

Genome Editing: Applications For GeneCopia CRISPR sgRNA Libraries

Ed Davis, Ph.D.

Biomedical researchers are enjoying a Renaissance in functional genomics, which aims to use a wealth of DNA sequence information—most notably, the complete sequence of the human genome—to determine the natural roles of the genes encoded by the genome. As a result, biochemical networks and pathways will be better understood, with the hope of leading to improved disease treatments.

A major approach of functional genomics is to ablate gene function, by either “knockdown” (reduction) or “knockout” (complete elimination). Since 2012, researchers have turned increasingly to CRISPR (clustered, regularly interspaced, short palindromic repeats) for functional genomics studies. CRISPR’s simple RNA-guided mechanism provides a quick, convenient, and relatively low-cost method for many applications, from gene knockout, in-frame fusion tagging, mutagenesis, and transgene knockin. Several groups recently adapted CRISPR for high-throughput knockout applications, by developing large-scale CRISPR sgRNA libraries. GeneCopia recently launched a number of smaller, pathway- and gene group-focused CRISPR sgRNA libraries, which offer several key advantages over the whole-genome libraries. In this Technical Note, we discuss the merits and applications for CRISPR sgRNA libraries, how to use CRISPR sgRNA libraries, the advantages of using small, pathway- and gene group-focused libraries, and how GeneCopia can help you with your high-throughput CRISPR knockout studies.

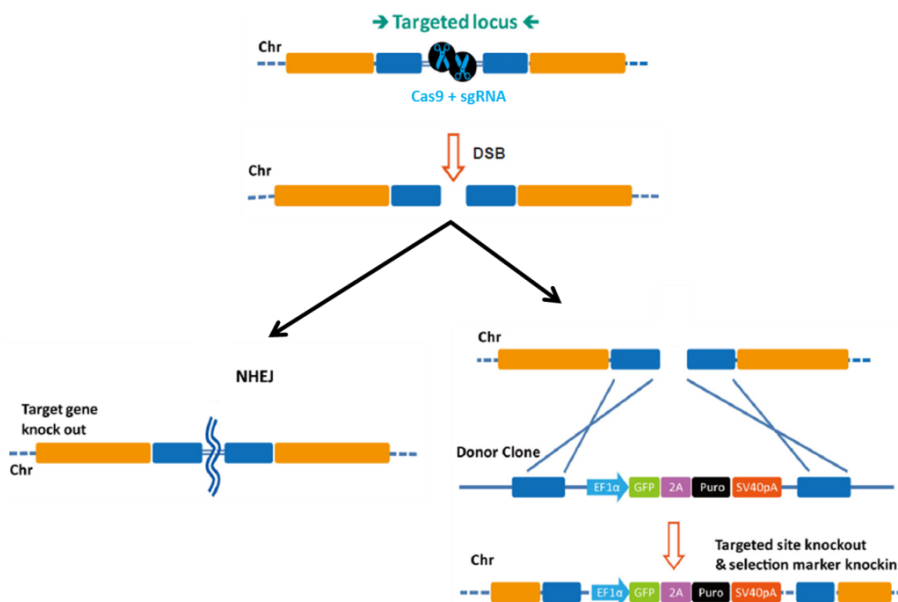


Figure 1. Pathways for repair of DSBs induced by genome CRISPR. Left: NHEJ. Right: HR in the presence of a donor construct.

What are CRISPR sgRNA libraries?

CRISPR sgRNA libraries are collections of hundreds-to-thousands of plasmids, each expressing a unique chromosome target-specific single guide RNA (sgRNA), which, in the presence of *Streptococcus pyogenes* Cas9 nuclease causes double-strand breaks (DSBs) in the chromosome (Figure 1). DSBs are repaired

either by homologous recombination (HR) in the presence of a homologous donor construct, or nonhomologous end joining (NHEJ), which leads to knockout mutations

CRISPR sgRNA libraries enable researchers to knock out many genes simultaneously in mammalian cells, opening the door to many applications, such as drug target identification, drug target validation, phenotypic changes, and reporter assays. Previously, these applications were achieved using RNA interference (RNAi). For example, Sethi, et al., (2012) used an shRNA library directed against 6,000 “druggable” genes to identify potential targets in ovarian cancer cells. Cooper and Brockdorff (2013) used an shRNA library to identify factors that activated a Gata6-neo promoter reporter. Finally, Zhang, et al, (2014) used fluorescence activated cell sorting (FACS) to isolate cells expressing the cell surface marker PSA-NCAM, which is expressed in human embryonic stem cells (hESCs).

While shRNA libraries are useful for high-throughput loss-of-function screens, RNAi has a number of disadvantages compared with CRISPR: 1) RNAi causes a knockdown of the level of gene expression. Therefore, false negatives, resulting from residual gene expression, can be missed; 2) RNAi acts only on cytoplasmic RNA, and so cannot silence nuclear RNAs like long non-coding RNAs. Conversely, CRISPR makes permanent changes to the genetic code. Thus, CRISPR can make a complete knockout of all alleles of a gene, whether their transcription products are localized to the nucleus or cytoplasm.

How are CRISPR sgRNA libraries used?

Using sgRNA libraries is straightforward. However, it is important to have a “readout”, (observable phenotype or assay) in order for the libraries to be useful (Figure 2). Briefly, sgRNA libraries are generally used as pools of lentiviral particles, although the plasmids can also transfect cells as DNA. Our library clones contain sgRNA target sites, but not Cas9, which is required. So you will need to either co-transfect /co-transduce your cells with sgRNAs and Cas9, or transfect/transduce a cell line already stably expressing Cas9. For lentiviral transduction, we recommend first establishing a cell line stably expressing Cas9; the large size of Cas9 limits the viral titer, so transduction is more efficient with the sgRNA-expressing lentiviruses alone. In any case, cells are transfected with DNA or infected with the lentiviral pools, followed by screening for the response of interest. Examples of some screening readouts include:

1. Cell death. Cells carrying lethal knockouts will be underrepresented in a pool of surviving cells.
2. Drug resistance: Cells resistant to a drug of interest will be overrepresented in the pool after applying drug treatment.
3. Reporter assay: Cells that lose the ability to express a fluorescent reporter (e.g. GFP) will be underrepresented in the population after FACS.

After the readout method or assay is applied, DNA sequencing is used to identify which sgRNAs are either under- or over-represented in the pools. For a very large, whole-genome scale library, it is necessary to use next-generation sequencing to identify these sgRNAs targeting candidate genes.

Development of CRISPR sgRNA libraries

In light of the advantages of CRISPR over RNAi for eliminating gene function, a few researchers have produced large CRISPR sgRNA libraries. In one study, Wang, et al. (2014) constructed pools of sgRNA-expressing lentivirus particles targeting 7,300 genes, which were used to infect a haploid cell line that was stably expressing Cas9 nuclease. In screening the transduced line for cells resistant to 6-thioguanosine, these researchers observed enrichment for cells carrying sgRNAs targeting genes in the mismatch repair pathway, but not in other DNA repair pathways. Similarly, Feng Zhang’s group at MIT

used a lentiviral sgRNA library targeting approximately 18,000 genes to identify melanoma and stem cells that were either sensitive or resistant to vemurafenib (Shalem, et al., 2014). The Zhang lab has since followed up on that initial study by producing second-generation libraries that contain several improvements and target a larger number of genes in human and mouse. (Sanjana, et al. 2014).

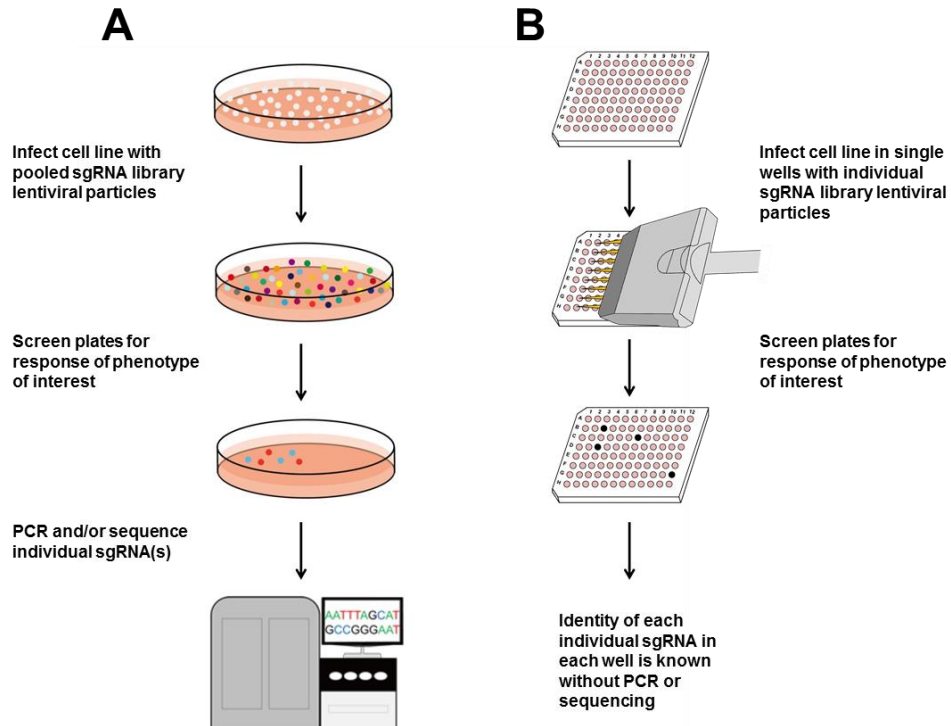


Figure 2. Basic workflow for using CRISPR sgRNA libraries. A. Pooled screen. Cells are infected with each sgRNA library pool. Infected pools are screened for the desired readout or phenotype. Pooled cells are subjected to Sanger sequencing for individual sgRNAs, or deep sequencing to look for over- or under-representation of individual sgRNAs. B. The GeneCopoeia Genome-CRISP™ CRISPR sgRNA libraries are available either as pools or individual sgRNAs. If desired, cells can be infected with lentiviruses expressing individual sgRNAs. Plates or wells are screened for the readout or phenotype of interest. Individual sgRNAs corresponding to the phenotype of interest are already known without sequencing.

GeneCopoeia Genome-CRISP™ CRISPR sgRNA libraries

Recently, GeneCopoeia launched a series of pre-made CRISPR sgRNA libraries, which are targeted to smaller gene groups and pathways (Table 1).

Catalog number	Library name	Number of sgRNAs	Number of genes
L01-LS03	Innate kinases & ubiquitin ligases	475	239
L02-LS03	Nuclear hormone receptors	236	118
L03-LS03	Tumor metastasis genes	114	57
L04-LS03	Oncogenes	576	288
L05-LS03	Tumor suppressor genes	462	231
L06-LS03	Protein kinases	1,316	658
L07-LS03	Key genes in 50 pathways	278	139

Table 1. GeneCopoeia Genome-CRISPR™ pre-made human CRISPR sgRNA libraries. Each library can be ordered as pools of either bacterial stock, transfection-ready DNA, or lentiviral particles from our website.

Each sgRNA is individually constructed, sequence-verified, and cultured in *E. coli* before pooling, in contrast to other CRISPR sgRNA libraries, which are constructed from the beginning as pools. GeneCopoeia's CRISPR sgRNA libraries have several advantages over other sgRNA libraries, such as:

1. Individual construction of each sgRNA makes it possible to purchase the libraries either as pools or as individual clones.
2. Sequence verification provides high quality of each sgRNA.
3. Individual construction and culturing of each sgRNA ensures the best representation in pools.
4. Small library sizes reduce the amount of work and time spent on analyzing your results. You can screen the library for candidates using either Sanger or next-generation sequencing.

Our library sgRNAs are cloned into a very versatile vector backbone (Figure 3), which can be used for either direct transfection of cells, packaging into lentiviral particles, or *in vitro* transcription.

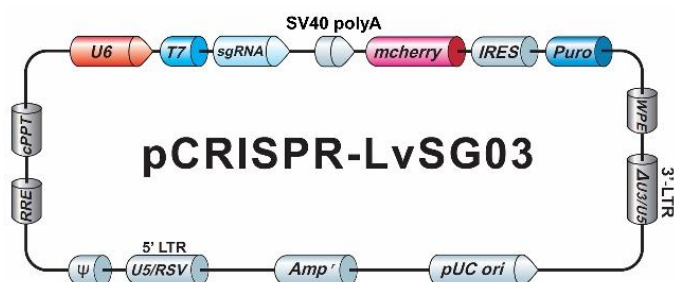


Figure 3. Plasmid backbone carrying individual sgRNAs in the GeneCopoeia Genome-CRISPR™ CRISPR sgRNA libraries.

Another significant advantage of the GeneCopoeia Genome-CRISPR™ CRISPR sgRNA libraries is that, because the sgRNA clones are individually constructed before pooling, researchers can order custom libraries in which we “mix and match” sgRNA clones from different libraries. In addition, we can also create custom CRISPR sgRNA libraries targeting new sets of genes.

Related products and services

In addition to our versatile, high-quality CRISPR sgRNA libraries, we also offer a number of companion products and services (Table 2). These include pre-made cell lines stably expressing Cas9 nuclease from the human AAVS1 Safe Harbor locus, and donor clones that you can use with our human AAVS1 Safe Harbor knockin kits to knock Cas9 in to AAVS1 in your cell line of choice. We also offer stable cell line services, in which we can construct a cell line carrying virtually any modification of interest, including expression of Cas9 from AAVS1 in a chosen cell line.

Conclusions

In this Technical Note, we highlighted a new and exciting technological development in the ever-evolving world of Genome Editing: CRISPR sgRNA libraries. As an alternative to whole genome libraries, GeneCopoeia provides several powerful, high-quality, and versatile CRISPR sgRNA libraries targeting smaller, focused pathways and gene groups. To find out more about how the GeneCopoeia Genome-

CRISPTM CRISPR sgRNA libraries can accelerate your research, visit our website: <http://www.genecopoeia.com>, or call us at 1-866-360-9531.

Catalog number	Product
Cas9 clones	
CP-C9LvNU-01	Cas9 nuclease lentiviral expression clone
CP-C9LvNU-02	Cas9 nuclease lentiviral expression clone
Cas9 lentiviral particles	
LPP-CP-LvC9NU-01-100	Cas9 Nuclease Purified Lentifact™ Lentiviral Particles (100 µl x 1 vial)
LPP-CP-LvC9NU-02-100	Cas9 Nuclease Purified Lentifact™ Lentiviral Particles (100 µl x 1 vial)
Pre-made Cas9-expressing stable cell lines	
SCL-01-CA1	Human cell line H1299 stably expressing CRISPR Cas9, single clone
SCL-02-CA2	Human cell line HEK293T stably expressing CRISPR Cas9, single clone
Cas9 AAVS1 knockin donor clones	
DC-C9NU-03	Cas9 AAVS1 knockin donor clone (Puro selection)
DC-C9NU-04	Cas9 AAVS1 knockin donor clone (Hygro selection)
DC-C9NU-05	Cas9 AAVS1 knockin donor clone (Neo selection)
Human AAVS1 Safe Harbor knockin kits	
SH-AVS-K100	Genome-TALER™ human AAVS1 safe harbor gene knock-in kit
SH-AVS-K000	Genome-TALER™ human AAVS1 safe harbor gene knockin kit (without donor)
SH-AVS-K200	Genome-CRISP™ human AAVS1 safe harbor gene knockin kit
SH-AVS-K002	Genome-CRISP™ human AAVS1 safe harbor gene knockin kit (without donor)

Custom stable cell line services available on request. [Contact us for a quote.](#)

Table 2. Companion products and services for the GeneCopoeia™ Genome-CRISP™ CRISPR sgRNA libraries.

References

- Cooper & Brockdorff (2013). Genome-wide shRNA screening to identify factors mediating Gata6 repression in mouse embryonic stem cells. *Development* **140**, 4110.
- Sethi, *et al.*, (2012). An RNA interference lethality screen of the human druggable genome to identify molecular vulnerabilities in epithelial ovarian cancer. *PLoS ONE* **7**.
- Shalem, *et al.* (2014). Genome-scale CRISPR-Cas9 knockout screening in human cells. *Science* **343**, 84.
- Sanjana, *et al.* (2014). Improved vectors and genome-wide libraries for CRISPR screening. *Nature Methods* **11**, 783.
- Wang, *et al.* (2014). Genetic screens in human cells using the CRISPR-Cas9 system. *Science* **343**, 80.
- Zhang, *et al.* (2014). Functional genomic screen of human stem cell differentiation reveals pathways involved in neurodevelopment and neurodegeneration. *Proc Natl Acad Sci U S A* **110**, 12361.