

CRISPR/Cas9-mediated Genome Editing of Human Primary T-cells using the Neon™ Transfection System

Pub. No. MAN0017851 Rev. A.0

Note: Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Overview

This protocol describes the isolation and activation conditions for human primary T-cells, as well as the procedure for the delivery of Cas9/gRNA (guide RNA) complexes via electroporation using the Neon™ Transfection System to T-cells.

Required materials not supplied

All materials are available through thermofisher.com.

Table 1 Materials required for T-cell isolation, activation, and CRISPR-Cas9 genome editing using the Neon™ Transfection System

Item	Catalog No.
Dynabeads™ Untouched™ Human T Cells Kit	11344D
CTS™ OpTmizer™ T Cell Expansion SFM	A1048501
CTS™ Immune Cell SR	A2596101
IL2 (Interleukin 2) CTS™ Recombinant Human Protein	CTP0023
GlutaMAX™ Supplement	35050061
Dynabeads™ Human T-Activator CD3/CD28	11131D
DPBS, no calcium, no magnesium	14190144
TrueCut™ Cas9 Protein v2	A36497
GeneArt™ Precision gRNA Synthesis Kit	A29377
Neon™ Transfection System	MPK5000
Neon™ Transfection System 10-µL Kit	MPK1025, MPK1096
<i>(Optional)</i> GeneArt™ Genomic Cleavage Detection Kit	A24372

Isolate and activate T-cells

1. Isolate the T-cells from peripheral blood mononuclear cells (PBMC) derived from healthy donors using the Dynabeads™ Untouched™ Human T Cells Kit.
2. Activate the T-cells (at 1×10^6 cells/mL) with Dynabeads™ Human T-Activator CD3/CD28 in CTS™ OpTmizer™ T Cell Expansion SFM medium containing 100 U/mL IL2 CTS™ Recombinant Human Protein, 6 mM GlutaMAX™ Supplement, and 2% CTS™ Immune Cell SR or human AB serum.
3. Culture the T-cells in a humidified 37°C, 5% CO₂ incubator for 3 days before electroporation.

Prepare CRISPR-Cas9/gRNA complex

1. Generate the gRNA using the GeneArt™ Precision gRNA Synthesis Kit, or order TrueGuide™ Synthetic gRNA at thermofisher.com/trueguide.

2. Add 1 µg of TrueCut™ Cas9 Protein v2 and 250 ng of gRNA to 5 µL of Resuspension Buffer R. Mix well gently.

Note: Use high concentration Cas9 protein (5 µg/µL) and gRNA to keep the volume of Cas9/gRNA complex at less than 10% of total reaction volume (e.g., 1 µL of Cas9 protein + gRNA in 10 µL total reaction volume).

Note: The amounts of reaction components are for a single electroporation using the 10-µL Neon™ tip. For 100 µL, you can linearly scale up the payload amount.

3. Incubate the Cas9/gRNA complex in Resuspension Buffer R at room temperature for 5–20 minutes.

Prepare T-cells for electroporation

1. To remove the Dynabeads™ Human T-Activator CD3/CD28 beads from the T-cells, place the tube on a magnetic rack for 1–2 minutes, then transfer the supernatant containing the T-cells to a new tube.

2. Count the T-cells, then collect 2×10^5 cells for each 10 µL Neon™ electroporation.

Note: The optimal T-cell concentration for both 10-µL and 100-µL electroporations is 2×10^7 – 3×10^7 cells/mL.

3. Wash the T-cells once with DPBS (no calcium, no magnesium) in 1.5-mL centrifuge tubes.

4. Resuspend the T-cells in 5 µL of Buffer R, then gently mix with 6 µL of Cas9/gRNA complex.

Note: We recommend preparing extra amount of cells needed to avoid pipette errors. For example, prepare 4×10^5 of T-cells in 10 µL Buffer R, then transfer 5 µL of cells for one reaction.

5. *(Optional)* For knock-in studies, add 0.5–1 µg of double-stranded or 50–100 pmol of single-stranded DNA (Shumman K, et al., 2015) into the mix, then electroporate. Alternatively, use AAV6 to deliver donor DNA after Cas9/gRNA complex delivery by the Neon™ Transfection System.

Electroporate using the Neon™ Transfection System

1. Pipette 10 µL of the T-cells mixed with Cas9/gRNA complexes into the Neon™ 10-µL tip.

IMPORTANT! Avoid creating bubbles, which can hinder electroporation.

2. Use program #24 (1600 V/10 ms/3 pulses) for electroporation.

Note: Programs #16 and #23 also work well for T-cells.

3. Immediately transfer the electroporated cells into a 24-well plate containing 0.5 mL of pre-warmed culture medium described in "Isolate and activate T-cells".

4. Transfer the plate to a humidified 37°C, 5% CO₂ incubator, then incubate the cells for 48 hours.

5. Verify the editing efficiency using flow cytometry or the GeneArt™ Genomic Cleavage Detection Kit.

Reference

Shumman K, Lin S, Boyer E, et al. (2015) Generation of knock-in primary human T cells using Cas9 ribonucleoproteins. *Proc Natl Acad Sci USA* 112:10437–10442. doi: 10.1073/pnas.1512503112.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at thermofisher.com/support.



Manufacturer: Life Technologies Corporation | 5791 Van Allen Way | Carlsbad, CA 92008

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

© 2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

thermofisher.com/support | thermofisher.com/askaquestion

thermofisher.com

01 May 2018

ThermoFisher
S C I E N T I F I C