# DH5a Competent Cells for Subcloning

#### Catalog Number EC0111

Pub. No. MAN0018593 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**.

# **Product description**

Thermo Scientific<sup> $\mathbb{T}$ </sup> DH5 $\alpha$  Competent Cells for Subcloning are recommended for routine subcloning into plasmid vectors. Subcloning efficiency cells are not suitable for the generation of cDNA libraries. The  $\varphi 80dlacZ\Delta M15$  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

 $F^- \varphi 80 lac Z \Delta M 15 \Delta (lac ZYA-argF) U169 \ rec A1 \ end A1 \ hs dR17 \ (r_k-, m_k+) \ phoA \ sup E44 \ \lambda^- \ thi^{-1} \ gyrA96 \ rel A1 \ rel A1 \ rel A1 \ rel A1 \ rel A2 \ rel A2 \ rel A2 \ rel A2 \ rel A3 \ rel A3 \ rel A3 \ rel A3 \ rel A4 \ rel$ 

### **Contents and storage**

Contents	Amount	Storage
DH5a Competent Cells for Subcloning	4 × 500 μL	–80°C (Do not store in liquid nitrogen)
pUC19 DNA (100 pg/µL)	20 µL	-80°C

### Guidelines for tranforming cells

- For best results, thaw each vial of cells only once. Subsequent freeze-thaw cycles significantly lower transformation efficiency.
- 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 250 pg (2.5  $\mu$ L) of the pUC19 DNA stock solution. Spread 100  $\mu$ L of the pUC19 DNA control reaction on a LB plate containing 100  $\mu$ g/mL of ampicillin. The cells should have a transformation efficiency of  $\geq$ 1 × 10<sup>6</sup> cfu/ $\mu$ g.

# Transform competent cells

- 1. Thaw competent cells on wet ice. Place the required number of 1.5-mL polypropylene microcentrifuge tubes on wet ice.
- 2. Gently mix the cells, then make  $50 \ \mu L$  aliquots of competent cells in the chilled 1.5-mL microcentrifuge tubes.
- 3. Add 1–5  $\mu$ L of sample DNA directly into a tube of competent cells. Mix well by gently flicking tube several times.
- **4.** Incubate the cells on ice for 30 minutes.
- Heat-shock the cells for exactly 20 seconds in a 42°C water bath. Do not mix or shake the tube.
- 6. Incubate the cells on ice for 2 minutes.
- 7. Add 950 µL of pre-warmed growth medium (e.g., S.O.C. or LB).
- **8**. Place the tube on its side in a shaking incubator. Use tape to secure the tube in place.
- 9. Shake the tube at 225 rpm for 1 hour at 37°C.
- 10. 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.
- 11. Invert the plates and incubate overnight at 37°C.

### **Calculate transformation efficiency**

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

CFU in plate	v	<u>1 × 10<sup>6</sup> pg</u>	v	dilution factor(s)
og of DNA used in transformation	^	hà	^	ullution lactor(s)

For example, if 250 pg of pUC19 DNA yields 100 colonies when 100  $\mu$ L of a 1:10 dilution is plated, then:

CFU/µg = -	100 CFU	x	1 x 10 <sup>6</sup> pg	х	1000 µL	$x 10 = 4 \times 10^6$
	250 pg		μg		100 µL plated	X 10 - 4X 10



# 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. If you have any questions, please contact Life Technologies at 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**.

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.