

USER GUIDE

invitrogen™
by *life* technologies™

GeneArt® *Synechococcus* Engineering Kits

For expression of recombinant proteins in
Synechococcus elongatus

Catalog Numbers A14259, A14263

Publication Number MAN0005339

Revision 2.0

For Research Use Only. Not for use in diagnostic procedures.

life
technologies™

Information in this document is subject to change without notice.

DISCLAIMER

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

NOTICE TO PURCHASER: LIMITED USE LABEL LICENSE: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008.

TRADEMARKS

The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate(s) or their respective owners.

© 2013 Life Technologies Corporation. All rights reserved.

Contents

Product Information	2
Contents and Storage.....	2
Description of the System	4
Experiment Outline.....	6
Methods	7
Cloning into pSyn_1.....	7
Transforming One Shot® TOP10 Competent <i>E. coli</i> Cells.....	9
Analyzing <i>E. coli</i> Transformants	10
Guidelines for Culturing <i>Synechococcus elongatus</i> PCC 7942.....	12
Thawing <i>Synechococcus elongatus</i>	13
Transforming <i>Synechococcus elongatus</i>	15
Screening for Integration by Colony PCR.....	17
Appendix A: Support Protocols	18
Media and Plates.....	18
Appendix B: Vectors	19
Map and Features of pSyn_1 Vector	19
Map of pSyn_2/Control Vector	21
Appendix C: Ordering Information	22
Accessory Products.....	22
Documentation and Support	24
Obtaining Support.....	24
References.....	25

Product Information

Contents and Storage

Types of kits

This manual is supplied with the products listed below. For a list of components supplied with each catalog number, see below.

Product	Catalog number
GeneArt® <i>Synechococcus</i> Engineering Kit	A14259
GeneArt® <i>Synechococcus</i> Engineering Kit with 6 L media	A14263

Kit components

Each GeneArt® *Synechococcus* Engineering Kit contains the components listed below. See the next page for a detailed description of each of the components.

Box	Component	Catalog number	
		A14259	A14263
1	GeneArt® <i>Synechococcus elongatus</i> Cells	✓	✓
2	GeneArt® <i>Synechococcus</i> Vector Set	✓	✓
3	Gibco® BG-11 Media		✓

Shipping/Storage

The GeneArt® *Synechococcus* Engineering Kits are shipped in separate boxes as described below. Upon receipt, store each box as detailed below. All reagents are guaranteed for six months if stored properly.

Box	Component	Shipping	Storage
1	GeneArt® <i>Synechococcus elongatus</i> Cells	Dry ice	-80°C
2	GeneArt® <i>Synechococcus</i> Vector Set	Dry ice	-20°C
3	Gibco® BG-11 Media	Gel ice	4°C

Continued on next page

Contents and Storage, continued

GeneArt® *Synechococcus* *elongatus* Cells

Each GeneArt® *Synechococcus* Engineering Kit is supplied with 10 vials of GeneArt® *Synechococcus elongatus* PCC 7942 cells, with each vial containing 100 µL of frozen cells. Store the cells at –80°C upon receipt.

GeneArt® *Synechococcus* Vector Set

The table below lists the components of the GeneArt® *Synechococcus* Vector Set (Box 2). Store the contents of Box 2 at –20°C.

Component	Concentration	Amount
pSyn_1 Vector	20 µL of vector at 0.5 µg/µL in TE buffer, pH 8.0*	10 µg
pSyn_2/Control Vector	20 µL of vector at 0.5 µg/µL in TE buffer, pH 8.0	10 µg

*TE buffer, pH 8.0: 10 mM Tris-HCl, 1 mM EDTA, pH 8.0

Gibco® BG-11 Media

Gibco® BG-11 Media, included in the GeneArt® *Synechococcus* Engineering Kit with 6 L media (Cat. no. A14265), is supplied in 6 × 1 L bottles and is optimized for the growth and maintenance of GeneArt® *Synechococcus elongatus* PCC 7942 cells. Store the Gibco® BG-11 Medium at 4°C.

Note: Gibco® BG-11 Media (Cat. nos. A1379901, A1379902) are also available separately from Life Technologies. See page 22 for ordering information.

Description of the System

Introduction

The GeneArt® *Synechococcus* Engineering Kit is a prokaryotic photosynthetic model system based on the unicellular cyanobacterium *Synechococcus elongatus* PCC 7942 (<http://genome.kazusa.or.jp/cyanobase/SYNPCC7942>), which offers a simplified approach for cloning and expressing genes of interest for the study of circadian rhythms, nutrient regulation, environmental response, lipid metabolism, and protein expression.

Synechococcus elongatus PCC 7942

Synechococcus elongatus PCC 7942 is a freshwater unicellular cyanobacterium. Cyanobacteria, sometimes referred to as blue-green algae, are prokaryotes that are able to obtain their energy through photosynthesis. *Synechococcus elongatus* has a rod-shaped appearance and is oligotrophic, having the ability to survive in freshwater environments with low nutrients. This organism has a circular chromosome of ~2.7 Mb (fully sequenced) with a GC content of 55.5%, which contains the genes for 2,612 proteins and 53 RNAs (Vijayan *et al.*, 2011).

The cyanobacterium *Synechococcus elongatus* PCC 7942 is an excellent synthetic biology chassis and a model system for studying prokaryotic circadian rhythms, nutrient regulation, environmental responses, and lipid metabolism because of its small genome size and the ease with which it can be genetically manipulated by natural transformation or conjugation from *E. coli* (Atsumi *et al.*, 2009; Ducat *et al.*, 2011; Lan & Liao, 2011; Min & Golden, 2000; Simkovsky *et al.*, 2012; Taniguchi *et al.*, 2012).

Transformation of *Synechococcus elongatus* PCC 7942

The transformation of *Synechococcus elongatus* PCC 7942 relies on homologous recombination between the cell's chromosome and exogenous DNA that is not autonomously replicating and containing sequences homologous to the chromosome. The location of integration into the chromosome (neutral site, NS1) has been developed as a cloning locus (Clerico *et al.*, 2007) as it can be disrupted without any aberrant phenotype, thus allowing the homologous recombination of ectopic sequences. When transformed with vectors containing an antibiotic resistance cassette and neutral site sequences, a double homologous recombination event occurs between the neutral site vector and the *Synechococcus elongatus* chromosome. The selective marker (Spectinomycin) and the gene of interest driven by a promoter are inserted into the neutral site and the vector backbone (pUC) is lost, allowing the expression of recombinant genes in *Synechococcus elongatus* PCC 7942.

Continued on next page

Description of the System, continued

pSyn_1 Vector

The pSyn_1 Vector is designed to facilitate rapid cloning of your gene of interest (GOI) for expression in *Synechococcus elongatus* PCC 7942. Some of the features of the vector are listed below. For a map of the vector, see page 19.

- A versatile multiple cloning site for simplified cloning of your gene of interest
- NS1 (neutral site 1) homologous recombination sites for the integration of the vector into the *Synechococcus elongatus* genome
- The weak constitutive promoter of solanesyl diphosphate synthase gene from *Synechocystis* sp. strain PCC 6803 driving the basal expression of your gene of interest

Note: Use of this weak constitutive promoter is a good fit for applications that are hindered by strong expression, such as pathway engineering or complementation of mutant genes normally expressed at low levels (Kaneko *et al.*, 1996; Simkovsky *et al.*, 2012).

- Spectinomycin resistance gene for selection in *E. coli* and *Synechococcus elongatus* PCC 7942
 - pUC origin for maintenance in *E. coli*
-

Experiment Outline

Workflow

The table below describes the major steps needed to clone and express your gene of interest (GOI) in *Synechococcus elongatus* PCC 7942. For details, refer to the pages indicated.

Step	Action	Page
1	Clone your gene of interest (GOI) directly into pSyn_1Vector	7
2	Transform One Shot® TOP10 <i>E. coli</i> with pSyn_1 Vector containing your GOI and select the transformants on LB plates containing spectinomycin	9
3	Analyze transformants by restriction digestion or PCR	10
4	Thaw and resuscitate <i>Synechococcus elongatus</i> cells	13
5	Transform <i>Synechococcus elongatus</i> cells and select transformants	15
7	Perform colony PCR to screen the transformed <i>Synechococcus elongatus</i> colonies for full integration of the GOI	17

Methods

Cloning into pSyn_1

General molecular biology techniques

For help with PCR amplification, DNA ligations, *E. coli* transformations, restriction enzyme analysis, DNA sequencing, and DNA biochemistry, refer to *Molecular Cloning: A Laboratory Manual* (Sambrook *et al.*, 1989) or *Current Protocols in Molecular Biology* (Ausubel *et al.*, 1994).

E. coli host

For cloning and transformation, we recommend using a recombination deficient (*recA*) and endonuclease A-deficient (*endA*) strain such as TOP10 (available separately; see page 22). Note that other *recA*, *endA* *E. coli* strains are also suitable.

Genotype of TOP10:

F *mcrA* Δ (*mrr-hsdRMS-mcrBC*) Φ 80*lacZ* M15 Δ *lac74* *recA1* *araD139* Δ (*ara-leu*)7697 *galU galK rpsL* (Str^R) *endA1 nupG*

Maintaining pSyn_1 vector

To propagate and maintain the pSyn_1 vector, use 10 ng of the vector to transform a *recA*, *endA* *E. coli* strain like TOP10, DH5 α TM, JM109, or equivalent. Select *E. coli* transformants on LB plates containing 100 μ g/mL spectinomycin. Be sure to prepare a glycerol stock of the plasmid for long-term storage (see page 11 for a protocol).

Cloning considerations

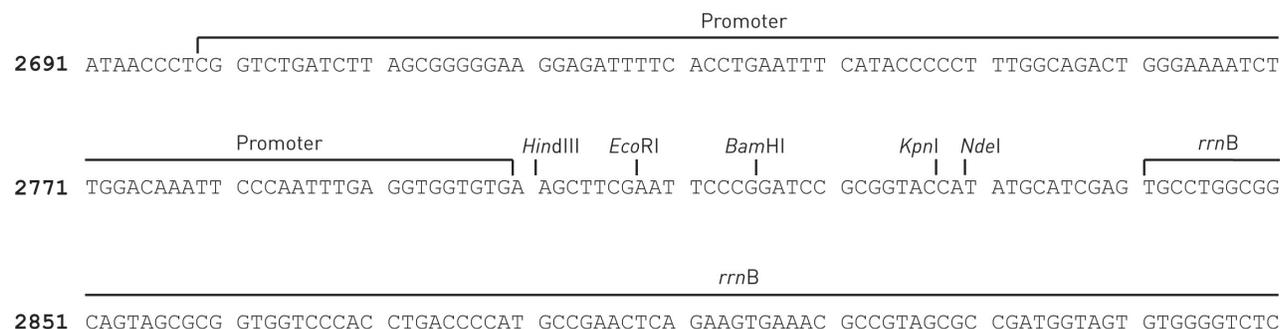
- pSyn_1 vector does not contain an ATG initiation codon for proper initiation of translation. Be sure to design your insert to contain an ATG initiation sequence. You can either use the native sequence containing the start codon or design your forward PCR primer so that your insert is amplified with the start codon in frame with your GOI.
 - pSyn_1 vector does not contain a ribosome binding site (RBS). To obtain maximal level of translation, you can include a strong RBS (such as GAAGGAG) 9–11 bp from the ATG initiation codon of your gene of interest. You can also use the RBS calculator programs from Life Technologies Corporation's Vector NTI[®] product lines or web based tools to design RBSs with various strengths.
 - Your insert must also contain a stop codon for proper termination of your mRNA. Although the pSyn_1 vector already contains a *rrnB* transcription termination region, you need to include a stop codon at the end of your gene of interest to guarantee the proper termination of your gene.
-

Continued on next page

Cloning into pSyn_1, continued

Multiple cloning site of pSyn_1

Below is the multiple cloning site for pSyn_1. Restriction sites are labeled to indicate the cleavage site. The multiple cloning site has been confirmed by sequencing and functional testing. Use the diagram below to design suitable PCR primers to clone and express your PCR product in pSyn_1. The vector sequence of pSyn_1 is available for downloading at www.lifetechnologies.com or by contacting Technical Support (page 24).



Ligation

Once you have determined a cloning strategy and PCR amplified your gene of interest, digest pSyn_1 with the appropriate restriction enzyme(s) and ligate your insert containing your gene of interest into pSyn_1 using standard molecular biology techniques.

E. coli transformation method

You may use any method of your choice for transformation. Chemical transformation is the most convenient for most researchers. Electroporation is the most efficient and the method of choice for large plasmids. For your convenience, a protocol for chemical transformation using One Shot[®] TOP10 Chemically Competent *E. coli* is provided on page 9; however, you may also transform electrocompetent cells.

Transforming One Shot[®] TOP10 Competent *E. coli* Cells

Introduction

Once you have performed the cloning reaction, you will transform your pSyn_1 construct into competent *E. coli*. The following protocol for transforming One Shot[®] TOP10 Chemically Competent *E. coli* (available separately; see page 22) is included for your convenience. Note that you may also transform electrocompetent cells using the protocol supplied with the electrocompetent cells.

Materials needed

- pSyn_1 construct containing your gene of interest
 - One Shot[®] TOP10 Chemically Competent *E. coli* (Cat. no. C4040; see page 22)
 - S.O.C. Medium (supplied with Cat. no. C4040)
 - pSyn_2 positive control (recommended for verifying transformation efficiency)
 - 42°C water bath
 - LB plates containing 100 µg/mL spectinomycin (two for each transformation)
 - 37°C shaking and non-shaking incubator
-

Preparing for transformation

For each transformation, you will need one vial of competent cells and two selective plates.

1. Equilibrate a water bath to 42°C.
 2. Warm the vial of S.O.C. medium to room temperature.
 3. Warm LB plates containing 100 µg/mL spectinomycin at 37°C for 30 minutes.
 4. Thaw **on ice** 1 vial of One Shot[®] TOP10 for each transformation.
-

One Shot[®] chemical transformation protocol

1. Add 1–5 µl of the DNA (10 pg to 100 ng) into a vial of One Shot[®] Chemically Competent *E. coli* and mix gently. **Do not mix by pipetting up and down.**
 2. Incubate on ice for 5 to 30 minutes.
Note: Longer incubations on ice seem to have a minimal effect on transformation efficiency.
 3. Heat-shock the cells for 30 seconds at 42°C without shaking.
 4. Immediately transfer the tubes to ice and incubate on ice for 2 minutes.
 5. Add 250 µL of room temperature S.O.C. Medium.
 6. Cap the tube tightly and shake the tube horizontally (200 rpm) at 37°C for 1 hour.
 7. Spread 50–200 µL from each transformation on a prewarmed selective plate and incubate overnight at 37°C. We recommend that you plate two different volumes to ensure that at least one plate will have well-spaced colonies.
 8. Proceed to **Analyzing *E. coli* Transformants**, page 10.
-

Analyzing *E. coli* Transformants

Picking positive clones

1. Pick 5–10 colonies and culture them overnight in LB medium containing 100 µg/mL spectinomycin.
 2. Isolate plasmid DNA using your method of choice. If you need ultra-pure plasmid DNA for automated or manual sequencing, we recommend using the PureLink® HQ Mini Plasmid Purification Kit (Cat. no. K2100-01, see page 22).
 3. Analyze the plasmids by restriction analysis or PCR (see below) to confirm the presence and correct orientation of the insert.
-

Analyzing transformants by PCR

Use the protocol below (or any other suitable protocol) to analyze positive transformants using PCR. You will have to determine the primer sequences and amplification conditions based on your gene of interest. Design one of the primers to hybridize to the vector backbone flanking your insert and the other one to hybridize within your insert. If you are using this technique for the first time, we recommend performing restriction analysis in parallel. Artifacts may be obtained because of mispriming or contaminating template.

Materials Needed:

- PCR Super Mix High Fidelity (Cat. no. 10790-020)
- Appropriate forward and reverse PCR primers (20 µM each)

Procedure:

1. For each sample, aliquot 48 µL of PCR SuperMix High Fidelity into a 0.2 mL PCR tube. Add 1 µL each of the forward and reverse PCR primer.
 2. Pick 5–10 colonies and resuspend them individually in 50 µL of the PCR SuperMix containing PCR primers (remember to make a patch plate to preserve the colonies for further analysis).
 3. Incubate reaction for 10 minutes at 94°C to lyse the cells and inactivate the nucleases before proceeding to the normal PCR cycling protocol.
 4. Amplify for 20 to 30 cycles.
 5. For the final extension, incubate at 72°C for 10 minutes. Store at 4°C.
 6. Visualize by agarose gel electrophoresis.
-

Continued on next page

Analyzing *E. coli* Transformants, continued

Analyzing transformants by sequencing

Once you have identified the correct clone(s), you may sequence your construct to confirm that your gene is cloned in the correct orientation. Use a pair of primers that hybridize to the vector backbone flanking your insert to help you sequence your insert. For the complete sequence of the pSyn_1 Vector, refer to our website (www.lifetechnologies.com) or contact Technical Support (see page 24).

Long-term storage of *E. coli* clones

Once you have identified the correct clone, purify the colony and make a glycerol stock for long-term storage. We recommend that you store a stock of plasmid DNA at -20°C .

1. Streak the original colony out on LB plates containing 100 $\mu\text{g}/\text{mL}$ spectinomycin.
 2. Isolate a single colony, inoculate into 1–2 mL of LB containing 100 $\mu\text{g}/\text{mL}$ spectinomycin, and grow until culture reaches stationary phase.
 3. Mix 0.85 mL of culture with 0.15 mL of sterile glycerol and transfer to a cryovial, and store at -80°C .
-

Guidelines for Culturing *Synechococcus elongatus* PCC 7942

General guidelines for *Synechococcus* culture

- All solutions and equipment that may contact cells must be sterile. Always use proper aseptic technique and work in a laminar flow hood.
 - *Synechococcus elongatus* liquid cultures should be grown at $34^{\circ}\text{C} \pm 1^{\circ}\text{C}$ with CO_2 (1–2% in air) under continuous illumination using moderate light intensities of cool fluorescent white ($50\text{--}100 \mu\text{E m}^{-2} \text{s}^{-1}$) with agitation on a gyrotary shaker set to 100 rpm.
Note: *Synechococcus elongatus* cultures can also be grown using light intensities of $50\text{--}400 \mu\text{E m}^{-2} \text{s}^{-1}$ using only atmospheric CO_2 (i.e., without the addition of CO_2).
 - The presence of CO_2 is needed to obtain optimal growth of *Synechococcus elongatus* in liquid culture as it is a photosynthetic organism; however, additional CO_2 is not necessary during transformation where the cells are grown on agar support and are exposed to atmospheric CO_2 .
 - If you are bubbling the culture with CO_2 enriched air or CO_2 gas, you need to prepare the BG-11 with 50 mM NaHCO_3 and adjust the pH to 7.5. The presence of NaHCO_3 in the medium prevents the medium from becoming acidic.
 - *Synechococcus elongatus* solid cultures on BG-11 agar plates can be grown at 34°C under continuous illumination using $100\text{--}200 \mu\text{E m}^{-2} \text{s}^{-1}$ of cool fluorescent white light.
Note: You can also incubate the cultures at room temperature if sufficient light is provided; however, the growth will slower.
 - The optimal equipment for culturing *Synechococcus elongatus* is an algal growth chamber (e.g., Percival Algal Chamber from Geneva Scientific) with regulatable light supply and a light meter to guide adjustments. If an algal growth chamber is not available, the cells can be grown in a standard cell culture incubator illuminated with cool fluorescent lights placed within 12 inches of the culture plates. Standard room lights and incubation at room temperature provide sub-optimal growth conditions.
 - Do not stack the culture plates to allow continuous uniform illumination.
 - Grow the cells using Gibco® BG-11 medium, which is specifically formulated for optimal growth and maintenance of *Synechococcus elongatus* cells.
 - Grow the cells until the culture reaches OD_{750} of ≥ 1 before transformation.
 - Take the OD measurements at 750 nm.
 - *Synechococcus elongatus* is classified as a GRAS (generally regarded as safe) organism with no known viral or bacterial pathogens. However, we recommend following general safety guidelines under Biosafety Level 1 (BL-1) containment, similar to working with *E. coli* or yeast. For more information on BL-1 guidelines, refer to *Biosafety in Microbiological and Biomedical Laboratories*, 5th ed., published by the Centers for Disease Control, which is available for downloading at: www.cdc.gov/od/ohs/biosfty/bmb15/bmb15toc.htm.
-

Thawing *Synechococcus elongatus* PCC 7942

Guidelines for thawing *Synechococcus elongatus* PCC 7942

- Do not thaw more than 3 vials at a time.
 - Frozen *Synechococcus elongatus* cells are very sensitive to temperature fluctuations.
 - Before the cells are thawed, the cells must be transferred from the -80°C freezer into a dry ice container as quickly as possible and the vials should be buried in dry ice.
 - To maximize the recovery of the cells when thawing, warm the cells very quickly by placing the tubes directly from the dry ice container into a 35°C water bath. Once the cells are completely thawed, immediately dilute them into Gibco® BG-11 medium, pre-warmed to room temperature.
-

Materials needed

- 35°C water bath with a dark lid
Note: If a dark lid is not available, cover the water bath with aluminum foil.
 - Algal growth chamber (e.g., Percival Algal Chamber from Geneva Scientific) set to $34^{\circ}\text{C} \pm 1^{\circ}\text{C}$ and 1% CO_2 , optimal, under continuous illumination with $50 \mu\text{E m}^{-2} \text{s}^{-1}$
Note: If an algal growth chamber is not available, you can use a standard cell culture incubator under continuous illumination using moderate light intensities of cool fluorescent white ($50 \mu\text{E m}^{-2} \text{s}^{-1}$).
 - Gyrotory shaking platform set to 100 rpm
 - 6-well clear-bottom culture plates
 - Gibco® BG-11 medium (Cat. no. A1379901 or A1379902), pre-warmed to room temperature
 - 70% ethanol
 - Dry ice
-

Continued on next page

Thawing *Synechococcus elongatus* PCC 7942, continued

- Thawing procedure
1. Remove the frozen cells from -80°C storage and immediately place them in a dry ice container. Bury the vial(s) containing the cells in dry ice to minimize temperature fluctuations before thawing.
 2. Add 6 mL of Gibco[®] BG-11 medium, pre-warmed to room temperature, into each well of a 6-well plate.
 3. Remove the cryovial containing the frozen cells from the dry ice storage and **immediately** place it into a 35°C water bath.
 4. Quickly thaw the cells by placing the vial containing the cells in the 35°C water bath until the last ice crystal has melted (~ 2 minutes). Do not agitate the cells while thawing (i.e., do not swirl the vial).
 5. After the cells have thawed, wipe the outside of the vial with 70% ethanol, and place the vial in a rack at room temperature. Proceed immediately to the next step
 6. Transfer 100 μL of thawed cells from the vial into each well of the 6-well plate containing 6 mL of Gibco[®] BG-11 medium.
 7. Place the 6-well plate(s) in the algal growth chamber set to set to $34^{\circ}\text{C} \pm 1^{\circ}\text{C}$ with 1% CO_2 and illuminated with constant light of $50 \mu\text{E m}^{-2} \text{s}^{-1}$. Do not incubate under light intensity of $>50 \mu\text{E m}^{-2} \text{s}^{-1}$ because the cells are sensitive to light immediately after resuscitation. **Do not stack the plates.**
 8. Incubate the cells with gentle agitation on a gyrotary shaker set to 100 rpm.
 9. On Day 2, transfer 400 μL of the cell suspension and into a disposable plastic cuvette containing 400 μL of Gibco[®] BG-11 medium to measure the optical density.
 10. Measure the cell density at 750 nm (OD_{750}). If the OD_{750} is greater than 1, proceed to the transformation step (page 15). If the culture has not yet reached $\text{OD}_{750} = 1$, return it to the algal growth chamber and continue the incubation.
-

Transforming *Synechococcus elongatus* PCC 7942

Guidelines for transforming *Synechococcus elongatus* PCC 7942

- *Synechococcus elongatus* is naturally transformable with highest efficiencies of transformation when the culture is in the log phase of growth (OD_{750} of 1–2).
 - Transform *Synechococcus elongatus* using circular, supercoiled DNA.
 - Incubate the transformation reaction at 34°C, in the dark.
Note: Darkness increases the transformation efficiency.
 - The quality and the concentration of DNA used play a central role for the efficiency of transformation. Use a commercial kit such as the PureLink® HQ Mini Plasmid Purification or the PureLink® HiPure Plasmid Miniprep kits that deliver pure DNA (see page 22 for ordering information).
 - Pre-warm the selective BG-11 + spectinomycin plates to room temperature for 1 hour before plating the transformants.
-

Materials needed

- pSyn_1 construct carrying your gene of interest
 - pSyn_2/Control Vector
 - Gibco® BG-11 medium (Cat. no. A1379901 or A1379902), pre-warmed to room temperature
 - BG-11 agar plates containing 10 µg/mL spectinomycin, pre-warmed to room temperature (see page 18 for recipe)
Note: You will need 2 plates per transformation.
 - 34°C water bath with a dark lid or covered with aluminum foil
 - Algal Growth Chamber (e.g., Percival Algal Chamber from Geneva Scientific) set to 34°C ± 1°C and 1–2% CO₂, under continuous illumination with 100 µE m⁻² s⁻¹
Note: If an Algal Chamber is not available, you can use a standard cell culture incubator under continuous illumination using moderate light intensities of cool fluorescent white (100 µE m⁻² s⁻¹).
 - 70% ethanol
 - Sterile microcentrifuge tubes
 - Disposable spreaders
-

Continued on next page

Transforming *Synechococcus elongatus* PCC 7942, continued

Transformation procedure

1. Measure the optical density of the *Synechococcus elongatus* cultures (from step 10, page 14) at 750 nm (i.e., OD₇₅₀).
Note: For best performance, the OD₇₅₀ of cultures should be greater than 1 and less than 2.
 2. Harvest 1.5 mL of the cells (per transformation) by centrifugation at 14,000 rpm for 3 minutes at room temperature.
 3. Remove the supernatant by pipetting.
 4. Resuspend the cells in 1 mL of Gibco® BG-11 medium by gently pipetting up and down.
 5. Centrifuge the cells at 14,000 rpm for 1 minute at room temperature, and remove the supernatant by pipetting.
 6. Resuspend the cells in 100 µL of Gibco® BG-11 medium by gently pipetting up and down.
 7. Add 100 ng of supercoiled plasmid DNA (i.e., pSyn_1 construct containing your gene of interest) into the resuspended cells. Mix the DNA-cell suspension gently by flicking the tube. In a separate tube, prepare a control transformation with the pSyn_2/Control Vector.
 8. Incubate the cell-DNA mixture(s) in the 34°C water bath with a dark lid for 4 hours. After the incubation is complete, remove the tube(s) from the water bath and wipe them with 70% ethanol.
 9. Plate 80 µL and 5 µL of each transformation mixture on separate BG-11 agar plates containing 10 µg/mL spectinomycin and pre-warmed to room temperature.
 10. Place the plates with agar side down on illuminated shelves at room temperature (25–30°C). Do not stack the plates to ensure continuous and even illumination.
 11. Incubate the plates for 5–7 days or until the colonies are ready to pick. Control vector should produce a minimum of 100 transformants per transformation. The results from the transformation with the pSyn_1 construct will depend on the nature of your gene of interest.
-

Screening for Integration by Colony PCR

Introduction

Use the protocols below to prepare cell lysates and perform colony PCR to screen the transformed *Synechococcus elongatus* colonies for full integration of the promoter and the gene of interest. You will have to design the forward and reverse PCR primers specific to your insert and determine the amplification conditions. We recommend using the AccuPrime™ *Pfx* Polymerase SuperMix or the PCR SuperMix High Fidelity for best results; however, other DNA polymerases may also work.

Materials needed

- AccuPrime™ *Pfx* SuperMix (Cat. no. 12344-040) or PCR SuperMix High Fidelity (Cat. no. 10790-020)
 - Appropriate forward and reverse primers (10 μM each)
-

Colony PCR procedure

1. Streak or patch colonies onto fresh BG-11 agar plates containing 10 μg/mL of spectinomycin and allow them to grow for 1–2 days or until you have sufficient growth before proceeding with the colony PCR protocol below.
2. Prepare the PCR reaction as shown in table below.
3. Pick up cells with a pipette tip from the re-streaked plates (lift enough material that is equivalent to about 1 small colony, too much material will inhibit the PCR reaction) and resuspend in the PCR reaction mix.

Reagent	Amount
AccuPrime™ <i>Pfx</i> SuperMix or PCR SuperMix High Fidelity	45 μL
Primer pre-mix (10 μM each of forward and reverse primers)	1 μL
Colony	1

4. Mix the contents of the tubes and load into a thermal cycler.
 5. Heat at 95°C for 5 minutes for the initial denaturation step before proceeding with the normal PCR cycling protocol.
 6. Maintain reaction at 4°C after cycling. Samples can be stored at –20°C.
 7. Analyze the results by agarose gel electrophoresis.
-

Scale-up

We recommend initiating a seed culture first before scaling up your clones. To initiate a seed culture, scrape up as much of the streak or patch as possible (from step 1, **Colony PCR procedure**, above) and transfer it into each well of a 6-well plate containing 6 mL of Gibco® BG-11 medium with antibiotic, and grow them at 34°C ± 1°C with 1–2% CO₂ and illuminated with constant light of 50 μE m⁻² s⁻¹. Do not stack the plates.

Incubate the cells with gentle agitation on a gyrotary shaker set to 100 rpm until they reach the log phase of growth (OD₇₅₀ of 1–2). Use these seed cultures to initiate larger cultures in shake flasks for further investigation by diluting them to 1/20 (v/v). We recommend the addition of antibiotic (10 μg/mL spectinomycin) to the culture, especially for genes that are toxic to the cells.

Appendix A: Support Protocols

Media and Plates

LB (Luria-Bertani) medium and plates

LB medium:

1. For 1 liter, dissolve 10 g tryptone, 5 g yeast extract, and 10 g NaCl in 950 mL of deionized water.
2. Adjust the pH of the solution to 7.0 with NaOH and bring the volume up to 1 liter.
3. Autoclave the solution on liquid cycle for 20 minutes at 15 psi. Allow the solution to cool to 55°C and add the appropriate antibiotics, if needed.
Note: Use spectinomycin at a final concentration of 100 µg/mL.
4. Store the medium at room temperature or at 4°C.

LB agar plates

1. Prepare LB medium as above, but add 15 g/L agar before autoclaving.
 2. Autoclave the medium plus agar on liquid cycle for 20 minutes at 15 psi.
 3. After autoclaving, cool the medium to ~55°C, add the appropriate antibiotics, and pour into 10 cm plates.
 4. Let the agar harden, then invert the plates and store them at 4°C, in the dark.
-

Preparing BG-11 agar plates

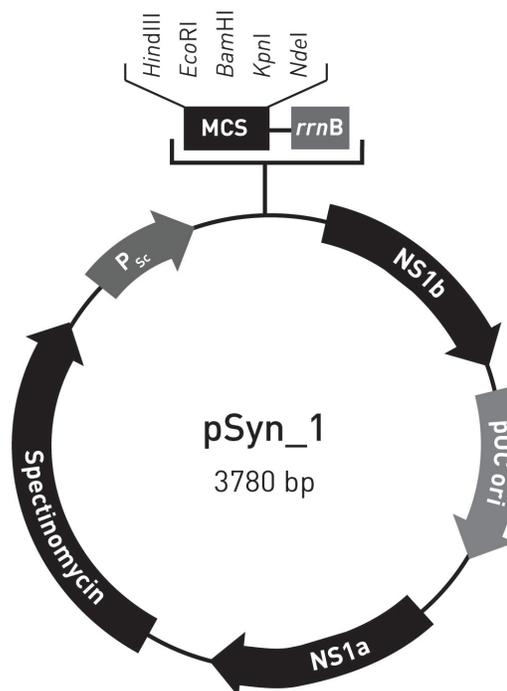
1. Add 15 g of agar to 200 mL of Gibco® BG-11 medium in an autoclaveable flask.
Note: Use high quality agar, such as Calbiochem, Cat no.12177, or Sigma, Cat. no. A1296.
 2. Autoclave on liquid cycle for 20 minutes.
 3. Warm 800 mL of Gibco® BG-11 medium to 55–60°C in a water bath.
 4. After autoclaving, cool the agar containing flask to ~55°C.
 5. Combine the agar containing flask with 800 mL of Gibco® BG-11 medium.
 6. Add spectinomycin to a final concentration of 10 µg/mL (if required), and pour into 10 cm plates.
 7. Let the plates harden (do **not** overdry), invert them, and store at 4°C in the dark. Final agar concentration will be 1.5%.
-

Appendix B: Vectors

Map and Features of pSyn_1 Vector

Map of pSyn_1 Vector

The map below shows the features of the pSyn_1 vector. The complete sequence of the vector is available for downloading at www.lifetechnologies.com or by contacting Technical Support (page 24).



Features of pSyn_1

3780 nucleotides

pUC origin:	58–673
NS1a (neutral site 1a):	745–1543
Spectinomycin resistance gene:	1550–2694
Promoter (P _{sc}):	2699–2799
Multiple cloning site (MCS):	2800–2834
<i>rrnB</i> transcriptional termination region:	2841–2998
NS1b (neutral site 1b):	2999–3780

Continued on next page

Map and Features of pSyn_1, continued

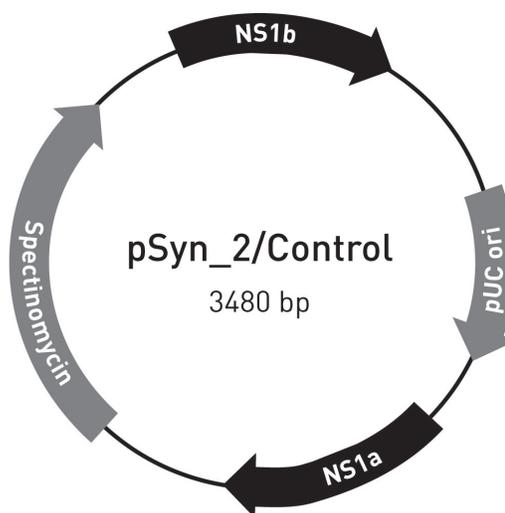
Features of pSyn_1 Vector The pSyn_1 Vector contains the following elements. All features have been functionally tested.

Feature	Benefit
pUC origin	Allows high-copy replication and growth in <i>E. coli</i> .
NS1a and NS1b (neutral site 1)	Sites also present on <i>Synechococcus elongatus</i> genome to guide double homologous recombination of DNA contained between the neutral sites in the vector (Clerico <i>et al.</i> , 2007).
Spectinomycin promoter (P_{Sp})	Allows expression of the spectinomycin resistance gene in <i>E. coli</i> and <i>Synechococcus elongatus</i> .
Spectinomycin resistance gene (<i>aadA1</i>)	Allows selection of the plasmid in <i>E. coli</i> and <i>Synechococcus elongatus</i> (Liebert <i>et al.</i> , 1999).
Promoter (P_{Sc})	Weak constitutive promoter from <i>Synechocystis</i> sp. strain PCC 6803 that allows a minimal level of expression of the GOI (Simkovsky <i>et al.</i> , 2012).
Multiple cloning site with 5 unique sites (<i>Hind</i> III, <i>Eco</i> RI, <i>Bam</i> HI, <i>Kpn</i> I, <i>Nde</i> I)	Allows insertion of your gene into pSyn_1.
<i>rrnB</i> transcription termination region	Strong transcription termination region.

Map of pSyn_2/Control Vector

pSyn_2/Control Vector

The map below shows the features of the pSyn_2/Control Vector. The control vector is used as a positive control for *Synechococcus elongatus* transformations and confers spectinomycin resistance to successfully transformed cells. When used as detailed on page 16, the control vector should produce a minimum of 100 transformants per transformation. The complete sequence of the vector is available for downloading at www.lifetechnologies.com or by contacting Technical Support (page 24).



Features of pSyn_2/Control 3480 nucleotides

pUC origin:	58–673
NS1a (neutral site 1a):	745–1543
Spectinomycin resistance gene:	1550–2694
NS1b (neutral site 1b):	2699–3480

Appendix C: Ordering Information

Accessory Products

Proofreading DNA polymerases

Life Technologies offers a variety of proofreading, thermostable DNA polymerases for generating blunt-end PCR products. Ordering information is provided below. For details, visit www.lifetechnologies.com.

Product	Quantity	Cat. no.
Platinum [®] Pfx DNA Polymerase	100 units	11708-013
AccuPrime [™] Pfx DNA Polymerase	200 reactions	12344-024
Pfx50 [™] DNA Polymerase	100 reactions	12355-012
PCR SuperMix High Fidelity	100 reactions	10790-020

Competent cells

Chemically competent and electrocompetent cells that can be used with GeneArt[®] *Synechococcus* Engineering Kits are also available separately from Life Technologies. Ordering information is provided below. For details, visit www.lifetechnologies.com.

Product	Quantity	Cat. no.
One Shot [®] TOP10 Chemically Competent Cells	10 reactions	C4040-10
	20 reactions	C4040-03
One Shot [®] TOP10 Electrocomp [™] <i>E. coli</i>	10 reactions	C4040-50
	20 reactions	C4040-52
TOP10 Electrocomp [™] Kits	20 reactions	C664-55
	40 reactions	C664-11
	120 reactions	C664-24

Additional products

The following reagents are recommended for use with the GeneArt[®] *Synechococcus* Engineering Kits. Ordering information for these reagents is provided below. For details, visit www.lifetechnologies.com.

Product	Quantity	Cat. no.
Gibco [®] BG-11 Media: Optimized for <i>Synechococcus</i>	1 L	A1379901
	6 × 1 L	A1379902
PureLink [®] Growth Block	50 blocks	12256-020
PureLink [®] HQ Mini Plasmid Purification Kit	100 preps	K2100-01
PureLink [®] HiPure Plasmid Miniprep Kit	25 preps	K2100-02
	100 preps	K2100-03
LB Broth (1X), liquid	500 mL	10855-021

Continued on next page

Accessory Products, continued

Other GeneArt® algal kits

In addition to the GeneArt® *Synechococcus* Engineering Kits, Life Technologies offers the following products as model algal hosts. Ordering information is provided below. For details, visit www.lifetechnologies.com.

Product	Quantity	Cat. no.
GeneArt® <i>Synechococcus</i> TOPO® Engineering Kit	1 kit	A14261
GeneArt® <i>Synechococcus</i> TOPO® Engineering Kit with 6 L media	1 kit	A14265
GeneArt® <i>Chlamydomonas</i> Engineering Kit	1 kit	A14258
GeneArt® <i>Chlamydomonas</i> Engineering Kit with 6 L media	1 kit	A14262
GeneArt® <i>Chlamydomonas</i> TOPO® Engineering Kit	1 kit	A14260
GeneArt® <i>Chlamydomonas</i> TOPO® Engineering Kit with 6 L media	1 kit	A14264

Documentation and Support

Obtaining Support

Technical Support

For the latest services and support information for all locations, go to www.lifetechnologies.com.

At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
 - Search through frequently asked questions (FAQs)
 - Submit a question directly to Technical Support (techsupport@lifetech.com)
 - Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents
 - Obtain information about customer training
 - Download software updates and patches
-

Safety Data Sheets (SDS)

Safety Data Sheets (SDSs) are available at www.lifetechnologies.com/sds.

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 www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

References

- Atsumi, S., Higashide, W., and Liao, J. C. (2009) Direct photosynthetic recycling of carbon dioxide to isobutyraldehyde. *Nat Biotechnol* 27, 1177-1180
- Ausubel, F. M., Brent, R., Kingston, R. E., Moore, D. D., Seidman, J. G., Smith, J. A., and Struhl, K. (1994) *Current Protocols in Molecular Biology*, Greene Publishing Associates and Wiley-Interscience, New York
- Clerico, E. M., Ditty, J. L., and Golden, S. S. (2007) Specialized techniques for site-directed mutagenesis in cyanobacteria. *Methods Mol Biol* 362, 155-171
- Ducat, D. C., Sachdeva, G., and Silver, P. A. (2011) Rewiring hydrogenase-dependent redox circuits in cyanobacteria. *Proc Natl Acad Sci U S A* 108, 3941-3946
- Kaneko, T., Sato, S., Kotani, H., Tanaka, A., Asamizu, E., Nakamura, Y., Miyajima, N., Hirose, M., Sugiura, M., Sasamoto, S., Kimura, T., Hosouchi, T., Matsuno, A., Muraki, A., Nakazaki, N., Naruo, K., Okumura, S., Shimpo, S., Takeuchi, C., Wada, T., Watanabe, A., Yamada, M., Yasuda, M., and Tabata, S. (1996) Sequence analysis of the genome of the unicellular cyanobacterium *Synechocystis* sp. strain PCC6803. II. Sequence determination of the entire genome and assignment of potential protein-coding regions. *DNA Res* 3, 109-136
- Lan, E. I., and Liao, J. C. (2011) Metabolic engineering of cyanobacteria for 1-butanol production from carbon dioxide. *Metab Eng* 13, 353-363
- Liebert, C. A., Watson, A. L., and Summers, A. O. (1999) Transposon Tn21, Flagship of the Floating Genome. *Microbiol. Mol. Biol. Rev.* 63, 507-522
- Min, H., and Golden, S. S. (2000) A new circadian class 2 gene, *opcA*, whose product is important for reductant production at night in *Synechococcus elongatus* PCC 7942. *J Bacteriol* 182, 6214-6221
- Sambrook, J., Fritsch, E. F., and Maniatis, T. (1989) *Molecular Cloning: A Laboratory Manual*, Second Ed., Cold Spring Harbor Laboratory Press, Plainview, New York
- Simkovsky, R., Daniels, E. F., Tang, K., Huynh, S. C., Golden, S. S., and Brahmsha, B. (2012) Impairment of O-antigen production confers resistance to grazing in a model amoeba-cyanobacterium predator-prey system. *Proc Natl Acad Sci U S A* 109, 16678-16683
- Taniguchi, Y., Nishikawa, T., Kondo, T., and Oyama, T. (2012) Overexpression of *lalA*, a paralog of *labA*, is capable of affecting both circadian gene expression and cell growth in the cyanobacterium *Synechococcus elongatus* PCC 7942. *FEBS Lett* 586, 753-759
- Vijayan, V., Jain, I. H., and O'Shea, E. K. (2011) A high resolution map of a cyanobacterial transcriptome. *Genome Biol* 12, R47
-

Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA | Phone +1 760 603 7200 | Toll Free in USA 800 955 6288

For support visit lifetechnologies.com/support or email techsupport@lifetech.com

lifetechnologies.com

18 March 2013

