Takara Bio USA, Inc.

Guide-it™ Recombinant Cas9 (3 µg/µl) User Manual

Cat. Nos. 632640, 632641 (021121)

 Takara Bio USA, Inc.

 1290 Terra Bella Avenue, Mountain View, CA 94043, USA

 U.S. Technical Support: technical support@takarabio.com

Guide-it™ Recombinant Cas9 (3 µg/µl) User Manual

Tab	e of Contents		
I. I	ntroduction	3	
A.	Summary	3	
II. I	ist of Components	3	
III.	Additional Materials Required	3	
А.	Electroporation Supplies	3	
B.	Mammalian Cell Culture Supplies	4	
C.	. General Supplies		
D.	sgRNA Development and Production	4	
E.	Detection and Characterization of Gene Editing	4	
IV.	Protocol Overview	5	
V. E	Electroporation Protocol for Neon Transfection System	5	
А.	Protocol: Preparation of Cells and Media	5	
В.	Protocol: Preparation of Cas9-sgRNA RNP Complex		
C.	Protocol: Electroporation		
VI.	Electroporation Protocol for 4D-Nucleofector System		
А.	Protocol: Preparation of Cells		
В.	3. Protocol: Preparation of Cas9-sgRNA RNP Complex		
C.	Protocol: Electroporation	8	
VII.	References	9	

Table of Figures

Figure 1. Protoco	ol overview for Gu	de-it Recombinant	: Cas9 (3 µg/µl).	
1.8				

I. Introduction

A. Summary

The CRISPR/Cas9 system has emerged as a powerful tool for gene editing because of its high targeting specificity, editing efficiency, and ease of use in virtually any organism. CRISPR/Cas9 technology consists of two key components that form a complex: Cas9 endonuclease and a single guide RNA (sgRNA) that directs Cas9 to cleave genomic DNA in a sequence-specific manner (Jinek et al. 2012). This RNA-programmable method exploits the error-prone nature of the non-homologous end joining DNA repair pathway (NHEJ) to generate gene knockouts (via insertion/deletion). The method can also be used to generate knockins via the homology-directed repair (HDR) pathway.

CRISPR/Cas9 system components have been delivered successfully into target cells through a variety of approaches, including vector-based expression systems, transfection of RNA, and more recently, introduction of Cas9-sgRNA ribonucleoprotein (RNP) complexes. Delivery of Cas9-sgRNA RNPs provides a fast turnaround for gene-editing experiments while minimizing the likelihood of off-target effects compared to vector-based approaches (Sander and Joung 2014), and this approach has been optimized for various cell types using microinjection, electroporation, and lipid-mediated transfection (Liang et al. 2015).

Guide-it Recombinant Cas9 (3 μ g/ μ l) is a recombinant wild-type *Streptococcus pyogenes* Cas9 nuclease expressed with a C-terminal nuclear-localization signal (NLS) and purified from *E. coli* for use in CRISPR/Cas9-mediated gene editing experiments. The Cas9 protein solution has been verified to be sterile and well-tolerated by mammalian cells when electroporated as a ribonucleoprotein complex (RNP) with a single guide RNA (sgRNA) for knockout experiments, or as an RNP with a donor repair template for knockin experiments.

II. List of Components

Guide-it Recombinant Cas9 (3 µg/µl) (Cat. No. 632641)

• 100 µg Guide-it Recombinant Cas9 (3 µg/µl)

Guide-it Recombinant Cas9 (3 µg/µl) (Cat. No. 632640)

- $3 \times 100 \ \mu g$ Guide-it Recombinant Cas9 ($3 \ \mu g/\mu l$)
- Store Guide-it Recombinant Cas9 (3 μ g/ μ l) at -70°C.
- Avoid repeated freeze/thaw cycles. We recommend preparing aliquots upon initial thawing of Guide-it Recombinant Cas9 (3 μg/μl).

III. Additional Materials Required

The following reagents/materials are required but not included.

A. Electroporation Supplies

Use of this product requires an electroporator, electroporation chamber (typically cuvettes or tips), and an electroporation buffer that is suitable for your target cells. Here we provide separate guidelines for the Neon Transfection System (Thermo Fisher Scientific, Cat. No. MPK5000) and the 4D-Nucleofector System (Lonza, Cat. No. AAF-1002B).

B. Mammalian Cell Culture Supplies

- Culture medium, supplies, and additives specific to your target cells
- Cell culture plates
- PBS without Ca²⁺ or Mg²⁺
- Trypsin/EDTA or equivalent
- Humidified incubator (set at 37° C, 5% CO₂)

C. General Supplies

- Single-channel pipettes
- Nuclease-free thin-wall PCR tubes or strips

D. sgRNA Development and Production

CRISPR/Cas9 gene editing requires a custom sgRNA with a user-designed targeting sequence that is homologous to the target gene or genomic region of interest. Selecting an appropriate DNA sequence at the target region is critical for maximizing the potential for efficient cleavage at the target site and for minimizing the likelihood of non-specific cleavage events. There are several freely available online tools that can be helpful for determining suitable sgRNA target sequences for a given organism and genomic target. For a list of these tools, please refer to:

http://www.takarabio.com/US/Products/Genome_Editing/CRISPR_Cas9/Resources/Online_tools_for_guide_RNA_design.

NOTE: For many applications, it is advisable to design and test several variant sgRNAs against the same genomic target region.

Candidate sgRNAs must ultimately be produced in sufficient quantity for the generation of functional Cas9-sgRNA RNPs. For development and production of user-designed sgRNAs, we recommend either of the following kits:

- For constructing and purifying sgRNAs: Guide-it sgRNA *In Vitro* Transcription Kit (Takara Bio, Cat. No. 632635).
- For constructing and purifying sgRNAs, and testing target cleavage efficiencies *in vitro*: Guide-it Complete sgRNA Screening System (Takara Bio, Cat. No. 632636).

E. Detection and Characterization of Gene Editing

These items are recommended for determining the efficiency of gene editing and the nature of the edits:

<u>Cat. No</u> .	<u>Product</u>	<u>Size</u>
631443	Guide-it Mutation Detection Kit	100 rxns
631448	Guide-it Mutation Detection Kit	25 rxns
632611	Guide-it Genotype Confirmation Kit	100 rxns
631444	Guide-it Indel Identification Kit	10 rxns

IV. Protocol Overview

Please read each relevant protocol completely before starting. Successful results depend on understanding and performing the following steps correctly.



Figure 1. Protocol overview for Guide-it Recombinant Cas9 (3 µg/µl).

V. Electroporation Protocol for Neon Transfection System

Here we provide protocols for performing knockout and knockin experiments in hiPS cells and CD34-positive stem cells using the Neon Transfection System. While these protocols may serve as a helpful starting point for electroporation of other cell types as well, further optimization will be required. Please refer to the Neon Transfection System. System User Manual and manufacturer's website for detailed operating instructions for the Neon Transfection System.

A. Protocol: Preparation of Cells and Media

Cultured target cells are harvested, washed, and resuspended in the appropriate buffer.

- Prepare a sufficient number of fresh cells for your experiment.
 NOTE: Each electroporation requires 1 x 10⁵ cells. However, due to the potential variation of pipette and tip volumes, we recommend preparing 1.5X the necessary volume of cell suspension (i.e., 1.5 x 10⁵ cells) for electroporation with a 10-μl Neon Tip to ensure that there is sufficient volume.
- 2. For hiPS cells (adherent cells), continue to Step 3. For CD34-positive stem cells (suspension cells), skip to Step 5.
- 3. Aspirate the medium, wash the cell layer once with PBS (without Ca²⁺ and Mg²⁺), and dissociate the cells using TrypLE Select Enzyme (1X) (Thermo Fisher Scientific, Cat. No. 12563011).
- 4. Harvest the cells in growth medium.
- 5. Take an aliquot of the cell suspension and measure the cell density using your preferred method.
- 6. Harvest the cells by centrifugation at 400g for 5 min in a 15-ml conical tube.
- Wash the cells once with PBS (without Ca²⁺ and Mg²⁺), and then resuspend hiPS cells in Buffer R and CD34-positive stem cells in Buffer T (included with Neon kits) at a concentration of 2 x 10⁷ cells/ml (i.e., 1.5 x 10⁵ cells in 7.5 μl).

NOTE: Use Resuspension Buffer R for established adherent and suspension cells as well as primary adherent cells, and use Resuspension Buffer T for primary blood-derived suspension cells.

8. Keep the cell suspension on ice until use.

B. Protocol: Preparation of Cas9-sgRNA RNP Complex

Cas9 and sgRNA components are combined to form RNP complexes for electroporation.

- Thaw Guide-it Recombinant Cas9 (3 μg/μl) and sgRNA solutions at room temperature.
 NOTE: We recommend preparing aliquots upon initial thawing of Guide-it Recombinant Cas9 (3 μg/μl) to avoid repeated freeze/thaw cycles.
- Combine the following components in a 200-μl PCR tube to mix the Cas9 protein and sgRNA at a 5:1 mass ratio. The molar ratio of Cas9 protein to sgRNA will be approximately 1:1 in this mixture, and the total volume will be 7.5 μl. Be sure to use the same buffer that was used to resuspend the cells.
 NOTE: The reaction volume indicated below is 1.5X the required volume.

Per reaction:

0.45 μl*	sgRNA (e.g., 1 µg/µl)
0.75 µl	Guide-it Recombinant Cas9 (3 µg/µl)
6.3 μl*	Resuspension Buffer R or T
7.5 ul	Total volume

*The added volume of sgRNA will vary depending on sgRNA concentration, and the added volume of Resuspension Buffer should be adjusted such that the total reaction volume is 7.5 μ l. The volumes indicated above are based on a sgRNA concentration of 1 μ g/ μ l.

NOTE: Make a master mix if you are performing multiple electroporations.

NOTE: The optimal amount of RNP complex may vary for different cell types.

NOTE: To maximize electroporation efficiency, the combined volume of the Cas9 and sgRNA solutions should be $\leq 20\%$ of the total volume of the Cas9-sgRNA RNP complex reaction (e.g., for the 7.5-µl reaction specified above, the combined volume of the sgRNA and Cas9 solutions should be ≤ 1.5 µl). **NOTE:** If you plan to use donor DNA to induce HDR-mediated knockin, add the DNA after the subsequent incubation step (Step 3). We recommend using ≤ 1 µg of DNA for knockin experiments. Adjust the volume of Resuspension Buffer R or T included in the reaction such that the final volume upon addition of donor DNA is 7.5 µl.

3. Mix the reaction well by gently pipetting up and down. Incubate using a thermal cycler preheated to 37°C with the following program:

37°C	5 min
4°C	hold

4. <u>OPTIONAL:</u> Add donor DNA and keep on ice until use.

C. Protocol: Electroporation

Cas9-sgRNA RNPs are electroporated into target cells.

- 1. Fill the Neon Tube with 3 ml of Buffer E (included with Neon kits) and insert the Neon Tube into the Neon Pipette Station.
- 2. Using the touchscreen on the Neon system, set up the electroporation parameters as follows:

Pulse voltage / Pulse width / Pulse number = 1100 v / 20 ms / 2 pulses

NOTE: We have used these parameters successfully for hiPS cells and CD34-positive stem cells using the Neon Transfection System. Optimization of electroporation parameters will be required for different target cell types. Suggested parameters for different cell types are included in the supplementary material for (Liang et al. 2015).

- 3. Gently resuspend the cells by tapping, and transfer 7.5 μl of the cell suspension into the PCR tube containing the 7.5 μl of Cas9-sgRNA RNP complex solution.
- 4. Mix well by gently pipetting up and down.
- 5. Insert the Neon Pipette into the Neon Tip and confirm that the pipette and tip are tightly connected.
- Using the Neon Pipette, aspirate the mixture slowly into the Neon Tip.
 NOTE: Avoid any air bubbles in the tip. If you notice air bubbles, place the sample back into the PCR tube and aspirate again into the tip without any air bubbles.
- 7. Insert the Neon Pipette into the Neon Tube placed in the Neon Pipette Station and run the program.
- 8. Remove the pipette very carefully and transfer the cells into a cell culture plate with pre-warmed medium.

NOTE: Use an appropriate well plate for your target-cell type. We had success using 24 and 48-well plates for CD34-positive stem cells and hiPS cells, respectively. hiPS cells typically require greater confluence than regular adherent cells.

9. Shake the plate appropriately to disperse the cells and incubate at 37°C in a humidified incubator with 5% CO₂ until the next necessary procedure.

VI. Electroporation Protocol for 4D-Nucleofector System

Here we provide protocols for performing knockout and knockin experiments in Jurkat and CD34-positive stem cells using the 4D-Nucleofector System with 16-well Nucleocuvette Strips. While these protocols may serve as a helpful starting point for electroporation of other cell types as well, further optimization will be required. Please refer to the 4D-Nucleofector System User Manual and manufacturer's website for more detailed information.

A. Protocol: Preparation of Cells

Cultured target cells are harvested, washed, and resuspended in the appropriate solution.

- 1. Prepare a sufficient number of fresh cells for your experiment. **NOTE:** Each electroporation requires 2×10^5 cells.
- 2. Take an aliquot of the cell suspension and measure the cell density using your preferred method.
- 3. Harvest the cells by centrifugation at 400g for 5 min in a 15-ml conical tube.
- 4. Wash once with PBS (without Ca²⁺ and Mg²⁺), and then resuspend Jurkat cells in SE Nucleofector Solution (with supplement) and CD34-positive stem cells in P3 Nucleofector Solution (with supplement) at a concentration of 1 x 10⁷ cells/ml (i.e., 2 x 10⁵ cells in 20 μl).
 NOTE: Please refer to the 4D-Nucleofector System User Manual and manufacturer's website for more information about working with other cell types.
 NOTE: 20 μl of cell suspension will be needed per well of the Nucleocuvette Strip.
- 5. Keep the cell suspension on ice until use.

B. Protocol: Preparation of Cas9-sgRNA RNP Complex

Cas9 and sgRNA components are combined to form RNP complexes for electroporation.

- Thaw Guide-it Recombinant Cas9 (3 μg/μl) and sgRNA solutions at room temperature.
 NOTE: We recommend preparing aliquots upon initial thawing of Guide-it Recombinant Cas9 (3 μg/μl) to avoid repeated freeze/thaw cycles.
- 2. Combine the following components in a 200-µl PCR tube to mix the Cas9 protein and sgRNA at a 5:1 mass ratio:

Per reaction (e.g. 5 µl):

2µl*	sgRNA (e.g., 1 µg/µl)
3.3 µl	Guide-it Recombinant Cas9 (3 µg/µl)
5.3 µl	Total volume

*The added volume of sgRNA will vary depending on sgRNA concentration. The volume indicated above is based on an sgRNA concentration of 1 μ g/ μ l.

NOTE: Make a master mix if you are performing multiple electroporations.

NOTE: The RNP volume required for efficient transfection needs to be optimized for different cell types. Usually $\leq 10 \ \mu$ l of the RNP solution will be tested for electroporation in each well of the 16-well Nucleocuvette Strip.

3. Mix the reaction well by gently pipetting up and down. Incubate using a thermal cycler preheated to 37°C with the following program:

37°C	5 min
4°C	Hold

C. Protocol: Electroporation

Cas9-sgRNA RNPs are electroporated into target cells.

- 1. Label wells of Nucleocuvette Strips to be used for electroporation.
- Combine 20 μl of cell suspension with 5 μl Cas9-sgRNA RNP complex solution in each well of the Nucleocuvette Strip, and mix well by gently pipetting up and down.
 NOTE: If you plan to use donor DNA to induce HDR-mediated knockin, add the donor DNA at this step.
- 3. Select the program CL-120 for Jurkat cells or the program D0-100 for CD34-positive stem cells.
- 4. Insert the Nucleocuvette Strip into the Nucleofector machine and run the program.
- 5. Add 80 μl of pre-warmed medium to each cuvette and allow electroporated cells to recover for 12 min post-transfection at room temperature.
- 6. Gently collect the cells along with the media from each well and transfer to individual wells of a 48well plate containing pre-warmed medium.
- 7. Shake the plate appropriately to disperse the cells and incubate at 37°C in a humidified incubator with 5% CO₂ until the next necessary procedure.

VII. References

- 1. Jinek, M., Chylinsky, K., Fonfara, I., Hauer, M., Doudna, J.A., & Charpentier, E. A programmable dual-RNA-guided DNA endonuclease in adaptive bacterial immunity. *Science*. **337(6096)**, 816–21 (2012).
- 2. Liang, X. *et al.* Rapid and highly efficient mammalian cell engineering via Cas9 protein transfection. *J. Biotechnol.* **208**, 44–53 (2015).
- 3. Sander, J.D. & Joung, J.K. CRISPR-Cas9 systems for genomic editing, regulation and targeting. *Nat. Biotechnol.* **32**, 347-55 (2014).

Contact Us		
Customer Service/Ordering	Technical Support	
tel: 800.662.2566 (toll-free)	tel: 800.662.2566 (toll-free)	
fax: 800.424.1350 (toll-free)	fax: 800.424.1350 (toll-free)	
web: <u>takarabio.com</u>	web: <u>takarabio.com</u>	
e-mail: ordersUS@takarabio.com	e-mail: techUS@takarabio.com	

Notice to Purchaser

Our products are to be used for **Research Use Only**. They may not be used for any other purpose, including, but not limited to, use in humans, therapeutic or diagnostic use, or commercial use of any kind. Our products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without our prior written approval.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product's web page at <u>takarabio.com</u>. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements

© 2021 Takara Bio Inc. All Rights Reserved.

All trademarks are the property of Takara Bio Inc. or its affiliate(s) in the U.S. and/or other countries or their respective owners. Certain trademarks may not be registered in all jurisdictions. Additional product, intellectual property, and restricted use information is available at <u>takarabio.com</u>.

This document has been reviewed and approved by the Quality Department.