

Optimizing non-viral genome editing workflows for induced pluripotent stem cells with advanced electroporation technology

Olga Cohen, Shahan Molla, Gregory B. Bonello, Arseny Smirnov, Namritha Ravinder

Thermo Fisher Scientific, 5781 Van Allen Way, Carlsbad, CA 92008, USA

Key Takeaways

- The development of versatile and efficient non-viral based cell engineering workflows using electroporation (EP) delivery method has significant implications in cell and gene therapy field.
- The Neon NxT Electroporation System with 8-Channel Pipette enables rapid and efficient assessment of electroporation condition to achieve high genome editing efficiency of cells of interest.
- This study highlights the importance of optimizing electroporation conditions to achieve optimal knock-in efficiency balanced with superior CAR-iPSC recovery with CultureCEPT™ supplement.
- Our findings demonstrate the potential of CultureCEPT™ supplement in cell engineering applications.

Material & Methods

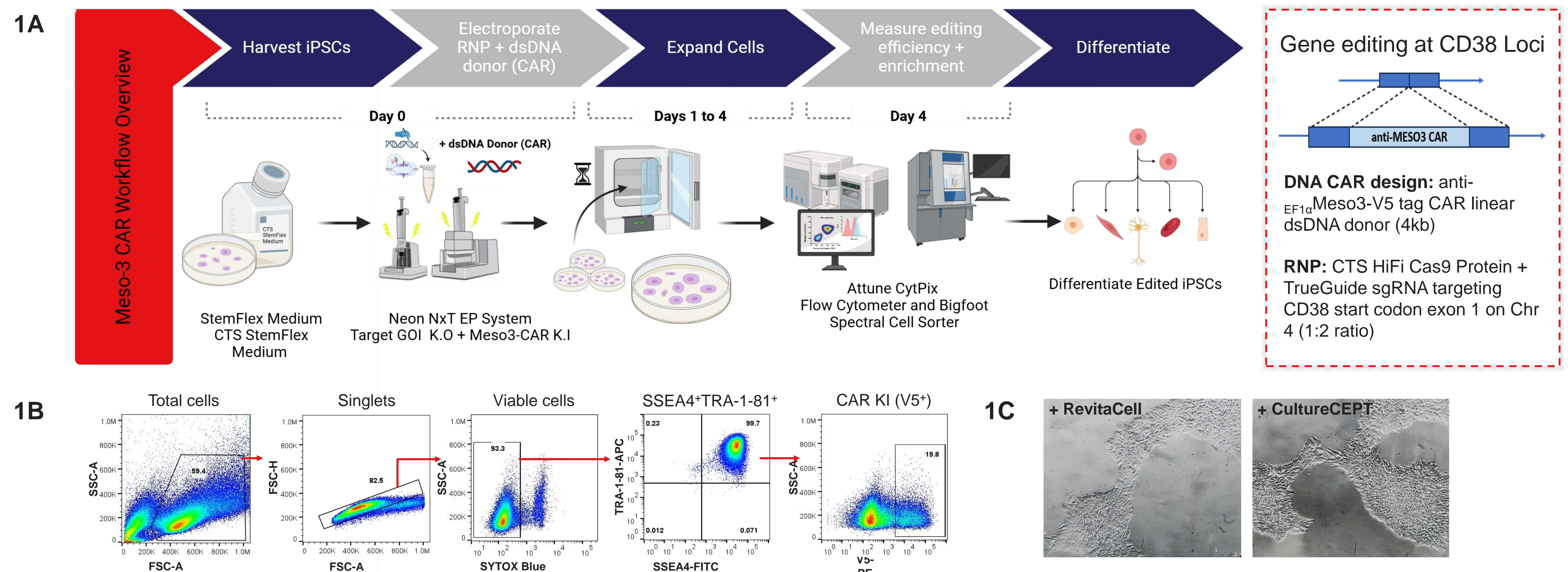
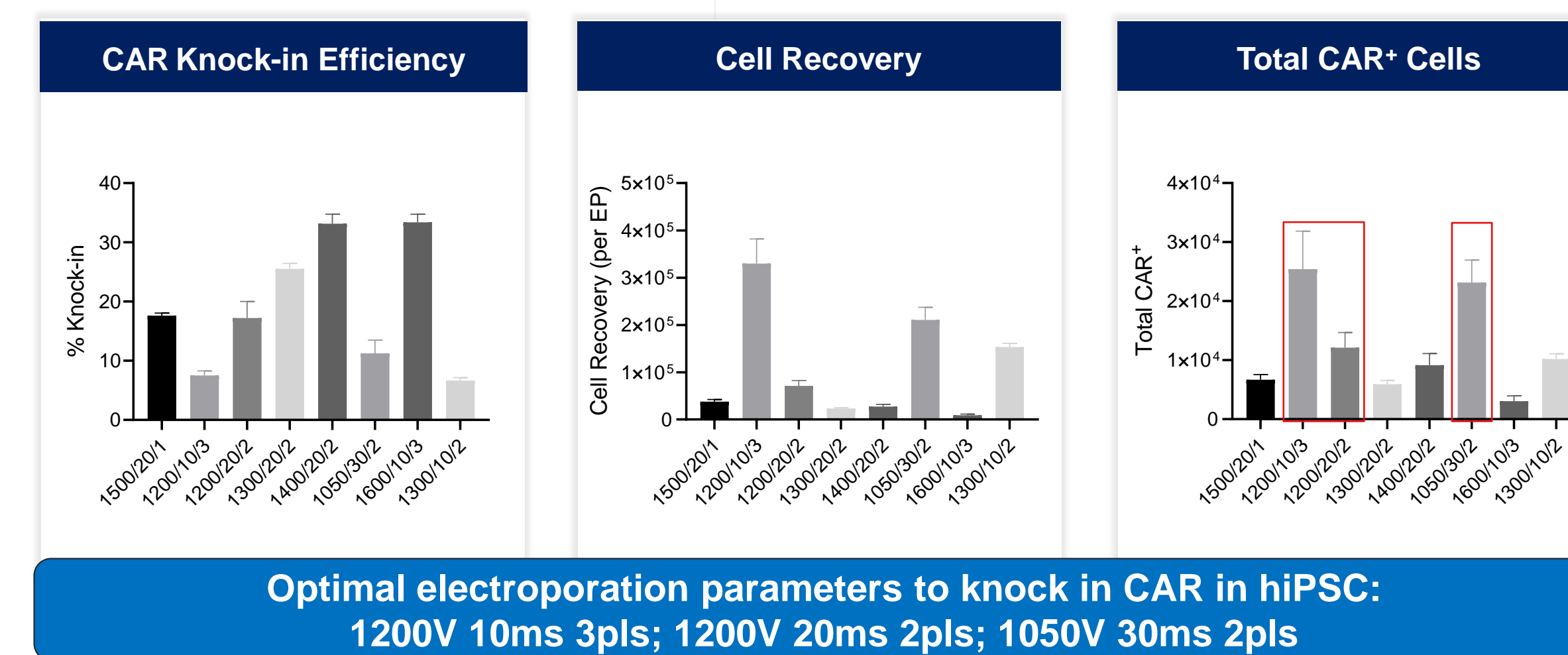


Figure 1. iPSC cell engineering protocol workflow diagram

A. Human iPSCs (TFSi001-A) were grown in Gibco™ CTS™ StemFlex™ Medium (cat#A5465001) in Gibco™ CTS™ Vitronectin (cat#CTS279S3)-coated culture dishes. At the confluency of 70-80%, the cells were harvested using CTS™ Versene™ Solution (cat#A4239101), washed once with Gibco CTS DPBS (cat#A1285801) to prepare for electroporation. Genome editing was performed using CTS HiFi Cas9, TrueGuide™ Synthetic Guide RNA target CD38 loci and anti- $EF1\alpha$ Meso3-CAR (V5-tagged) dsDNA payloads. Neon™ NxT 8 channel EP system and 10 μ l tips were used to deliver gene editing payloads to 10e6 cells/mL in R buffer. After electroporation, cells were cultured in the same condition medium supplemented with Gibco™ CTS™ RevitaCell™ Supplement (cat#A4238401) for 24 hours. Next day the medium was replaced without RevitaCell™ Supplement and cells were maintained in culture for three more days. **B.** 4 days post EP cells were analyzed for viability, recovery and anti-Meso3-CAR knock-in efficiency (%V5 positive cells) using Attune™ CytPix™ Flow Cytometer. SYTOX™ Blue Dead Cell Stain (cat#S34857) was used to measure cell viability. The iPSCs were stained with anti-V5 tag-PE (cat#12-6796-42) to quantify CAR knock-in efficiency and with anti-TRA-1-81-APC (cat#17-8883-42) and anti-SSEA4-FITC (cat#53-8843-42) antibodies to confirm pluripotency. The cells were stained in eBioscience™ Flow Cytometry Staining Buffer (cat#00-4222-26) at 4°C for 40 minutes prior to flow cytometry. **C.** Representative images of CAR-iPSC recovered in presence of CultureCEPT or RevitaCell. Images were taken on day 4 post EP using EVOS™ M5000 Imaging System.

Results

2A. Neon NxT electroporation program optimization



2B. Neon NxT electroporation buffer comparison

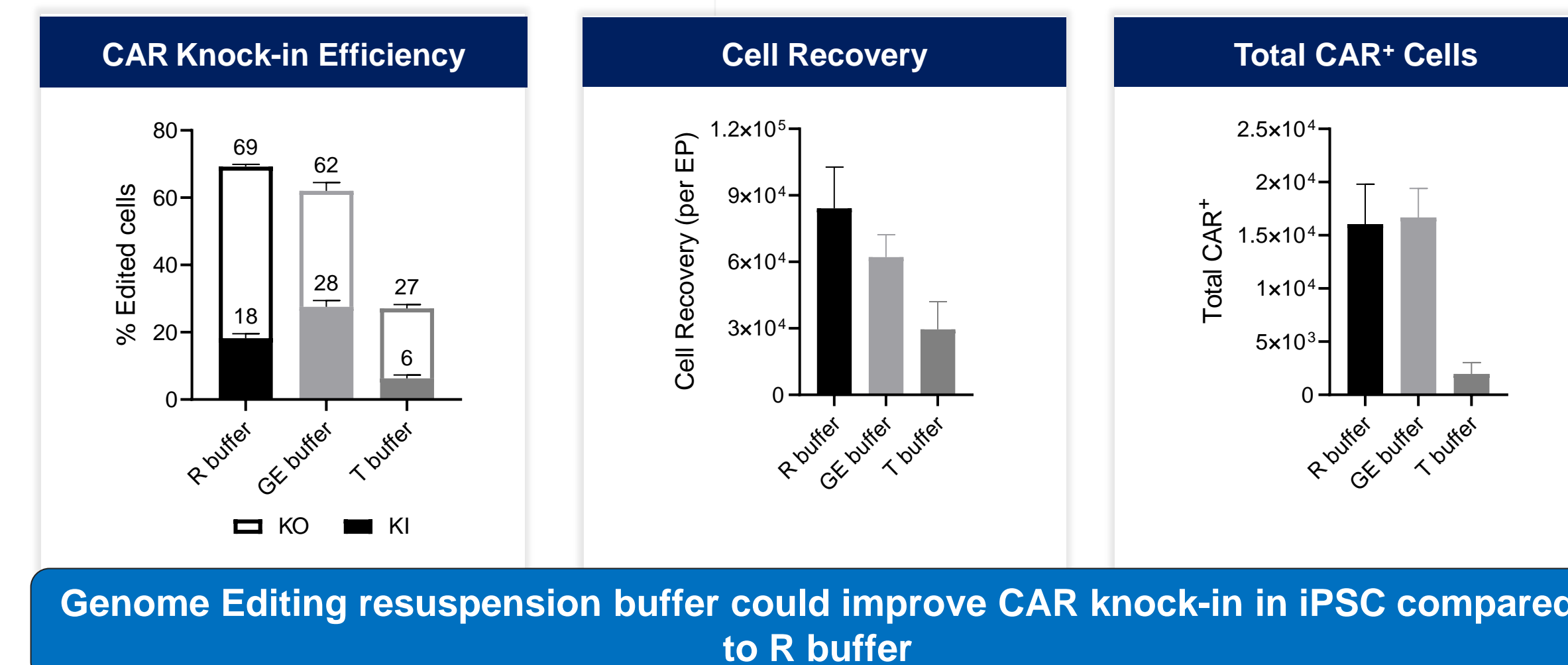


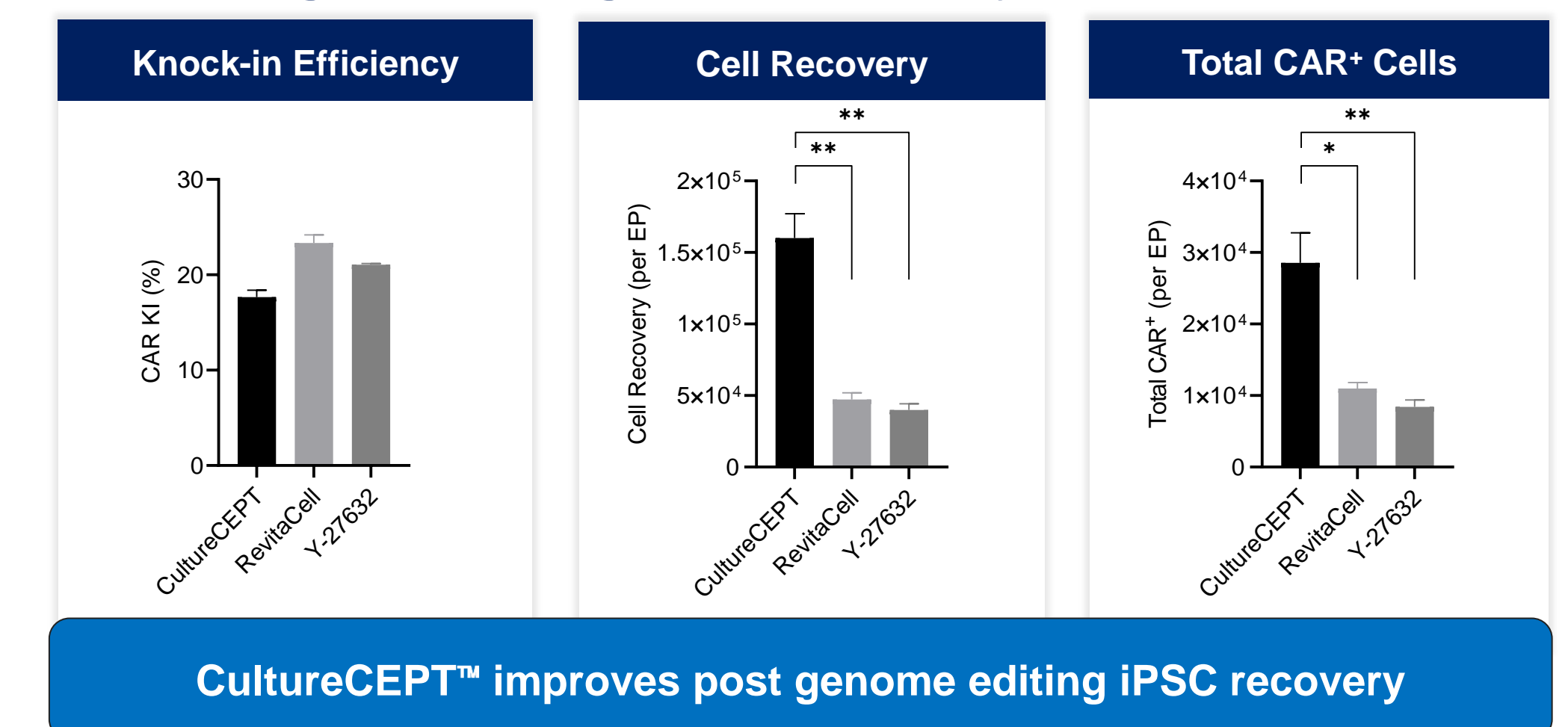
Figure 2. Optimization of electroporation conditions for CAR knock-in in hiPSC

A. Several preselected electroporation parameters were assessed to evaluate CAR knock-in efficiency, total viable cell number and total CAR-iPSC positive cells recovered four days post electroporation. 10e6/mL cells were electroporated with Cas9 RNP + 50 μ g/mL anti-Meso3 CAR payload using 10 μ l tips and R buffer. **B.** Different Neon NxT buffers (R, GE and T) were evaluated at selected EP condition (1200V 20ms 2pls) and post EP, % editing efficiency, total viable cell number and total CAR-iPSC positive cells were analyzed ($n \geq 3$). **C.** Post genome editing cell recovery in presence of CultureCEPT™ vs Y-27632 or RevitaCell. The analysis represents CAR knock-in efficiency, total viable cell number and total CAR-iPSC positive cells recovered on day 4 post electroporation. CAR knock-in was performed using 1200V 20ms 2pls electroporation parameters in GE buffer. Electroporated cells were cultured in Gibco™ CTS™ StemFlex™ Medium supplemented with either Gibco™ CultureCEPT™, CTS™ RevitaCell™ Supplement or ROCK inhibitor Y-27632 from another vendor for 24 hours ($n \geq 3$). Cells recovered using media supplemented with CultureCEPT showed better recovery and Total CAR positive cells. **D.** Characterization of CAR-iPSCs after clonal selection with KaryoStat assay to assess genomic stability and by PluriTest assay to confirm CAR-iPSC pluripotency and genetic background.

Conclusions

The Neon NxT Electroporation System with 8-channel pipette demonstrates remarkable versatility in rapidly optimizing electroporation conditions ensuring consistent and efficient workflows. The ability of Neon NxT Electroporation System to achieve high editing efficiency along with efficient cell recovery in presence of CultureCEPT™ supplement underscores the system's robustness and reliability for non-viral CAR-iPSC engineering applications.

2C. Post gene editing iPSC recovery with CultureCEPT



2D. Characterization of CAR-iPSCs

