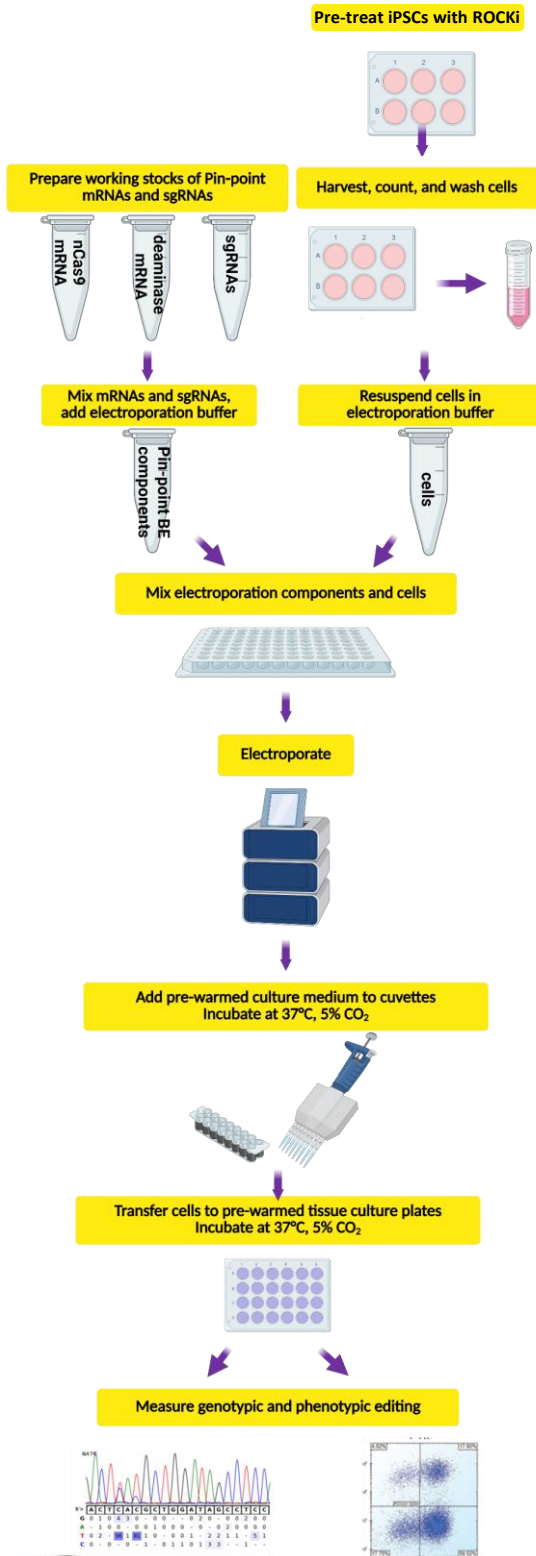


# Pin-point™ cytidine base editing (CBE) platform

## Short protocol for electroporation of human induced pluripotent stem cells (iPSCs) using the Lonza 4D-Nucleofector® System

### Experimental workflow:



Created with BioRender

The following is a protocol for delivering unmodified [Pin-point CBE nCas9 mRNA](#) (Cat # PNP12744, PNP12746, and PNP12748), [Pin-point CBE rAPOBEC deaminase mRNA](#) (Cat # PNP12745, PNP12747, and PNP12749), and Pin-point CBE sgRNAs ([validated controls](#) or [custom](#)) to stimulated human iPSCs using the Lonza 4D-Nucleofector System with the P3 Primary Cell 96-well Nucleofector® Kit. For more details, please refer to the [Pin-point platform CBE technical manual](#). (EP = electroporation)

Day -3	<b>Plate cells</b>	Coat plasticware with appropriate iPSC culture matrix (e.g Vitronectin XF). Seed cells at appropriate density to achieve ~70% confluence at Day 0. <i>NOTE: We recommend mTeSR™ PLUS culture medium (STEMCELL Technologies)</i>			
	<b>Prepare post-EP plates &amp; treat iPSCs with ROCKi</b>	2 hours prior to electroporation replace culture media with iPSC culture medium + Y-27632 (ROCKi) (10 µM). Coat 12-well plates with iPSC culture matrix. Add iPSC cell culture medium + ROCKi (10 µM) to matrix-coated plates: 1 mL per well of 12-well plate. Incubate at 37°C and 5% CO <sub>2</sub> .			
Day 0	<b>Prepare Pin-point base editing components</b>	Prepare working stock solutions of mRNAs and sgRNAs on ice according to the table below.			
	<b>Prepare the cells</b>	Remove iPSC culture media and rinse cells twice with PBS. Dissociate iPSC colonies by incubation at 37°C for 6-10 mins with Accutase (Gibco). Add 1 mL iPSC culture medium to dissociated iPSC colonies and gently triturate to single cells by pipetting. Count cells and transfer the desired number into a 15 mL conical tube: 2 x 10 <sup>5</sup> cells/ 20 µL electroporation. Centrifuge at 200 x g for 5 mins. Resuspend cell pellet in P3 nucleofector solution at a final concentration of 1 x 10 <sup>7</sup> cells/mL.			
	<b>Mix EP components</b>	Gently mix mRNAs + sgRNAs + cells:			
Day 1	<b>Change media</b>	Remove iPSC culture media containing ROCKi and add fresh iPSC culture medium: 2 mL per well of 12-well plate. Incubate at 37°C and 5% CO <sub>2</sub> .			
	<b>Post-EP analysis</b>	Proceed with desired genotypic (Sanger sequencing) and/or phenotypic (flow cytometry) analyses of base editing levels.			

	Pin-point component	Working stock*	Final amount per 20 µL EP	Volume per 20 µL EP
<b>mRNA and sgRNA mix</b>	CBE nCas9 mRNA	2 µg/µL	2.56 µg	1.28 µL
	CBE rAPOBEC deaminase mRNA	2 µg/µL	0.74 µg	0.37 µL
<b>Cells</b>	Synthetic sgRNA	200 µM	40 pmol	0.2 µL
	P3 nucleofector solution	-	-	to 10 µL
	Cells in P3 nucleofector solution	-	2 x 10 <sup>5</sup> cells	10 µL

\* CBE nCas9 mRNA and CBE deaminase mRNA shipped at 2 µg/µL

If you have questions or comments, please reach out to [Scientific Support](#).

This product contains material of biological origin. Use for research purposes only. Do not use in humans or for diagnostic purposes. The purchaser assumes all risk and responsibility concerning reception, handling and storage.