

Pin-point™ Base Editing System: A Versatile Editing Platform Driving Cell Therapies

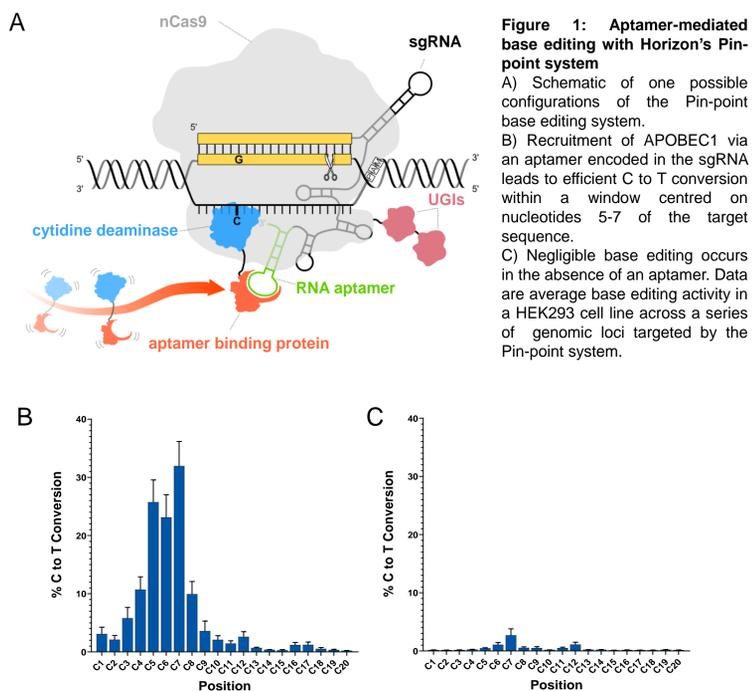
Bronwyn Joubert, Robert Blassberg, Jennifer Harbottle, Olga Mielczarek, Paul Russell, Jesse Stombaugh, Anastasia Lomova Kaufman, John Lambourne, Kevin Hemphill, Matthew Perket, Žaklina Strezoska, Anja van Brabant Smith, Immacolata Porreca

Horizon Discovery, 2650 Crescent Drive, Suite 100, Lafayette, CO 80026, United States | Email address: basediting@horizondiscovery.com

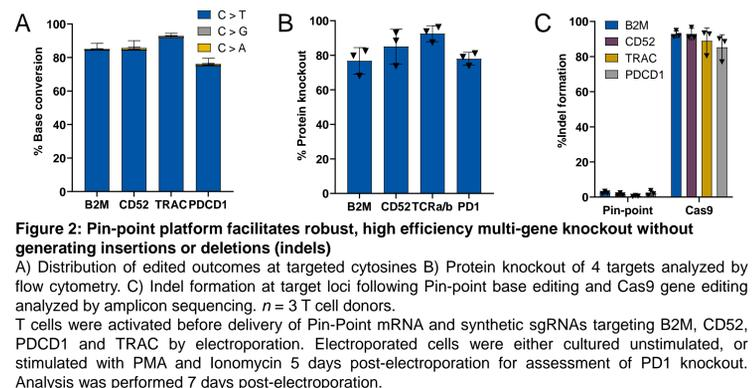
Introduction

Gene editing technologies have successfully been used to improve the next generation of cell and gene therapies. However, standard gene editing platforms such as CRISPR-Cas9 trigger concerning cell toxicity and off-target effects due to the formation of DNA double-strand breaks (DSBs) that could hinder wider clinical applications. Base editing offers an ability to correct disease-causing mutations or knock out genes in a multiplexed manner, without introduction of DSBs. The first cohort of base editing therapeutics entered clinical trials only a few years after the development of the technology. Horizon's modular Pin-point base editing system can use a nickase Cas9 with an aptameric guide RNA to recruit a deaminase to the site of interest, facilitating highly efficient and precise nucleotide conversion. We optimized design and delivery conditions of chemically modified synthetic guide RNAs and enzymatic mRNA to apply multiplex editing with our Pin-point base editing system to the development of engineered CAR-T cells. We target a set of therapeutically relevant loci for the development of allogeneic CAR-T cells achieving high knockout efficiency and editing purity at all sites simultaneously with a safer editing profile and enhanced cell viability compared to traditional nuclease systems. Our technology also enables robust simultaneous targeted knock-in and multiplex knockout without the requirement of additional sequence-targeting components. The ability to perform complex genome editing in multiple cell types (such as T cells, iPSCs, HSCs) safely, efficiently, and precisely opens the door to the application of the Pin-point system in a range of advanced cell therapies.

Horizon's Pin-point base editing system



Multiplex gene editing in T cells



Safety profile of multiplex edited T cells

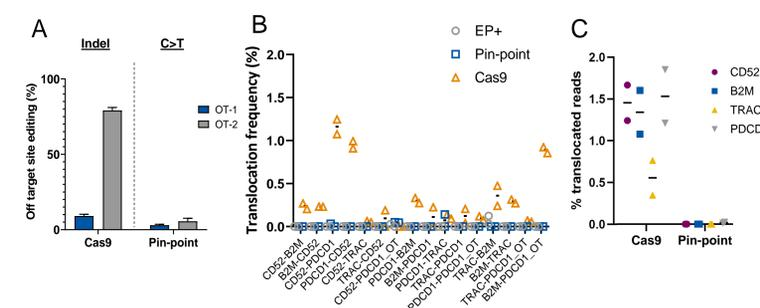
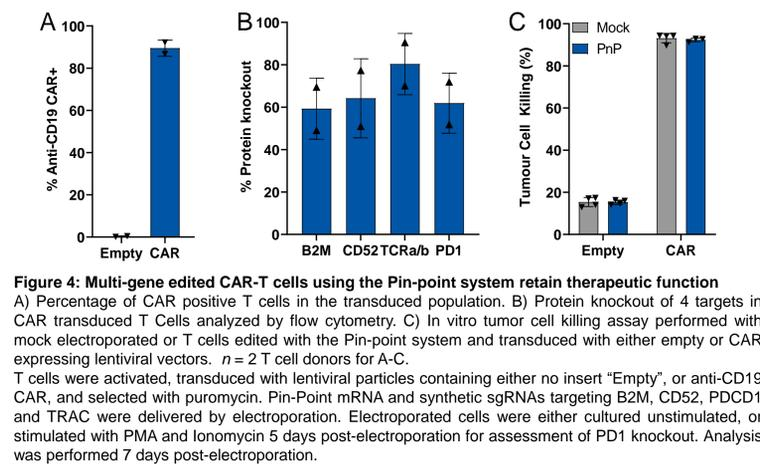
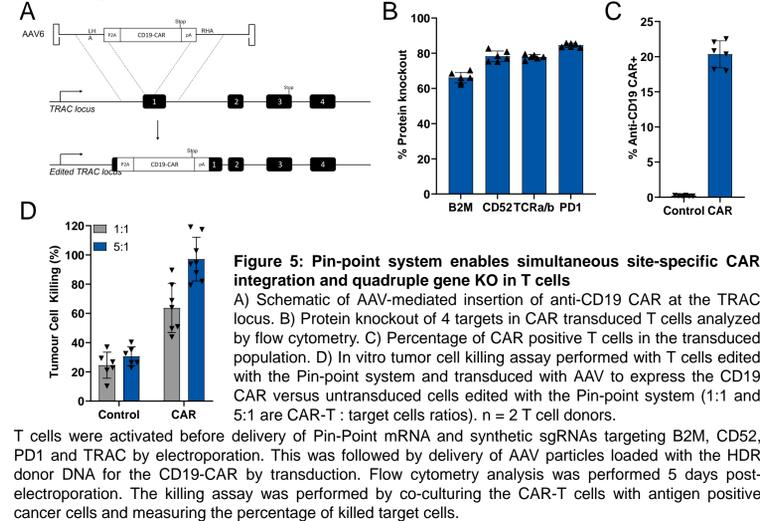


Figure 3: Minimal unintended editing outcomes following multi-gene knockout with the Pin-point system
 A) Off-target editing at two in silico predicted sites following base editing with the Pin-point system and Cas9 gene editing analyzed by amplicon sequencing. B) Frequency of translocations between 4 target sites and an off-target site analyzed by digital droplet PCR. C) Frequency of translocations at each of 4 target sites analyzed by target enrichment sequencing. $n = 3$ T cell donors for A; $n = 2$ T cell donors for B, C.

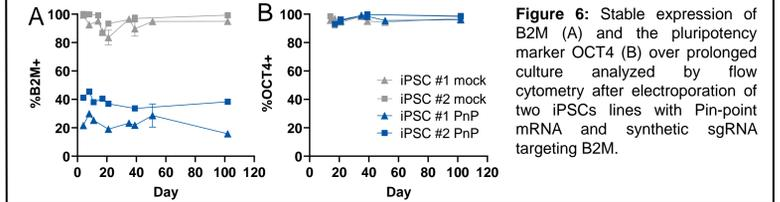
Chimeric Antigen Receptor (CAR)-T cell engineering



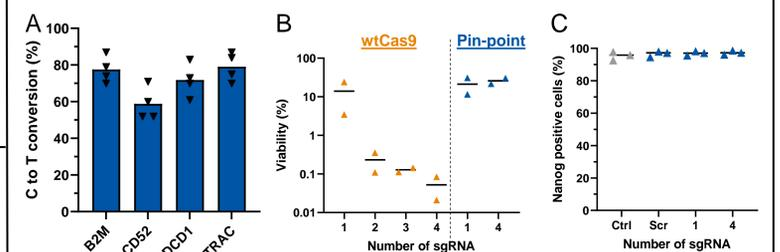
One-step simultaneous knock-in and multiplex knock-out in T cells



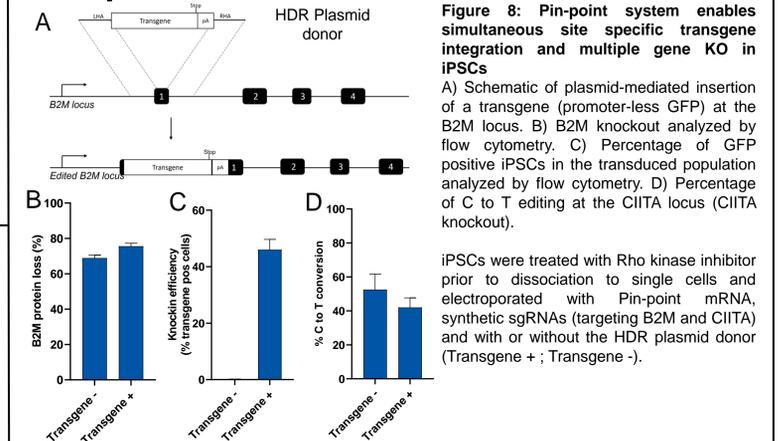
iPSCs edited with the Pin-point system retain pluripotency



Multiplex base editing in iPSCs



One-step simultaneous knock-in and multiplex knock-out in iPSC



Conclusions

- We applied multiplex base editing with the Pin-point system for the development of engineered CAR-T cells and hypoimmunogenic iPSCs.
- Base editing with the Pin-point system achieved greater than 70% knockout efficiency and high purity at therapeutically relevant target sites (B2M, CD52, TRAC and PDCD1) in T cells and iPSCs.
- Multiplex gene editing with the Pin-point platform substantially reduces guide-dependent off target editing and chromosomal translocations compared to Cas9-mediated knockout.
- CAR-T cells retain their cytotoxic activity in vitro following multiplex gene editing with the Pin-point system.
- The Pin-point platform enables simultaneous site-specific transgene knock-in and multi-gene knockout in T cells and iPSCs.
- iPSCs edited with the Pin-point platform retain pluripotency.
- We observed greatly improved viability of iPSCs following multiplex gene editing with the Pin-point system compared to Cas9.

References

Collantes et al. The CRISPR Journal. Feb 2021.58-68