

TOPO® TA Cloning® Kits

Catalog Numbers K4500-01SC, K4550-01SC, K4600-01SC

Doc. Part no. 25-0647

Pub. no. MAN0005242

Rev. A.0

Introduction

Follow these instructions to use the TOPO® TA Cloning® Kit to clone your PCR product into the TOPO® vector and transform the reaction into chemically competent *E. coli* cells. For transformation of electrocompetent cells, a map of the TOPO® vector, a diagram of the multiple cloning site, and additional detailed instructions, refer to the TOPO® TA Cloning® Kit manual available from www.lifetechnologies.com/manuals.

Produce PCR products

Produce PCR products using *Taq* polymerase and your own protocol. End the PCR reaction with a final 7–30 minutes extension step.

TOPO® kit cloning reaction

1. Set up the following 6 µL reaction:

Reagent	Amount*
Fresh PCR Product	0.5–4 µL
Salt Solution	1 µL
Sterile Water	add to a total volume of 5 µL
TOPO® Vector	1 µL
Final Volume	6 µL

- * For transformation of chemically competent *E. coli* only.
2. Mix gently and incubate for 5 minutes at room temperature.
 3. Place tubes on ice. Proceed to
Transformation and Analysis.

Transformation and Analysis

Follow the protocol in this section to transform chemically competent cells and to analyze positive clones. Use the **Rapid One Shot® chemical transformation** protocol if you are using *ampicillin selection only* and if you wish to obtain transformants as quickly as possible. For instructions to transform electrocompetent cells, refer to the TOPO® TA Cloning® manual.

One Shot® chemical transformation

1. Thaw 1 vial of One Shot® *E. coli* cells *on ice* for each transformation.
2. Add 2 µL of the TOPO® cloning reaction to each vial of One Shot® cells to be transformed, and mix gently.
3. Incubate the vials on ice for 5–30 minutes.
4. Heat-shock the cells for 30 seconds at 42°C without shaking.
5. Add 250 µL of room temperature S.O.C. medium to the cells.
6. Cap the tubes and shake them at 37°C for 1 hour.
7. Spread 10–50 µL from each transformation on pre-warmed LB plates containing X-gal and 50–100 µg/mL ampicillin or 50 µg/mL kanamycin. For TOP10F' *E. coli*, add both IPTG and X-gal to LB plates.
8. Incubate plates overnight at 37°C.
9. An efficient TOPO® cloning reaction should produce several hundred colonies. Pick ~10 white or light blue colonies for analysis. Do not pick dark blue colonies. Proceed to **Analyze positive clones**.

Transformation and Analysis, Continued

Rapid One Shot® chemical transformation (for ampicillin selection only)

1. Thaw 1 vial of One Shot® *E. coli* cells *on ice* for each transformation.
2. Add 4 µL of the TOPO® Cloning reaction to each vial of One Shot® cells to be transformed, and mix gently.
3. Incubate the vials on ice for 5 minutes.
4. Spread 50 µL from each transformation on pre-warmed LB plates containing X-gal and 50–100 µg/mL ampicillin. For TOP10F' *E. coli*, add both IPTG and X-gal to LB plates.
5. Incubate plates overnight at 37°C.
6. An efficient TOPO® cloning reaction should produce several hundred colonies. Pick ~10 white or light blue colonies for analysis. Do not pick dark blue colonies. Proceed to **Analyze positive clones**.

Analyze positive clones

1. Culture the 10 colonies overnight in LB medium containing 50–100 µg/mL ampicillin or 50 µg/mL kanamycin.
2. Isolate plasmid DNA using your method of choice. For ultra-pure plasmid DNA, we recommend the PureLink® HQ Mini Plasmid Purification Kit (Cat. no. K2100-01).
3. Analyze the plasmid by restriction analysis or by sequencing to confirm the presence and orientation of the insert.

Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

DISCLAIMER

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

Life Technologies is a Thermo Fisher Scientific brand. © 2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries.

Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA

Phone +1 760 603 7200 | Toll Free in USA 800 955 6288

For support visit lifetechnologies.com/support or
email techsupport@lifetech.com

lifetechnologies.com

25 February 2014

