



Complete MicroRNA Solutions

• Target Validation, Overexpression, Inhibition, Quantitation & Profiling •

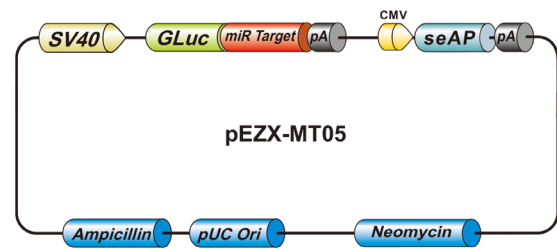
Functional Study Tools

miTarget™ miRNA 3' UTR Target Clones

Designed for microRNA (miRNA) target identification and functional validation of predicted targets, or to study the regulatory effect of miRNA on target genes.

Advantages:

- ✓ No lysis of the cells is necessary
- ✓ Dual-reporter vector system
- ✓ High-throughput compatible
- ✓ High sensitivity
- ✓ Convenience



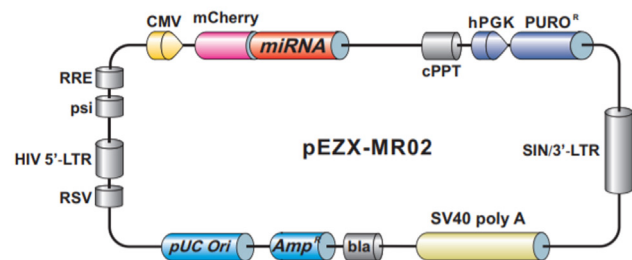
Vector	Reporter Gene	Tracking Gene	Compatible Assay Kit	Feature
pEZX-MT05	<i>Gaussia</i> luciferase	Secreted alkaline phosphatase (SEAP)	Secrete-Pair™ Dual Luminescence Assay Kit	No lysis of the cells is necessary
pEZX-MT06	Firefly luciferase	<i>Renilla</i> luciferase	Luc-Pair™ Luciferase Assay Kit	Cell lysis required

miExpress™ Precursor miRNA Expression Clones

Hairpin precursor miRNAs of approximately 150 nucleotides are cloned into lentiviral or non-viral vectors for delivery in virtually all cell types. Synthetic miRNA mimics are also available.

Advantages:

- ✓ Full coverage
- ✓ Flexible delivery systems
- ✓ Optimized expression cassettes
- ✓ Selection markers and reporters



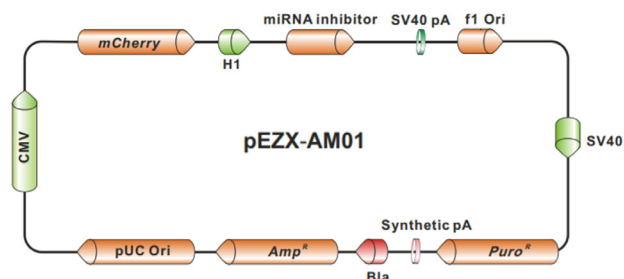
Vector	Promoter	Selection Marker	Reporter Gene	Type
pEZX-MR02	CMV	Puromycin	mCherry	HIV Lentiviral
pEZX-MR03	CMV	Puromycin	eGFP	HIV Lentiviral
pEZX-MR04	CMV	Puromycin	eGFP	Mammalian

miArrest™ miRNA Inhibitor Expression Clones

GeneCopoeia offers miRNA inhibitors as vector-based clones or synthetic oligonucleotides. Vector-based expression clones are available in lentiviral and non-viral vectors.

Advantages :

- √ Full coverage
- √ Flexible delivery systems
- √ Advanced and reliable design
- √ Superior performance



Vector	Promoter	Selection Marker	Reporter Gene	Type
pEZX-AM01	H1	Puromycin	mCherry	Mammalian
pEZX-AM02	U6	Puromycin	mCherry	Mammalian
pEZX-AM03	H1	Hygromycin	mCherry	HIV Lentiviral
pEZX-AM04	U6	Hygromycin	mCherry	HIV Lentiviral
pEZX-AM04p	U6	Puromycin	mCherry	HIV Lentiviral

miRNA Clone Sets

GeneCopoeia also offers pre-made or custom-selected clone sets. Uniform, high-quality clone sets are ideal for large-scale high content and high-throughput screening studies in functional genomics, proteomics and systems biology. We also provide custom clone set services.

Advantages :

- √ Large collection
- √ Quick delivery
- √ High Quality

Product type	Promoter	Selection Marker	Reporter Gene	Species	Vector Type
miRNA Precursor clone sets	CMV	Puromycin	eGFP	<ul style="list-style-type: none"> ● Human ● Mouse ● Rat 	<ul style="list-style-type: none"> ● Lentiviral ● Non-viral
miRNA Inhibitor clone sets	H1	Hygromycin	mCherry	<ul style="list-style-type: none"> ● Human ● Mouse ● Rat 	<ul style="list-style-type: none"> ● Lentiviral

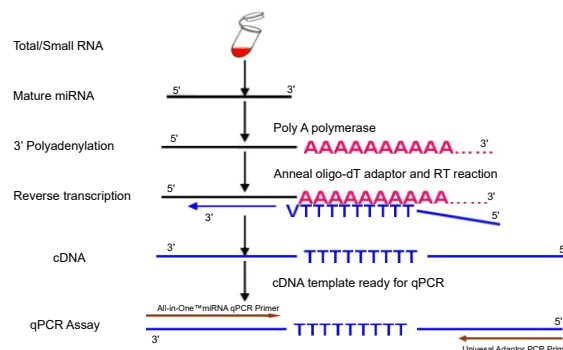
Profiling Tools

All-in-One™ qRT-PCR Reagent Kits and Validated Primers

The All-in-One™ miRNA qRT-PCR reagent kits include the miRNA qRT-PCR detection kits and the miRNA first-strand cDNA synthesis kits. When used with All-in-One™ Validated miRNA Primers or miProfile™ miRNA qPCR Arrays, the result is high-performance expression-profiling of all miRNAs.

Advantages :

- ✓ Detect miRNAs from as little as 10 pg of small RNA or 20 pg of total RNA
- ✓ Detect only mature miRNA
- ✓ Compatible with GeneCopoeia miRNA qPCR arrays or validated miRNA-specific primers



Overview of steps involved in the All-in-One™ miRNA qRT-PCR Detection Kit. During 3' polyadenylation, M-MLV RTase and a unique Oligo dT Adaptor primer, reverse transcribes the poly A miRNAs.

Catalog #	Product	Description
QP015	All-in-One™ miRNA qRT-PCR Detection Kit (20 RT and 200 qPCR reactions)	Poly A Polymerase, RTase Mix, qPCR Mix, ROX Reference Dye, Universal Adaptor PCR Primer and other buffers (for use with miRNA qPCR primers)
QP016	All-in-One™ miRNA qRT-PCR Detection Kit (60 RT and 600 qPCR reactions)	Poly A Polymerase, RTase Mix, qPCR Mix, ROX Reference Dye, Universal Adaptor PCR Primer and other buffers (for use with miRNA qPCR primers)
QP017	All-in-One™ miRNA First-Strand cDNA Synthesis Kit for miRNA qPCR array (20 RT reactions)	Poly A Polymerase, RTase Mix, PAP/RT buffer, spike-in control (for use with miRNA qPCR arrays)
QP018	All-in-One™ miRNA First-Strand cDNA Synthesis Kit for miRNA qPCR array (60 RT reactions)	Poly A Polymerase, RTase Mix, PAP/RT buffer, spike-in control (for use with miRNA qPCR arrays)
QP029	All-in-One™ miRNA Universal Adaptor PCR Primer (50 µM ; 20 µl)	miRNA Universal Adaptor PCR Primer (use with All-in-One™ miRNA First-Strand cDNA Synthesis kit QP017 or QP018)
Variable	All-in-One™ miRNA qPCR Primer (200 reactions)	Validated miRNA human, mouse and rat primers

miProfile™ miRNA qPCR Arrays

The miProfile™ miRNA PCR Arrays are designed for profiling the expression of pre-defined or customized sets of miRNAs in various tissues or cells. The resulting differential expression of profiled miRNAs help researchers to identify those miRNAs that are of biological significance and importance relevant to their research. GeneCopoeia can also make custom miRNA qPCR arrays of the miRNAs of your interest.

Advantages:

- √ Validated miRNA primers
- √ Robust performance
- √ Genome-wide coverage, pre-arranged groups, or customized groups

Product	Description	Species
miRNome qPCR Arrays	For profiling of miRNA whole genome expression	Human/Mouse
Cancer miRNA qPCR Arrays	For profiling of different types of cancer-related miRNA expression	Human/Mouse
Disease and Focus-Group miRNA qPCR Arrays	For profiling of disease or biological process-related miRNA expression	Human/Mouse
Exosome miRNA qPCR Arrays	For focused group profiling of exosome-related miRNA expression	Human
Custom miRNA qPCR Arrays	We will help you build custom qPCR arrays for your specific instrument type	Human/Mouse

Contact us:

GeneCopoeia, Inc.

9620 Medical Center Dr. Suite 101,
Rockville, MD 20850 USA

Email inquiry@genecopoeia.com
Tel +1 (301) 762-0888
Toll free +1 (866) 360-9531
Fax +1 (301) 762-3888
Website www.genecopoeia.com