

GeneCopoeia Introduces Genome-Wide MicroRNA (miRNA) Functional Analysis Tools

GERMANTOWN, Md., March 5 /PRNewswire/ -- GeneCopoeia announces the launch of a suite of miRNA functional analysis reagents and tools. They include 1) miExpress(TM) vector based precursor miRNA expression clones; 2) miTarget(TM) miRNA target sequence (3' UTR) expression clones; and 3) All-in-One(TM) miRNA RT-qPCR system.

Constructed in a lentiviral vector system, [miExpress\(TM\) precursor miRNA expression clone collections](#) cover all known miRNAs for human (750), mouse (450), and rat (270). These lentiviral vector based expression constructs allow efficient transduction of miRNAs into dividing and non-dividing cells transiently or stably, and can be used for *in vitro* assays and *in vivo* animal studies. The transduction efficiencies are monitored by the expression level of an indicator/reporter gene.

[miTarget\(TM\) miRNA target sequence \(3' UTR\) expression clones](#) cover genome-wide human, mouse and rat genes. These expression ready constructs are made in a mammalian expression vector system with dual reporter genes, which provides a ready-for-use and single vector solution for both transcriptional regulation reporting and an internal control for normalization reporting.

In addition to the two types of expression constructs, GeneCopoeia plans to launch All-in-One(TM) miRNA RT-qPCR system in the next few weeks. This system consists of a reverse transcription kit, a SYBR green RT-qPCR master mix kit, and a complete set of validated primers for human, mouse, and rat miRNAs, which allows specific and sensitive quantification of miRNAs individually or in an array format.

GeneCopoeia also plans to release soon a set of vector based miRNA suppressor construct products currently in development. Together with other miRNA products aforementioned, this product line will provide a complete set of tools and solutions for miRNA functional analysis.

For more information, please visit: <http://www.genecopoeia.com/product/mirna/>

About GeneCopoeia

[GeneCopoeia, Inc](#) is a Maryland, US, based manufacturer and provider of a wide range of biological and biomedical research reagent products as well as a full spectrum of custom services. GeneCopoeia, together with its joint venture, [FulenGen](#), has produced the largest number of full-sequence-verified human and mouse [full length protein coding ORF cDNA clones](#), which are offered to customers in more than 80 vectors with a choice of over 50 different fusion tags.

Sun Lu, Ph. D.
19520 Amaranth Drive
Germantown, Maryland 20874
USA
Tel: 301-515-6982
Fax: 301-515-6983
marketing@genecopoeia.com
www.genecopoeia.com