



QuikChange HT Protein Engineering System

Protocol

Version B.0, June 2015

SureSelect platform manufactured with Agilent SurePrint Technology

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



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2015

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

Manual Part Number

G5900-90000

Edition

Version B.0, June 2015

Printed in USA

Agilent Technologies, Inc.
5301 Stevens Creek Rd
Santa Clara, CA 95051 USA

Technical Support

For technical product support, contact your local Agilent Support Services representative.

For US and Canada, call (800) 227-9770 (option 3,4,4). For other countries, find your support center telephone numbers at www.agilent.com/chem/contactus.

Or send an e-mail to:
SureSelect.Support@agilent.com

Warranty

The material contained in this document is provided “as is,” and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

Notice to Purchaser

Use of this product is covered by one or more of the following U.S. Patent Nos. 5,789,166, 5,932,419, 6,391,548, 6,713,285, 7,132,265, and 7,176,004.

This product is provided under an agreement between Bio-Rad Laboratories and Agilent Technologies, Inc., and the manufacture, use, sale or import of this product is subject to US. Pat. No. 6,627,424 and EP Pat. No. 1 283 875 B1, owned by Bio-Rad Laboratories, Inc. Purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in PCR (but not real-time PCR) in the Research Field including all Applied Research Fields (including but not limited to forensics, animal testing, and food testing).

Safety Notices

CAUTION

A **CAUTION** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a **CAUTION** notice until the indicated conditions are fully understood and met.

WARNING

A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

In this Guide...

This document describes how to use the QuikChange HT Protein Engineering System to construct mutant plasmid libraries.

1 Before You Begin

Make sure you read and understand the information in this chapter and have the necessary equipment and reagents listed before you start an experiment.

2 Procedures

This chapter contains protocols for Agilent's QuikChange HT Protein Engineering System.

3 Reference

This chapter contains reference information related to the protocol.

What's New in Version B.0

- Updated product labeling statement

Content

1 Before You Begin

| | |
|--------------------|----|
| Required Reagents | 10 |
| Required Equipment | 11 |
| Procedural Notes | 12 |

2 Procedures

| | |
|--|----|
| Outline of the Protocol and Positive Control | 14 |
| Overview of the Workflow | 14 |
| Outline of the Protocol | 15 |
| Use of the Positive Control | 16 |
| Step 1. PCR Amplify the Custom Mutagenesis Library | 17 |
| Step 2. Purify the Amplified Mutagenesis Library | 19 |
| Step 3. Construct the Mutant Plasmid Library using Thermal Cycling | 21 |
| Step 4. Design the Library Recovery Strategy using a Pilot Transformation Reaction | 23 |
| Troubleshooting | 26 |

3 Reference

| | |
|-----------------------------------|----|
| Kit Contents | 30 |
| Preparation of Media and Reagents | 32 |



1 Before You Begin

| | |
|--------------------|----|
| Required Reagents | 10 |
| Required Equipment | 11 |
| Procedural Notes | 12 |

Make sure you read and understand the information in this chapter and have the necessary equipment and reagents listed before you start an experiment.



Required Reagents

Table 1 Required Reagents for the QuikChange HT Protein Engineering System protocol

| Description | Vendor and part number |
|--|--|
| QuikChange HT Protein Engineering System Kit* | Agilent |
| 150 nt, 10 sites, Non-academic Users | p/n G5900A |
| 150 nt, 20 sites, Non-academic Users | p/n G5900B |
| 200 nt, 10 sites, Non-academic Users | p/n G5901A |
| 200 nt, 20 sites, Non-academic Users | p/n G5901B |
| 150 nt, 10 sites, Academic Users | p/n G5902A |
| 150 nt, 20 sites, Academic Users | p/n G5902B |
| 200 nt, 10 sites, Academic Users | p/n G5903A |
| 200 nt, 20 sites, Academic Users | p/n G5903B |
| PCR-grade water | Ambion p/n AM9930, or equivalent |
| Ethanol, molecular biology grade | Sigma-Aldrich p/n E7023, or equivalent |
| 5 mM Tris-HCl, pH 8.0, 0.1 mM EDTA | General laboratory supplier |
| NZY ⁺ broth | Prepare as directed on page 32 |
| LB–ampicillin agar plates | Prepare as directed on page 32 |
| isopropyl-1-thio- β -D-galactopyranoside (IPTG) | Agilent p/n 300127 |
| 5-bromo-4-chloro-3-indolyl- β -D-galactopyranoside (X-gal) | Agilent p/n 300201 |

* The custom oligonucleotide libraries and custom primers provided in each kit are designed using the Agilent eArray design system. You must complete the design process in eArray before ordering the kit.

Required Equipment

Table 2 Required Equipment for the QuikChange HT Protein Engineering System protocol

| Description | Vendor and part number |
|--|---|
| Thermal Cycler | Agilent SureCycler 8800, p/n G8800A or equivalent |
| 8-well PCR strip tubes | Agilent, p/n 410092 |
| Tube cap strips (8 domed caps per strip) | Agilent, p/n 410096 |
| P20 and P200 pipettes | Pipetman or equivalent |
| Ice bucket | General laboratory supplier |
| Vortex mixer | General laboratory supplier |
| Microcentrifuge | Eppendorf Centrifuge model 5804 or equivalent |

Procedural Notes

- The plasmid to be mutagenized must be isolated from a dam^+ *E. coli* strain. The majority of commonly-used *E. coli* strains are dam^+ . Plasmid DNA isolated from dam^- strains, such as the JM110 and SCS110 strains, is not a suitable target for mutagenesis using this system.
- Efficiency of the mutagenesis reaction and of mutant plasmid recovery by transformation will vary for each custom library. The protocol in this manual includes a pilot transformation experiment to determine the number of transformation reactions required to reach the target library size. Additional competent cells may be required to recover the targeted number of mutagenized plasmids.
- The provided SoloPack Gold Supercompetent cells are resistant to tetracycline and chloramphenicol. If the plasmid to be mutagenized contains only the tet^r and chl^r resistance markers, an alternative tetracycline-sensitive strain of competent cells must be used.
- The custom Mutagenesis Library is used for both control and protein engineering mutagenesis reactions. The oligonucleotide required to mutate the positive control pWS4.5 plasmid in the positive control reaction is present in each custom oligonucleotide library mixture.
- The Positive Control Primer Mix contains primers with the following sequences:
5'-CCACTAGTTCTAGAGCGGC-3'
5'-ACAGCTATGACCATGATTACG-3'



2 Procedures

| | |
|--|----|
| Outline of the Protocol and Positive Control | 14 |
| Step 1. PCR Amplify the Custom Mutagenesis Library | 17 |
| Step 2. Purify the Amplified Mutagenesis Library | 19 |
| Step 3. Construct the Mutant Plasmid Library using Thermal Cycling | 21 |
| Step 4. Design the Library Recovery Strategy using a Pilot Transformation Reaction | 23 |
| Troubleshooting | 26 |

This chapter contains protocols for Agilent's QuikChange HT Protein Engineering System.



Outline of the Protocol and Positive Control

Overview of the Workflow

The QuikChange HT Protein Engineering System allows you to construct a diverse collection of engineered mutant clones using Agilent's QuikChange technology, using the workflow summarized in [Figure 1](#).

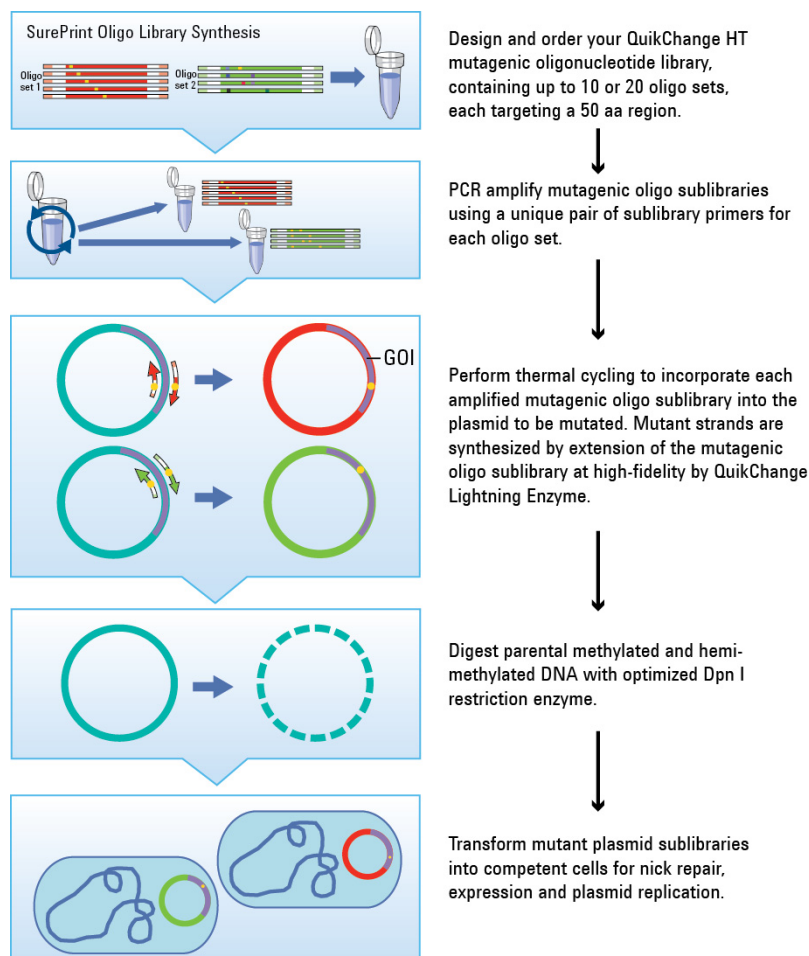


Figure 1 Overall QuikChange HT workflow.

Outline of the Protocol

The QuikChange HT Protein Engineering System protocol uses a library of custom mutagenic oligonucleotides and custom primers to produce an engineered mutant plasmid library. The mutagenic oligonucleotide library and custom primers must be designed using Agilent's eArray website before ordering a custom QuikChange HT Protein Engineering System kit.

The protocol steps for mutant plasmid library construction are summarized below. Detailed instructions are provided starting on [page 17](#).

- In step 1 (see [page 17](#)) the provided custom mutagenic oligonucleotide library is PCR amplified using the appropriate pair of provided custom primers.
- In step 2 (see [page 19](#)) the amplified mutagenesis library is purified using kit-provided components.
- In Step 3 (see [page 21](#)) the mutagenic oligonucleotide library is incorporated into the plasmid to be mutagenized by a linear amplification reaction using a thermocycler. Components of the thermal cycling reaction include the supercoiled, double-stranded DNA template, the amplified oligonucleotide library containing the desired mutations, and the kit-provided enzyme blend featuring a *Pfu* Fusion DNA polymerase, which extends the mutagenic primers with high fidelity. After thermal cycling, the parental DNA template is removed by treating the reaction products with the restriction endonuclease *Dpn* I (target sequence: 5´-Gm6ATC-3´). *Dpn* I acts specifically on methylated and hemimethylated DNA, and does not cleave the newly synthesized mutagenized DNA library. DNA isolated from almost all *E. coli* strains is dam methylated and therefore susceptible to digestion.
- In step 4 (see [page 23](#)) the mutagenized plasmid library is transformed into the provided SoloPack Gold supercompetent cells. The instructions provided are for the initial pilot transformation to determine the appropriate clone recovery strategy for the custom library.

Use of the Positive Control

The kit includes a *LacZ*-based control template and control primers to verify performance of the QuikChange HT Protein Engineering System custom library, reagents, and protocol.

The supplied 4.5 kb pWS4.5 control plasmid contains the *LacZ* coding sequence with a stop codon mutation that prevents expression of full-length *LacZ*. The *LacZ* gene product is responsible for the production of blue colonies when appropriate *E. coli* transformants are grown on media containing X-gal in the presence of the inducer IPTG. The provided SoloPack Gold supercompetent cells transformed with the pWS4.5 control plasmid appear white on LB-ampicillin agar plates containing IPTG and X-gal, because the stop codon in the control plasmid prevents production of active beta-galactosidase.

In addition to the custom-designed oligonucleotides, each QuikChange HT Mutagenesis library also contains mutagenic oligonucleotides that direct reversion of the mutant stop codon in the pWS4.5 control plasmid. After completing the QuikChange HT Protein Engineering System protocol using the provided Positive Control Primer Mix for amplification from the custom library, a mutagenized control plasmid library is constructed. Following transformation of the control library into SoloPack Gold supercompetent cells, colonies are scored for the blue color production phenotype, where a blue colony indicates the production of a mutagenized control plasmid expressing the full length *LacZ* gene product. Each custom mutagenesis library should be used to construct a positive control plasmid library in parallel with the custom mutant plasmid library to verify efficient amplification and transformation of the custom library.

Step 1. PCR Amplify the Custom Mutagenesis Library

- 1 For each Forward Custom Primer and Reverse Custom Primer to be used in the experiment, dilute 1 μL of the supplied primer solution with 19 μL of PCR-grade water, for a final concentration of 5 μM .
- 2 Prepare 50- μL custom library and control PCR amplification reactions by combining the reagents in [Table 3](#). Mix thoroughly by vortexing.

Table 3 Amplification of Mutagenesis Library

| Component | Custom Library Amplification | Positive Control Amplification |
|--|------------------------------|--------------------------------|
| 2 \times PfuUltra II HS Master Mix AD | 25 μL | 25 μL |
| QuikChange HT Mutagenesis Library solution | 1 μL | 1 μL |
| Forward Custom Primer dilution (5 μM , prepared in step 1) | 5 μL | — |
| Reverse Custom Primer dilution (5 μM , prepared in step 1) | 5 μL | — |
| 10 \times Positive Control Primer Mix (5 μM each primer) | — | 5 μL |
| PCR-grade H ₂ O | 14 μL | 19 μL |

- 3 Cycle the amplification reactions according to [Table 4](#) below.

Table 4 Thermal cycler program for Mutagenic Oligo Library amplification

| Segment | Cycles | Temperature | Time |
|---------|--------|-----------------------------|------------|
| 1 | 1 | 95°C | 30 seconds |
| 2 | 30 | 95°C | 20 seconds |
| | | primer T _m –5°C* | 10 seconds |
| | | 72°C | 30 seconds |
| 3 | 1 | 72°C | 2 minutes |
| 4 | 1 | 4°C | Hold |

* For custom library amplification, obtain custom primer T_m values from the supplied Custom Primer tube labels. For the positive control reaction, use an annealing temperature of 55°C.

2 Procedures

Step 1. PCR Amplify the Custom Mutagenesis Library

- 4 Analyze 5- μ L samples of the custom library amplicons and of the positive control amplicon by electrophoresis on a 2% (w/v) agarose gel. Load 5 μ L of the provided 1.1 kb Gel Standard marker DNA solution in an adjacent lane.
- 5 After staining the gel with a suitable DNA-staining agent, the intensity of each amplicon band should be comparable to the intensity of the band in the 5- μ L 1.1 kb Gel Standard lane.

If the library amplicon band (from 5 μ L sample) is substantially less intense than the 1.1 kb Gel Standard band (from 5 μ L of the standard solution), troubleshoot the amplification reaction steps before proceeding to purification and plasmid library construction steps.

Step 2. Purify the Amplified Mutagenesis Library

- 1 Before starting the purification procedure with a new set of reagents, prepare 1× buffer solutions according the instructions in [Table 5](#).

After adding the ethanol, be sure to mark the ethanol-added checkbox on the label for reference by later users.

Table 5 Preparation of 1× Buffer Solutions

| 1× Buffer to be Prepared | Preparation Instructions |
|--------------------------|--|
| 1× DNA Binding Buffer | Add 2.5 mL of 80% ethanol to the provided DNA Binding Buffer bottle (shipped containing 2.5 mL of 2× buffer) |
| 1× PCR Wash Buffer | Add 20 mL of 100% ethanol to the vial of 5× PCR Wash Buffer (shipped containing 5.0 mL of 5× buffer) |

- 2 Purify the amplicons from each library amplification reaction using the provided microspin cups and receptacle tubes according to the following protocol.
 - a To the remaining 45 μ L of amplification reaction, add 100 μ L of 1× DNA Binding Buffer and mix well by vortexing.
 - b For each library amplification reaction, seat a microspin cup (provided) in a receptacle tube (provided).
 - c Transfer each amplicon plus binding buffer mixture (approximately 145 μ L) into a seated microspin cup. Avoid touching the fiber matrix of the microspin cup with the pipet tip.
 - d Cap the spin cup with the provided receptacle tube cap, then spin in a microcentrifuge for 30 seconds at maximum speed (13,000 to 14,000 rpm).
 - e Retain the microspin cup, with bound DNA, and discard the flow-through solution in the receptacle tube (retaining the receptacle tube).
 - f Re-seat the spin cup in the receptacle tube, and add 500 μ L of 1× Wash Buffer.
 - g Cap the spin cup, then spin in a microcentrifuge for 30 seconds at maximum speed.

2 Procedures

Step 2. Purify the Amplified Mutagenesis Library

- h** Retain the microspin cup, and discard the flow-through solution in the receptacle tube (retaining the receptacle tube).
 - i** Re-seat and cap the spin cup, then spin in a microcentrifuge for an additional 2 minutes at maximum speed.
 - j** Transfer the microspin cup to a fresh 1.5-mL microcentrifuge tube, then add 40 μ L of TE elution buffer (5 mM Tris-HCl, 0.1 mM EDTA, pH 8.0) to the spin cup.
 - k** Incubate the spin cup at room temperature for 1 minute.
 - l** Cap the spin cup with the tube cap, then spin in a microcentrifuge for 1 minute at maximum speed.
 - m** Retain the 40- μ L purified amplicon solution in the microcentrifuge tube, and discard the microspin cup.
- 3** Keep the amplified library solutions on ice until use on [page 21](#).

Stopping Point

If you do not continue immediately to the next step of the protocol, store the samples at -20°C until use.

Step 3. Construct the Mutant Plasmid Library using Thermal Cycling

Prepare the custom mutant plasmid library using the linear amplification mutant strand synthesis protocol detailed below. In this step, the plasmid to be mutagenized serves as template and the mutagenesis sublibrary amplicons from [page 20](#) serve as primer in the mutant strand synthesis reaction.

For the positive control reaction, use the provided Positive Control Plasmid pWS4.5 as template and the control mutagenesis library amplicons as primer.

After amplification, the products are digested with *Dpn* I restriction enzyme to remove the parental (non-mutated) supercoiled dsDNA.

- 1 Prepare the 25- μ L mutant plasmid library synthesis reactions by combining the reagents listed in [Table 6](#).

Table 6 Mutant plasmid library synthesis

| Component | Custom Library Preparation | Positive Control |
|---|----------------------------|------------------|
| 10 \times QuikChange Lightning Buffer | 2.5 μ L | 2.5 μ L |
| dNTP Mix | 1 μ L | 1 μ L |
| QuikSolution | 0.75 μ L | 0.75 μ L |
| Mutagenesis target plasmid (25 ng/ μ L) | 1 μ L | — |
| Positive Control Plasmid pWS4.5 (25 ng/ μ L) | — | 1 μ L |
| Purified custom mutagenesis library amplicon (from step m on page 20) | 15 μ L | — |
| Purified control mutagenesis library amplicon (from step m on page 20) | — | 5 μ L |
| QuikChange Lightning Enzyme | 1 μ L | 1 μ L |
| PCR-grade H ₂ O | 3.75 μ L | 13.75 μ L |

2 Procedures

Step 3. Construct the Mutant Plasmid Library using Thermal Cycling

- 2 Cycle the reactions using the cycling parameters in [Table 7](#) below.

Table 7 Thermal cycler program for mutant plasmid library synthesis

| Segment | Cycles | Temperature | Time |
|---------|--------|-------------|--------------------------------------|
| 1 | 1 | 95°C | 2 minutes |
| 2 | 18 | 95°C | 20 seconds |
| | | 60°C | 10 seconds |
| | | 68°C | 30 seconds/kb target plasmid length* |
| 3 | 1 | 68°C | 5 minutes |

* For the pWS4.5 positive control reaction, use a 2.5-minute extension time.

- 3 Add 1 μL of the provided *Dpn* I Enzyme directly to each amplification reaction.
- 4 Gently and thoroughly mix each reaction mixture by pipetting the solution up and down several times. Briefly centrifuge the reaction tube and then immediately incubate at 37°C for 5 minutes.

Dpn I-digestion is required to remove the non-mutated dsDNA supplied as template in the amplification reaction.

NOTE

The protocol detailed on the following pages is designed to determine the number of transformants obtained from 1.5 μL of the mutagenesis library. Store the remaining library solution at -20°C for additional rounds of transformation following the pilot transformation experiment.

Step 4. Design the Library Recovery Strategy using a Pilot Transformation Reaction

Efficiency of the mutagenesis reaction and of mutant plasmid recovery by transformation may vary for each custom library. The purpose of the pilot transformation is to determine the number of transformation reactions required to reach the target library size and to determine the optimal plating strategy for the transformation reactions.

If desired, the transformation efficiency of the competent cells may be verified by transforming a tube of cells with 0.01 ng of pUC18 DNA. (Dilute the provided pUC18 DNA prior to use.) Using the conditions specified below, the expected efficiency is $\geq 1 \times 10^9$ cfu/ μ g pUC18 DNA.

1 Preheat NZY⁺ broth to 42°C for use as the outgrowth medium in [step 8](#).

NOTE

The optimized transformation protocol uses NZY⁺ broth for cell outgrowth (see [page 32](#) for a recipe). SOC medium may also be used for cell outgrowth, but transformation efficiency will be reduced.

- 2** For each custom sublibrary and positive control sample to be transformed, thaw one tube of SoloPack Gold supercompetent cells on ice.
- 3** When the cells have thawed, swirl the tube gently to mix the cells.
- 4** Transfer 1.5 μ L of the *Dpn* I-treated DNA from each mutant sublibrary construction reaction and control reaction to a separate tube of cells.
- 5** Swirl the transformation reaction gently to mix and then incubate the tube on ice for 30 minutes.
- 6** After removing any ice from the outside of the tube, heat-pulse the tube in a 42°C water bath for 30 seconds. The temperature and duration of the heat pulse are critical for maximum efficiency.
- 7** Incubate the tube on ice for 2 minutes.
- 8** Add 250 μ L of preheated (42°C) NZY⁺ broth and incubate the tube at 37°C for 1 hour with shaking at 225-250 rpm.

2 Procedures

Step 4. Design the Library Recovery Strategy using a Pilot Transformation Reaction

- 9 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. Prior to plating the cells, pipet 100 μL of NZY⁺ broth onto the surface of the agar. Pipet the transformation mixture into the pool of liquid medium and then spread the mixture with a sterile spreader.

Table 8 Transformation reaction plating volumes and media

| Reaction Type | Volume to Plate | Plating Medium |
|-------------------------------------|---|--|
| pWS4.5 positive control mutagenesis | 2 μL and 10 μL | LB-ampicillin agar containing 20 mM IPTG and 80 $\mu\text{g}/\text{mL}$ X-gal |
| Custom mutant plasmid library | 10 μL and 50 μL on each of two plates | LB agar containing the appropriate antibiotic (and mutagenesis indicators, where applicable) |

- 10 Incubate the plates at 37°C overnight. For the pWS4.5 positive control plates and any other plates used for blue-white color screening, incubate the plates at 37°C for at least 16 hours to allow full color development. Blue color may also be enhanced by subsequent incubation of the plates at 4°C for 2 hours.

- 11 Evaluate the transformation results and plan the optimal large-scale library transformation plating strategy.

- For the pWS4.5 positive control mutagenesis library, $\geq 80\%$ of the total colony forming units (cfu) should be blue-colored, indicating a mutagenesis efficiency $\geq 80\%$.
- For the custom mutagenesis library, colony number and mutagenesis efficiency will vary, depending on the plasmid size and sequence, library design and sublibrary primer design.

Determine the number of transformation reactions required to reach the target library size (target number of mutagenized plasmid clones) and to determine the optimal plating strategy for each transformation reaction using the steps below:

- a Count and record the number of cfu on each of the custom mutagenesis library transformation plates.

Step 4. Design the Library Recovery Strategy using a Pilot Transformation Reaction

- b** Determine the total number of cfu expected from transformation of each 1.5- μ L sample of the *Dpn* I-digested mutagenesis reaction. A sample calculation is shown below.

Table 9 Sample calculation of total cfu expected per transformation

| Volume of transformation reaction plated | Average cfu obtained | Average cfu obtained per μ l of transformation reaction | Calculated cfu per 300- μ l transformation reaction | Average calculated cfu per 300- μ l transformation reaction containing 1.5 μ l of mutagenesis library |
|--|----------------------|---|---|---|
| 10 μ L | 50 | 5 | 1500 | 1350 |
| 50 μ L | 200 | 4 | 1200 | |

- c** Determine the number of library transformation reactions required to generate the target library size. Using the example plating data in [Table 9](#), if your target library size is 1×10^4 clones, you would need to complete eight transformation reactions.
- d** Determine the optimal plating volume for your custom mutagenesis library. Typically, the optimal plating volume will yield approximately 200 well-isolated cfu per 100-mm plate. Using the example plating data in [Table 9](#), plating the full-scale transformation reactions at 50 μ L per plate is expected to produce a suitable cfu density for clone recovery.

Once the appropriate plating strategy is determined for your custom library, transform additional 1.5- μ L aliquots of each *Dpn* I-treated mutagenized plasmid sublibrary into additional tubes of SoloPack Gold supercompetent cells, using the transformation protocol starting on [page 23](#). Additional tubes of SoloPack Gold supercompetent cells may be purchased separately (Agilent p/n 230350).

Troubleshooting

| Observation | Suggestions |
|---|--|
| No product or low yield of mutagenesis library amplicons (from step 5 on page 18) | <p>Amplification yield may be improved for some custom library and custom primer pair combinations by substitution of the 2× PfuUltra II HS Master Mix AD with QuikChange Lightning components. Prepare the amplification reaction using the following reagents:</p> <ul style="list-style-type: none">32 μL PCR-grade water5 μL 10× QuikChange Lightning Buffer1 μL QuikChange HT Mutagenesis Library solution5 μL Forward Custom Primer dilution (5 μM)5 μL Reverse Custom Primer dilution (5 μM)1 μL dNTP Mix1 μL QuikChange Lightning Enzyme <p>Use the cycling conditions described in Table 4 on page 17 and continue with the protocol according to the instructions provided.</p> <hr/> <p>Titrate the concentration of the Forward and Reverse Custom Primers between 0.5 μM and 2.5 μM in the final reaction.</p> <hr/> <p>For library designs with high GC content or secondary structures, it may be beneficial to include a cosolvent in the PCR reaction. Titrate addition of QuikSolution in the reaction at 1% to 10% (v/v) final concentration or glycerol at 5% to 20% (v/v) final concentration.</p> <hr/> <p>Increase the cycle number to greater than 30.</p> |
| Multiple bands produced in the mutagenesis library amplification reaction (from step 5 on page 18) | <p>Ensure that the primer annealing temperature is optimal. Increase the annealing temperature in 5°C increments or try an annealing temperature gradient from 55°C to 70°C.</p> |

| Observation | Suggestions |
|--|--|
| <p>Low colony number or low mutagenesis efficiency after transformation of the LacZ control mutant plasmid library synthesis reaction (results from step 11 on page 24)</p> | <p>Verify the transformation efficiency of the SoloPack Gold Supercompetent Cells by transforming a tube of cells with 0.01 ng of the provided pUC18 DNA. (Dilute the provided pUC18 DNA prior to use.) Using the provided transformation protocol, the expected efficiency is $\geq 1 \times 10^9$ cfu/μg pUC18 DNA. Ensure that supercompetent cells are stored at the bottom of a -80°C freezer immediately upon receipt.</p> <hr/> <p>Verify that the agar plates were prepared correctly. See Preparation of Media and Reagents on page 33 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 and stored at 4°C overnight.</p> <hr/> <p>Increase amount of control mutagenesis library amplicon added in Table 6 on page 21 from 5 μL to 15 μL.</p> <hr/> <p>Allow sufficient time for the <i>Dpn</i> I enzyme to completely digest the parental template; increase the digestion time to 30 minutes.</p> <hr/> <p>Avoid multiple freeze-thaw cycles for the dNTP mix. If necessary, prepare single-use aliquots of the provided dNTP mix, and store the aliquots at -20°C.</p> |
| <p>Low colony number or low mutagenesis efficiency after transformation of the custom mutant plasmid library synthesis reaction (results from step 11 on page 24 to page 25)</p> | <p>Analyze the plasmid DNA used in the reaction by gel electrophoresis to verify the quantity and quality of the template. Nicked or linearized plasmid DNA will not generate complete circular products; verify that the template DNA is at least 80% supercoiled.</p> <hr/> <p>Repeat the strand synthesis reaction using 50 to 100 ng plasmid DNA.</p> <hr/> <p>Titrate addition of QuikSolution in the strand synthesis reaction at 0.5 μL increments from 0 to 2.5 μL.</p> <hr/> <p>Increase the amount of the mutagenesis library amplicon added to the strand synthesis reaction by eluting the amplicon in 20 μL of elution buffer (step j on page 20). Alternatively, concentrate the 40-μL eluate before use using a method of choice.</p> <hr/> <p>Increase extension time in the strand synthesis reaction from 30 seconds/kb to 45 second/kb.</p> <hr/> <p>Increase the amount of <i>Dpn</i> I-treated DNA used in the transformation reaction to 4 μL. If needed, you can purify and precipitate DNA from the entire strand synthesis reaction. After resuspending the DNA in a small volume, transform competent cells using the entire volume of DNA.</p> <hr/> <p>Allow sufficient time for the <i>Dpn</i> I enzyme to completely digest the parental template. Increase the digestion time to 30 minutes or repeat the <i>Dpn</i> I digestion if too much DNA was present.</p> <hr/> <p>Avoid multiple freeze-thaw cycles for the dNTP mix. If necessary, prepare single-use aliquots of the provided dNTP mix, and store the aliquots at -20°C.</p> |

2 **Procedures**

Troubleshooting



3 Reference

Kit Contents 30
Preparation of Media and Reagents 32

This chapter contains reference information related to the protocol.



Kit Contents

QuikChange HT Protein Engineering System Kits contain the following component kits:

Table 10 QuikChange HT Protein Engineering System Kit Contents

| Component Kits | Storage Condition | 10 Sites | 20 Sites |
|--|--------------------|-----------|------------|
| QuikChange HT Mutagenesis Library and Primers [*] | -20°C | 5190-7454 | 5190-7454 |
| QuikChange HT Mutagenesis Reagents | -20°C [†] | 5190-7452 | 5190-7455 |
| QuikChange HT DNA Cleanup Kit | Room Temperature | 5190-7456 | 5190-7456 |
| SoloPack Gold Supercompetent Cells | -80°C | 230350 | 230350 × 2 |

* Kits may also be provided with the QuikChange HT Mutagenesis Library and the Sublibrary Primers in separate boxes (p/n 5190-7451 and 5190-7454, respectively).

† Store at -20°C upon receipt. After thawing the 2× PfuUltra II HS Master Mix AD, store the master mix at 4°C; do not subject to multiple freeze-thaw cycles. Store the remainder of the QuikChange HT Mutagenesis Reagents kits at -20°C throughout use of the product.

The contents of the multi-component kits listed in [Table 10](#) are described in [Table 11](#) through [Table 14](#) below.

Table 11 QuikChange HT Mutagenesis Library and Primers Content

| Reagent | Format |
|--|---|
| QuikChange HT Mutagenesis Library | single tube with clear cap |
| QuikChange HT Mutagenesis Sublibrary Primers | variable number of primer-containing tubes (one pair of primer tubes per sublibrary in custom design) |

Table 12 QuikChange HT Mutagenesis Reagents Content

| Reagent | Format |
|--|----------------------|
| 2× PfuUltra II HS Master Mix AD | tube with clear cap |
| 10× QuikChange Lightning Buffer | tube with clear cap |
| QuikChange Lightning Enzyme | tube with red cap |
| 40 mM dNTP | tube with yellow cap |
| QuikChange Solution | tube with green cap |
| Positive Control Plasmid pWS4.5 (25 ng/μL) | tube with green cap |
| 10× Positive Control Primer Mix | tube with green cap |
| 1.1 kb Gel Standard | tube with blue cap |
| Dpn I Enzyme | tube with black cap |

Table 13 QuikChange HT DNA Cleanup Kit Content

| Kit Component | Format |
|-------------------------------------|--|
| DNA Binding Solution | bottle (contains 2× solution; see page 19 for dilution instructions) |
| PCR Wash Buffer | bottle (contains 5× solution; see page 19 for dilution instructions) |
| Microspin cups and receptacle tubes | bag containing 24 cups and tubes |

Table 14 SoloPack Gold Supercompetent Cells Kit Content

| Reagent | Format |
|--|---|
| SoloPack Gold Supercompetent Cells | 15 single-use transformation tubes (yellow caps) [*] |
| pUC18 control plasmid (0.1 ng/μl in TE buffer) | tube with blue cap [*] |

^{*} Two sets of 15 transformation tubes and two tubes of pUC18 are provided with 20 Site QuikChange HT Protein Engineering System Kits.

Preparation of Media and Reagents

NZY+ Broth (per Liter)

10 g NZ amine (casein hydrolysate)

5 g yeast extract

5 g NaCl

Add deionized H₂O to a final volume of 1 liter

Adjust pH to 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

LB Agar (per Liter)

10 g tryptone

5 g yeast extract

10 g NaCl

20 g 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 (approximately 25 mL per 100-mm plate)

LB-Ampicillin Agar (per Liter)

Prepare 1 liter of LB agar

Autoclave

Cool to 55°C

Add 10 ml of 10 mg/mL filter-sterilized ampicillin

Pour into petri dishes (approximately 25 mL per 100-mm plate)

LB-Ampicillin Agar with IPTG and X-gal for blue-white screening (per Liter)

Prepare 1 liter of LB agar

Autoclave

Cool to 55°C, then add the following components:

10 ml of 10 mg/mL filter-sterilized ampicillin

X-gal to a final concentration of 80µg/mL

IPTG to a final concentration of 20 mM

Pour into petri dishes (approximately 25 mL per 100-mm plate)

NOTE

Alternatively, IPTG and X-gal may be spread on pre-poured agar plates using the guidelines provided below.

10 mM IPTG (per 10 mL) for spreading on pre-poured agar plates

24 mg isopropyl-1-thio-β-d-galactopyranoside (IPTG)

10 mL sterile dH₂O

Store at -20°C

Spread 100 µL per 100-mm LB-agar plate for blue-white color screening
(do not mix the IPTG and X-gal solutions prior to spreading)

2% X-gal (per 10 mL) for spreading on pre-poured agar plates

0.2 g 5-bromo-4-chloro-3-indolyl-β-D-galactopyranoside (X-gal)

10 mL dimethylformamide (DMF)

Store at -20°C

Spread 100 µL per 100-mm LB-agar plate for blue-white color screening
(do not mix the X-gal and IPTG solutions prior to spreading)

www.agilent.com

In This Book

This guide contains information to run the QuikChange HT Protein Engineering System protocol.

© Agilent Technologies, Inc. 2015

Version B.0, June 2015



G5900-90000



Agilent Technologies