

Protocol for Q5[®] Site-Directed Mutagenesis Kit (NEB #E0554)

Materials Required but not Supplied

Q5[®] Hot Start High-Fidelity 2X Master Mix

- Template DNA and associated forward and reverse primers
- Nuclease-free Water (NEB #B1500)

KLD Enzyme Mix

- Nuclease-free Water (NEB #B1500)

Overview

Site-directed mutagenesis (SDM) is a method to confer a desired mutation in a double-stranded DNA plasmid using custom-designed oligonucleotide primers. This protocol for the Q5[®] Site-Directed Mutagenesis Kit can be completed in less than two hours for moderately sized plasmids. We recommend using the NEB online primer design software [NEBaseChanger[®]](#).

Protocol

Step I: Exponential Amplification (PCR)

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

	25 µl RXN	FINAL CONC.
Q5 Hot Start High-Fidelity 2X Master Mix	12.5 µl	1X
10 µM Forward Primer	1.25 µl	0.5 µM
10 µM Reverse Primer	1.25 µl	0.5 µM
Template DNA (1–25 ng/µl)	1 µl	1-25 ng
Nuclease-free water	9.0 µl	

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

STEP	TEMP	TIME
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. For pre-designed, back-to-back primer sets, a $T_a = T_m + 3$ rule can be applied, but optimization may be necessary. When using pre-designed primer sets, the T_m should be calculated based on the homologous binding region of the primer, not including any 5' non-homologous tails.

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

1. Assemble the following reagents:

	VOLUME	FINAL CONC.
PCR Product	1 μ l	
2X KLD Reaction Buffer	5 μ l	1X
10X KLD Enzyme Mix	1 μ l	1X
Nuclease-free Water	3 μ l	

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

Step III: Transformation

1. Thaw a tube of NEB 5-alpha Competent *E. coli* (High Efficiency) cells on ice. Add 5 μ 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.
2. Place the mixture on ice for 30 minutes.
3. Heat shock at 42°C for 30 seconds.
4. Place on ice for 5 minutes.
5. Pipette 950 μ l of room temperature SOC Outgrowth Medium into the mixture.
6. Incubate at 37°C for 60 minutes with shaking (250 rpm).
7. Mix the cells thoroughly by flicking the tube and inverting, then spread 50-100 μ 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

Resources:

- [NEBaseChanger](#)