

# DH5α Competent Cells

Catalog Number EC0112

Pub. No. MAN0018520 Rev. A.0

 **WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](http://thermofisher.com/support).

## Product description

Thermo Scientific™ DH5α Competent Cells are ideal for construction of gene banks or generation of cDNA libraries using plasmid-derived vectors. The  $\phi 80\text{dlacZ}\Delta\text{M15}$  marker provides  $\alpha$ -complementation of the  $\beta$ -galactosidase gene from pUC or similar vectors to allow blue/white colony screening on bacterial agar plates containing Bluo-Gal or X-Gal.

## Genotype

F<sup>-</sup>  $\phi 80\text{dlacZ}\Delta\text{M15}$   $\Delta(\text{lacZYA-argF})$  U169 *recA1 endA1 hsdR17* (*r<sub>k</sub>*<sup>-</sup>, *m<sub>k</sub>*<sup>+</sup>) *phoA supE44*  $\lambda^-$  *thi-1 gyrA96 relA1*

## Contents and storage

Contents	Amount	Storage
DH5α Competent Cells	10 × 100 μL	-80°C (Do not store in liquid nitrogen)
pUC19 DNA (10 pg/μL)	50 μL	-80°C
S.O.C. Medium	6 mL	4°C or room temperature

## Guidelines for transforming cells

- For best results, thaw each vial of cells only once. Subsequent freeze-thaw cycles significantly lower transformation efficiency.
- Use S.O.C. Medium for the cell recovery procedures after heat shock. Transformation efficiency is reduced if other types of media are used.
- Maximum transformation efficiency is obtained with plasmid DNA that is free of phenol, ethanol, protein, and detergents. Transformation of unpurified sample DNA or ligation reactions will result in slightly lower transformation efficiencies.
- To determine the transformation efficiency of the cells, perform a control reaction using 10 pg (1 μL) of the pUC19 DNA stock solution. Dilute pUC19 DNA control reaction 1:10 with S.O.C. Medium immediately before plating. Spread 30 μL of the pUC19 DNA control reaction on a LB plate containing 100 μg/mL of ampicillin. The cells should have a transformation efficiency of  $\geq 1 \times 10^9$  cfu/μg.

## Transform competent cells

- Thaw competent cells on wet ice. Place the required number of 1.5-mL polypropylene microcentrifuge tubes on wet ice.
- Gently mix the cells, then make 50 μL aliquots of competent cells in the chilled 1.5-mL microcentrifuge tubes.
- Add 1–5 μL of sample DNA directly into a tube of competent cells. Mix well by gently flicking tube several times.
- Incubate the cells on ice for 30 minutes.
- Heat-shock the cells for exactly 30 seconds in a 42°C water bath. Do not mix or shake the tube.
- Incubate the cells on ice for 2 minutes.
- Add 250 μL of room-temperature S.O.C. Medium.
- Place the tube on its side in a shaking incubator. Use tape to secure the tube in place.
- Shake the tube at 225 rpm for 1 hour at 37°C.
- If necessary, dilute the cells 1:10 with S.O.C. Medium.
- Spread at least two different volumes (20–200 μL) of cells from each transformation reaction on separate LB plates containing the appropriate selective antibiotic. Label the plates with the plating volume so that the amount providing the best colony density can be identified.
- Invert the plates and incubate overnight at 37°C.

## Calculate transformation efficiency

Transformation efficiency can be 10- to 100-fold lower for transformation of a ligation reaction mixture than for transformation of an intact plasmid such as pUC19 DNA.

Calculate the transformation efficiency (CFU/μg) as follows:

$$\frac{\text{CFU in plate}}{\text{pg of DNA used in transformation}} \times \frac{1 \times 10^6 \text{ pg}}{\mu\text{g}} \times \text{dilution factor(s)}$$

For example, if 10 pg of pUC19 DNA yields 200 colonies when 30 μL of a 1:10 dilution is plated, then:

$$\text{CFU}/\mu\text{g} = \frac{200 \text{ CFU}}{10 \text{ pg}} \times \frac{1 \times 10^6 \text{ pg}}{\mu\text{g}} \times \frac{300 \mu\text{L}}{30 \mu\text{L plated}} \times 10 = 2 \times 10^9$$

## 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 at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).

---



Life Technologies Corporation | 5781 Van Allen Way | Carlsbad, CA 92008

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](http://thermofisher.com/symbols-definition).

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Important Licensing Information:** This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.