

## Q5<sup>®</sup> Site-Directed Mutagenesis Kit (Without Competent Cells)

NEB #E0552S

10 reactions

Version 2.0\_7/20

### Table of Contents

Quick Protocol .....	2
Introduction .....	3
General Guidelines .....	4
Protocol.....	6
Protocol for Control Reaction.....	7
Expected Results.....	7
Troubleshooting.....	8
Quality Controls .....	9
Ordering Information.....	9
Revision History .....	10

### The Q5 Site-Directed Mutagenesis Kit (Without Competent Cells) Includes:

*The Q5 Site-Directed Mutagenesis Kit (Without Competent Cells) is stable at –20°C for two years. For flexibility, the mutagenesis reagents and control reactions are supplied without competent cells so that any chemically-competent E. coli cells suitable for cloning may be used.*

Q5 Hot Start High-Fidelity 2X Master Mix (2X)

KLD Enzyme Mix (10X)

KLD Reaction Buffer (2X)

Control SDM Primer Mix (10 μM each)

Control SDM Plasmid (5 μg/ml)

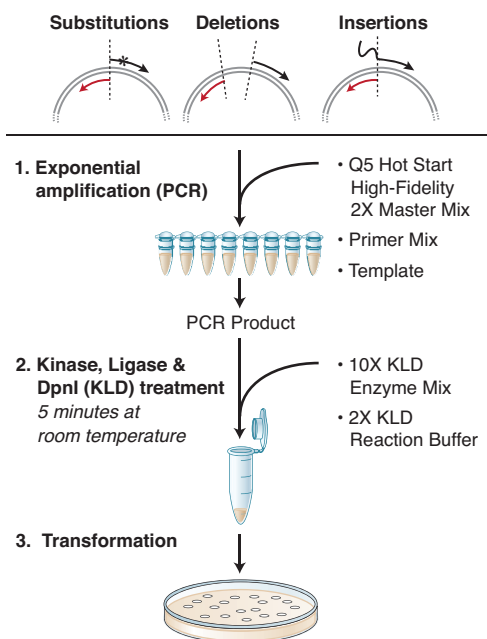
### Required Materials Not Included:

Chemically-competent *E. coli* cells

SOC Outgrowth Media

Selection Plates

## Q5 Site-Directed Mutagenesis Kit (Without Competent Cells) Quick Protocol



Primers should be designed with 5' ends annealing back-to-back. We recommend using the NEB online design software, NEBaseChanger<sup>®</sup>, which can be found at: [NEBaseChanger.neb.com](http://NEBaseChanger.neb.com).

### Step 1: Exponential Amplification

	25 $\mu$ l REACTION	FINAL CONCENTRATION
Q5 Hot Start High-Fidelity 2X Master Mix	12.5 $\mu$ l	1X
10 $\mu$ M Forward Primer	1.25 $\mu$ l	0.5 $\mu$ M
10 $\mu$ M Reverse Primer	1.25 $\mu$ l	0.5 $\mu$ M
Template DNA (1–25 ng/ $\mu$ l)	1 $\mu$ l	1–25 ng
Nuclease-free Water	9.0 $\mu$ l	

Cycling Conditions:

STEP	TEMP	TIME
Initial Denaturation	98°C	30 seconds
25 Cycles	98°C	10 seconds
	50–72°C*	10–30 seconds
	72°C	20–30 seconds/kb
Final Extension	72°C	2 minutes
Hold	4–10°C	

\*For mutagenic primers, please use the Ta provided by the online NEB primer design software (NEBaseChanger.neb.com).

### Step 2: KLD Reaction

	VOLUME	FINAL CONCENTRATION
PCR Product	1 $\mu$ l	
2X KLD Reaction Buffer	5 $\mu$ l	1X
10X KLD Enzyme Mix	1 $\mu$ l	1X
Nuclease-free Water	3 $\mu$ l	

Incubate for 5 minutes at room temperature.

### Step 3: Transformation

1. Add 5  $\mu$ l of KLD mix to 50  $\mu$ l of chemically-competent *E. coli* cells.\*
2. Incubate on ice for 30 minutes.
3. Heat shock at 42°C for 30 seconds.
4. Incubate on ice for 5 minutes.
5. Add 950  $\mu$ l SOC, gently shake at 37°C for 1 hour.
6. Spread 40–100  $\mu$ l onto appropriate selection plate, incubate overnight at 37°C.

\*see page 7 for recommendations on competent cells.

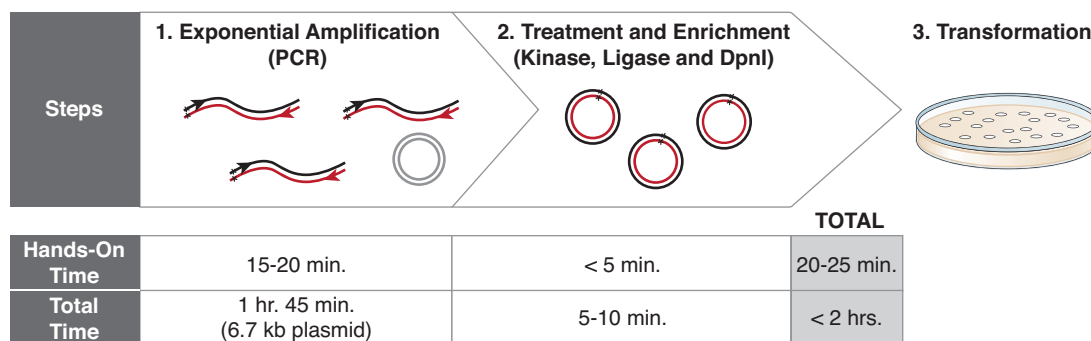
### Introduction

The Q5 Site-Directed Mutagenesis Kit (Without Competent Cells) enables rapid, site-specific mutagenesis of double-stranded plasmid DNA in less than 2 hours (Figure 1). The kit utilizes the robust Q5 Hot Start High-Fidelity DNA Polymerase along with custom mutagenic primers to create insertions, deletions and substitutions in a wide variety of plasmids. After PCR, the amplified material is added directly to a unique Kinase-Ligase-DpnI (KLD) enzyme mix for rapid (5 minutes), room temperature circularization and template removal (Figure 2). Transformation into high-efficiency chemically-competent *E. coli*, not supplied, ensures robust results with plasmids up to at least 14 kb in length.

### Advantages

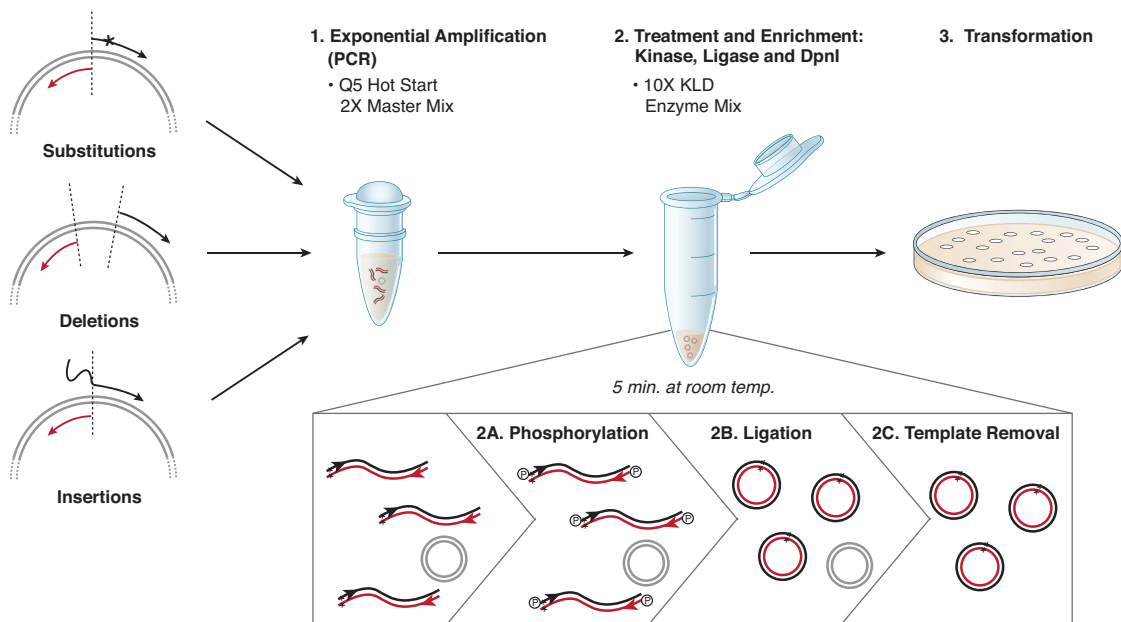
- Non-overlapping primer design ensures robust, exponential amplification, generating a high percentage of desired mutations from a wide range of templates.
- Intramolecular ligation and transformation into NEB high-efficiency competent cells results in high colony yield.
- Extremely low error rate of Q5 Hot Start High-Fidelity DNA Polymerase reduces screening time .
- Hot start polymerase enables room temperature reaction set-up
- DpnI background reduction permits a wide range of starting template concentrations
- Use of standard primers eliminates additional expenses from phosphorylated or purified oligos
- Easy-to-use PCR master mix and unique multi-enzyme KLD mix offer convenience and quality
- Rapid and direct treatment step proceeds at room temperature in 5 minutes
- Allows the use of any chemically-competent *E. coli* cells suitable for cloning

**Figure 1: Site-specific mutagenesis proceeds in less than 2 hours.**



The use of a master mix, a unique multi-enzyme KLD enzyme mix, and a fast polymerase ensures that, for most plasmids, the mutagenesis reaction is complete in less than two hours.

**Figure 2: Q5 Site-Directed Mutagenesis Overview.**



This kit is designed for rapid and efficient incorporation of insertions, deletions and substitutions into double-stranded plasmid DNA. The first step is an exponential amplification using standard primers and a master mix formulation of Q5 Hot Start High-Fidelity DNA Polymerase. The second step involves incubation with a unique enzyme mix containing a kinase, a ligase and DpnI. Together, these enzymes allow for rapid circularization of the PCR product and removal of the template DNA. The last step is a high-efficiency transformation into chemically-competent cells (not provided).

## General Guidelines

### Primer Quality/Modification

Standard desalting of the oligonucleotides is recommended. Phosphorylation and purification are not required.

### Primer Design

Although general recommendations are presented here, for best results, primers should be designed using NEBaseChanger, the NEB online primer design software, which can be found at: [NEBaseChanger.neb.com](http://NEBaseChanger.neb.com).

#### Substitutions:

Substitutions are created by designing a mismatch in the center of the mutagenic primer (Figure 3A). Include at least 10 nucleotides that are complementary to your plasmid at the 3' end of the primer. To accommodate many mutations (from 7 to 50 per primer), changes should be incorporated at the 5' end of the mutagenic primer.

The 5' end of the second primer will begin at the base next to the 5' end of the first primer, and proceed in the opposite direction on the complementary strand. This second primer can be 100% complementary to the plasmid sequence or can contain mismatches, if desired. The absence of any overlap ensures that exponential (rather than linear) amplification will take place. The maximum size of the substitution is largely dictated by oligo synthesis limitations.

#### Deletions:

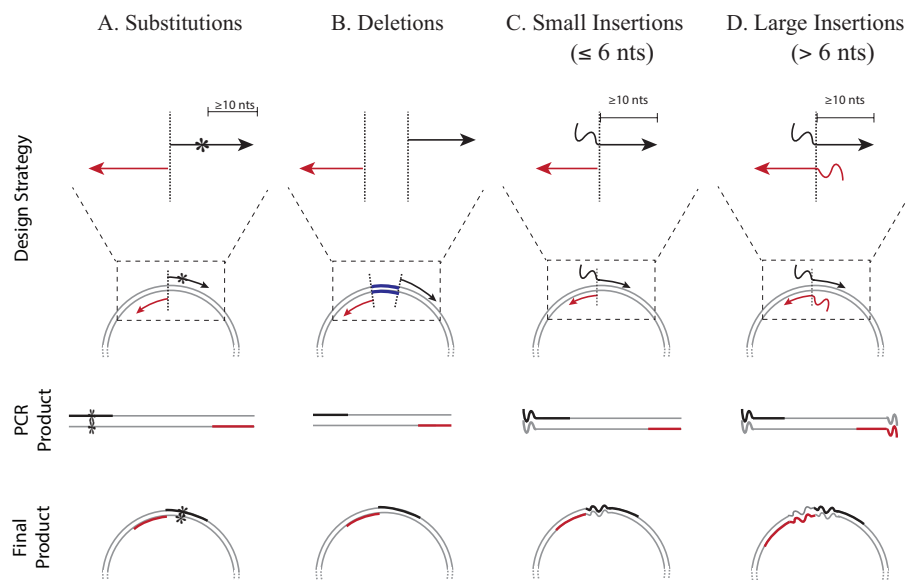
Deletions are created by designing primers that flank the region to be deleted (Figure 3B). The two primers should be designed in opposite directions with their 5' ends adjacent to the area to be deleted. The primers can be 100% complementary to the plasmid sequence or can contain mismatches and/or insertions if desired.

#### Insertions:

The sequence to be inserted should be added to the 5' end of the mutagenic primer (Figure 3C). For insertions of more than 6 nucleotides, the insertion sequence can be split between the two primers. Half of the new sequence should be added to the 5' end of the forward primer and the other half should be added to the 5' end of the reverse primer (Figure 3D). As described for substitutions, there should be at least 10 nucleotides that are complementary to your plasmid on the 3' end of each primer.

The maximum size of the insertion is largely dictated by oligo synthesis limitations. Insertions of up to 100 nucleotides (50 nucleotides at the 5' end of each primer) can routinely be accommodated using this kit.

**Figure 3: Primer Design for Q5 Site-Directed Mutagenesis Kit.**



Substitutions, deletions and insertions are incorporated into plasmid DNA through the use of specifically designed forward (black) and reverse (red) primers. **Unlike kits that rely on linear amplification, primers designed for the Q5 Site-Directed Mutagenesis Kit should not overlap to ensure that the benefits of exponential amplification are realized.** A) Substitutions are created by incorporating the desired nucleotide change(s) (denoted by \*) in the center of the forward primer, including at least 10 complementary nucleotides on the 3' side of the mutation(s). The reverse primer is designed so that the 5' ends of the two primers anneal back-to-back. B) Deletions are engineered by designing standard, non-mutagenic forward and reverse primers that flank the region to be deleted. C) Insertions less than or equal to 6 nucleotides are incorporated into the 5' end of the forward primer while the reverse primer anneals back-to-back with the 5' end of the complementary region of the forward primer. D) Larger insertions can be created by incorporating half of the desired insertion into the 5' ends of both primers. The maximum size of the insertion is largely dictated by oligonucleotide synthesis limitations.

### PCR Master Mix

Q5 Hot Start High-Fidelity 2X Master Mix contains the polymerase buffer, dNTPs, Mg<sup>++</sup> and Q5 Hot Start High-Fidelity DNA Polymerase.

For added convenience, Q5 Hot Start High-Fidelity DNA Polymerase is inactive at room temperature, allowing room temperature reaction set up. Due to the unique aptamer-based hot start format, there is no requirement for a separate activation step in the cycling protocol.

*Note: After thawing Q5 Hot Start High-Fidelity 2X Master Mix completely, a precipitate (most noticeable after the first 1–2 freeze/thaw cycles) is not uncommon. To ensure optimal performance, the master mix should be thawed completely and resuspended prior to use. Stability testing using up to 15 freeze/thaw cycles has shown no negative effect on master mix performance.*

### Template

We recommend starting with 1–25 ng of plasmid DNA; however, due to the robust nature of Q5 DNA Polymerase, experiments have also been successful using as little as 10 pg and as much as 100 ng of starting plasmid.

### Volume

The kit is designed for 25 µl PCR final reaction volumes, however, 50 µl reaction volumes work equally well. For either reaction volume, 1 µl of PCR product should be added to Step II (see page 6).

### Enzyme Mix

The 10X Kinase-Ligase-DpnI Mix (10X KLD) contains a blend of kinase, ligase and DpnI enzymes in a buffer that preserves the activity of the enzymes. This formulation allows efficient phosphorylation, intramolecular ligation/circularization and template removal in a single 5 minute reaction step at room temperature.

### Competent Cells

Q5 Site-Directed Mutagenesis protocols are optimized for use with NEB 5-alpha Competent *E. coli* (High Efficiency) Cells (NEB #C2987). Other chemically-competent *E. coli* strains suitable for cloning may be substituted, but results will vary depending upon the quality and efficiency of the cells.

## Protocol

### Step 1: Exponential Amplification (PCR)

1. Assemble the following reagents in a thin-walled PCR tube.

	25 $\mu$ l REACTION	FINAL CONCENTRATION
Q5 Hot Start High-Fidelity 2X Master Mix	12.5 $\mu$ l	1X
10 $\mu$ M Forward Primer	1.25 $\mu$ l	0.5 $\mu$ M
10 $\mu$ M Reverse Primer	1.25 $\mu$ l	0.5 $\mu$ M
Template DNA (1–25 ng/ $\mu$ l)	1 $\mu$ l	1–25 ng
Nuclease-free Water	9.0 $\mu$ l	

2. Mix reagents completely, then transfer to a thermocycler.
3. Perform the following cycling conditions:

Thermocycling Conditions for a Routine PCR:

STEP	TEMP	TIME
Initial Denaturation	98°C	30 seconds
25 Cycles	98°C	10 seconds
	50–72°C*	10–30 seconds
	72°C	20–30 seconds/kb
Final Extension	72°C	2 minutes
Hold	4–10°C	

\* For a Q5-optimized annealing temperature of mutagenic primers, please use NEBaseChanger, the online NEB primer design software ([www.NEBaseChanger.neb.com](http://www.NEBaseChanger.neb.com)). For pre-designed, back-to-back primer sets, a  $T_a = T_m + 3$  rule can be applied, but optimization may be necessary.

### Step II: Kinase, Ligase & DpnI (KLD) Treatment

1. Assemble the following reagents:

	VOLUME	FINAL CONCENTRATION
PCR Product	1 $\mu$ l	
2X KLD Reaction Buffer	5 $\mu$ l	1X
10X KLD Enzyme Mix	1 $\mu$ l	1X
Nuclease-free Water	3 $\mu$ l	

2. Mix well by pipetting up and down and incubate at room temperature for 5 minutes.

### Step III: Transformation

1. Thaw a 50  $\mu$ l aliquot of chemically competent *E. coli* cells [NEB 5-alpha Competent *E. coli* (High Efficiency), NEB #C2987, are recommended] on ice.
2. Add 5  $\mu$ l of the KLD mix from Step II to the tube of thawed cells. Carefully flick the tube 4–5 times to mix. Do not vortex.
3. Place the mixture on ice for 30 minutes.
4. Heat shock at 42°C for 30 seconds.
5. Place on ice for 5 minutes.
6. Pipette 950  $\mu$ l of room temperature SOC into the mixture.
7. Incubate at 37°C for 60 minutes with shaking (250 rpm).
8. Mix the cells thoroughly by flicking the tube and inverting, then spread 50–100  $\mu$ l onto a selection plate and incubate overnight at 37°C. It may be necessary (particularly for simple substitution and deletion experiments) to make a 10- to 40-fold dilution of the transformation mix in SOC prior to plating, to avoid a lawn of colonies.

## Protocol for Control Reaction

A control mutagenesis reaction is included in this kit, allowing blue/white screening of mutagenesis efficiency. The Control SDM Plasmid is 6.7 kb and contains a premature stop-codon at amino acid #23 in the *lacZα* open-reading frame (1), as well as two BglII sites at positions #1948 and #2841. The Control SDM Primer Mix contains one primer that is 100% complementary to the Control SDM Plasmid and a second mutagenic primer. The mutagenic primer is designed to revert the aforementioned stop-codon to glutamine, thus allowing *lacZα* expression and blue colony formation upon transformation and plating on media containing X-gal and IPTG into *E. coli* containing the *lacZ* deletion mutant, Δ (*lacZ*)M15. The mutagenic primer also removes the BglII site at position #2841, allowing for restriction endonuclease screening of positive colonies. A BglII digestion reaction of the control plasmid will yield two DNA fragments of 893 bp and 5.8 kb. After a successful mutagenesis reaction, only a single DNA fragment of 6.7 kb will be present after BglII digestion.

CONTROL SDM PRIMER MIX	
Forward	AACCCTGGCGtAcCcAACTTAATCG
Reverse	TTCCCAGTCACGACGTTGTA AAA

### Step 1: Exponential Amplification (PCR)

For the control mutagenesis reaction, assemble PCR components as described in the following table:

	25 µl REACTION	FINAL CONCENTRATION
Q5 Hot Start High-Fidelity 2X Master Mix	12.5 µl	1X
Control SDM Primer Mix	1.25 µl	0.5 µM
Control SDM Plasmid	2 µl	10 ng
Nuclease-free Water	9.25 µl	

Thermocycling Conditions for the Control PCR:

STEP	TEMP	TIME
Initial Denaturation	98°C	30 seconds
25 Cycles	98°C	10 seconds
	60°C*	20 seconds
	72°C	1 min 45 sec
Final Extension	72°C	2 minutes
Hold	4–10°C	

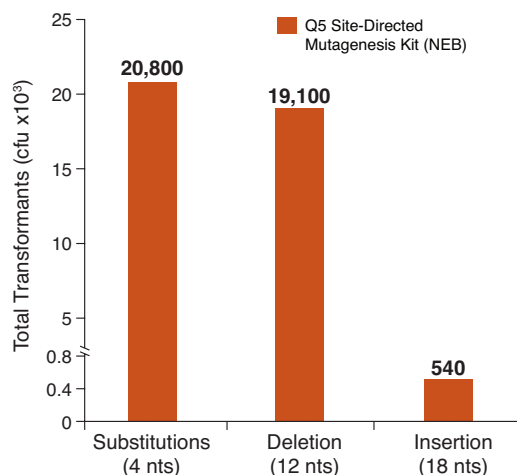
\* Although the control reaction works optimally with a 60°C annealing temperature, it can be successful with an annealing temperature from 54–68°C, allowing the control reaction to be cycled with similar experimental samples.

For Step II (KLD reaction) and Step III (Transformation), follow the protocol found on page 6. Prior to plating, make a 40-fold dilution of cells in room temperature SOC (25 µl of cells and 975 µl of SOC) to avoid a lawn of colonies. For blue/white screening, spread cells on room temperature LB plates containing 100 µg/ml of ampicillin, 40 µg/ml X-gal and 50 µg/ml IPTG (stock solutions of X-gal and IPTG are made in dimethylformamide to 40 mg/ml and 50 mg/ml, respectively). Allow the plates to grow overnight at 37°C. For optimal color development, let plates sit at room temperature or 30°C for an additional 24 hours..

## Expected Results

Exponential amplification with Q5 Hot Start High-Fidelity DNA Polymerase, combined with the intramolecular ligation reaction that precedes high-efficiency transformation, enables high numbers of transformants for simple mutagenesis experiments (substitutions and deletions), and more complex strategies (insertions). In contrast, linear amplification methods typically generate fewer colonies in substitution experiments and are unable to perform large deletions or insertions. Typical results from three different mutagenesis experiments with three different Q5 amplified site-directed templates are shown in Figure 4. For substitutions and deletions, thousands of colonies were observed with > 90% of the colonies having incorporated the desired mutation. An insertion of 18 nucleotides (6 amino acids for the addition of a 6X His tag) at the end of an open-reading frame still generated over 500 colonies. All isolates screened for the proper insertion sequences were found to be correct. As with all PCR experiments, success rate is critically dependent on proper primer design. We highly recommend designing primers using NEBaseChanger, the NEB online primer design software, which can be found at: [NEBaseChanger.neb.com](http://NEBaseChanger.neb.com).

**Figure 4: Typical Results from Substitution, Deletion and Insertion Mutagenesis Reactions using Q5 Site-Directed Mutagenesis Kit.**



Typical results from a substitution reaction (4 nt) using the Control SDM Primer Mix, Control SDM Plasmid (6.7 kb) and NEB 5-alpha Competent *E. coli* (High Efficiency) Cells (NEB #C2987) are shown. For comparison, results from a 12 nt deletion experiment (5.8 kb plasmid) and an 18 nt insertion experiment (7.0 kb plasmid) are also shown. In all three cases, over 90% of the resultant colonies contained the desired mutation(s). Results are normalized to the total number of transformants if cells were not diluted prior to plating.

## Troubleshooting

### No/Low Colonies

- Be sure to use high-efficiency chemically-competent *E. coli* cells. The following competent *E. coli* cells have been shown to work with this kit:

NEB #C2987, NEB 5-alpha (High Efficiency) (standard recommendation)

NEB #C2992, NEB 5-alpha F'Iq (High Efficiency)

NEB #C3019, NEB 10-beta (High Efficiency)

NEB #C2984, NEB Turbo

NEB #C2566, T7 Express

NEB #C3029, Shuffle<sup>®</sup> T7

For convenience, we offer a version of the Q5 Site Directed Mutagenesis Kit prepackaged with NEB #C2987 NEB 5-alpha Competent *E. coli* (High Efficiency) cells in product number NEB #E0554.

Other chemically competent *E. coli* strains suitable for cloning can be substituted. Results will vary according to the quality and efficiency of the cells.

- Check that the transformation efficiency of the competent cells is  $\sim 1 \times 10^9$  colony forming units (cfu) per  $\mu\text{g}$ . To calculate transformation efficiency, transform 2  $\mu\text{l}$  of pUC19 DNA (NEB #N3041) (100 pg) into 50  $\mu\text{l}$  of cells. Follow the transformation protocol on page 6. Prior to plating, dilute 10  $\mu\text{l}$  of cells up to 1 ml in SOC. Plate 100  $\mu\text{l}$  of this dilution. In this case, 150 colonies will yield a transformation efficiency of  $1.5 \times 10^9$  cfu/ $\mu\text{g}$  ( $\mu\text{g DNA}=0.0001$ , dilution= $10/1000 \times 100/1000$ ).
- Ensure that your primers are designed properly. To take advantage of the exponential nature of the amplification reaction, the 5' ends of the two primers should align back-to-back unless deletions are being made (see Figure 3). For best results, primers should be designed and annealing temperatures calculated using NEBaseChanger, the NEB online primer design software, which can be found at: [NEBaseChanger.neb.com](http://NEBaseChanger.neb.com).
- Ensure there is a clean PCR product by visualizing 2–5  $\mu\text{l}$  of the reaction on an agarose gel. Follow the suggestions below for low or impure PCR products.
- Only use 1  $\mu\text{l}$  of PCR product in the KLD reaction. Carrying too much PCR product forward can decrease transformation efficiency. If the PCR yield is low, more product can be added to the KLD reaction, however a buffer exchange step, such as PCR purification, must be included prior to transformation.
- Only use 5  $\mu\text{l}$  of the KLD reaction in the transformation. If more KLD reaction is added, a buffer exchange step, such as PCR purification, should be included prior to transformation.
- Ensure that the selectable marker in the plasmid matches the selection agent used in the plates
- Ensure the *E. coli* cells have been stored at  $-80^\circ\text{C}$ .



### No/Low PCR Product

- Ensure that the optimal annealing temperature ( $T_a$ ) is used. High-Fidelity polymerases benefit from a  $T_m+3$  annealing temp. Use NEBaseChanger, the NEB online primer design software, to calculate  $T_a$  ([NEBaseChanger.neb.com](http://NEBaseChanger.neb.com)). Alternatively, the optimal annealing temperature could be determined using a gradient PCR followed by agarose gel analysis.
- Ensure that the elongation time is adequate for the plasmid length. We recommend 10–20 seconds per kb of plasmid.
- Ensure that the final concentration of each primer is 0.5  $\mu\text{M}$ .
- Purify the primers with polyacrylamide gel electrophoresis (PAGE).

### Resulting Plasmids Do Not Contain the Desired Mutation

- Ensure proper design of the mutagenic primers
- Optimize the PCR conditions (see above)
- Use 1–25 ng of template in the PCR step. A small increase in the number of clones with no/incorrect mutation incorporated can occur if less than 1 ng or more than 25 ng of template is used.

### Quality Controls

Product specifications for individual components in the Q5 Site-Directed Mutagenesis Kit (Without Competent Cells) are available separately. The Q5 Site-Directed Mutagenesis Kit (Without Competent Cells) has been tested using the control plasmid and primer mix provided in the kit.

### Reference

1. Kalnins et al., (1983) *The EMBO Journal* 2, 593-597.

### Ordering Information

NEB #	PRODUCT	SIZE
E0552S	Q5 Site-Directed Mutagenesis Kit (Without Competent Cells)	10 reactions

#### COMPANION PRODUCTS

NEB #	PRODUCT	SIZE
E0554S	Q5 Site-Directed Mutagenesis Kit	10 reactions
M0494S/L	Q5 Hot Start High-Fidelity 2X Master Mix	100/500 reactions
C29871/H	NEB 5-alpha Competent <i>E. coli</i> (High Efficiency)	6 x 0.2 ml/ 20 x 0.05 ml
B9020S	SOC Outgrowth Medium	4 x 25 ml

## Revision History

REVISION #	DESCRIPTION	DATE
1.0	N/A	3/14
2.0	New Format applied	7/20

This product is intended for research purposes only. This product is not intended to be used for therapeutic or diagnostic purposes in humans or animals.

This product is covered by one or more patents, trademarks and/or copyrights owned or controlled by New England Biolabs, Inc. For more information about commercial rights, please email us at [busdev@neb.com](mailto:busdev@neb.com) While NEB develops and validates its products for various applications, the use of this product may require the buyer to obtain additional third party intellectual property rights for certain applications.

This product is licensed for research and commercial use from Bio-Rad Laboratories, Inc., under U.S. Pat. Nos. 6,627,424, 7,541,170, 7,670,808, 7,666,645, and corresponding patents in other countries. No rights are granted for use of the product for Digital PCR or real-time PCR applications, with the exception of quantification in Next Generation Sequencing workflows.

B CORPORATION™ is a trademark of B Lab, Inc.

© Copyright 2020, New England Biolabs, Inc.; all rights reserved



be INSPIRED  
drive DISCOVERY  
stay GENUINE