sample

Insight





## Cells, fresh tissue and frozen tissue

- miRNeasy Mini Kit
- miRNeasy Micro Kit
- miRNeasy 96 Kit

## FFPE tissue

- miRNeasy FFPE Kit

## Fluids (serum, plasma, urine, CSF and saliva)

- miRNeasy Serum/Plasma Kit

## Exosome enrichment/isolation from serum/plasma

- exoRNeasy Serum/Plasma Midi Kit
- exoRNeasy Serum/Plasma Maxi Kit

## PAXgene Blood miRNA Kit

## PAXgene Tissue miRNA Kit

## QIAseq miRNA Library Kit

- 12 rxn: 331502
- 96 rxn: 331505
- What is included? 3' ligation, 5' ligation, reverse-transcription, cDNA cleanup, library amplification, library cleanup reagents and quality control primers

## QIAseq miRNA NGS 12 Index

- 12 rxn: 331592
- What is included? Sequencing adapters, primers and indices compatible with Illumina platforms – 12 indices for 12 samples

## QIAseq miRNA NGS 48 Index

- 96 rxn: 331595
- What's included? Sequencing adapters, primers and indices compatible with Illumina platforms – two 48 indices for 96 samples





## What is the kit?

miRNA-focused next-generation sequencing library prep kit and integrated bioinformatics/data analysis solution

- Compatible with Illumina sequencers

## What can be done with the sequencing data?

- Differential expression calculations of miRNA from highly multiplexed samples
- Novel miRNA discovery

## What are the distinguishing features of the prep kit?

- Gel-free, rapid workflow
- Broad RNA input: 1–500 ng
  - No adapter dimers at any RNA input amount
- Library prep from serum, plasma, biofluids, cells and tissues (any species)
- Integrated Unique Molecular Index (UMI) technology
- Highly optimized chemistry
- All-in-one-box solution

● QIAseq miRNA Library Kit: Unparalleled miRNA-focused sequencing for robust miRNA quantification and discovery.

## miRNA mapping rates routinely observed

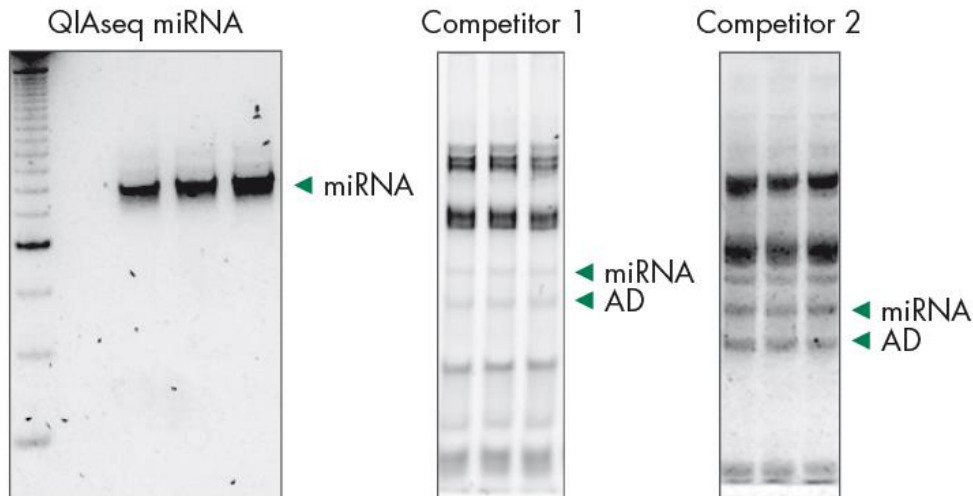
	QIAseq miRNA Kits	Competitors' kits
<b>Cell lines</b>	50–60% or greater	20–30% average post gel
<b>Tissues</b>	75% or greater	30–40% average post gel
<b>Serum/plasma</b>	15–30% or greater	1–15% average post gel

## QIAseq miRNA Kit: Specifications

- **Sample type:**
  - Cells, fresh or frozen tissue, FFPE tissue, serum/plasma and other biofluids
  - Animal and plant samples
  - Any species
- **Total RNA input range (cells/tissues):** 1–500 ng
- **Total RNA input recommendation (serum/plasma):** 5 µl when RNA has been isolated from 200 µl of sample
- **What RNAs are included in library prep?**
  - Highly optimized for miRNA
  - piRNAs will also be efficiently sequenced
- **Multiplex capability:** 48 samples
- **Sequencer compatibility:** Illumina
- **Total library construction time:** 8 hours (3 hours hands-on)

- QIAseq miRNA vs. two competitor (C) kits
  - RNA amounts: 100 ng (QIAseq miRNA), 1  $\mu$ g (C1) and 100 ng (C2)

## PAGE gel after standard library prep protocol



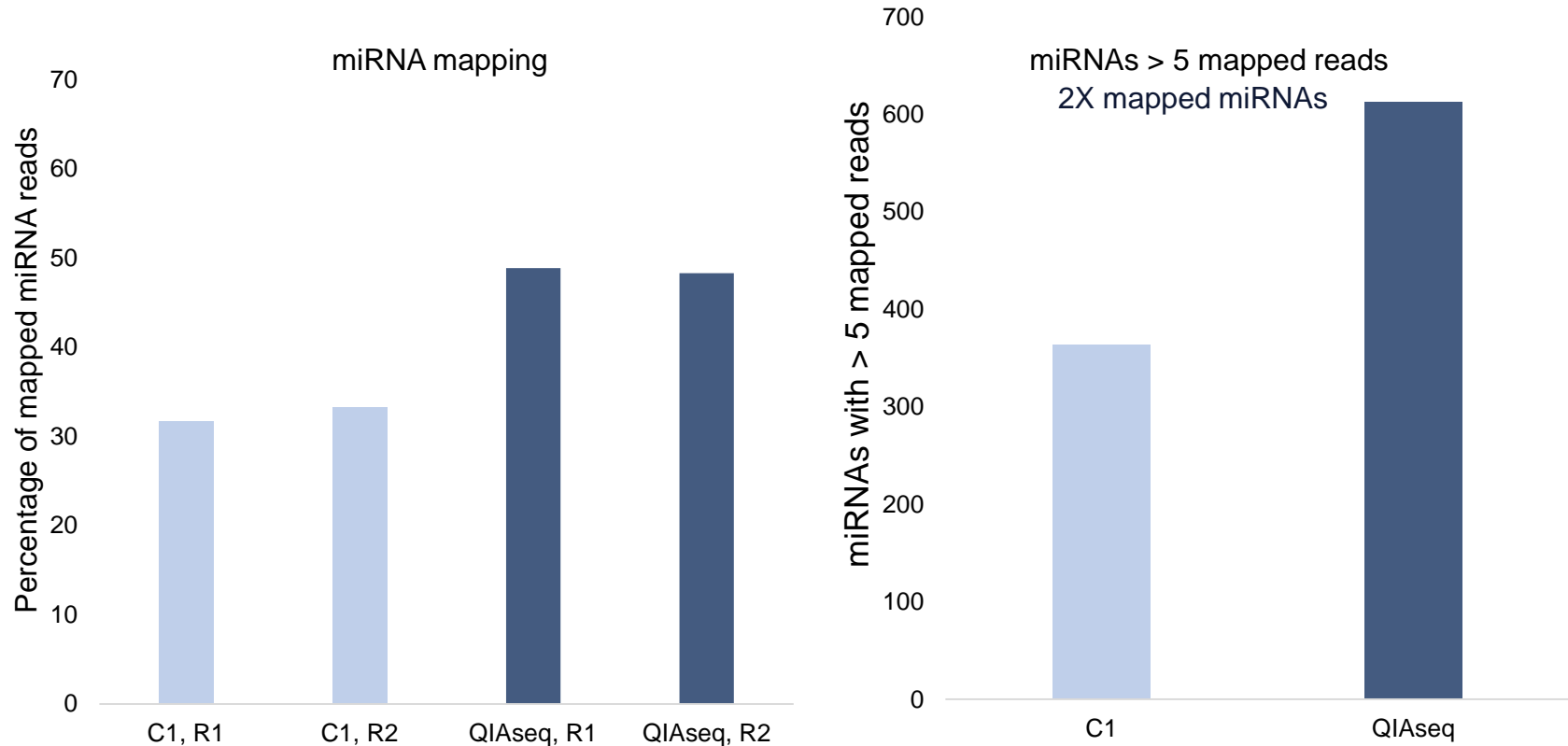
AD: Adapter dimer

- With the QIAseq miRNA standard protocol, a robust, specific miRNA library is generated with negligible background. Other commercial options are fraught with side-products, including adapter dimers.

# QIAseq miRNA Library Kit outperforms the competition

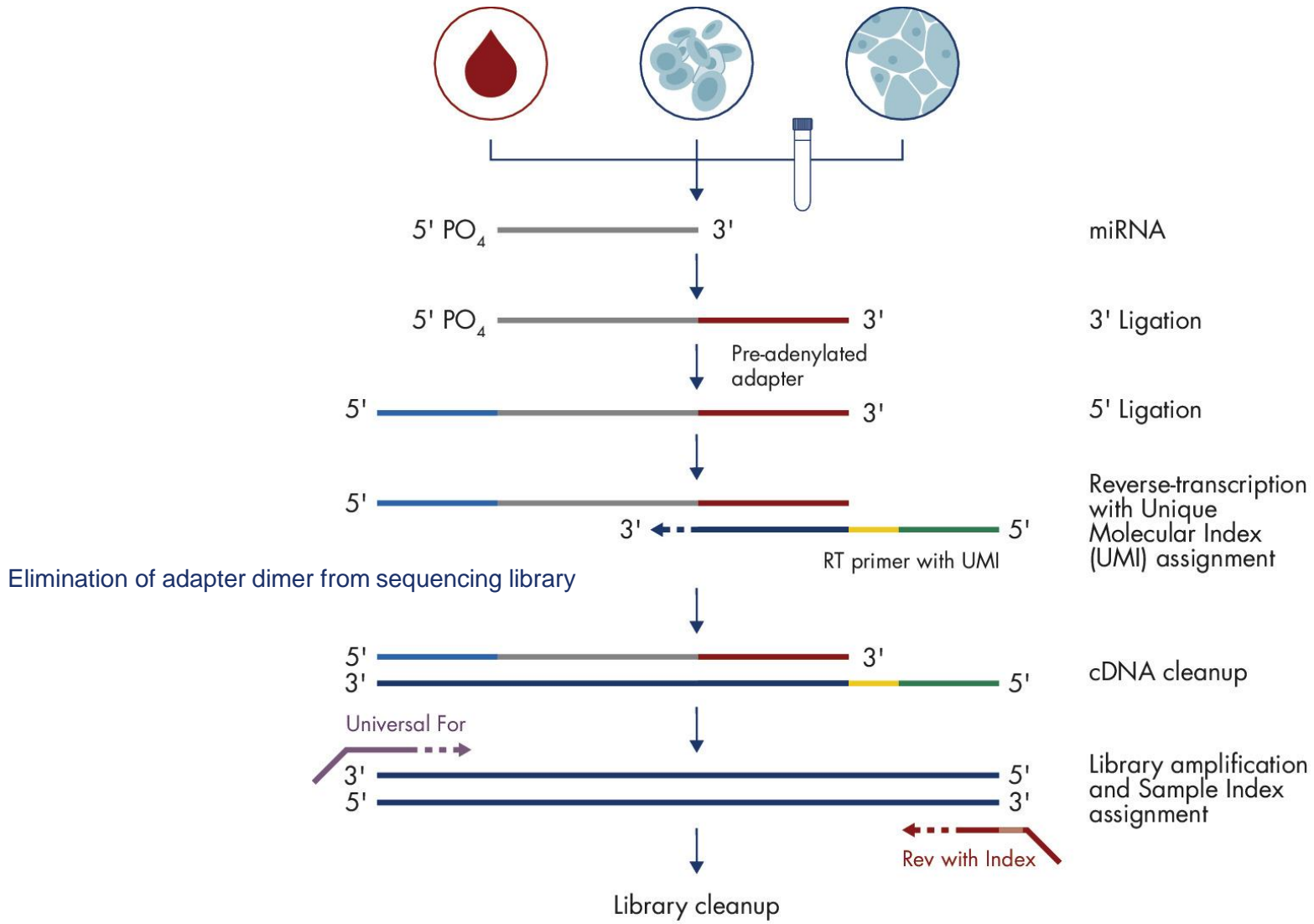
## Next-generation sequencing: QIAseq miRNA and competitor 1 (C1)

- For C1, prior to sequencing, miRNA library was excised and purified from a PAGE gel
- MiSeq: 75 bp Single-Read (QIAseq miRNA) and 50 bp Single-Read (C1)



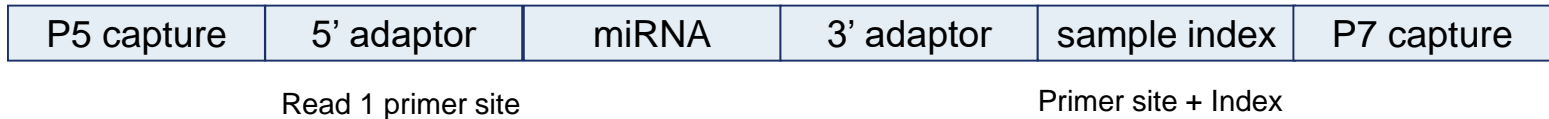
With QIAseq miRNA, increase your mapped miRNA reads (as a result of reduced bias and improved sensitivity) while reducing your workflow time.

# QIAseq miRNA: Library construction



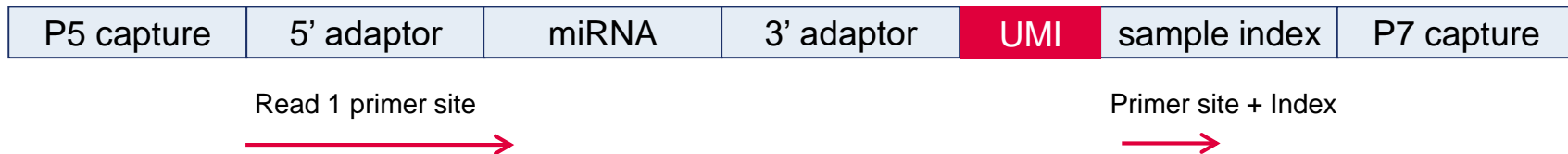
Library pre-seq QC, determining library concentration, preparation for sequencing and data analysis

## Other kits



- 50 bp read
- 48 possible sample indices

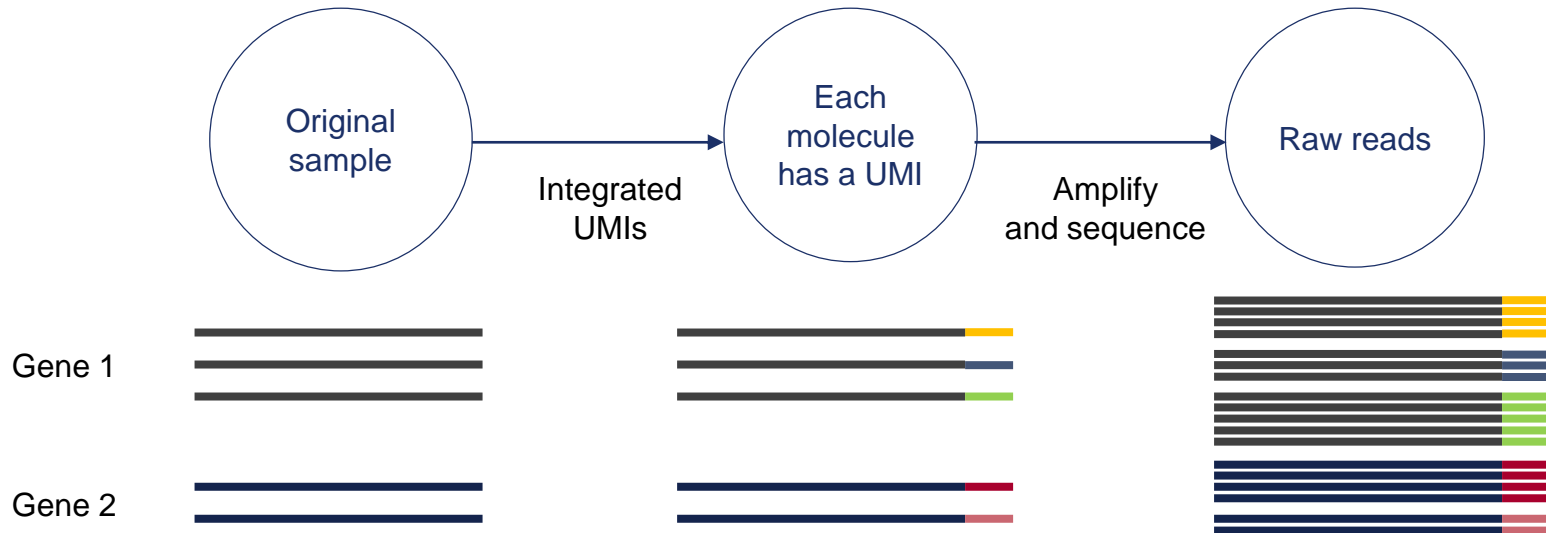
## QIAseq miRNA Library Kit



- 75 bp read to make use of the UMI
  - Note: 50 bp read is possible, but the UMI will not be sequenced
- 48 possible sample indices

● UMI enables precision quantification of miRNA molecules following next-generation sequencing.

# The principle of Unique Molecular Indices (UMIs)



Original sample (3:2 ratio of gene 1 to gene 2)

- Gene 1: 3 molecules
- Gene 2: 2 molecules

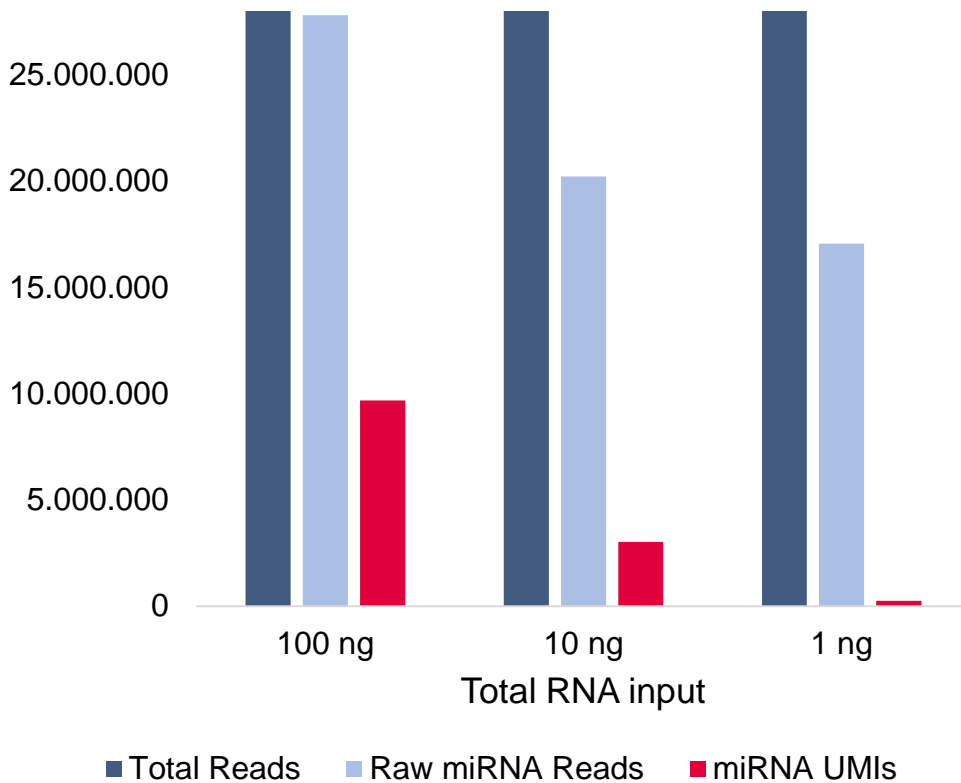
Interpretation of **raw reads** (2:1 of gene 1 to gene 2)

- Gene 1: 12 reads
- Gene 2: 6 reads

Interpretation of **UMIs** (3:2 ratio of gene 1 to gene 2)

- Reads are collapsed based on molecule counts
- Gene 1: 12 reads but 3 molecules are identified due to UMIs
- Gene 2: 6 reads but 2 molecules are identified due to UMIs

● Quantification based on UMIs reflects quantities of original RNA molecules.



## Assessment of raw miRNA reads

- Sequencing of the same miRNA molecule over and over results in an overestimation of miRNA expression
- The lower the RNA input, the lower the effect

## Assessment of miRNA UMIs

- Individual miRNA molecules are being counted, resulting in a true assessment of miRNA expression
- The lower the RNA input, the more powerful the UMIs

● UMIs give a true readout of miRNA expression.



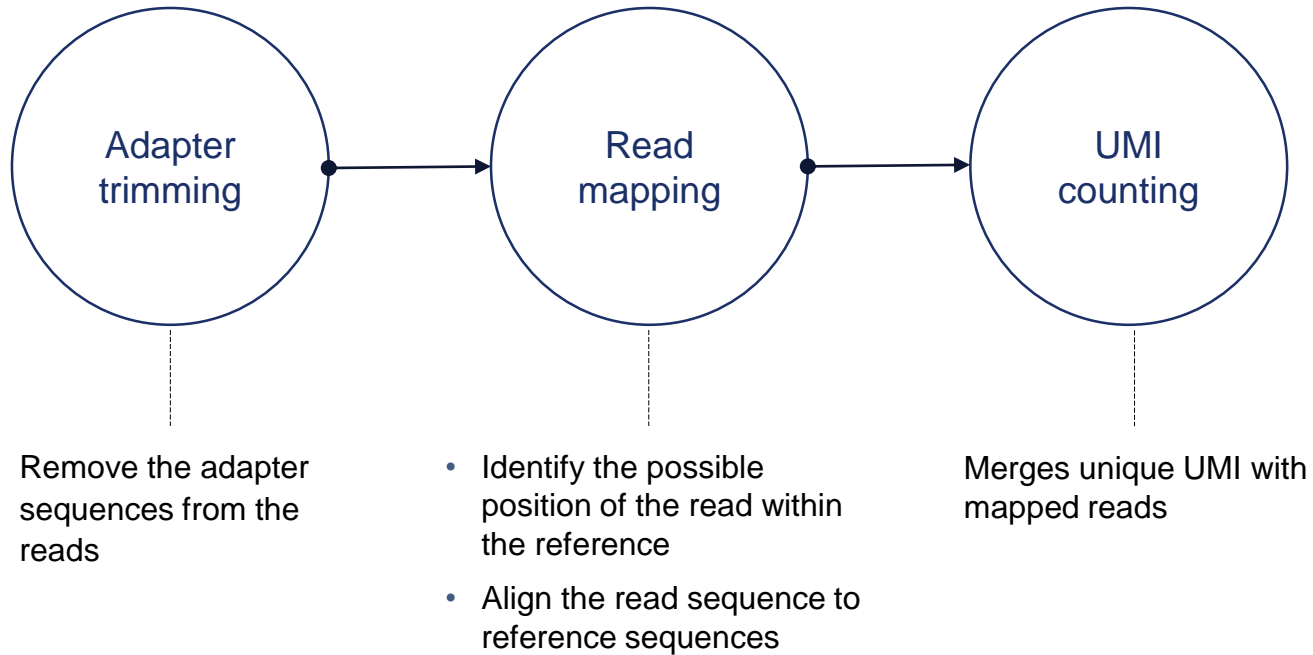


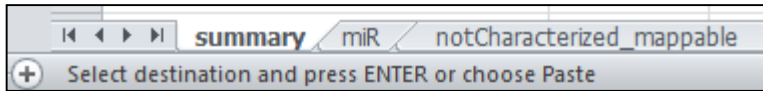
- 1 miRNA background
- 2 QIAsseq miRNA sequencing: Product overview
- 3 Data analysis overview**
- 4 Performance and application data
- 5 Summary and questions

- Free, easy to use primary data analysis: [www.qiagen.com/geneglobe](http://www.qiagen.com/geneglobe)
- Species: Human, mouse, rat or other species (all of miRBase)

[Shop](#)[Products](#)[Resources](#)[Support](#)[About QIAGEN](#)[File Upload](#)[File Management](#)[miRNA Quantification Jobs](#)[Add files...](#)[Start upload](#)[Cancel upload](#)

**Note:** Acceptable file extensions are ".fastq" or ".fastq.gz" for Illumina reads, and ".basecaller.bam" for Ion reads. Please submit only unaligned base-caller bam files generated by Torrent Server 3.4.1 or higher. Please do not submit aligned bam files. Please **DO NOT** refresh the browser or navigate to other pages while uploading files.





## Summary tab

	A	B	C
1	read set	Sample 1, Rep 2	Sample 2, Rep 1
2	total_reads	3,485,926	3,454,577
3	no_adapter_reads	245,252	321,093
4	too_short_reads	916,139	737,228
5	UMI_defective_reads	143,102	192,311
6	miRNA_Reads	1,250,002	1,333,379
7	hairpin_Reads	3,319	2,787
8	piRNA_Reads	26,771	29,049
9	rRNA_Reads	99,637	93,880
10	tRNA_Reads	22,192	18,248
11	mRNA_Reads	12,305	12,127
12	otherRNA_Reads	115,790	149,227
13	notCharacterized_Mappable	143,611	135,715
14	notCharacterized_notMappable	507,806	429,533

## miR tab

	A	B	C	D	E
1	miRNA	Sample 1, Rep 2 - UMI	Sample 2, Rep 1 - UMI	Sample 1, Rep 2 - Reads	Sample 2, Rep 1 - Reads
2	hsa-miR-16-5p	112,095	96,568	224,770	196,617
3	hsa-miR-126-3p	80,983	77,107	164,836	160,661
4	hsa-let-7a-5p	43,816	39,426	84,078	76,639
5	hsa-miR-223-3p	36,192	50,904	71,624	104,480

## Features of the primary analysis report

- **Summary tab:** Important mapping metrics for the sequencing run
- **miR tab:** Mapped reads and UMI counts for each miRNA found in miRBase
- **notCharacterized mappable tab:** Mapped reads and UMI counts for each sequence that aligns to the genome

QIaseq miRNA primary analysis output gives a complete summary of important sequencing metrics and UMI counts for each mapped miRNA.



The screenshot shows the QIAGEN Data Analysis Center website. At the top, there is a navigation bar with links for Shop, Products, Resources, Support, About QIAGEN, and Careers, along with a search bar and a user greeting for Jonathan Shaffer. A left sidebar contains a 'GeneGlobe' menu with options like 'Browse for Targeted Sequencing' and 'Browse for qPCR Arrays/Assays'. The main content area features a 'Data Analysis Center' header with a large image of a scientist and a line graph. Below the image is a text block describing the center's capabilities for analyzing real-time PCR and NGS data. A central navigation bar includes 'Analysis', 'Demo', and 'Resources'. Under 'Analysis', there is a 'Choose format' section with buttons for 'NGS', 'qPCR', and 'Microbial'. Below this are several dropdown menus for 'Choose Array / Panel', 'Choose Instrument / Pack Size', 'Choose Data Analysis Type', and 'Specify CatNo'. On the right side, there is a 'GeneGlobe Lists' section with counts for 'My Gene Lists', 'My miRNA Lists', 'My Pathway Lists', and 'My Publication Lists', and a 'Contact QIAGEN' section with links for 'Global contacts', 'Technical Service', and 'Customer Care'. At the bottom right, there is a section for 'QIAGEN Apps' with an image of a tablet and a link to 'See apps'.

Free, easy-to-use secondary data analysis: [qiagen.com/geneglobe](http://qiagen.com/geneglobe)

What can you do?

- Perform gene expression analysis

How can you normalize the data?

- geNorm
- Total Molecular Tag Count
- DESeq2
- Trimmed Mean of M (edgeR)

What is the output?


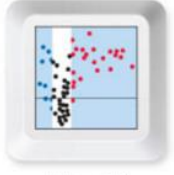

- Fold-regulation results
- Scatter plots, volcano plots and clustergrams
- Ingenuity Pathway Analysis (IPA) upload file

## Visual representations of data

▶ Upload data
▶ Analysis setup
▶ Analysis
▼ Plots & charts
▶ Export data

Plot Home
Scatter Plot
Volcano Plot
Clustergram

Select a Plot or Chart Icon below to launch a separate window.

<p><b>Notes:</b></p> <ol style="list-style-type: none"> <li>1. Before launching a plot or chart, please upload your readout data or take a test run.</li> <li>2. Please disable any Pop-Up Window Blockers in your web browser.</li> <li>3. To save the figure:             <ol style="list-style-type: none"> <li>a. Using Windows: Right-click on the figure and save the image/picture</li> <li>b. Using OS X: Hold down the Control key &amp; Click and save the image/picture</li> </ol> </li> </ol>	 <p><b>Scatter Plot</b></p>	<p>The scatter plot compares the normalized expression of every gene on the array between two groups by plotting them against one another to quickly visualize large gene expression changes. The central line indicates unchanged gene expression. Set the boundary (fold regulation cut-off) and the experimental groups to compare. Then, export the lists of genes whose expression changes are greater than the selected boundary.</p>
 <p><b>Volcano Plot</b></p>	 <p><b>Clustergram</b></p>	<p>The clustergram performs non-supervised hierarchical clustering of the entire dataset to display a heat map with dendrograms indicating co-regulated genes across groups or individual samples.</p>

\* NOTE: This plot requires three or more replicates in each group.



Scatter plot



Volcano plot



Clustergram



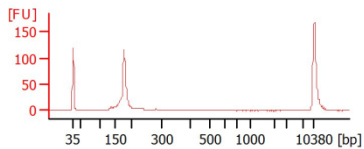
- 1 miRNA background
- 2 QIAseq miRNA sequencing: Product overview
- 3 Data analysis overview
- 4 Performance and application data**
- 5 Summary and questions

# QIAseq miRNA: Input as low as 1 ng without adapter dimers

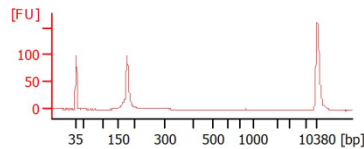
## QIAseq miRNA Library Kit workflow on kidney total RNA

- RNA amounts: 500 ng, 100 ng, 10 ng and 1 ng
- Sample QC: Bioanalyzer (BA)
- Sequencing: NextSeq®, 75 bp Single Read

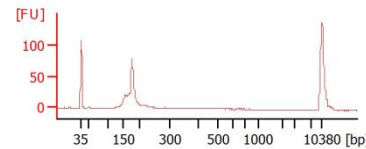
BA: 500 ng



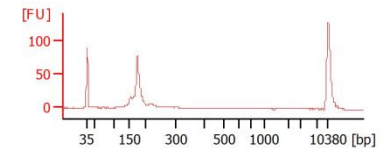
BA: 100 ng



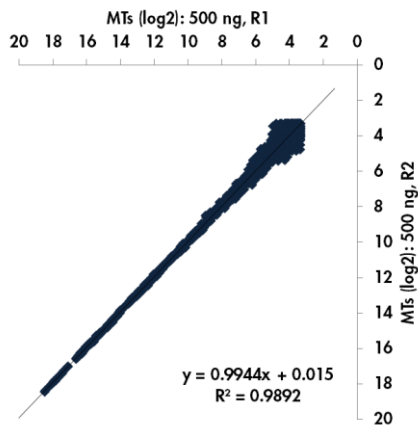
BA: 10 ng



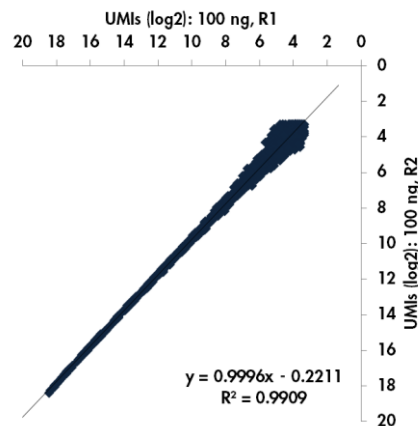
BA: 1 ng



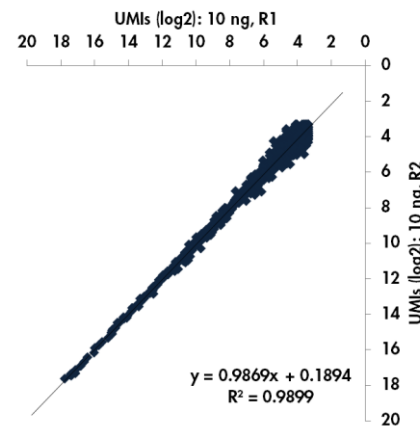
UMIs: 500 ng



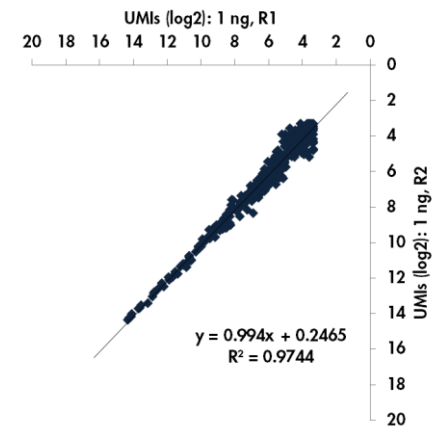
UMIs: 100 ng



UMIs: 10 ng



UMIs: 1 ng

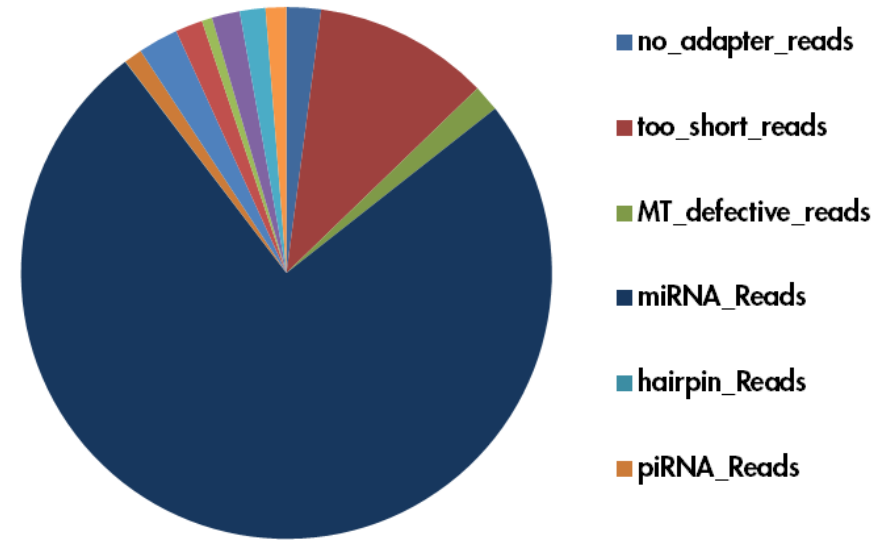


The QIAseq miRNA workflow enables robust, reproducible results from 1–500 ng.

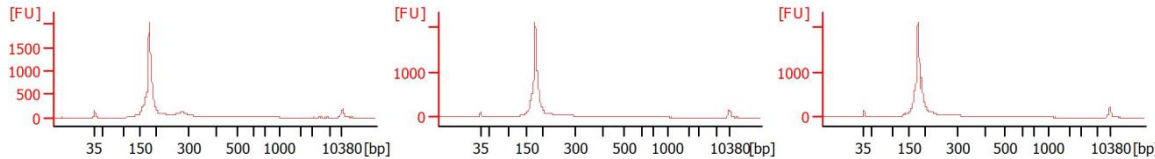


## QIAseq miRNA Library Kit workflow on kidney (K) and lung (L) total RNA

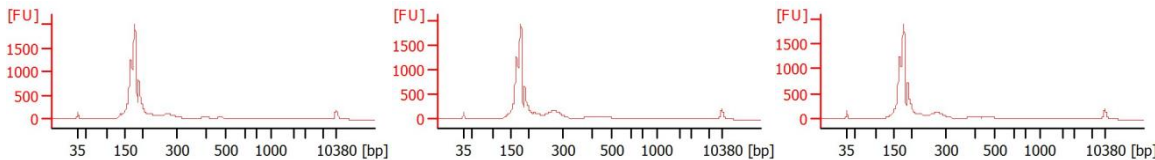
- RNA amount: 100 ng (n = 3 of each tissue)
- Sample QC: Bioanalyzer (BA)
- Sequencing: NextSeq, 75 bp Single Read



### BA: Kidney libraries



### BA: Lung libraries

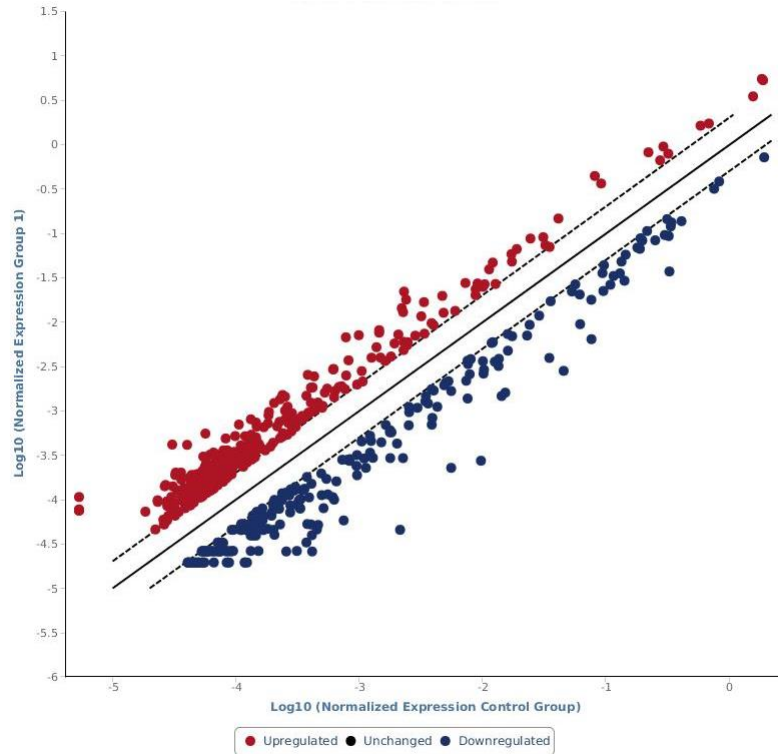


### Sequencing summary

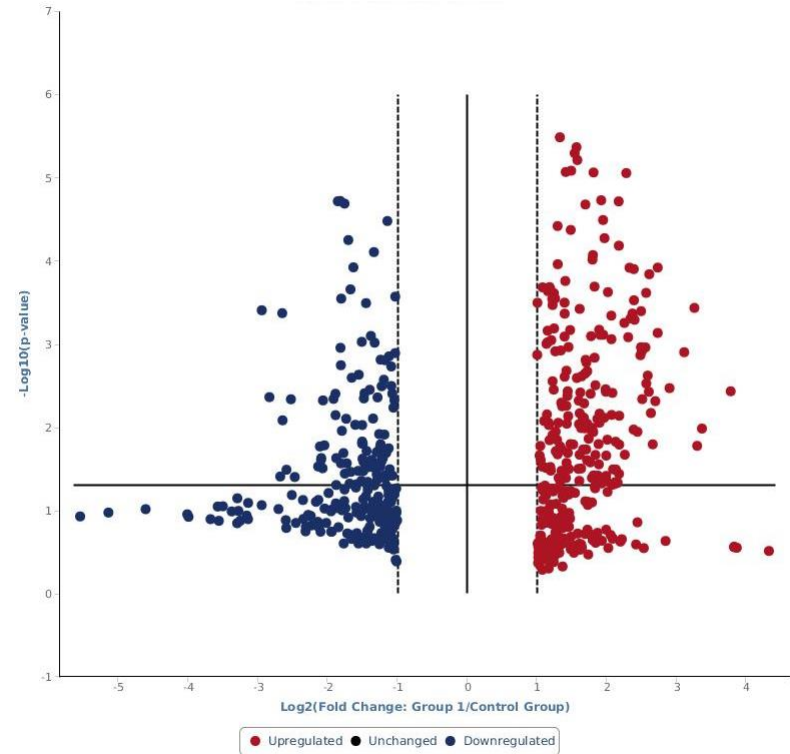
Sample	Mapped miRNA
K1	75%
K2	67%
K3	67%
L1	64%
L2	68%
L3	67%

The QIAseq miRNA workflow prepares robust libraries from tissues, enabling a high percentage of miRNA reads without gel excision.

Scatter plot (lung vs. kidney)



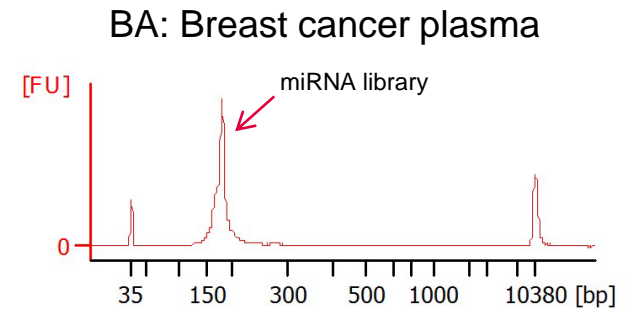
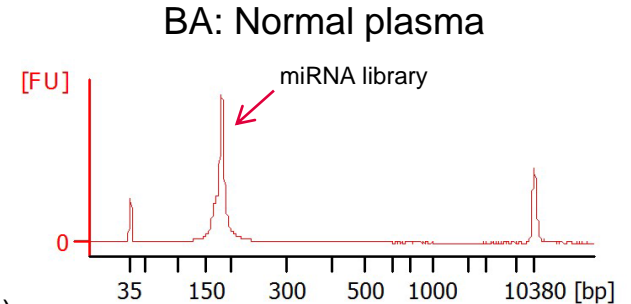
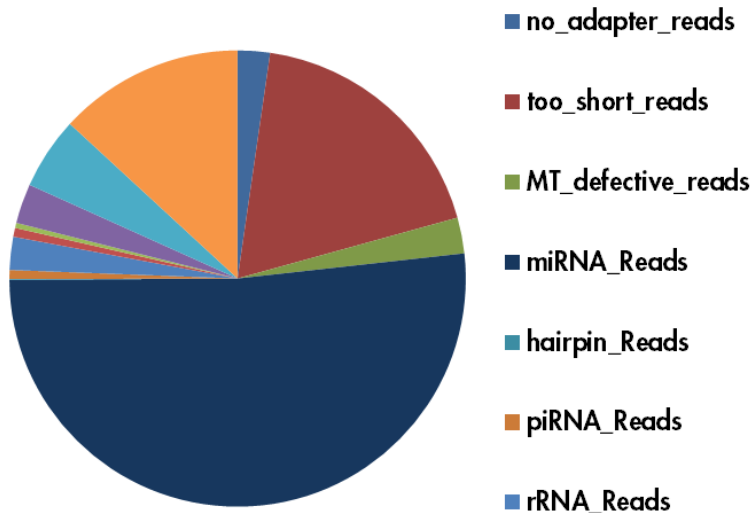
Volcano plot (fold-regulation vs. p-value)



Use the GeneGlobe Data Analysis Center to easily identify differentially expressed miRNAs.

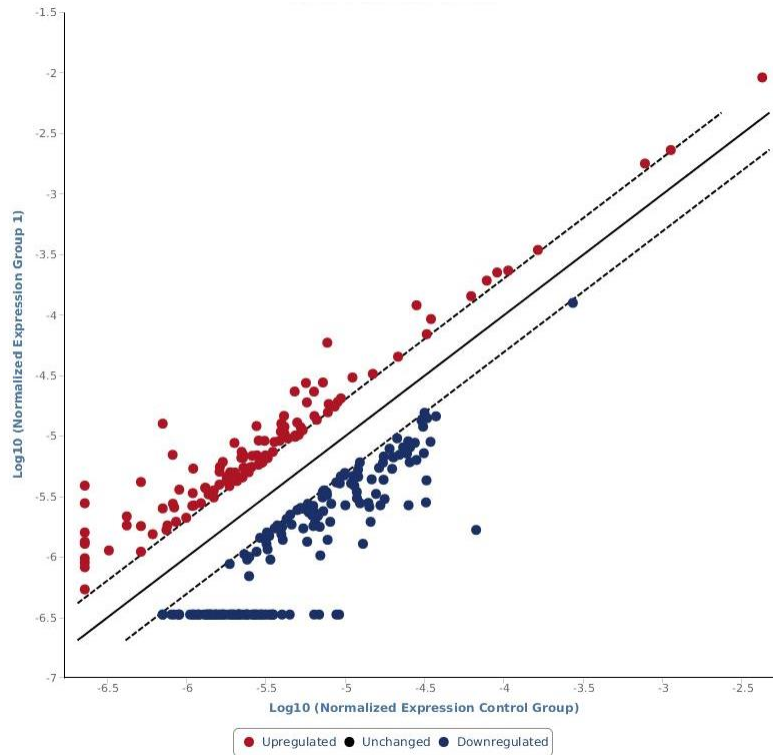
## QIAseq miRNA Library Kit workflow on plasma total RNA

- miRNeasy Serum/Plasma Kit: 200 µl input
- Total RNA input: 5 µl RNA eluate (80 µl of serum equivalents)
  - Normal (N) plasma (n = 3)
  - Breast cancer plasma (n = 3)
- Sequencing: NextSeq, 75 bp Single Read
- miRNA mapping %: 52 (N1), 54 (N2), 47 (N3), 49 (BC1), 41 (BC2), 57 (BC3)

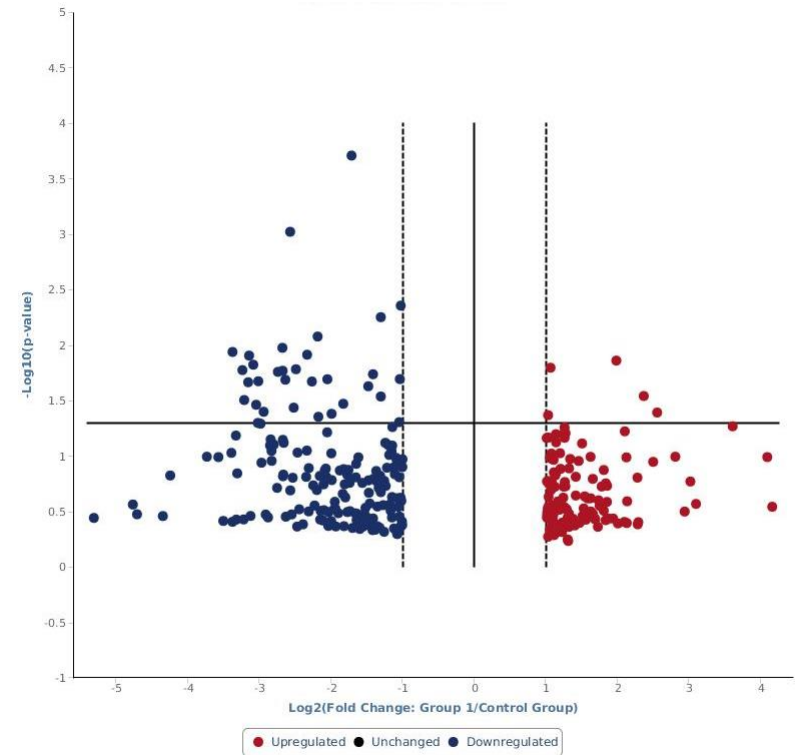


The QIAseq miRNA workflow prepares robust libraries from plasma, enabling a high percentage of miRNA reads without gel excision.

Scatter plot (breast cancer vs. normal)



Volcano plot (fold-regulation vs. p-value)

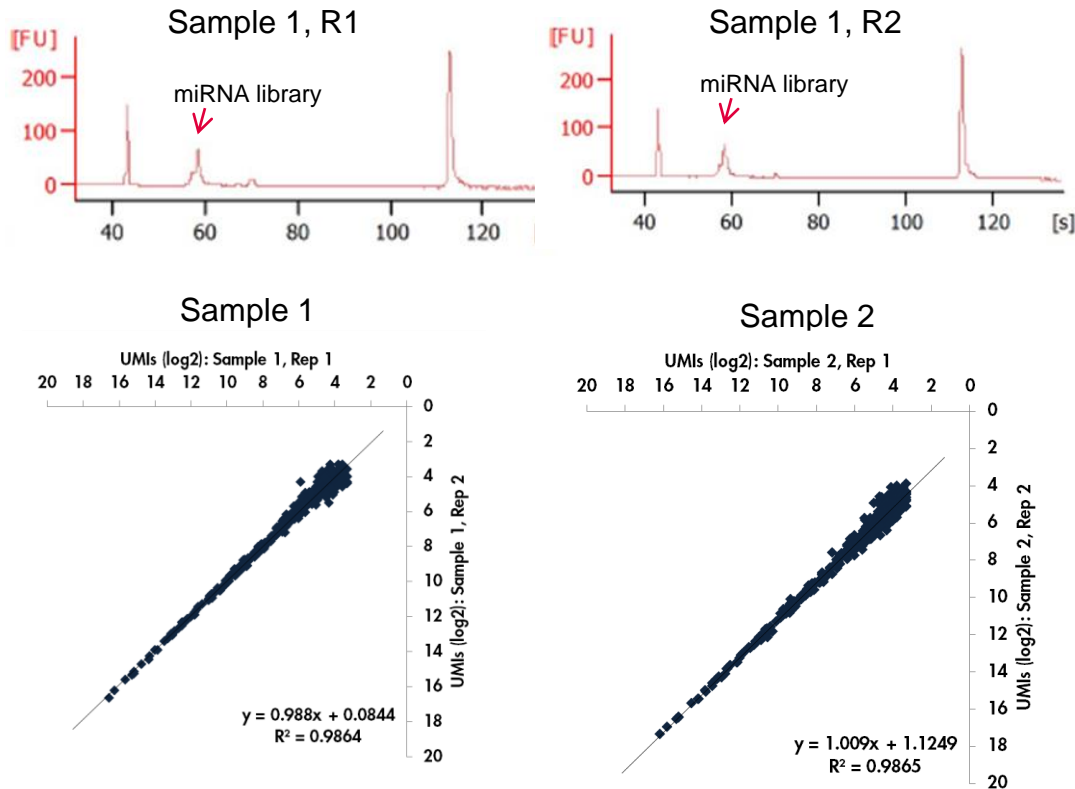


- 123 miRNAs upregulated (6 significantly)
  - miR-520g-3p: Associated with important prognostic factors in breast cancer patients
- 179 miRNAs downregulated (31 significantly)

● Use the GeneGlobe Data Analysis Center to easily identify differentially expressed miRNAs.

# Exosome samples: High mapped miRNA percentages

- Isolation: exoRNeasy (1 ml plasma processed)
- Samples: Four total RNA samples (2 donors, 2 replicates)
- RNA input: 5 µl of RNA eluate



read set	Sample 1, Rep 1	Sample 1, Rep 2	Sample 2, Rep 1	Sample 2, Rep 2
total_reads	3,454,577	3,539,076	2,531,228	6,230,468
no_adapter_reads	321,093	276,815	345,974	562,212
too_short_reads	737,228	799,712	461,630	1,240,571
UMI_defective_reads	192,311	154,736	194,358	398,158
miRNA_Reads	1,333,379	1,424,014	913,946	2,413,667
hairpin_Reads	2,787	2,851	2,078	6,554
piRNA_Reads	29,049	30,768	23,773	62,763
rRNA_Reads	93,880	92,517	83,114	201,736
tRNA_Reads	18,248	18,566	16,370	41,872
mRNA_Reads	12,127	12,383	9,533	24,309
otherRNA_Reads	149,227	152,885	88,306	239,613
notCharacterized_Mappable	135,715	139,635	120,375	326,720
notCharacterized_notMappable	429,533	434,194	271,771	712,293
miRNA Mapping %	38.6	40.2	36.1	38.7

Outcome: High mapping percentage to miRNAs; low mapping percentage to OtherRNA reads (often observed with other commercial kits).



- 1 miRNA background
- 2 QIAseq miRNA sequencing: Product overview
- 3 Data analysis overview
- 4 Performance & application data
- 5 Summary and questions**

## QIAseq miRNA

miRNA-focused next-generation sequencing library prep kit and integrated bioinformatics/data analysis solution

### What can be done with the sequencing data?

- Differential expression calculations of miRNA from highly multiplexed samples
- Novel miRNA discovery

### What are distinguishing features of the QIAseq miRNA Kit?

- Gel-free, rapid workflow
- Broad RNA input: 1–500 ng
- Library prep from serum, plasma, biofluids, cells and tissues
- Integrated Unique Molecular Index (UMI) technology
- Highly optimized chemistry
- All-in-one-box solution



Questions?

Jonathan Shaffer, Ph.D.

[Jonathan.Shaffer@qiagen.com](mailto:Jonathan.Shaffer@qiagen.com)