

Applications For CRISPR-Cas9 Stable Cell Lines

Presenter:

**Ed Davis, Ph.D.
Senior Application Scientist
GeneCopoeia, Inc.**

March 22, 2017

GeneCopoeia products & services

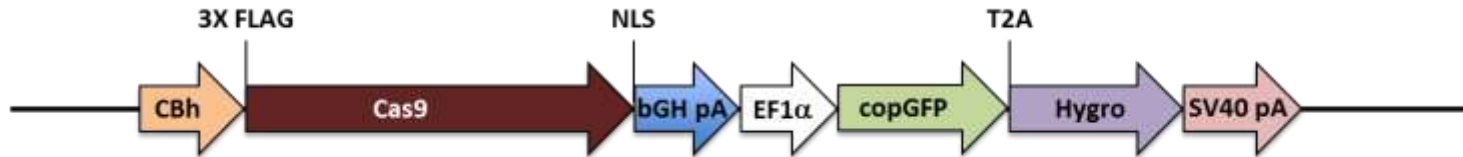
Functional Genomics & Cell Biology

<i>Clones</i>	<i>Viral systems</i>	<i>Kits & reagents</i>	<i>Fluorescent detection</i>
ORF	Lentivirus	Transfection	Cell function assays
Promoter	AAV	Luciferase	Nucleic acid detection
miRNA		FISH probes	Cell structure probes
CRISPR		Indel detection	Fluorescent dyes
siRNA		Cloning	



GeneCopoeia genome editing services

Cas9-expressing stable cell lines



- ❖ Cell lines with stably expressing Cas9
- ❖ Have pre-made lines, or can have us integrate Cas9 in your cell line
- ❖ Plasmids are available for DIY stable cell line creation
- ❖ Ideal for sgRNA library screening, validation, inducible CRISPR, and more

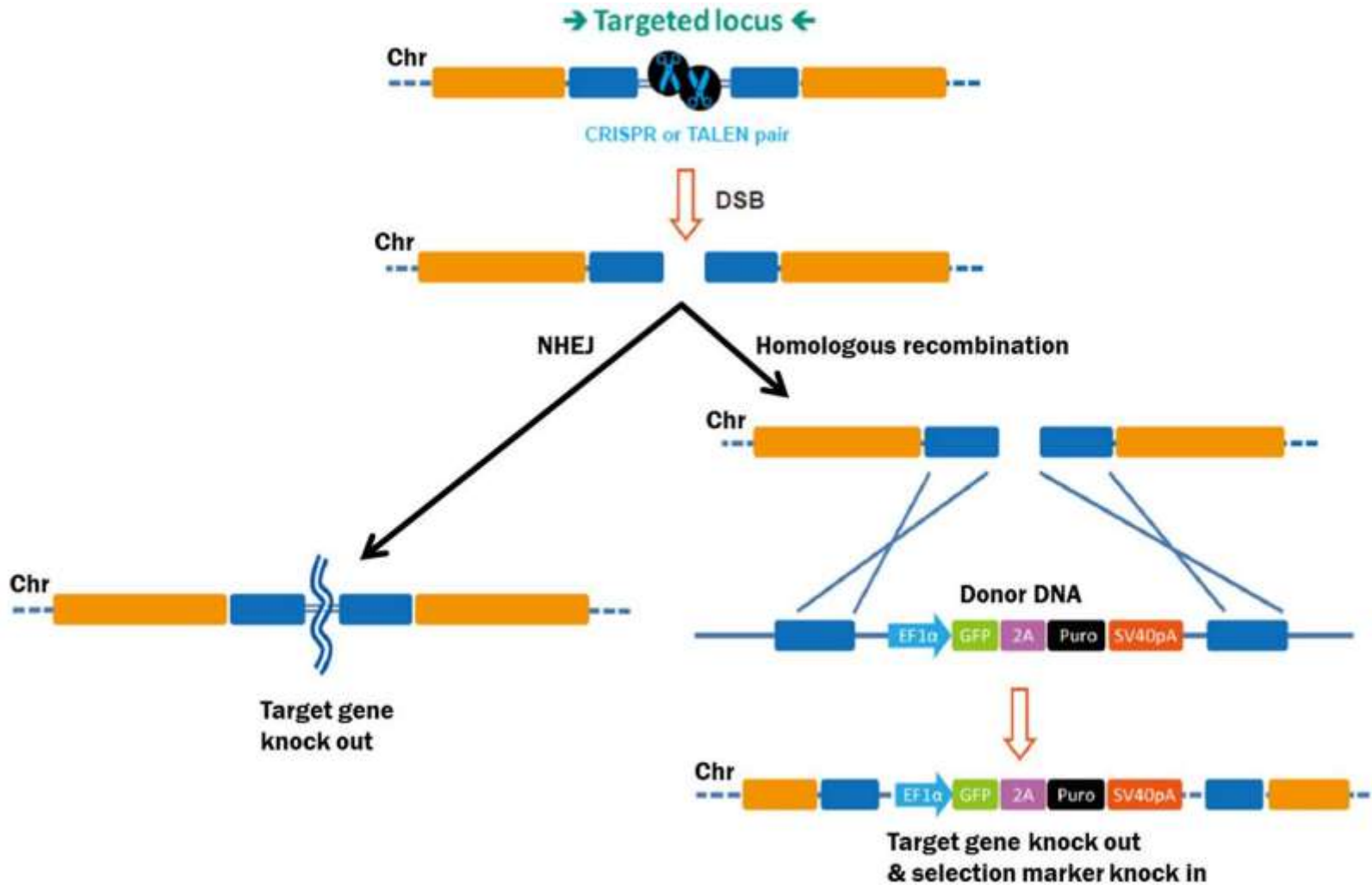
Outline

- ❖ Genome editing: Technologies
- ❖ Applications for genome editing
- ❖ sgRNA validation
- ❖ Lentiviral CRISPR
- ❖ CRISPR library screening
- ❖ Inducible CRISPR

Outline

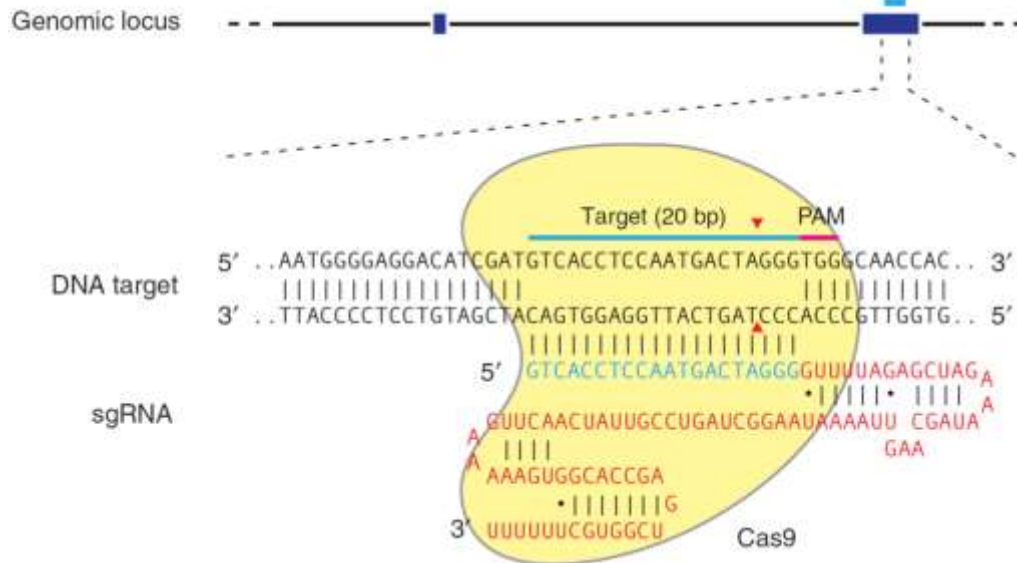
- ❖ **Genome editing: Technologies**
- ❖ Applications for genome editing
- ❖ sgRNA validation
- ❖ Lentiviral CRISPR
- ❖ CRISPR library screening
- ❖ Inducible CRISPR

Targeted DNA editing by DSB induction



CRISPR genome editing technology

CRISPR-Cas9: RNA-guided endonuclease

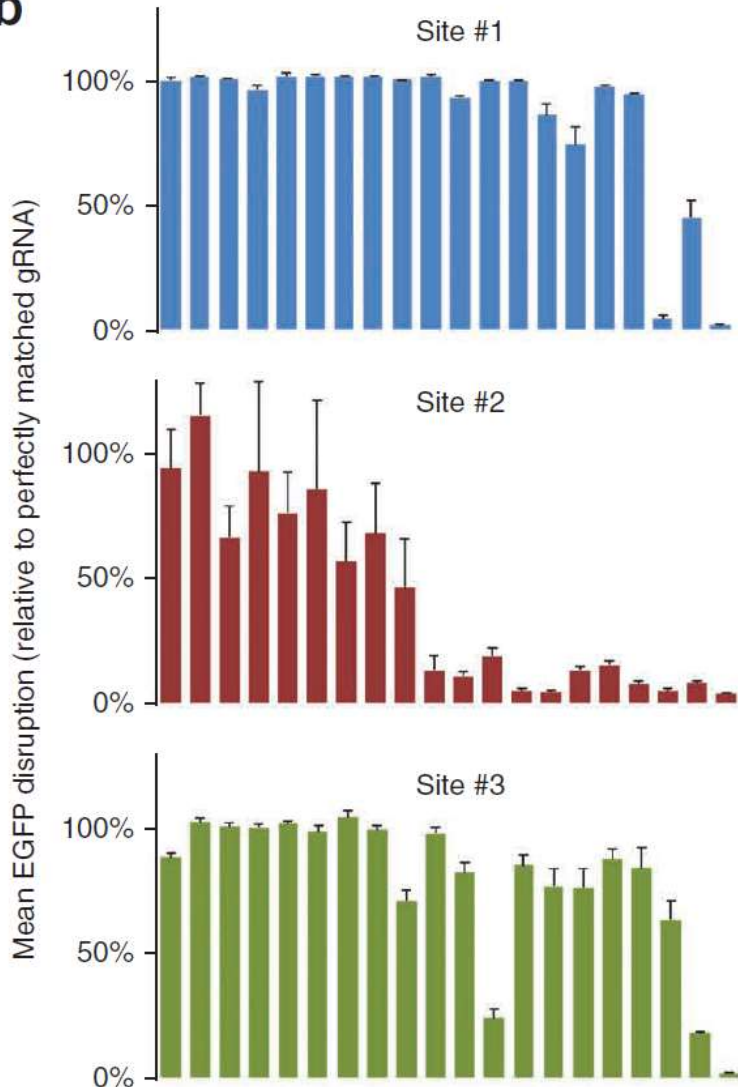


Ran, et al. (2013). Nature Protocols 8, 2281

- ❖ 20 nt single guide RNA (sgRNA) guides Cas9 nuclease to target site.
- ❖ Requires NGG “PAM” site immediately downstream of sgRNA target sequence.
- ❖ Cas9-RNA complex makes DSB 3-4 nt upstream of PAM.
- ❖ Target almost any gene in any cell

CRISPR-Cas9: Specificity

b



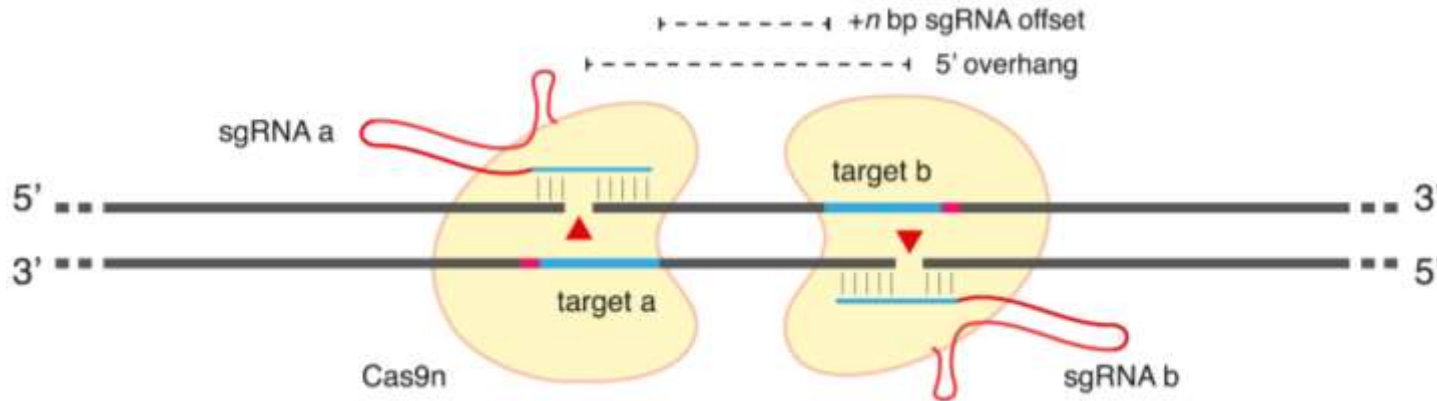
High-frequency off-target mutagenesis induced by CRISPR-Cas nucleases in human cells

Yanfeng Fu¹⁻⁴, Jennifer A Foden¹⁻³, Cyd Khayter¹⁻³, Morgan L Maeder^{1-3,5}, Deepak Reyon¹⁻⁴, J Keith Joung¹⁻⁵ & Jeffrey D Sander¹⁻⁴

- ❖ Showed that some sgRNAs with single, double, and even up to 5 transversion mismatches could still direct Cas9 to mutate EGFP.
- ❖ Found that for 4 of 6 tested sgRNAs, evidence of off-target mutagenesis (5.6% -125% of the intended targets).

CRISPR-Cas9: Specificity

Double Nicking by RNA-Guided CRISPR Cas9 for Enhanced Genome Editing Specificity

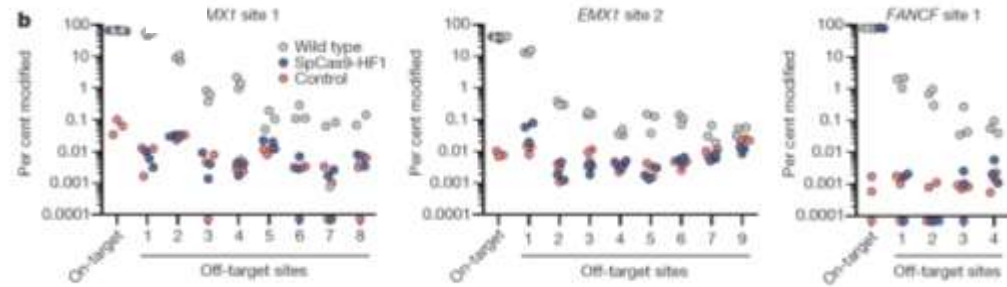
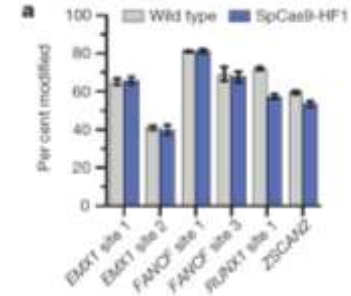
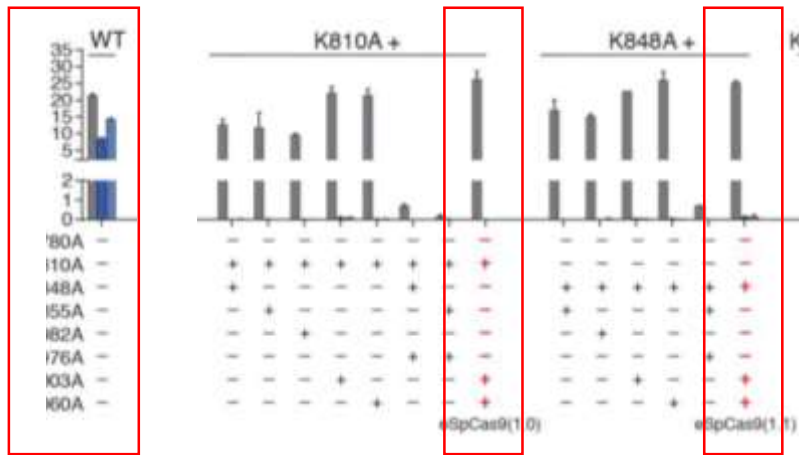


Ran, et al. (2013). Cell 154, 1380

- Cas9 D10A “nickase” mutant creates single-strand nicks instead of DSBs
- Off-target nicks repaired by high-fidelity base excision repair
- Permits ability to generate dimer-like chimeric endonuclease, similar to TALEN. 2 nicks will create a DSB
- Dramatically (50x-1,500x) reduces incidence of off-target effects

CRISPR-Cas9: Specificity

Cas9 point mutants



Slymaker, et al. (2015). Science

Kleinstiver, et al. (2016). Nature **529**, 490

Outline

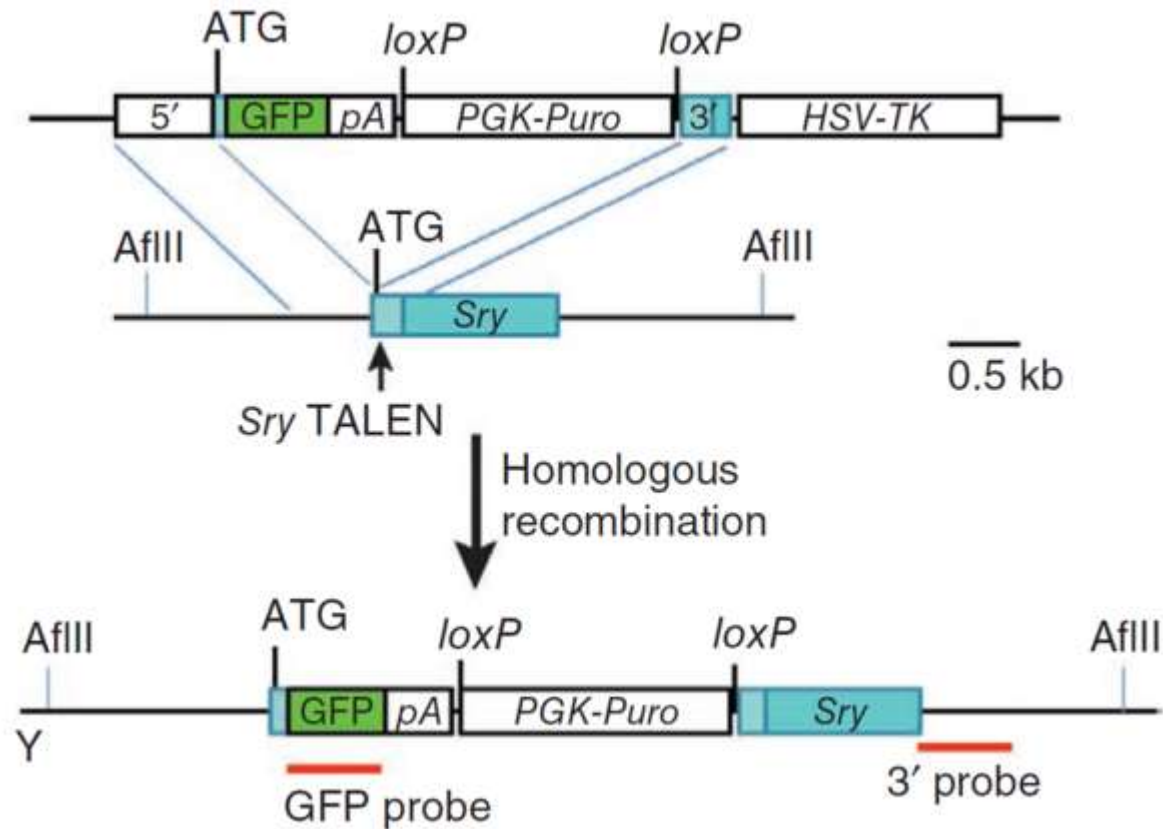
- ❖ Genome editing: Technologies
- ❖ Applications for genome editing
- ❖ sgRNA validation
- ❖ Lentiviral CRISPR
- ❖ CRISPR library screening
- ❖ Inducible CRISPR

Applications for genome editing

Application	Example	Technology	Reference
Gene knockout	Knockout SRY, UTY genes in cultured cells	TALEN via NHEJ and plasmid donor	Wang, et al. (2013). Nature Biotech. 31, 530
Gene knockout	Knockout of tet genes in transgenic rats	CRISPR via NHEJ	Li, et al. (2013). Nature Biotech. 8, 684
Correction of disease mutations	Cured heritable cataracts in transgenic mice	CRISPR via oligonucleotide donor	Wu, et al. (2013). Cell Stem Cell 13, 659
Engineered disease resistance	Knockout CCR5 gene to cure patients of HIV	ZFN via NHEJ	Perez, et al. (2008). Nat Biotech 26, 808
Forward mutagenesis screens	Identified genes in mismatch repair pathway by selecting for 6-thioguanosine resistance	CRISPR via NHEJ	Wang, et al. (2014). Science 343, 80
In-frame fusion tagging	C-terminal tag Sox2p with V5	CRISPR via oligonucleotide donor	Yang, et al. (2013). Cell 154, 1370
Transgene knock-in	Sox2, Oct4 ORFs KI into human AAVS1 "Safe Harbor"	TALEN via plasmid donor	GeneCopoeia internal data

Applications for genome editing

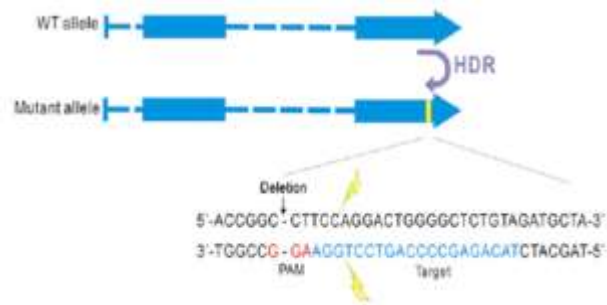
Knockout *via* HDR: Donor plasmid



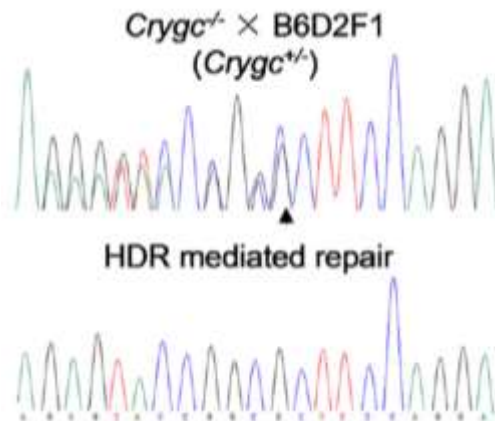
Wang, et al. (2013). Nature Biotech. 31, 530

Applications for genome editing

Mutagenesis *via* HDR: Oligo donor

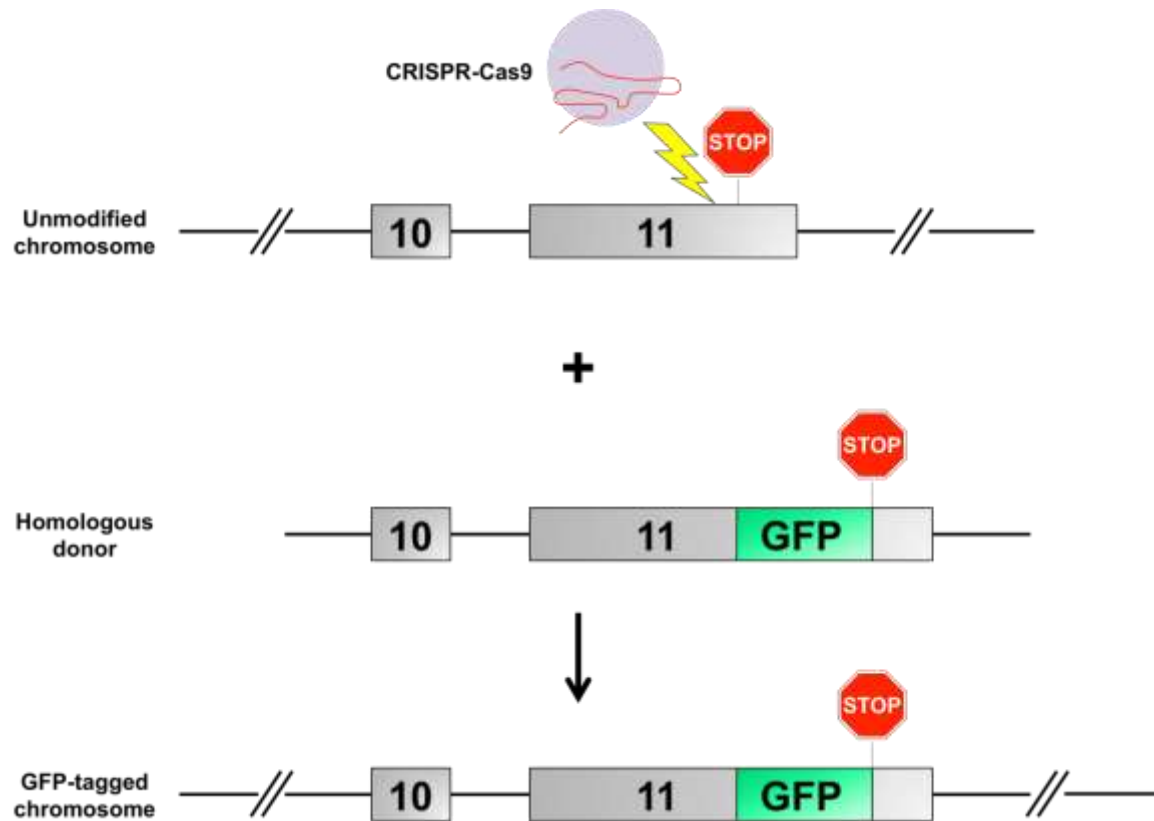


- ❖ Use single strand oligonucleotide (ssODN) to introduce base changes or small deletions.
- ❖ Use for mutagenesis or disease correction.
- ❖ Wu, et al.: Used CRISPR + ssODN to cure heritable cataract disease in mice



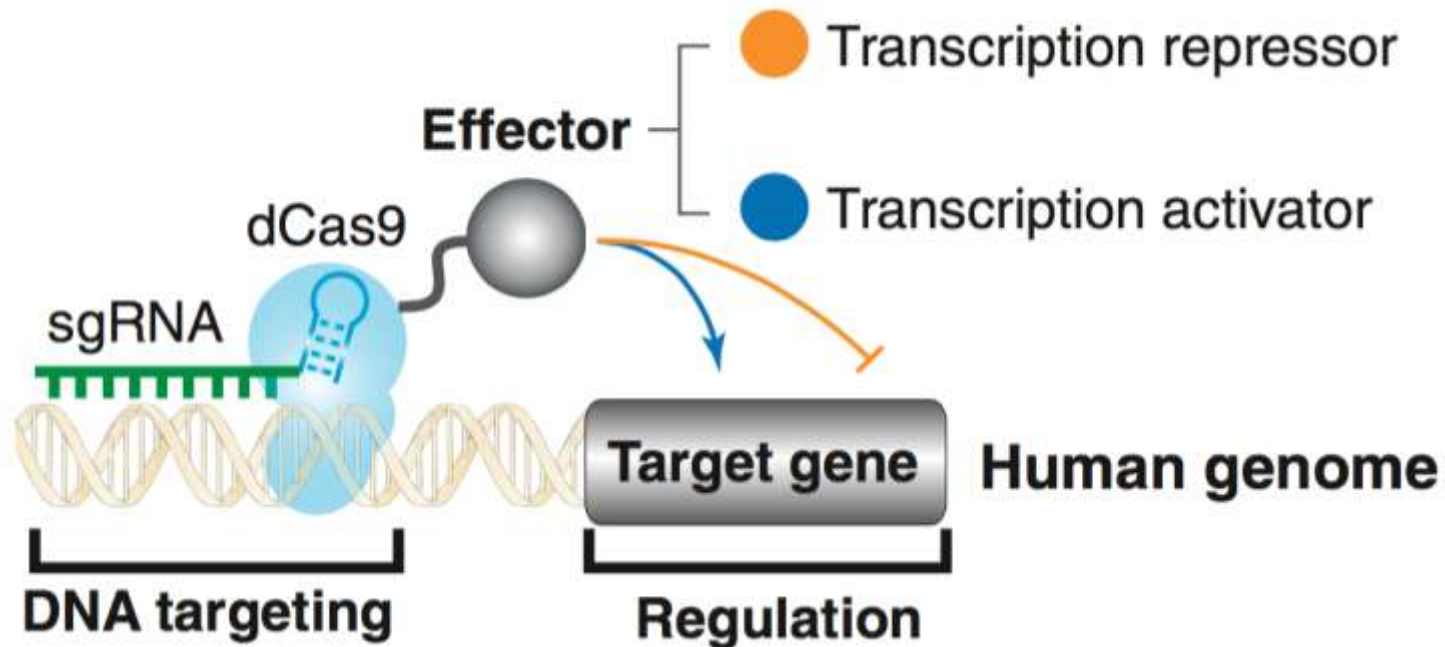
Applications for genome editing

In-frame fusion tagging



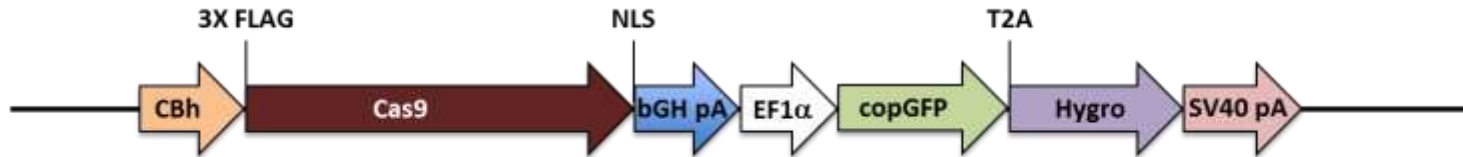
Applications for genome editing

Targeted gene activation/repression



GeneCopoeia genome editing services

Cas9-expressing stable cell lines



- ❖ Cell lines with stably expressing Cas9
- ❖ Have pre-made lines, or can have us integrate Cas9 in your cell line
- ❖ Plasmids are available for DIY stable cell line creation
- ❖ Ideal for sgRNA library screening, validation, inducible CRISPR, and more

Outline

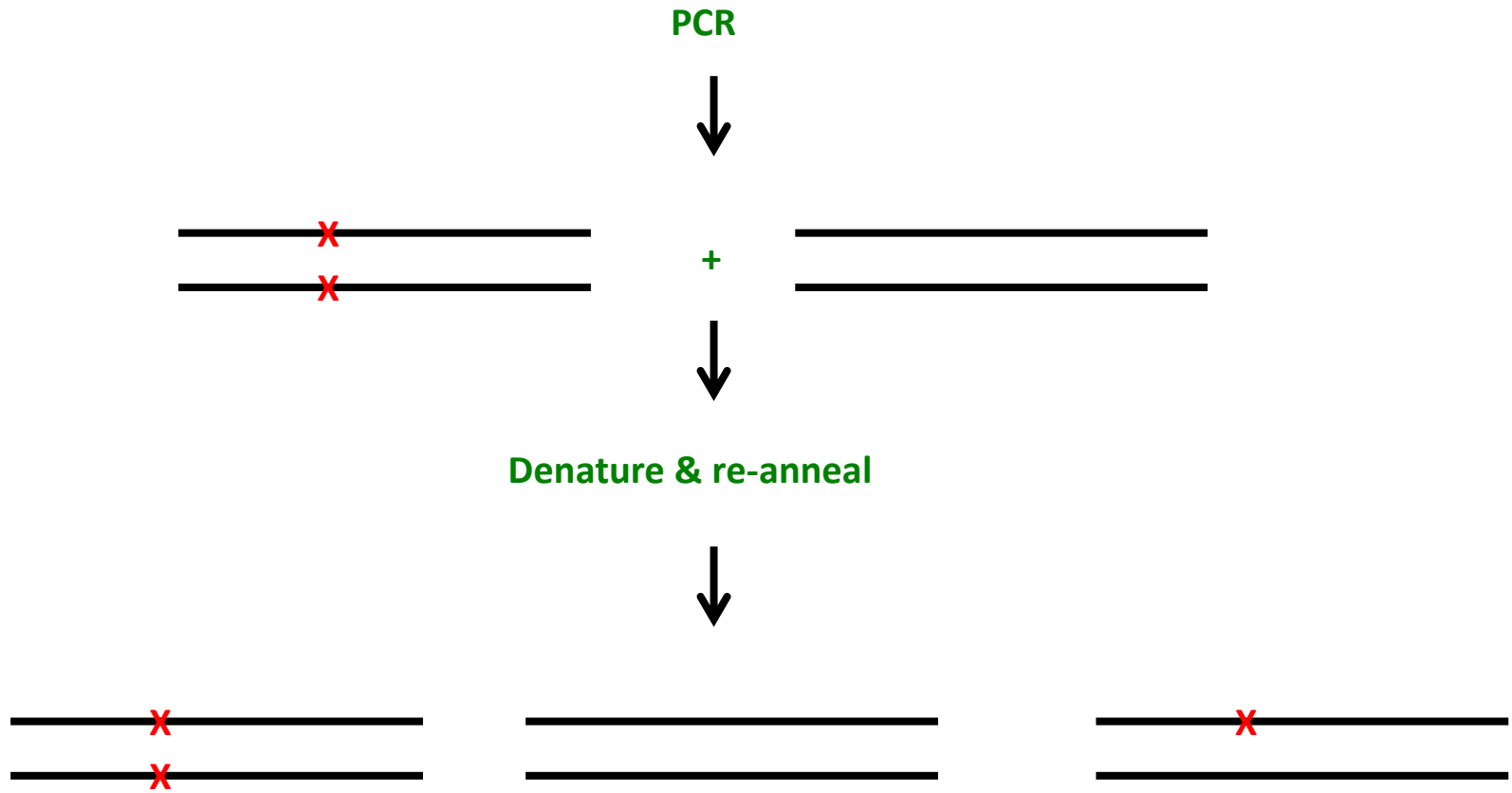
- ❖ Genome editing: Technologies
- ❖ Applications for genome editing
- ❖ **sgRNA validation**
- ❖ Lentiviral CRISPR
- ❖ CRISPR library screening
- ❖ Inducible CRISPR

Functional validation

Why do functional validation?

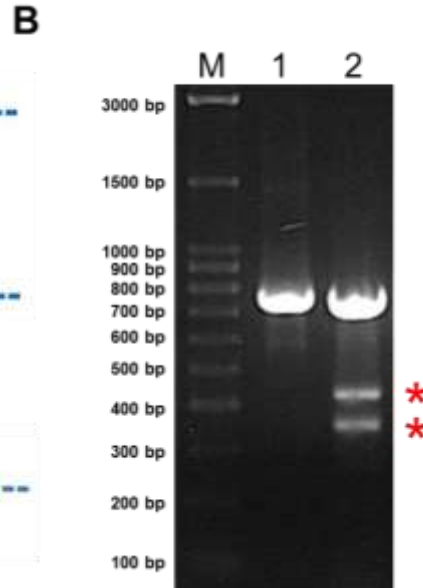
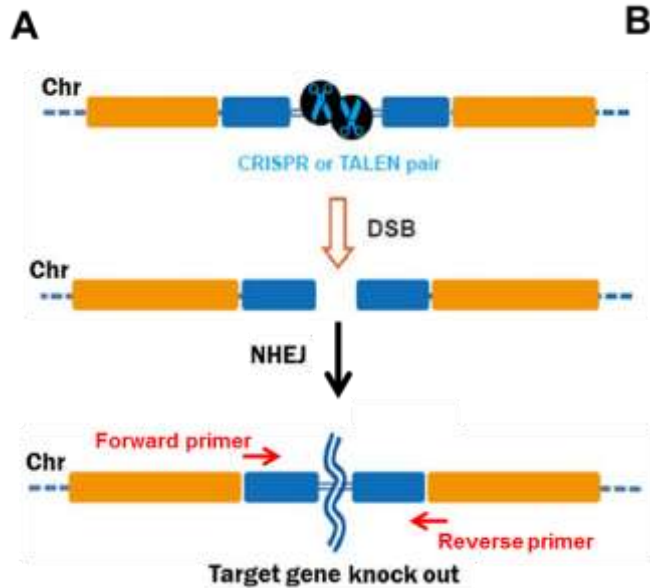
- ❖ Cell culture-based genome editing projects can take 3 or more months to complete
- ❖ Not all CRISPR sgRNAs are created equal! Test first before undergoing long genome editing project to avoid wasting time and expense.
- ❖ Recommend cell culture based strategy using T7 Endonuclease I assay.

Functional validation



Functional validation

IndelCheck™ T7 Endonuclease System

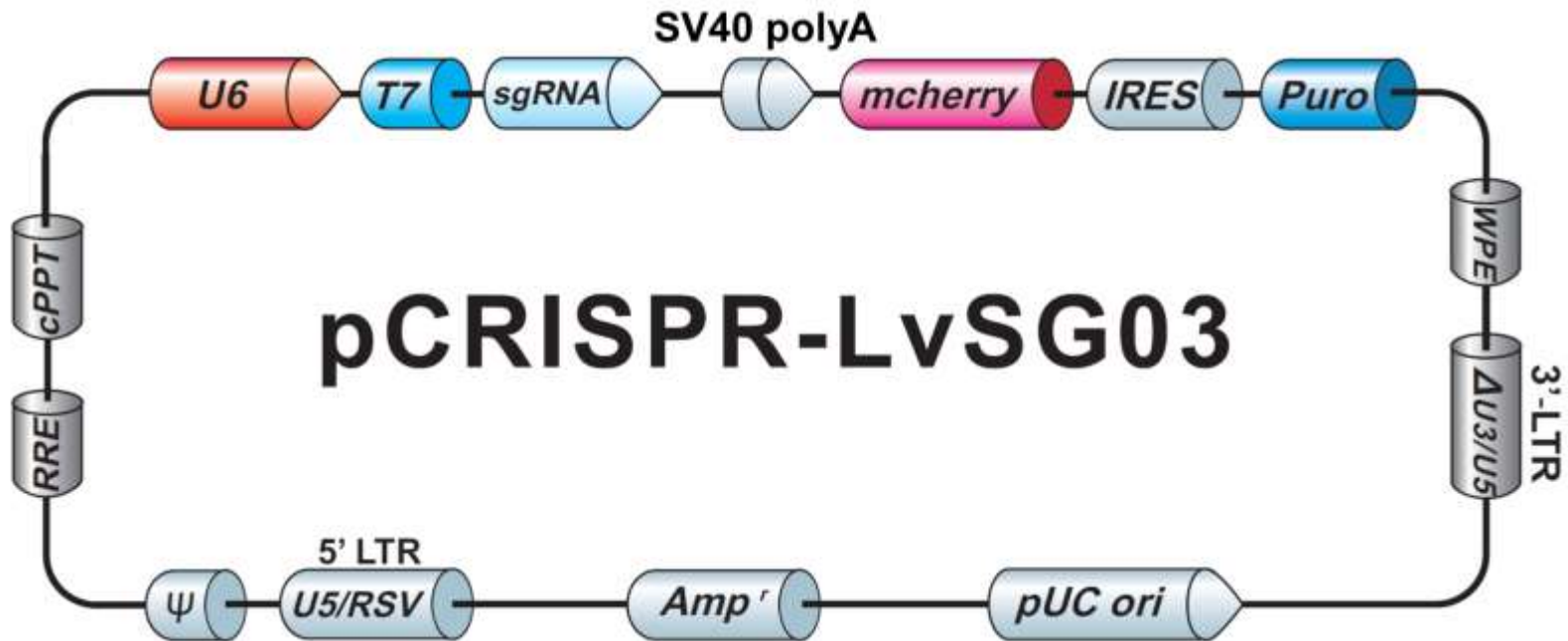


- ❖ Contains PCR and T7 endonuclease I reagents
- ❖ Purchase with or without target-specific primers

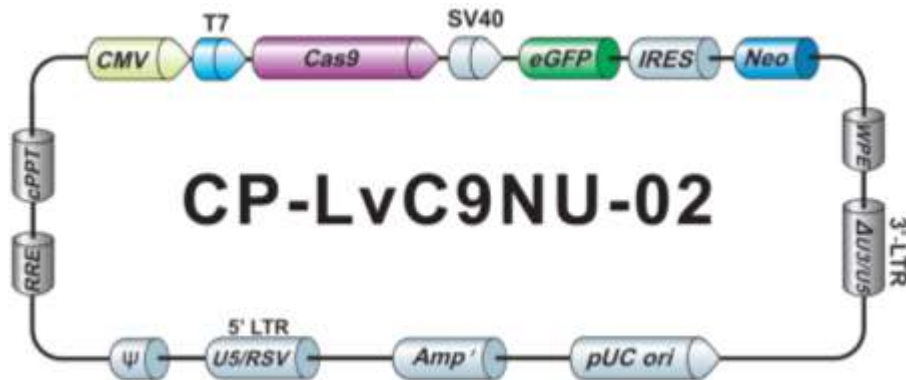
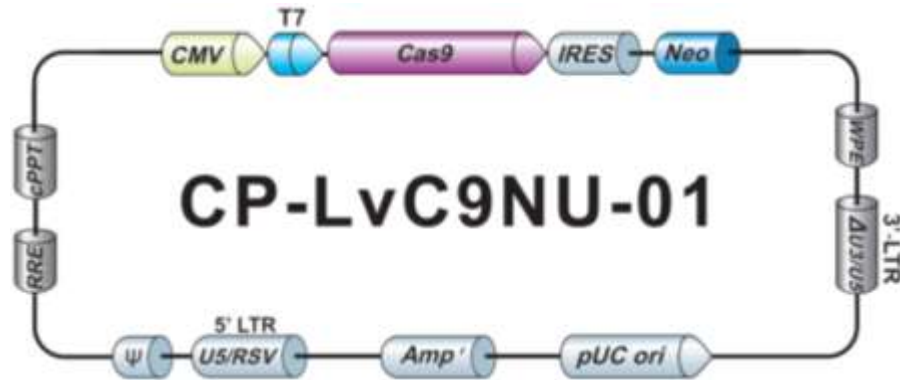
Outline

- ❖ Genome editing: Technologies
- ❖ Applications for genome editing
- ❖ sgRNA validation
- ❖ **Lentiviral CRISPR**
- ❖ CRISPR library screening
- ❖ Inducible CRISPR

Lentiviral CRISPR



Lentiviral CRISPR



Outline

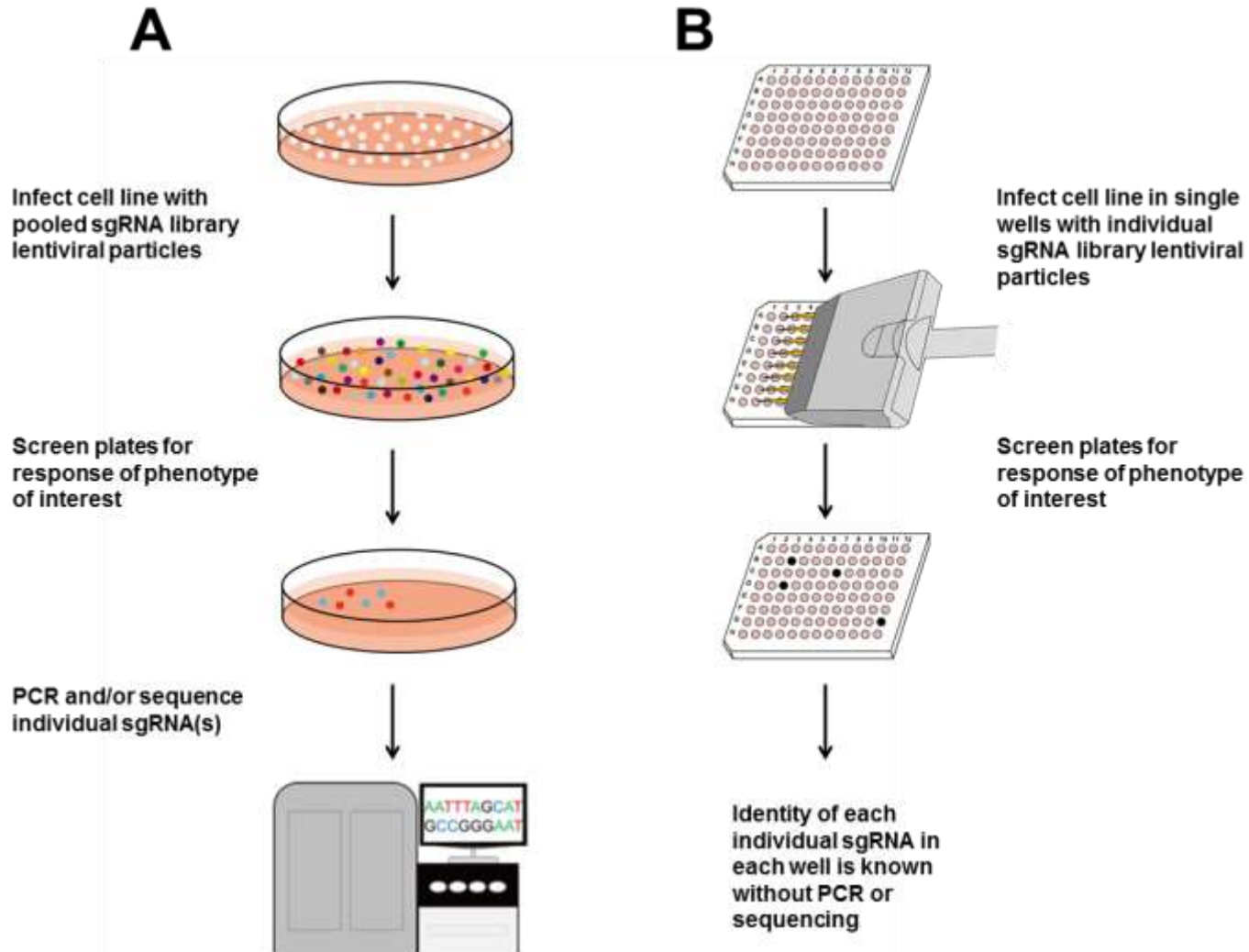
- ❖ Genome editing: Technologies
- ❖ Applications for genome editing
- ❖ sgRNA validation
- ❖ Lentiviral CRISPR
- ❖ **CRISPR library screening**
- ❖ Inducible CRISPR

CRISPR or RNAi?

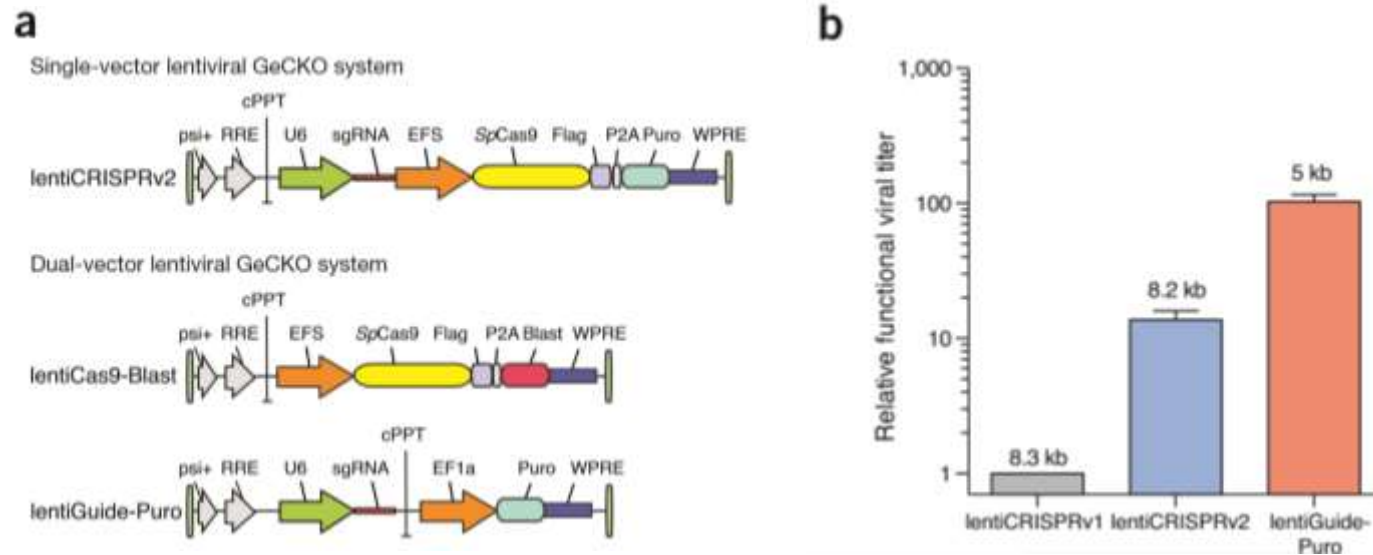
Knock down vs. Knock out

Method	Change expression level	Knock down	Knock out	Change genetic code
CRISPR	✓		✓	✓
RNAi	✓	✓		

CRISPR sgRNA libraries



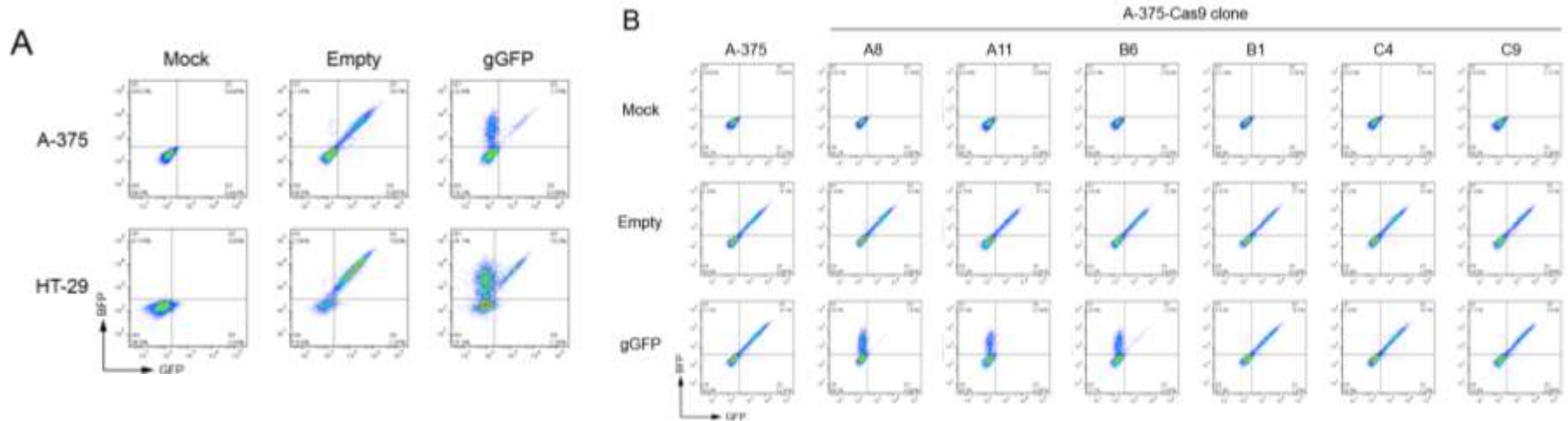
CRISPR sgRNA libraries



Sanjana, et al. (2014). *Nature Methods* 11, 783

- ❖ CRISPR libraries often used for high-throughput knockout or expression screening
- ❖ Viral titers of sgRNA-alone constructs is much higher than “all-in-one” (Cas9 + sgRNA) constructs

CRISPR sgRNA libraries



Tzelepis, et al. (2016). Cell Reports 17, 1193

- ❖ Transduced cells with Cas9-expressing lentivirus
- ❖ Found that bulk population had some cells that did not express Cas9
- ❖ Subcloned bulk cells to get clones with uniformly-expressing Cas9

GeneCopoeia sgRNA libraries

Pathway & gene group sgRNA libraries

Library name	Number of genes
Innate kinases & ubiquitin ligases	239
Nuclear hormone receptors	118
Tumor metastasis genes	57
Oncogenes	288
Tumor suppressor genes	231
Protein kinases	658
Key genes in 50 pathways	139
Custom	Made-to-order

Available as bacteria, DNA, or lentiviral particles either:

- A. Pooled**
- B. Individually arrayed**

GeneCopoeia sgRNA libraries

Advantages

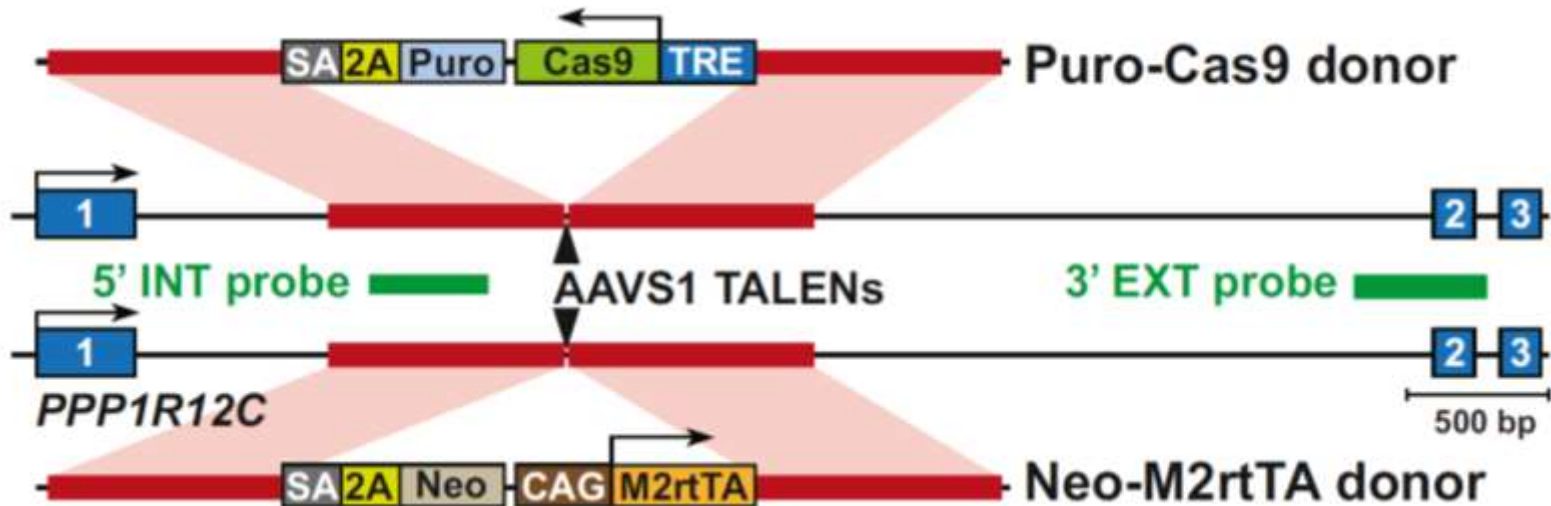
- ❖ Individually constructed and cultured in *E. coli* before pooling. Makes possible use in pools or as individual sgRNAs.
- ❖ Pools limited to 150 sgRNAs, ensuring excellent representation of each sgRNA
- ❖ Sequence verification provides high quality of each sgRNA
- ❖ Small library sizes: Reduces time and cost of screening

Outline

- ❖ Genome editing: Technologies
- ❖ Applications for genome editing
- ❖ sgRNA validation
- ❖ Lentiviral CRISPR
- ❖ CRISPR library screening
- ❖ **Inducible CRISPR**

Inducible CRISPR

Tet system

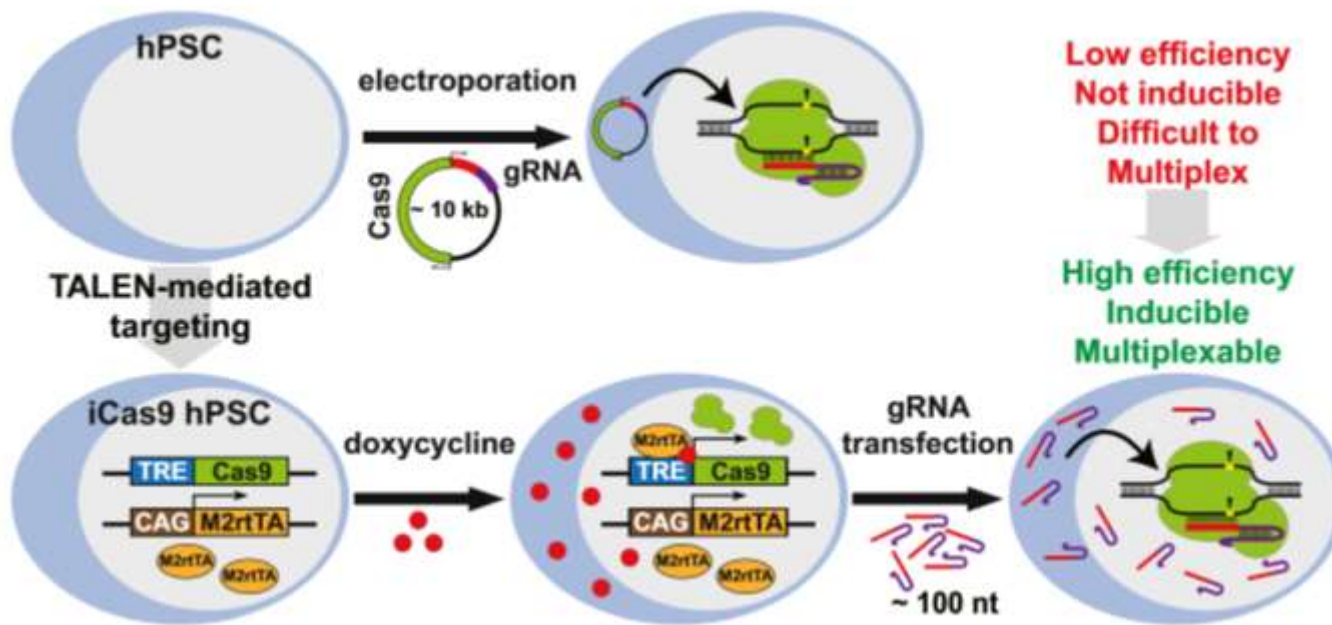


Gonzalez, et al. (2014). Cell Stem Cell 15, 1

- ❖ One chromosome: Construct donor with Tet-inducible promoter at AAVS1. Marked with puro.
- ❖ Other chromosome: Construct donor with CMV-driven tet transactivator (M2rtTA) at AAVS1. Marked with neo.
- ❖ Used TALEN with puro + G418 selection in one transfection for knockin at AAVS1.

Inducible CRISPR

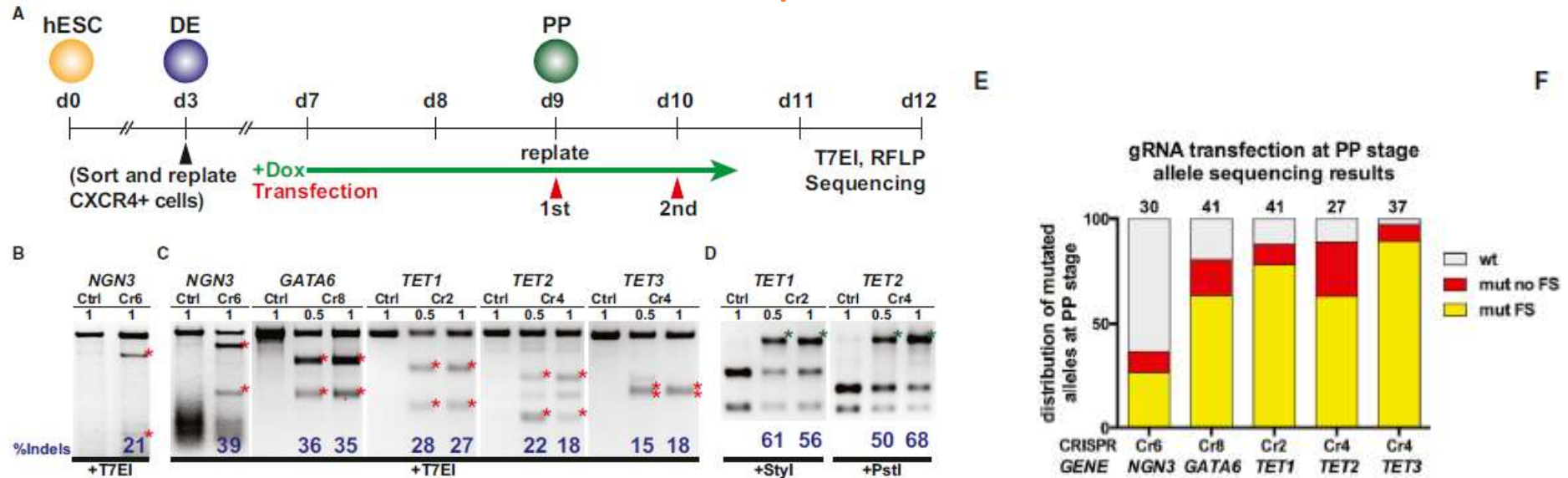
Tet system



Gonxalez, et al. (2014). Cell Stem Cell 15, 1

Inducible CRISPR

Tet system



- ❖ 2 cell lines. Differentiated to pancreatic progenitor (PP) stage in absence of dox and sgRNA (A).
- ❖ At PP stage, transfected sgRNAs and added dox
- ❖ T7 (B, C, and D) shows stage-specific induction of indels at 5 genes.
- ❖ Sequencing (E) shows that mutation frequencies are very high and most are indels.

Cas9-expressing stable cell lines

Human lung cancer cell lines

Catalog #	Cell line	Selection marker	Cas9 integration site
SL501	NCI-1299	Puro	AAVS1
SL533	NCI-H1299	Hygro	Random
SL504	A549	Hygro	AAVS1
SL529	H1975	Hygro	Random
SL536	NCI-H1437	Hygro	Random

Cas9-expressing stable cell lines

Human cervical cancer cell line

Catalog #	Cell line	Selection marker	Cas9 integration site
SL503	HeLa	Hygro	AAVS1

Cas9-expressing stable cell lines

Human breast cancer cell lines

Catalog #	Cell line	Selection marker	Cas9 integration site
SL514	MCF-7	Hygro	AAVS1
SL524	MCF-7	Hygro	Random
SL515	MDA-MB-231	Hygro	Random
SL516	MDA-MB-468	Hygro	Random
SL517	T47D	Hygro	Random
SL526	SNU-1	Hygro	Random
SL535	SK-BR-3	Hygro	Random
SL544	HCC70	Hygro	Random
SL547	DU4475	Hygro	Random

Cas9-expressing stable cell lines

Human liver cancer cell lines

Catalog #	Cell line	Selection marker	Cas9 integration site
SL518	HepG2	Hygro	AAVS1
SL522	SNU-475	Hygro	Random
SL530	SNU-449	Hygro	Random
SL531	PLC/PRF/5	Hygro	Random
SL537	SNU-387	Hygro	Random
SL538	SNU-423	Hygro	Random
SL543	C3A	Hygro	Random

Cas9-expressing stable cell lines

Human stomach/gastric cancer cell lines

Catalog #	Cell line	Selection marker	Cas9 integration site
SL520	AGS	Hygro	Random
SL527	KATO111	Hygro	Random
SL545	SNU-16	Hygro	Random

Cas9-expressing stable cell lines

Human pancreatic cancer cell lines

Catalog #	Cell line	Selection marker	Cas9 integration site
SL521	BXPC-3	Hygro	Random
SL540	CFPAC-1*	Hygro	Random

*derived from metastatic liver

Cas9-expressing stable cell lines

Human colon cancer cell lines

Catalog #	Cell line	Selection marker	Cas9 integration site
SL523	HT-29	Hygro	Random
SL525	LoVo	Hygro	Random
SL539	RKO	Hygro	Random
SL541	T84	Hygro	Random
SL542	COLO 205*	Hygro	Random

*derived from metastatic site

Cas9-expressing stable cell lines

Human caecum cancer cell line

Catalog #	Cell line	Selection marker	Cas9 integration site
SL532	LS411N	Hygro	Random

Cas9-expressing stable cell lines

Human bone/bone marrow/blood cancer cell lines

Catalog #	Cell line	Selection marker	Cas9 integration site
SL546	U-2 OS	Hygro	Random
SL552	K562	Hygro	Random
SL555	Jurkat	Hygro	Random

Cas9-expressing stable cell lines

Human embryonic kidney cell line

Catalog #	Cell line	Selection marker	Cas9 integration site
SL502	HEK293T	Puro	AAVS1
SL553	HEK293T*	Hygro	AAVS1

*Expresses high-fidelity Cas9

Cas9-expressing stable cell lines

Human neuroblastoma cancer cell line

Catalog #	Cell line	Selection marker	Cas9 integration site
SL550	SH-SY5Y	Hygro	Random

Cas9-expressing stable cell lines

Mouse neuroblastoma cancer cell lines

Catalog #	Cell line	Selection marker	Cas9 integration site
SL509	Neuro2a	Hygro	ROSA26
SL510	Neuro2a	Puro	ROSA26
SL511	Neuro2a	Neo	ROSA26

Cas9-expressing stable cell lines

Rat glioma cancer cell line

Catalog #	Cell line	Selection marker	Cas9 integration site
SL534	C6	Hygro	Random

Summary

- ❖ CRISPR is a highly effective method for many applications, from knockout, knockin, activation, & more
- ❖ GeneCopoeia provides a large number of cell lines stably expressing the CRISPR-Cas9 nuclease
- ❖ GeneCopoeia's CRISPR-Cas9-expressing stable cell lines provide great utility for several genome editing applications, such as:
 - ❖ sgRNA functional validation
 - ❖ Lentiviral CRISPR
 - ❖ sgRNA library screening
 - ❖ Inducible CRISPR



Cas9-expressing stable cell lines

GeneCopoeia Technical Note: Cas9 stable cell lines



TECHNICAL NOTE

Genome Editing: Cas9 Stable Cell Lines for CRISPR sgRNA Validation, Library Screening, and More

Ed Davis, Ph.D.

Introduction

The CRISPR-Cas9 system has become greatly popular for genome editing in recent years, due to its ease-of-design, efficiency, specificity, and relatively low cost (Wang, et al., 2016). In mammalian cell culture systems, most genome editing is achieved using transient transfection or lentiviral transduction, which works well for routine, low-throughput applications. However, for other applications, it would be beneficial to have a system in which one component, namely the CRISPR-Cas9 nuclease, was stably integrated into the genome. In this Technical Note, we introduce GeneCopoeia's suite of Cas9 stable cell lines, and discuss the great utility that these cell lines provide for genome editing applications.

The CRISPR-Cas9 system for genome editing has been adapted from anti-viral immunity mechanisms present in many bacterial species. The most commonly used CRISPR-Cas9 system to date is from *Streptococcus pyogenes* (SpCas9). SpCas9 is a DNA endonuclease that, in the presence of a chimeric

Download from:

<http://www.genecopoeia.com/wp-content/uploads/2017/03/GeneCopoeia-Technical-Note-Cas9-stable-cell-lines-03-2017.pdf>

Upcoming webinar!

**How To Choose Between
Adeno-associated Virus and
Lentivirus For DNA Delivery**

Wednesday, April 12, 2017 12:00 pm ET

Register here:

[https://attendee.gotowebinar.com/register/14390983
74222287363](https://attendee.gotowebinar.com/register/1439098374222287363)

Upcoming webinar!

**Applications For Safe Harbor
Transgenesis in Genome
Editing**

Wednesday, April 19, 2017 12:00 pm ET

Register here:

[https://attendee.gotowebinar.com/register/73880445
19385274881](https://attendee.gotowebinar.com/register/7388044519385274881)

Thank you!

If you have any additional
questions, please call

1-866-360-9531 x227

Email: edavis@genecopoeia.com

Or visit us on the web:

www.genecopoeia.com

GeneCopoeia, Inc.

9260 Medical Center Drive Suite 101

Rockville, Maryland USA 20850

