# MultiShot<sup>™</sup> StripWell Competent Cell Kits

**Catalog Numbers** C8696-01, C7396-01, C6096-01, C4496-01, C4096-01, C4096-05, and C4096-10

Publication Number MAN0017711 Revision 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**

The Invitrogen  $^{\text{\tiny{TM}}}$  MultiShot  $^{\text{\tiny{TM}}}$  StripWell Competent Cell Kits contain chemically competent  $E.\ coli$  cells in StripWell tubes to simplify medium- and high-volume bacterial transformations. The StripWell format provides the flexibility to set up the exact number of transformations required for your particular experiment.

## Contents and storage

Kit type	Cat. No.
MultiShot <sup>™</sup> StripWell BL21 Star <sup>™</sup> (DE3) Competent Cells, 1 rack	C6096-01
MultiShot <sup>™</sup> StripWell DH5a <sup>™</sup> T1 <sup>R</sup> Competent Cells, 1 rack	C4496-01
MultiShot <sup>™</sup> StripWell Mach1 <sup>™</sup> T1 <sup>R</sup> Competent Cells, 1 rack	C8696-01
MultiShot <sup>™</sup> StripWell Stbl3 <sup>™</sup> Competent Cells, 1 rack	C7396-01
MultiShot <sup>™</sup> StripWell TOP10 Competent Cells, 1 rack	C4096-01
MultiShot <sup>™</sup> StripWell TOP10 Competent Cells, 5 racks	C4096-05
MultiShot <sup>™</sup> StripWell TOP10 Competent Cells, 10 racks	C4096-10

Contents	1 rack kit	5 rack kit	10 rack kit	Storage
Chemically competent cells (50 µL/tube)	96 StripWell tubes	5 × 96 StripWell tubes	10 × 96 StripWell tubes	-80°C to -70°C
pUC19 vector, supercoiled (10 pg/ μL) [1]	50 μL	50 μL	2 × 50 μL	-60 0 10 -70 0
S.O.C. medium	2 × 15 mL	10 × 15 mL	20 × 15 mL	Room temperature or 4°C

 $<sup>^{[1]}</sup>$  In 5 mM Tris-HCl, 0.5 mM EDTA, pH 8



## Genotype

Strain	Genotype	
BL21 Star <sup>™</sup> (DE3)	F-ompT hsdSB (r <sub>B</sub> - m <sub>B</sub> -) gal dcm rne131 (DE3)	
DH5a <sup>™</sup> T1 <sup>R</sup>	$F^-$ Φ80 $lac$ ZΔM15 Δ( $lac$ ZYA- $arg$ F)U169 $rec$ A1 $end$ A1 $hsd$ R17 ( $r_{K}^ m_{K}^+$ ) $pho$ A $sup$ E44 $thi$ -1 $gyr$ A96 $rel$ A1 $ton$ A	
Mach1 <sup>™</sup> T1 <sup>R</sup>	F- Φ80 $lac$ ZΔM15 Δ $lac$ X74 $hsd$ ( $r_{K}^-m_{K}^-$ ) Δ $rec$ A1398 $end$ A1 $ton$ A	
Stbl3 <sup>™</sup>	F^ $mrrhsd$ S20(r <sub>B</sub> $^-$ m <sub>B</sub> $^-$ ) $rec$ A13 $sup$ E44 $ara$ -14 $gal$ K2 $lac$ Y1 $pro$ A2 $rps$ L20(Str $^R$ ) $xyl$ -5 $\lambda^ leu$ $mtl$ -1	
T0P10	F- mcrA Δ(mrr-hsdRMS-mcrBC) Φ80lacZΔM15 ΔlacX74 recA1 araD139 Δ(araleu)7697 galU galK rpsL20(Str <sup>R</sup> ) endA1 nupG	

## Required materials not provided

- 42°C water bath (High Efficiency Transformation Protocol only)
- 37°C shaking incubator
- LB agar plates with appropriate antibiotic (e.g., ampicillin for transformation controls)
- Ice bucket with ice

## Procedural guidelines

- **IMPORTANT!** Do not transform chemically competent cells by electroporation. The salt content of the buffer will cause arcing and kill the cells.
- Be extremely gentle when working with competent cells. Competent cells are highly sensitive to changes in temperature or mechanical lysis caused by pipetting. Start transformation immediately after thawing the cells on ice. Mix cells by gently tapping the tubes, or stirring with a pipette tip; do not mix by pipetting up and down.
- Practice sterile technique when handling S.O.C. medium to avoid contamination.
- Avoid freeze-thaw cycles for pUC19 control DNA. Prepare aliquots of pUC19 control DNA and store at -80°C.
- Minimize handling of StripWell tubes. Remove the required number of tubes for an experiment and return the unused tubes to the freezer.
- Handle StripWell tubes carefully when using a water bath to avoid accidental contamination of cells.
- Perform a transformation control to test the transformation efficiency of the competent cells in the kit. See "Calculate transformation efficiency" on page 5 for details.

## **Transformation procedures**

Two protocols are provided to transform MultiShot<sup>TM</sup> StripWell Competent Cells. Consider the following factors when choosing which protocol to use.

If you wish to	Then use the	
maximize the number of transformants obtained	"High Efficiency Transformation Protocol"	
use an antibiotic other than ampicillin to select for your DNA		
obtain transformants as quickly as possible	"Rapid Transformation Protocol" on page 4	

#### High Efficiency Transformation Protocol

The High Efficiency Transformation Protocol is used to achieve the highest efficiency when transforming MultiShot<sup>™</sup> competent cells with your DNA of interest.

#### Before you begin

- Warm S.O.C. medium to room temperature if necessary.
- Pre-warm LB selection plates in a 37°C incubator for at least 30 minutes.
- Pre-heat water bath or heat block to 42°C.

#### Prepare cells

- 1. Thaw tube(s) MultiShot<sup>™</sup> competent cells on ice for 2–5 minutes.
- **2.** Carefully remove the caps from the tube(s) and set them aside for later use.

#### Transform cells

1. Add  $\leq 5 \mu L$  of DNA directly into each tube.

**Note:** For optimal results, do not add more than 5  $\mu$ L (10% of the cell volume) of DNA.

- **2.** Mix the cells by gently stirring the mixture with a pipette tip three times. **Do not** mix by pipetting up and down.
- **3.** Cap the tubes and incubate the cells on ice for 30 minutes.
- **4.** Incubate the tube(s) of cells at 42°C in a water bath. Do not mix or shake.

Strain	42°C incubation
BL21 Star <sup>™</sup> (DE3) , DH5a <sup>™</sup> T1 <sup>R</sup> , Mach1 <sup>™</sup> T1 <sup>R</sup> , and T0P10	30 seconds
Stbl3 <sup>™</sup>	45 seconds

**5.** Incubate the tube(s) on ice for 2 minutes.

**6.** Place the tube(s) on the benchtop, then add 250  $\mu L$  of S.O.C. medium to each tube.

**Note:** S.O.C. is a rich medium; good sterile technique must be practiced to avoid contamination.

- 7. Cap the tube(s) and secure them at a 45° angle in a shaking incubator for optimal agitation and aeration.
- **8.** Incubate the tube(s) at 37°C with shaking at 225 rpm for 1 hour.
- **9.** Dilute transformation reaction with S.O.C. medium as needed.

Strain	Dilution factor
DH5a <sup>™</sup> T1 <sup>R</sup> , Mach1 <sup>™</sup> T1 <sup>R</sup> , and T0P10	1:10
BL21 Star <sup>™</sup> (DE3) and Stbl3 <sup>™</sup>	No dilution

**Note:** Ensure the transformation reaction is thoroughly mixed before dilution.

#### Plate transformed cells

1. Plate 50–100  $\mu$ L of cells from each transformation reaction on pre-warmed LB selection plates, then invert the plates and incubate overnight at 37°C.

**Note:** Ensure the transformation reaction is thoroughly mixed before plating.

2. Select colonies for further analysis by plasmid isolation, PCR, or sequencing.

#### Rapid Transformation Protocol

The Rapid Transformation Protocol is used to obtain transformants as quickly as possible, and is only compatible with vectors with ampicillin resistance. Using this protocol to transform  $MultiShot^{TM}$  competent cells can result in a 1-log reduction in transformation efficiency.

#### Before you begin

Pre-warm LB selection plates in a 37°C incubator long enough to ensure the plates reach 37°C (a minimum of 1 hour). Prepare one plate for each transformation.

#### Transform competent cells

- 1. Thaw tube(s) of MultiShot<sup>™</sup> competent cells on ice for 2–5 minutes.
- **2.** Add  $\leq 5 \mu L$  of DNA directly into each tube.
- **3.** Mix the cells by gently stirring the mixture with a pipette tip three times. **Do not** mix by pipetting up and down.
- 4. Cap the tubes and incubate the cells on ice for 5 minutes.
- 5. Immediately plate 5–20  $\mu$ L of cells from each transformation reaction on prewarmed LB selection plates, then invert the plates and incubate overnight at 37°C.

#### Perform transformation control

- 1. Prepare competent cells as described in "Prepare cells".
- 2. Add 1  $\mu$ L of the pUC19 control DNA to thawed cells, then follow the instructions in "Transform cells" starting from step 2.
- 3. Plate 30–50  $\mu$ L of cells on pre-warmed LB agar plates containing 100  $\mu$ L/mL ampicillin (see "Plate transformed cells" for detailed instructions).

## Calculate transformation efficiency

Use the following formula to calculate transformation efficiency in terms of colony forming units per  $\mu g$  of DNA (CFU/ $\mu g$ ).

$$\frac{\text{Transformation efficiency}}{(\text{CFU/}\mu\text{g})} = \frac{\text{\# of colonies (CFU)}}{\mu\text{g transformed DNA}} \times \frac{\text{Transformation volume (}\mu\text{L})}{\text{Plating volume (}\mu\text{L})} \times \frac{\text{Dilution factor of transformed cells}}{\text{Transformation volume (}\mu\text{L})}$$

Use the following formula to determine the amount of DNA used for transformation for ligation reactions.

$$\mu g$$
 transformed DNA =  $\frac{DNA \text{ in ligation reaction } (\mu g)}{Ligation \text{ volume } (\mu L)}$   $\times$   $\frac{Dilution factor of}{ligation product}$   $\times$   $\frac{Volume \text{ of ligation product}}{used \text{ for transformation}}$ 

#### Example of transformation efficiency calculation

50 ng of DNA is ligated in a 20  $\mu$ L reaction. After ligation, the reaction is diluted 2-fold and 5  $\mu$ L of the diluted ligation mixture is added to 100  $\mu$ L of competent cells for transformation.

$$\mu$$
g transformed DNA =  $\frac{0.5 \mu g}{20 \mu L}$  × 2 × 5  $\mu$ L = 0.025  $\mu$ g

After transformation, the cell suspension is diluted 5-fold and 200  $\mu$ L of the diluted cells are plated. 300 colonies are formed after overnight incubation.

Transformation efficiency (CFU/
$$\mu$$
g) =  $\frac{300 \text{ CFU}}{0.025 \, \mu \text{g}}$  ×  $\frac{100 \, \mu \text{L}}{200 \, \mu \text{L}}$  × 5 =  $3 \times 10^4 \, \text{CFU/} \mu \text{g}$ 

## **Troubleshooting**

For troubleshooting help on competent cells, go to **thermofisher.com/ compcellsupport** 

Observation	Possible cause	Recommended action
Few or no colonies	Problem with competent cells.	Carry out the pUC19 control DNA transformation to obtain information about the performance of the cells.
	Problem with the antibiotic in the plate.	Confirm that the correct selection antibiotic and concentration was used in the LB agar plate.
	Using a vector with a selection marker that is not ampicillin when performing the Rapid Transformation Protocol.	If the antibiotic selection marker for the vector being transformed is not ampicillin, use the High Efficiency Transformation Protocol.
	LB plate used for growing colonies insufficiently warmed.	Ensure that the LB agar plates are prewarmed to 37°C prior to plating. Incubate LB agar plates for at least 30 minutes if using the High Efficiency Transformation Protocol, and for at least 1 hour if using the Rapid Transformation Protocol.
Transformants contain incorrect or truncated DNA	Instability in DNA being cloned.	Use Stbl3™ competent cells for transformation to prevent plasmid recombination.
inserts	DNA mutation.	Mutations may have occurred during plasmid propagation in transformed cells. Pick a sufficient number of colonies for representative screening by sequencing. If all colonies carry the same mutation, it may have originated from the original template.

## **Accessory products**

Product	Cat. No.
LB Broth, 500 mL	10855021
LB Broth Base, powder 500 g	12780052
LB Agar, powder 500 g	22700025
X-gal, 100 mg	15520034
S.O.C. Medium, 10 × 10 mL	15544034
StripWell Caps	15088

## For European customers

The Mach1<sup>™</sup> T1 Phage-Resistant Chemically Competent *E. coli* strain is genetically modified to carry the  $lacZ\Delta M15$  hsdR lacX74 recA endA tonA genotype. As a condition of sale, this product must be in accordance with all applicable local legislation and guidelines, including EC Directive 90/219/EEC on the contained use of genetically modified organisms.

## **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 <a href="https://www.thermofisher.com/us/en/home/global/terms-and-conditions.html">www.thermofisher.com/us/en/home/global/terms-and-conditions.html</a>. If you have any questions, please contact Life Technologies at <a href="https://www.thermofisher.com/support">www.thermofisher.com/support</a>.



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

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES 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.

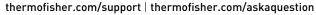
#### Revision history: Pub. No. MAN0017711

······································			
Revision	Date	Description	
Α	11 April 2018	New document	

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

TRADEMARKS: All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2018 Thermo Fisher Scientific Inc. All rights reserved.



thermofisher.com

