

**BL21-AI<sup>™</sup> One Shot<sup>®</sup>**  
**Chemically Competent**  
***E. coli***

Catalog nos. C6070-03

**Rev. Date: 7 June 2010**  
Manual part no. 25-0468

MAN0000259



# Table of Contents

About the Kit..... 1

Basic Transformation Procedure..... 4

Expression Guidelines..... 6

Testing Transformation Efficiency ..... 10

Regulation by L-Arabinose..... 11

Technical Support..... 12

Purchaser Notification ..... 13

References ..... 14



# About the Kit

## Shipping/ Storage

The BL21-AI™ One Shot® Chemically Competent *E. coli* Kit is shipped on dry ice. Upon receipt, store at -80°C.

## Kit Contents

The table below describes the items included in the BL21-AI™ One Shot® Chemically Competent *E. coli* Kit. The transformation efficiency of BL21-AI™ cells is  $1 \times 10^8$  cfu/μg of pUC18 plasmid DNA.

**Store at -80°C.**

Item	Composition	Amount
SOC Medium (store at room temperature or 4°C)	2% Tryptone, 0.5% Yeast Extract, 10 mM NaCl, 2.5 mM KCl, 10 mM MgCl <sub>2</sub> , 10 mM MgSO <sub>4</sub> , 20 mM glucose	6 mL
20% L-arabinose	20% L-arabinose in sterile water	1 mL
BL21-AI™ chemically competent cells	—	21 × 50 μL
pUC19 Control DNA	10 pg/μL in 5 mM Tris-HCl, 0.5 mM EDTA, pH 8	50 μL

## Genotype

**BL21-AI™:** F<sup>-</sup> *ompT hsdS<sub>B</sub> (r<sub>B</sub><sup>-</sup>m<sub>B</sub><sup>-</sup>) gal dcm araB::T7RNAP-tetA*

The BL21-AI™ strain is an *E. coli* B/r strain and does not contain the *lon* protease. It is also deficient in the outer membrane protease, OmpT. The lack of these proteases reduces degradation of heterologous proteins expressed in this strain.

The strain carries a chromosomal insertion of a cassette containing the T7 RNA polymerase (T7 RNAP) gene in the *araB* locus, allowing expression of T7 RNAP to be regulated by the *araBAD* promoter (see page 1 for more information).

The presence of the *tetA* gene confers resistance to tetracycline and permits verification of strain identity using tetracycline.

*Continued on next page*

# About the Kit, Continued

---

## The BL21-AI™ Strain

The BL21-AI™ *E. coli* strain is derived from the BL21 strain (Grodberg and Dunn, 1988; Studier and Moffatt, 1986) and is intended for use as an expression host. The BL21-AI™ strain contains a chromosomal insertion of the gene encoding T7 RNA polymerase (T7 RNAP) into the *araB* locus of the *araBAD* operon, placing regulation of T7 RNAP under the control of the *araBAD* promoter. The *araB* gene is deleted in this strain.

---

## Regulating Expression of T7 RNA Polymerase

Because the T7 RNAP gene is inserted into the *araB* locus of the *araBAD* operon, expression of T7 RNA polymerase can be regulated by the sugars, L-arabinose and glucose.

- To induce expression from the *araBAD* promoter, use L-arabinose (Lee, 1980; Lee *et al.*, 1987). To modulate expression, simply vary the concentration of L-arabinose added.
- To repress basal expression from the *araBAD* promoter, use glucose.

**Note:** In the absence of glucose, basal expression from the *araBAD* promoter is generally low (Lee, 1980; Lee *et al.*, 1987). Adding glucose further represses expression from the *araBAD* promoter by reducing the levels of 3', 5'-cyclic AMP (Miyada *et al.*, 1984).

For more information on the mechanism of expression and repression of the *ara* regulon, see page 11 or refer to Schleif, 1992.

---

## Expression of Heterologous Genes

**The BL21-AI™ strain is suitable for high-level recombinant protein expression from any T7-based expression vector.**

Because T7 RNA polymerase levels can be tightly regulated, the BL21-AI™ strain is especially useful to express genes that may be toxic to other BL21 strains where basal expression of T7 RNA polymerase is leakier (*e.g.* BL21 Star™(DE3) or BL21(DE3)). The yield of recombinant protein obtained from BL21-AI™ is generally similar to that obtained from other BL21 strains.

---

*Continued on next page*

# About the Kit, Continued

---

## **When to Use BL21-AI™**

In general, we recommend using the BL21-AI™ strain to express your gene of interest if:

- You are using a T7-based expression vector (either high-copy or low-copy)
- You observe growth inhibitory effects (i.e. toxicity) when using other BL21 strains (see page 8 for more information about indications of toxicity in bacterial cells)

You are expressing a known toxic gene

---

## **Intended Use**

**For research use only.** Not intended for human or animal diagnostic or therapeutic uses.

---

# Basic Transformation Procedure

---

## Introduction

A basic transformation protocol for BL21-AI™ cells is provided below. Once you have selected transformants, we recommend proceeding directly to expression using your own protocol. Note that the BL21-AI™ strain is designed for expression, not cloning or subcloning.

---

## Materials Supplied by the User

- Plasmid DNA (ready for transformation)
  - 42°C water bath
  - 37°C shaking and non-shaking incubator
  - Ice bucket with ice
  - Spectrophotometer to measure optical density of the cell cultures
  - Microcentrifuge tube rack (optional)
- 

## Before Starting

- Prepare LB agar plates containing the appropriate concentration of antibiotic for plasmid selection
  - Equilibrate a water bath to 42°C
  - Warm the vial of SOC medium to room temperature
  - Place the plates in a 37°C incubator to remove excess moisture (use two plates for each transformation)
- 

*Continued on next page*

# Basic Transformation Procedure, Continued

---

## Basic Transformation Procedure

1. Thaw one vial of BL21-AI™ One Shot® cells on ice per transformation.
2. Add 5–10 ng of DNA, in a volume of 1–5  $\mu\text{L}$  to the cells and mix by tapping gently. **Do not mix cells by pipetting.**
3. Incubate the vial(s) on ice for 30 minutes.
4. Heat shock the cells by incubating the vial(s) for exactly 30 seconds in the 42°C water bath. **Do not mix or shake.**
5. Remove the vial(s) from the 42°C bath and quickly place on ice.
6. Add **250  $\mu\text{L}$**  of pre-warmed SOC medium to the vial(s). (SOC is a rich medium; use proper sterile technique to avoid contamination.)
7. Secure the vial(s) in a microcentrifuge rack with tape. Place the rack in a shaking incubator, and shake the vial(s) at 37°C for 1 hour at 225 rpm.
8. Plate two different volumes of the transformation reaction onto LB plates containing the appropriate antibiotic for plasmid selection. Include 34  $\mu\text{g}/\text{mL}$  chloramphenicol if using BL21(DE3)pLysS or BL21(DE3)pLysE cells. Select two volumes ranging from 20–200  $\mu\text{L}$  to ensure well-spaced colonies on at least one plate. The remaining transformation reaction may be stored at 4°C and plated out the next day, if needed.
9. Invert the plates and incubate at 37°C overnight.
10. Select transformants from the plates and culture as described on page 7.

**Note:** Clones may exhibit differences in expression of heterologous genes. We recommend choosing 3–4 transformants when characterizing clones for protein expression.

---

# Expression Guidelines

---

## Introduction

If you have an expression protocol for the plasmid that you are working with, we recommend that you use your own protocol. This section provides some general guidelines for the use of T7 RNA polymerase-based expression plasmids in BL21-AI™ cells.

---



Use BL21-AI™ cells **for expression only**. To maintain your construct, transform your expression plasmid into a strain that does not bear the gene for T7 RNA polymerase (i.e. TOP10, DH5α™).

---

## Modulating Gene Expression

To induce expression of your gene of interest in BL21-AI™ cells, you will need the following reagents:

- L-arabinose to induce expression of T7 RNA polymerase. L-arabinose is supplied with the kit, but is also available from Sigma (Catalog no. A3256).
- Isopropyl β-D-thiogalactoside (IPTG; only required for T7-based expression vectors containing the *lacI* gene). IPTG is available from Invitrogen (Catalog no. 15529-019).

**Note:** You may repress **basal** transcription of T7 RNA polymerase and thereby, your gene of interest by adding glucose to the plates and/or media (see pages 8 and 11 for more information). By varying the concentration of L-arabinose and glucose, you may modulate expression of your gene of interest.

---

## Materials Needed

You should have the following stock solutions on hand before beginning. Prepare all solutions in sterile, deionized water.

- 20% L-arabinose (supplied with the kit)
  - 1 M IPTG (if needed)
  - 20% glucose (if needed)
- 

*Continued on next page*

# Expression Guidelines, Continued

---

## Expression Guidelines

General guidelines are provided below to express your gene of interest in BL21-AI™.

- Pick 3–4 transformants for overnight culture in 5 mL LB medium containing antibiotic to select for your expression plasmid. Grow overnight at 37°C with shaking until the OD<sub>600</sub> reaches 0.6–1.0.
- Use the overnight cultures to inoculate fresh LB medium containing antibiotic to an OD<sub>600</sub> of 0.05–0.1 (~1:20 dilution of the overnight culture). This dilution allows the cells to quickly return to logarithmic growth and reach the appropriate cell density. Use a volume appropriate for taking time points, if desired.
- Use the remainder of each overnight culture to create glycerol stocks. Once you have identified the clone that best expresses your protein, you can use the glycerol stock to perform additional expression experiments.
- Grow the cultures until they reach mid-log phase (OD<sub>600</sub> ~0.4; 2 to 3 hours).
- Induce the cultures (see below), and culture for an additional 2–3 hours. You may also take time points to analyze for optimal expression of your protein.

**For T7 expression vector containing the *lacI* gene** (e.g. Invitrogen's pET vectors), induce by adding L-arabinose to a final concentration of 0.2% **AND** IPTG to a final concentration of 1 mM.

**For T7 expression vector with no *lacI* gene** (e.g. Invitrogen's pCR®T7 vectors), induce by adding L-arabinose to a final concentration of 0.2%. Culture for an additional 2–3 hours.

- Analyze clones by western blot or enzymatic assay to determine which clone best expresses your protein of interest. Use the glycerol stock created from this clone for expression experiments. If you find that expression levels in subsequent inductions decrease, or you find that you lose your plasmid, your protein may be toxic to *E. coli* (see page 8 for additional information).

---

*Continued on next page*

# Expression Guidelines, Continued

---

## Indications of Toxicity

When expressing recombinant proteins in BL21-AI™ strains, one can generally assume that the recombinant protein is toxic to bacterial cells when any of the following occurs:

- No transformants are obtained after following Steps 1-9 of the **Basic Transformation Protocol** on page 5 **OR** a combination of large and small, irregular colonies appears on the plate
- The initial culture does not grow (see previous page)
- It takes longer than 5 hours after a 1:20 dilution of the initial culture for the fresh culture to reach an  $OD_{600}=0.4$  (see previous page)
- The cells lyse after induction with L-arabinose (or L-arabinose and IPTG) (see previous page)

---

## Precautions

Review the guidelines below when basal level expression of a gene of interest is toxic. These guidelines assume that the T7 expression plasmid has been correctly designed and created.

- Propagate and maintain your expression plasmid in a strain that does not contain T7 RNA polymerase (i.e. TOP10, DH5 $\alpha$ ™, etc.).
- Perform a fresh transformation of BL21-AI™ cells before each induction experiment.
- Minimize the amount of time that the cells bearing the gene of interest are cultured before induction.
- Supplement LB plates in **Basic Transformation Protocol** (page 5 Step 7), and growth medium for expression experiments (page 7) with 0.1% glucose. The presence of glucose represses basal expression of T7 RNA polymerase.
- Following transformation of BL21-AI™ cells, grow cells in SOC medium for 1 hour and go directly to protein expression. Do not plate the transformation mixture to select for individual clones. See next page for details.

---

*Continued on next page*

# Expression Guidelines, Continued

---

## Transformation/ Expression Protocol for Toxic Genes

This protocol is used with BL21-AI™ cells. Please note that other protocols are possible, depending on your needs.

### Transformation

1. Follow the basic transformation protocol on page 5 through Step 7.
2. After growing the transformation reaction in SOC for 1 hour (page 5, Step 7), add the **entire** transformation reaction (300  $\mu$ L) to 50–200 mL of LB medium pre-warmed to 37°C containing the appropriate selective antibiotic for your expression plasmid (and 0.1% glucose, if desired).

### Induction

3. Incubate the vial(s) with shaking at 37°C until the cells reach mid-log phase ( $OD_{600} = 0.4$ ). **Note:** Doubling times may vary (30 to 90 minutes) depending on the protein expressed.
  4. Induce expression of the recombinant protein by adding L-arabinose to a final concentration of 0.2% (and IPTG to a final concentration of 1 mM, if needed) and grow for 2–3 more hours. You may take time points, if desired.
  5. Harvest cells by centrifugation and use immediately for analysis, or store the cell pellet at –80°C.
-

# Testing Transformation Efficiency

---

## Introduction

To test the transformation efficiency of the competent cells contained in the One Shot<sup>®</sup> kit, use the supercoiled pUC19 plasmid supplied with the kit as described below. An extra vial of cells is included for this purpose.

---

## Before Starting

- Prepare LB agar plates containing 50 µg/mL ampicillin.
  - Equilibrate a water bath to 42°C.
  - Warm the vial of SOC medium to room temperature.
  - Place the plates in a 37°C incubator to remove excess moisture (use two plates for each transformation).
- 

## Transformation

Follow the transformation protocol on page 5 to transform pUC19 into BL21-AI<sup>™</sup>. Use the specific modifications below.

- Transform cells with 1 µL (10 pg) of pUC19
  - Plate 50 µL each onto two LB plates containing 50 µg/mL ampicillin
  - Calculate the transformation efficiency as transformants per µg of plasmid (see below). The cells should have an efficiency of  $1 \times 10^8$  transformants/µg of supercoiled plasmid
- 

## Calculation

Use the formula below to calculate transformation efficiency.

$$\frac{\# \text{ of colonies}}{10 \text{ pg transformed DNA}} \times \frac{10^6 \text{ pg}}{\mu\text{g}} \times \frac{300 \text{ }\mu\text{L transformed cells}}{X \text{ }\mu\text{L plated}} = \frac{\# \text{ transformants}}{\mu\text{g plasmid DNA}}$$

---

# Regulation by L-Arabinose

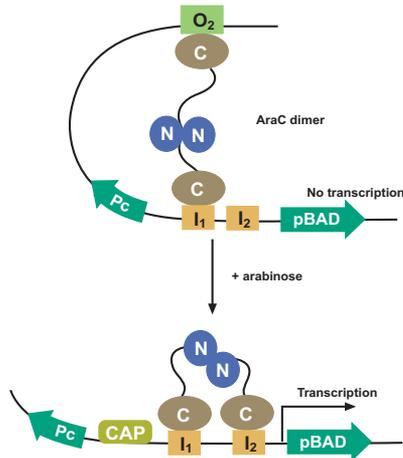
## Introduction

The L-arabinose regulatory circuit is briefly described below.

## Regulation of the *araBAD* ( $P_{BAD}$ ) Promoter

The *araBAD* promoter ( $P_{BAD}$ ) used to control expression of T7 RNA polymerase in BL21-AI™ is both positively and negatively regulated by the product of the *araC* gene (Ogden *et al.*, 1980; Schleif, 1992). AraC is a transcriptional regulator that forms a complex with L-arabinose. In the absence of L-arabinose the AraC dimer contacts the O<sub>2</sub> and I<sub>1</sub> half sites of the *araBAD* operon, forming a 210 bp DNA loop (see figure below). For maximum transcriptional activation two events are required.

- L-Arabinose binds to AraC and causes the protein to release the O<sub>2</sub> site and bind the I<sub>2</sub> site which is adjacent to the I<sub>1</sub> site. This releases the DNA loop and allows transcription to begin.
- The cAMP activator protein (CAP)-cAMP complex binds to the DNA and stimulates binding of AraC to I<sub>1</sub> and I<sub>2</sub>.



## Glucose Repression

Basal expression levels can be repressed by adding glucose to the growth medium. Glucose lowers cAMP levels, which in turn decreases the binding of CAP. As cAMP levels are lowered, transcriptional activation is decreased.

# Technical Support

---

## Web Resources



Visit the Invitrogen website at [www.invitrogen.com](http://www.invitrogen.com) for:

- Technical resources including manuals, vector maps and sequences, application notes, SDSs, FAQs, formulations, citations, handbooks, etc.
  - Complete technical support contact information.
  - Access to the Invitrogen Online Catalog.
  - Additional product information and special offers.
- 

## Contact Us

For more information or technical assistance, call, write, fax, or email. Additional international offices are listed on [www.invitrogen.com](http://www.invitrogen.com).

### Corporate Headquarters:

5791 Van Allen Way  
Carlsbad, CA 92008 USA  
Tel: 1 760 603 7200  
Tel (Toll Free): 1 800 955 6288  
Fax: 1 760 602 6500  
E-mail: [tech\\_support@invitrogen.cc](mailto:tech_support@invitrogen.cc)

### Japanese Headquarters:

LOOP-X Bldg. 6F  
3-9-15, Kaigan  
Minato-ku, Tokyo 108-0022  
Tel: 81 3 5730 6509  
Fax: 81 3 5730 6519  
E-mail: [jpinfo@invitrogen.cc](mailto:jpinfo@invitrogen.cc)

### European Headquarters:

Inchinnan Business Park  
3 Fountain Drive  
Paisley PA4 9RF, UK  
Tel: +44 (0) 141 814 6100  
Tech Fax: +44 (0) 141 814 6117  
E-mail: [eurotech@invitrogen.com](mailto:eurotech@invitrogen.com)

---

## SDS and Certificate of Analysis

Safety Data Sheets (SDSs) are available at [www.invitrogen.com/sds](http://www.invitrogen.com/sds). The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to [www.invitrogen.com/support](http://www.invitrogen.com/support) and search for the Certificate of Analysis by product lot number, which is printed on the box.

---

# Purchaser Notification

---

## Information for European Customers

The BL21-AI™ *E. coli* strain is genetically modified and carries a chromosomal insertion of a cassette containing the T7 RNA polymerase (T7 RNAP) gene. 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 Use Label License No: 30 T7 Expression System

The composition and/or use of this product may be claimed in U.S. Patent No. 5,693,489 licensed to Life Technologies Corporation by Brookhaven Science Associates, LLC. The T7 expression system is based on technology developed at Brookhaven National Laboratory under contract with the U.S. Department of Energy, and is the subject of patents and patent applications assigned to Brookhaven Science Associates, LLC (BSA.). By provisions of the Distribution License Agreement granted to Invitrogen covering said patents and patent applications, Invitrogen grants you a non-exclusive sub-license under patents assigned to BSA for the use of this technology, including the enclosed materials, based upon the following conditions: 1 – these materials are to be used for non-commercial research purposes only. A separate license under patents owned by BSA is required for any commercial use, including the use of these materials for research purposes or production purposes by any commercial entity. Information about commercial license may be obtained from The Office of Technology Transfer, Brookhaven National Laboratory, Bldg. 475D, P.O. Box 5000, Upton, New York 11973-5000. Phone (516) 344-7134. 2 - No materials that contain the cloned copy of the T7 gene 1, the gene for T7 RNA polymerase, may be distributed further to third parties outside of your laboratory, unless the recipient receives a copy of this sub-license and agrees to be bound by its terms. This limitation applies to strains BL21(DE3), BL21(DE3)pLysS and BL21(DE3)pLysE, CE6, BL21-SI Competent Cells and any derivatives that are made of them. You may refuse this sub-license by returning this product unused in which case Invitrogen accept return of the product with a full refund. By keeping or using this product, you agree to be bound by the terms of this license.

---

## Limited Warranty

Invitrogen (a part of Life Technologies Corporation) is committed to providing our customers with high-quality goods and services. Our goal is to ensure that every customer is 100% satisfied with our products and our service. If you should have any questions or concerns about an Invitrogen product or service, contact our Technical Support Representatives. All Invitrogen products are warranted to perform according to specifications stated on the certificate of analysis. The Company will replace, free of charge, any product that does not meet those specifications. This warranty limits the Company's liability to only the price of the product. No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. The Company reserves the right to select the method(s) used to analyze a product unless the Company agrees to a specified method in writing prior to acceptance of the order. Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore the Company makes no warranty of any kind regarding the contents of any publications or documentation. If you discover an error in any of our publications, report it to our Technical Support Representatives. **Life Technologies Corporation shall have no responsibility or liability for any special, incidental, indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive. No other warranty is made, whether expressed or implied, including any warranty of merchantability or fitness for a particular purpose.**

---

## References

---

- Grodberg, J., and Dunn, J. J. (1988). *ompT* Encodes the *Escherichia coli* Outer Membrane Protease that Cleaves T7 RNA Polymerase During Purification. *J. Bacteriol.* *170*, 1245-1253.
- Lee, N. (1980) Molecular Aspects of *ara* Regulation. In *The Operon*, J. H. Miller and W. S. Reznikoff, eds. (Cold Spring Harbor, N.Y.: Cold Spring Harbor Laboratory), pp. 389-410.
- Lee, N., Francklyn, C., and Hamilton, E. P. (1987). Arabinose-Induced Binding of AraC Protein to *araI*<sub>2</sub> Activates the *araBAD* Operon Promoter. *Proc. Natl. Acad. Sci. USA* *84*, 8814-8818.
- Miyada, C. G., Stoltzfus, L., and Wilcox, G. (1984). Regulation of the *araC* Gene of *Escherichia coli*: Catabolite Repression, Autoregulation, and Effect on *araBAD* Expression. *Proc. Natl. Acad. Sci. USA* *81*, 4120-4124.
- Ogden, S., Haggerty, D., Stoner, C. M., Kolodrubetz, D., and Schleif, R. (1980). The *Escherichia coli* L-Arabinose Operon: Binding Sites of the Regulatory Proteins and a Mechanism of Positive and Negative Regulation. *Proc. Natl. Acad. Sci. USA* *77*, 3346-3350.
- Schleif, R. S. (1992). DNA Looping. *Ann. Rev. Biochem.* *61*, 199-223.
- Studier, F. W., and Moffatt, B. A. (1986). Use of Bacteriophage T7 RNA Polymerase to Direct Selective High-Level Expression of Cloned Genes. *J. Mol. Biol.* *189*, 113-130.

---

©2010 Life Technologies Corporation. All rights reserved.

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.





**Corporate Headquarters**

5791 Van Allen Way  
Carlsbad, CA 92008

T: 1 760 603 7200

F: 1 760 602 6500

E: [tech\\_support@invitrogen.com](mailto:tech_support@invitrogen.com)

For country-specific contact information visit our web site at [www.invitrogen.com](http://www.invitrogen.com)