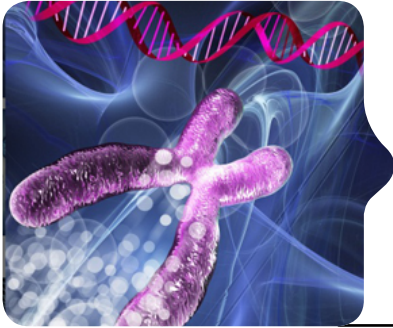


CRISPR Genome Editing Tools

Quick delivery • Sequence guaranteed • Complete solutions



Genome-CRISP™ CRISPR sgRNA Cloning Services

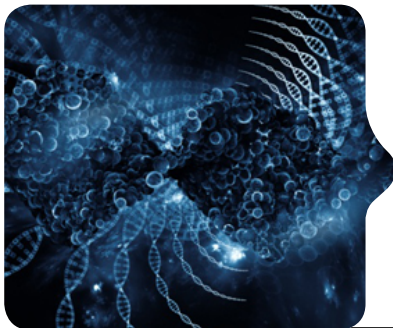
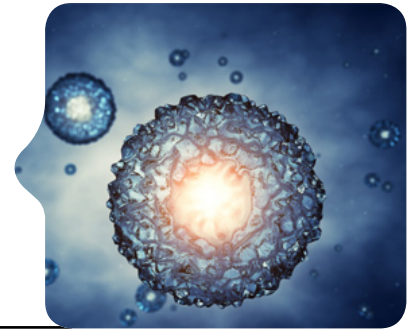
For precision RNA-guided CRISPR genome editing

We provide genome-wide, sequence-validated sgRNA clones for human, mouse and rat. Available in non-viral and lentiviral vectors. Lentiviral and AAV particles also available.

Genome-CRISP™ Premade Cas9 Stable Cell Lines

For sgRNA library screening and functional genomics

Our unique collection of Cas9-expressing stable cell lines comprises more than 70 human, mouse and rat background cell lines. A Cas9 integration kit is available to generate your own stable cell line.



Genome-CRISP™ CRISPR Human sgRNA Libraries

For high-throughput knockout of human genes in specific or custom gene groups or pathways

Our sgRNA libraries are available in lentiviral particles ready for transduction. Pathway-specific sgRNA libraries available for pathway screening in combination with Cas9 stable cell lines.

Safe Harbor Gene Knock-in Kits and Clones

For the expression of transgenes in cells and animals

Integration of a gene of interest in well-characterized safe harbor sites via CRISPR-mediated recombination. Pre-made Cas9 donor clones available to express Cas9 nuclease from the safe harbor sites.



CRISPR Genome Editing Tools

Additional Products and Services

Product/Service	Feature	Description
IndelCheck™ insertion/deletion detection system	Pre-validation of sgRNA and screening for genome modification	T7 endonuclease I assay with optimized conditions and positive control.
CRISPR Knock-In Donor construction services	HDR donor cloning vectors	Standard or customized donor vectors with multiple features for DIY construction.
	Custom HDR donor construction	Designed and constructed for CRISPR HR-mediated applications, such as knockout or knock-in.
Stable cell line services	Customized CRISPR-modified monoclonal stable cell lines.	
Transgenic mouse services	Mouse model generation	Customized CRISPR-modified transgenic mice.
CRISPR lentiviral packaging services	Cas9 lentivirus and customized lentiviral sgRNA production	Cas9 and sgRNA lentiviral particles for difficult-to-transfect cell lines.
Chromosome FISH Probes	FISH probes for chromosomal copy number determination	Chromosome enumeration probes for determining target chromosome copy number in CRISPR-mediated genome editing projects.

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