

## HB101 Chemically Competent Cells Transformation Protocol

### Introduction

GoldBio's HB101 chemically competent *E. coli* cells allow you to obtain high transformation efficiency in applications such as cloning and sub-cloning. Our *E. coli* HB101 proceeds from the K12 x B hybrid strain. HB101 strain also contains the *recA13* mutation useful in the insert stability and minimization of recombination. *E. coli* HB101 also has the *hsdS20(rB-mB-)* genotype that prevents the cleavage of cloned DNA by endogenous restriction enzymes. Here, we present a detailed protocol for transformation using HB101 Chemically Competent *E. coli* cells.

### Materials

- HB101 Chemically Competent *E. coli* cells (GoldBio Catalog # CC-150)
- pUC19 Control DNA, 10 pg/μl
- Recovery medium (GoldBio Catalog # CC-300)

### Storage and Handling

- This product may be shipped on dry ice. HB101 Chemically Competent *E. coli* cells should be stored at -80°C, pUC19 Control DNA should be stored at -20°C and recovery medium should be stored at 4°C immediately upon arrival. When stored under the recommended conditions and handled correctly, these products should be stable for at least 1 year from the date of receipt.
- Thaw HB101 Chemically Competent *E. coli* cells and pUC19 Control DNA ice and mix by gentle vortexing. After thawing, these products should be kept on ice before use. These products can be refrozen for storage.

**Note:** The genotype of HB101 Chemically Competent *E. coli* cells *F- Lambda- araC14 leuB6(Am) DE(gpt-proA)62 lacY1 glnX44(AS) galK2(Oc) recA13 rpsL20(strR) xylA5 mtl-1 thiE1 hsdS20(rB-, mB-)*.

**Note:** Transformation efficiency is tested by using the pUC19 control DNA supplied with the kit and using given below. Transformation efficiency should be  $\geq 1 \times 10^8$  CFU/μg pUC19 DNA. Untransformed cells are tested for appropriate antibiotic sensitivity.

## Method

### Transformation protocol

Use this procedure to transform HB101 Chemically Competent *E. coli* cells. We recommend verifying the transformation efficiency of the cells using the pUC19 control DNA supplied with the kit. Do not use these cells for electroporation.

**Note: Handle the competent cells gently as they are highly sensitive to changes in temperature or mechanical lysis caused by pipetting.**

**Note: Thaw competent cells on ice and transform cells immediately following thawing. After adding DNA, mix by tapping the tube gently. Do not mix cells by pipetting or vortexing.**

1. Remove competent cells from the -80°C freezer and thaw completely on ice (10-15 minutes).
2. Aliquot 1-5 µl (1 pg-100 ng) of DNA into the chilled microcentrifuge tubes on ice.
3. When the cells are thawed, add 50 µl of cells to each DNA tube on ice and mix gently by tapping 4-5 times. For the pUC19 control, add 1 µl of (10 pg/µl) DNA to the 50 µl of cells on ice. Mix well by tapping. Do not pipette up and down or vortex to mix, this can harm cells and decrease transformation efficiency.
4. Incubate the cells with DNA on ice for 30 minutes.
5. After a 30-minute incubation on ice, heat shock the cells at 42°C for 45 seconds.
6. Transfer the tubes to ice for 2 minutes.
7. Add 950 µl of Recovery Medium or any other medium of choice to each tube.
8. Incubate tubes at 37°C for 1 hour at 210 rpm.
9. Spread 50 µl to 200 µl from each transformation on pre-warmed selection plates. We recommend plating two different volumes to ensure that at least one plate will have well-spaced colonies. For the pUC19 control, plate 50 µl on an LB plate containing 100 µg/ml ampicillin. Use a sterilized spreader or autoclaved plating beads to spread evenly.
10. Incubate the plates overnight at 37°C.

### 5 Minute Transformation Protocol

The following procedure results in only ~10% of the transformation efficiency as the protocol listed above.

1. Remove competent cells from the -80°C freezer and thaw the tube in your hand.
2. Aliquot 1-5  $\mu$ l (1 pg – 100 ng) of DNA to the microcentrifuge tubes. Do not pipette up and down or vortex to mix, this can harm cells and decrease transformation efficiency.
3. Incubate the cells with DNA on ice for 2 minutes.
4. After the 2-minute ice incubation, heat shock the cells at 42°C for 45 seconds.
5. Transfer the tubes to ice for 2 minutes.
6. Add 950  $\mu$ l of Recovery Medium at room temperature or any other medium of choice to each tube. Immediately spread 50  $\mu$ l to 200  $\mu$ l from each transformation onto pre-warmed selection plates. We recommend plating two different volumes to ensure that at least once plate will have well-spaced colonies. For the pUC19 control, plate 50  $\mu$ l on an LB plate containing 100  $\mu$ g/ml ampicillin. Use a sterilized spreader or autoclaved plating beads to spread evenly.
7. Incubate the plates overnight at 37°C.

## Calculations

Transformation efficiency (TE) is defined as the number of colony forming units (cfu) produced by transforming 1 µg of plasmid into a given volume of competent cells.

$$TE = \text{Colonies}/\mu\text{g}/\text{Dilution}$$

Where:

Colonies = the number of colonies counted

µg = amount of DNA transformed in µg

Dilution = total dilution of the DNA before plating

*Transform 1 µl of (10 pg/µl) pUC19 control plasmid into 50 µl of cells, add 950 µl of Recovery Medium. Dilute 10 µl of this in 990 µl of Recovery Medium and plate 50 µl. Count the colonies on the plate the next day. If you count 100 colonies, the TE is calculated as follows:*

$$\text{Colonies} = 100$$

$$\mu\text{g of DNA in } 10\text{pg} = 0.00001$$

$$\text{Dilution} = 50 \mu\text{l}/1000 \times 10 \mu\text{l}/1000 = 0.0005$$

$$TE = 100/0.00001/0.0005 = 2.0 \times 10^{10}$$

## Associated Products

- DH10B Chemically Competent *E. coli* Cells (GoldBio Catalog # CC-100)
- DH5-alpha Chemically Competent *E. coli* Cells (GoldBio Catalog # CC-101)
- Competent Cell Recovery Medium (GoldBio Catalog # CC-300)
- Ampicillin (GoldBio Catalog # A-301)