

QuikChange XL Site-Directed Mutagenesis Kit

Instruction Manual

Catalog # 200516 (30 reactions) and 200517 (10 reactions)

Revision F1

For Research Use Only. Not for use in diagnostic procedures.

200516-12



Agilent Technologies

LIMITED PRODUCT WARRANTY

This warranty limits our liability to replacement of this product. No other warranties of any kind, express or implied, including without limitation, implied warranties of merchantability or fitness for a particular purpose, are provided by Agilent. Agilent shall have no liability for any direct, indirect, consequential, or incidental damages arising out of the use, the results of use, or the inability to use this product.

ORDERING INFORMATION

Please visit www.agilent.com.

TECHNICAL SERVICES

United States and Canada

Email: genomics@agilent.com

Telephone: 800 227 9770 (option 3,4,3)

All Other Locations

Please visit www.agilent.com/en/contact-us/page.

QuikChange XL Site-Directed Mutagenesis Kit

CONTENTS

Materials Provided	1
Storage Conditions	1
Additional Materials Required	1
Introduction	2
QuikChange XL Mutagenesis Control	4
Mutagenic Primer Design	5
XL10-Gold Ultracompetent Cells	7
Protocol	8
QuikSolution	8
Mutant Strand Synthesis Reaction (Thermal Cycling).....	8
<i>Dpn</i> I Digestion of the Amplification Products	10
Transformation of XL10-Gold Ultracompetent Cells	10
Transformation Guidelines	13
Storage Conditions	13
Aliquoting Cells	13
Use of 14-ml Falcon Round-Bottom Polypropylene Tubes.....	13
Use of β -Mercaptoethanol	13
Quantity of DNA Added	13
Length and Temperature of the Heat Pulse.....	13
Preparing the Agar Plates for Color Screening.....	13
Troubleshooting	14
Preparation of Media and Reagents	15
References	16
Endnotes	16
MSDS Information	16
Quick-Reference Protocol	17

QuikChange XL Site-Directed Mutagenesis Kit

MATERIALS PROVIDED

Materials provided	Quantity	
	Catalog #200516 ^a 30 reactions	Catalog #200517 ^b 10 reactions
<i>PfuTurbo</i> DNA polymerase (2.5 U/μl)	80 U	25 U
10× reaction buffer ^c	500 μl	500 μl
<i>Dpn</i> I restriction enzyme (10 U/μl)	300 U	100 U
Oligonucleotide control primer #1 [34-mer (100 ng/μl)] 5' CCA TGA TTA CGC CAA GCG CGC AAT TAA CCC TCA C 3'	750 ng	750 ng
Oligonucleotide control primer #2 [34-mer (100 ng/μl)] 5' GTG AGG GTT AAT TGC GCG CTT GGC GTA ATC ATG G 3'	750 ng	750 ng
pWhitescript 4.5-kb control plasmid (5 ng/ μl)	50 ng	50 ng
QuikSolution	500 μl	500 μl
dNTP mix ^{d,e}	30 μl	10 μl
XL10-Gold ultracompetent cells ^f (yellow tubes)	10 × 135 μl	4 × 135 μl
XL10-Gold β-mercaptoethanol mix (β-ME)	2 × 50 μl	50 μl
pUC18 control plasmid (0.1 ng/μl in TE buffer ^c)	10 μl	10 μl

^a The QuikChange XL Site-Directed Mutagenesis Kit (Catalog #200516) contains enough reagents for 30 reactions total (control and experimental reactions combined).

^b The QuikChange XL Site-Directed Mutagenesis Kit (Catalog #200517) contains enough reagents for 10 reactions total (control and experimental reactions combined).

^c See Preparation of Media and Reagents.

^d Thaw the dNTP mix once, prepare single-use aliquots, and store the aliquots at –20°C. **Do not subject the dNTP mix to multiple freeze-thaw cycles.**

^e **The composition of the dNTP mix is proprietary.** This reagent has been optimized for the QuikChange XL site-directed mutagenesis protocols and has been qualified for use in conjunction with the other kit components. Do not substitute with dNTP mixes provided with other Agilent kits.

^f Genotype: Tet^rΔ (*mcrA*)183 Δ(*mcrCB-hsdSMR-mrr*)173 *endA1 supE44 thi-1 recA1 gyrA96 relA1 lac* Hte [F' *proAB lac*^qZΔM15 Tn10 (Tet^r) Amy Cam^r]. Available separately as catalog # 200314 and # 200315.

STORAGE CONDITIONS

XL10-Gold Ultracompetent cells, XL10-Gold β-ME, and pUC18 Control Plasmid: –80°C
All Other Components: –20°C

ADDITIONAL MATERIALS REQUIRED

14-ml Falcon round-bottom polypropylene tubes (Thermo Fisher Scientific p/n 352059)
 5-Bromo-4-chloro-3-indolyl-β-D-galactopyranoside (X-gal)
 Isopropyl-1-thio-β-D-galactopyranoside (IPTG)

INTRODUCTION

In vitro site-directed mutagenesis is an invaluable technique for studying protein structure-function relationships and gene expression, and for carrying out vector modification. Several approaches to this technique have been published, but these methods generally require single-stranded DNA (ssDNA) as the template¹⁻⁴ and are labor intensive or technically difficult. The QuikChange XL Site-Directed Mutagenesis Kit allows site-specific mutation in virtually any double-stranded plasmid, thus eliminating the need for subcloning into M13-based bacteriophage vectors and for ssDNA rescue.⁵ In addition, the QuikChange XL system requires no specialized vectors, unique restriction sites, or multiple transformations. This rapid four-step procedure generates mutants with greater than 80% efficiency. The protocol is simple and uses either miniprep plasmid DNA or cesium-chloride-purified DNA.

The QuikChange XL system is used to make point mutations, switch amino acids, and delete or insert single or multiple amino acids. The QuikChange XL method is performed using *PfuTurbo* DNA Polymerase and a thermal temperature cycler. *PfuTurbo* DNA polymerase replicates both plasmid strands with high fidelity and without displacing the mutant oligonucleotide primers. The basic procedure utilizes a supercoiled double-stranded DNA (dsDNA) vector with an insert of interest and two synthetic oligonucleotide primers containing the desired mutation (see Figure 1). The oligonucleotide primers, each complementary to opposite strands of the vector, are extended during temperature cycling by using *PfuTurbo* DNA polymerase. Incorporation of the oligonucleotide primers generates a mutated plasmid containing staggered nicks. Following temperature cycling, the product is treated with *Dpn* I. The *Dpn* I endonuclease (target sequence: 5'-Gm⁶ATC-3') is specific for methylated and hemimethylated DNA and is used to digest the parental DNA template and to select for mutation-containing synthesized DNA.⁶ DNA isolated from almost all *E. coli* strains is *dam* methylated and therefore susceptible to *Dpn* I digestion. The nicked vector DNA incorporating the desired mutations is then transformed into XL10-Gold Ultracompetent Cells. The small amount of starting DNA template required to perform this method, the high-fidelity of the *PfuTurbo* DNA polymerase, and the low number of thermal cycles all contribute to the high mutation efficiency and decreased potential for random mutations.

Note *While plasmid DNA isolated from almost all of the commonly used E. coli strains (dam⁺) is methylated and is a suitable template for mutagenesis, plasmid DNA isolated from the exceptional dam⁻ E. coli strains, including JM110 and SCS110, is not suitable.*

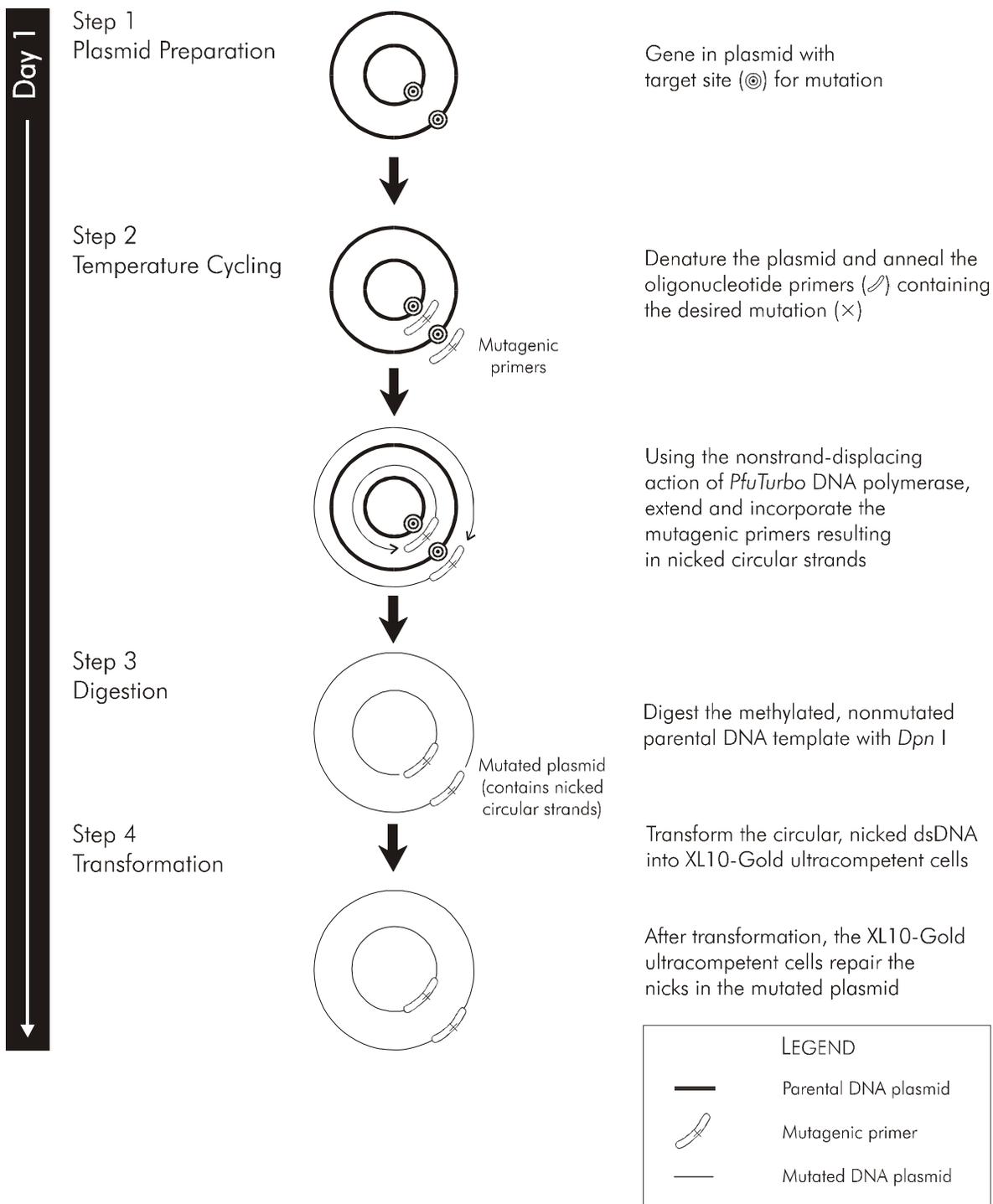


FIGURE 1 Overview of the QuikChange XL site-directed mutagenesis method.

The QuikChange XL site-directed mutagenesis kit is a specialized version of our popular QuikChange site-directed mutagenesis kit, created for efficient mutagenesis of large or otherwise difficult-to-mutagenize plasmid templates. The QuikChange XL kit features components specifically designed for more efficient DNA replication and bacterial transformation. The QuikChange solution is provided to facilitate replication of large plasmids, while XL10-Gold ultracompetent cells have been included to ensure the highest transformation efficiencies possible. The transformation efficiency of XL10-Gold cells is 5-fold higher than the transformation efficiency of XL1-Blue cells employed in the original QuikChange kit.⁷ Moreover, XL10-Gold cells contain the Hte phenotype, which increases the transformation efficiency of larger DNA plasmids.

QUIKCHANGE XL MUTAGENESIS CONTROL

To demonstrate the effectiveness of the QuikChange XL method, the pWhitescript 4.5-kb control plasmid is used to test the efficiency of mutant plasmid generation. The pWhitescript 4.5-kb control plasmid contains a stop codon (TAA) at the position where a glutamine codon (CAA) would normally appear in the β -galactosidase gene of the pBluescript II SK(-) phagemid (corresponding to amino acid 9 of the protein). XL10-Gold ultracompetent cells transformed with this control plasmid appear white on LB-ampicillin agar plates (see *Preparation of Media and Reagents*), containing IPTG and X-gal, because β -galactosidase activity has been obliterated. The oligonucleotide control primers create a point mutation that reverts the T residue of the stop codon (TAA) in the β -galactosidase gene encoded on the pWhitescript 4.5-kb control template to a C residue to produce a glutamine codon (Gln, CAA). Following transformation, colonies can be screened for β -galactosidase production (β -gal⁺) by virtue of a blue colony phenotype.

MUTAGENIC PRIMER DESIGN

Consider the following guidelines when designing mutagenic primers:

1. Agilent recommends the use of our web-based primer design tool, (<https://www.agilent.com/store/primerDesignProgram.jsp>) which was optimized to design mutagenic primer sequences specifically for the QuikChange kits.

Please note that the tool only designs one pair of primers. If the primer sequences need to be modified (e.g., due to predicted secondary structure formation), then you may need to use an alternate primer design tool or manually manipulate the sequences.

2. Both mutagenic primers must contain the desired mutation and anneal to the same sequence on opposite strands of the plasmid.
3. Primers should be between 25 and 45 bases in length with a melting temperature (T_m) of $\geq 78^\circ\text{C}$. Optimum primer sets for simultaneous mutagenesis should have similar melting temperatures. If necessary, primers can be longer than 45 bases to achieve a $T_m \geq 78^\circ\text{C}$. However, using longer primers increases the likelihood of secondary structure formation, which may affect the efficiency of the mutagenesis reaction. Use the following formula to estimate the T_m of primers:

$$T_m = 81.5 + 0.41(\%GC) - (675/N) - \% \text{ mismatch}$$

For calculating T_m :

- N is the primer length in bases.
- values for **%GC** and **% mismatch** are whole numbers

For calculating T_m for primers intended to introduce insertions or deletions, use this modified version of the above formula:

$$T_m = 81.5 + 0.41(\%GC) - (675/N)$$

where N does not include the bases which are being inserted or deleted.

4. The desired mutation should be located near the center of the primer, at least 10 bases from either end. The maximum number of mismatches is 3 bases in a row or 2 bases with a maximum of 9 bases in between.
5. Ideally, the primers have a minimum GC content of 40% and terminate in one or more C or G bases at the 3' end.
6. An online oligo sequence analysis tool can be used to check the primer sequences for potential secondary structure (i.e., hairpin) formation at or above the annealing temperature of 55°C . Primers predicted to form stable hairpin structures at 55°C or above may need to be redesigned or modified (manually or using a primer design tool).

7. To avoid primer dimer formation, the ΔG of the primer must be greater (i.e., closer to zero) than -9 kcal/mole (-9 kJ). The formation of primer dimers (hetero dimers and self dimers) can affect the efficiency of linear amplification during the mutant strand synthesis reaction.
8. Similarly, binding between the primer and off-target sites on the plasmid can also affect the mutant strand synthesis reaction. The NCBI blastn alignment tool (or similar alignment tool) can identify potential off-target binding between the primer and plasmid.
9. Primers should not be methylated.
10. Primers do not require 5' phosphorylation, but they must be purified by HPLC, FPLC, or PAGE. Failure to adequately purify the primers significantly decreases mutation efficiency.

XL10-GOLD ULTRACOMPETENT CELLS

XL10-Gold ultracompetent cells are derived from the highest-efficiency Agilent competent cell line, XL2-Blue MRF'. These strains possess the Hte phenotype, which increases transformation efficiency of ligated DNA.⁷ XL10-Gold cells are both endonuclease deficient (*endA1*) and recombination deficient (*recA*). The *endA1* mutation greatly improves the quality of plasmid miniprep DNA,⁸ and the *recA* mutation helps ensure insert stability. In addition, the McrA, McrCB, McrF, Mrr, and HsdR systems have been removed from XL10-Gold ultracompetent cells. The *mcrA*, *mcrCB* and *mrr* mutations prevent cleavage of cloned DNA that carries cytosine and/or adenine methylation, which is often present in eukaryotic DNA and cDNA.^{9, 10, 11} The McrA and McrCB systems recognize and restrict methylated cytosine DNA sequences, and the Mrr system recognizes and restricts methylated adenine DNA sequences. The Mrr system also restricts methylated cytosine DNA sequences with a specificity differing from that of McrA and McrCB. This activity has been named McrF. This McrF activity against methylated cytosines has been shown to be equal to or greater than the restriction activity of the McrA and McrCB systems.¹² The *hsdR* mutation prevents the cleavage of cloned DNA by the *EcoK* (*hsdR*) endonuclease system. XL10-Gold cells grow faster than XL1 or XL2-Blue cells, resulting in larger colonies. To permit blue-white color screening, the XL10-Gold ultracompetent cells contain the *lacI^qZΔM15* gene on the F' episome.

Host strain	References	Genotype
XL10-Gold ultracompetent cells	7, 13, 14	Tet ^R Δ(<i>mcrA</i>)183 Δ(<i>mcrCB</i> - <i>hsdSMR</i> - <i>mrr</i>)173 <i>endA1 supE44 thi-1 recA1 gyrA96 relA1 lac Hte</i> [F' <i>proAB lacI^qZΔM15</i> Tn10 (Tet ^R) Amy Cam ^R]

It is important to store the XL10-Gold ultracompetent cells at -80°C to prevent a loss of efficiency. For best results, please follow the directions outlined in the following sections.

PROTOCOL

QuikSolution

QuikSolution has been shown to improve linear amplification. Enhanced amplification efficiencies are observed when using between 2.5–3.5 μl QuikSolution/50 μl reaction, with 3 μl being optimal for most targets.

Mutant Strand Synthesis Reaction (Thermal Cycling)

Notes *Ensure that the plasmid DNA template is isolated from a dam^+ E. coli strain. The majority of the commonly used E. coli strains are dam^+ . Plasmid DNA isolated from dam^- strains (e.g. JM110 and SCS110) is not suitable.*

*To maximize temperature-cycling performance, use **thin-walled tubes**, which ensure ideal contact with the heat block. The following protocols were optimized using thin-walled tubes.*

Keep the primer concentration in excess relative to the ds-DNA template. When optimizing reaction conditions, vary the amount of template while keeping the primer constantly in excess.

1. Synthesize two complimentary oligonucleotides containing the desired mutation, flanked by unmodified nucleotide sequence. Purify these oligonucleotide primers prior to use in the following steps (see *Mutagenic Primer Design*).

2. Prepare the control reaction as indicated below:

5 μl of 10 \times reaction buffer (see *Preparation of Media and Reagents*)

2 μl (10 ng) of pWhitescript 4.5-kb control plasmid (5 ng/ μl)

1.25 μl (125 ng) of oligonucleotide control primer #1
[34-mer (100 ng/ μl)]

1.25 μl (125 ng) of oligonucleotide control primer #2
[34-mer (100 ng/ μl)]

1 μl of dNTP mix

3 μl of QuikSolution

36.5 μl of double-distilled water (ddH_2O) to a final volume of 50 μl

Then add

1 μl of *PfuTurbo* DNA polymerase (2.5 U/ μl)

3. Prepare the sample reaction(s) as indicated below:

Note Set up an initial sample reaction using 10 ng of dsDNA template. If this initial reaction is unsuccessful, set up a series of sample reactions using various concentrations of dsDNA template ranging from 5 to 50 ng (e.g., 5, 10, 20, and 50 ng of dsDNA template) while keeping the primer concentration constant.

5 μ l of 10 \times reaction buffer
 X μ l (10 ng) of dsDNA template
 X μ l (125 ng) of oligonucleotide primer #1
 X μ l (125 ng) of oligonucleotide primer #2
 1 μ l of dNTP mix
 3 μ l of QuikSolution
 ddH₂O to a final volume of 50 μ l

Then add

1 μ l of *PfuTurbo* DNA polymerase (2.5 U/ μ l)

4. Cycle each reaction using the cycling parameters outlined in Table I.

Note It is important to adhere to the 18-cycle limit when cycling the mutagenesis reactions. More than 18 cycles can have deleterious effects on the reaction efficiency.

5. Following temperature cycling, place the reaction tubes on ice for 2 minutes to cool the reactions to $\leq 37^\circ\text{C}$.

Note If desired, amplification may be checked by electrophoresis of 10 μ l of the product on a 1% agarose gel. A band may or may not be visualized at this stage. In either case proceed with Dpn I digestion and transformation.

TABLE I

Cycling Parameters for the QuikChange XL Method

Segment	Cycles	Temperature	Time
1	1	95°C	1 minute
2	18	95°C	50 seconds
		60°C	50 seconds
		68°C	1 minute/kb of plasmid length
3	1	68°C	7 minutes

* For example, a 5-kb plasmid requires 5 minutes at 68°C per cycle.

Dpn I Digestion of the Amplification Products

1. Add 1 μl of the *Dpn* I restriction enzyme (10 U/ μl) to each amplification reaction.
2. Gently and thoroughly mix each reaction mixture by pipetting the solution up and down several times. Spin down the reaction mixtures in a microcentrifuge for 1 minute, then immediately incubate the reactions at 37°C for 1 hour to digest the parental (i.e., the nonmutated) supercoiled dsDNA.

Transformation of XL10-Gold Ultracompetent Cells

Notes *Please read the Transformation Guidelines before proceeding with the transformation protocol.*

XL10-Gold cells are resistant to tetracycline and chloramphenicol. If the mutagenized plasmid contains only the tet^R or cam^R resistance marker, an alternative strain of competent cells must be used.

1. Gently thaw the XL10-Gold ultracompetent cells on ice. For each control and sample reaction to be transformed, aliquot 45 μl of the ultracompetent cells to a *prechilled* 14-ml Falcon round-bottom polypropylene tube.
2. Add 2 μl of the β -ME mix provided with the kit to the 45 μl of cells. (Using an alternative source of β -ME may reduce transformation efficiency.)
3. Swirl the contents of the tube gently. Incubate the cells on ice for 10 minutes, swirling gently every 2 minutes.
4. Transfer 2 μl of the *Dpn* I-treated DNA from each control and sample reaction to separate aliquots of the ultracompetent cells.

As an optional control, verify the transformation efficiency of the XL10-Gold ultracompetent cells by adding 1 μl of 0.01 ng/ μl pUC18 control plasmid (dilute the control provided 1:10 in high-quality water) to another 45- μl aliquot of cells.

Swirl the transformation reactions gently to mix and incubate the reactions on ice for 30 minutes.

5. Preheat NZY⁺ broth (see *Preparation of Media and Reagents*) in a 42°C water bath for use in step 8.

Note *Transformation of XL10-Gold ultracompetent cells has been optimized using NZY⁺ broth.*

- Heat-pulse the tubes in a 42°C water bath for 30 seconds. The duration of the heat pulse is *critical* for obtaining the highest efficiencies. Do not exceed 42°C.

Note *This heat pulse has been optimized for transformation in 14-ml Falcon round-bottom polypropylene tubes.*

- Incubate the tubes on ice for 2 minutes.
- Add 0.5 ml of preheated (42°C) NZY⁺ broth to each tube, then incubate the tubes at 37°C for 1 hour with shaking at 225–250 rpm.
- Plate the appropriate volume of each transformation reaction, as indicated in the table below, on agar plates containing the appropriate antibiotic for the plasmid vector.

For the mutagenesis and transformation controls, spread cells on LB–ampicillin agar plates containing 80 µg/ml X-gal and 20 mM IPTG (see *Preparing the Agar Plates for Color Screening*).

Transformation reaction plating volumes

Reaction Type	Volume to Plate
pWhitescript mutagenesis control	250 µl
pUC18 transformation control	5 µl (in 200 µl of NZY+ broth)*
Sample mutagenesis	250 µl on each of two plates (entire transformation reaction)

* Place a 200-µl pool of NZY⁺ broth on the agar plate, pipet the 5 µl of the transformation reaction into the pool, then spread the mixture.

- Incubate the transformation plates at 37°C for >16 hours.

Expected Results for the Control Transformations

The expected colony number from the transformation of the pWhitescript 4.5 kb control mutagenesis reaction is between 50 and 800 colonies. Greater than 80% of the colonies should contain the mutation and appear as blue colonies on agar plates containing IPTG and X-gal.

Note *The mutagenesis efficiency (ME) for the pWhitescript 4.5-kb control plasmid is calculated by the following formula:*

$$ME = \frac{\text{Number of blue colony forming units (cfu)}}{\text{Total number of colony forming units (cfu)}} \times 100\%$$

If transformation of the pUC18 control plasmid was performed, >100 colonies (>10⁹ cfu/µg) should be observed, with >98% having the blue phenotype.

Expected Results for Sample Transformations

The expected colony number is between 10 and 1000 colonies, depending upon the base composition and length of the DNA template employed. For suggestions on increasing colony number, see *Troubleshooting*. The insert of interest should be sequenced to verify that selected clones contain the desired mutation(s).

TRANSFORMATION GUIDELINES

Storage Conditions

Ultracompetent cells are sensitive to even small variations in temperature and must be stored at the bottom of a -80°C freezer. Transferring tubes from one freezer to another may result in a loss of efficiency. Ultracompetent cells should be placed at -80°C directly from the dry ice shipping container.

Aliquoting Cells

When aliquoting, keep ultracompetent cells on ice at all times. It is essential that the Falcon polypropylene tubes are placed on ice before the cells are thawed and that the cells are aliquoted directly into the prechilled tubes.

Use of 14-ml Falcon Round-Bottom Polypropylene Tubes

It is important that 14-ml Falcon round-bottom polypropylene tubes (Thermo Fisher Scientific p/n 352059) are used for the transformation protocol, since other tubes may be degraded by the β -mercaptoethanol used in the *Transformation Protocol*. In addition, the duration of the heat-pulse step is critical and has been optimized specifically for the thickness and shape of these tubes.

Use of β -Mercaptoethanol

β -Mercaptoethanol (β -ME) has been shown to increase transformation efficiency. The XL10-Gold β -mercaptoethanol mix provided in this kit is diluted and ready to use.

Quantity of DNA Added

Greatest efficiencies are observed when adding 2 μl of the ligation mixture. A greater number of colonies will be obtained when adding up to 50 ng, although the overall efficiency may be lower.

Length and Temperature of the Heat Pulse

There is a defined window of highest efficiency resulting from the heat pulse during transformation. Optimal efficiencies are observed when cells are heat-pulsed for 30 seconds. Heat-pulsing for at least 30 seconds is recommended to allow for slight variations in the length of incubation. Efficiencies decrease when incubating for <30 seconds or for >40 seconds. Do not exceed 42°C .

Preparing the Agar Plates for Color Screening

To prepare the LB agar plates for blue-white color screening, add 80 $\mu\text{g/ml}$ of 5-bromo-4-chloro-3-indolyl- β -D-galactopyranoside (X-gal), 20 mM isopropyl-1-thio- β -D-galactopyranoside (IPTG), and the appropriate antibiotic to the LB agar. Alternatively, 100 μl of 10 mM IPTG and 100 μl of 2% X-gal can be spread on the LB agar plates 30 minutes prior to plating the transformations. Prepare the IPTG in sterile dH_2O ; prepare the X-gal in dimethylformamide (DMF). Do not mix the IPTG and X-gal before pipetting them onto the plates because these chemicals may precipitate.

TROUBLESHOOTING

When used according to the guidelines outlined in this instruction manual, this kit provides a reliable means to conduct site-directed mutagenesis using dsDNA templates. Variations in the base composition and length of the DNA template and in thermal cycler performance may contribute to differences in mutagenesis efficiency. We provide the following guidelines for troubleshooting these variations.

Observation	Suggestion(s)
Low transformation efficiency or low colony number	Ensure that sufficient DNA template is used in the reaction. Visualize the DNA template on a gel to verify the quantity and quality. Repeat reaction using higher amounts of plasmid DNA (100 ng, 200 ng, 500 ng).
	Ensure that sufficient mutant DNA is synthesized in the reaction. <ul style="list-style-type: none"> • Titrate QuikSolution in 1-μl increments from 0 to 5 μl • Increase the amount of the <i>Dpn</i> I-treated DNA used in the transformation reaction to 4 μl • Increase the extension time to 2.5 min/kb • Precipitate the entire reaction and use all of it in the transformation
	Ensure sufficient mutant DNA is synthesized by adjusting the cycling parameters for the sample reaction to overcome differences in ramping efficiencies of thermal cyclers. Increase initial denaturation step (segment 1) to 1–2 minutes and denaturation cycles (segment 2) to 1 minute.
	Ethanol precipitate the <i>Dpn</i> I digested PCR product, and resuspend in a decreased volume of water before transformation.
Low mutagenesis efficiency or low colony number with the control reaction	Different thermal cyclers contribute to variations in cycling efficiencies. Optimize the cycling parameters (including ramp rates) for the control reaction then repeat the protocol for the sample reactions using the optimized conditions.
	Ensure that ultracompetent cells are stored at the bottom of a –80°C freezer immediately upon arrival; use XL10-Gold β -ME in the transformation reactions (see also <i>Transformation Guidelines</i>).
	Verify that the agar plates were prepared correctly. See <i>Preparing the Agar Plates for Color Screening</i> , and follow the recommendations for IPTG and X-Gal concentrations carefully.
	For best visualization of the blue (β -gal ⁺) phenotype, the control plates must be incubated for at least 16 hours at 37°C.
	Avoid multiple freeze-thaw cycles for the dNTP mix. Thaw the dNTP mix once, prepare single-use aliquots, and store the aliquots at –20°C. Do not subject the dNTP mix to multiple freeze-thaw cycles.
Low mutagenesis efficiency with the sample reaction(s)	Allow sufficient time for the <i>Dpn</i> I to completely digest the parental template; repeat the digestion if too much DNA template was present. Increase digestion time to 1.5–2.0 hours.
	Avoid multiple freeze-thaw cycles for the dNTP mix. Thaw the dNTP mix once, prepare single-use aliquots, and store the aliquots at –20°C. Do not subject the dNTP mix to multiple freeze-thaw cycles.
	The formation of secondary structures may be inhibiting the mutagenesis reaction. Increasing the annealing temperature up to 68°C may help to alleviate secondary structure formation and improve mutagenesis efficiency.

Table continues on following page

Table continues from the previous page

False positives	Poor quality primers can lead to false positives. Radiolabel the primers and check for degradation on an acrylamide gel or resynthesize the primers.
	False priming can lead to false positives. Increase the stringency of the reaction by increasing the annealing temperature up to 68°C.
Unwanted deletion or recombination of plasmid DNA following mutagenesis and transformation	Transform the mutagenesis reaction into competent cells that are designed to prevent recombination events, such as Agilent's SURE 2 Supercompetent Cells (Catalog #200152). Note that SURE 2 competent cells are not recommended for use with mutagenized plasmids greater than 10 kb in size; note also that SURE 2 cells are Kan ^r , Tet ^r , and Chl ^r , and are not compatible with plasmid selection using kanamycin, tetracycline, or chloramphenicol resistance markers.

PREPARATION OF MEDIA AND REAGENTS

<p>LB Agar (per Liter) 10 g of NaCl 10 g of tryptone 5 g of yeast extract 20 g of agar Add deionized H₂O to a final volume of 1 liter Adjust pH to 7.0 with 5 N NaOH Autoclave Pour into petri dishes (~25 ml/100-mm plate)</p>	<p>LB-Ampicillin Agar (per Liter) 1 liter of LB agar, autoclaved Cool to 55°C Add 10 ml of 10-mg/ml filter-sterilized ampicillin Pour into petri dishes (~25 ml/100-mm plate)</p>
<p>NZY⁺ Broth (per Liter) 10 g of NZ amine (casein hydrolysate) 5 g of yeast extract 5 g of NaCl Add deionized H₂O to a final volume of 1 liter Adjust to pH 7.5 using NaOH Autoclave Add the following filter-sterilized supplements prior to use: 12.5 ml of 1 M MgCl₂ 12.5 ml of 1 M MgSO₄ 20 ml of 20% (w/v) glucose (or 10 ml of 2 M glucose)</p>	<p>10× Reaction Buffer 100 mM KCl 100 mM (NH₄)₂SO₄ 200 mM Tris-HCl (pH 8.8) 20 mM MgSO₄ 1% Triton[®] X-100 1 mg/ml nuclease-free bovine serum albumin (BSA)</p> <p>TE Buffer 10 mM Tris-HCl (pH 7.5) 1 mM EDTA</p>

REFERENCES

1. Kunkel, T. A. (1985) *Proc Natl Acad Sci U S A* 82(2):488-92.
2. Sugimoto, M., Esaki, N., Tanaka, H. and Soda, K. (1989) *Anal Biochem* 179(2):309-11.
3. Taylor, J. W., Ott, J. and Eckstein, F. (1985) *Nucleic Acids Res* 13(24):8765-85.
4. Vandeyar, M. A., Weiner, M. P., Hutton, C. J. and Batt, C. A. (1988) *Gene* 65(1):129-33.
5. Papworth, C., Bauer, J. C., Braman, J. and Wright, D. A. (1996) *Strategies* 9(3):3-4.
6. Nelson, M. and McClelland, M. (1992) *Methods Enzymol* 216:279-303.
7. Jerpseth, B., Callahan, M. and Greener, A. (1997) *Strategies* 10(2):37-38.
8. Wnendt, S. (1994) *Biotechniques* 17(2):270, 272.
9. Kohler, S. W., Provost, G. S., Kretz, P. L., Dyaico, M. J., Sorge, J. A. *et al.* (1990) *Nucleic Acids Res* 18(10):3007-13.
10. Kretz, P. L., Kohler, S. W. and Short, J. M. (1991) *J Bacteriol* 173(15):4707-16.
11. Raleigh, E. A. and Wilson, G. (1986) *Proc Natl Acad Sci U S A* 83(23):9070-4.
12. Jerpseth, B., Greener, A., Short, J. M., Viola, J. and Kretz, P. L. (1992) *Strategies* 5(3):81-83.
13. Bullock, W. O., Fernandez, J. M. and Short, J. M. (1987) *Biotechniques* 5(4):376-378.
14. Greener, A. and Jerpseth, B. (1993) *Strategies* 6(2):57.
15. Choi, J. H., Oh, E. S., & Oh, M. H. (2021) *Genes & genomics*, 43(11), 1269-1276. <https://doi.org/10.1007/s13258-021-01148-2>

ENDNOTES

Triton® is a registered trademark of Rohm and Haas Co.

MSDS INFORMATION

Material Safety Data Sheets (MSDSs) are provided online at www.agilent.com. MSDS documents are not included with product shipments.

QuikChange XL Site-Directed Mutagenesis Kit

Catalog #200516 and #200517

QUICK-REFERENCE PROTOCOL

- Prepare the control and sample reaction(s) as indicated below:

Note Set up an initial sample reaction using 10 ng of dsDNA template. If this initial sample reaction is unsuccessful, set up a series of reactions using various concentrations of dsDNA template ranging from 5 to 50 ng (e.g., 5, 10, 20, and 50 ng of dsDNA template) while keeping the primer concentration constant.

Control Reaction

5 μ l of 10 \times reaction buffer
2 μ l (10 ng) of pWhitescript 4.5-kb control template (5 ng/ μ l)
1.25 μ l (125 ng) of oligonucleotide control primer #1 [34-mer (100 ng/ μ l)]
1.25 μ l (125 ng) of oligonucleotide control primer #2 [34-mer (100 ng/ μ l)]
1 μ l of dNTP mix
3 μ l of QuikSolution
36.5 μ l ddH₂O to a final volume of 50 μ l

Sample Reaction

5 μ l of 10 \times reaction buffer
X μ l (10 ng) of dsDNA template
X μ l (125 ng) of oligonucleotide primer #1
X μ l (125 ng) of oligonucleotide primer #2
1 μ l of dNTP mix
3 μ l of QuikSolution
ddH₂O to a final volume of 50 μ l

- Then add 1 μ l of *PfuTurbo* DNA polymerase (2.5 U/ μ l) to each control and sample reaction.
- Cycle each reaction using the cycling parameters outlined in the following table:

Segment	Cycles	Temperature	Time
1	1	95°C	1 minute
2	18	95°C	50 seconds
		60°C	50 seconds
		68°C	1 minute/kb of plasmid length
3	1	68°C	7 minutes

- Add 1 μ l of *Dpn* I restriction enzyme (10 U/ μ l).
- Gently and thoroughly mix each reaction, spin down in a microcentrifuge for 1 minute, and immediately incubate at 37°C for 1 hour to digest the parental supercoiled dsDNA.
- Transform 2 μ l of the *Dpn* I-treated DNA from each control and sample reaction into separate 45- μ l aliquots of XL10-Gold ultracompetent cells (see *Transformation of XL10-Gold Ultracompetent Cells* in the instruction manual).