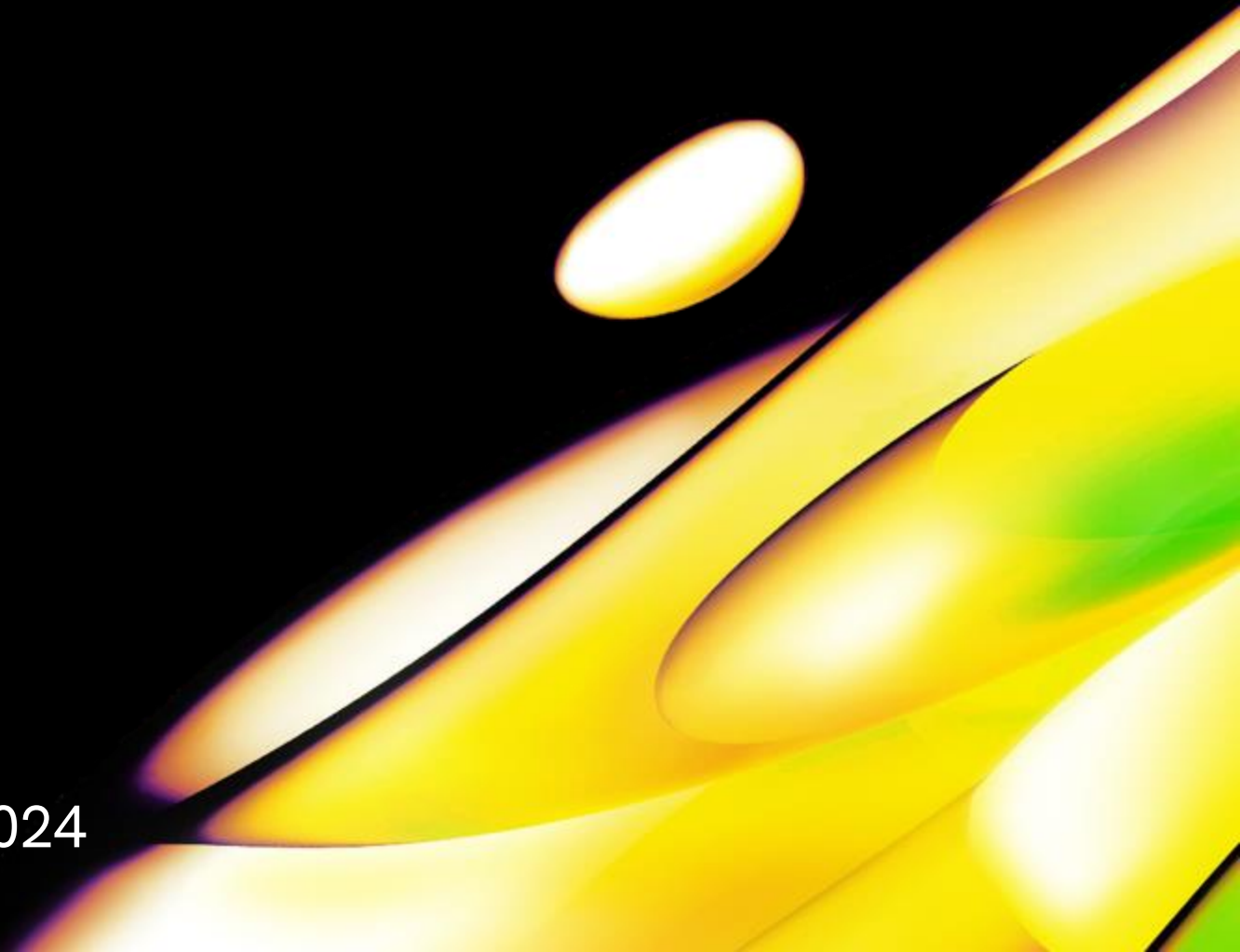


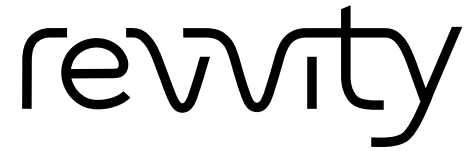
# The Pin-point™ platform

A novel modular base editing system

revvity

Updated 16 July 2024





## Outline

1. [Introduction](#)
2. [Data Overview](#)
3. [Summary](#)

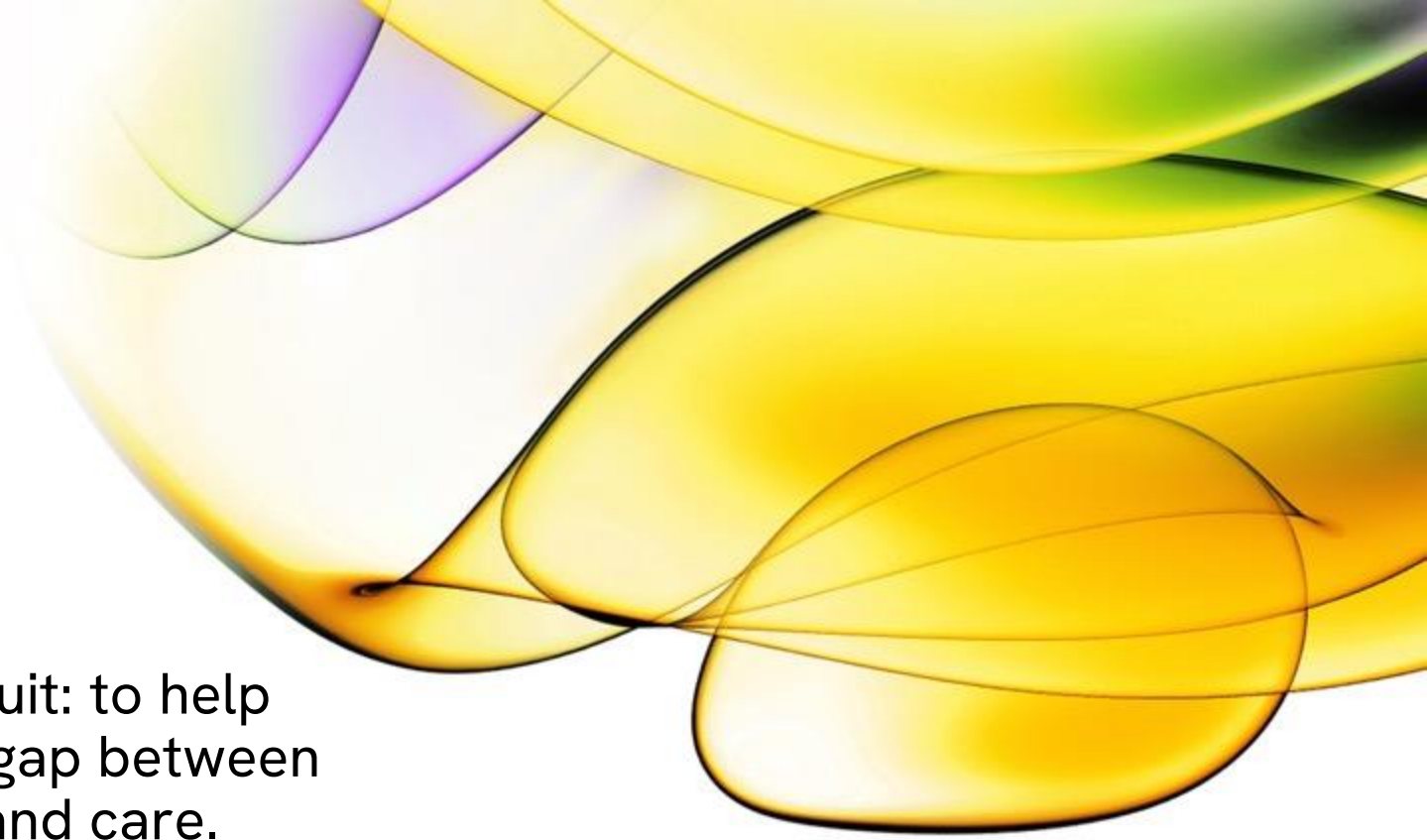
## Supplemental Sections

1. [Knockout explained](#)
2. [T cell datapack](#)
3. [iPSC datapack](#)
4. [HSPC datapack](#)
5. [Complex Engineering](#)
6. [Flexible Modularity](#)
7. [Screening - Arrayed & Pooled](#)

# revvity

Revvity is born of a single-minded pursuit: to help improve human health by bridging the gap between science and people through precision and care.

We innovate and collaborate to empower our partners to see science in unexpected ways that deliver breakthrough results.





Targeted knock-in

Programmable

Dividing, non-dividing,  
& primary cells

Efficient

Flexible

Predictable

Precise

Modular

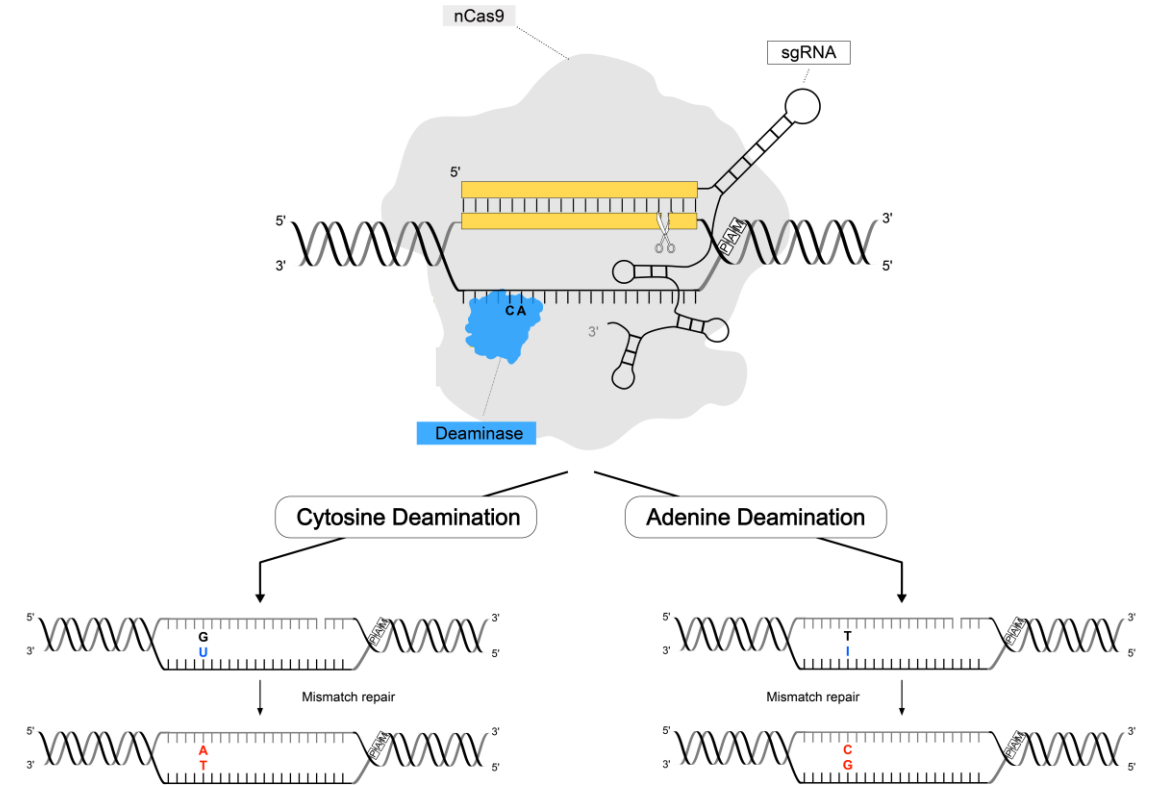
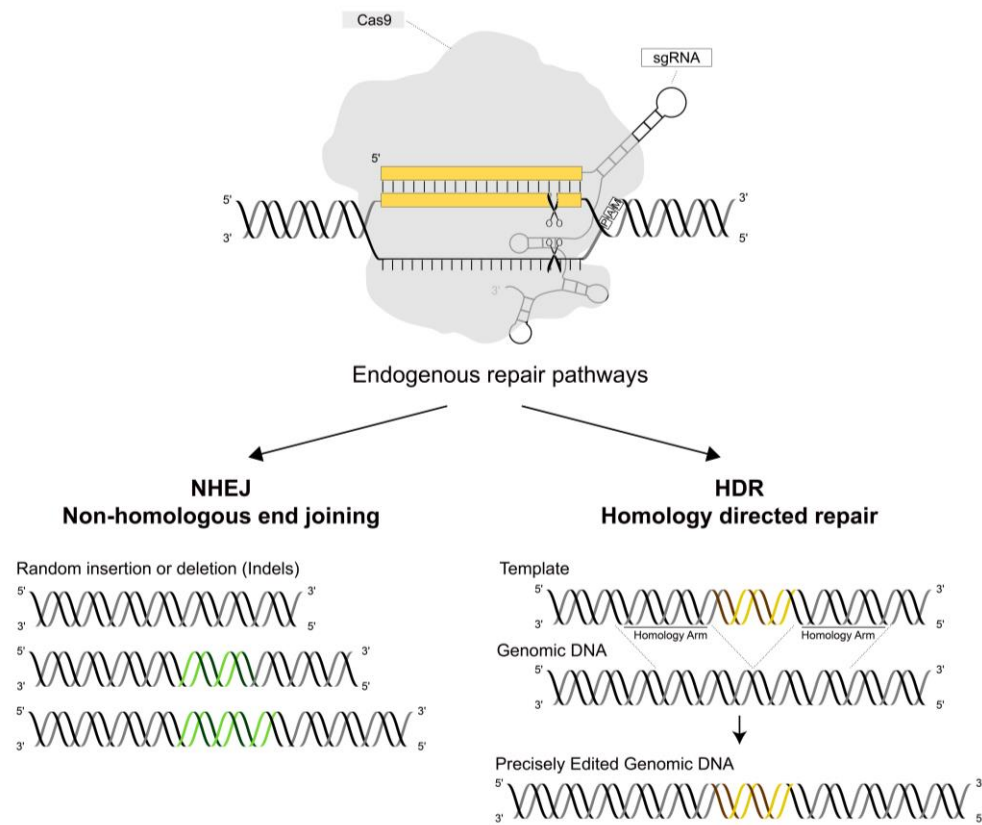
Multiplexing

Safe

Avoid DSBs

# CRISPR gene editing

# Base editing



## GENE DISRUPTION BY A DsDNA BREAK

- Indel formation to disrupt gene sequence
- complex population of indels

## GENE MODIFICATION BY POINT MUTATIONS

- Creation of stop codons or splice site disruption for knockout
- Introduction of single base conversion

# Competitive advantages of base editing

## 1<sup>st</sup> generation Cas enzymes

Gene disruption by a dsDNA break

## 2<sup>nd</sup> generation base editing

Gene modification by point mutation

- ✓ Avoidance of double strand DNA breaks for **reduced cytotoxicity** and **high viability**
- ✓ **Permanent change** to the DNA
- ✓ Flexible creation of stop codons or splice site disruption for **knockout** or introduction of **single base conversion**
- ✓ **Predictable, precise, and efficient** single or multi-gene editing
- ✓ Components are easy to **design, synthesize, and deliver**

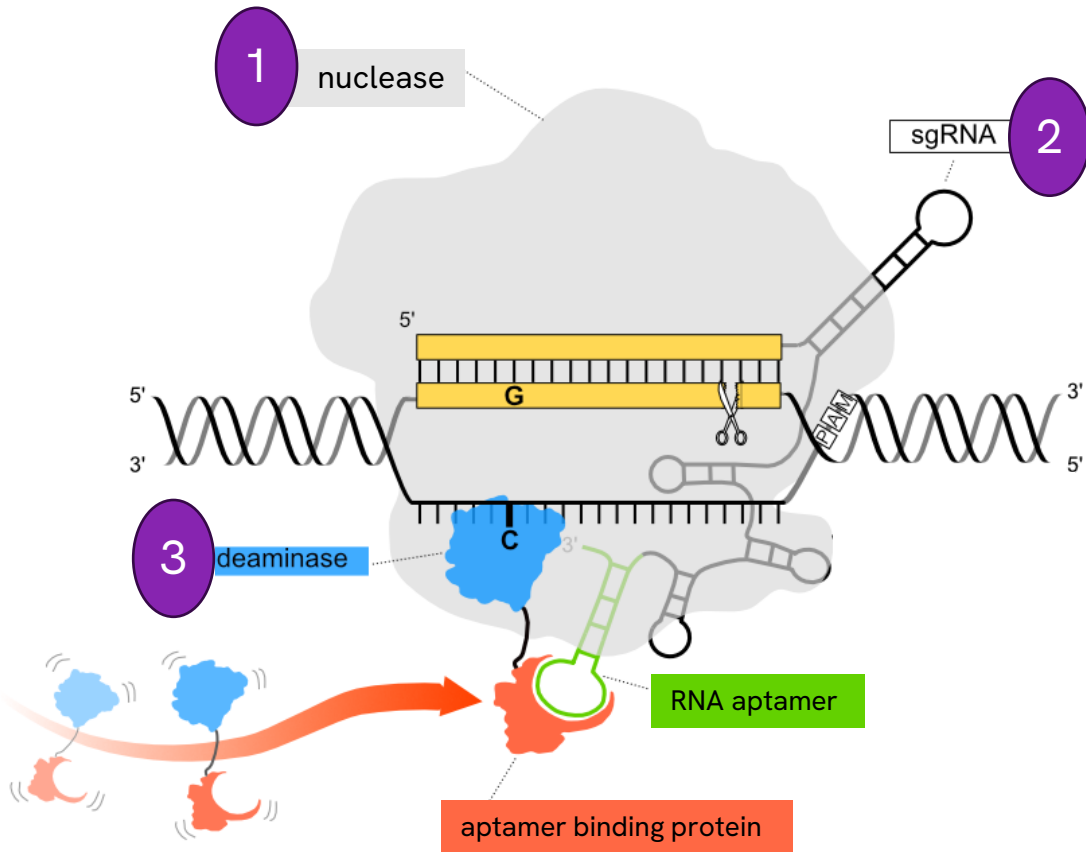


## New generation Pin-point™ base editing platform

- ✓ Simultaneous knock-in and knockout in a single reaction
- ✓ Nuclease and deaminase flexible
- ✓ Modular control over target and editing window to specifically reach your gene of interest
- ✓ A single novel, patented aptamer-recruited base editing platform for your therapeutic development

# What is the Pin-point™ platform?

Based on a patented aptamer-recruited base editing arrangement



\*Schematic depicts nCas9 configuration

## 3 component system

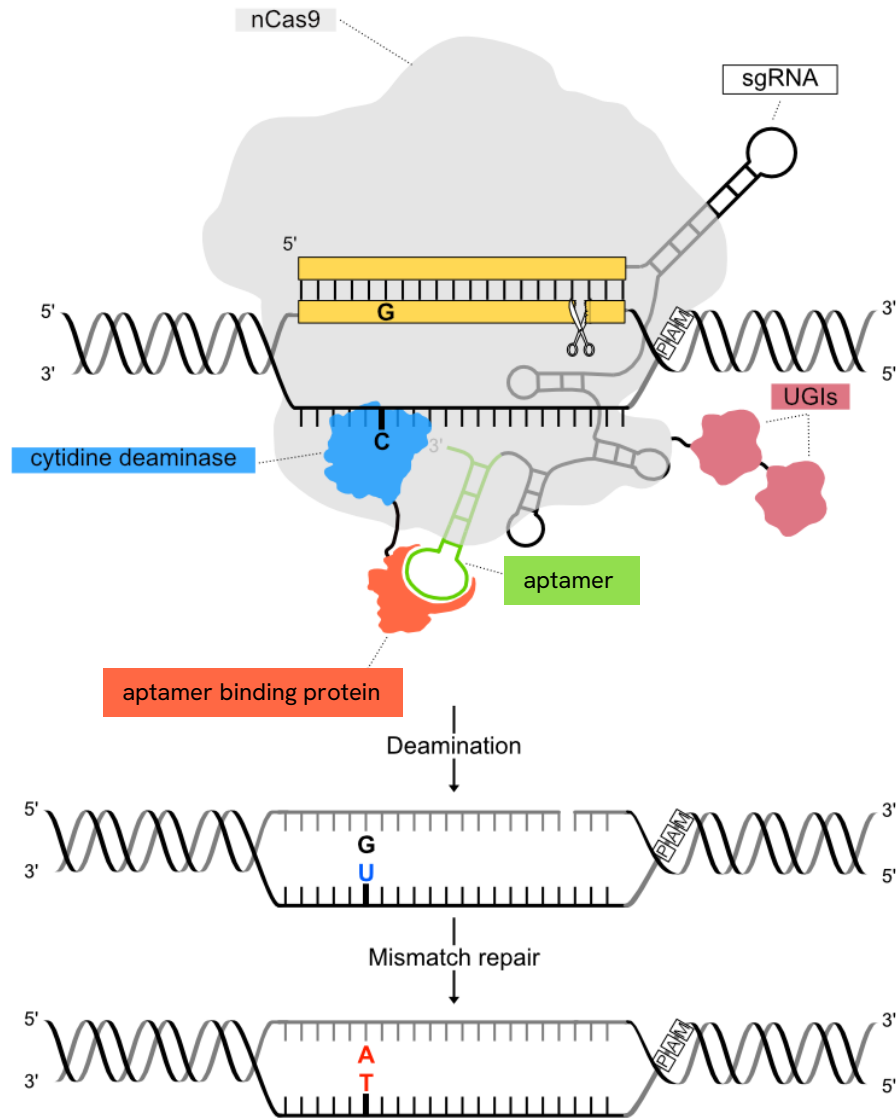
1. RNA-guided enzyme
2. Guide RNA with aptamer
3. Deaminase and recruitment protein

## Demonstrated advantages

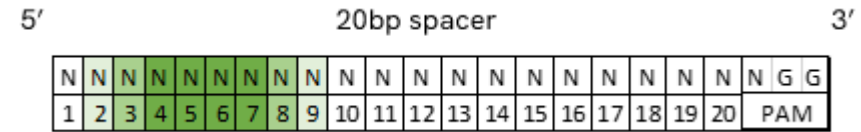
- ✓ Multiplex gene editing including knock-in and knockout with high efficiency and safety
- ✓ Validated performance in T cells, iPSCs, and HSPCs
- ✓ Mix-and-match for target specificity and efficiency



# Base editing terminology



## "Base editing window"



*nCas9/rat APOBEC is most likely to edit C's in positions 4-7*

"Bystander editing" is any editing other than the target base of interest

"Off-target editing" is any editing other than at the locus that is targeted

# The Pin-point™ base editing platform is accelerating therapeutic development research

Molecular Therapy  
Original Article



## An aptamer-mediated base editing platform for simultaneous knockin and multiple gene knockout for allogeneic CAR-T cells generation

Immacolata Porreca,<sup>1</sup> Robert Blassberg,<sup>1</sup> Jennifer Harbottle,<sup>1,7</sup> Bronwyn Joubert,<sup>1</sup> Olga Mielczarek,<sup>1</sup> Jesse Stombaugh,<sup>2</sup> Kevin Hemphill,<sup>2</sup> Jonathan Sumner,<sup>2</sup> Devidas Pazeraitis,<sup>3</sup> Julia Liz Touza,<sup>4</sup> Margherita Francescato,<sup>4</sup> Mike Firth,<sup>2</sup> Tommaso Selmi,<sup>1,9</sup> Juan Carlos Collantes,<sup>5</sup> Zaklina Strezoska,<sup>2</sup> Benjamin Taylor,<sup>3</sup> Shengkan Jin,<sup>6</sup> Ceri M. Wiggins,<sup>1,8</sup> Anja van Brabant Smith,<sup>2</sup> and John J. Lambourne<sup>1,10</sup>

<sup>1</sup>Revvity, 8100 Cambridge Research Park, Cambridge CB25 9TL, UK; <sup>2</sup>Revvity, 2650 Crescent Drive, Lafayette, CO 80026, USA; <sup>3</sup>AstraZeneca, Discovery Sciences, R&D, 1 Francis Crick Avenue, Cambridge Biomedical Campus, Cambridge CB2 0AA, UK; <sup>4</sup>AstraZeneca, Discovery Sciences, BioPharmaceuticals R&D Unit, AstraZeneca, Pepparedsleden 1, 431 83 Mölndal, Sweden; <sup>5</sup>Departamento de Biotecnología, Colegio de Ciencias Biológicas y Ambientales, Universidad San Francisco de Quito, Campus Cumbayá, Casilla Postal 17-1200-841, Quito 170901, Ecuador; <sup>6</sup>Pharmacology Department, Rutgers, The State University of New Jersey, Robert Wood Johnson Medical School, 675 Hoes Lane West, Piscataway, NJ 08854, USA

Gene editing technologies hold promise for enabling the next generation of adoptive cellular therapies. In conventional gene editing platforms that rely on nuclease activity, such as clustered regularly interspaced short palindromic repeats CRISPR-associated protein 9 (CRISPR-Cas9), allow efficient introduction of genetic modifications; however, these modifications occur via the generation of DNA double-strand breaks (DSBs) and can lead to unwanted genomic alterations and genotoxicity. Here, we apply a novel modular RNA aptamer-mediated Pin-point base editing platform to simultaneously introduce multiple gene knockouts and site-specific integration of a transgene in human primary T cells. We demonstrate high editing efficiency and purity at all target sites and significantly reduced frequency of chromosomal translocations compared with the conventional CRISPR-Cas9 system. Site-specific knockin of a chimeric antigen receptor and multiplex gene knockout are achieved within a single intervention and without the requirement for additional sequence-targeting components. The ability to perform complex genome editing efficiently and precisely highlights the potential of the Pin-point platform for application in a range of advanced cell therapies.

cell product. To expand the scope of these innovative off-the-shelf therapies to solid tumors, further edits will also be required to ensure therapeutic cells retain their efficacy in the refractory and heterogeneous tumor microenvironment.<sup>8</sup> These factors, together with the need to provide new functions to the cells to make effective and safe therapies that offer wider patient accessibility and therapy deployment, ultimately demand increasingly refined genome editing strategies.

Gene editing technologies such as zinc-finger nucleases, transcription activator-like effector nucleases and CRISPR-Cas9 have all been employed to successfully perform targeted editing at genomic loci for effective knockout and knockin applications. However, the generation of double-strand breaks (DSBs) inherent to their mechanism of conferring a DNA edit can lead to chromosomal loss or structural variation.<sup>9-11</sup> The occurrence of chromosomal aberrations is enhanced in the context of multi-gene editing as more concurrent DSBs are generated, and the extent of this damage is expanded if DNA breaks also occur at off-target sites. Although many structural aberrations in a cell may not be viable, it has been reported that

### INTRODUCTION

Gene editing technologies have entered the clinic and show significant potential for advancing next-generation therapies, particularly in the development of more efficient chimeric antigen receptor (CAR)-T cell therapies to address hematological malignancies.<sup>1-3</sup> To overcome the logistical and infrastructure-related challenges and product variability barriers of the autologous cell therapy paradigm, recent focus has shifted to realizing the potential of allogeneic cell therapies. The manufacture of allogeneic cell products requires multiple edits to prevent both graft-versus-host disease and immune rejection by the host, which would otherwise limit efficacy, persistence, and safety of the

Received 21 June 2023; accepted 24 June 2024;  
<https://doi.org/10.1016/j.jmthe.2024.06.033>.

<sup>1</sup>Present address: AstraZeneca, Safety Sciences, Cell Therapy Oncology team, R&D, CB2 0QQ Cambridge, UK

<sup>2</sup>Present address: AstraZeneca, Discovery Sciences, R&D, 1 Francis Crick Avenue, Cambridge Biomedical Campus, CB2 0AA Cambridge, UK

<sup>3</sup>Present address: Consiglio Nazionale delle Ricerche, Istituto di Tecnologia Biomedica, Via Fratelli Cervi 93, 20054 Segrate (MI), Italy

<sup>4</sup>Present address: Penick Biosciences Ltd, 21G1, Alderley Park, Macclesfield, Cheshire SK11 9ATG, UK

Correspondence: Immacolata Porreca, Revvity, 8100 Cambridge Research Park, Cambridge CB25 9TL, UK.

E-mail: [immacolata.porreca@revvity.com](mailto:immacolata.porreca@revvity.com)

Molecular Therapy Vol. 32 No 8 August 2024 © 2024 The Authors.  
Published by Elsevier Inc. on behalf of The American Society of Gene and Cell Therapy.  
This is an open access article under the CC BY-NC-ND license (<http://creativecommons.org/licenses/by-nc-nd/4.0/>).

GEN Biotechnology  
Volume 1, Number 1, 2022  
Mary Ann Liebert, Inc.  
DOI: 10.1089/genbio.2021.0010



GEN  
Biotechnology

### REVIEW

## Genome Editing of Pluripotent Stem Cells for Adoptive and Regenerative Cell Therapies

Robert Blassberg<sup>1</sup>

### Abstract

The explosion of interest in CRISPR-Cas systems over recent years has delivered rapid advances in precision genome engineering technologies that hold enormous potential for application across the biotechnology sector. Within the cell therapy space, adoptive immunotherapies continue to lead the translation of increasingly sophisticated genetic engineering strategies toward clinical application. Although early focus was placed on engineering primary immune cells collected from patient blood, the democratization of these therapeutics will require the development of off-the-shelf allogeneic cell therapy products that can be manufactured at scale. Pluripotent stem cell (PSC) represent an attractive substrate for such products, and its adoptive cell therapies are common cures how genome editing can overcome generation technologies in the des

The CRISPR Journal  
Volume 4, Number 1, 2021  
© Mary Ann Liebert, Inc.  
DOI: 10.1089/crispr.2020.0035

### RESEARCH ARTICLE

## Development and Characterization of a Modular CRISPR and RNA Aptamer Mediated Base Editing System

Juan Carlos Collantes,<sup>1</sup> Victor M. Tan,<sup>1</sup> Huting Xu,<sup>1</sup> Melany Ruiz-Uriguen,<sup>1</sup> Amer Alasadi,<sup>1</sup> Jingjing Guo,<sup>1</sup> Hanlin Tao,<sup>1</sup> Chi Su,<sup>1</sup> Katarzyna M. Tyc,<sup>2</sup> Tommaso Selmi,<sup>3</sup> John J. Lambourne,<sup>4</sup> Jennifer A. Harbottle,<sup>5</sup> Jesse Stombaugh,<sup>6</sup> Jinchuan Xiao,<sup>7</sup> Ceri M. Wiggins,<sup>8</sup> and Shengkan Jin<sup>1\*</sup>

### Abstract

Conventional CRISPR approaches for precision genome editing rely on the introduction of DNA breaks (DSB) and activation of homology-directed repair (HDR), which is inherently genotoxic in somatic cells. The development of base editing (BE) systems that edit a target base without requirement of DSB or HDR offers an alternative. Here, we describe a novel BE system called Pin-point™, its base-modifying enzyme through an RNA aptamer within the gRNA molecule. Pin-point is capable of modifying base pairs in the human genome with precision and low on-target indel formation, potentially be applied for correcting pathogenic mutations, installing premature stop codons, and introducing other types of genetic changes for basic research and therapeutic development.

The CRISPR  
Journal



Press  
Release  
Details  
News Details

VIEW ALL NEWS

Revvity Announces New License Agreement for Next-Generation Base Editing Technology

05/18/2023

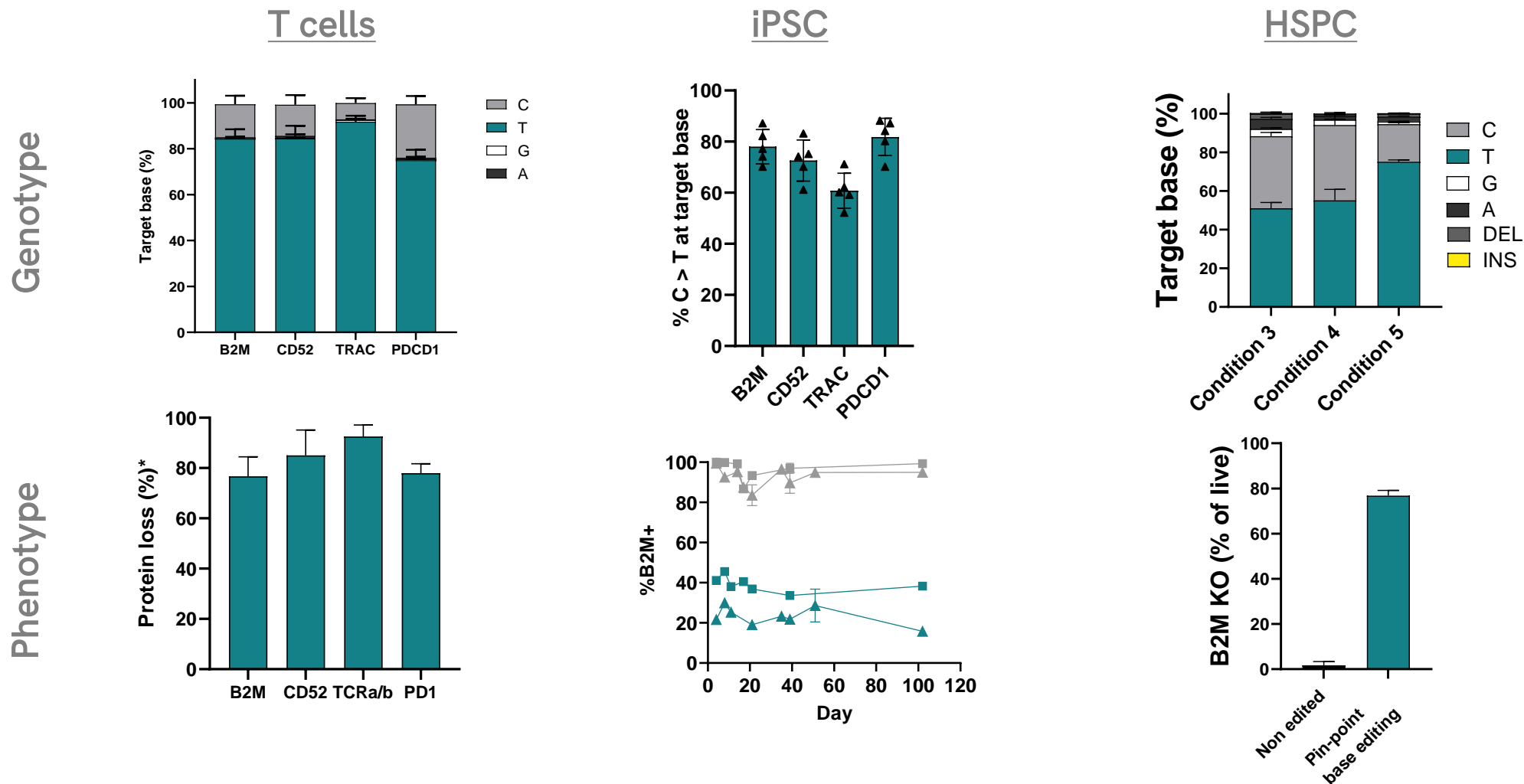
<https://www.sciencedirect.com/science/article/pii/S1525001624004234>

<https://doi.org/10.1089/genbio.2021.0010>

<https://www.biorxiv.org/content/10.1101/2023.06.20.545315v1>

Press release

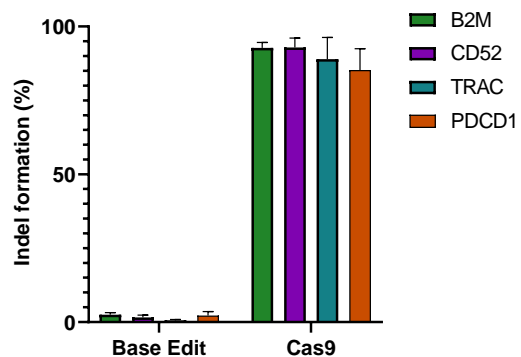
# Highly efficient and precise single and multiplex editing in T cells, iPSCs, and HSPCs



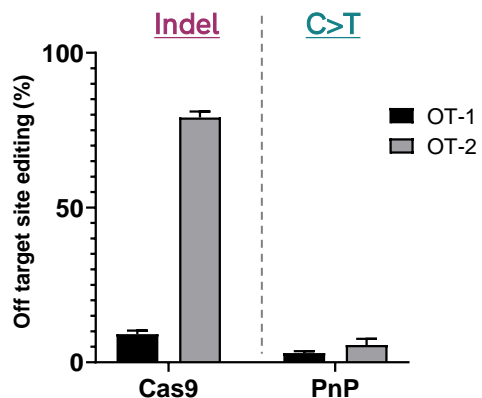
The Pin-point™ base editing platform is highly efficient in multiple cell types

# Strong safety profile compared to CRISPR/Cas9 editing

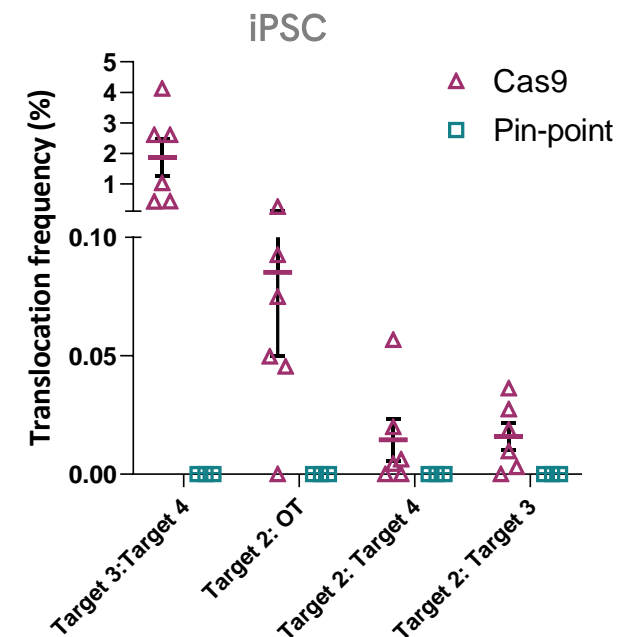
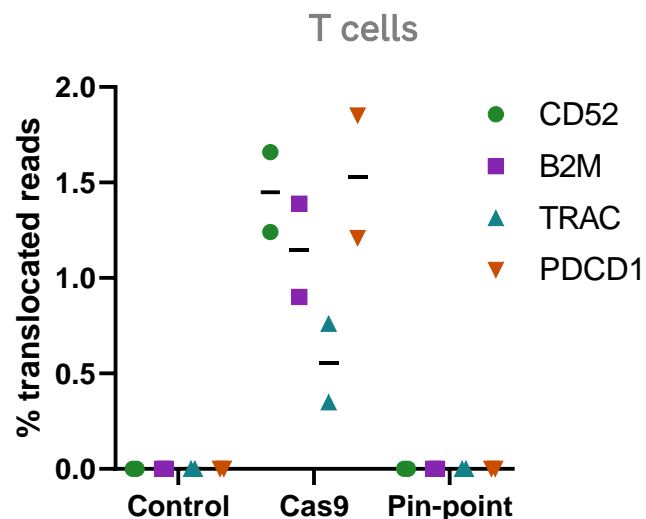
Reduced indel formation  
(T cells)



Negligible off-target edits (T cells)



No detected translocations with the Pin-point platform

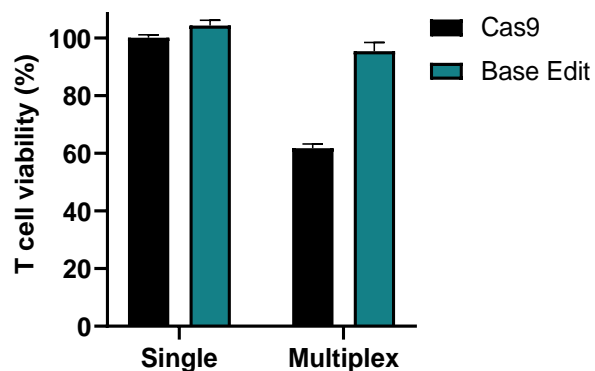


The Pin-point™ base editing platform is precise and avoids translocations

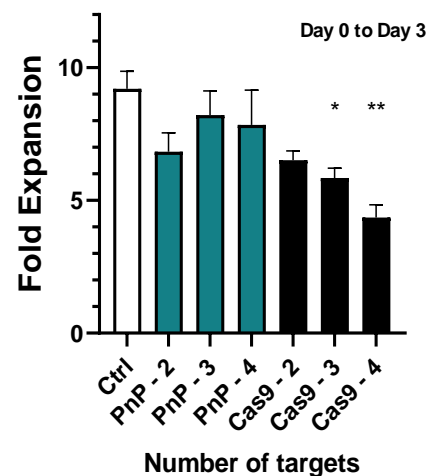
# High viability and cell health, even when editing multiple targets

## T cells

Cell viability maintained in multiplex

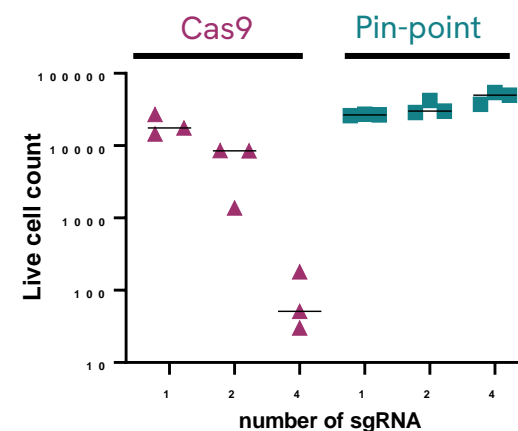


Rate of cell expansion unaffected

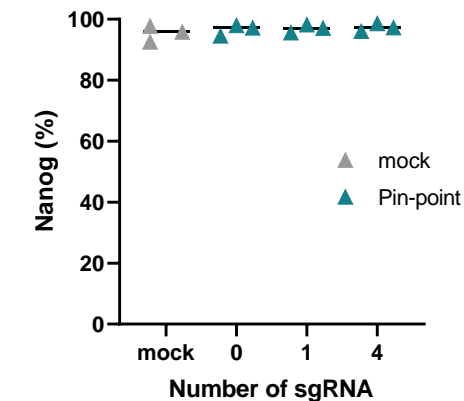


## iPSCs

High viability in multiplex

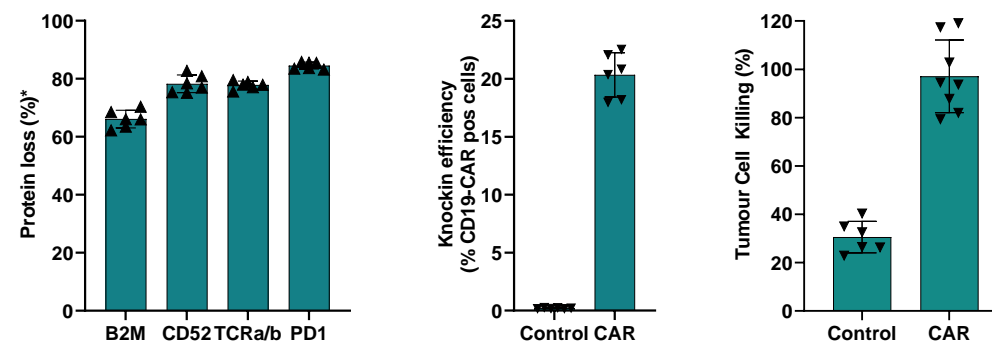
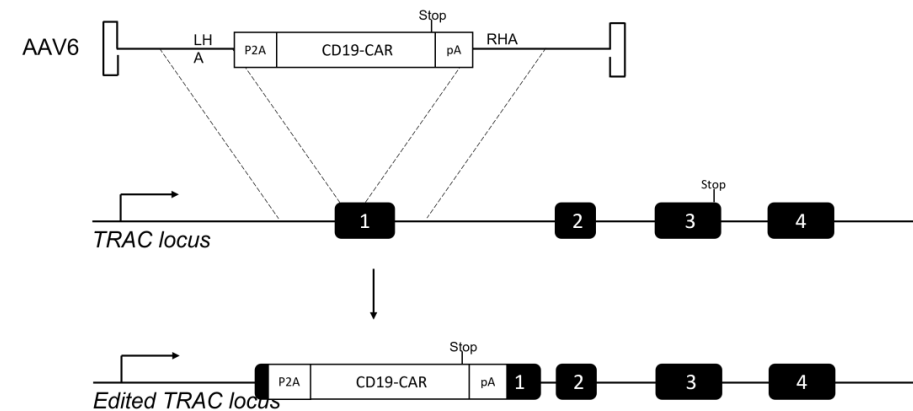
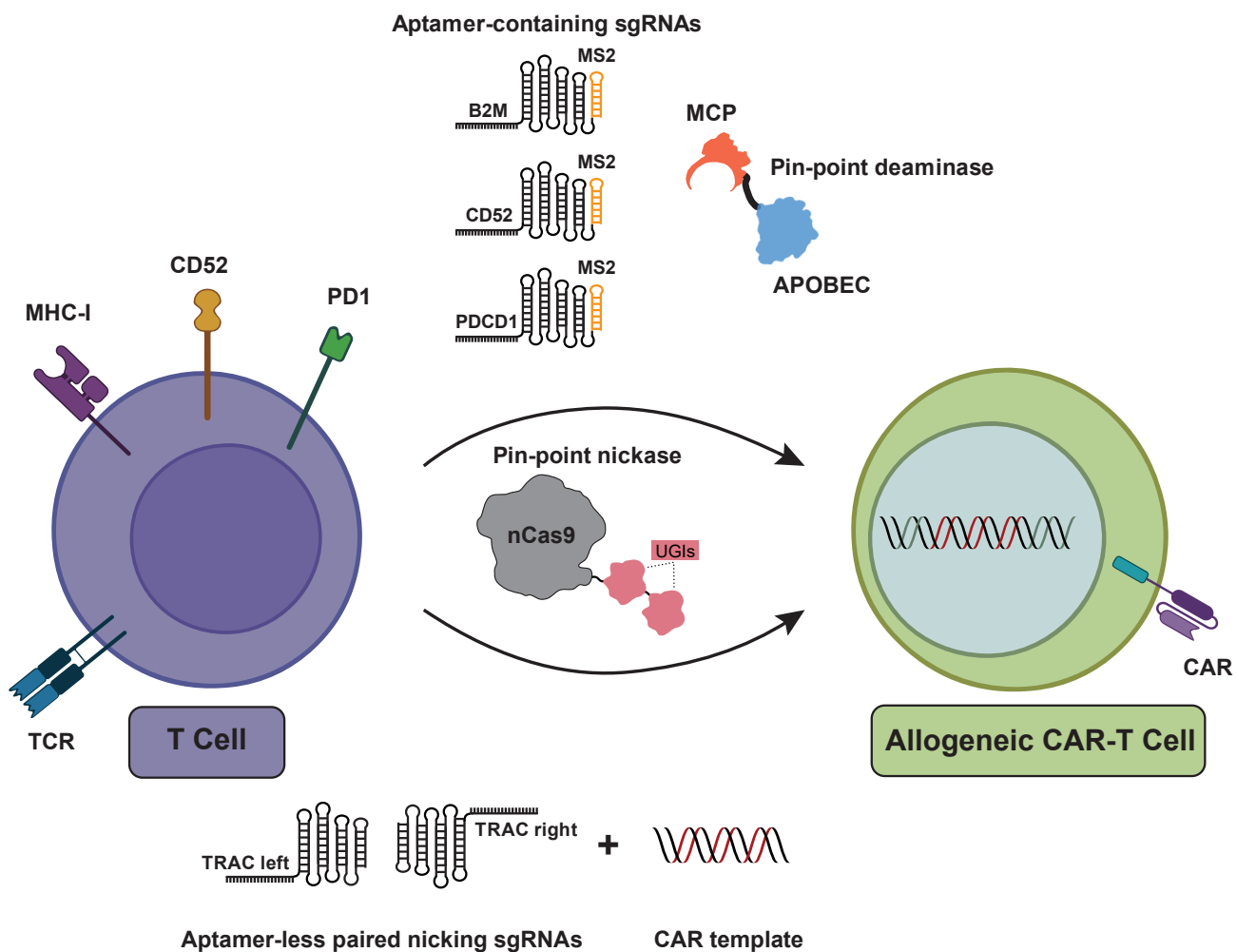


Cells retain pluripotency



The Pin-point™ base editing platform is gentle on sensitive cell types

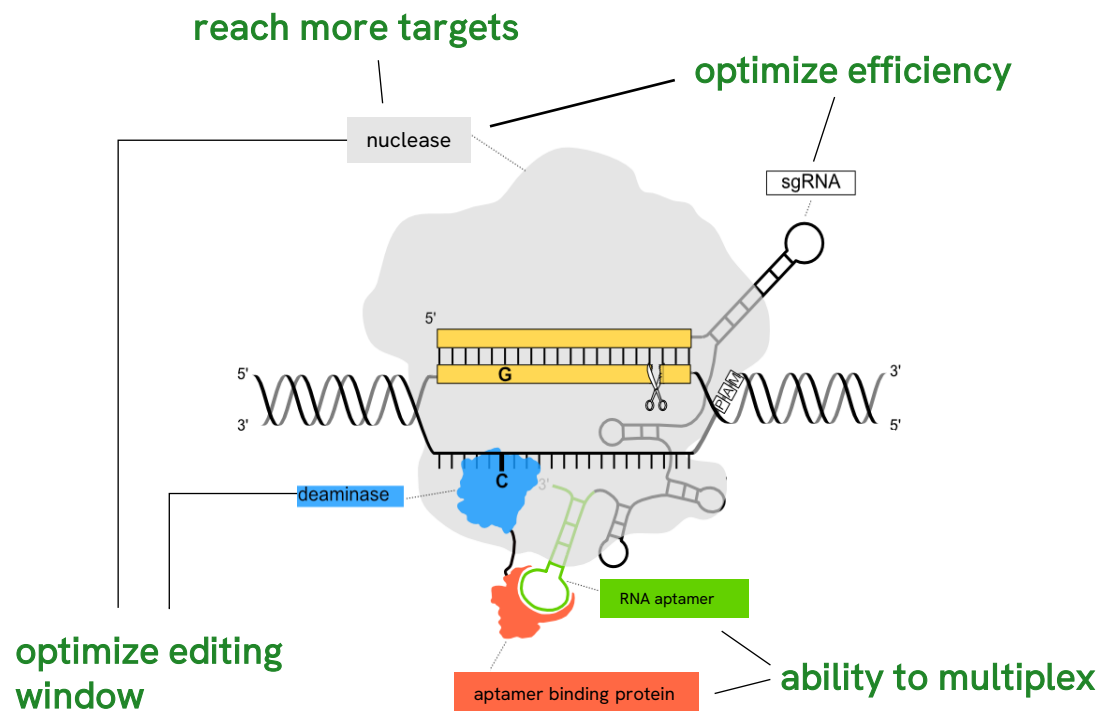
# A solution for complex engineering: One-step simultaneous knock-in and knockout in T cells



Efficient and accurate for concurrent transgene insertion and multiplex base editing

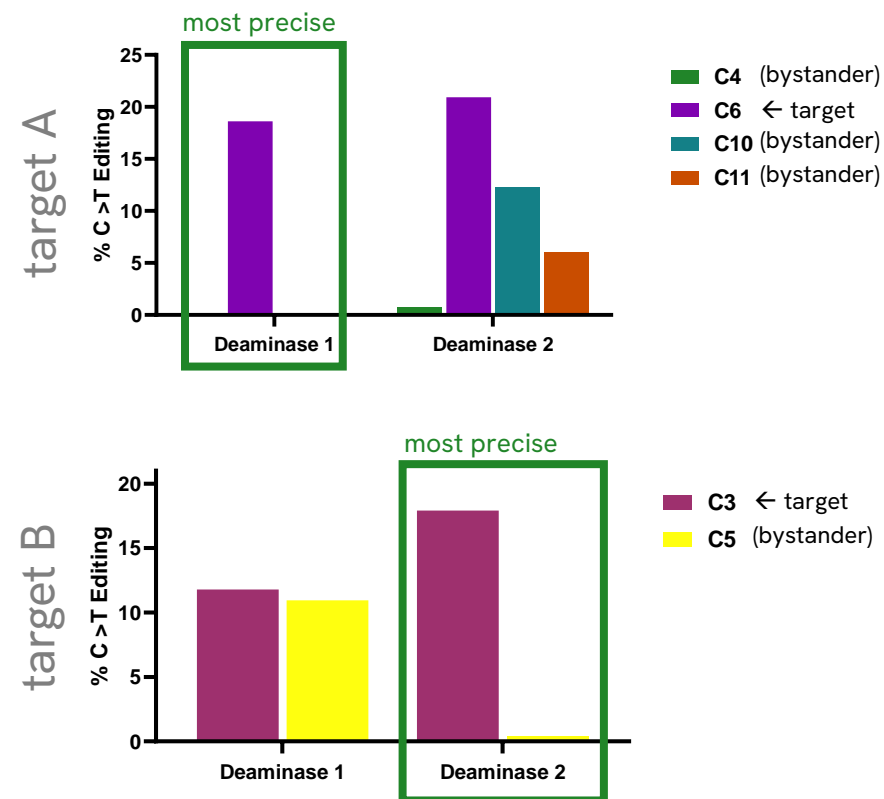
# Choose components for locus-specific optimization

Increased potential to correct more pathogenic SNVs that are not reachable with existing published systems\*



Schematic depicts nCas9 configuration

Example of optimization of the editing window by selecting the best guide RNA and deaminase pairs

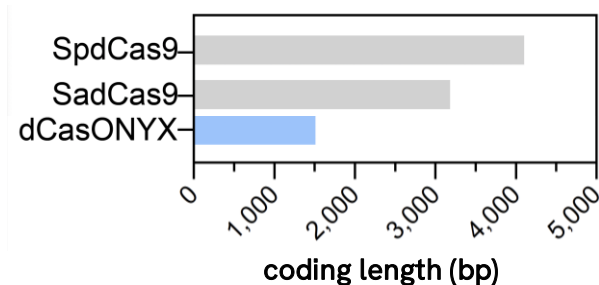


The modular Pin-point™ platform can be customized to combine optimal components for a wide range of base editing applications

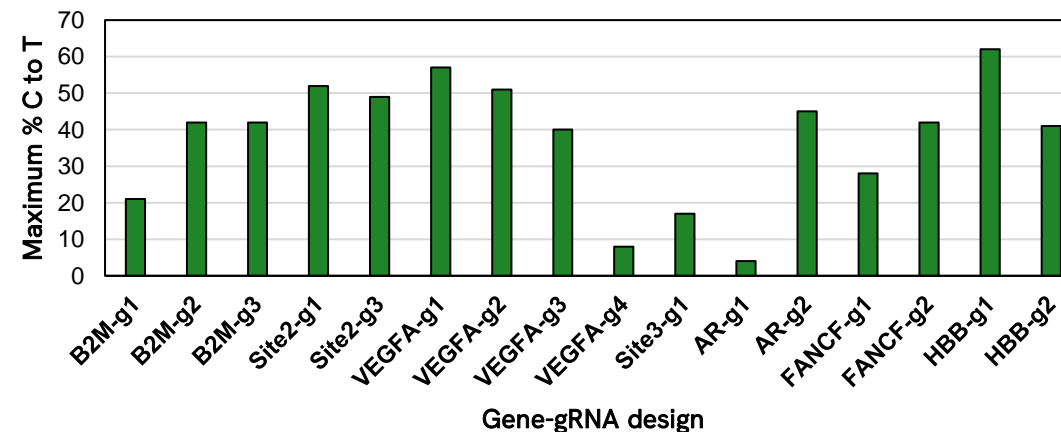
# The Pin-point™ platform configured with Epic Bio's ultracompact Type V effector protein, dCasONYX



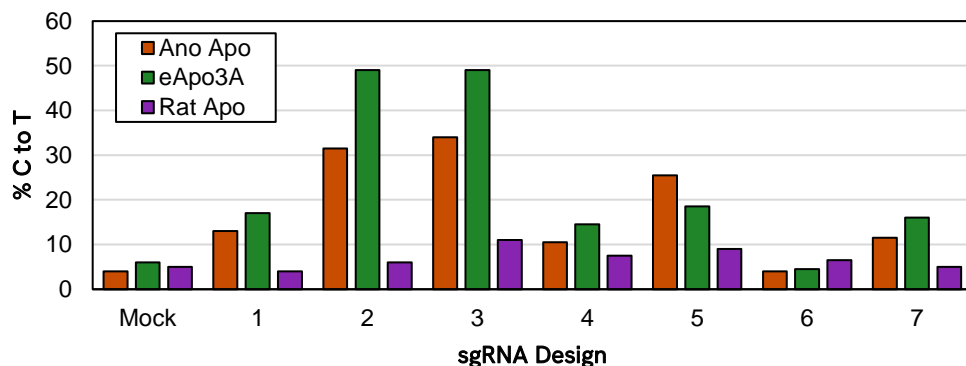
## Ultra-compact engineered Cas12f variant dCasONYX is under 1.5 kb



## Robust targeting capability



## Efficient editing with multiple deaminases when optimal gRNA scaffold is used



## Additional benefits of dCasONYX

- ✓ **Rapidly advancing to the clinic:** Epic Bio's asset EPI-321 for the potential cure for FSHD
- ✓ **Superior off-target profile:** described in Xin et al. Nature Communications<sup>1</sup>
- ✓ **Fully deactivated nuclease:** base editing without the risk of cutting the DNA
- ✓ **Low immunogenicity:** no prior exposure to dCasONYX in 10 human T cell donors, while 80% of human population have prior exposure to Cas9<sup>2</sup>
- ✓ **Small size:** Coding length less than 1.5 kb ideal for AAV packaging

A dCasONYX Pin-point configuration is one potential alternate to Cas9 for therapeutic applications



# The Pin-point™ platform is a transformational next-generation gene editing technology



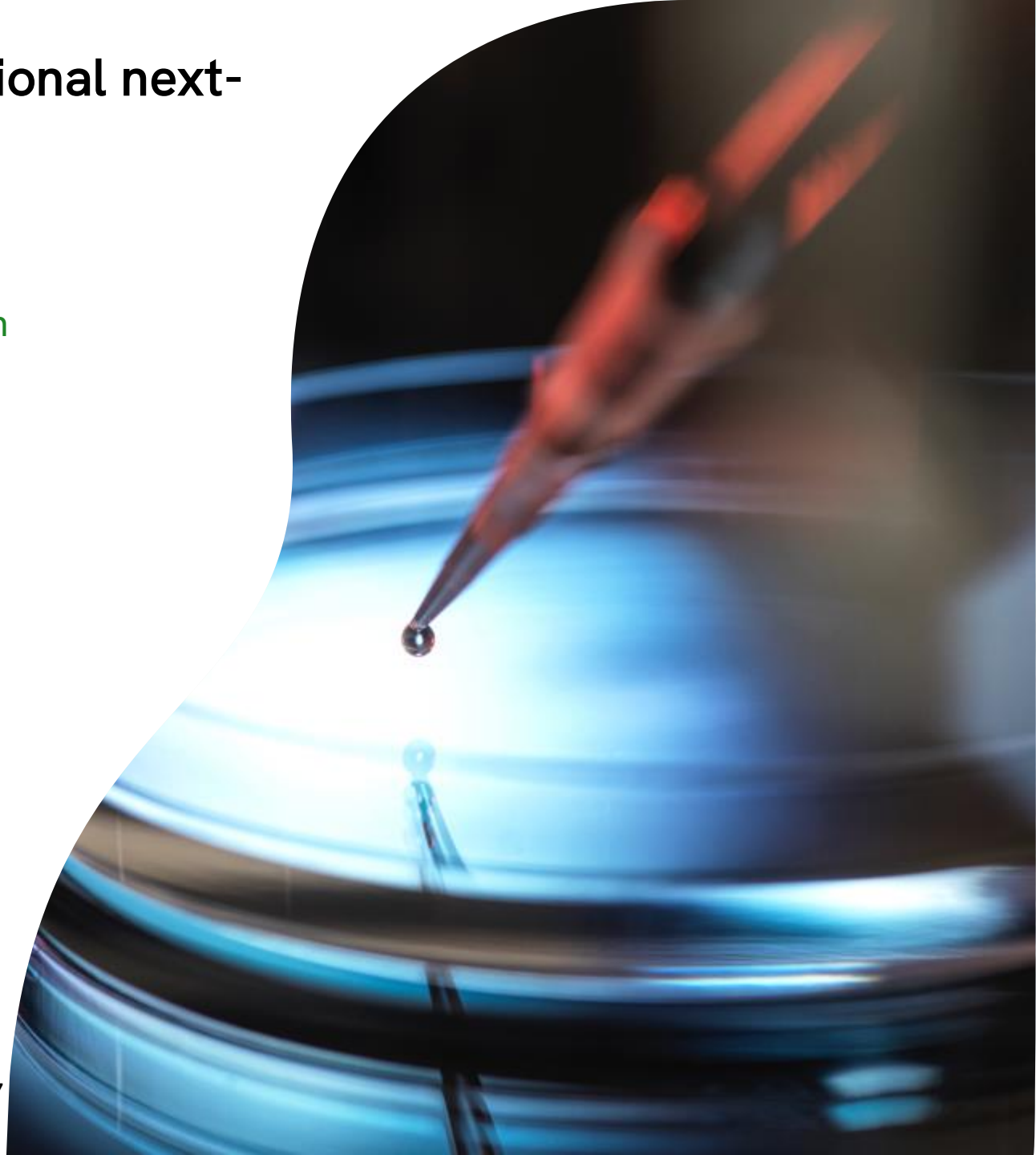
Highly effective editing platform, even for complex edits



Versatile technology modular and capable of generating locus-specific effects for novel therapies



Improved safety compared to standard CRISPR-Cas9 systems



# Access the Pin-point™ base editing platform



## Licensing

- Licenses for therapeutic development
- Comprehensive support
- Collaboration opportunities



## Research Reagents

- Synthetic off-the-shelf reagents
- Validated controls
- Custom guide RNAs



## Services

- Tiled pooled screening
- Functional genomics
- Cell models

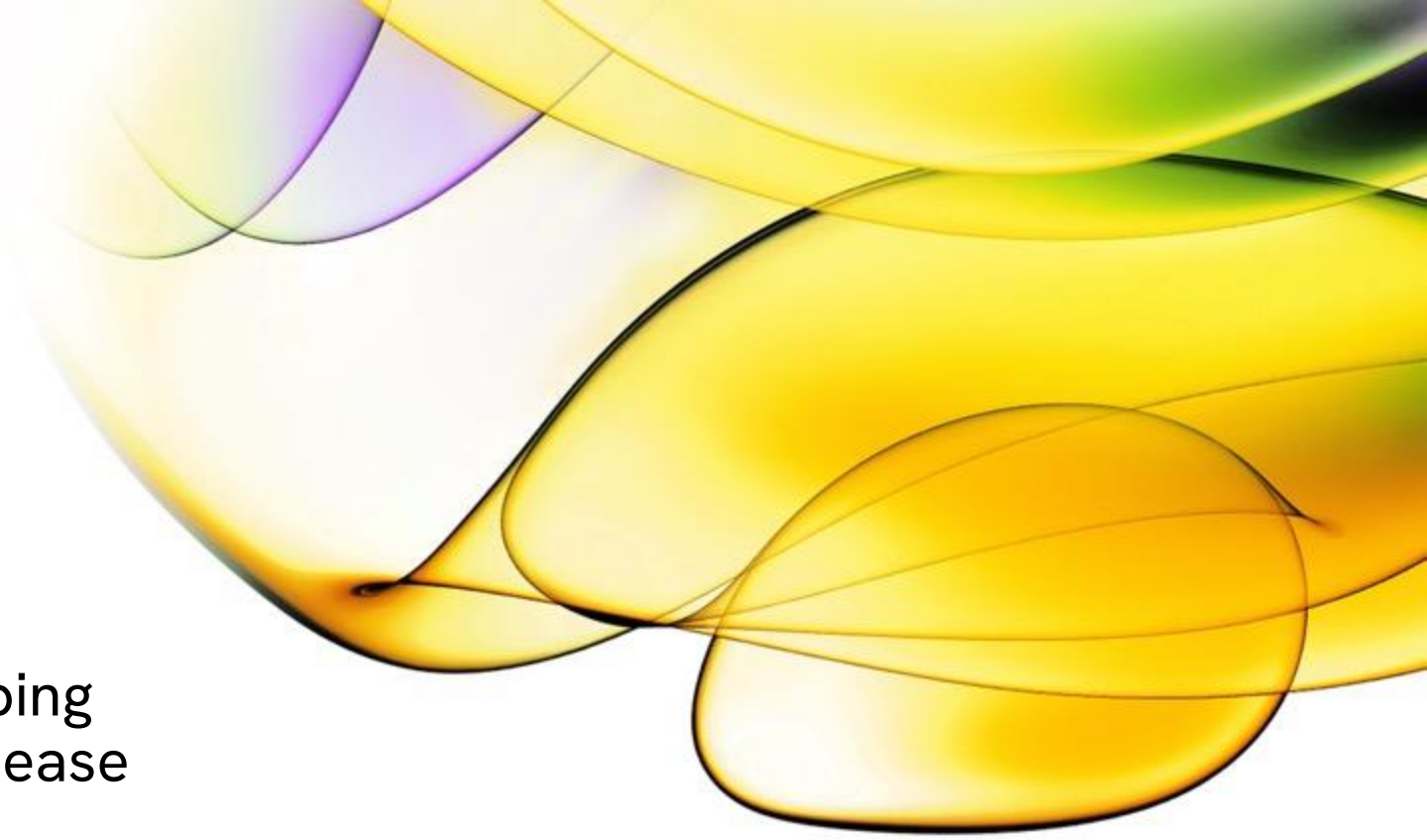
<https://horizondiscovery.com/en/gene-editing/pin-point-base-editing-platform>

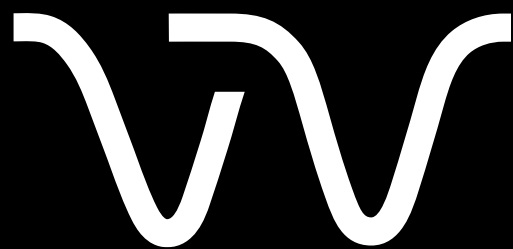
[BaseEditing@HorizonDiscovery.com](mailto:BaseEditing@HorizonDiscovery.com)

# revvity

We are a visionary partner in developing technologies and solutions across disease research pathways.

Here for a healthier humankind.





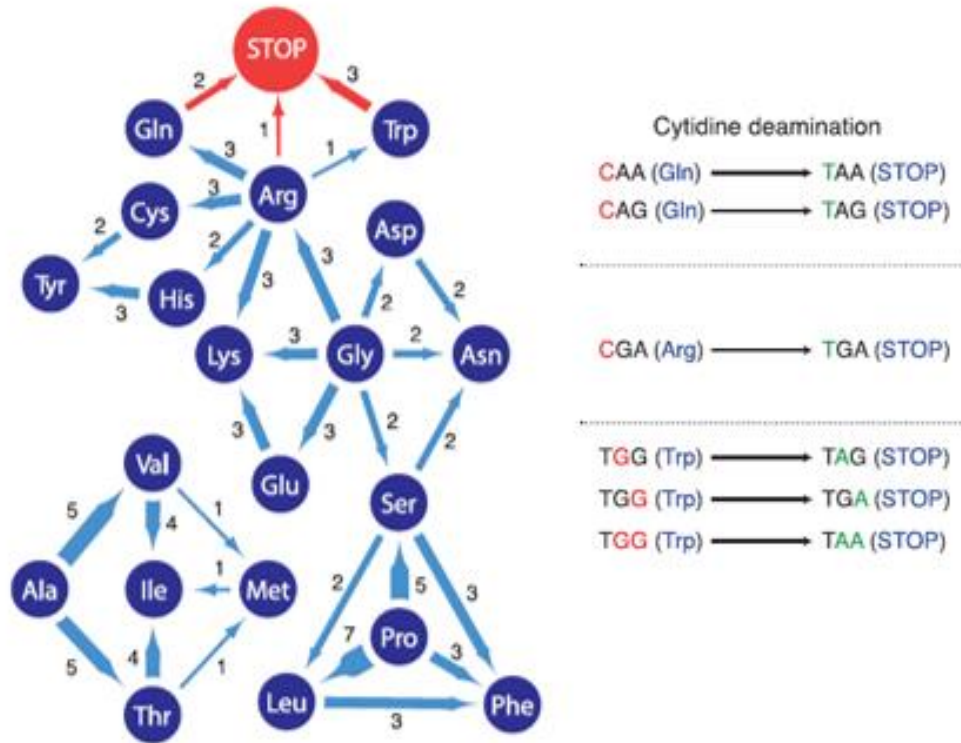
Revvity.com



Supplemental  
slides

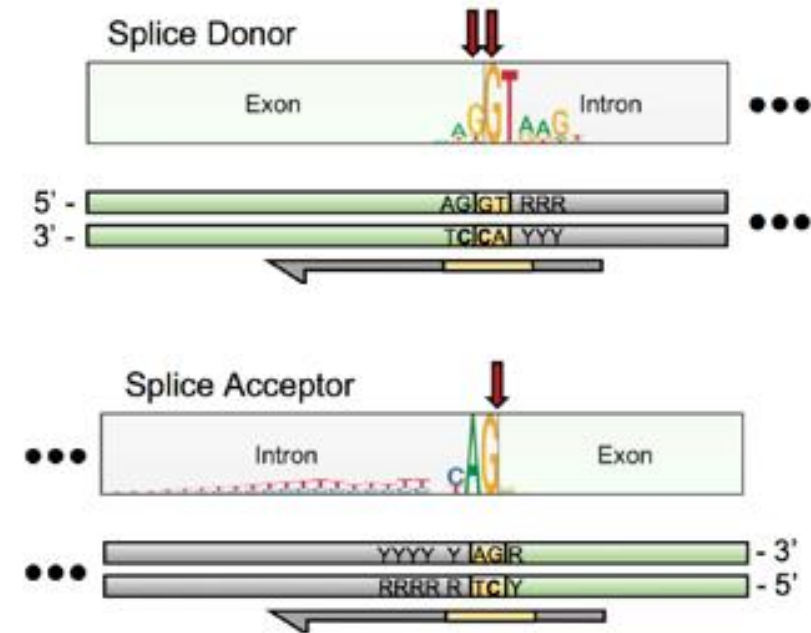
# How can base editing be used to create gene knockout?

## Premature Stop Codons



Billon et al. 2017. doi.org/10.1016/j.molcel.2017.08.008

## Splice Sites Disruption



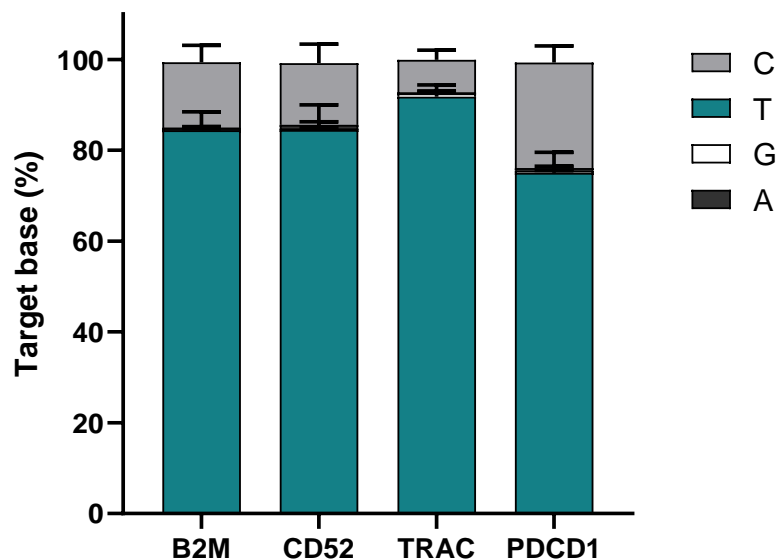
Webber et al 2019. doi.org/10.1038/s41467-019-13007-6



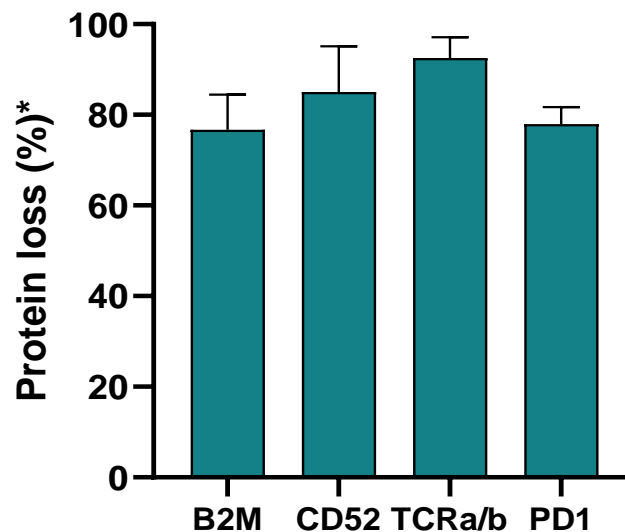
Validated  
performance in  
primary T cells

# Highly efficient and precise multiplex T cell editing

>75% editing in each target base without enrichment

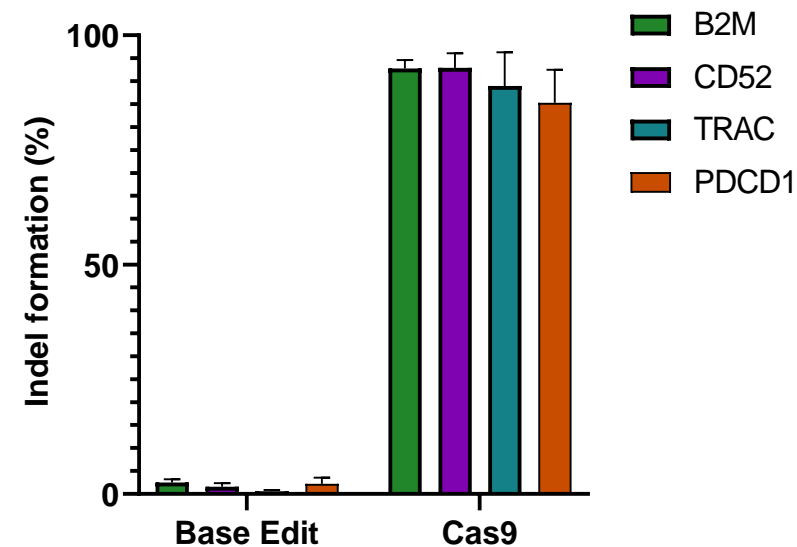


~50% of all 4 proteins knocked out (no enrichment)



\*Normalized to controls

No indel formation

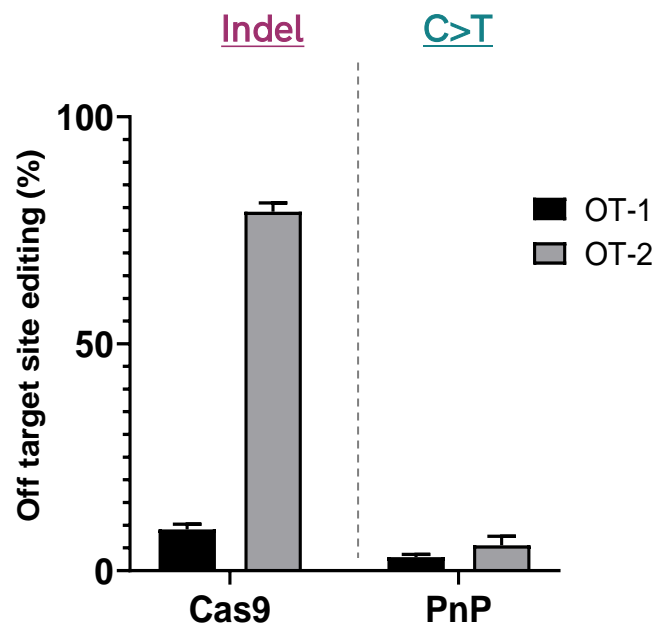


Pin-point™ base editing system is highly efficient and avoids potentially catastrophic DNA damage

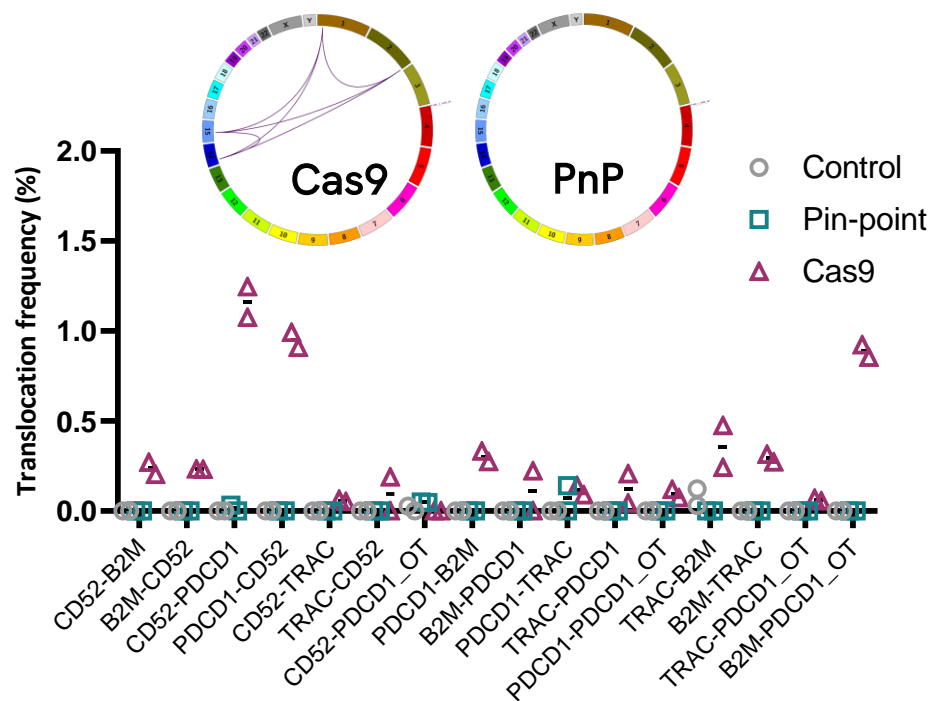


# Strong safety profile in T cells

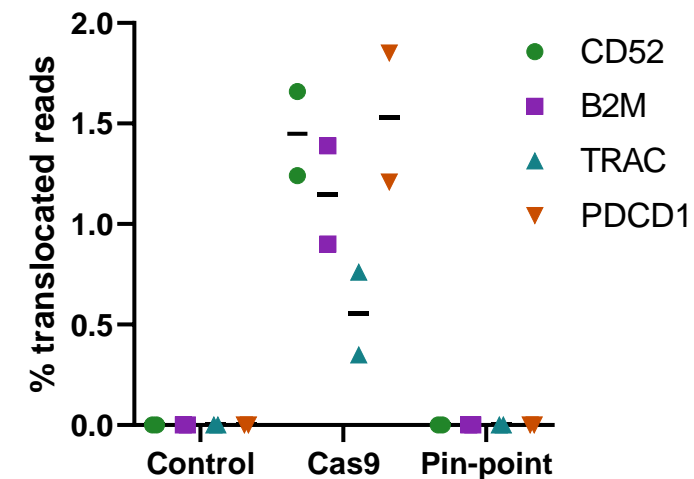
Negligible off-target edits



No detected translocations:  
targets and known off-target sites



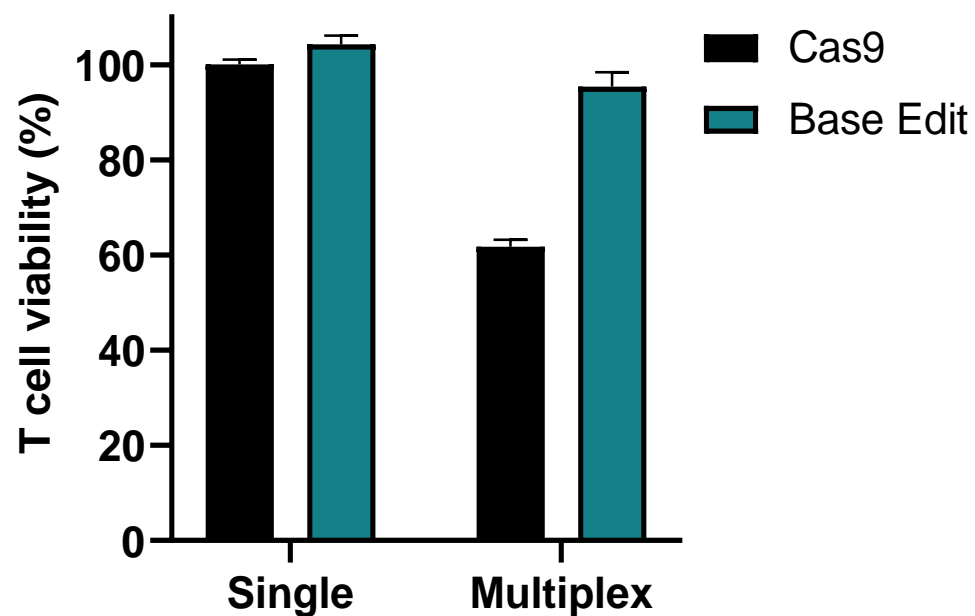
No detected translocations:  
fusions



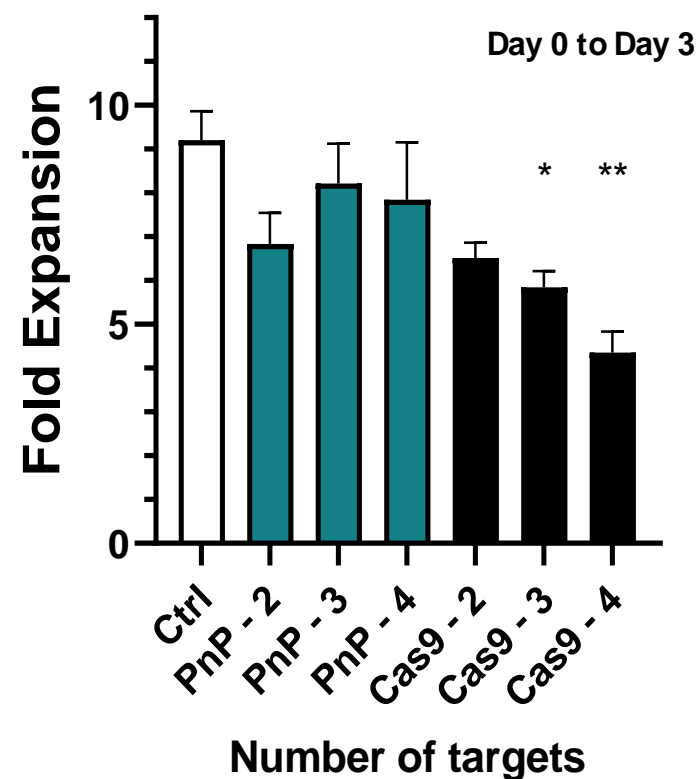
A cleaner and safer approach to multiplex gene editing in T cells

# No impact on T cell health

Cell viability maintained

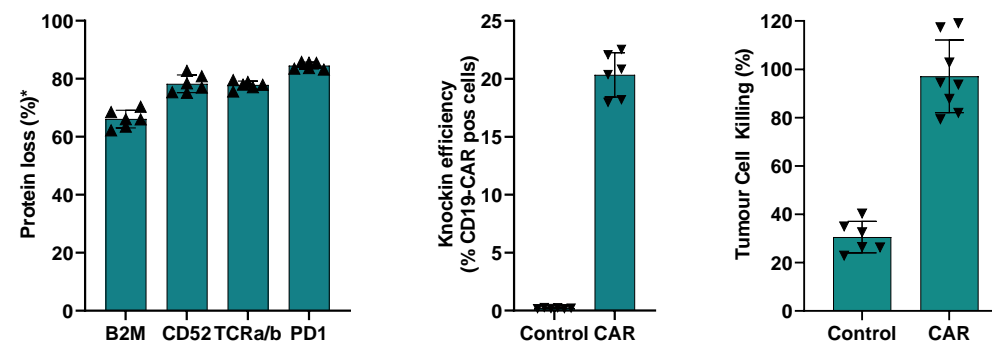
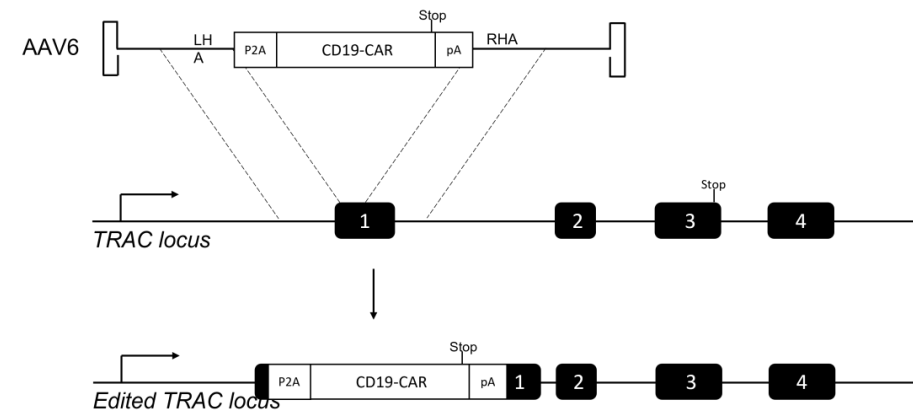
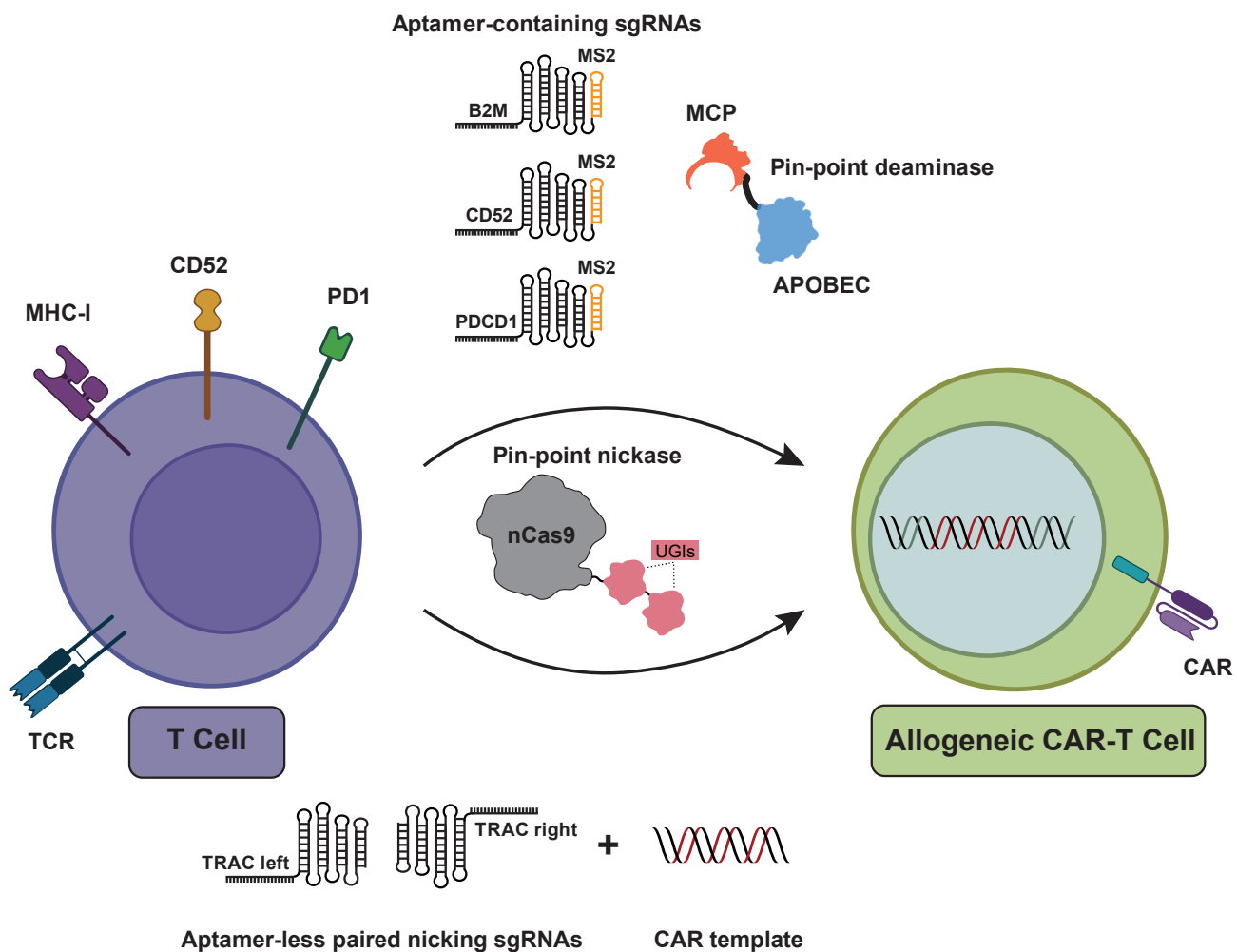


Rate of cell expansion unaffected

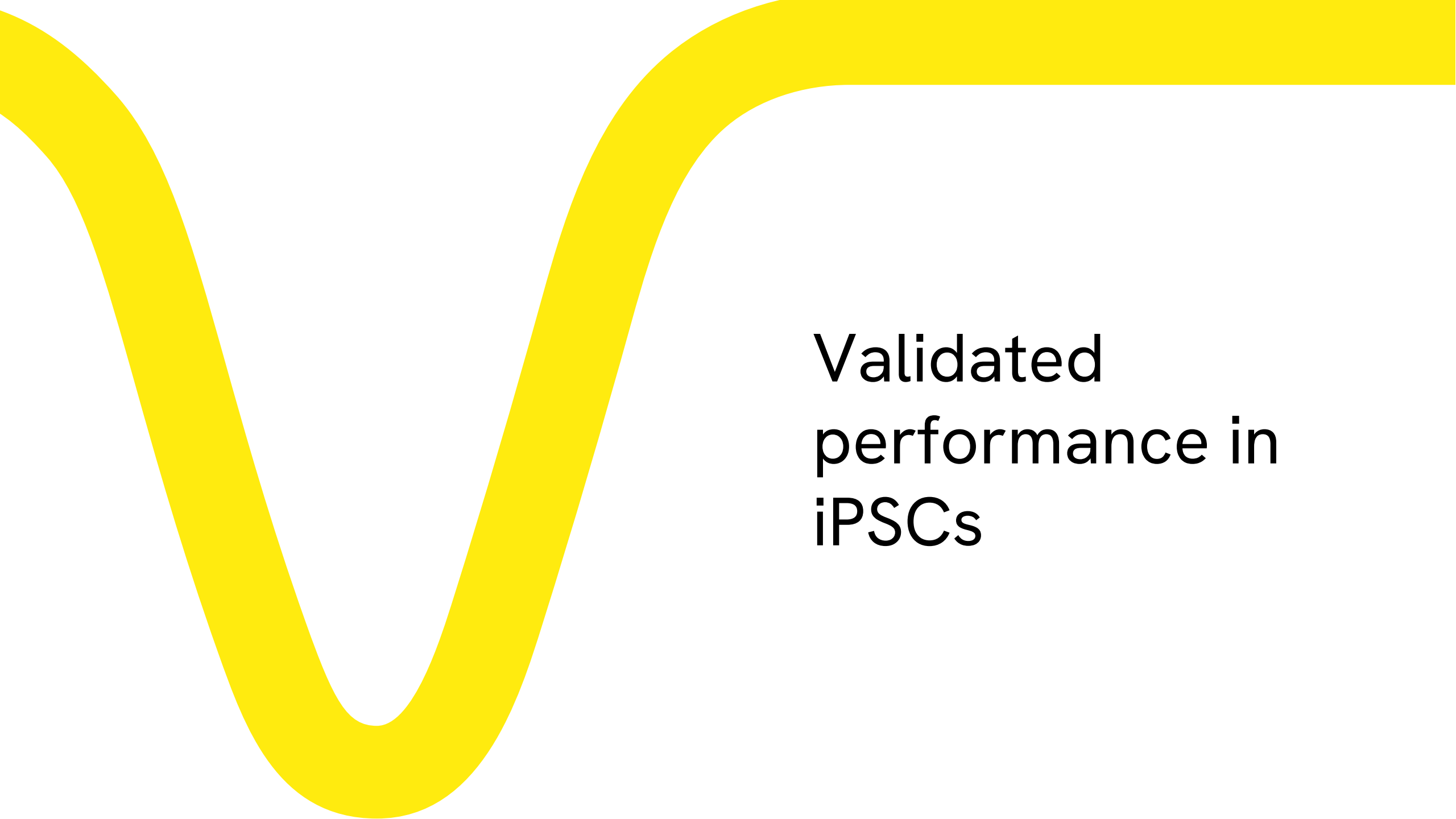


High multiplexing does not compromise cellular health or yield

# A solution for complex engineering: One-step simultaneous knock-in and knockout in T cells



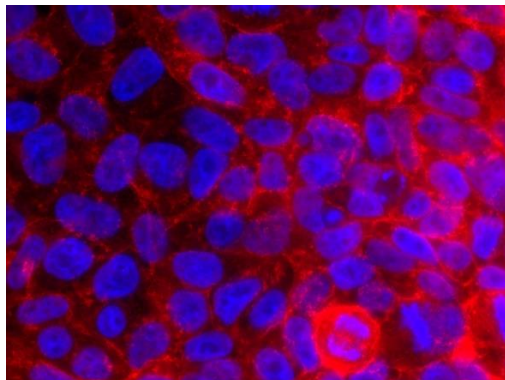
Efficient and accurate for concurrent transgene insertion and multiplex base editing



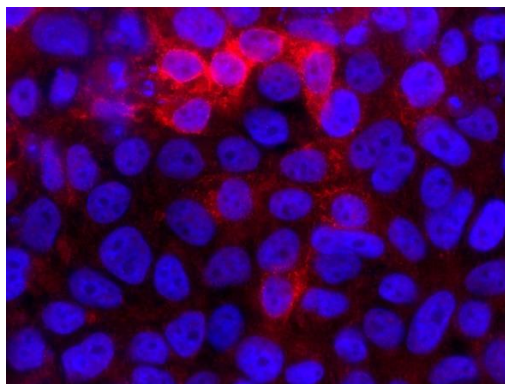
Validated  
performance in  
iPSCs

# Base editing with a Pin-point™ platform in iPSCs

Mock

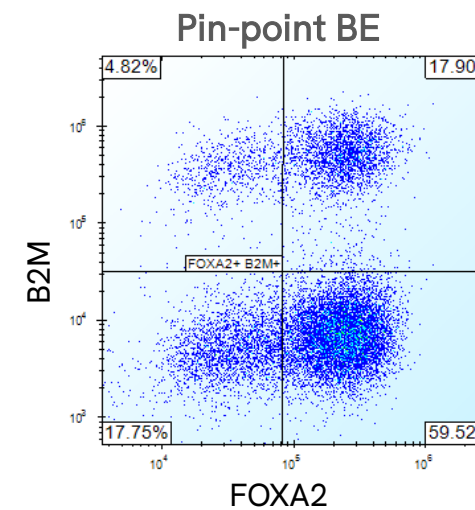
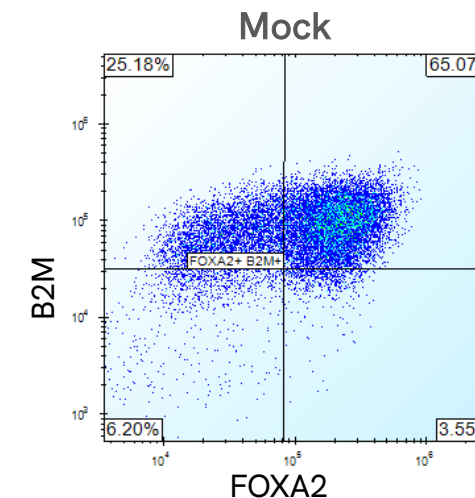
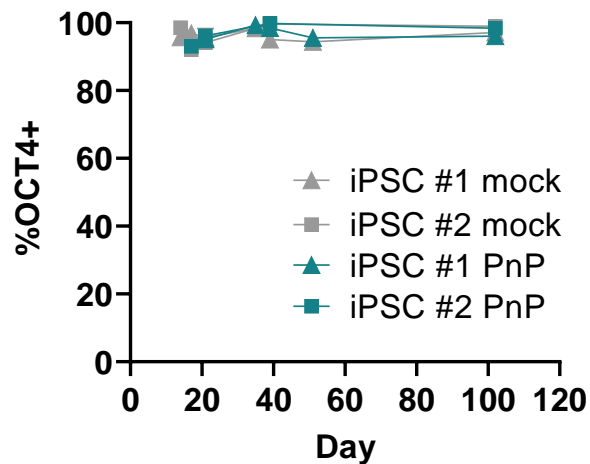
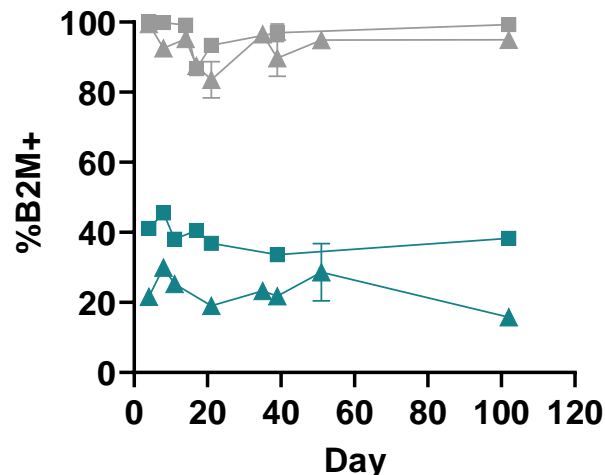


Pin-point BE



DAPI (nucleus)

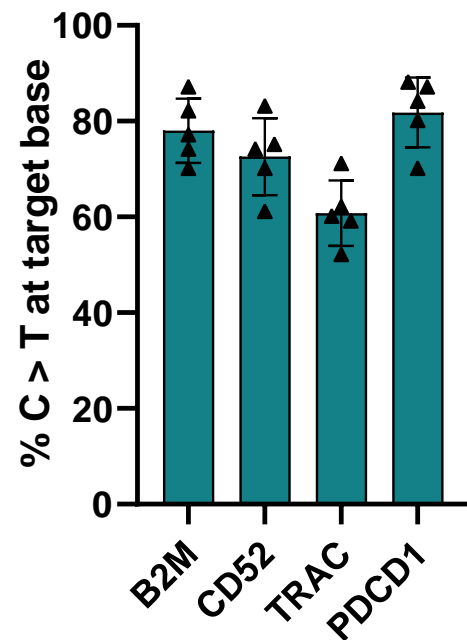
B2M



Edited iPSCs are stable with no growth defects when cultured up to 100 days and retain differentiation potential

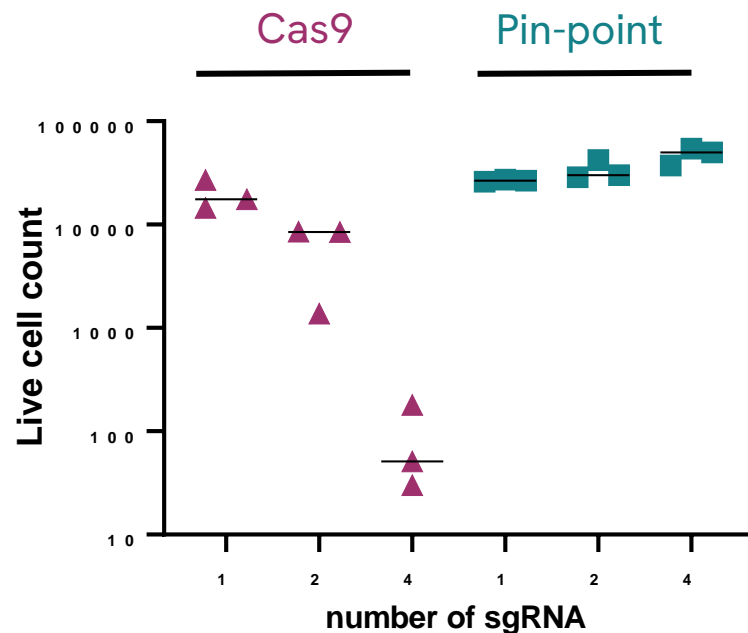
# Multi-gene editing in iPSCs

### Effective multiplex base editing



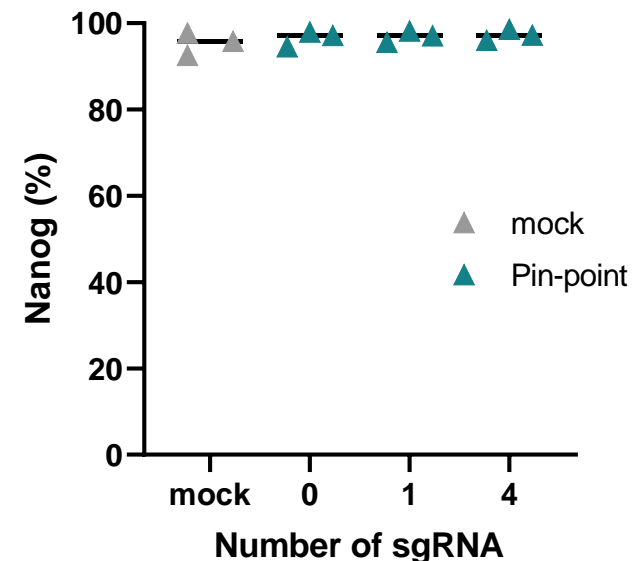
High base editing efficiency at target loci in a multiplex setting

### Edited cells are viable



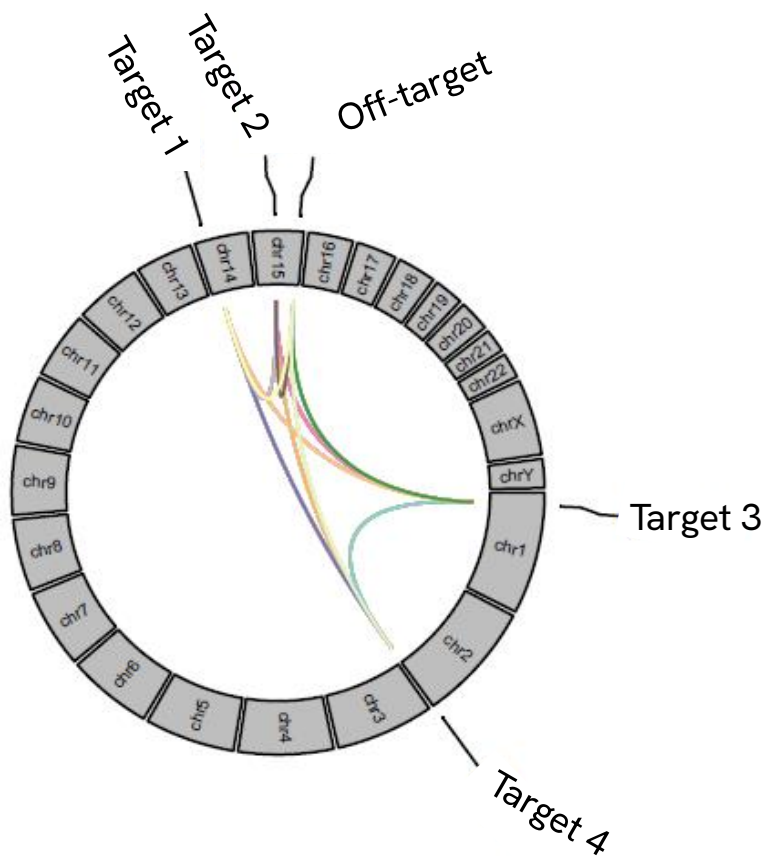
High survival of multi-edited iPSCs with a Pin-point™ system

### Edited cells retain their pluripotency



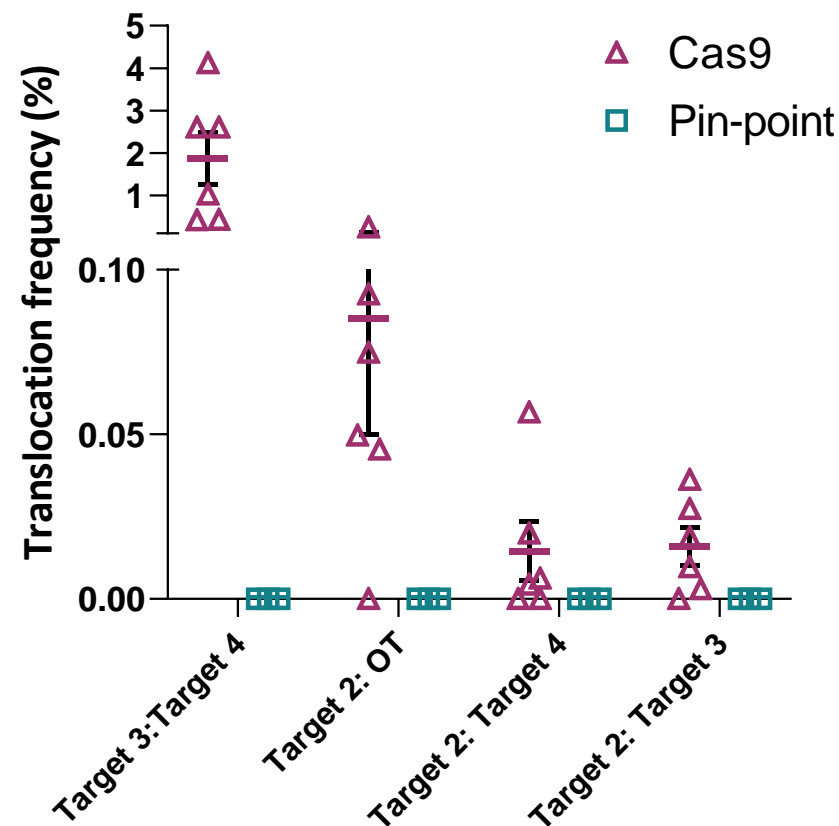
Pluripotency is retained in iPSCs edited with a Pin-point system

# Strong safety profile in iPSCs



in-silico predicted translocations  
previously validated in T cells

## Undetectable translocations after multiplex base editing with a Pin-point™ system



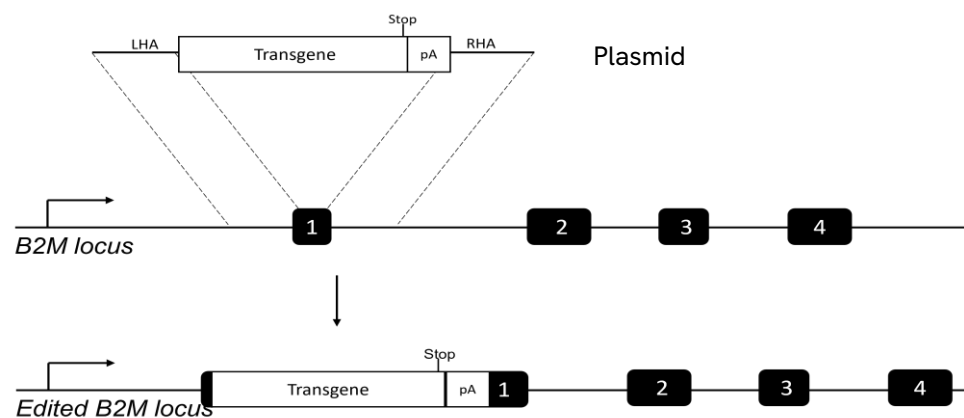
A cleaner and safer approach to multiplex gene editing in iPSCs

# Demonstrated simultaneous knock-in and multiple knockout in iPSCs

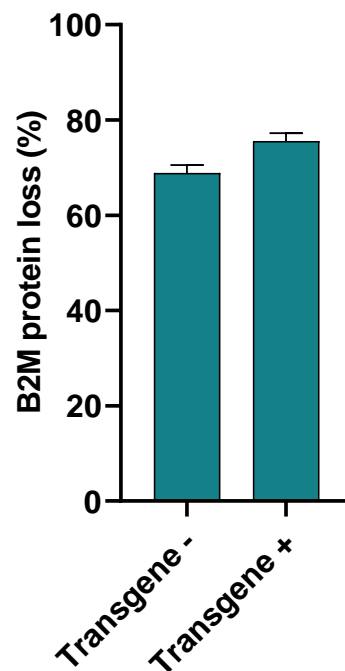
Base editing  
Knockout CIITA

&

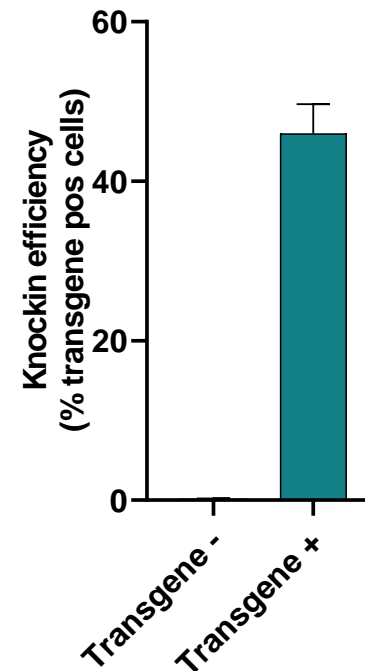
Insertion of a transgene  
in B2M (promoter-less GFP)



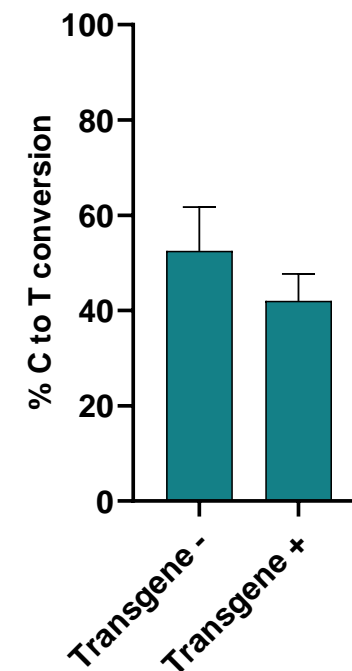
High level  
of B2M  
knockout



High level of  
transgene  
knock-in

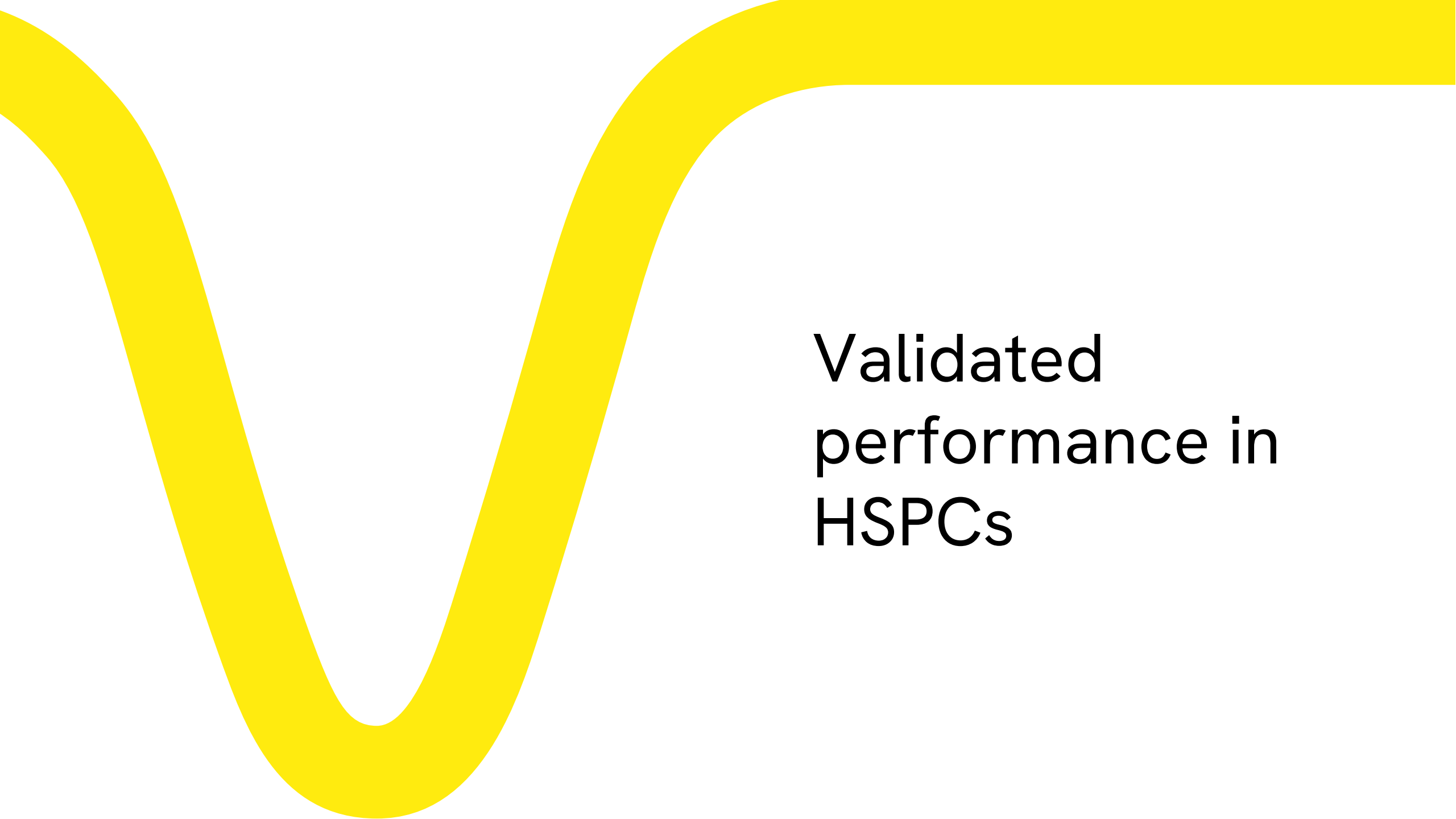


... and base  
editing  
(CIITA KO)



The Pin-point™ platform enables one-step simultaneous knock-in and multiple knockout in iPSCs

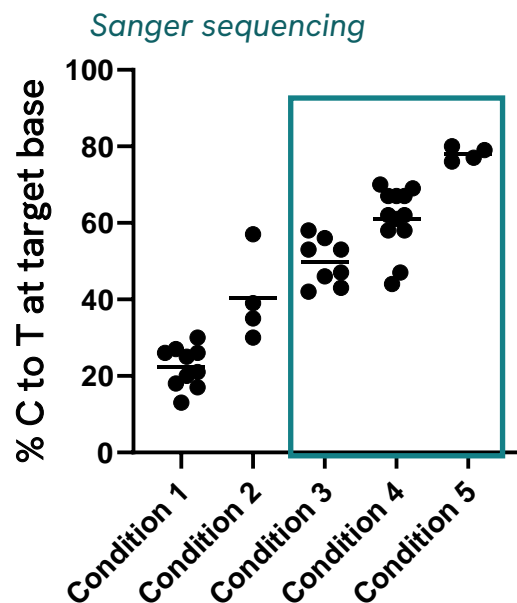




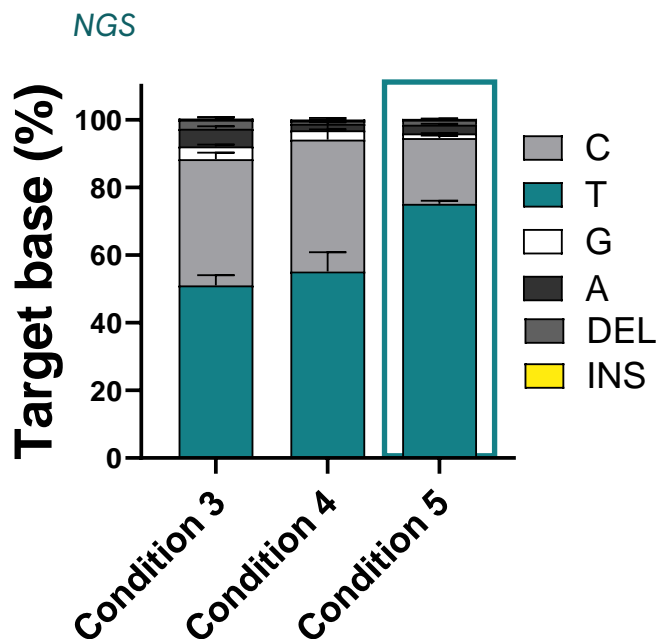
Validated  
performance in  
HSPCs

# Highly efficient base editing in HSPCs

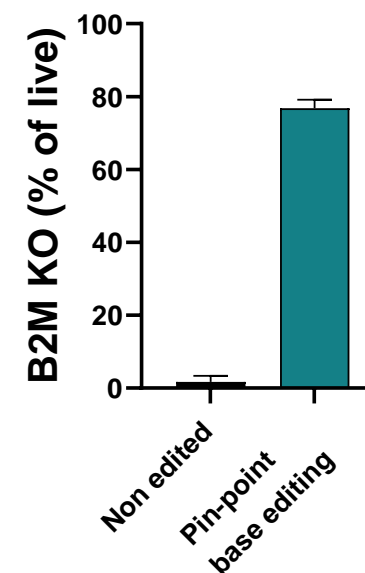
High level of editing achieved with optimised conditions



High level of editing and purity achieved at the target site



High level of B2M phenotypic knock-out

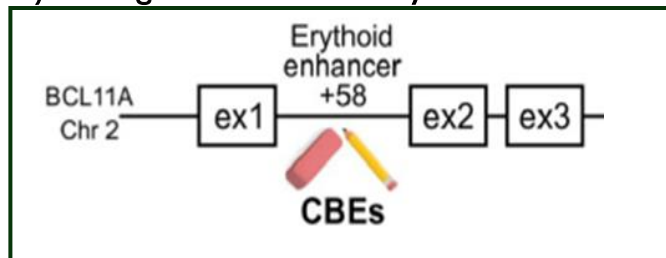


Optimized application of the Pin-point™ platform achieves high levels of editing in HSPCs with high purity of C to T conversion

# Therapeutic editing of HSPCs with the Pin-point™ platform

## Reactivating Fetal Haemoglobin

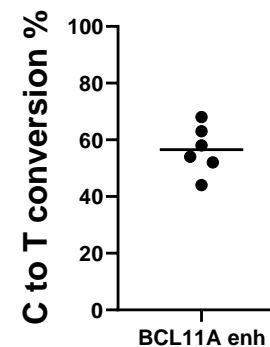
### 1) Editing of the BCL11A erythroid enhancer



## Editing Efficiency

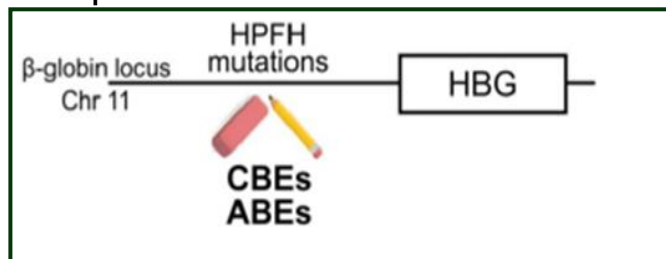
GATA1 binding site

| Base | T | T | T | A | T | C  | A | C | A | G | G | C | T | C | C | A | G | G | A | A |
|------|---|---|---|---|---|----|---|---|---|---|---|---|---|---|---|---|---|---|---|---|
| G    | 0 | 0 | 0 | 0 | 0 | 4  | 0 | 0 | 6 | - | - | 0 | 0 | 0 | 0 | 4 | - | - | 0 | 5 |
| A    | 0 | 0 | 0 | - | 0 | 0  | - | 0 | 0 | 0 | 0 | 0 | 0 | 0 | 0 | - | 0 | 0 | - | - |
| T    | - | - | - | 0 | - | 68 | 0 | 5 | 0 | 0 | 0 | 0 | - | 5 | 0 | 0 | 0 | 0 | 0 | 0 |
| C    | 0 | 0 | 0 | 0 | 0 | -  | 0 | - | 0 | 0 | 0 | - | 6 | - | - | 0 | 0 | 0 | 0 | 0 |



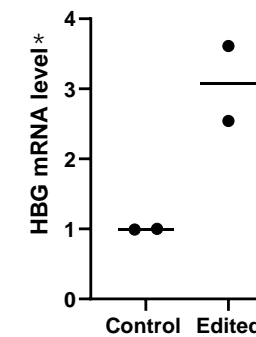
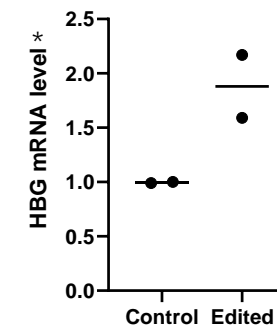
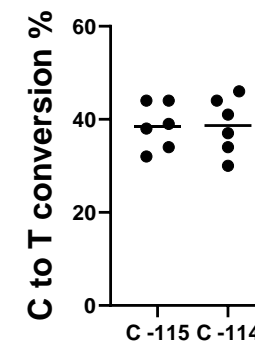
## Induction of Fetal Haemoglobin

### 2) Editing of the BCL11A binding site in the HBG promoter



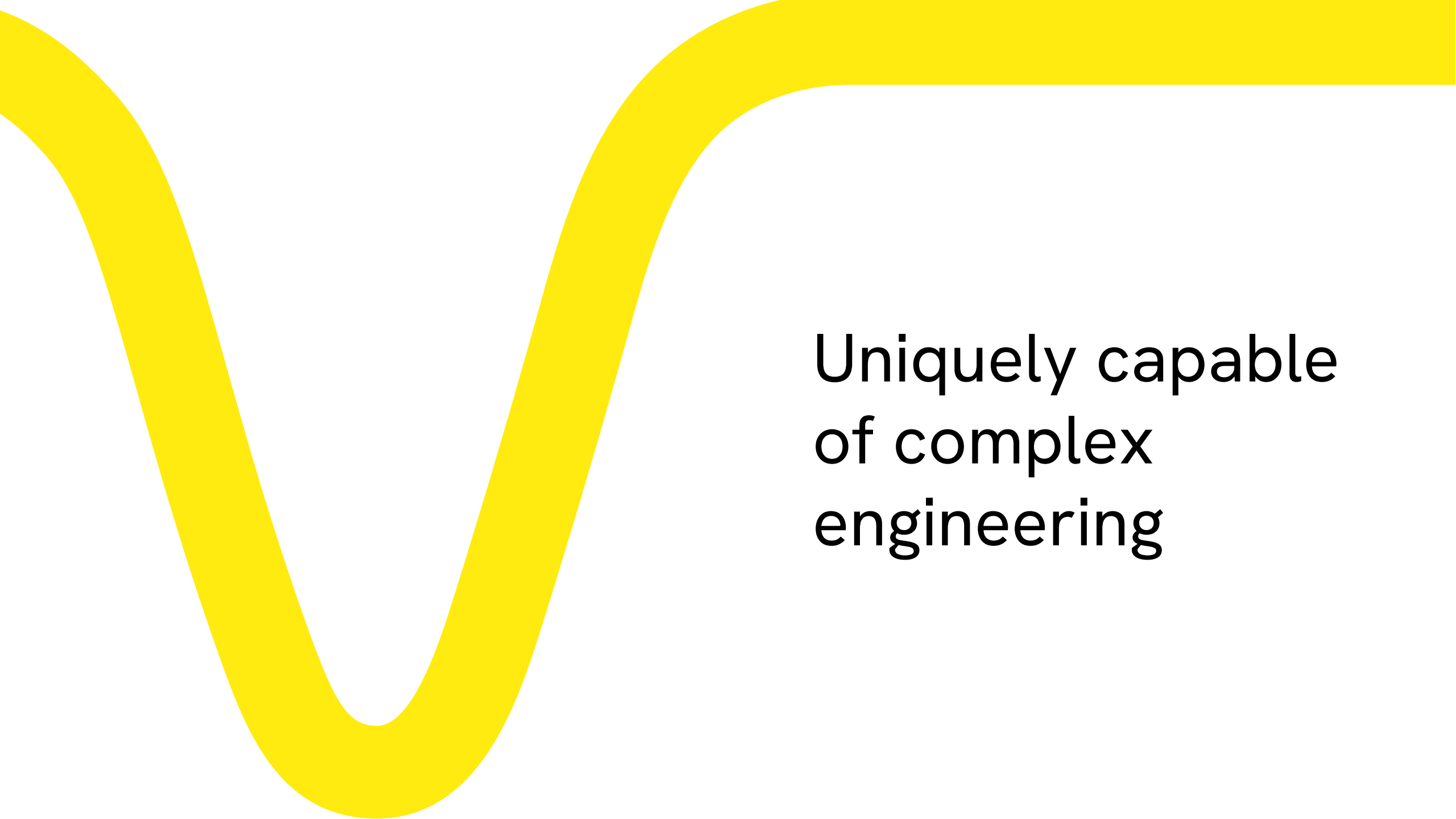
BCL11A binding site

| Base | C | T | T | G | A | C  | C  | A | A | T | A | G | C | C | T | T | G | A | C | A |
|------|---|---|---|---|---|----|----|---|---|---|---|---|---|---|---|---|---|---|---|---|
| G    | 0 | 0 | 0 | - | 0 | 0  | 0  | 0 | 0 | 0 | 0 | - | 0 | 0 | 0 | 0 | - | 0 | 0 | 0 |
| A    | 0 | 0 | 0 | 0 | - | 0  | 0  | - | - | 0 | - | 0 | 0 | 0 | 0 | 0 | 0 | - | 0 | - |
| T    | 0 | - | - | 0 | 0 | 38 | 37 | 0 | 0 | 0 | - | 0 | 5 | 0 | 0 | - | - | 0 | 0 | 0 |
| C    | - | 0 | 0 | 0 | 0 | -  | -  | 0 | 0 | 0 | 0 | 0 | - | - | 0 | 0 | 0 | 0 | 0 | 0 |



\*qPCR data expressed as HBG/HBA and normalised on expression in control samples

The Pin-point base editing platform achieves therapeutic editing in HSPCs



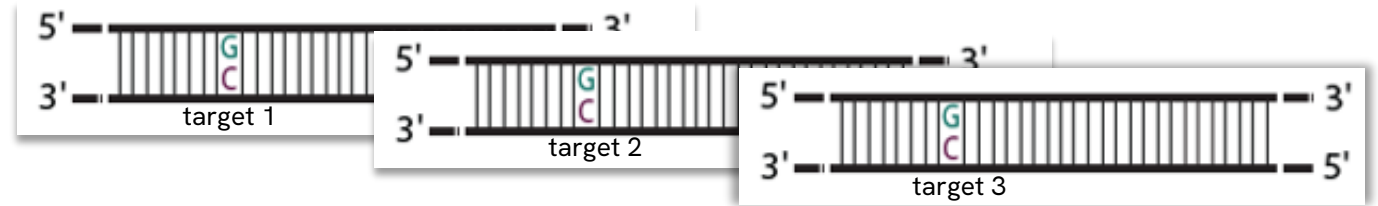
Uniquely capable  
of complex  
engineering

# A solution for complex engineering

One-step simultaneous knock-in and multiple knockout in T cells

Base Editing with aptamer gRNAs

Knockout B2M, CD52, PDCD1

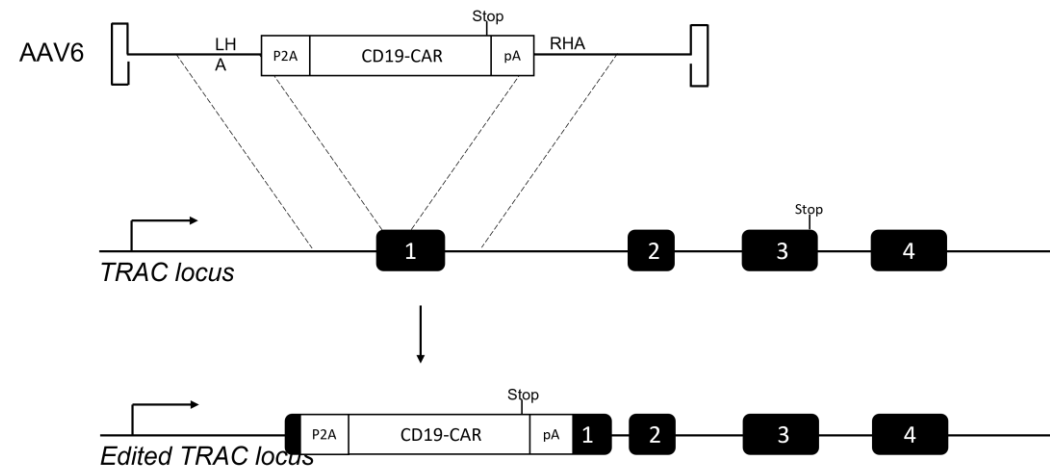


## One transfection

- Pin-point deaminase
- Pin-point aptamer gRNA (3)
- Nickase Cas
- Nicking gRNAs (2)
- Donor insert template

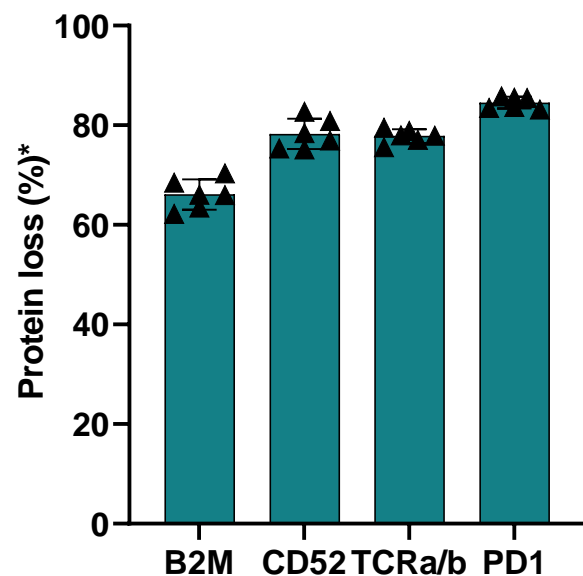
Insertion of a transgene by non-aptamer nicking gRNAs

CAR in TRAC

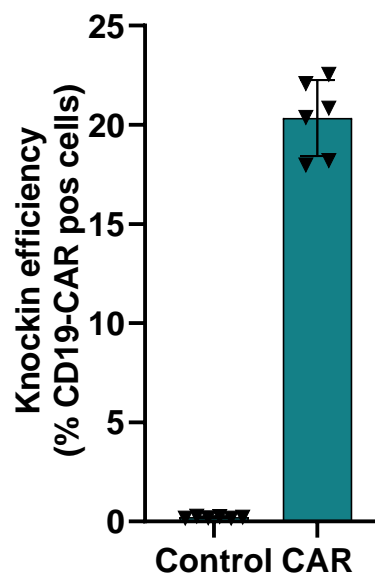


# Streamlined creation of CAR-T cells is enabled with the Pin-point™ platform

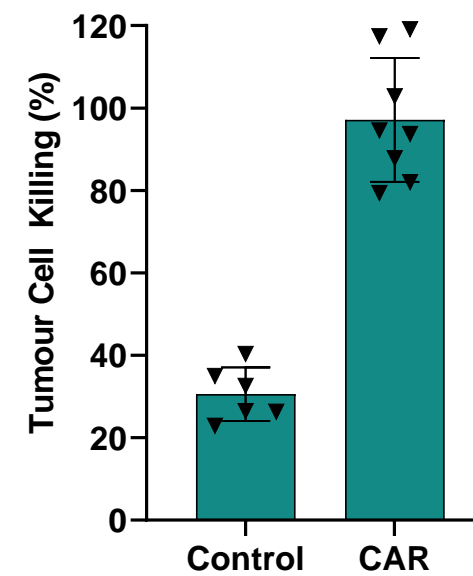
Multiple proteins are knocked out



... while enabling protein knock-in



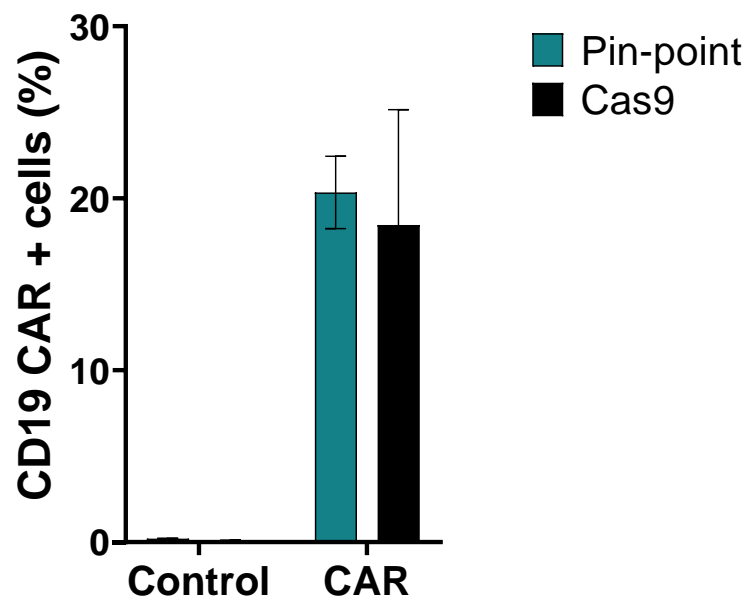
... and weaponizing T cells against cancer cells



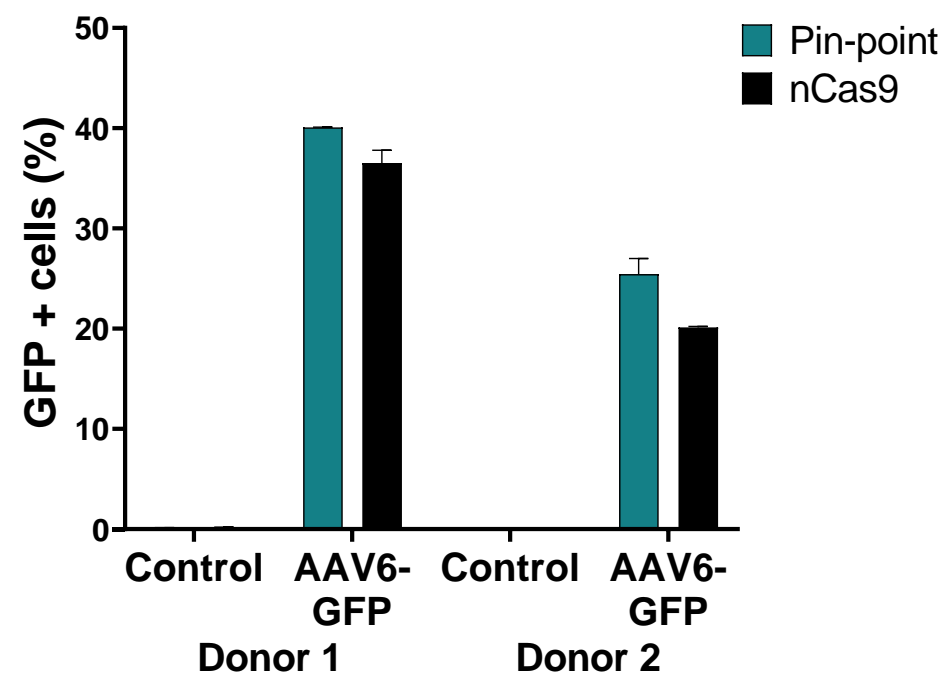
The Pin-point platform is efficient and accurate for concurrent transgene insertion and multiplex base editing

# No loss of efficiency in payload deliveries

Equivalent to dsDNA knock-in



Presence of modular deaminase has no impact on knock-in



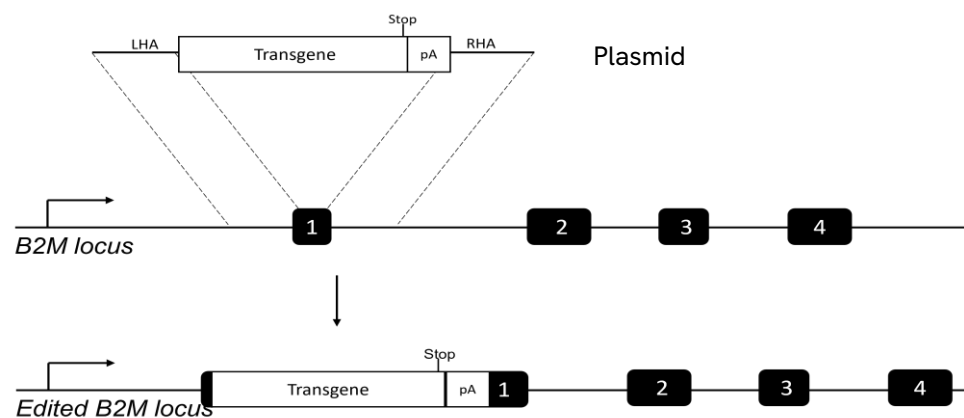
The Pin-point™ platform can deliver payloads equivalently to standard Cas9 or nCas9 knock-in strategies

# Demonstrated simultaneous knock-in and multiple knockout in iPSCs

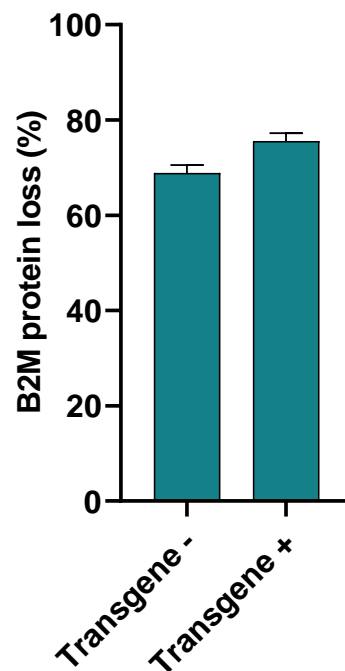
Base editing  
Knockout CIITA

&

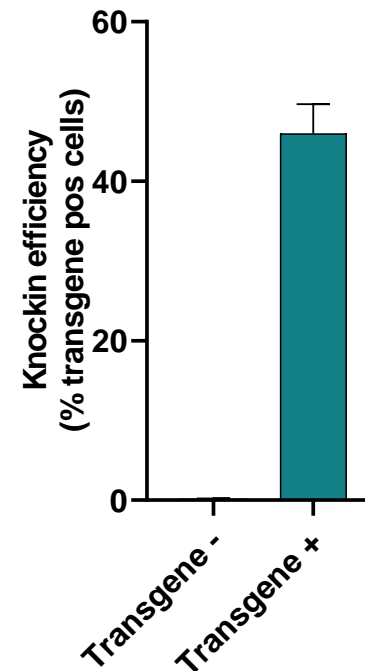
Insertion of a transgene  
in B2M (promoter-less GFP)



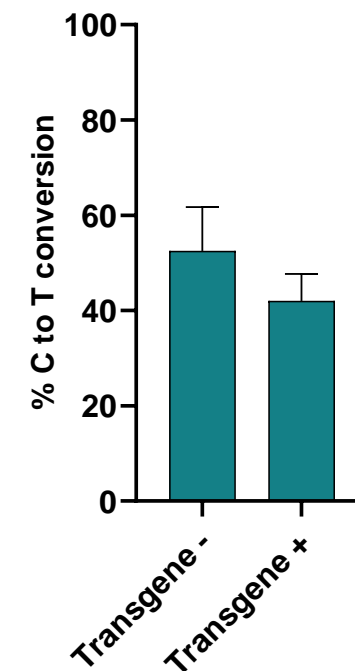
High level  
of B2M  
knockout



High level of  
transgene  
knock-in



... and base  
editing  
(CIITA KO)



The Pin-point™ platform enables one-step simultaneous knock-in and multiple knockout in iPSCs

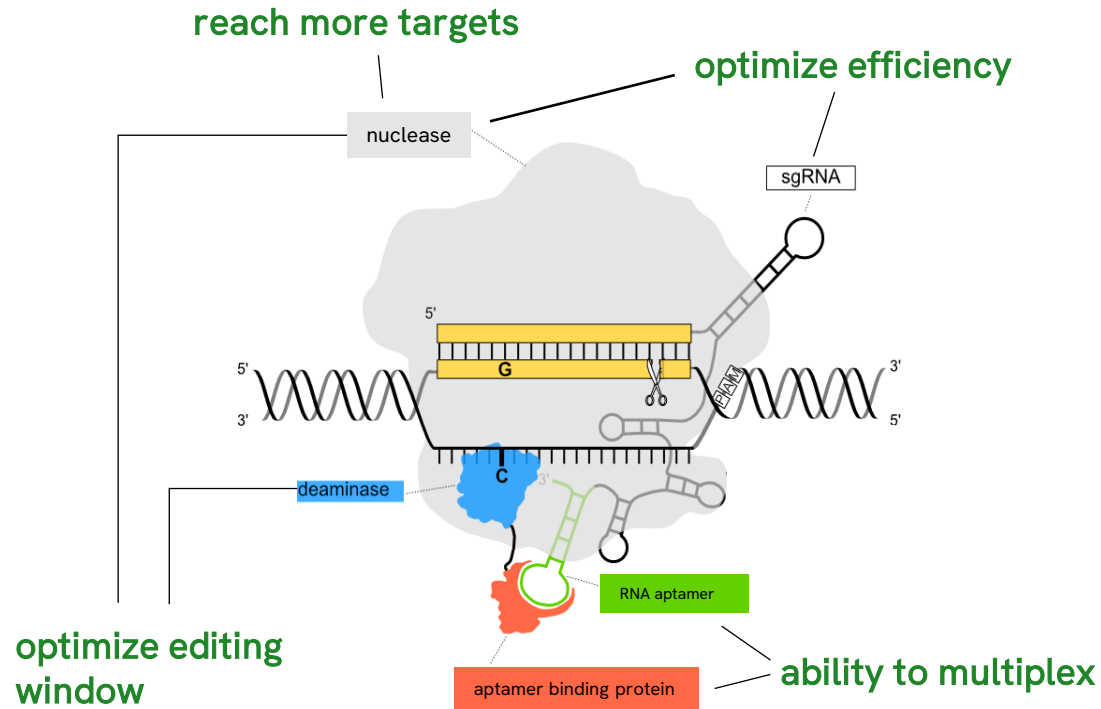




Flexibility for  
target optimization

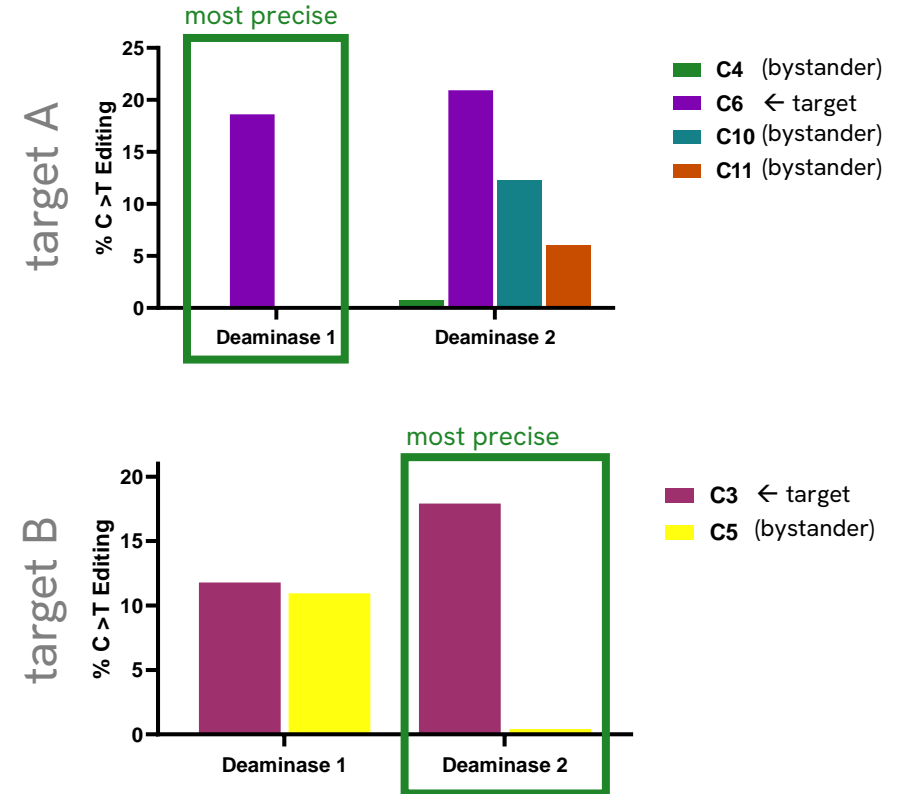
# Choose components for locus-specific optimization

Increased potential to correct more pathogenic SNVs that are not reachable with existing published systems\*



Schematic depicts nCas9 configuration

Example of optimization of the editing window by selecting the best guide RNA and deaminase pairs



The modular Pin-point™ platform can be customized to combine optimal components for a wide range of base editing applications

# A benefit of modularity of the Pin-point™ platform

## Demonstrated compatibility with numerous nucleases

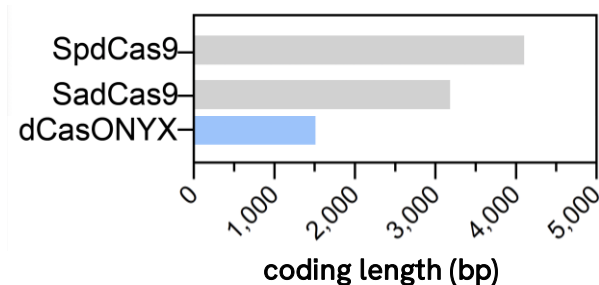
|   | Type II |             |             | Type V      |             |             |             |             |             |             |  |
|---|---------|-------------|-------------|-------------|-------------|-------------|-------------|-------------|-------------|-------------|--|
|   | A       | B           | C           | D           | E           | F           | G           | H           | I           | J           |  |
| Enzyme activity                                   | nickase | nickase     | nickase     | deactivated | deactivated | deactivated | deactivated | deactivated | deactivated | deactivated |  |
| Demonstrated nuclease activity in mammalian cells | ✓       | ✓           | ✓           | ✓           | ✓           | ✓           | ✓           | ✓           | ✓           | ✓           |  |
| Demonstrated with the Pin-point system            | ✓       | ✓           | In progress | ✓           | In progress | ✓           | ✓           | ✓           | ✓           | In progress |  |
| sgRNA optimized                                   | ✓       | In progress |             | In progress |             | ✓           | ✓           | ✓           | ✓           |             |  |
| Enzyme optimized                                  | ✓       |             |             |             |             | ✓           |             |             |             |             |  |
| Confirmed at multiple targets (2+)                | ✓       | In progress |             |             |             | ✓           | ✓           | ✓           | ✓           |             |  |
| Demonstrated in multiple cell types (2+)          | ✓       | In progress |             |             |             | ✓           | In progress | ✓           |             |             |  |
| Demonstrated with multiple deaminases (2+)        | ✓       |             |             |             |             |             | In progress | ✓           | ✓           |             |  |

The Pin-point platform enables utilization of a variety of RNA-guided nucleases, which can be further optimized for editing efficiency

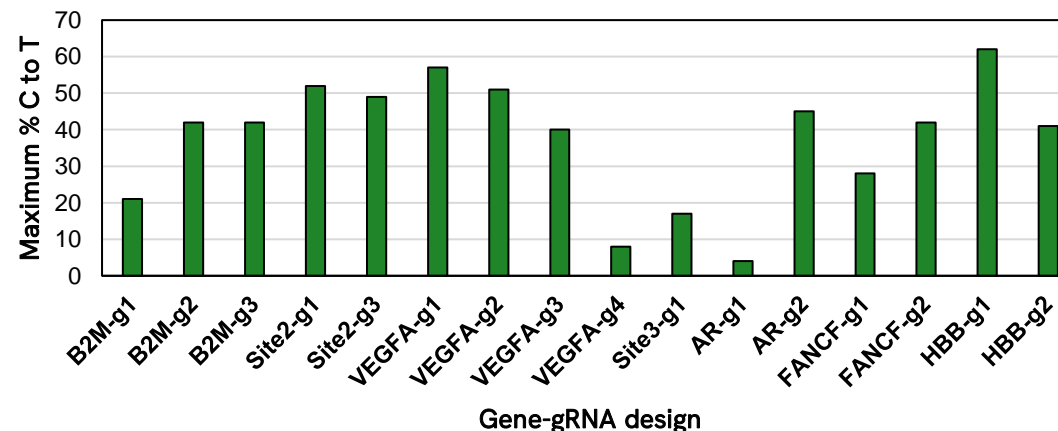
# The Pin-point™ platform configured with Epic Bio's ultracompact Type V effector protein, dCasONYX



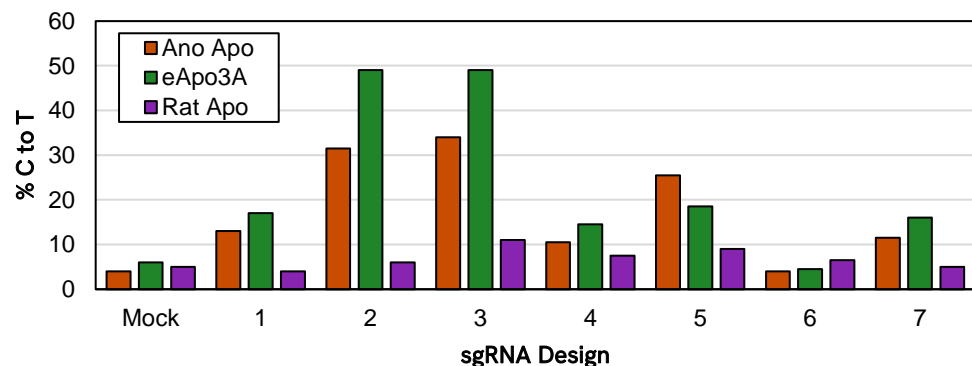
## Ultra-compact engineered Cas12f variant dCasONYX is under 1.5 kb



## Robust targeting capability



## Efficient editing with multiple deaminases when optimal gRNA scaffold is used

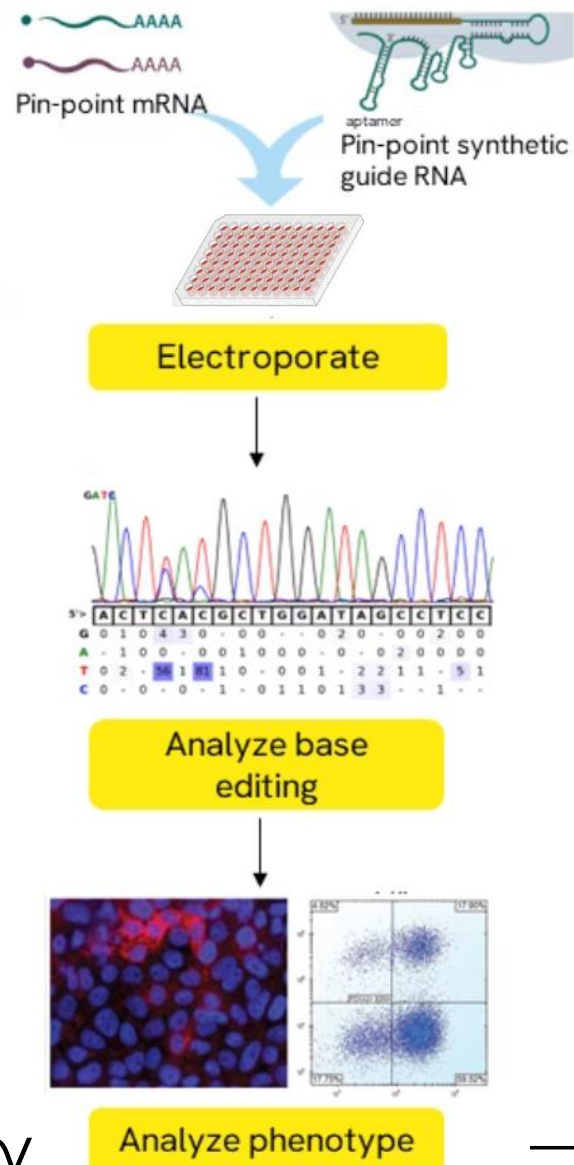


## Additional benefits of dCasONYX

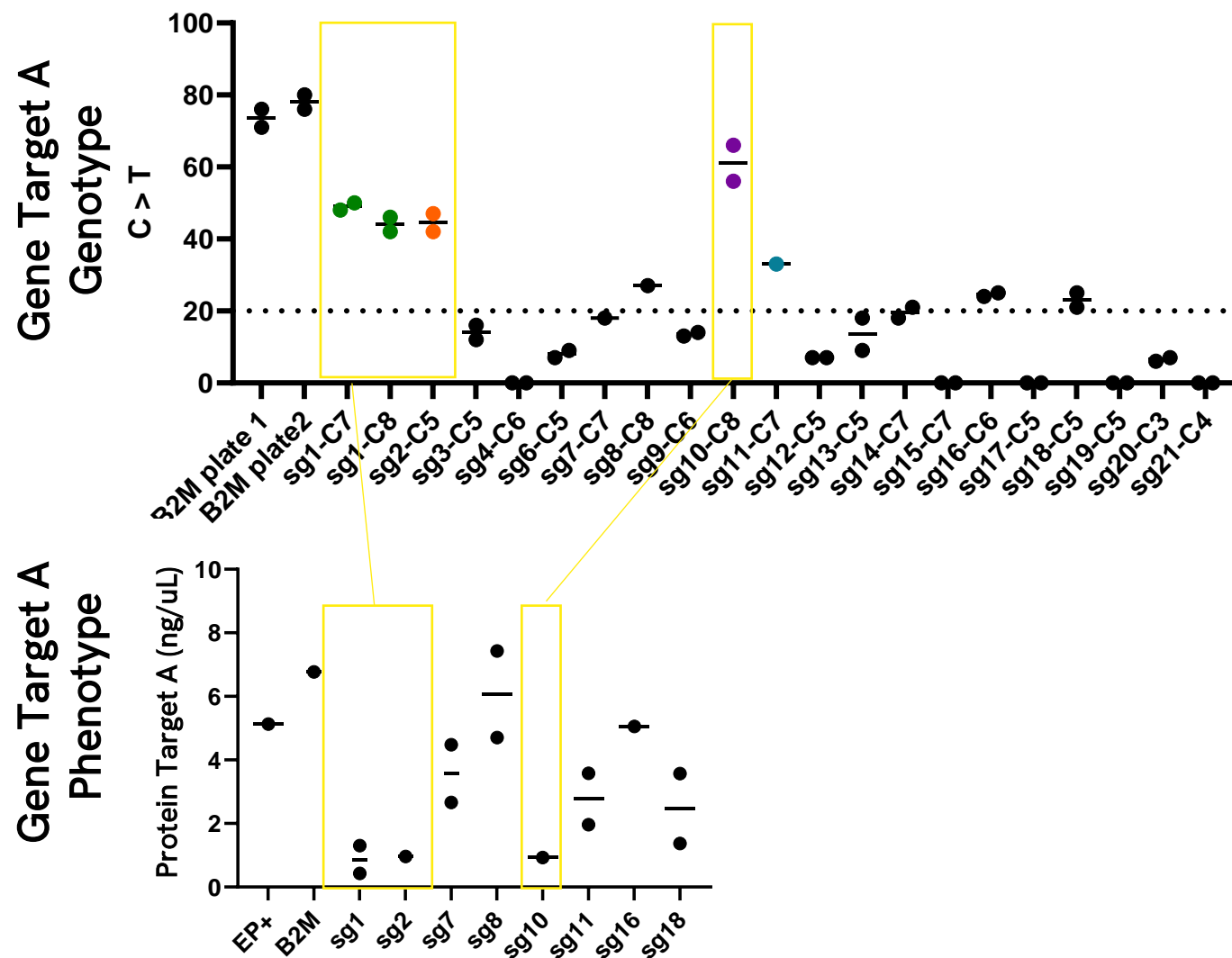
- ✓ **Rapidly advancing to the clinic:** Epic Bio's asset EPI-321 for the potential cure for FSHD
- ✓ **Superior off-target profile:** described in Xin et al. Nature Communications<sup>1</sup>
- ✓ **Fully deactivated nuclease:** base editing without the risk of cutting the DNA
- ✓ **Low immunogenicity:** no prior exposure to dCasONYX in 10 human T cell donors, while 80% of human population have prior exposure to Cas9<sup>2</sup>
- ✓ **Small size:** Coding length less than 1.5 kb ideal for AAV packaging

A dCasONYX Pin-point configuration is one potential alternate to Cas9 for therapeutic applications

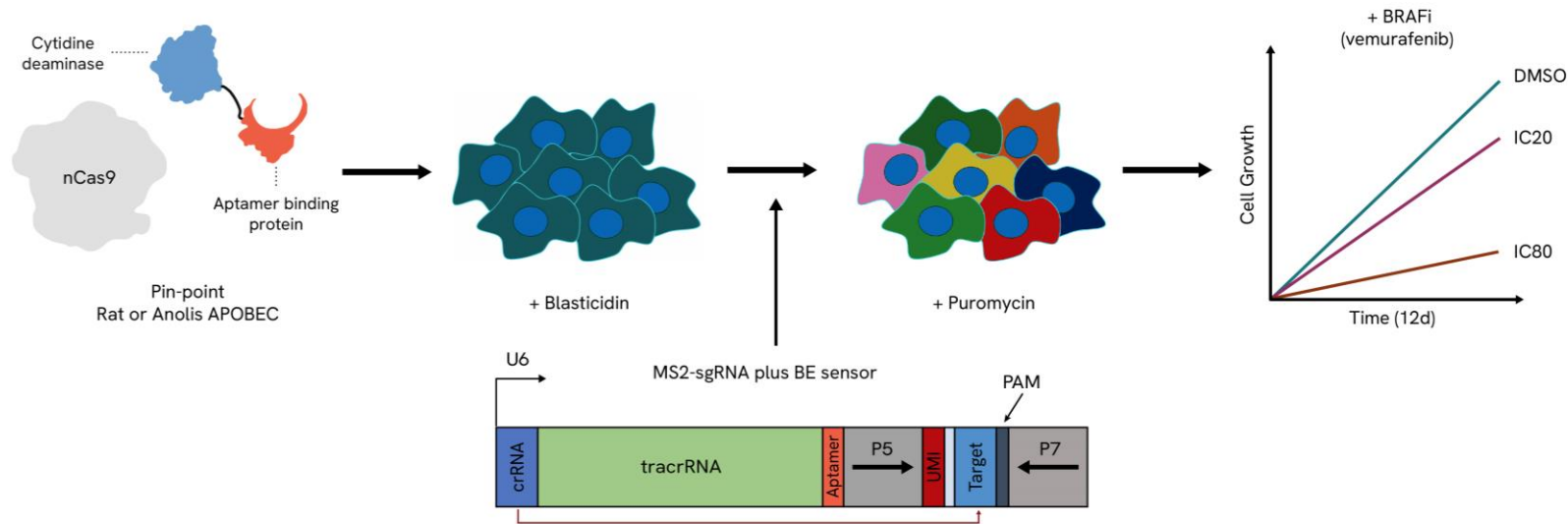
# Synthetic reagents for arrayed screening applications



Confirm results of genotype with phenotype



# Pooled tiled base editing screening services

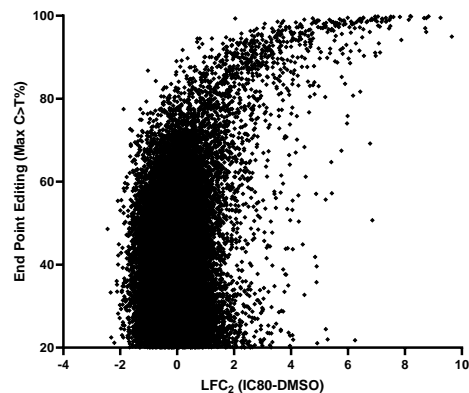


Modify up to every PAM-accessible cytidine in the gene of interest to gain unparalleled understanding of the genotype-phenotype relationship.

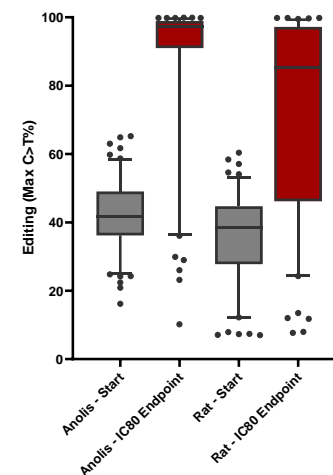
- Target splice sites, introduce premature stop codons, and introduce all possible missense mutations to recapitulate and then go way beyond CRISPRko and CRISPRi screens.
- Include **loss of function** and **gain of function** mutations in a single screen to elucidate possible mechanisms of drug sensitivity and resistance, or protein function.
- Utilize **different deaminases** to maximize the editing window and evaluate more of the genetic sequence.
- Data analysis to generate hit lists and potentially **identify putative causative mutations**.

# Pooled tiled base editing screening services | example BRAFi resistance screen

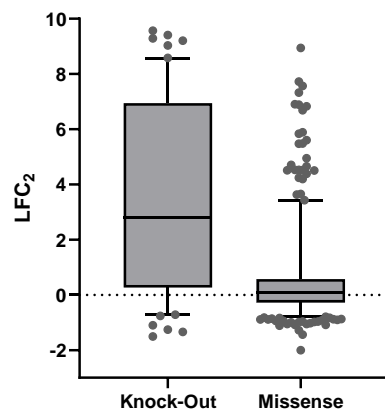
Enrichment of guides that confer resistance to the drug correlate with increased levels of editing



Strong enrichment for edited genomes



NF1 knockout by targeting either the mRNA slice sites or introducing nonsense mutations are more likely to confer drug resistance



resistance hits show no hotspot regions in the NF1 gene, agrees with clinical observations. Multiple guides introducing the same resistance conferring mutation are shown to demonstrate similar screen phenotypes.

