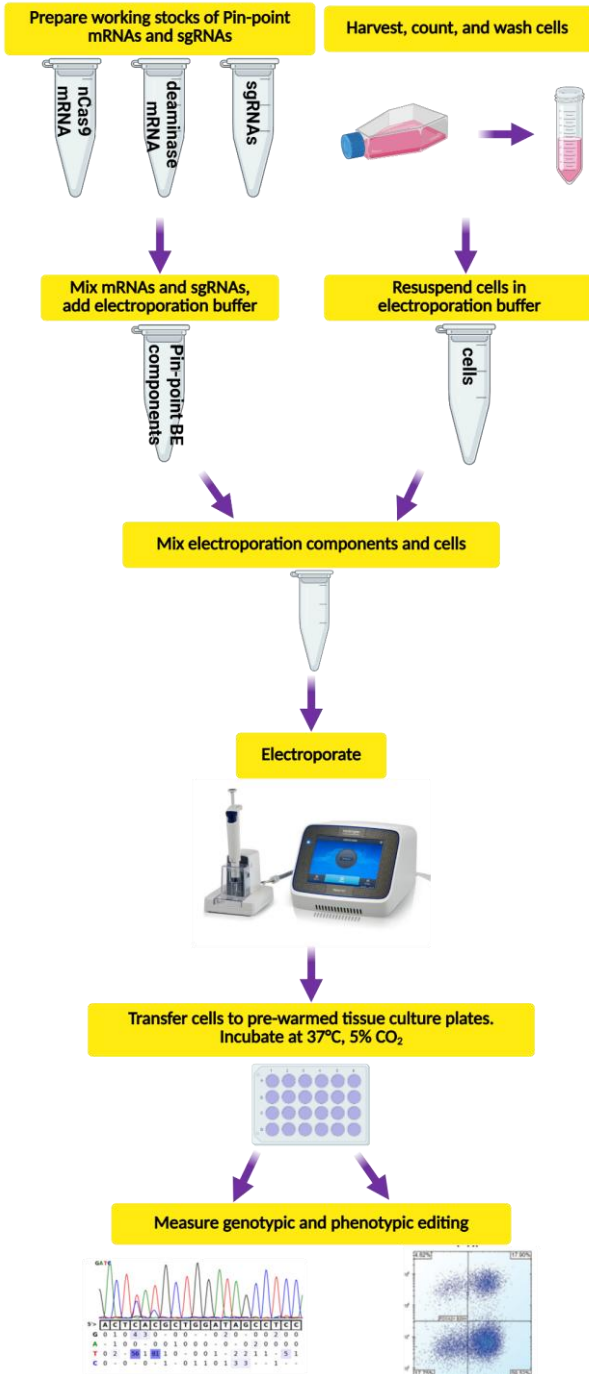


# Pin-point™ adenine base editing (ABE) platform

## Short protocol for electroporation of human hematopoietic stem and progenitor cells (HSPCs) using the Neon™ Nxt Electroporation System

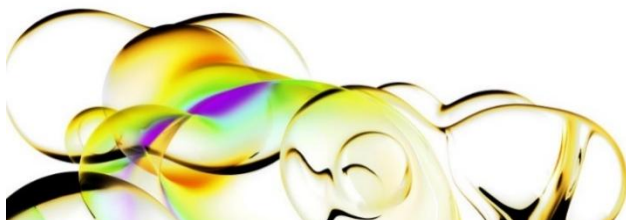
### Experimental workflow:



Created with BioRender

The following is a protocol for delivering 5moU-modified [Pin-point ABE nCas9 mRNA](#) (Cat # PNP12744, PNP12746, PNP12748), [Pin-point ABE-exact deaminase mRNA](#) (Cat # PNP13785, PNP13786, PNP13787), or [Pin-point ABE-flex deaminase mRNA](#) (Cat # PNP13777, PNP13778, PNP13779), and Pin-point sgRNAs ([validated controls](#) or [custom](#)) to stimulated human HSPCs using the Neon Nxt Electroporation System. For more details, please refer to the [Pin-point platform ABE technical manual](#). (EP = electroporation)

Day -1	<b>Plate cells</b>	Thaw cryopreserved cells and seed at 200,000 cells/mL in appropriate culture medium. We recommend StemSpan™ SFEM II medium supplemented with StemSpan CD34+ expansion supplement, 1 $\mu$ M SR1, 1000 nM UM729 and Pen/Step.			
	<b>Prepare post-EP plates</b>	Add complete antibiotic-free cell culture medium to plates: 200 $\mu$ L in 96-well plate or 500 $\mu$ L in 24-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>	Harvest and count the cells. Transfer the desired number of cells for electroporation into a centrifuge tube. Wash with PBS, centrifuge at 200 x g for 10 minutes. Resuspend cell pellet in buffer R at a final concentration of 1 x 10 <sup>7</sup> cells/mL.			
	<b>Mix EP components</b>	Gently mix mRNAs + sgRNAs + cells:			
			<b>Pin-point component</b>	<b>Working stock*</b>	<b>Final amount per 10 <math>\mu</math>L EP</b>
<b>mRNA and sgRNA mix</b>		ABE nCas9 mRNA	2 $\mu$ g/ $\mu$ L	1.5 $\mu$ g	0.75 $\mu$ L
		ABE deaminase mRNA	2 $\mu$ g/ $\mu$ L ABE-exact, 0.8 $\mu$ g/ $\mu$ L ABE-flex*	1 $\mu$ g ABE-exact, 0.4 $\mu$ g ABE-flex	0.5 $\mu$ L
		synthetic sgRNA	200 $\mu$ M	6.25 $\mu$ M	0.312 $\mu$ L
	Buffer R	-	-	to 5 $\mu$ L	
<b>Cells</b>	Cells in buffer R	-	50,000 cells	5 $\mu$ L	
* ABE nCas9 mRNA and ABE deaminase mRNA shipped at 2 $\mu$ g/ $\mu$ L					
	<b>EP</b>	Electroporate 10 $\mu$ L on Neon Nxt Electroporation System using the following conditions: 1400V, 10ms, 3 pulses. Pipette cells into prepared plates. Incubate at 37°C, 5% CO <sub>2</sub> for 48-72 hours.			
Days 3 - 7	<b>Post-EP analysis</b>	Proceed with desired genotypic (Sanger sequencing) and/or phenotypic (flow cytometry or qPCR) analyses of base editing and protein expression levels.			



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.