



The following is a protocol for delivering unmodified Pin-point nCas9 mRNA (Cat # PNP12744, PNP12746, PNP12748), Pin-point rat APOBEC mRNA (Cat # PNP12745, PNP12747, PNP12749), and Pin-point sgRNAs (Cat # PNP-02000-01, PNP-02200-01, PNP-02300-01) to cultured mammalian cells using the Neon™ NxT Electroporation System. For more details, please see [full protocol](#).

This protocol is written for electroporation of HEK293T cells and activated primary human T cells.

Day -2/-1	Plate cells	Seed cells at appropriate density. HEK293T cells: 3 x 10 ⁶ cells/10cm dish; Activated T cells: 1 x 10 ⁶ cells/mL																																															
Day 0	Prepare post-electroporation plates	Add appropriate cell culture medium to plates HEK293T cells: 96-well plates with 100µL medium per well; Activated T cells: 24-well plates with 500µL medium per well. Incubate at 37°C and 5% CO ₂																																															
	Prepare Pin-point base editing components	Prepare working stock solutions of mRNAs and sgRNAs according to the table below																																															
	Prepare the cells	Harvest and count the cells. Transfer the desired number of cells for electroporation into a centrifuge tube. Wash with PBS (avoid centrifuge speeds > 400 x g) Resuspend cell pellet in R buffer: HEK293T cells: 1 x 10 ⁷ cells/mL; Activated T cells: 5 x 10 ⁷ cells/mL.																																															
	Mix electroporation components	Gently mix mRNAs + sgRNAs + cells HEK293T cells: <table border="1" style="display: inline-table; margin-right: 20px;"> <thead> <tr> <th>Pin-point component</th> <th>Working stock*</th> <th>Final amount per reaction</th> <th>Volume per reaction</th> </tr> </thead> <tbody> <tr> <td>nCas9 mRNA</td> <td>2 µg/µL</td> <td>1 µg</td> <td>0.5 µL</td> </tr> <tr> <td>Rat APOBEC mRNA</td> <td>200 ng/µL</td> <td>100 ng</td> <td>0.5 µL</td> </tr> <tr> <td>synthetic sgRNA</td> <td>100 µM</td> <td>6 µM</td> <td>0.6 µL</td> </tr> <tr> <td>Buffer R</td> <td>-</td> <td>-</td> <td>3.4 µL</td> </tr> <tr> <td>Cells</td> <td></td> <td>50,000 cells</td> <td>5 µL</td> </tr> </tbody> </table> Activated T cells: <table border="1" style="display: inline-table;"> <thead> <tr> <th>Pin-point component</th> <th>Working stock*</th> <th>Final amount per reaction</th> <th>Volume per reaction</th> </tr> </thead> <tbody> <tr> <td>nCas9 mRNA</td> <td>2 µg/µL</td> <td>1.56 µg</td> <td>0.78 µL</td> </tr> <tr> <td>Rat APOBEC mRNA</td> <td>1 µg/µL</td> <td>0.222 µg</td> <td>0.222 µL</td> </tr> <tr> <td>synthetic sgRNA</td> <td>200 µM</td> <td>2 - 6µM</td> <td>0.1 µL of each sgRNA (0.3µL total when using 3 sgRNAs)</td> </tr> <tr> <td>Buffer R</td> <td>-</td> <td>-</td> <td>to 10 µL</td> </tr> <tr> <td>Cells</td> <td></td> <td>250,000 cells</td> <td>5 µL</td> </tr> </tbody> </table>	Pin-point component	Working stock*	Final amount per reaction	Volume per reaction	nCas9 mRNA	2 µg/µL	1 µg	0.5 µL	Rat APOBEC mRNA	200 ng/µL	100 ng	0.5 µL	synthetic sgRNA	100 µM	6 µM	0.6 µL	Buffer R	-	-	3.4 µL	Cells		50,000 cells	5 µL	Pin-point component	Working stock*	Final amount per reaction	Volume per reaction	nCas9 mRNA	2 µg/µL	1.56 µg	0.78 µL	Rat APOBEC mRNA	1 µg/µL	0.222 µg	0.222 µL	synthetic sgRNA	200 µM	2 - 6µM	0.1 µL of each sgRNA (0.3µL total when using 3 sgRNAs)	Buffer R	-	-	to 10 µL	Cells		250,000 cells
Pin-point component	Working stock*	Final amount per reaction	Volume per reaction																																														
nCas9 mRNA	2 µg/µL	1 µg	0.5 µL																																														
Rat APOBEC mRNA	200 ng/µL	100 ng	0.5 µL																																														
synthetic sgRNA	100 µM	6 µM	0.6 µL																																														
Buffer R	-	-	3.4 µL																																														
Cells		50,000 cells	5 µL																																														
Pin-point component	Working stock*	Final amount per reaction	Volume per reaction																																														
nCas9 mRNA	2 µg/µL	1.56 µg	0.78 µL																																														
Rat APOBEC mRNA	1 µg/µL	0.222 µg	0.222 µL																																														
synthetic sgRNA	200 µM	2 - 6µM	0.1 µL of each sgRNA (0.3µL total when using 3 sgRNAs)																																														
Buffer R	-	-	to 10 µL																																														
Cells		250,000 cells	5 µL																																														
	Electroporate	Electroporate on Neon NxT Electroporation System with the following conditions: HEK293T cells: 1150V, 20ms, 2 pulses; Activated T cells: 1600V, 10ms, 3 pulses. Pipette cells into prepared plates. Incubate at 37°C, 5% CO ₂ for 48-72 hours.																																															
Days 3-7	Post-electroporation analysis	Proceed with desired genotypic (Sanger sequencing) and phenotypic (flow cytometry) analyses of base editing 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.

