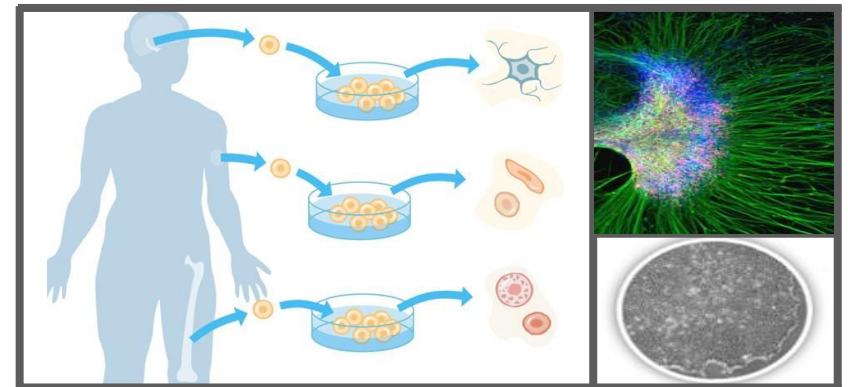


Clontech **Takara** cellartis

Efficient, Footprint-Free Gene Editing and Single-Cell Cloning of iPS Cells Using CRISPR/Cas9

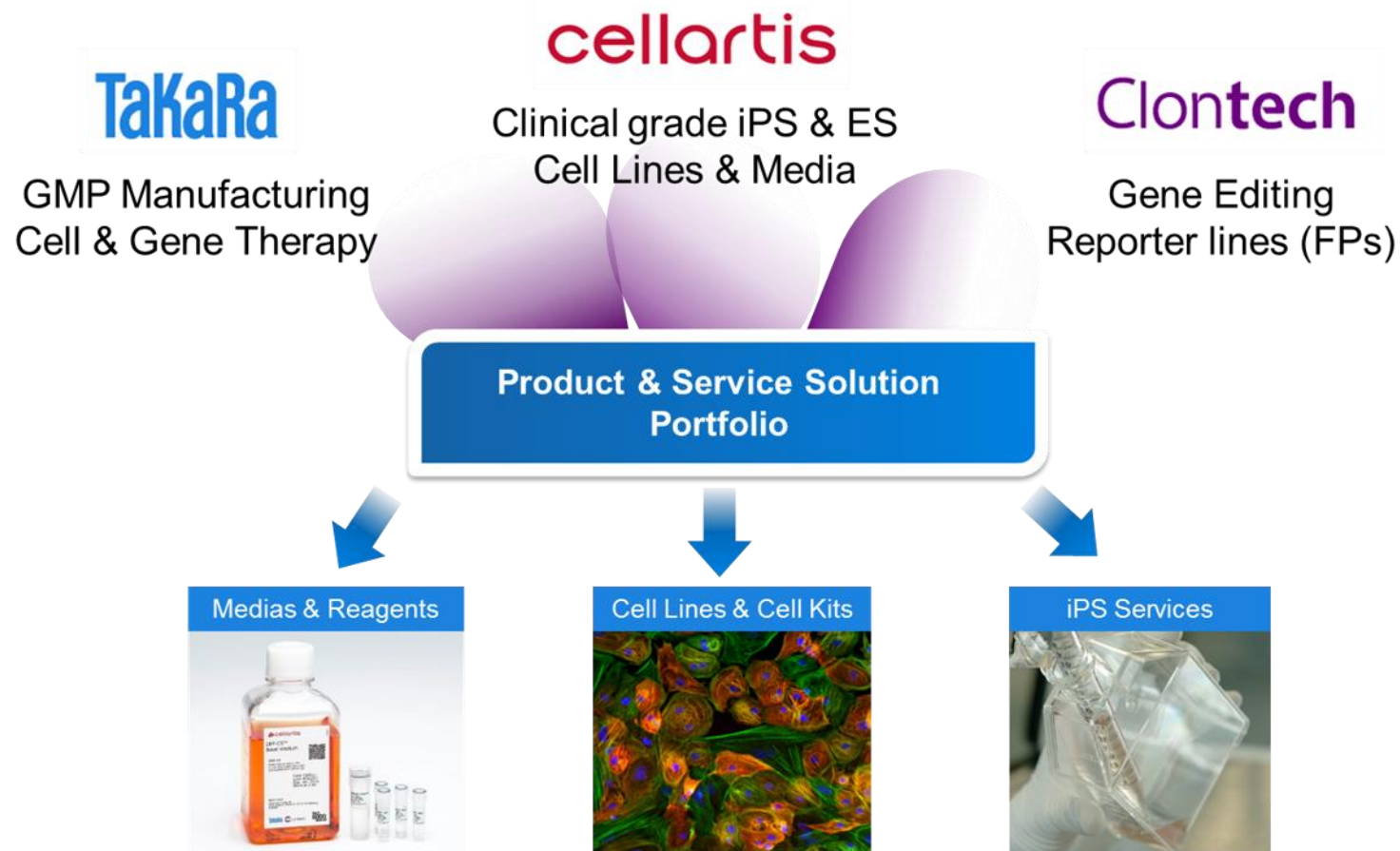
Liz Quinn, PhD

June 15, 2017



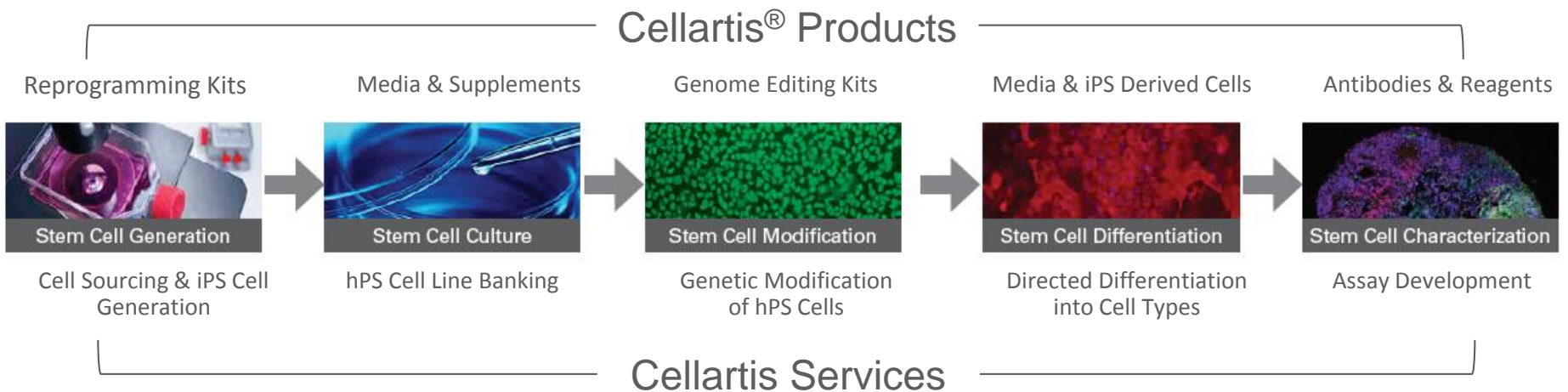
that's
GOOD
science!®

Stem Cell Research Expertise



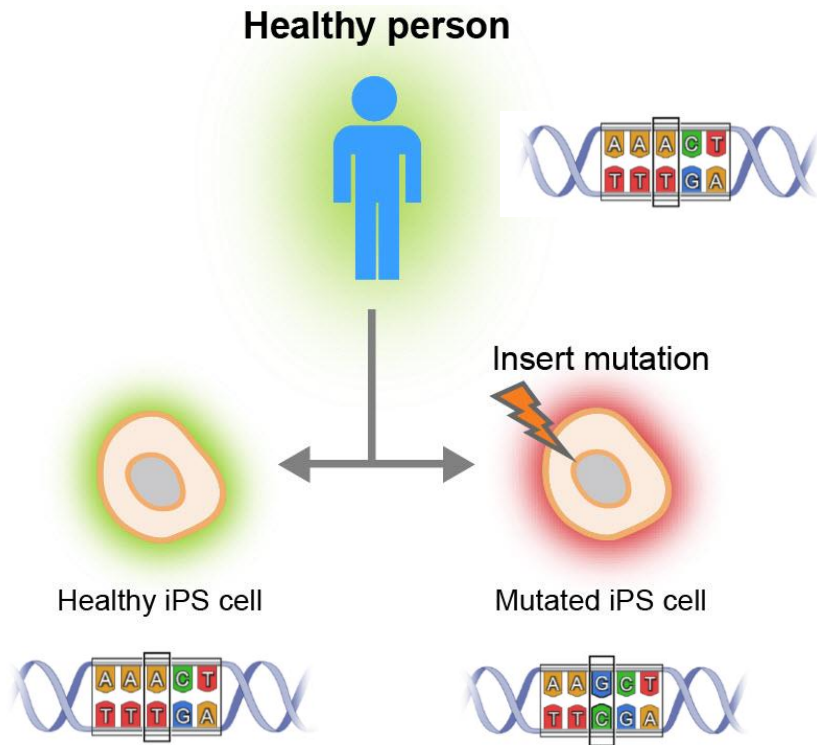
We focus on innovative research & clinical grade stem cell products & services

Comprehensive Products & Services



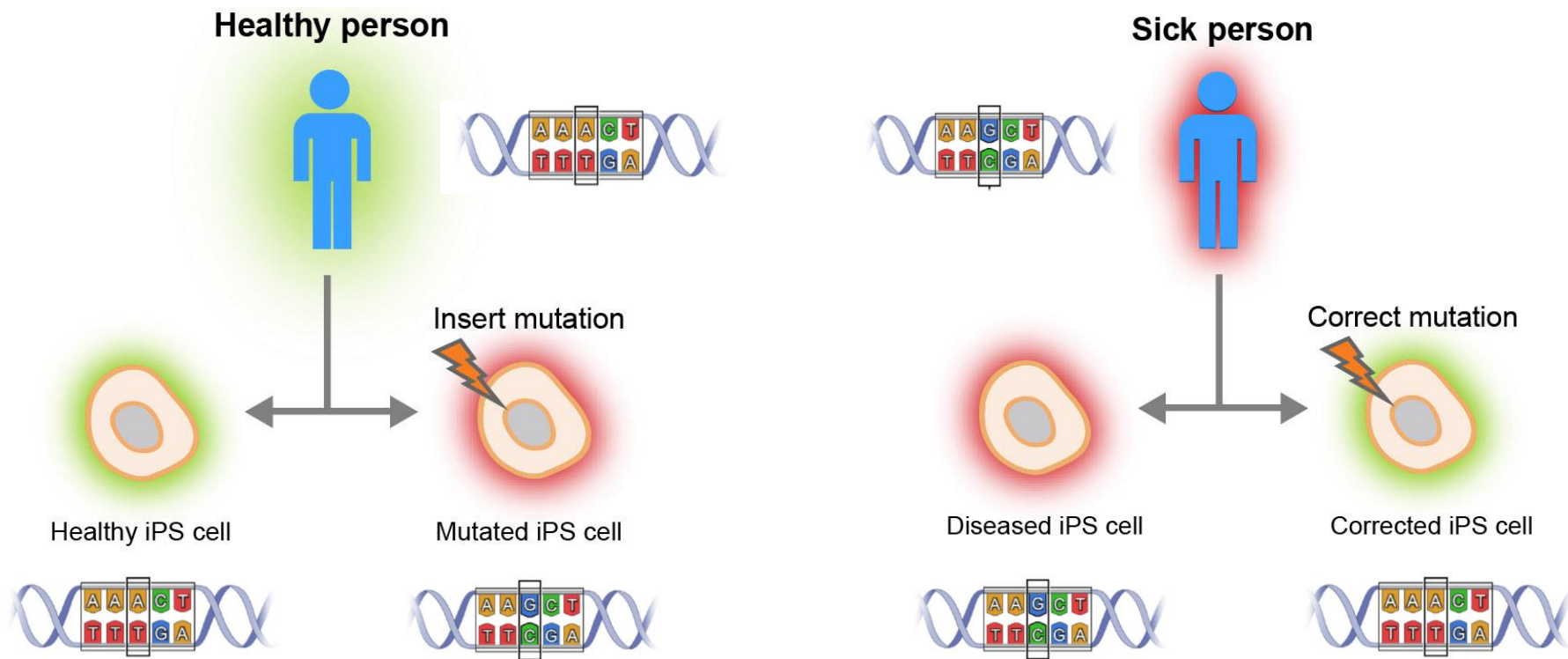
Utility of Stem Cells and Gene Editing

- Stem cells offer a renewable, expandable source of edited cells that can become multiple cell types



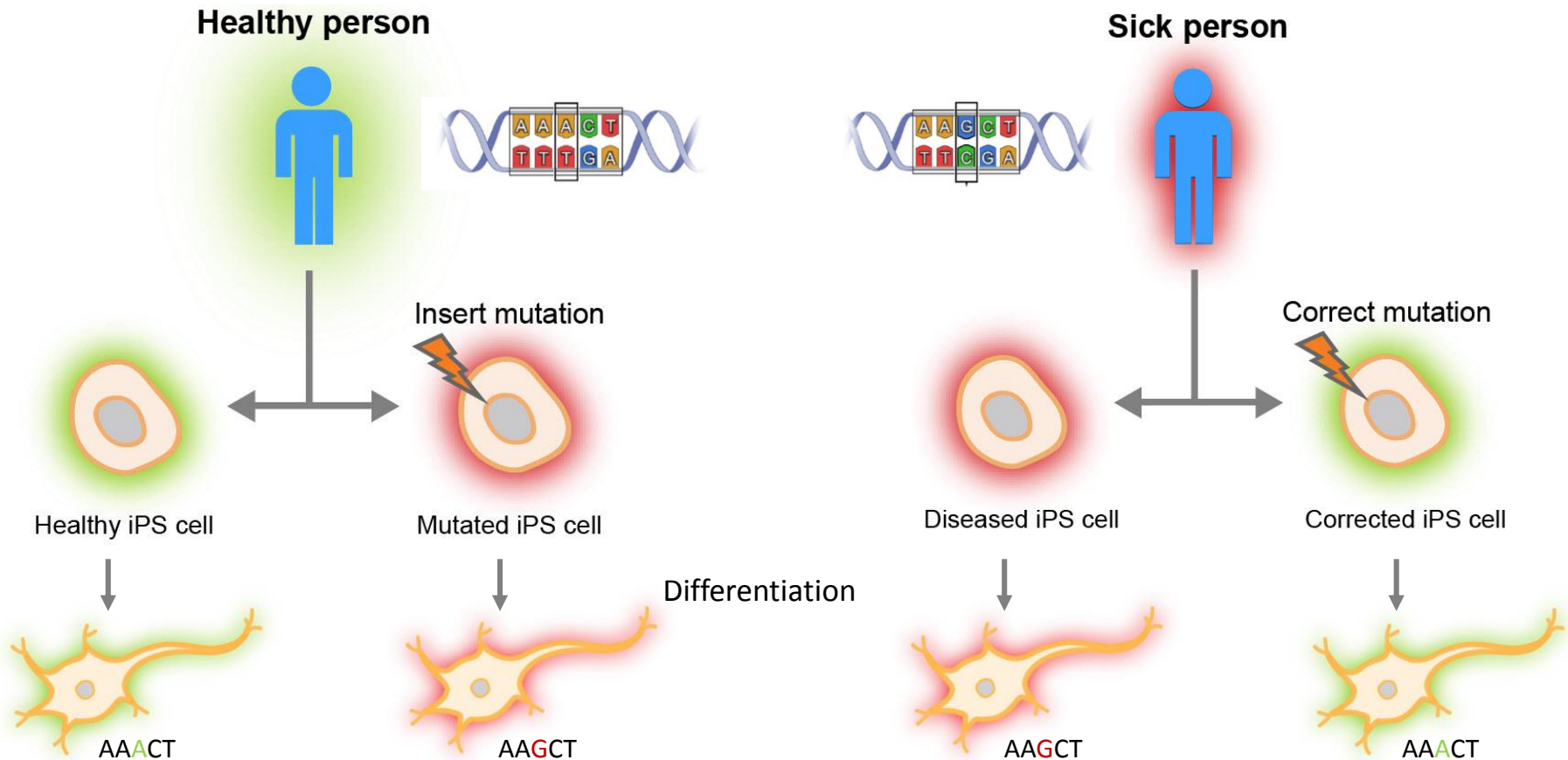
Utility of Stem Cells and Gene Editing

- Stem cells offer a renewable, expandable source of edited cells that can become multiple cell types

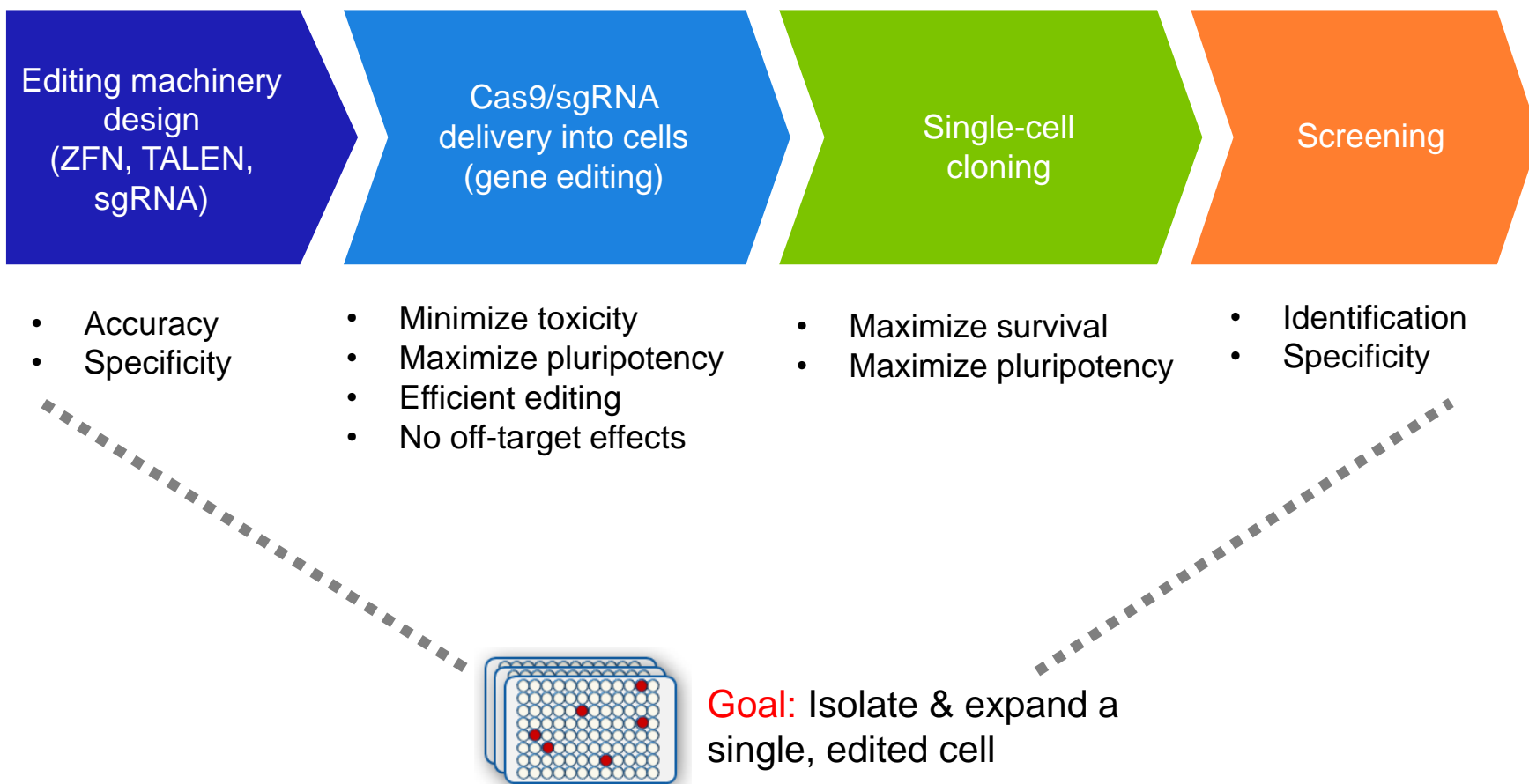


Utility of Stem Cells and Gene Editing

- Stem cells offer a renewable, expandable source of edited cells that can become multiple cell types

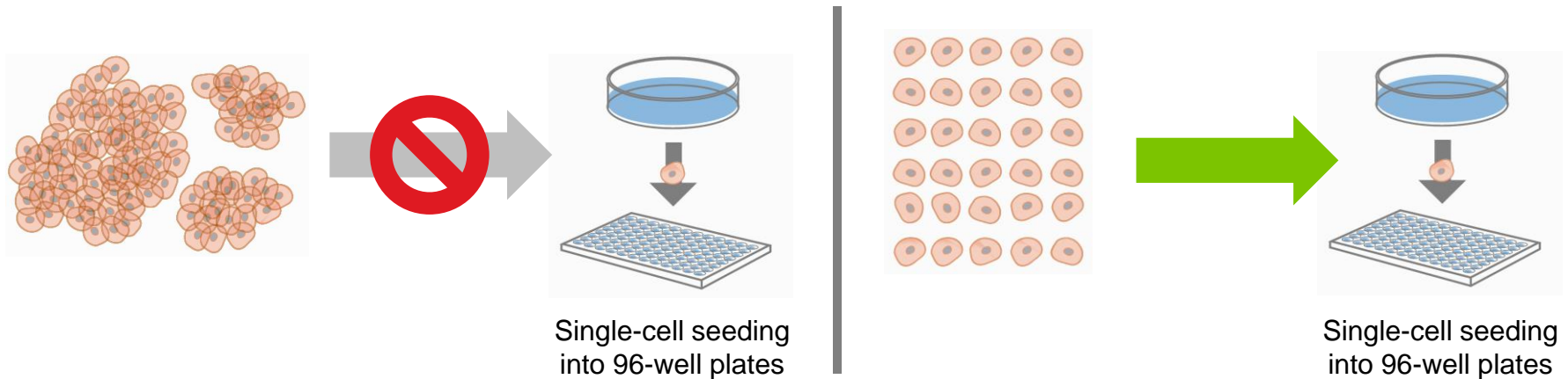


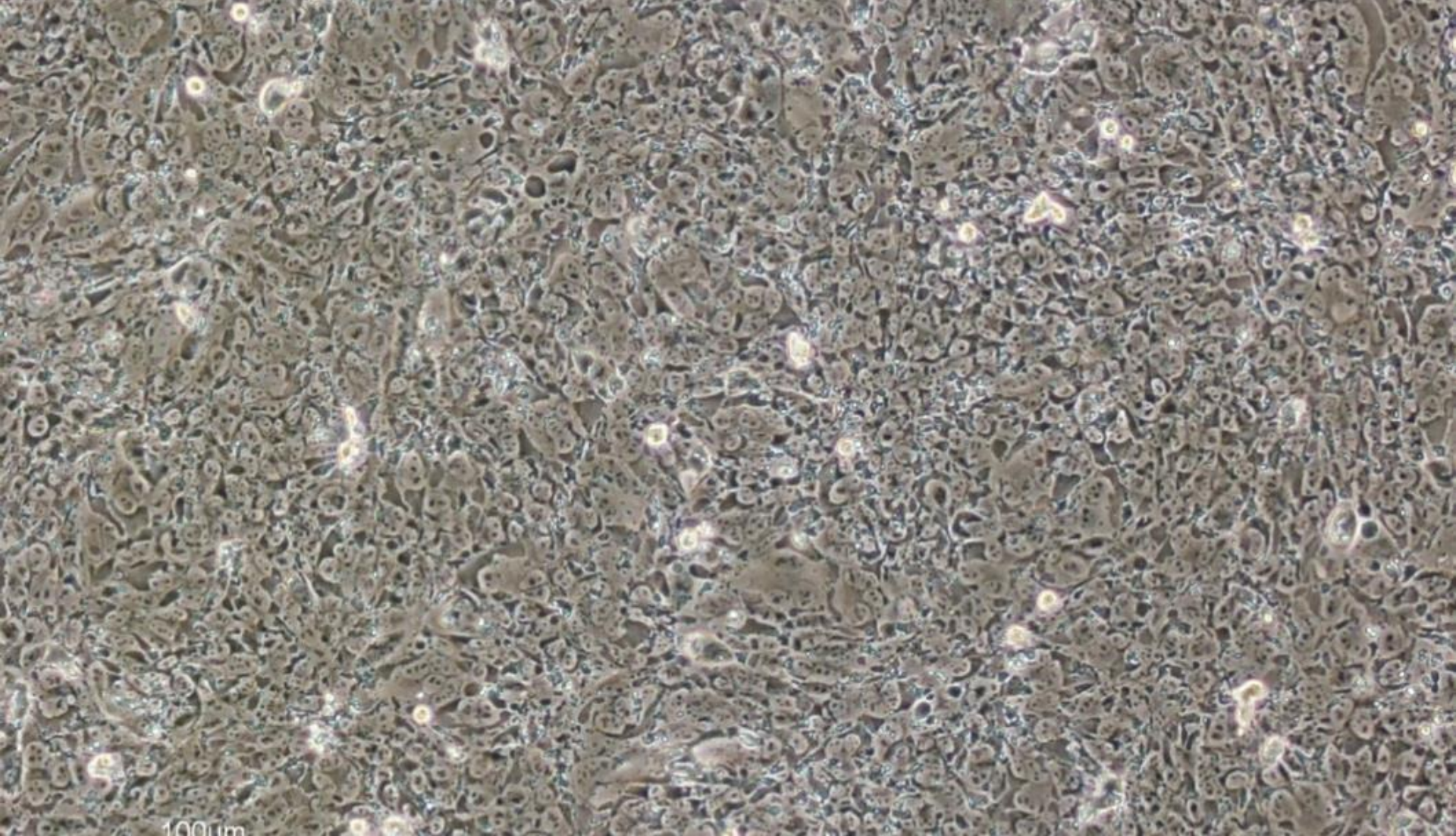
Challenges in Editing Stem Cells



Clonal Expansion of Single Pluripotent Stem Cells is a Bottleneck

- Isolating and clonally expanding edited cells
 - Pluripotent stem cells traditionally grow in colonies
 - Screening a colony is time-consuming and challenging
 - Single pluripotent cells die or differentiate
 - Need for single-cell culture of pluripotent stem cells

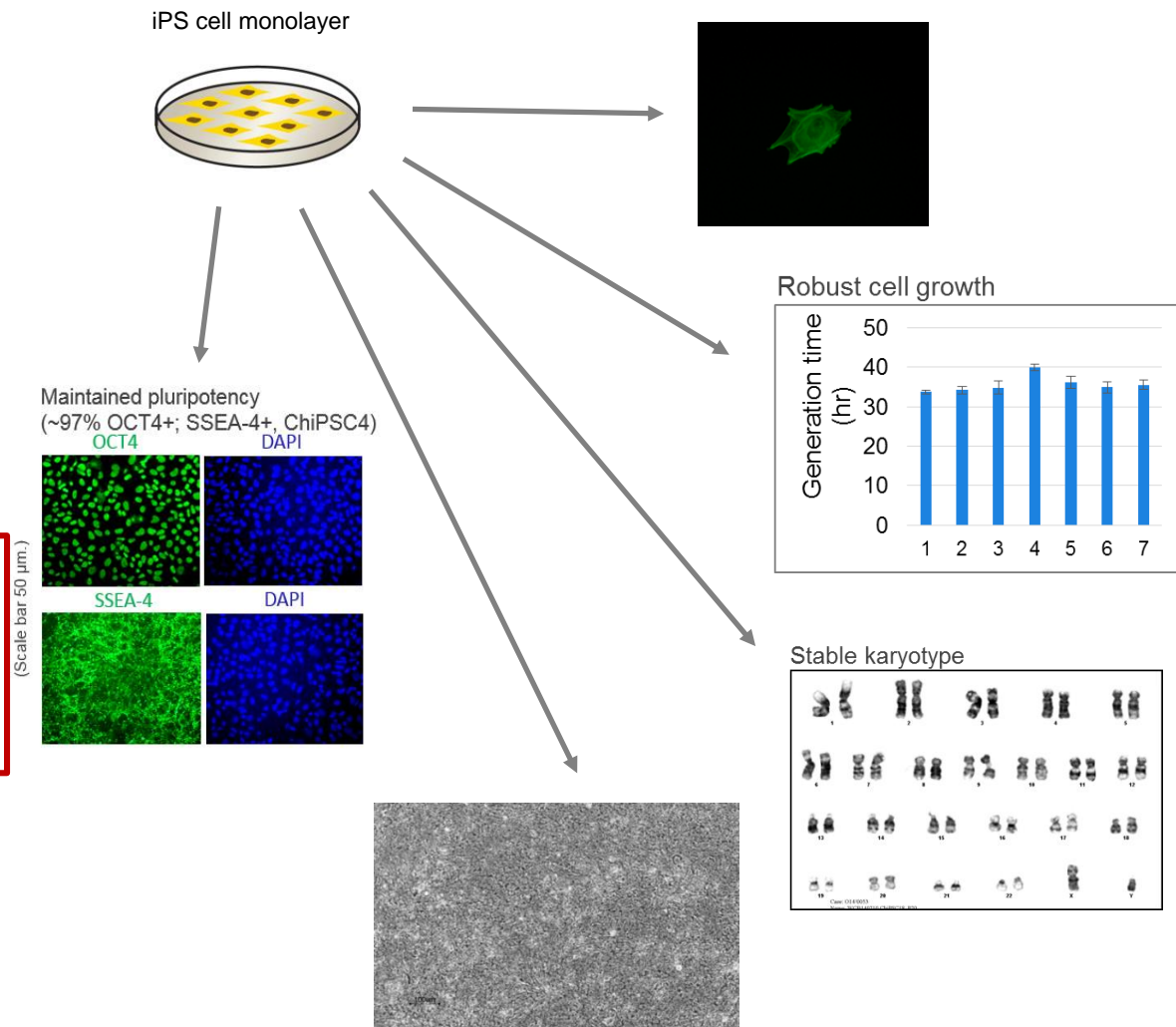




Human iPSC Culture System

Cellartis DEF-CS™ System & Single-Cell Cloning of iPSC

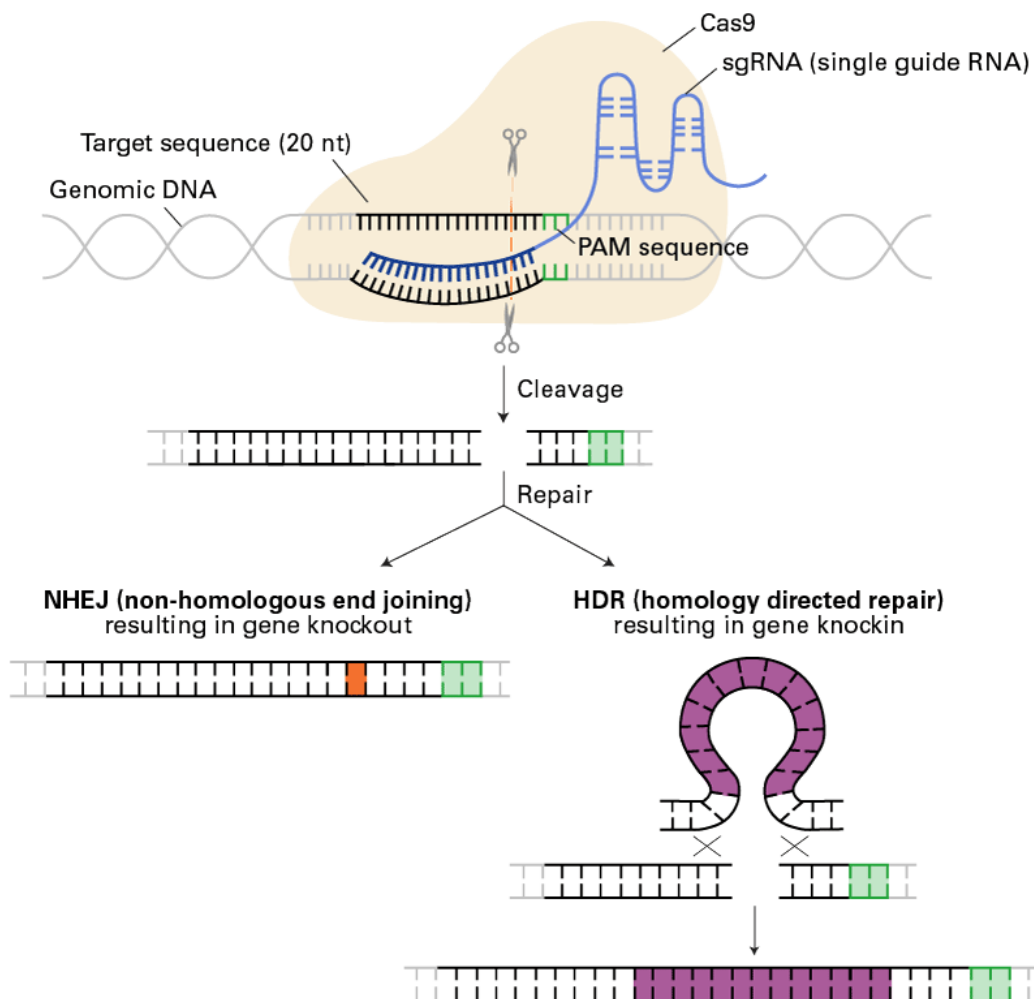
- Maintains cells in a highly undifferentiated state
- Allows for culturing iPSC cells in a monolayer
- Feeder-free — no contamination, less time consuming, increased consistency
- Enables survival and expansion of single cells
- Maintains normal karyotype
- Allows rapid expansion for further downstream applications and analysis





Gene Editing Tools

Gene Editing with Cas9



Repair template (HDR template) production methods

- AAV
- PCR
- Plasmid
- Synthetic oligos
- Takara Long ssDNA Kit

3 Components required:

- Cas9 nuclease
- Guide RNA
- Repair template

CRISPR/Cas9 Genome Editing

Target Gene



↓ Artificially create a Double-strand break to allow editing

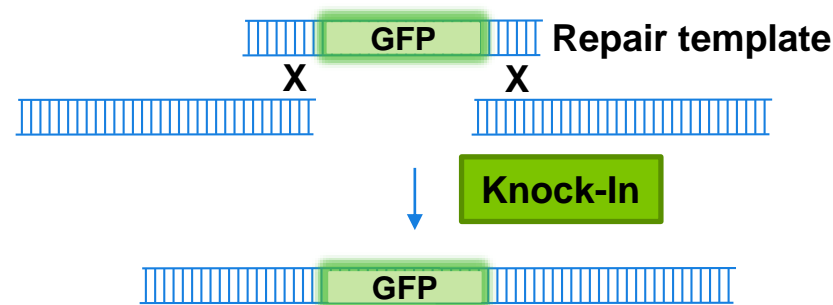
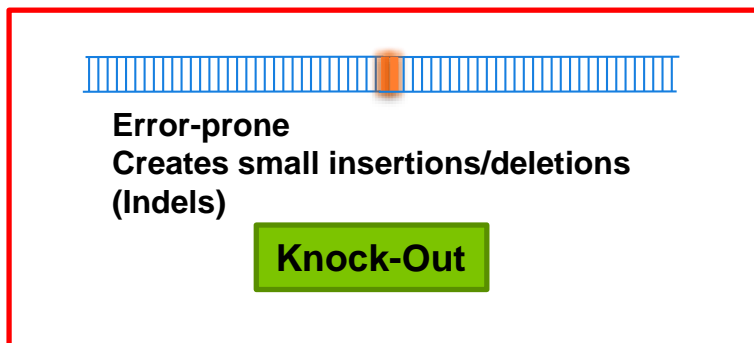


Non-Homologous End Joining (NHEJ)

“Quick repair”

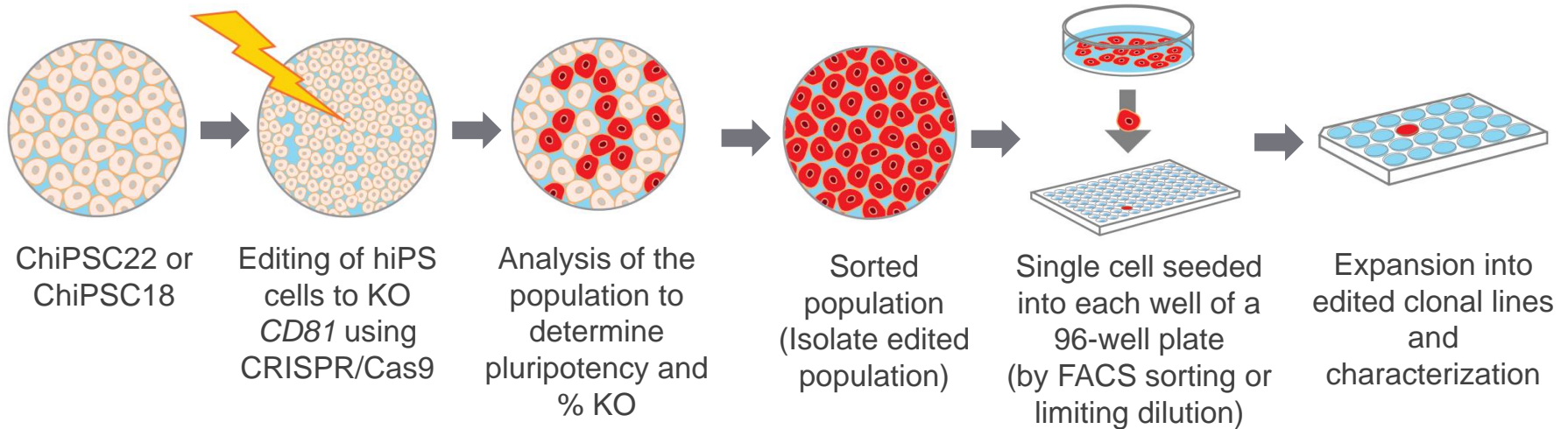
Homologous recombination (HR)

“Slower, more accurate repair”



- Delivery of two components
 - Cas9/sgRNA (RNP)
- Knockout efficiency via RNP acceptable/high even in primary cells or hiPS cells

CRISPR/Cas9 Gene Editing Workflow



Two options for Cas9 Delivery

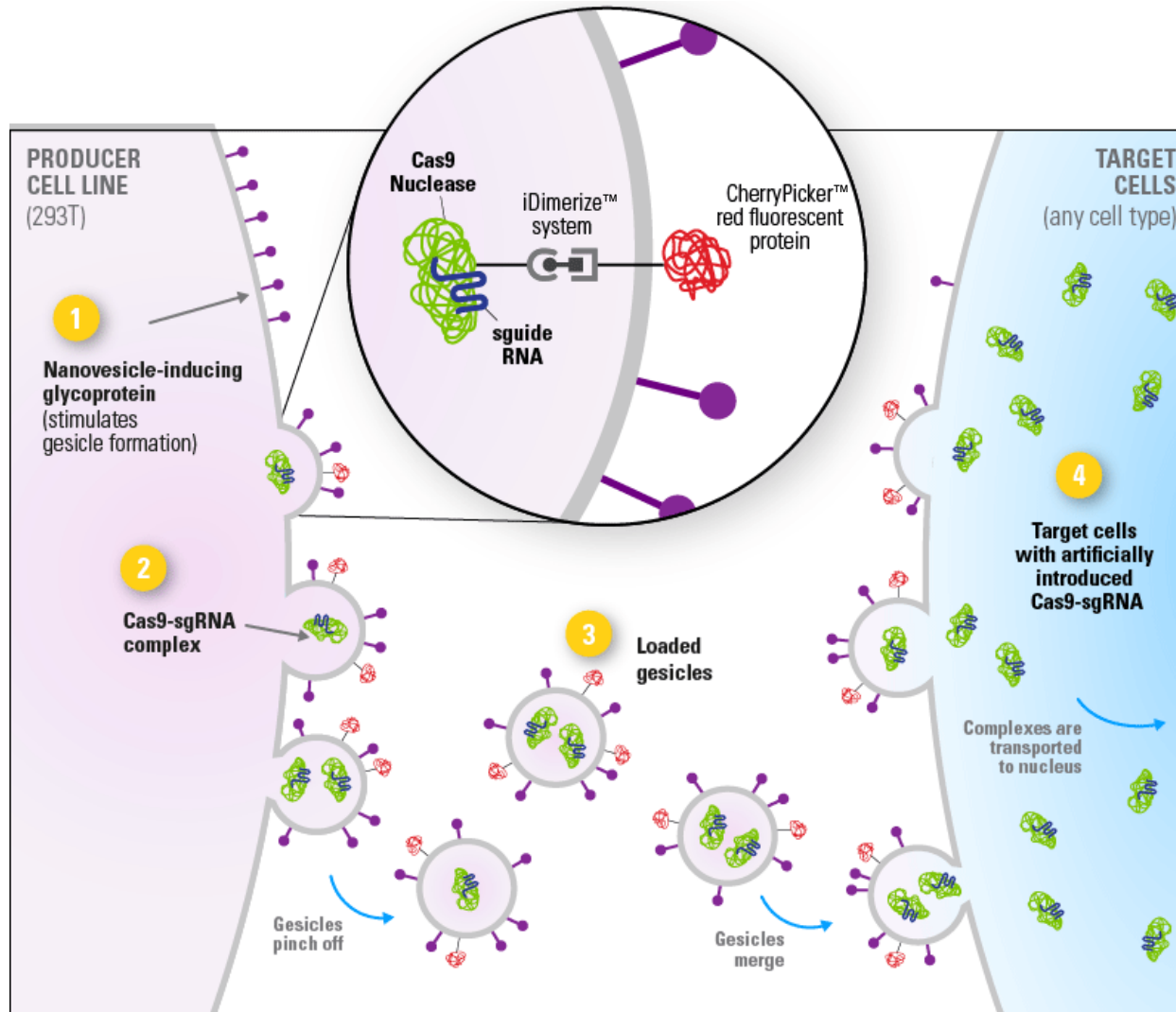
Guide-it™ rCas9 for electroporation:

- ~2 µg rCas9
- ~0.5 µg sgRNA (*in vitro transcribed*)
- ~0.5–1µg of donor DNA
- Using a Neon Electroporator

Guide-it CRISPR/Cas9 vesicle production kit:

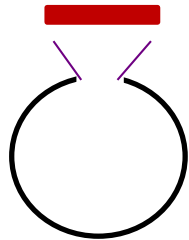
- Delivery of a Cas9/sgRNA RNP complex via cell derived nanovesicles

Guide-it CRISPR/Cas9 Gesicle Production Kit

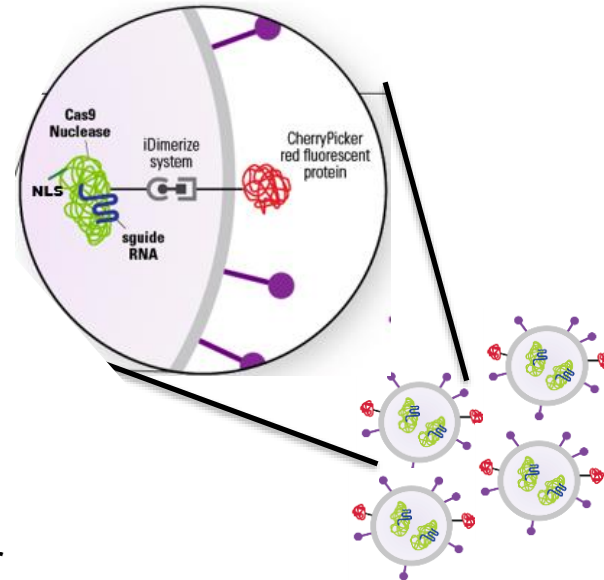


Guide-it CRISPR/Cas9 Gesicle Production Kit

sgRNA of interest



Pre-linearized sgRNA expression vector



Add 10µg of Plasmid encoding for sgRNA (600µl to final) to Packaging Mix1



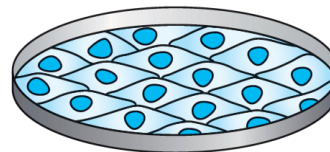
Add 600µl from Mix1 to Mix2



Vortex and incubate nanoparticles complexes for 10min

Apply nanoparticle complexes to sub-confluent (50-60%) cell culture

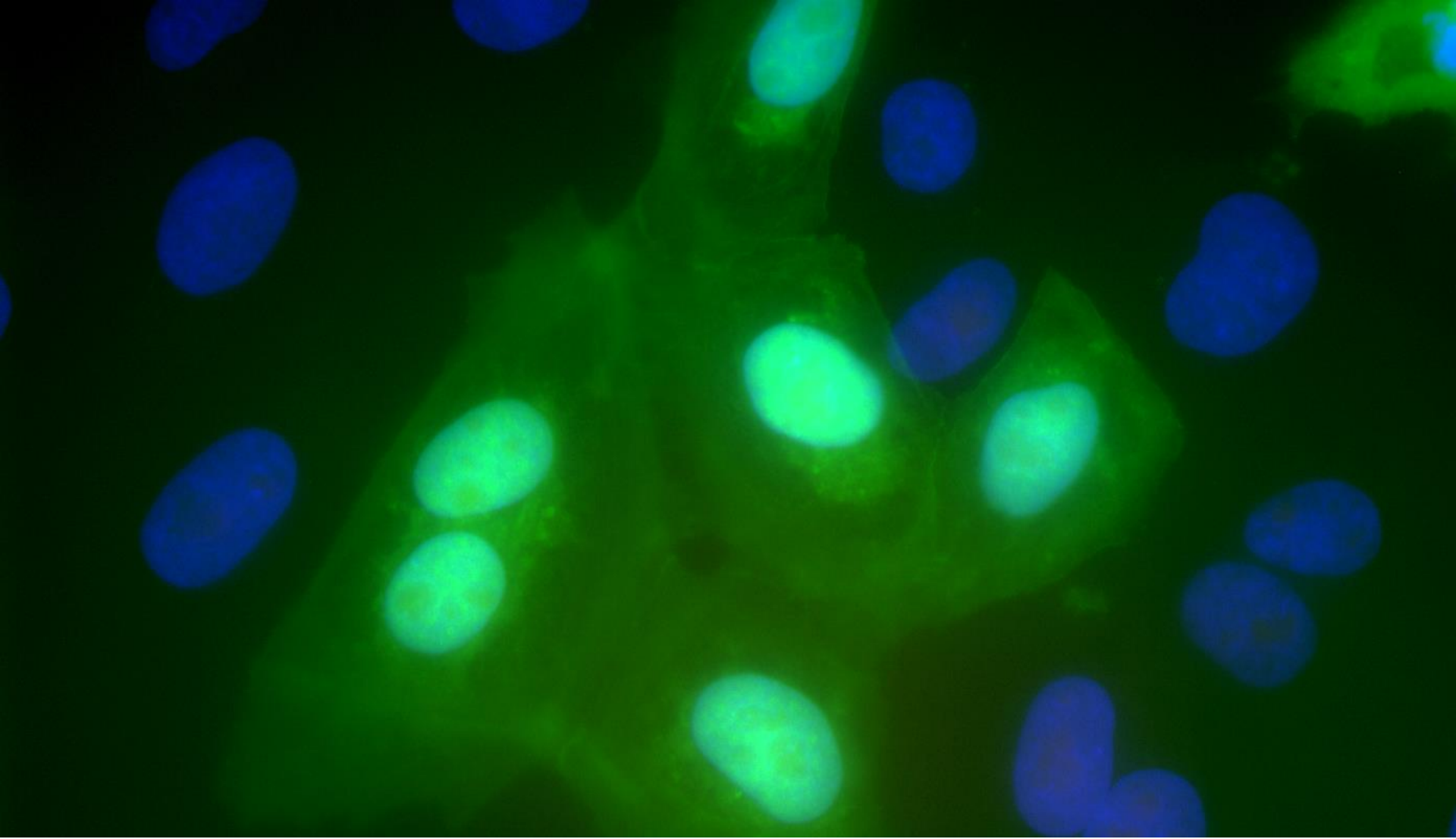
Add A/C Heterodimerizer



Incubate overnight at 37°C

Add fresh media

Collect Cas9 gesicle supernatant 48-72h post-transfection



Case Studies - Knock Out

Maintenance of Pluripotency after Gene Editing: *AcGFP1* KO Test Case

Cellartis Human iPS
Cell Line 22 (ChiPSC22)
stably expressing *AcGFP1*



CRISPR/Cas9
gene editing
targeting *AcGFP1*



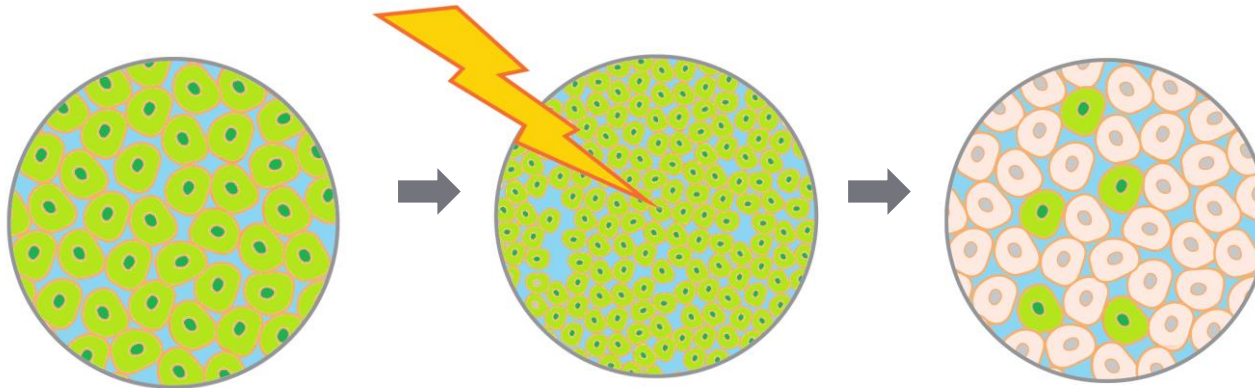
Cells grown
for 9 days



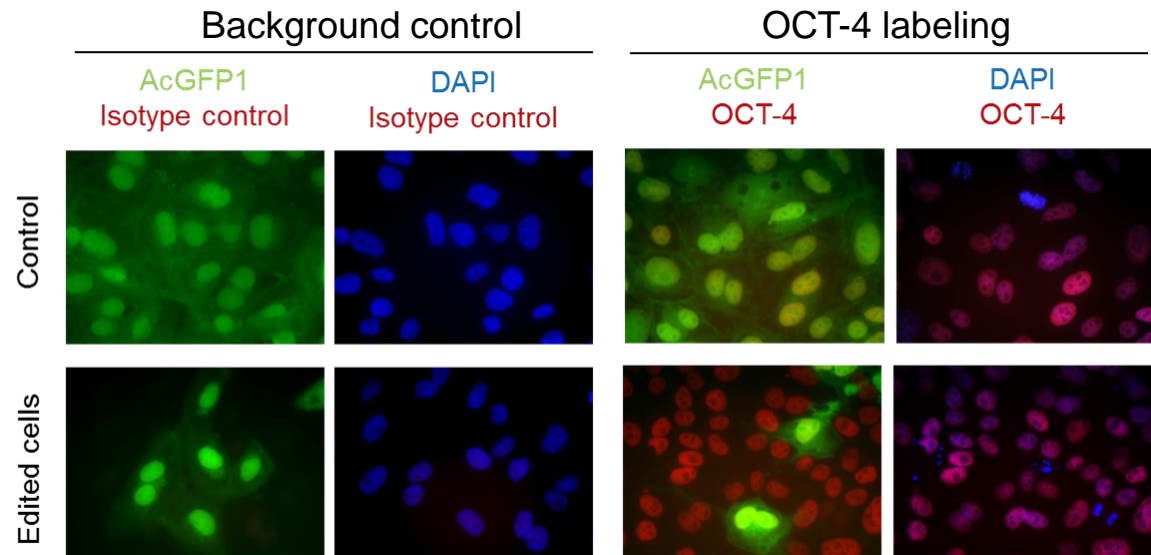
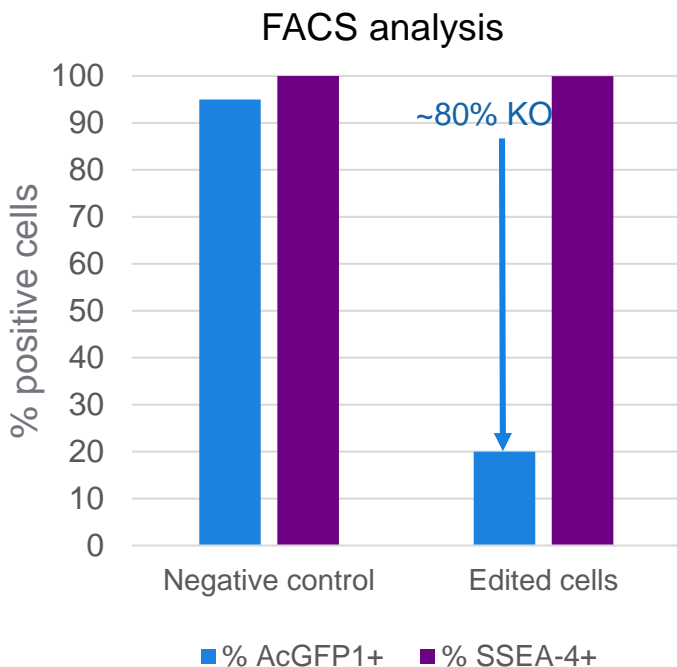
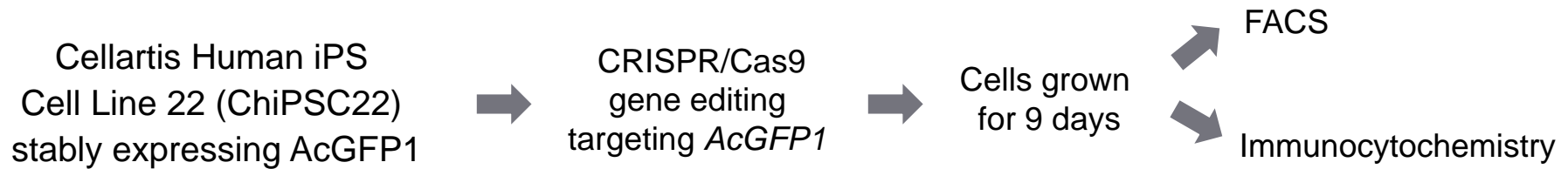
FACS



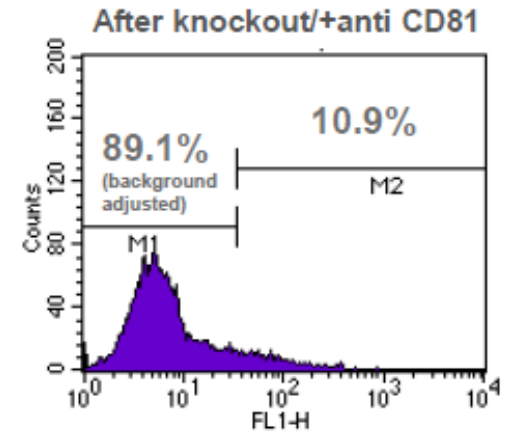
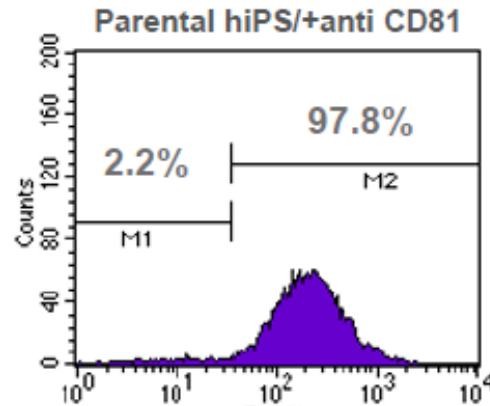
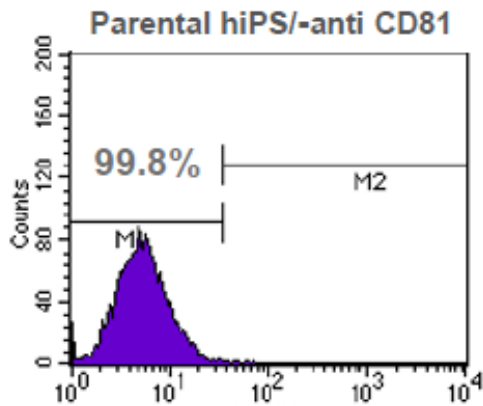
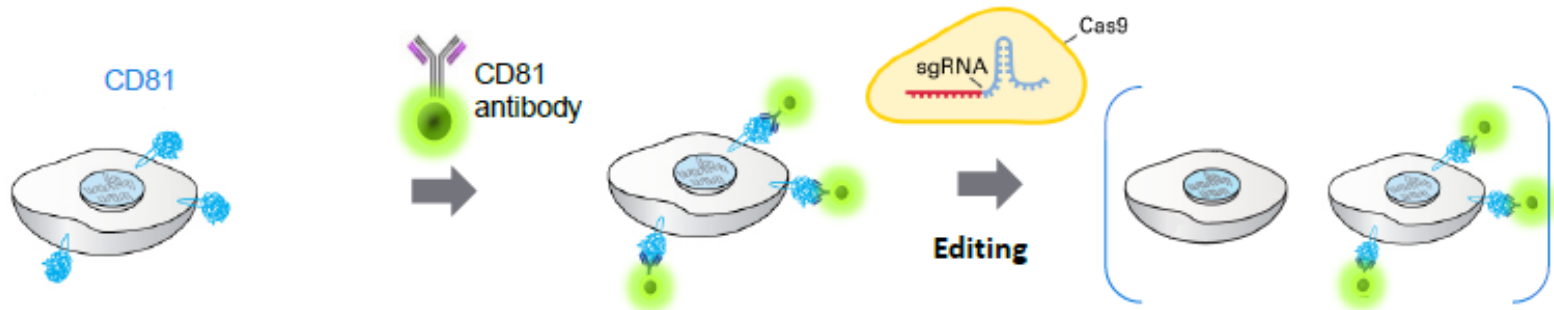
Immunocytochemistry



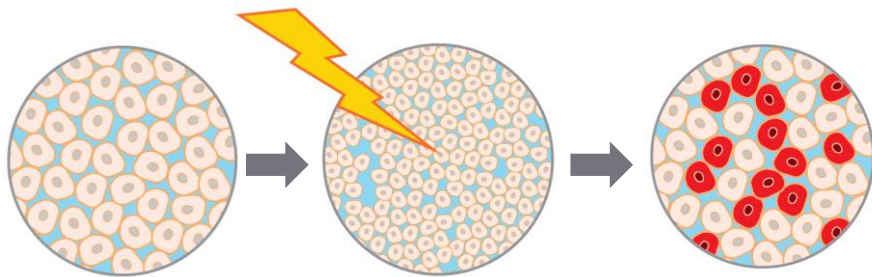
Maintenance of Pluripotency after Gene Editing: *AcGFP1* KO Test Case



Model System - CD81 KO in hiPSC



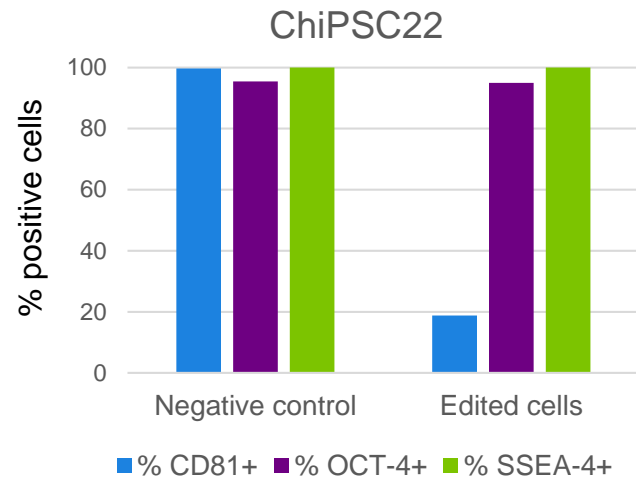
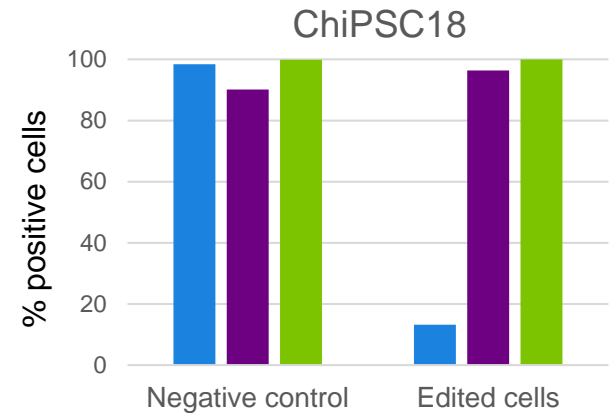
Pluripotency Maintained after *CD81* Knockout



ChiPSC22 or
ChiPSC18

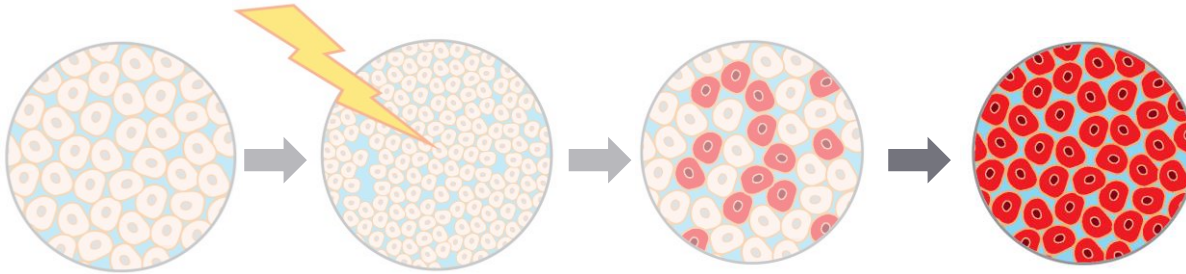
Editing of
hiPS cells to
KO *CD81*

FACS analysis
of the population
to determine
pluripotency and
% KO



Sorting the *CD81* Negative Cell Population

CD81/SSEA-4 double labeling

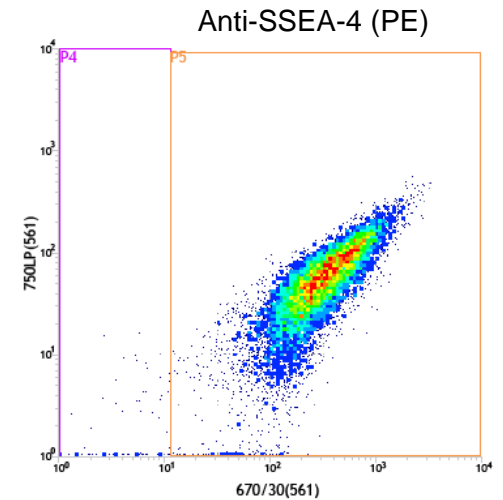
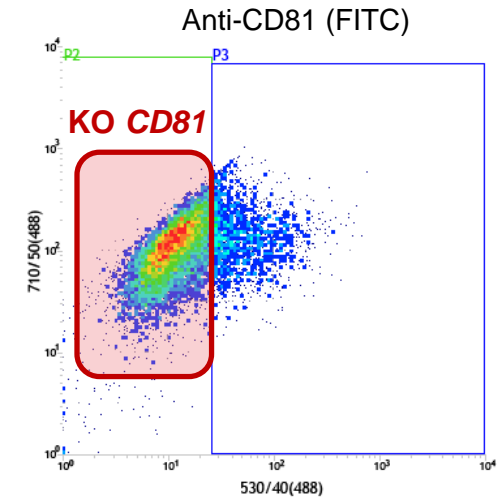


ChiPSC22 or
ChiPSC18

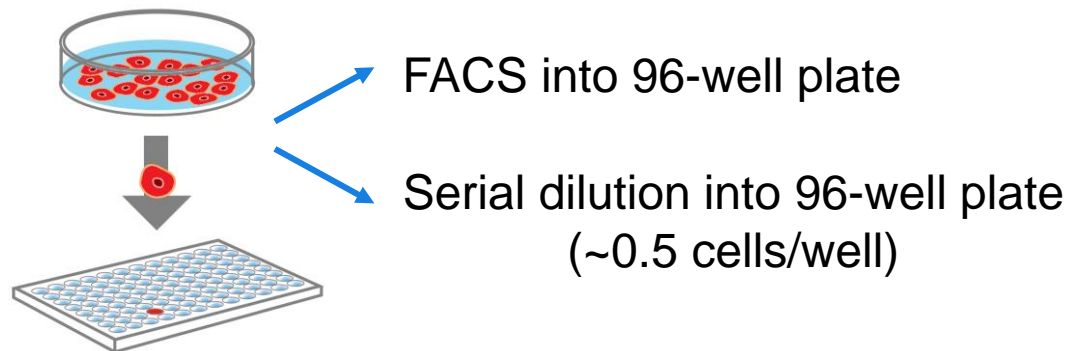
Editing of
hiPS cells to
KO *CD81*

FACS analysis
of the population
to determine
pluripotency and
% KO

Sorted
population

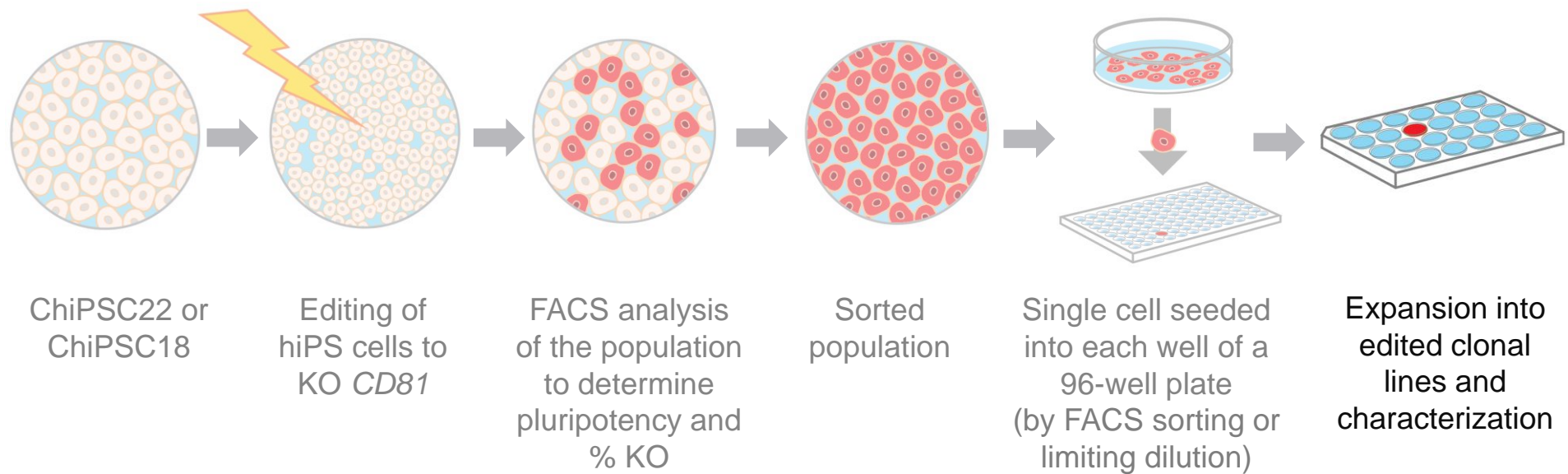


Cloning of Edited hiPSCs via FACS or Limiting Dilution



Cell line	Isolation method	Single clones	Double clones	Total clones (proportion)	Total clones (%)
ChiPSC22	FACS	8	0	8/96	8.5%
ChiPSC22	Limiting dilution	39	15	54/55	98%
ChiPSC18	FACS	52	0	52/96	54%
ChiPSC18	Limiting dilution	46	12	58/55	105%

Expansion of Edited Clonal Lines



Robust Expansion of Edited Clones

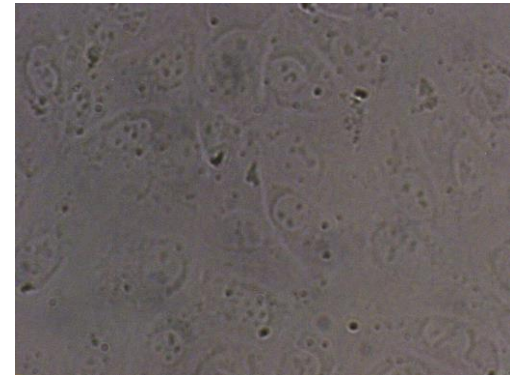
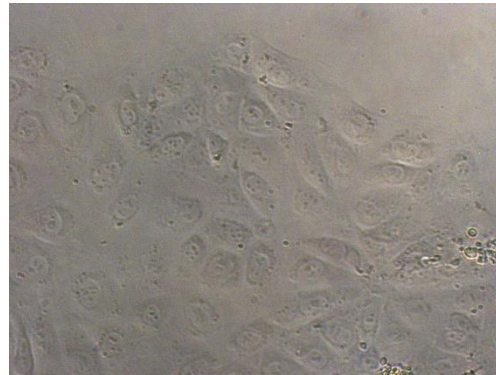
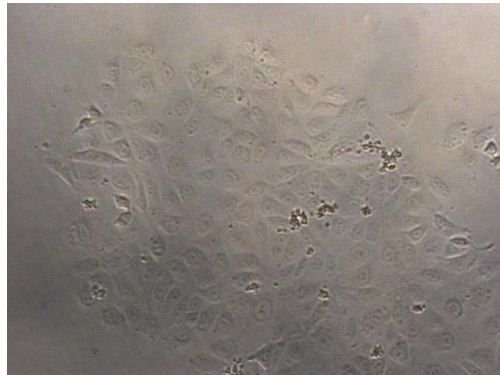
Images of colonies that originated from a single cell

10X

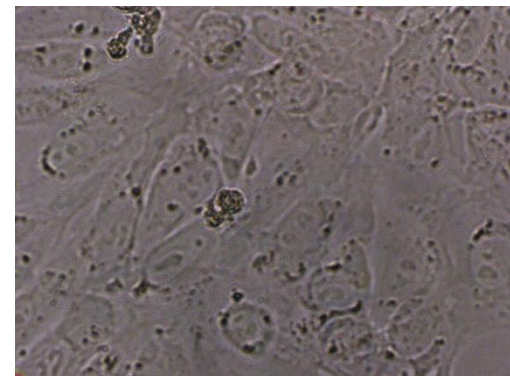
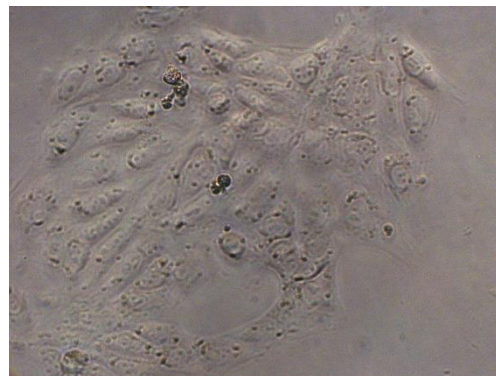
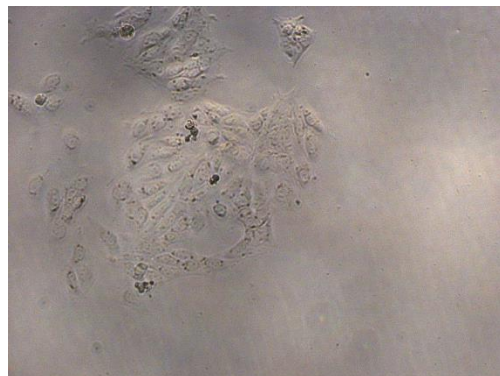
20X

40X

ChiPSC18

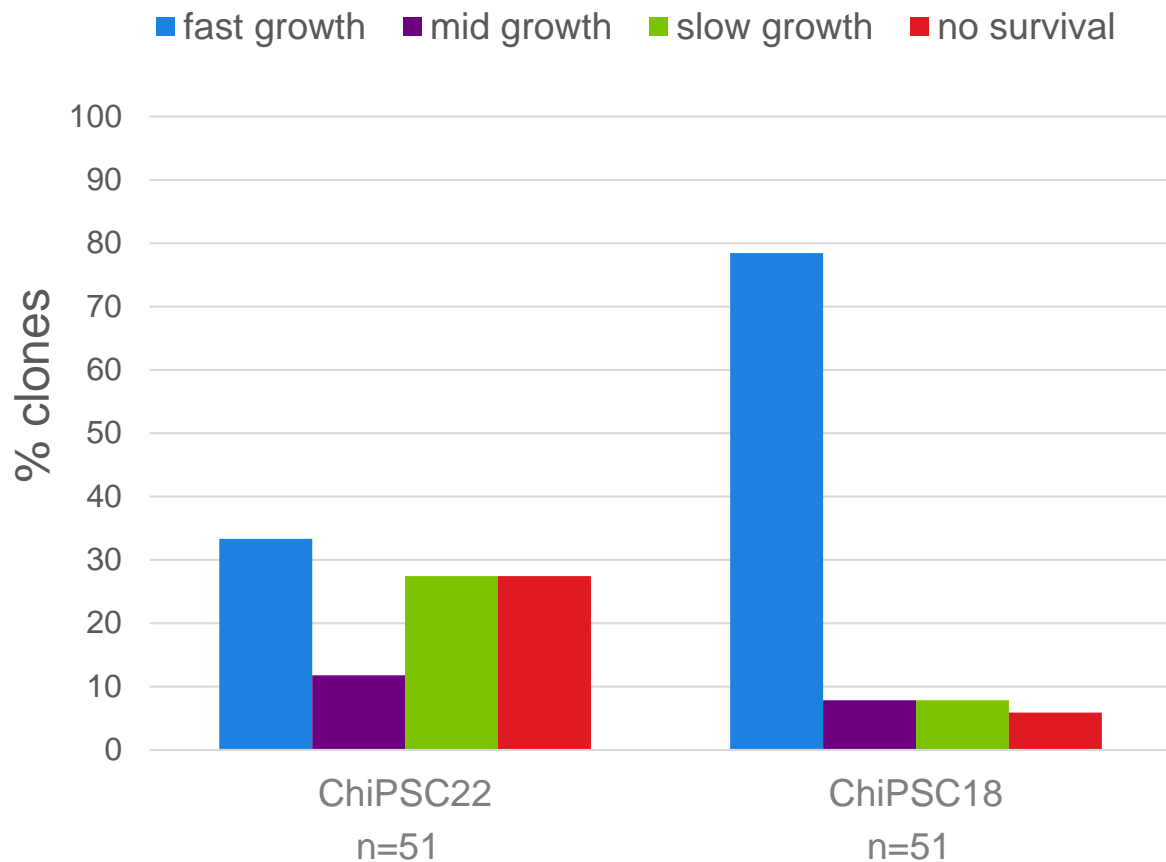


ChiPSC22



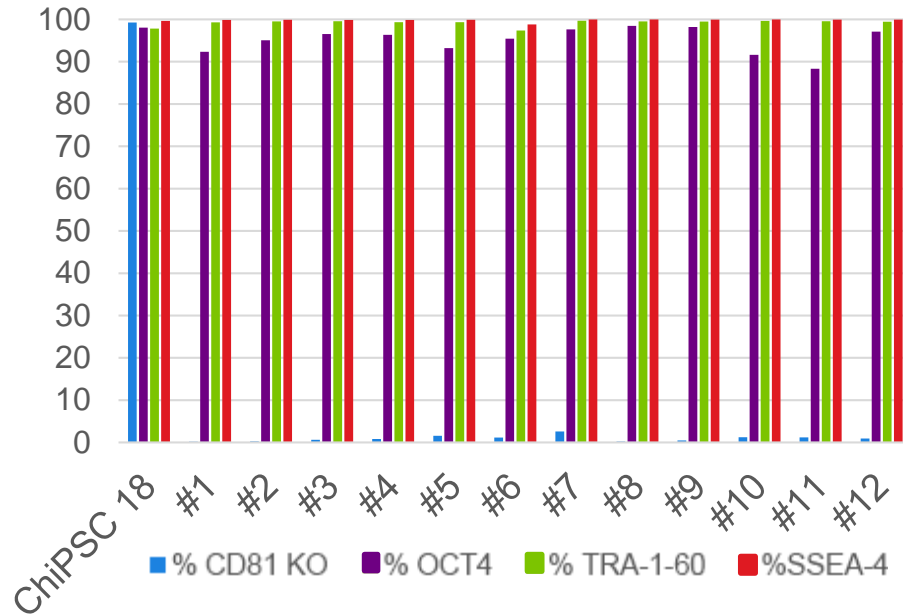
Robust Expansion of Edited Clones

Growth characteristics of clonal lines



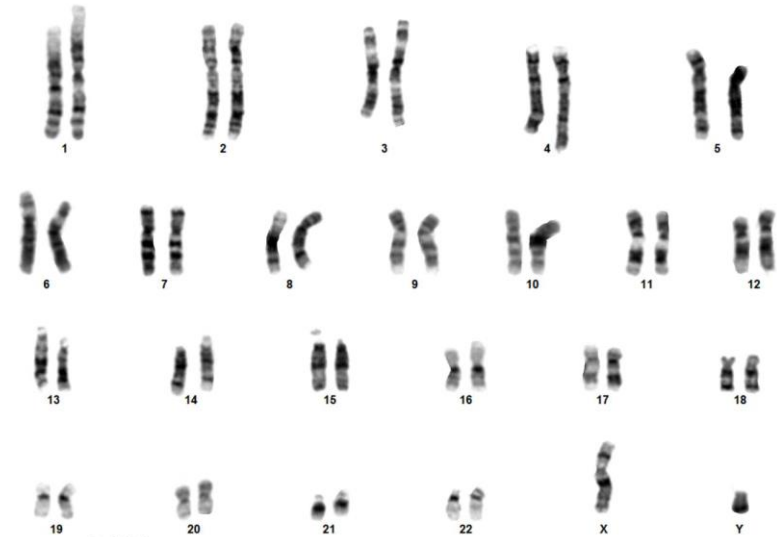
Edited Single-Cell Clones in DEF-CS Maintain Pluripotency & Stable Karyotype

Clone Analysis



Pluripotency maintained following Limiting Dilution

Karyotype Analysis

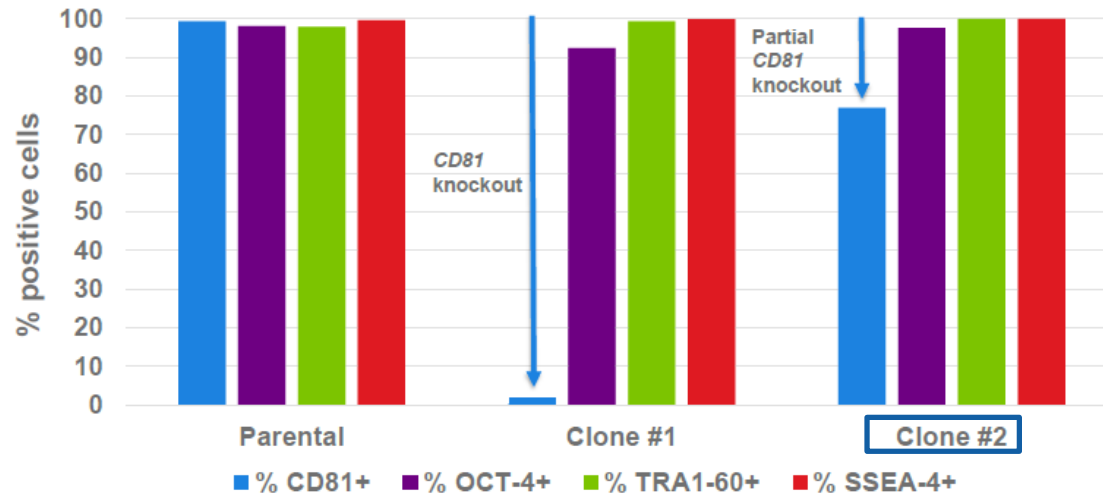


ChiPSC18, c.8, Passage 14

Stable karyotype observed in all edited clones tested after 14 to 15 passages (21 doublings)

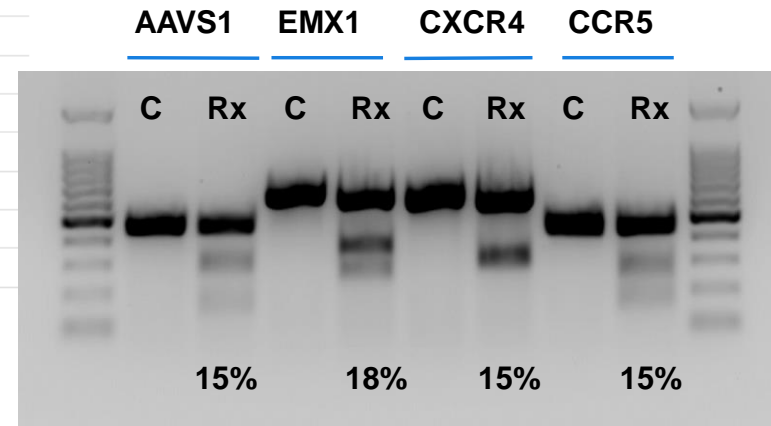
Genomic Characterization of Edited Clones in ChiPSC18

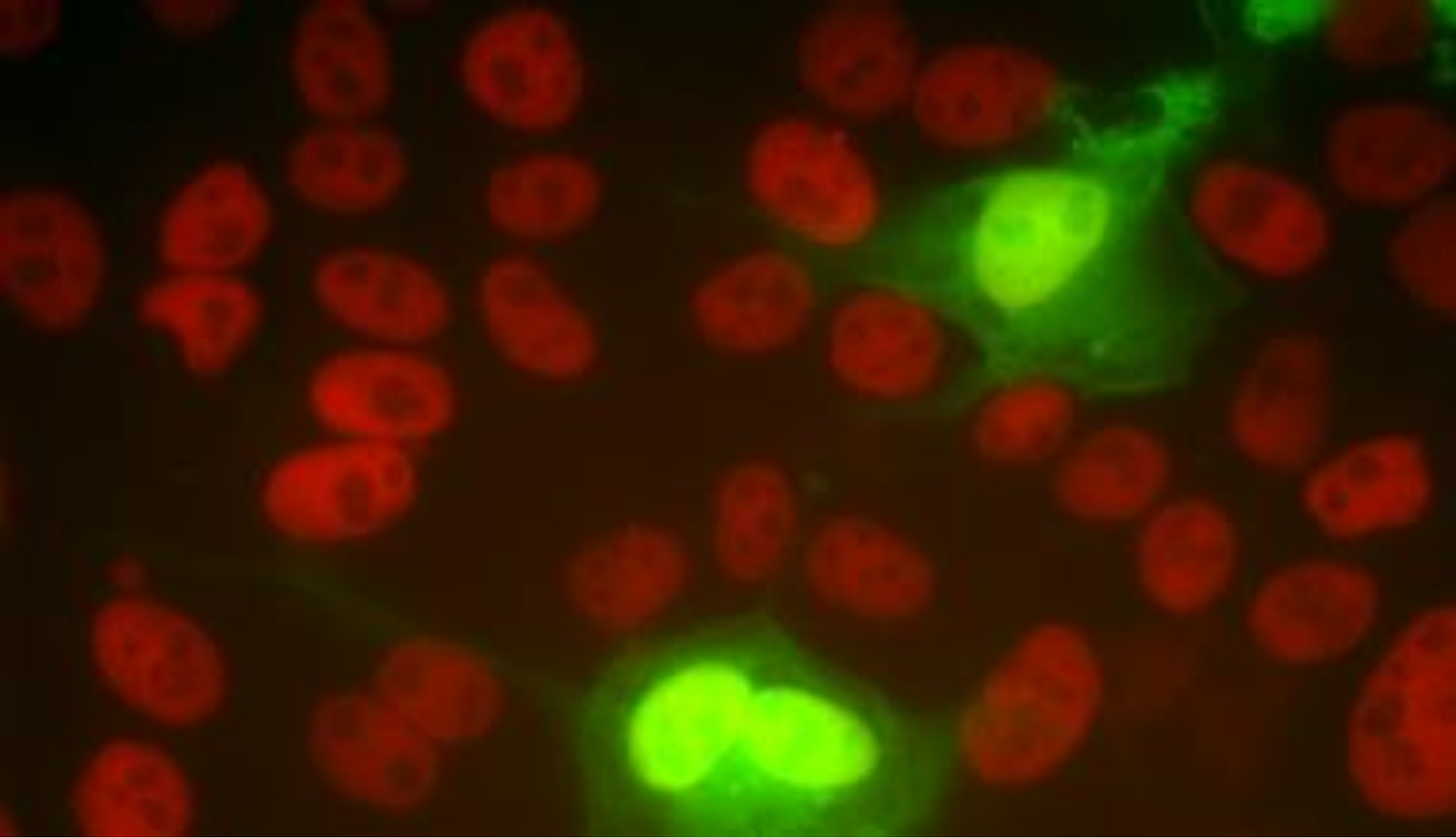
CD81



Other Genes

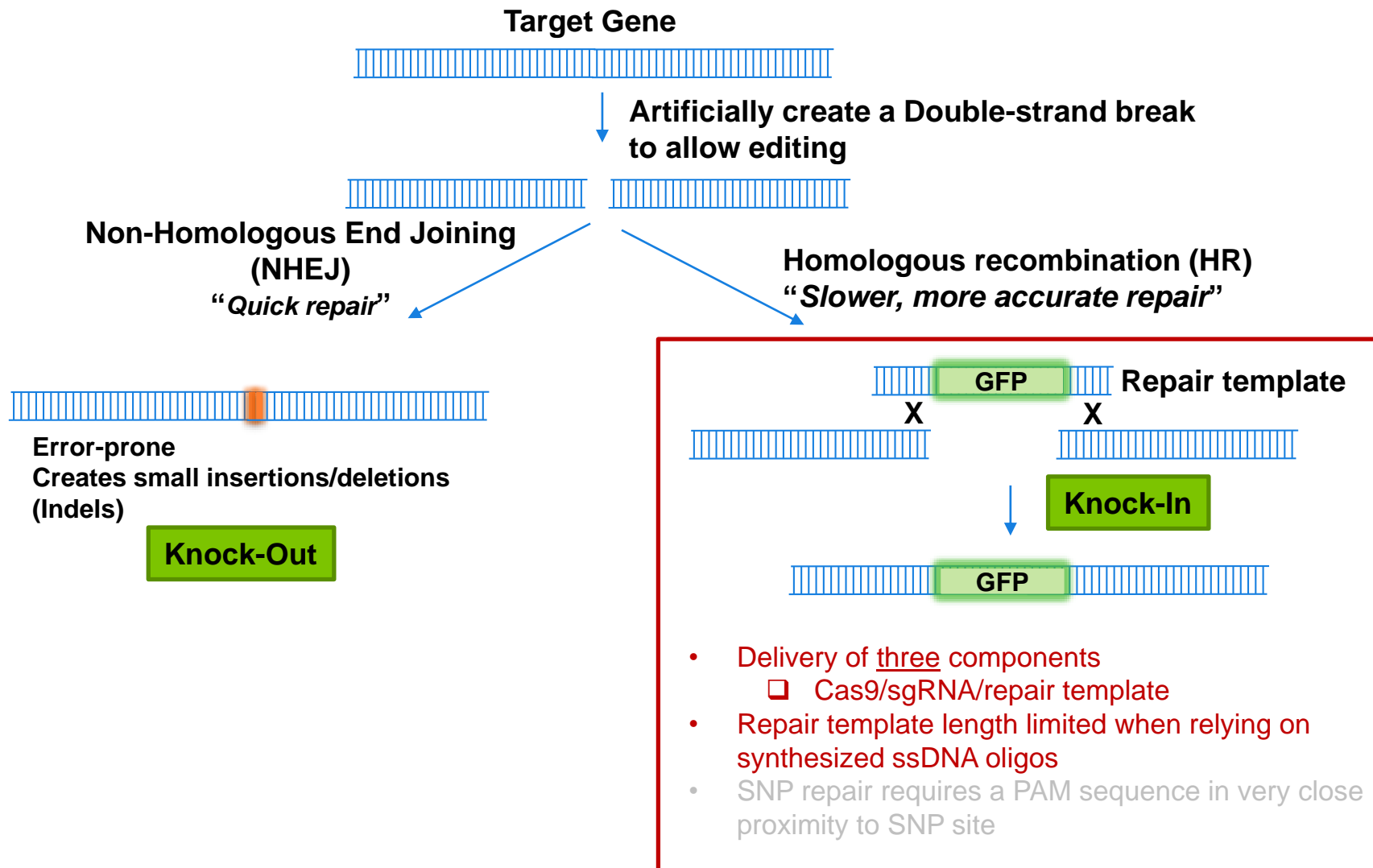
“Mutation detection assay”





Case Study - Knockin

Bottleneck: Homologous Recombination



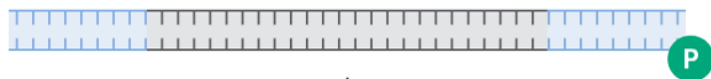
Guide-it Long ssDNA Production System – How it Works

Prepare an HDR template using a method such as cloning or fusion PCR

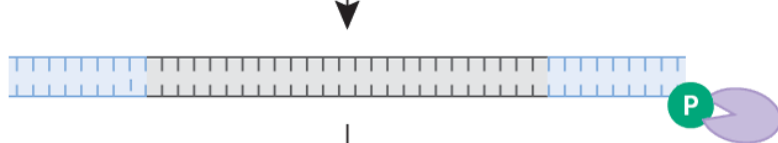


Use PCR to generate starting dsDNA material for the strandase reaction

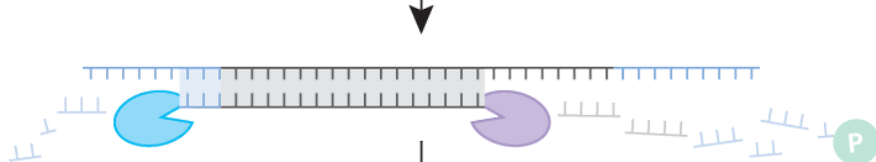
Sense strand generation



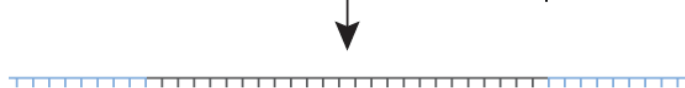
Add Strandase Mix A to begin digesting the phosphorylated strand



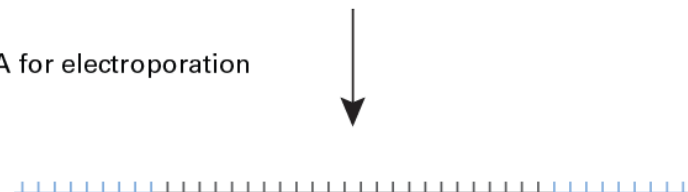
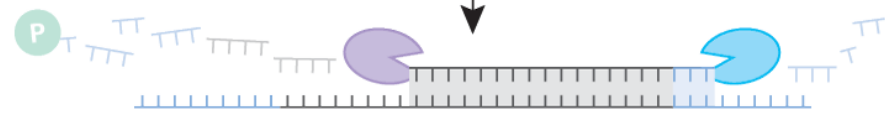
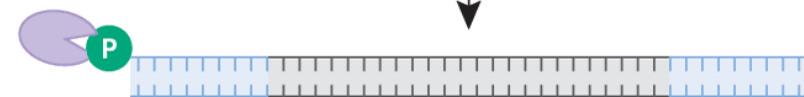
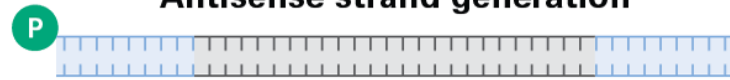
Add Strandase Mix B to finish digesting the strand



Clean up strandase reaction to prepare ssDNA for electroporation

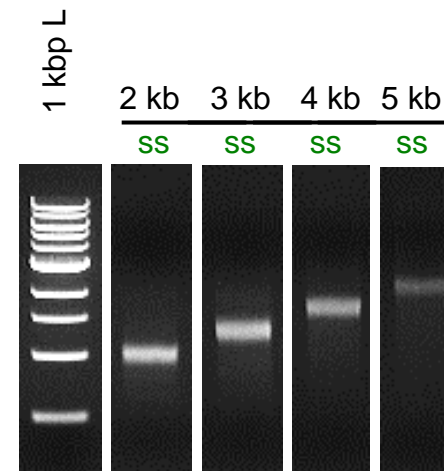
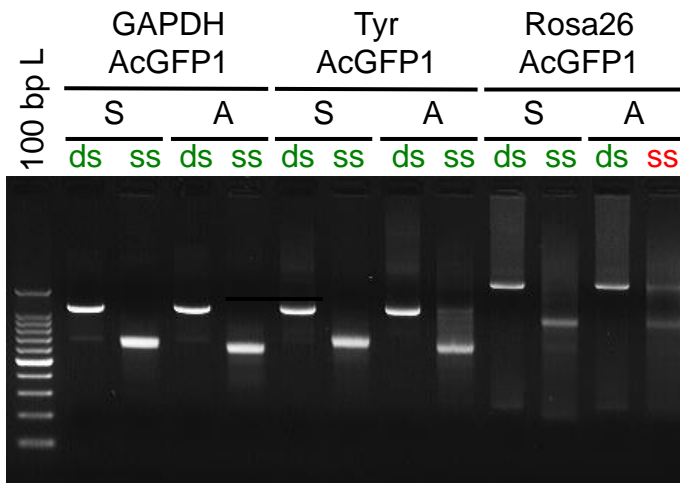


Antisense strand generation



Long ssDNA Prep: *Agarose gel*

* ssDNA is much less sensitive to Et-Br



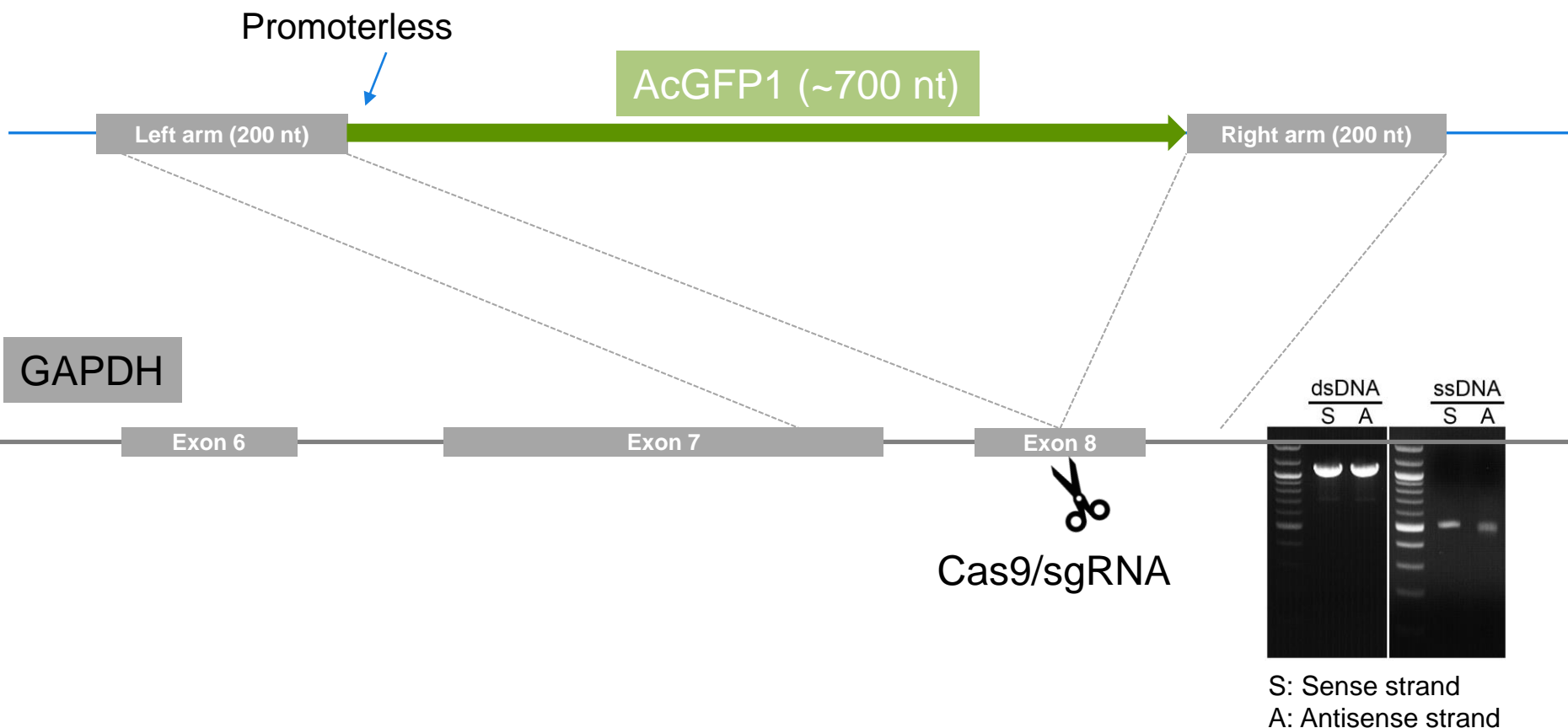
Gene Editing Data

Key things to note:

- dsDNA can result in random integration in the absence of Cas9
- ssDNA integrates seamlessly with a very low error rate at the junctions
- dsDNA will have significant background expression when delivering expression cassettes (e.g. CMV→GFP), making selection for edited cells difficult
- dsDNA can cause significant cellular toxicity

Proof of Concept in HEK293 cells

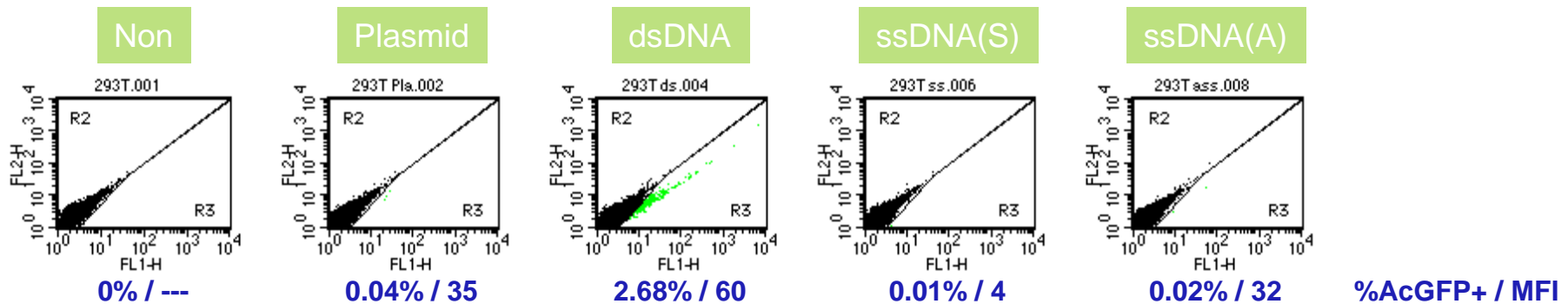
Knock-in of promoterless AcGFP1 at the C-terminus of GAPDH



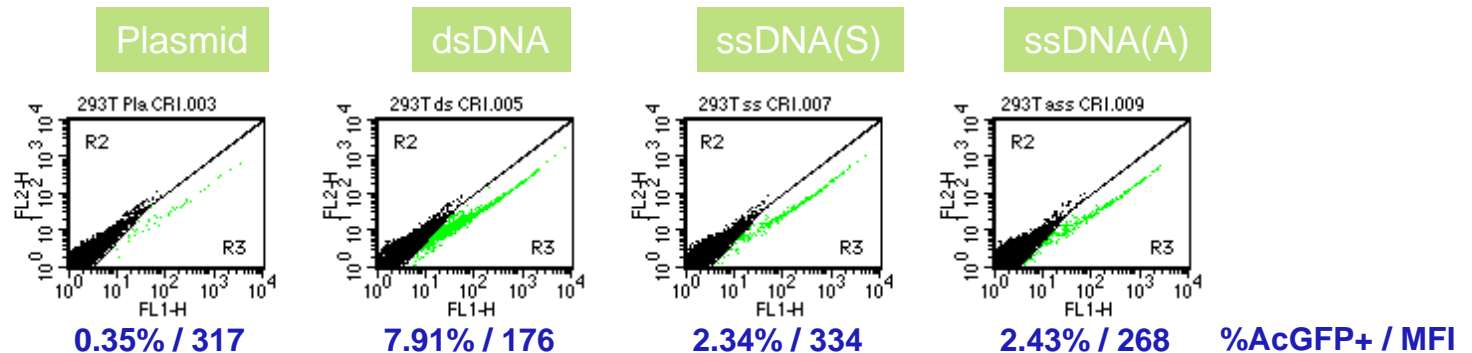
dsDNA Integrates Randomly

HEK293 (GAPDH/AcGFP1 donor)

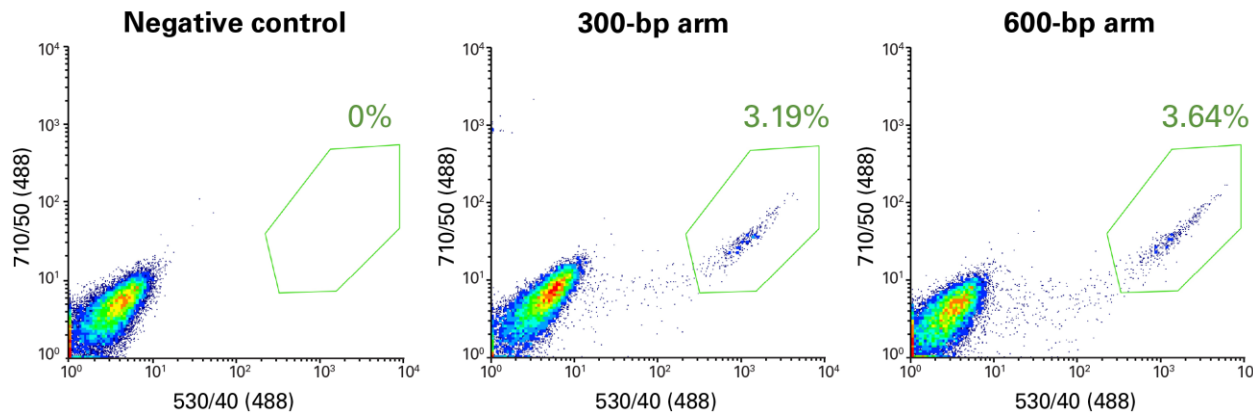
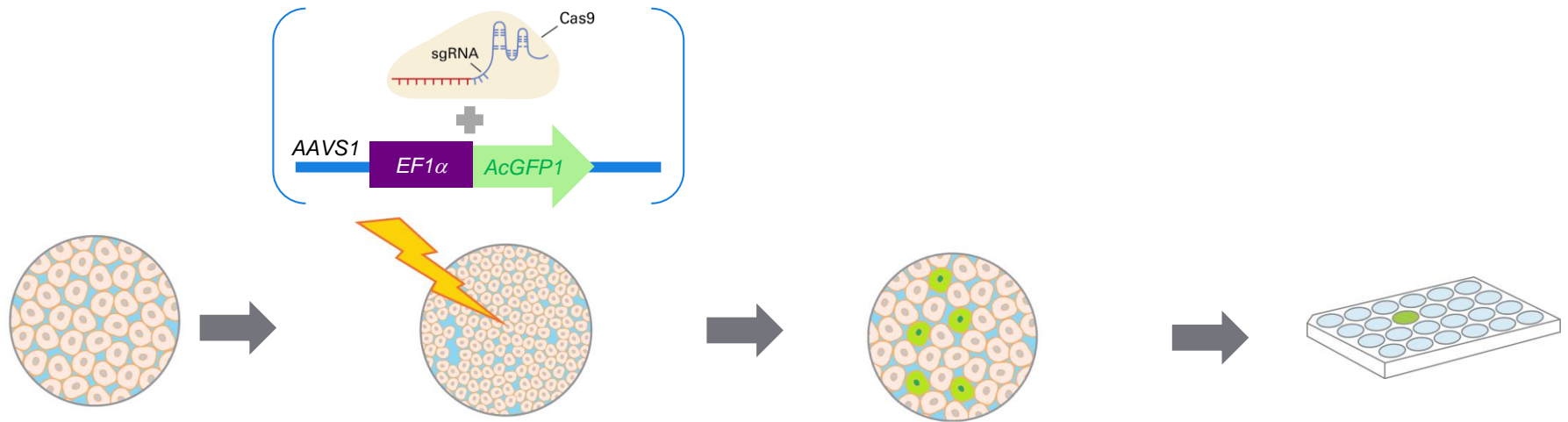
Without Cas9/sgrNA



With Cas9/sgrNA

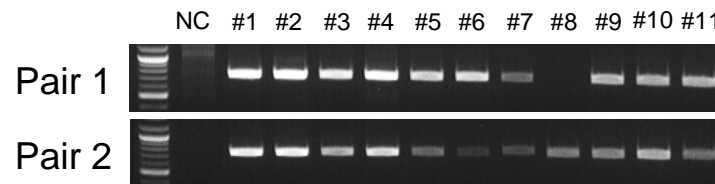
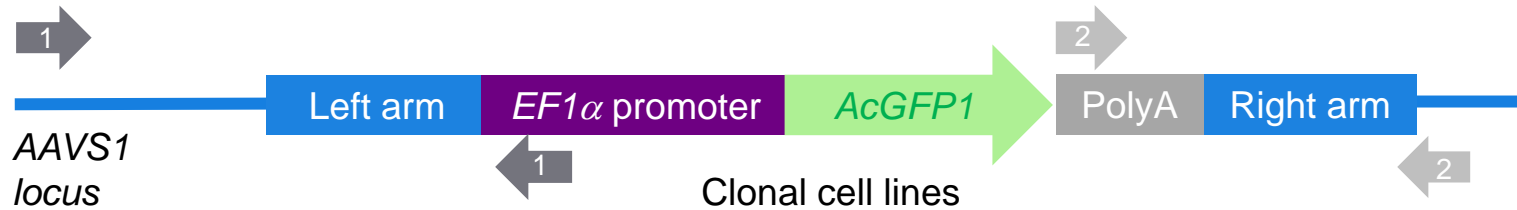


Knock-in of EF1 α -AcGFP1 at AAVS1 Site of hiPS cells (Clone 18)



Isolated AcGFP1 positive single cells

Isolated Clonal Lines Do Not Contain Mutations



Left arm

Right arm

tttccggagcacttccttctcggcgcctgaccacgtagtgctctgagcggatcctcccgtgctctgggtcctctcgg
 TTTCGGAGCAC TTCCTCTCGGCCTGCACCACGTGATGTCTCTGAGCGGATCCTCCCCTGCTG6GTCTCTCCGG
 TTTCGGAGCAC TTCCTCTCGGCCTGCACCACGTGATGTCTCTGAGCGGATCCTCCCCTGCTG6GTCTCTCCGG
 TTTCGGAGCAC TTCCTCTCGGCCTGCACCACGTGATGTCTCTGAGCGGATCCTCCCCTGCTG6GTCTCTCCGG
 TTTCGGAGCAC TTCCTCTCGGCCTGCACCACGTGATGTCTCTGAGCGGATCCTCCCCTGCTG6GTCTCTCCGG
 TTTCGGAGCAC TTCCTCTCGGCCTGCACCACGTGATGTCTCTGAGCGGATCCTCCCCTGCTG6GTCTCTCCGG
 TTTCGGAGCAC TTCCTCTCGGCCTGCACCACGTGATGTCTCTGAGCGGATCCTCCCCTGCTG6GTCTCTCCGG
 TTTCGGAGCAC TTCCTCTCGGCCTGCACCACGTGATGTCTCTGAGCGGATCCTCCCCTGCTG6GTCTCTCCGG
 TTTCGGAGCAC TTCCTCTCGGCCTGCACCACGTGATGTCTCTGAGCGGATCCTCCCCTGCTG6GTCTCTCCGG
 TTTCGGAGCAC TTCCTCTCGGCCTGCACCACGTGATGTCTCTGAGCGGATCCTCCCCTGCTG6GTCTCTCCGG
 TTTCGGAGCAC TTCCTCTCGGCCTGCACCACGTGATGTCTCTGAGCGGATCCTCCCCTGCTG6GTCTCTCCGG
 TTTCGGAGCAC TTCCTCTCGGCCTGCACCACGTGATGTCTCTGAGCGGATCCTCCCCTGCTG6GTCTCTCCGG

taggtgttcaccaggctggtggccgctctactcccttctctttctccatccttcttcccttaaagagtcctccagtgcta
 TAGGTGTTACCAAGTCGTG6CCGCCTCTACTCCCTTCTCTTTCTCCATCCTTCTTCCCTAAAGAGTCCCCAGTGCTA
 TAGGTGTTACCAAGTCGTG6CCGCCTCTACTCCCTTCTCTTTCTCCATCCTTCTTCCCTAAAGAGTCCCCAGTGCTA
 TAGGTGTTACCAAGTCGTG6CCGCCTCTACTCCCTTCTCTTTCTCCATCCTTCTTCCCTAAAGAGTCCCCAGTGCTA
 TAGGTGTTACCAAGTCGTG6CCGCCTCTACTCCCTTCTCTTTCTCCATCCTTCTTCCCTAAAGAGTCCCCAGTGCTA
 TAGGTGTTACCAAGTCGTG6CCGCCTCTACTCCCTTCTCTTTCTCCATCCTTCTTCCCTAAAGAGTCCCCAGTGCTA
 TAGGTGTTACCAAGTCGTG6CCGCCTCTACTCCCTTCTCTTTCTCCATCCTTCTTCCCTAAAGAGTCCCCAGTGCTA
 TAGGTGTTACCAAGTCGTG6CCGCCTCTACTCCCTTCTCTTTCTCCATCCTTCTTCCCTAAAGAGTCCCCAGTGCTA
 TAGGTGTTACCAAGTCGTG6CCGCCTCTACTCCCTTCTCTTTCTCCATCCTTCTTCCCTAAAGAGTCCCCAGTGCTA
 TAGGTGTTACCAAGTCGTG6CCGCCTCTACTCCCTTCTCTTTCTCCATCCTTCTTCCCTAAAGAGTCCCCAGTGCTA
 TAGGTGTTACCAAGTCGTG6CCGCCTCTACTCCCTTCTCTTTCTCCATCCTTCTTCCCTAAAGAGTCCCCAGTGCTA

Clonal
cell
lines

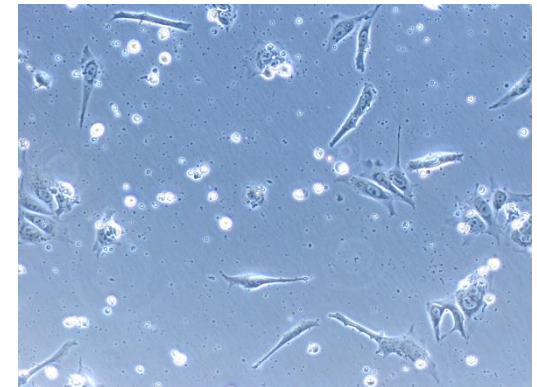
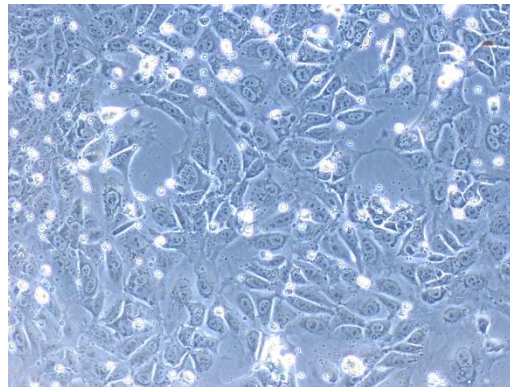
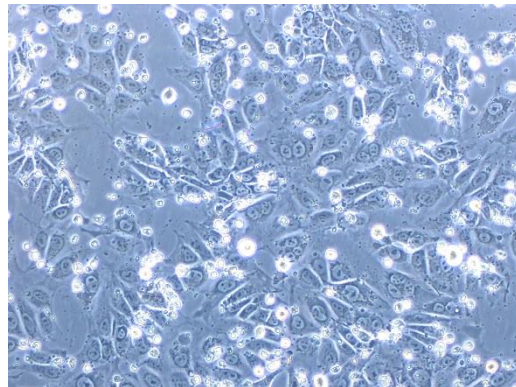
dsDNA-induced toxicity in ChiPSC18

No donor template

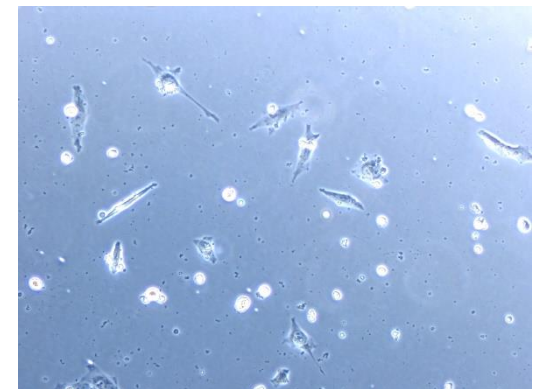
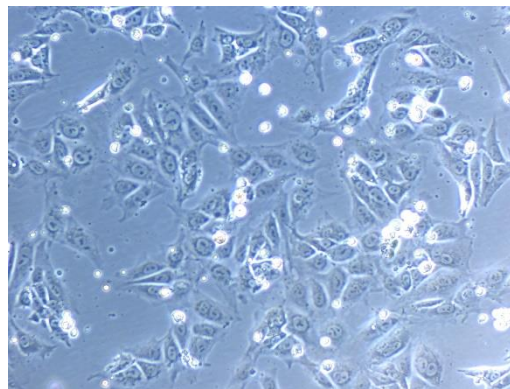
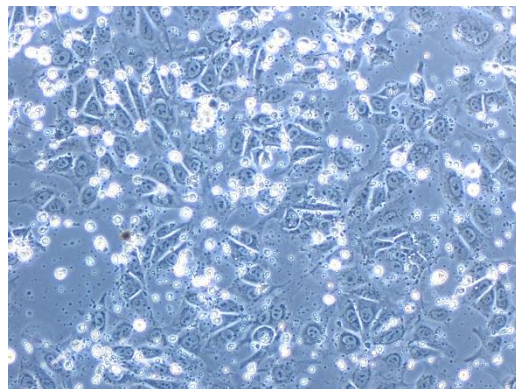
ssDNA

dsDNA

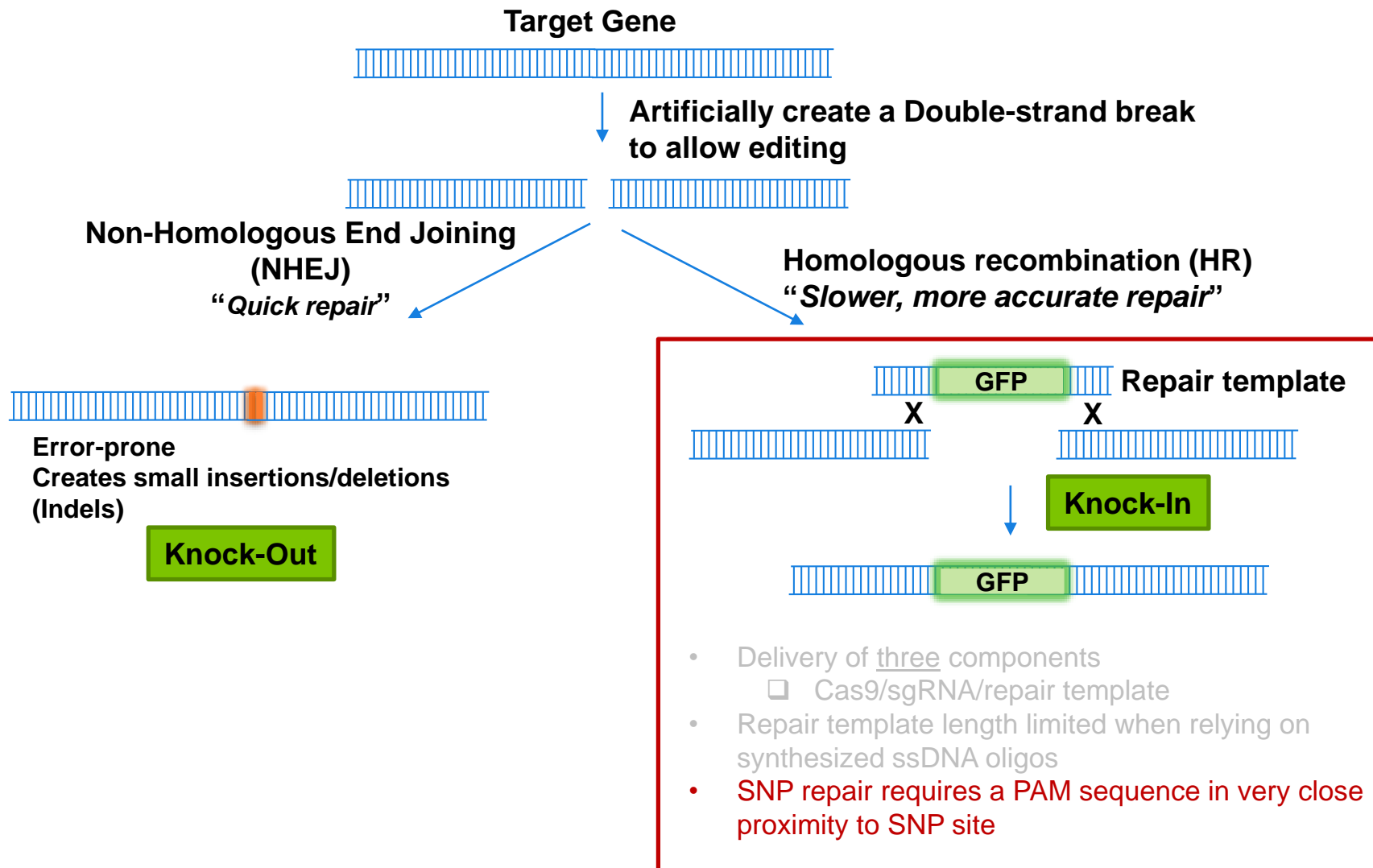
-Cas9
+sgRNA



+Cas9
+sgRNA



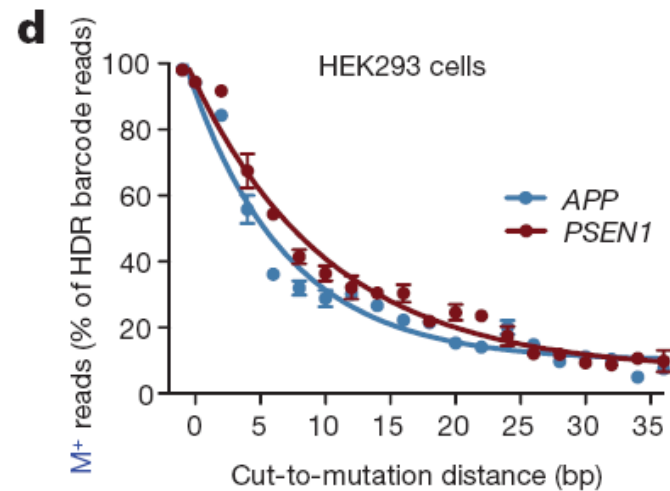
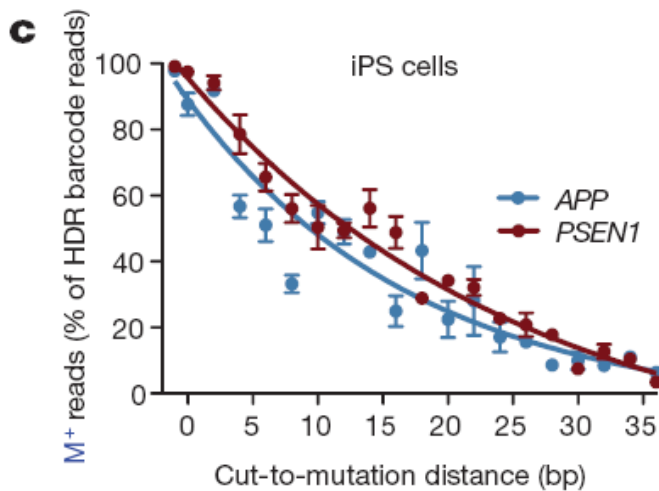
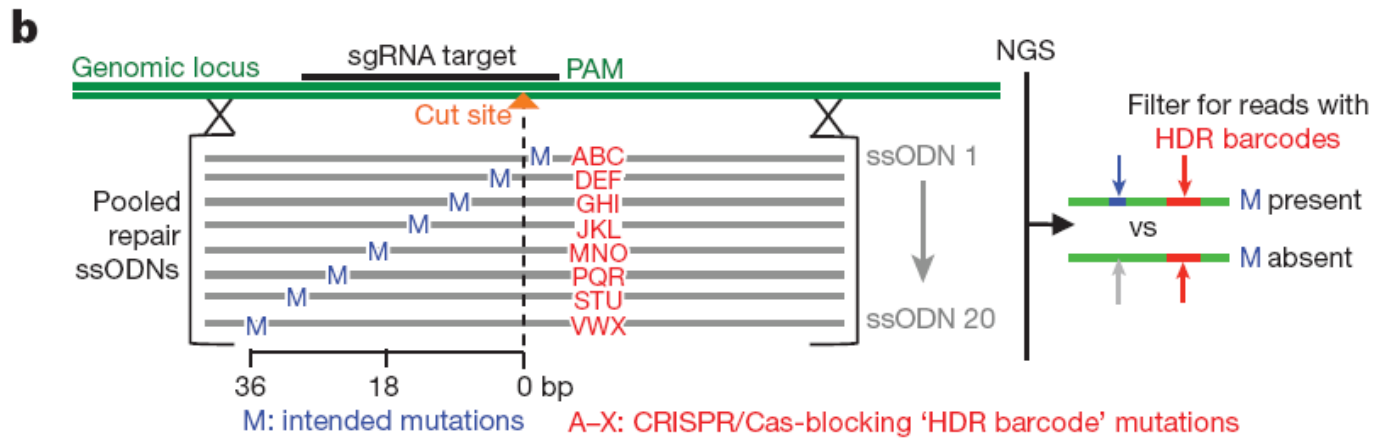
Bottleneck: Homologous Recombination



Why is SNP Editing/Screening Important?

- Majority of genetic diseases are “caused” by (or can be correlated to) SNPs
- 46,589 SNPs and 6,356 diseases and phenotypes (source: <http://www.disgenet.org/>)
- Essential for disease model development
- “Personalized medicine” by predicting efficacy of a new drug dependent on a SNP “fingerprint”
- Identifying how a complex set of different SNPs can cause similar disease phenotypes

Efficiency of SNP Repair Relies on Close Proximity to PAM Site



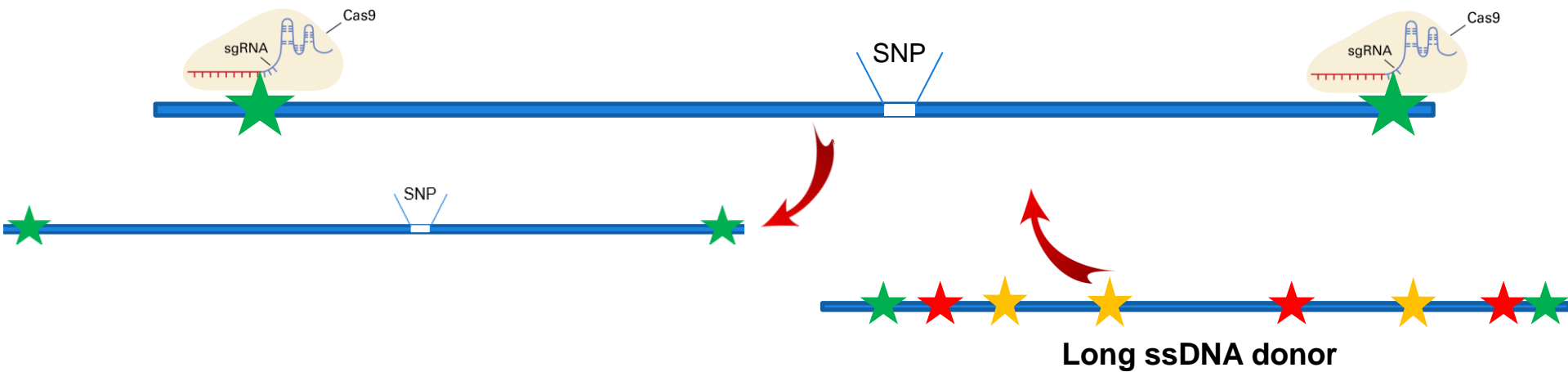
Nature 533, 125–129 (05 May 2016)

Advantage of Using Long ssDNA for SNP Repair via Homologous Recombination

SNP allele



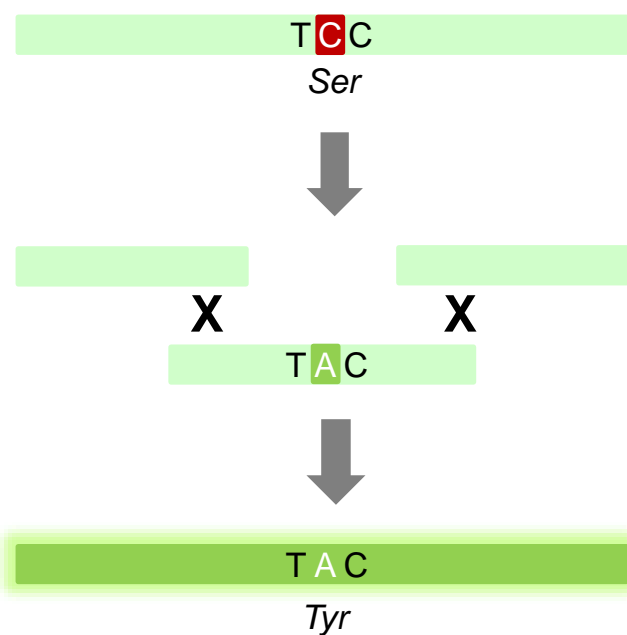
- ★ lousy CRISPR/Cas9 PAM/cut site
- ★ mediocre CRISPR/Cas9 PAM/cut site
- ★ great CRISPR/Cas9 PAM/cut site



Repaired allele



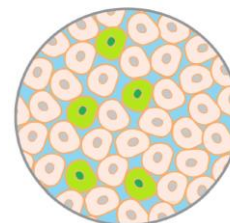
Homology-Directed Knockin of Point Mutations in hiPSCs



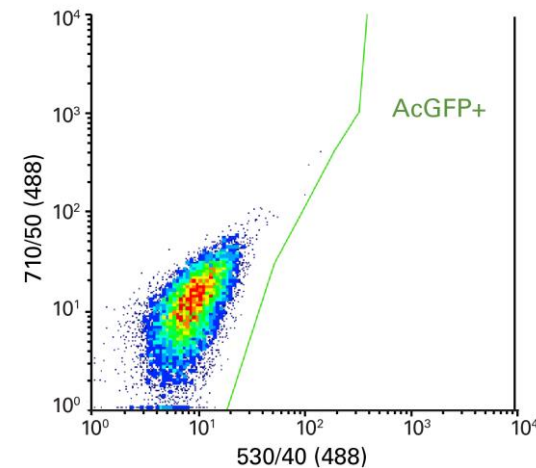
Non-fluorescent mutant AcGFP1



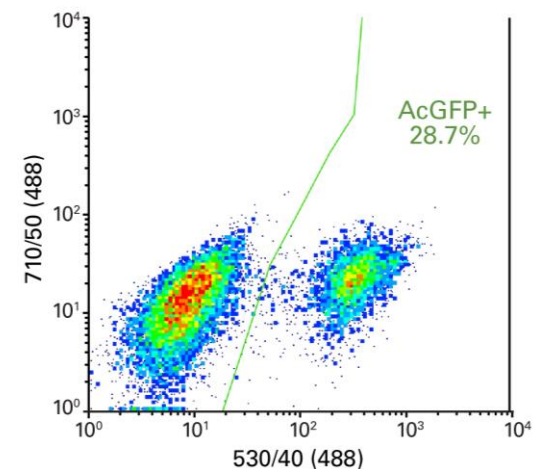
Fluorescent AcGFP1



Negative control



Cas9/sgRNA + 1 µg ssDNA oligo



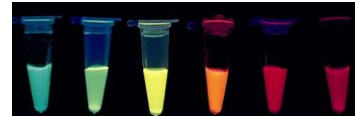
Cell Biology Portfolio Enabling Gene Function Analysis

Cellartis:
-Products
-Service business

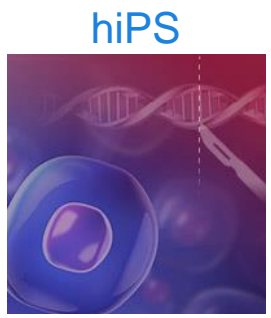
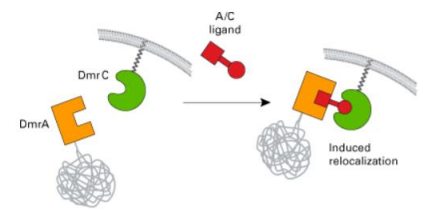
Genome editing



Fluorescent proteins and reporters

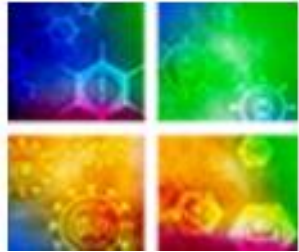


Ligand induced protein interaction

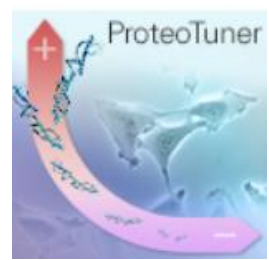


Gene Function

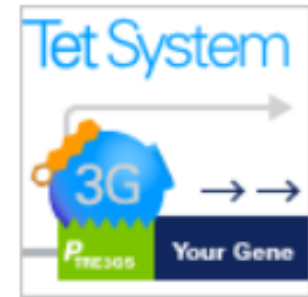
Viral delivery



Ligand induced protein stabilization

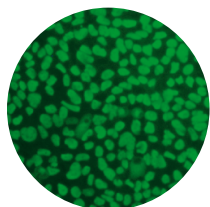


Inducible gene expression systems



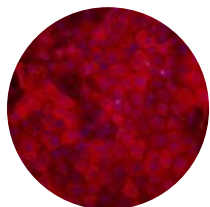
Stem Cell Research

Stem cell innovations for today and the future



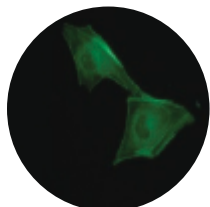
Cell Culture

- Expansion & maintenance
- Research-grade, Xeno-Free & GMP
- Differentiation



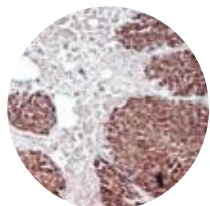
ES/iPS-derived Cells

- Hepatocytes & cardiomyocytes
- DE cells
- Beta cells
- Neural progenitors



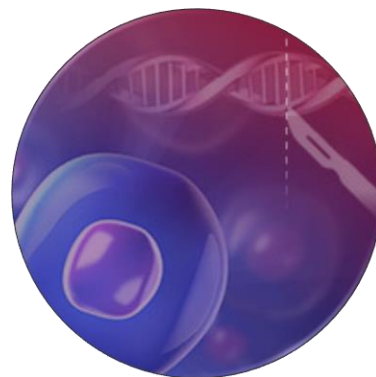
Stem Cell Services

- Sourcing & reprogramming
- Clinical-grade cell line generation & banking
- Genome modification



Characterization and Detection

- Antibodies to verify pluripotency, differentiation, etc.
- qPCR primer sets



New products: Human iPS Gene Editing Systems

- Optimized delivery for clonal expansion
- Flexible gene methods



that's
GOOD
science!®

Clontech **TAKARA** cellartis