

# LyoComp GT116

Lyophilized competent *E. coli* GT116 cells

Catalog # lyo-116-11, lyo-116-21

For research use only

Version # 09G07-MM

## PRODUCT INFORMATION

### Content:

LyoComp GT116 cells are provided lyophilized and are available in two sizes;

- lyo-116-11: 4 x 0.5 ml (5 - 10 transformations)
- lyo-116-21: 4 x 1 ml (10-20 transformation)

With each product, 8 ml of sterile reconstitutive solution is provided.

**GT116 Genotype:** *F* *mcrA*  $\Delta$ (*mrr-hsdRMS-mcrBC*)  $\phi$ 80*lacZ*M15  $\Delta$ *lacX74* *recA1* *rspL* (*StrA*) *endA1*  $\Delta$ *sbcC-sbcD*

### Storage and stability:

- LyoComp cells are shipped at room temperature.
- Upon receipt, store LyoComp cells at -20°C.
- LyoComp cells are stable for 6 months when stored at -20°C.

### Quality control:

The transformation efficiency of LyoComp cells are evaluated periodically and are guaranteed to be stable for 6 months when properly stored (-20°C).

## DESCRIPTION / PROPERTIES

GT116 is a *sbcCD* deletion strain specifically engineered to support the growth of plasmid DNAs carrying hairpin structures, such as psiRNA a vector expressing small interfering RNA. Hairpin structures are known to be unstable in *E. coli* due to their elimination by a protein complex called SbcCD that recognizes and cleaves hairpins<sup>1</sup>. To increase their stability in *E. coli*, InvivoGen has developed GT116 by deleting the *sbcC* and *sbcD* genes. This modification significantly improves the number of recombinant clones harboring a plasmid with hairpin structures. This strain contains the *rpsL* (*StrA*) gene which confers resistance to streptomycin.

LyoComp GT116 was specifically designed for transformation with psiRNA ligation products or recombinant plasmids. LyoComp GT116 cells exhibit a lower competency than standard chemically competent cells but the competency is sufficient to allow successful cloning into psiRNA plasmids.

Transformation efficiency:  $1 \times 10^6$  cfu/ $\mu$ g

1. Connelly JC. et al., 1998. The SbcCD nuclease of escherichia coli is a structural maintenance of chromosomes (SMC) family protein that cleaves hairpin DNA. Proc. Natl. Acad. Sci. USA 95:7969-7974

## TRANSFORMATION

The following protocol describes a method used to introduce DNA into bacterial host for efficient and convenient construction or maintenance of plasmid recombinants, and blue/white screening.

### Additional required materials to be supplied by user:

- LB agar plates with appropriate antibiotic. For optimal results we recommend the use of InvivoGen's selective Fast-Media®.
- 37°C shaking Incubator
- Ice bucket
- 42°C water bath
- LB or SOC medium

### Method:

#### **Before starting:**

- Prepare LB agar plates containing the appropriate antibiotic.
- Set water bath to 42°C.
- Pre-chill appropriate number of 1.5 ml sterile tubes in ice.

- 1- Place the appropriate number of LyoComp GT116 cell vials (100  $\mu$ l per ligation or transformation reaction) in ice for 5 minutes.
- 2- Rehydrate LyoComp GT116 using cold reconstitutive solution
  - for lyo-116-11, add 0.5 ml per vial
  - for lyo-116-21, add 1 ml per vialand store in ice for 5 minutes.
- 3- Gently homogenize and allow the cells to completely rehydrate in ice for 25-30 minutes.
- 4- Introduce 10  $\mu$ l of ligation product (or 1  $\mu$ g supercoiled plasmid DNA in pre chilled 1.5 ml tubes and return tubes to ice.
- 5- Gently flick the cells twice to homogenize and add 100  $\mu$ l of cells to each DNA-containing tubes.
- 6- Mix by tapping gently and place in ice immediately.
- 7- Incubate the tubes in ice for 30 minutes.
- 8- Incubate the tubes in a 42°C water bath for exactly 30 seconds, then place the tubes back in ice for 1-2 minutes.
- 9- Add 900  $\mu$ l of room temperature SOC (or LB) medium to each reaction. (Practice sterile techniques to avoid contamination.)
- 10- Incubate tubes at 37°C for 1h30 with shaking at 250 rpm.
- 11- Spread each transformation reaction (4 x 150  $\mu$ l if using a ligation product or 100  $\mu$ l of  $10^{-1}$  and  $10^{-2}$  dilutions if using a supercoiled plasmid) onto LB agar plates containing the appropriate antibiotic.
- 12- Incubate plates at 37°C overnight.

## TECHNICAL SUPPORT

Toll free (US): 888-457-5873  
Outside US: (+1) 858-457-5873  
Europe: +33 562-71-69-39  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)  
Website: [www.invivogen.com](http://www.invivogen.com)



3950 Sorrento Valley Blvd. Suite A  
San Diego, CA 92121 - USA