

Product Information

Cas9 Lambda Red and gRNA Plasmids For CRISPR-Mediated Recombineering in *E. coli*

Cas9 Lambda Red Homologous Recombination Plasmid for *E. coli*
 Catalog Number **CAS9BAC1P**

Custom CRISPR HR gRNA Plasmid
 Catalog Number **CRISPRBACD**

Storage Temperature -20 °C

CRISPR Non-Target Negative Control Plasmid for Bacteria
 Catalog Number **CRISPR31**

CRISPR *LacZ* Positive Control Plasmid for Bacteria
 Catalog Number **CRISPR30**

TECHNICAL BULLETIN

Product Description

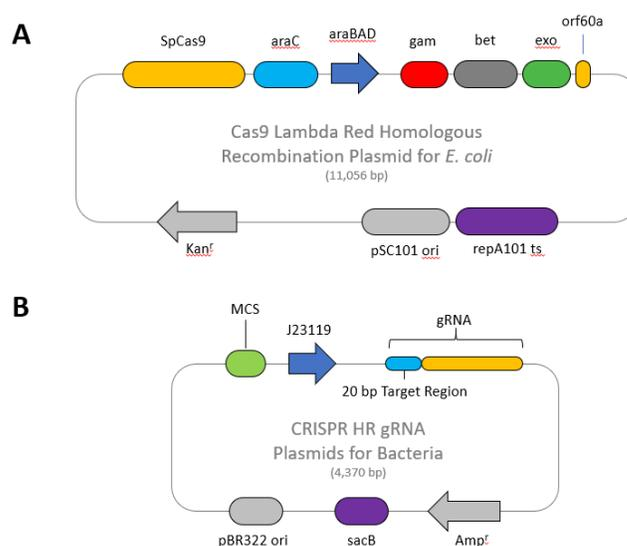
These plasmids constitute a two-part system for CRISPR-mediated recombineering in *Escherichia coli* (*E. coli*).

The Cas9 Lambda Red Homologous Recombination Plasmid for *E. coli* (Catalog Number CAS9BAC1P) contains the gene for Cas9 from *Streptococcus pyogenes* (spCas9) expressed from its native promoter, as well as the genes for λ-red recombinases *exo*, *beta*, and *gam* under the control of the arabinose-inducible ParaB promoter. This plasmid confers kanamycin resistance and possesses the repA101ts temperature-sensitive origin of replication, allowing for easy plasmid maintenance and curing (see Figure 1A).

- The Custom CRISPR HR gRNA Plasmid (Catalog Number CRISPRBACD) is a custom plasmid containing the spCas9 gRNA with a custom-designed spacer sequence specific to the gene of interest.
- The CRISPR *LacZ* Positive Control Plasmid for Bacteria (Catalog Number CRISPR30) contains a gRNA spacer targeting the *lacZ* gene in wild-type *E. coli*.
- The CRISPR Non-Target Negative Control Plasmid for Bacteria (Catalog Number CRISPR31) contains a non-targeting spacer.

All gRNAs are expressed constitutively from a J23119 promoter. These plasmids also contain an ampicillin resistance marker, a pBR322 origin of replication, and a *sacB* gene from *Bacillus subtilis* for counter-selection-based curing (see Figure 1B).

Figure 1.
 Bacterial CRISPR HR Vector Constructs



(A) Cas9 Lambda Red Homologous Recombination Plasmid for *E. coli* Vector Map. (B) CRISPR HR gRNA Plasmids for Bacteria. The vector map shown in B illustrates the structure of all described gRNA plasmids, including the Custom CRISPR HR gRNA Plasmid, CRISPR Non-Target Negative Control Plasmid for Bacteria, and CRISPR *LacZ* Positive Control Plasmid for Bacteria. These gRNA plasmids vary only in the sequence of the 20 bp target region of the gRNA.

Background Information

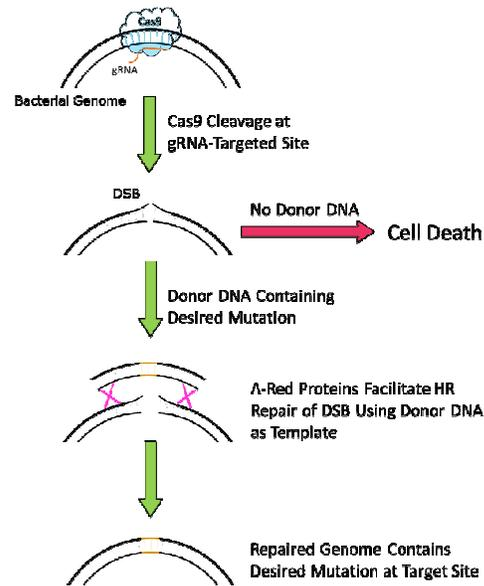
Microbial genome engineering is used for many biotechnological and biomedical purposes, including metabolic engineering, strain optimization, and general molecular biology research. One of the most commonly engineered genomes is that of *Escherichia coli* (*E. coli*).¹ Our CRISPR/Cas9-mediated λ -Red vector system offers a simple, quick method for markerless and scarless gene editing that is significantly more efficient and amenable to multiplexing than traditional methods.

E. coli has an active endogenous homologous recombination (HR) repair system, so chromosomal integration can be achieved simply by donating DNAs with long (~1,000 bp) homology arms to recombine at the target insertion site. Unfortunately, this method is very inefficient.² Recombination-mediated genetic engineering, or recombineering, is a popular method that utilizes phage-derived proteins to assist HR and increase the efficiency of chromosomal integration.³

The CRISPR system is a bacterial immune system that has been exploited by researchers as a revolutionary gene editing tool. In this system, a CRISPR-associated (Cas) endonuclease is guided by a gRNA to a specific DNA sequence, where it creates a double-stranded break (DSB).⁶

Until now, this nuclease-based method has been mostly overlooked as a microbial gene editing tool because most bacteria lack DSB repair mechanisms, rendering the technology toxic.⁷ However, when CRISPR/Cas9 is used to mediate recombineering, this cytotoxic quality offers an advantage: Cas9-induced DSBs kill those cells that fail to recombine with the donor DNA, avoiding the need for selectable markers to select edited cells⁵ (see Figure 2).

Figure 2. CRISPR-Mediated Recombineering Eliminates the Need for Selection.



Recent publications using CRISPR/Cas9-mediated recombineering in *E. coli* tout editing efficiencies near 100%,⁵ making CRISPR/Cas9-mediated recombineering the most powerful bacterial genome engineering method to date. In addition, Cas9-mediated recombineering overcomes the dependence on a second recombination step, avoids the creation of destabilizing scar sites, can be used in multiplexing, and is less time-consuming than previous procedures.⁸

Here we present a novel dual-vector CRISPR/Cas-mediated λ -Red system for improved recombineering in *E. coli*. This system is shown to facilitate homology-directed repair of DSBs created by the Cas9 endonuclease, enabling genetic alterations through chromosomal integration of a donor DNA. These vectors have been tested in multiple strains of *E. coli*, and have been shown to yield insertions, deletions, and point mutations, as well as multi-gene manipulations, all with high efficiency. This donor can be either ssDNA or dsDNA with homology arms of 45–50 or 150–500 nucleotides, respectively. For more information about donor design, please see the corresponding donor design sections in the procedure.

Procedure Overview

In general, the steps required for successful engineering of *E. coli* using this system are as follows:

1. Design and order custom gRNA plasmid.
2. Design custom donor DNA template(s); order or PCR amplify design(s).
3. Prepare *E. coli* strain of interest via electroporation of the Cas9 Lambda Red Homologous Recombination Plasmid for *E. coli*.
4. Electroporate *E. coli* strain from step 3 with gRNA plasmid and donor DNA.
5. Harvest genomic DNA or perform colony PCR and assay mutations.

Note: The provided procedure uses electroporation for transformation of *E. coli*. Other methods of transformation may be used, if preferred.

Components

Product Format

One vial containing 1 µg of plasmid DNA in Tris-EDTA (TE) buffer at a concentration of 20 ng/µl.

Note: The Custom CRISPR HR gRNA Plasmid (Catalog Number CRISPRBACD) is configurable and may be ordered in quantities larger than 1 µg.

Reagents and Equipment Recommended but Not Provided.

- Custom oligonucleotides from Sigma-Aldrich for PCR primers and ssDNA donor templates
- Electroporation System for Bacterial Cells
- SOC medium (Catalog Number S1797)
- LB Agar Plates with Kanamycin
- LB Agar Plates with 100 µg/ml Carbenicillin, 60 µg/ml X-Gal, and 0.3 mM IPTG
- L-(+)-Arabinose (Catalog Number A3256)
- LB Agar Plates with 5% Sucrose
- GenElute™ Bacterial Genomic DNA Miniprep Kit (Catalog Number G1N70)

- JumpStart™ REDTaq® ReadyMix™ (Catalog Number P0982) for colony PCR
- JumpStart™ REDAccuTaq® LA DNA Polymerase (Catalog Number D1313) for amplification of dsDNA donor templates
- Water, PCR Reagent (Catalog Number W1754)
- Gel Loading Buffer (Catalog Number G2526)
- Tris-Borate EDTA Buffer, 5× concentrate, powder blend (Catalog Number T3913)
- Ethidium Bromide Solution, 10 mg/ml in water (Catalog Number E1510)

Precautions and Disclaimer

For Research Use Only. Not for use in diagnostic procedures. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

Storage/Stability

These plasmids are shipped as frozen stocks on dry ice. Once received, it is recommended the plasmids are stored at -20 °C. Avoid repeated freezing and thawing.

Procedures

The following procedures provide general guidelines for designing and ordering custom gRNA plasmids and donor template DNAs for the target of interest, as well as a detailed account of the *LacZ* control experiment for Cas9-mediated recombineering using this system.

Note: For the *LacZ* positive control experiment, one will need to order the recommended ssDNA donor template (see Figure 3).

Custom gRNA Plasmid Design

To design and order custom gRNA plasmid(s) for the target(s) of interest, please go to the CRISPR Design Tools page at milliporesigmabioinfo.com.

1. If you need to design a gRNA based on your target gene, use the CRISPR design tool and follow the prompts to design your gRNA sequence(s). If you already know your gRNA sequence, you can input this sequence directly via the same tool.
2. When prompted, select the CRISPRBACD J23119-gRNA (bacterial plasmid). Once you have configured your custom gRNA vectors(s), submit them to be redirected to the ordering page.

- For the positive and negative control experiments, one will also need to order the corresponding positive control gRNA plasmid (Catalog Number CRISPR30) and negative control gRNA plasmid (Catalog Number CRISPR31).

Single-Stranded DNA (ssDNA) Donor Template Design

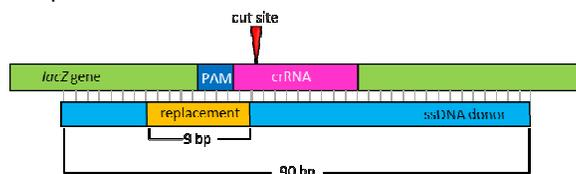
We recommend using single-stranded DNA donors when the desired mutation is a small deletion (<1 kb) or replacement (<20 bp).⁵ For insertions, as well as larger deletions and replacements, we recommend using dsDNA donors, which are covered in the next section of this procedure.

- Once you have designed your gRNA and have determined the cut site, design your ssDNA donor template using the following guidelines:
 - The donor should consist of ~90 nucleotides surrounding the cut site, altering bases as necessary to incorporate your desired mutation.
 - The resulting mutation must remove or mutate the PAM sequence or the protospacer sequence proximal to the PAM, as this prevents toxicity from repeated Cas9 cutting.⁷

Note: We recommend trying donors that are homologous to both the sense and antisense strands, as one is typically more efficient than the other, depending on the location of the target in the chromosome.

- For the positive control experiment, you will need to order the control ssDNA donor template. This 90-nt oligonucleotide contains a 9 bp replacement that introduces two premature stop codons in *lacZ* in wild-type *E. coli*, effectively knocking out *lacZ* function upon recombination at the target site.⁵ The donor design scheme and sequence are shown in Figure 3, top right.

Figure 3.
LacZ 9 bp Replacement ssDNA Donor Design and Sequence



LacZ Positive Control ssDNA Donor Sequence

```
C*C*CATTACGGTCAATCCGCCGTTTGTTCACGGA
GAATCCGTGAACCTGATACTCGCTCACATTTAATGTT
GATGAAAGCTGGCTA*C*A
```

*denotes phosphorothioate linkage

- Order your synthetic ssDNA donor design(s) through Sigma-Aldrich at sigmaaldrich.com/oligos. We recommend adding phosphorothioate linkages to the 5' and 3' termini to protect the ssDNA from nuclease degradation upon transformation.

Double-Stranded DNA (dsDNA) Donor Template Design

For larger deletions (>1 kb), replacements (>20 bp), and insertions, we recommend using a double-stranded DNA donor. Cas9-mediated λ -red recombineering has been shown to successfully delete regions up to 12 kb using 500-bp homology arms.⁵

- Once you have designed your gRNA and have determined the cut site, design your dsDNA donor template using the following guidelines:
 - The donor should consist of the desired mutation surrounded by homology arms on each side. These arms are homologous to the target site, which enables recombination between the genomic target and the dsDNA donor.
 - For most experiments, we recommend 300–500-bp homology arms; though, arms as short as 150 bp may be sufficient in many cases. We recommend trying donors of varying sizes for each of your target mutations and strains of interest.
 - The resulting mutation must remove or mutate the PAM sequence or the protospacer sequence proximal to the PAM, as this prevents toxicity from repeated Cas9 cutting.⁷
- You may order these dsDNA donors as synthetic gene fragments. Alternatively, you can construct your dsDNA donors using fusion PCR. This involves separately PCR-amplifying each homology arm (and the sequence to be inserted, if applicable) from your target strain's genomic DNA using primers that create overlaps between the two fragments, then fusing these arms together as one product in a subsequent round of PCR.⁹ For this step, we recommend using KOD Hot Start DNA Polymerase (Catalog Number 71086).

Preparation of *E. coli* Strain

We recommend transforming your *E. coli* strain of interest in two steps. In this first step, you will transform your strain with the Cas9 plasmid. We recommend saving this strain as a glycerol stock for future experiments. In a later step, you will transform this strain with your gRNA plasmids and corresponding donor DNAs.

1. Transform the Cas9 Lambda Red Homologous Recombination Plasmid for *E. coli* via electroporation, using a procedure that is optimal for your strain(s) of interest.
2. To select for positive colonies, spread 100 μ l on LB agar plate(s) with 50 μ g/ml kanamycin. Incubate for 8–12 hours at 30 °C.
3. Once colonies appear, pick one and inoculate 3 ml of liquid LB with 50 μ g/ml kanamycin in a 15-ml culture tube. Incubate for 8–12 hours at 30 °C with shaking (250 rpm).
4. Save this culture as a 15% glycerol stock and store at –80 °C for future experiments.

Note: For the *lacZ* control experiment, you will need to use a strain of *E. coli* with a wild-type *lacZ* gene to enable gene knockout and blue-white screening. This typically excludes cloning strains. We recommend wild-type *E. coli* MG1655 or *E. coli* W3110.

Gene Editing of *E. coli* (*LacZ* Positive Control Experiment)

In this second step, you will induce λ -red activity and transform your strain with your custom gRNA plasmid(s) and corresponding donor DNA(s). These steps detail this procedure for the *lacZ* control experiment.

1. The evening (or 8–12 hours) before the experiment, inoculate 5 ml of liquid LB with 50 μ g/ml kanamycin in a 15-ml culture tube using the glycerol stock from the previous step.
2. The next morning, dilute the overnight culture 1:100 in 15-ml culture tubes with 3 ml of fresh liquid LB with 50 μ g/ml kanamycin. Be sure to inoculate enough cultures to dedicate 3 ml to each experimental condition. Grow at 30 °C with shaking (250 rpm) until the OD₆₀₀ reaches 0.3–0.5 (~3 hours).

3. Once the OD₆₀₀ reaches 0.3–0.5, induce λ -red expression by adding a sterile arabinose solution to give a final arabinose concentration between 3 mg/ml (0.3% w/v) and 20 mg/ml (2% w/v). Incubate for one more hour at 30 °C with shaking.
4. For the *lacZ* control experiment, prepare 3 vials of electrocompetent cells, concentrating each final cell pellet into 40 μ l of sterile 10% glycerol. Label these 3 tubes as L, K, and E for Lawn (negative control), Killing, and Editting, respectively.

Note: Use your own optimized electroporation procedure for your strain of interest. This procedure is recommended for wild-type *E. coli* MG1655 or W3110 strains using 1 mm cuvettes. The final total volume should be 50 μ l.
5. Electroporate the following components into each corresponding vial of cells (see Table 1), using an optimized procedure for your strain of interest:

Table 1.
LacZ Control Experimental Conditions

Rxn	gRNA Plasmid	Donor
L	60 - 100 ng of CRISPR Non-Target Negative Control Plasmid for Bacteria	None
K	60 - 100 ng of CRISPR <i>LacZ</i> Positive Control Plasmid for Bacteria	None
E	60 - 100 ng of CRISPR <i>LacZ</i> Positive Control Plasmid for Bacteria	5 μ l of 10 μ M <i>lacZ</i> Positive Control ssDNA Donor (1 μ M final concentration)

6. Add 950 μ l of pre-warmed SOC medium immediately after electroporation and incubate the outgrowth at 30 °C with shaking (250 rpm) for at least 1 hour.
7. Spread outgrowth cultures onto individual LB agar plates with 100 μ g/ml carbenicillin, X-Gal, and IPTG. Initially, spread 100 μ l of outgrowth directly. For L, spreading 100 μ l of 10⁻³ dilution may be required.
8. Incubate at 30 °C for 16–24 hours to observe colony formation. Expected results are provided in Table 2.

Table 2.
LacZ Control Experimental Conditions

Rxn	Expected Result	Appearance
L	Non-targeting gRNA does not target the genome; therefore, Cas9 does not cut. <i>lacZ</i> gene is functional.	Lawn of blue colonies.
K	Guided by the <i>lacZ</i> -targeting positive control gRNA, Cas9 cuts at <i>lacZ</i> . Lack of donor prevents repair by HR, leading to cell death.	Killing; significantly fewer colonies than plate L. Colonies can be blue or white.
E	Guided by the <i>lacZ</i> -targeting positive control gRNA, Cas9 cuts at <i>lacZ</i> . HR with positive control ssDNA donor knocks out <i>lacZ</i> function via 9-bp replacement.	Editing; more colonies than plate K. Colonies should be mostly white.

Mutation Detection

There are many methods to detect mutations produced by CRISPR systems. Mutations resulting in phenotypic changes are the most easily screened (e.g., mutations in *lacZ* can be detected via blue-white screening). For genotyping, we recommend designing primers around the target site in the organism's genomic DNA for PCR amplification of the mutated region. It is important to make sure these primers fall outside the donor DNA sequence to prevent false positives. We recommend performing this PCR on isolated genomic DNA, though colony PCR may be sufficient in many cases. If the mutation is an insertion or deletion of significant size, gel electrophoresis of the PCR amplicon may be used to observe this change. For easy colony PCR, we recommend using JumpStart REDTaq ReadyMix (Catalog Number P0982), which enables samples to be loaded directly into an agarose gel without supplementary loading dye. If the mutation does not result in an appropriate size difference in the amplicon, we recommend sequencing the PCR product.

Note: For the positive control experiment, the following primers may be used to verify the 9-bp replacement at *lacZ* following blue-white screening. We recommend performing colony PCR and subsequent Sanger sequencing. The same primers may be used for both steps. These primer sequences are provided in Table 3.

Table 3.
LacZ Control Experiment Sequencing Primers

Primer	Sequence
<i>lacZ</i> -Forward	GTGAAAAGAAAAACCCACCTGGC
<i>lacZ</i> -Reverse	TGTTACCCGTAGGTAGTCACGCAACT

Plasmid Curing

Upon obtaining a correct clone, the gRNA plasmids can be cured by SacB-sucrose based counter-selection and the Cas9 Lambda Red Homologous Recombination Plasmid for *E. coli* by incubating the clone at 37 °C. To remove both simultaneously, follow these steps:

1. Spread a 200 µl dilution (10^{-1} , 10^{-2} , or 10^{-3}) of the selected clone on an LB agar plate with 5% sucrose (Teknova, Catalog Number L5110) and incubate at 37 °C overnight (or for 8–12 hours).
2. The next day, pick 8–16 large colonies and streak each onto 3 different plates: LB agar, LB agar with 100 µg/ml ampicillin or carbenicillin, and LB agar with 50 µg/ml kanamycin. Incubate these plates at 37 °C overnight.
3. Observe the plates. Those clones that are sensitive to both ampicillin (or carbenicillin) and kanamycin have been cured of both editing plasmids. Save the successfully cured plasmids from the LB agar plate as glycerol stocks, if desired.

References

1. Jiang, Y., et al., Multigene Editing in the Escherichia coli Genome via the CRISPR-Cas9 System. *Appl Environ Microbiol.*, 81(7), 2506-2515 (2015). doi:10.1128/AEM.04023-14.
2. Jeong, J., et al., Genome-Scale Genetic Engineering in *Escherichia Coli*. *Biotechnol Adv.*, 31(6), 804-810 (2013).
3. Sharan, S.K., et al., Recombineering: a homologous recombination-based method of genetic engineering. *Nat Protoc.*, 4(2), 206-223 (2009). doi:10.1038/nprot.2008.227 .
4. Datsenko, K.A. and Wanner, B.L. One-step inactivation of chromosomal genes in *Escherichia coli* K-12 using PCR products. *PNAS*, 97(12), 6640-6645 (2000).
5. Li, Y., et al., Metabolic engineering of *Escherichia coli* using CRISPR-Cas9 mediated genome editing. *Metab Eng.*, 31, 13-21 (2015). doi:10.1016/j.ymben.2015.06.006.

6. Jinek, M., et al., A Programmable Dual-RNA-Guided DNA Endonuclease in Adaptive Bacterial Immunity. *Science* (80-), 337, 816-822 (2012).
7. Jiang, W., et al., RNA-guided editing of bacterial genomes using CRISPR-Cas systems. *Nat Biotechnol.*, 31(3), 233-239 (2013). doi:10.1038/nbt.2508.
8. Pyne, M.E., et al, Coupling the CRISPR/Cas9 system to lambda red recombineering enables simplified chromosomal gene replacement in *Escherichia coli*. *Appl Environ Microbiol.* (2015). doi:10.1128/AEM.01248-15.
9. Shevchuk, N.A., et al., Construction of long DNA molecules using long PCR-based fusion of several fragments simultaneously. *Nucleic Acids Res.*, 32(2), e19-e19 (2004). doi:10.1093/nar/gnh014.
10. Chayot, R., et al., An end-joining repair mechanism in *Escherichia coli*. *Proc Natl Acad Sci.*, 107(5), 141 LP-2146 (2010). doi:10.1073/pnas.0906355107.

GenElute, JumpStart, and ReadyMix are trademarks of Sigma-Aldrich Co. LLC.
 REDTaq is a registered trademark of Sigma-Aldrich Co. LLC.
 Pulser is a registered trademark of Bio-Rad.
 XCell is a trademark of Bio-Rad.

CHM,MAM, EE 06/20-1

Troubleshooting

If no cutting is observed and there is reason to suspect an experimental flaw is at fault, the following considerations may aid the researcher in troubleshooting the experiment.

Suspected Issue	Solution
One of the plasmids has been thawed and refrozen too many times.	Our plasmid DNA has been shown to withstand some freezing and thawing, but aliquoting each into smaller quantities will allow this potential issue to be avoided.
The electroporation or other transformation method is not working or is too toxic.	For any transformation method, the procedure should be optimized for each strain used. For troubled electroporations, refer to the manufacturer's procedure for further assistance. For the positive control experiment in <i>E. coli</i> MG1655 or <i>E. coli</i> W3110, we used the Gene Pulser® XCell™ device (BioRad, Catalog Number 165266) with Gene Pulser Electroporation Cuvettes, 0.1 cm gap (BioRad, Catalog Number 1652083). To electroporate, we used the pre-set program for <i>E. coli</i> in 0.1 cm cuvettes found on this device.
Blue and white colonies are seen on the "killing" control plate in the <i>lacZ</i> control experiment.	While most Cas9-generated DSBs will lead to cell death without a donor template for repair, some escaper colonies are to be expected, and are not necessarily an indicator of a larger problem. Blue colonies are often escapers of the CRISPR system, most commonly due to spontaneous mutations in the colony's plasmid copy of Cas9. White colonies often indicate mutations in the target sequence (<i>lacZ</i>), possibly due to very low levels of naturally occurring NHEJ repair.
The sequencing primers generate no PCR product after gene editing.	If no PCR product is observed, it is possible that the corresponding genomic binding sites are no longer present. This may be due to low levels of repair via the <i>E. coli</i> alternative end-joining pathway (A-EJ), which creates large genomic rearrangements/deletions (rather than precise changes like those generated via HR). ¹⁰ We recommend double-checking your primer sequences and optimizing your PCR reagents and methods; however, if a deletion is still suspected, you may choose to design alternate primers that bind farther from the target site.

US Patents Pending

License Agreements

CRISPR Use License Agreement

This Product and its use are the subject of one or more of the following patents controlled by The Broad Institute, Inc. (BROAD), the Massachusetts Institute of Technology (MIT), or the President and Fellows of Harvard College (HARVARD): U.S. Patent Nos. 8,697,357; 8,771,945, and any substitutions, divisions, continuations, reissues, renewals, re-examinations or extensions, and corresponding foreign patent applications and patents.

BEFORE OPENING OR USING THIS PRODUCT, PLEASE READ THE TERMS AND CONDITIONS SET FORTH IN THIS LICENSE AGREEMENT. YOUR USE OF THIS PRODUCT SHALL CONSTITUTE ACKNOWLEDGMENT AND ACCEPTANCE OF THESE TERMS AND CONDITIONS. If you do not agree to use this Product pursuant to the terms and conditions set out in this License Agreement, please contact Sigma Technical Services within ten days of receipt to return the unused and unopened Product for a full refund; provided, however, that custom-made Products may not be returned for a refund.

The purchase of this Product conveys to you, the buyer, the non-transferable right to use the purchased Product for Licensed Research Use (see definition below) subject to the conditions set out in this License Agreement. If you wish to use this Product for any purpose other than Licensed Research Use, you must first obtain an appropriate license (see information set out below).

This Product may not be used for any purpose other than Licensed Research Use. Your right to use this Product for Licensed Research Use is subject to the following conditions and restrictions:

"Licensed Research Use" means any use for research purposes, with the following exceptions: (i) you may not sell or otherwise transfer Licensed Products (including without limitation any material that contains a Licensed Product in whole or part) or any Related Material to any other person or entity, or use Licensed Products or any Related Material to perform services for the benefit of any other person or entity, (ii) You may use only the purchased amount of the Licensed Products and components of the Licensed Products, and shall use any Related Material, only for your internal research as a research tool for research purposes; provided, however, that notwithstanding the foregoing, you shall not use the Licensed Products for: (a) any clinical use, including, without limitation, diagnostic and prognostic use, (b) any human, veterinary, livestock or commercial agricultural use, (iii) modification or reverse-engineering of the Product in any way or creating any derivatives or sequence variants thereof, or (iv) the manufacture, distribution, importation, exportation, transportation, sale, offer for sale, marketing, promotion or other exploitation or use of, or as, a testing service, therapeutic or diagnostic for humans or animals, and not for any Commercial Purposes, (v) you shall use Licensed Products and any Related Material in compliance with all applicable laws and regulations, including without limitation applicable human health and animal welfare laws and regulations.

"Commercial Purposes" means (a) the practice, performance or provision of any method, process or service, or (b) the manufacture, sale, use, distribution, disposition or importing of any product, in each case (a) or (b) for consideration, e.g., a fee, or on any other commercial basis.

Your right to use the Product will terminate immediately if you fail to comply with these terms and conditions. You shall, upon such termination of your rights, destroy all Product, Modified Animals, and components thereof in your control, and notify Sigma of such in writing.

Broad, MIT and Harvard make no representations or warranties of any kind concerning the patent rights and hereby disclaim all representations and warranties, express or implied, including without limitation warranties of merchantability, fitness for a particular purpose, noninfringement of intellectual property rights of Broad, MIT, Harvard or third parties, validity, enforceability and scope of patent rights, validity of any claims, whether issued or pending, and the absence of latent or other defects, whether or not discoverable. In no event shall Broad, MIT, Harvard or their directors, trustees, officers, employees, agents, faculty, affiliated investigators or students, be liable for incidental or consequential damages of any kind, including without limitation economic damages or injury to property and lost profits, regardless of whether Broad, MIT or Harvard shall be advised, shall have other reason to know, or in fact shall know of the possibility of the foregoing.

For information on purchasing a license to this Product for purposes other than Licensed Research Use, contact your local Sigma Sales representative, who will refer you to the proper licensing representative, or in the USA call 800-325-3010.