

# RapidTrans™

## 96-Tube Chemically Competent *E. coli*

(version C2)

Catalog Nos. 11096, 11596

### Active Motif North America

1914 Palomar Oaks, Suite 150  
Carlsbad, California 92008, USA

Toll free: 877 222 9543

Telephone: 760 431 1263

Fax: 760 431 1351

### Active Motif Europe

1104 Avenue Franklin Roosevelt

B-1330 Rixensart, Belgium

UK Free Phone: 0800 169 31 47

France Free Phone: 0800 90 99 79

Germany Free Phone: 0800 181 99 10

Telephone: +32 (0)2 653 0001

Fax: +32 (0)2 653 0050

### Active Motif Japan

Azuma Bldg, 7th Floor

2-21 Ageba-Cho, Shinjuku-Ku

Tokyo, 162-0824, Japan

Telephone: +81 3 5225 3638

Fax: +81 3 5261 8733

Information in this manual is subject to change without notice and does not constitute a commitment on the part of Active Motif, Inc. It is supplied on an “as is” basis without any warranty of any kind, either explicit or implied. Information may be changed or updated in this manual at any time.

This documentation may not be copied, transferred, reproduced, disclosed, or duplicated, in whole or in part, without the prior written consent of Active Motif, Inc. This documentation is proprietary information and protected by the copyright laws of the United States and international treaties.

The manufacturer of this documentation is Active Motif, Inc.

© 2007 Active Motif, Inc., 1914 Palomar Oaks Way, Suite 150; Carlsbad, CA 92008. All rights reserved.

All trademarks, trade names, service marks, or logos referenced herein belong to their respective companies.

<b>TABLE OF CONTENTS</b>	<b>Page</b>
<b>Introduction</b> .....	1
<b>Kit Contents</b> .....	1
Additional Materials Required .....	1
<b>Protocol</b>	
Preparation .....	2
General Transformation Procedure .....	2
Control Transformation .....	3
Calculation of Efficiency .....	3
<b>Appendix</b>	
Section 1: Strain Information .....	4
Section 2: Media Information .....	5
Section 3: References .....	6
Section 4: Technical Services .....	7

## Introduction

---

RapidTrans™ is a 96-tube tray of competent *E. coli* giving the flexibility of using 1 to 96 tubes at your convenience. This eliminates the waste associated with freeze/thawing and aliquoting the competent cells. Each tube contains 50 µl of competent *E. coli*, which is sufficient for one reaction.

## Kit Contents

---

Component	Quantity	Store at
Chemically competent <i>E. coli</i>	96 rxns (50 µl per tube)	-80°C
pUC19 plasmid DNA	100 ng (10 ng/µl)	-20°C
SOC medium	30 ml	Room Temperature

- Important:**
- To maintain high competency level, please store cells at -80°C.
  - For strain genotypes and media recipes please see Appendix.

## Additional Materials Required

- Ice bucket with crushed ice
- 37°C incubators (shaking and non-shaking)
- 42°C water bath
- LB agar plates containing appropriate additives (antibiotics, IPTG, X-gal, etc.)

## Protocol

---

### Preparation

1. Ensure water bath is at 42°C.
2. Pre-warm LB agar plates to 37°C.
3. The box that holds the 96 transformation tubes can be used as a rack for all stages of the procedure. For heat-shock at 42°C, remove the bottom panel prior to immersion in the water bath.

### General Transformation Procedure

1. Remove required number of transformation reaction tubes from –80°C storage and place on ice to thaw.  
**Note:** Individual tubes can be removed by cutting cap strips with a sharp blade.
2. Add 1-5 µl of plasmid DNA/ligation to thawed cells. Mix by tapping tubes gently and replace on ice immediately.  
**Do not** mix by vortexing or pipetting.  
**Do not** add more than 5 µl (10% of competent cell volume) to reaction.
3. Incubate transformation reactions on ice for 30 minutes.
4. Heat-shock the tubes by immersing in a 42°C water bath for exactly 30 seconds.  
**Note:** The bottom of the 96-tube storage rack can be removed to allow for easier immersion of the tubes into the water bath.
5. Replace transformation reactions on ice for 2 minutes.
6. Aseptically add 250 µl SOC medium to each reaction.  
**Note:** A special reservoir is supplied with the kit to facilitate ease of addition of SOC medium when carrying out multiple transformations. Aseptically transfer SOC medium into the sterile reservoir and use a multi-channel pipette to add medium to transformation reactions.
7. Incubate tubes at 37°C for 1 hour with shaking at 225-250 rpm.
8. Using a sterile spreader, plate out 20-200 µl of each transformation on pre-warmed LB agar plates.
9. Allow plates to completely absorb any excess media.
10. Incubate inverted plates overnight at 37°C.

## Control Transformation

It is recommended that you test the transformation efficiency of the cells supplied in this kit. An aliquot of pUC19 supercoiled plasmid DNA is included with the kit for this purpose. The control procedure is similar to the general transformation procedure described above, with the following additional steps.

1. Prepare a 10 pg/μl solution of pUC19 in sterile H<sub>2</sub>O from the supplied 10 ng/μl stock solution by serial dilution. Add 1 μl of the 10ng/μl stock to 99 μl of sterile H<sub>2</sub>O to make a 100 pg/μl solution.
2. Add 1 μl of the 100 pg/μl solution to 9 μl of sterile H<sub>2</sub>O to make a 10 pg/μl solution. Store on ice until ready to use.
3. Transform 1 μl of the 10 pg/μl solution according to the general transformation procedure described above.
4. Plate 10 & 100 μl of the transformation reaction on two separate LB Amp plates (for easier plating add 90 μl SOC to the 10 μl transformation reaction).
5. Incubate overnight at 37°C.

## Calculation of Efficiency

$$\begin{array}{ccccccc} \text{No. of colonies} & \times & 10^6 \text{pg} & \times & \text{total transformation vol (300 } \mu\text{l)} & = & \text{No. of transformants} \\ 10 \text{ pg transformed pUC19} & & \mu\text{g} & & \text{transformation vol plated} & & \mu\text{g pUC19} \end{array}$$

## Appendix

---

### Section 1: Strain Information

#### For use in cloning, plasmid preparation and library construction

TAMI: *mcrA*  $\Delta$ (*mrr-hsdRMS-mcrBC*)  $\Phi$ 80*lacZ* $\Delta$ M15  $\Delta$ *lacX74* *recA1* *ara* $\Delta$ 139 (*ara-leu*)7697  
*galU galK rpsL endA1 nupG*

1. The *mcrA*, *mcrBC* and *mrr* mutations prevent cleavage of methylated DNA. The McrA and McrBC restriction systems<sup>1-4</sup> cleave DNA at methylated cytosines contained in the target sequences 5'-CG-3' for McrA and 5'-PuC-3' for McrBC. DNA from some sources (including human DNA) may be methylated at these sites and will therefore be cloned inefficiently in strains expressing McrA and/or McrBC. The Mrr restriction system<sup>1, 3, 5</sup> cleaves DNA at methylated adenines, although the precise recognition sequence is not known. DNA from some sources may be methylated at Mrr recognition sites and will therefore be cloned inefficiently in a *mrr+* strain.
2. The EcoK system (*hsdM* DNA methylase, *hsdR* endonuclease and *hsdS* specificity determinant) recognizes the sequence 5'-AACNNNNNGTT-3'<sup>1, 6, 7</sup>. The modification component protects the host DNA by methylation of the second A in each strand of the target sequence. DNA cloned in a *hsdM* host will be restricted if subsequently transferred into a *hsdR+* host.
3. The *lacZ* $\Delta$ M15 deletion removes the amino-terminal  $\alpha$  peptide (amino acids 11-41) of  $\beta$ -galactosidase. Cloning vectors that employ Lac selection carry a gene that codes for the  $\alpha$  peptide and rescues the *lacZ* $\Delta$ M15 mutation by  $\alpha$  complementation<sup>8</sup>.
4. RecA is responsible for general recombination, DNA repair and phage  $\lambda$  induction. Mutation in *recA* helps stabilize sequences with direct repeats.
5. EndA is a DNA-specific endonuclease I. Mutation in *endA* results in improved quality of plasmid miniprep DNA.

## Section 2: Media Information

### LB Agar (per liter)

10 g NaCl

10 g tryptone

5 g yeast extract

Dissolve in 950 ml of deionized H<sub>2</sub>O

Adjust pH to 7.0 with 5 N NaOH

15 g agar

Add deionized H<sub>2</sub>O to 1 L final volume

Autoclave and let cool to ~55°C before pouring or addition of antibiotics, etc.

### SOB Medium (per liter)

20 g tryptone

5 g yeast extract

0.5 g NaCl

Dissolve in 950 ml of deionized H<sub>2</sub>O

Add 10 ml of 250 mM KCl stock solution (1.86 g KCl/100 ml deionized H<sub>2</sub>O)

Adjust pH to 7.0 with 5 N NaOH

Add deionized H<sub>2</sub>O to 1 L final volume

Autoclave and let cool to ~55°C

Aseptically add 10 ml of sterile 1 M MgCl<sub>2</sub>.

### SOC Medium (per liter)

Aseptically add 7.2 ml of sterile 50% glucose to SOB medium.

### Blue/White Screening

#### X-gal Stock Solution

Make a 40 mg/ml solution of X-gal in dimethylformamide (DMF). Use a glass or polypropylene tube and wrap in aluminum foil to prevent damage by light. Store at -20°C.

Plate 40 µl of X-gal stock solution on top of the LB agar and let stand for approximately 1 hour. Alternatively, add the X-gal to the molten LB agar prior to pouring plates.



### Section 3: References

1. Raleigh, E.A., Leach, K. and Brent, R. (1989) in *Current Protocols in Molecular Biology* (F.M. Ausubel, R. Brent, R.E. Kingston, D.D. Moore, J.G. Seidman, J.A. Smith and K. Struhl, eds). John Wiley, New York, p. 1.4.6-1.4.10.
2. Raleigh, E.A. (1992) *Mol. Microbiol.*, 6, 1079.
3. Kelleher, J. and Raleigh, E.A. (1991) *J. Bacteriol.*, 173, 5220.
4. Sutherland, E., Coe, L. and Raleigh, E.A. (1992) *J. Mol. Biol.*, 225, 327.
5. Waite-Rees, P.A., Keating, C.J., Moran, L.S., Stalko, B.E., Hornstra, L.J. and Benner, J.S. (1991) *J. Bacteriol.*, 173, 5207.
6. Raleigh, E.A. (1987) *Meth. Enzymol.*, 152, 130.
7. Bickle, T. (1993) in *Nucleases* (S.M. Linn, R.S. Lloyd and R.J. Roberts, eds). Cold Spring Harbor Laboratory Press, New York, p. 89-109.
8. Ullman, A. and Perrin, D. (1970) in *The Lactose Operon* (J.R. Beckwith and D. Zipser, eds). Cold Spring Harbour Laboratory Press, New York, p. 143.

## Section 4: Technical Services

If you need assistance at any time, please call Active Motif Technical Service at one of the numbers listed below.

### Active Motif North America

1914 Palomar Oaks Way, Suite 150

Carlsbad, CA 92008

USA

Toll Free Tel: 877 222 9543

Telephone: 760 431 1263

Fax: 760 431 1351

E-mail: [tech\\_service@activemotif.com](mailto:tech_service@activemotif.com)

### Active Motif Europe

104 Avenue Franklin Roosevelt

B-1330 Rixensart, Belgium

UK Free Phone: 0800 169 31 47

France Free Phone: 0800 90 99 79

Germany Free Phone: 0800 181 99 10

Telephone: +32 (0)2 653 0001

Fax: +32 (0)2 653 0050

E-mail: [eurotech@activemotif.com](mailto:eurotech@activemotif.com)

### Active Motif Japan

Azuma Bldg, 7th Floor

2-21 Ageba-Cho, Shinjuku-Ku

Tokyo, 162-0824, Japan

Telephone: +81 3 5225 3638

Fax: +81 3 5261 8733

E-mail: [japantech@activemotif.com](mailto:japantech@activemotif.com)

Visit Active Motif on the worldwide web at <http://www.activemotif.com>

At this site:

- Read about who we are, where we are, and what we do
- Review data supporting our products and the latest updates
- Enter your name into our mailing list to receive our catalog, *Motivations* newsletter and notification of our upcoming products
- Share your ideas and results with us
- View our job opportunities

**Don't forget to bookmark our site for easy reference!**