

# pGLO<sup>™</sup> Transformation and Inquiry Kit

A ThINQ!<sup>™</sup> Investigation

**Student Manual**

**BIO-RAD**

## Dear Students

Mice that glow fluorescent green. Plants that turn red when grown near a land mine. Goats that make milk that can be spun into parachute fabric. Virus-resistant papayas. Cheese puff snacks. Insulin. What do all these things have in common? They are all either genetically modified organisms (GMOs), or produced by GMOs.

GMO stands for genetically modified organism, an organism that has undergone genetic transformation (literally “change caused by genes”). Whether you call it genetic modification, genetic transformation, or genetic engineering, it’s all one and the same — these terms all refer to the process of manipulating the genes of an organism to cause a change to traits of that organism.

Genetic engineering and GMOs are currently debated topics. Opponents of the technology claim, among other things, potential harm to the health of humans and the environment. They contend that the technology needs more testing and much more regulation before it is deemed safe. Meanwhile, supporters of the technology tout its many useful applications: increasing the food supply to accommodate a growing human population, producing medicines and other products **more efficiently, treating or eradicating disease, to name a few.** They assert the technology has no demonstrated ill effects on humans, and its utility cannot be ignored.

This instruction manual, and the experiments outlined within it, are designed to help you better understand genetic engineering by creating a GMO of your own — a bacterium that, like the mouse mentioned above, will glow bright green under ultraviolet (UV) light.

The experiments take you through the process of bacterial genetic transformation via a standard protocol used in molecular and cell biology laboratories worldwide. You will begin with a plate of bacteria that appear off-white under normal or UV light conditions, and a tube of DNA containing a gene that codes for green **fluorescent protein (GFP)**. **The source of this gene is the bioluminescent jellyfish *Aequorea victoria*.** You will learn all the steps needed to move the DNA into the bacteria, to have the **bacteria incorporate the DNA, and to induce the bacteria to produce the fluorescent protein**, which causes them to glow a brilliant bright green under UV light. You will then have the opportunity to expand and explore the process further by designing and performing your own experiments to explore transformation, antibiotic dosage, gene expression, or satellite colony formation. Finally, you’ll end the hands-on experience with a real-world science case study to understand how genetic engineering and bacterial transformation might one day play a major **role in the fight against malaria.**

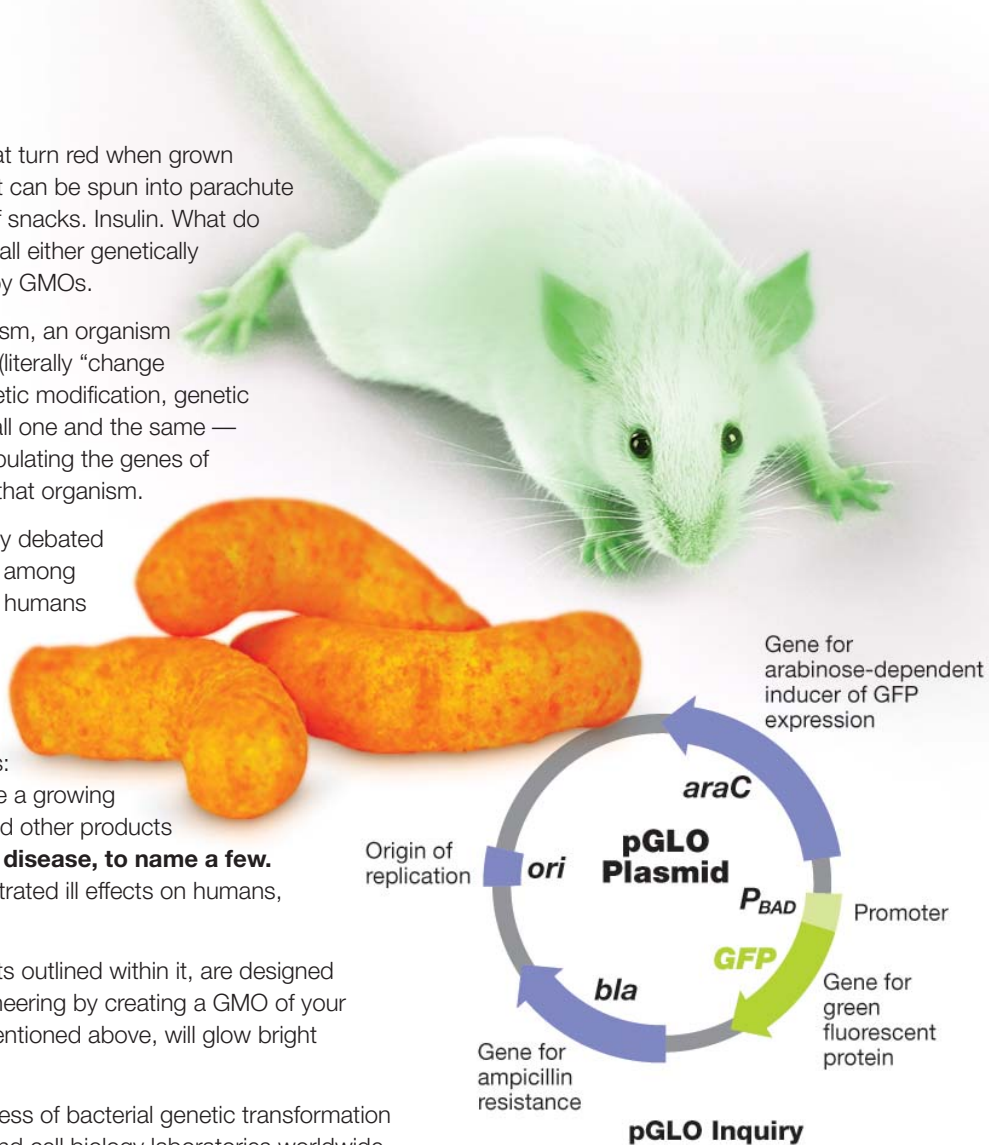
Throughout these experiments, we hope you ask a lot of questions so that you can widen your understanding of genetic engineering — its promises, its limitations, and the types of careful considerations that should be made before creating or using GMOs for any application. We hope the knowledge you gain in the following investigations will help you develop your own opinions on the complex ethical and practical debates surrounding GMOs and genetic engineering.

Bio-Rad’s Explorer Team

**Bio-Rad Laboratories**

6000 James Watson Drive, Hercules, CA 94547

***Biotechnology\_Explorer@bio-rad.com***



# Contents

<b>I. Introduction and Background .....</b>	<b>1</b>
Introduction to Bacterial Transformation.....	1
<b>II. Investigation #1: pGLO Bacterial Transformation Laboratory (Structured Inquiry).....</b>	<b>4</b>
Workstation Checklist .....	4
Protocol.....	5
Data Collection and Analysis of Results .....	9
The Interaction between Genes and Environment .....	11
Calculation of Transformation Efficiency .....	12
<b>III. Investigations #2–5: Guided and Open Inquiry .....</b>	<b>16</b>
Investigation #2: Transformation Efficiency Lab (Guided Inquiry).....	17
Investigation #3: Effect of Ampicillin on Bacterial Growth Lab (Open Inquiry) .....	22
Investigation #4: Effect of Arabinose on GFP Expression Lab (Open Inquiry) .....	27
Investigation #5: Satellite Colonies Lab (Open Inquiry).....	33
Post-Lab Assessment .....	38
<b>IV. Class Discussion.....</b>	<b>40</b>
Science Case Study: Can Bacterial Transformation Stop the Spread of Malaria? .....	40
<b>Appendix A: Serial Dilution and Dilution Instructions.....</b>	<b>48</b>
<b>Appendix B: Gene Regulation .....</b>	<b>52</b>

# I. Introduction to Bacterial Transformation

## Background

In this lab, you will perform a genetic transformation, which literally means “change caused by genes.” A gene is a piece of DNA that provides the instructions for making (it codes for, or encodes) a protein. This protein usually serves some cellular function and gives an organism a particular trait, also known as a phenotype. Genetic transformation involves inserting a gene into an organism in order to change a trait of that organism.

Genetic transformation is used in many areas of biotechnology. In the lab, bacterial transformation is a common method for producing more copies of genes (cloning) for insertion of that gene into a variety of model organisms. Expressing genes in model organisms can help determine the function of those genes in more complex organisms. In agriculture, genes coding for traits such as frost, pest, or spoilage resistance can be genetically transformed into plants. In bioremediation, bacteria can be transformed with genes that enable them to digest oil spills. In medicine, diseases caused by defective genes are beginning to be treated by gene therapy; that is, by genetically transforming a sick person’s cells with healthy copies of the defective gene that causes the disease.

Genes can be obtained from human, animal, or plant DNA and placed inside bacteria for a variety of purposes. One purpose is to use bacterial cells as microscopic factories, either to clone more copies of DNA or to produce a gene product. For example, the healthy human gene for the hormone insulin can be put into bacteria. Under the right conditions, these bacteria can make authentic human insulin, which can then be used to treat patients with diabetes, in whom insulin-producing cells do not function normally.

You will transform bacteria with a gene that codes for green fluorescent protein (GFP). The source of this gene is the bioluminescent jellyfish *Aequorea victoria*. When exposed to ultraviolet (UV) light, the GFP protein causes the jellyfish to fluoresce and glow green. Following the transformation procedure, your bacteria will express their newly acquired jellyfish gene and produce the fluorescent protein, which will cause them to glow a brilliant green under UV light.

In this activity, you will learn about the process of moving genes from one organism to another with the aid of a **plasmid**. In addition to one large chromosome, bacteria naturally contain one or more small circular pieces of DNA called plasmids. Plasmid DNA usually contains genes for one or more traits that may be beneficial to bacterial survival. In nature, bacteria can transfer plasmids back and forth among themselves, allowing them to share genes, which may permit them to adapt to new environments. The development of bacterial resistance to antibiotics is often due to such transmission of plasmids.

Bio-Rad’s unique pGLO plasmid contains several elements critical to its function:

- *Aequorea victoria* GFP gene, which confers green fluorescence in the presence of UV light when the gene is expressed
- *bla* gene, whose gene product confers ampicillin resistance to bacteria when it is expressed
- *araC* gene, a component of the arabinose operon whose encoded protein stimulates GFP gene expression from the  $P_{BAD}$  promoter when **arabinose** is present

Following bacterial transformation, selection for cells that have been transformed with pGLO DNA is accomplished by growth on **ampicillin** containing plates. When pGLO transformed bacterial cells are grown in nutrient medium containing **arabinose** as well, GFP expression is stimulated and the bacteria will glow brilliant green upon exposure to UV light. When **arabinose** is absent from the nutrient medium, the GFP gene remains turned off and the colonies appear white.

This is an excellent example of the central dogma of biology in action; that is, **DNA > RNA > Protein > Trait**. For more detailed information on the **arabinose** operon and how it is used in the pGLO plasmid to regulate GFP expression, see Appendix B.

## THINQ! Exercises

What are other ways to use genetic transformation? Search the Internet and explain other real-world applications of genetic transformation that are used in biotechnology.

Define the following:

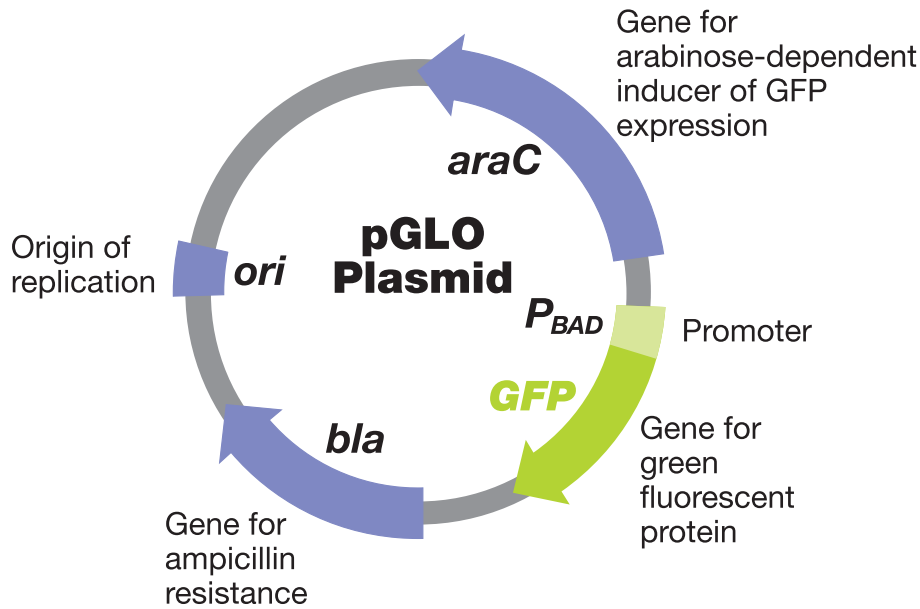
Plasmid:

LB nutrient medium:

Ampicillin:

Arabinose:

## pGLO Inquiry



**The pGLO plasmid.** *GFP*, green fluorescent protein; *araC*, gene for arabinose-dependent inducer of GFP expression; *bla*, gene for ampicillin resistance; *ori*, origin of replication.

You will be provided with tools and a protocol for performing genetic transformation. Your task will be to:

1. **Perform the genetic transformation.** The procedure involves three main steps: two to move the pGLO plasmid DNA through the *E. coli* cell membrane and one to provide an environment for the cells to express the newly acquired genes.
  - a) Incubate actively growing bacteria with plasmid DNA in CaCl<sub>2</sub> solution.
  - b) Rapidly heat and cool the bacteria in a process known as heat shock during which the plasmid DNA enters the bacteria.
  - c) Transfer the bacteria to growth media to recover and express their newly acquired genes.
2. **Assess the degree of success.** You will determine transformation efficiency.

Along the way, you will learn about the reasons for each step and how and to what extent each step influences gene transfer.

## THINQ! Exercises

Steps for designing a scientific investigation:

- Observe the natural world
- Ask questions about your observations
- Formulate a reasonable hypothesis to explain the observations
- Create and execute experiments testing the hypothesis and generating data
- Analyze the data, compare to the hypothesis, and communicate your findings

Define the following:

Phenotype:

Genotype:

### **Focus Questions**

Scientific investigations begin with an observation about the natural world and the formulation of questions about that observation. Below are a few questions for you to ponder as you take on the challenge of performing a genetic transformation lab.

#### **Question 1: Which organism should I choose, and why?**

1. To genetically transform an entire organism, the new gene must be in every cell in the organism. Considering this, which organism would be the simplest to work with for total genetic transformation: one composed of many cells, or one composed of a single cell?
2. Scientists often want to know if the genetically transformed organism can pass its new traits on to its offspring and future generations. To get this information, which would be a better candidate for your investigation: an organism that develops each new generation quickly, or one that reproduces more slowly?
3. Safety is another important consideration in choosing an experimental organism. What traits or characteristics should the organism have (or not have) to be sure it will not harm you or the environment?
4. Based on the above considerations, which would be the best choice for a genetic transformation: bacterium, earthworm, fish, or mouse? Describe your reasoning.

#### **Question 2: How can I tell if cells have been genetically transformed?**

The goal of genetic transformation is to change an organism's traits, also known as its **phenotype**. Before a change in the phenotype can be detected, however, a thorough examination of its natural (pretransformation) phenotype must be made.

1. Describe how you could use two LB/agar plates, some *E. coli*, and some ampicillin to determine how *E. coli* cells are affected by ampicillin.
2. What would you expect your experimental results to indicate about the effect of ampicillin on the *E. coli* cells?

## II. Investigation #1: pGLO Bacterial Transformation Laboratory (Structured Inquiry)

In this activity, you will perform a bacterial transformation, transforming a stock *E. coli* culture with the pGLO plasmid.

<b>Student workstations</b>	<b>Quantity</b>
<input type="checkbox"/> <i>E. coli</i> starter plate	1
<input type="checkbox"/> Poured nutrient agar plates (1 LB, 2 LB/amp, 1 LB/amp/ara)	4
<input type="checkbox"/> Transformation solution (1 ml)	1
<input type="checkbox"/> LB nutrient broth (1 ml)	1
<input type="checkbox"/> Inoculation loops (1 pk of 10)	7
<input type="checkbox"/> Disposable plastic transfer pipets (DPTPs)	4
<input type="checkbox"/> Foam microcentrifuge tube holder/float	1
<input type="checkbox"/> Container full of crushed ice (foam/paper cups)	1
<input type="checkbox"/> Microcentrifuge tubes	2
<input type="checkbox"/> Marking pen	1

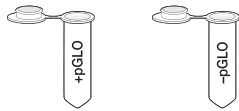
<b>Common workstation</b>	<b>Quantity</b>
<input type="checkbox"/> Rehydrated pGLO plasmid, vial	1
<input type="checkbox"/> 42°C water bath and thermometer	1
<input type="checkbox"/> UV pen light	1
<input type="checkbox"/> 37°C incubator	1
<input type="checkbox"/> Clock or timer for counting seconds	1

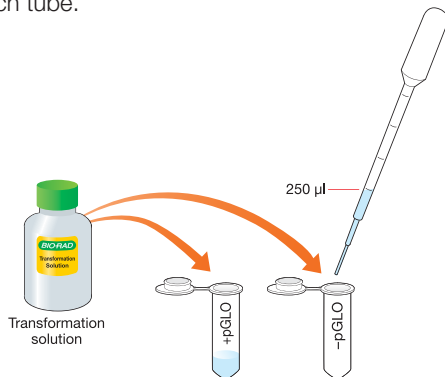
<b>Optional</b>	<b>Quantity</b>
<input type="checkbox"/> 2–20 µl adjustable volume micropipet	1
<input type="checkbox"/> 2–20 µl micropipet tips, box	1

### Protocol

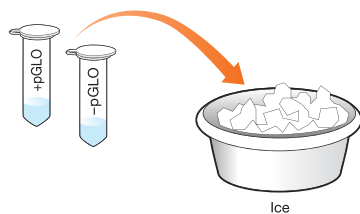
1. Label one microcentrifuge tube **+pGLO** and another **-pGLO**. Label both tubes with your group's name. Place them in the foam tube rack.



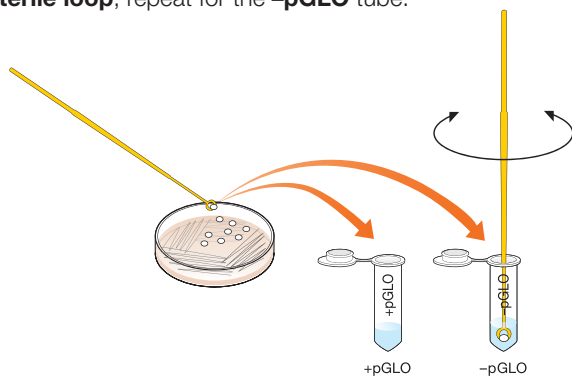
2. Open the tubes and use a sterile DPTP to transfer 250  $\mu$ l of transformation solution (50 mM  $\text{CaCl}_2$ ) into each tube.



3. Place the tubes on ice.



4. Use a sterile loop to pick **2-4 large colonies of bacteria** from the starter plate. Select colonies that are "fat" (1-2 mm in diameter). It is important to take individual colonies (not a swab of bacteria from the dense portion of the plate), since the bacteria must be actively growing to achieve high transformation efficiency. Pick up the **+pGLO** tube and immerse the loop into the transformation solution in the tube. Spin the loop between your index finger and thumb until the colonies are dispersed in the transformation solution (there are no floating chunks). Place the tube back in the tube rack in the ice. Using a **new sterile loop**, repeat for the **-pGLO** tube.



## THINQ! Exercises

**Collaborate and use outside resources to answer the following questions:**

Examine the bottle of pGLO plasmid DNA solution with the UV lamp. What do you see? Note your observations:

Look at the individual colonies of *E. coli* on your starter plates. On a separate piece of paper list all observable traits or characteristics that can be described. For example:

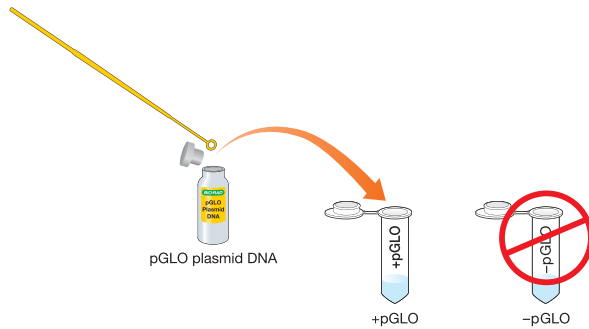
- Color of colonies
- Size of: 1) the largest colony; 2) the smallest colony; 3) the majority of colonies
- Shape of colonies (both 2-D and 3-D)
- Appearance of the colonies under regular and UV light

Why do you add the  $\text{CaCl}_2$  transformation solution?

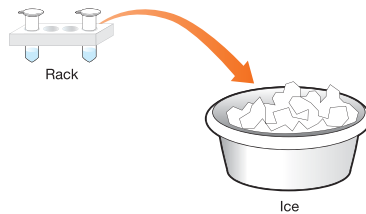
Why do you place the tubes on ice?



5. Immerse a **new sterile loop** into the tube of  $0.8 \times 10^{-1} \mu\text{g}/\mu\text{l}$  solution of pGLO plasmid DNA stock tube. Withdraw a loopful (10  $\mu\text{l}$ ). You should see a film of plasmid solution across the ring, similar to the soapy film across a ring for blowing soap bubbles. Mix the loop into the cell suspension of the **+pGLO** tube. **Do not add plasmid DNA to the -pGLO tube.** Close both the **+pGLO** and **-pGLO** tubes and return them to the rack on ice.



6. Incubate the tubes on ice for 10 min. Push the tubes all the way down in the rack so the bottoms of the tubes make contact with the ice.



7. While the tubes are on ice, label the four LB nutrient agar plates on the bottom (not the lid):
- Label one **LB/amp** plate: **+pGLO**
  - Label the **LB/amp/ara** plate: **+pGLO**
  - Label the **LB** plate: **-pGLO**
  - Label the other **LB/amp** plate: **-pGLO**



## THINQ! Exercises

**Collaborate and use outside resources to answer the following:**

Approximately how much volume is picked up by the loop if the solution is  $0.8 \times 10^{-1} \mu\text{g}/\mu\text{l}$  and a loopful of solution contains  $0.8 \mu\text{g}$  of pGLO plasmid?

Why do you incubate the tubes on ice for 10 minutes?

Why do you use only four LB nutrient agar plates, as opposed to 6 plates (2 LB, 2 LB/amp, and 2 LB/amp/ara)?

What would you expect to grow on the following LB nutrient agar plates?

**+pGLO on LB:**

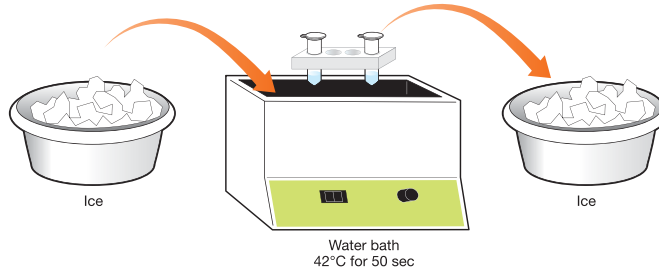
**-pGLO on LB/amp/ara:**

Why do you heat shock the cells?

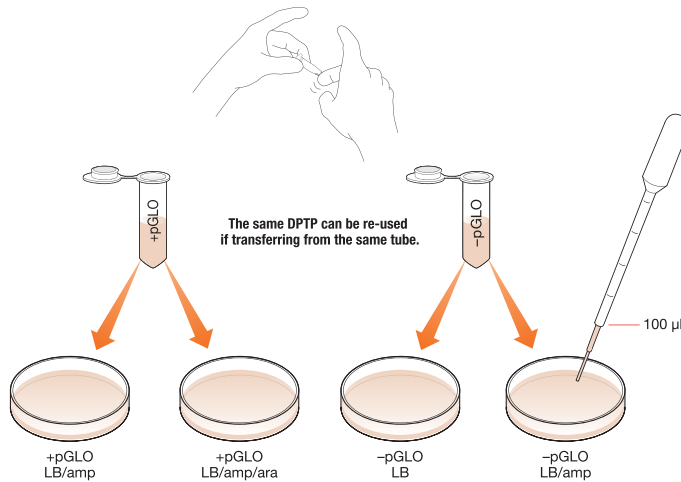
Why do you incubate on ice for 2 minutes?

8. **Heat shock.** Using the foam holder as a rack, transfer both the **+pGLO** and **-pGLO** tubes into the **42°C** water bath for **exactly 50 sec**. Push the tubes all the way down in the rack so the bottoms of the tubes make contact with the warm water.

When the 50 sec has passed, place both tubes back on ice. For best results, the transfer from the ice (**0°C**) to the **42°C** water and back to the ice must be rapid. Incubate tubes on ice for **2 min**.



9. Remove the rack of tubes from the ice and place it on the benchtop. Open a tube and use a new sterile DPTP to add 250  $\mu$ l of LB nutrient broth. Close the tube. Use the same DPTP for the other tube. Incubate the tubes for 10 min at room temperature.
10. Gently flick the closed tubes with your finger to mix and resuspend the bacteria. Using a new sterile DPTP for each tube, pipet 100  $\mu$ l of the transformation and control suspensions onto the appropriate agar plates.

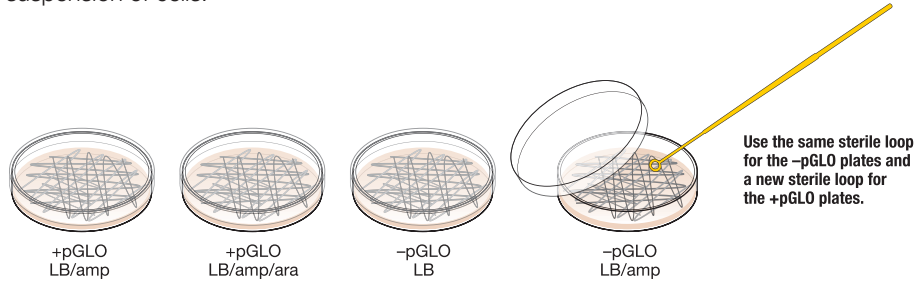


## THINQ! Exercises

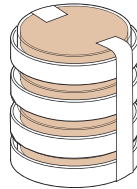
**Collaborate and use outside resources to answer the following.**

Which is your control plate? Why?

11. **Spread the suspensions evenly using one new sterile loop for the +pGLO plates and one new sterile loop for the -pGLO plates.** Spread the +pGLO LB/amp plate first, then the +pGLO LB/amp/ara plate. Then, using a new sterile loop for the -pGLO plates, spread the -pGLO LB plate first, then the -pGLO LB/amp plate. On the surface of each LB nutrient agar plate quickly skate the flat surface of the sterile loop back and forth across the plate surface. Do not press into the agar. Minimize contamination by uncovering one plate at a time and re-covering it immediately after spreading the suspension of cells.



12. Stack the plates and tape them together. Write your group name and class period on the bottom of the bottom plate in the stack and place the stack of plates **upside down** in the 37°C incubator until the next day.



**IMPORTANT:** After analyzing the results, **save transformed pGLO plates** for the inquiry labs that follow.

## THINQ! Exercises

**Collaborate and use outside resources to answer the following questions:**

Why do you need to place the stack of plates upside down?

Alternatively, you could incubate your plates at room temperature. What difference would you expect if plates were incubated at 37°C vs. room temperature (22°C)?

### Review Questions:

On which of the plates would you expect to find bacteria most like the original untransformed *E. coli* colonies you initially observed? Explain your prediction.

If there are any transformed bacterial cells, on which plate(s) would they most likely be located? Explain your prediction.

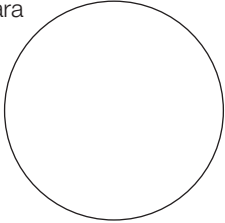
Which plates should be compared to determine if any genetic transformation has occurred? Why?

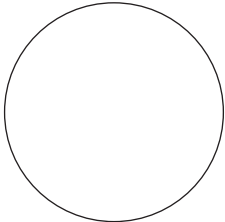
### Data Collection

Observe the results you obtained from the transformation lab under normal room lighting. Then turn out the lights and hold the ultraviolet (UV) light over the plates.

Carefully observe and draw what you see on each of the four plates. Put your drawings in the data table below. Record your data to allow you to compare your observations of the **+pGLO** cells with your observations of the non-transformed *E. coli*. Write down the following observations for each plate.

1. How much bacterial growth do you see on each plate, relatively speaking?
2. What color are the bacteria under normal light and UV light conditions?
3. How many bacterial colonies are on each plate (count the colonies you see).

Observations	
Transformation plates	+pGLO LB/amp 
	+pGLO LB/amp/ara 

Observations	
Control plates	-pGLO LB/amp 
	-pGLO LB 

### **Analysis of Results**

The goal of this analysis is to determine whether genetic transformation has occurred.

1. How did the traits you originally observed for *E. coli* alter?

2. If the transformed cells have acquired the ability to grow in the presence of ampicillin, then what might be inferred about their ability to glow bright green under UV light?

3. From the results that you obtained, how could you provide evidence to support your hypothesis/argument that the changes that occurred were due to the procedure that you performed?

### **The Interaction between Genes and Environment**

Look again at the four plates. Do you observe some *E. coli* growing on the LB plate that does not contain ampicillin or arabinose?

1. From your results, can you tell if these bacteria are ampicillin resistant by looking at them on the LB plate? Explain your answer.
  
2. What might happen to these bacteria if you moved them to plates containing ampicillin?
  
3. Often an organism's traits are caused by a combination of its genes and its environment. Think about the green color you saw in the genetically transformed bacteria as you consider these questions:
  - a) What two factors must be present in the bacteria's environment for you to see the green color?
  
  
  
  
  
  
  
  
  
  
  - b) Provide another example of a change in the environment causing expression of a different trait.
  
  
  
  
  
  
  
  
  
  
  - c) What advantage would there be for an organism to be able to turn on or off particular genes in response to certain conditions?

### Calculation of Transformation Efficiency

The next task in this investigation will be to determine how many of the *E. coli* cells were transformed. This quantitative measurement is referred to as transformation efficiency.

In many experiments, it is important to transform as many cells as possible. For example, in some types of gene therapy, cells are collected from a patient, transformed in the laboratory, and then put back into the patient. The more cells that are transformed to produce the needed protein, the more likely the therapy will work. Transformation efficiency helps scientists determine how well transformation is working.

### The Task

You are about to calculate the transformation efficiency for this experiment, which indicates how effective you were in getting new DNA molecules into bacterial cells.

Transformation efficiency is a number: the total number of colonies growing on the plate divided by the amount of DNA spread on the plate. It represents the total number of bacterial cells transformed using one microgram of DNA. Each colony on the plate can be assumed to derive from a single cell. As individual cells reproduce, more and more cells accumulate, developing into a colony. The most direct way to determine the total number of bacteria that were transformed with the pGLO plasmid is to count the number of colonies on the plate. Transformation efficiency is calculated using the following formula:

$$\text{Transformation efficiency} = \frac{\text{Total number of colonies growing on the agar plate}}{\text{Amount of DNA spread on the agar plate (in } \mu\text{g)}}$$

Therefore, before you can calculate the efficiency of your transformation, you will need two pieces of information:

1. The total number of green fluorescent colonies growing on your LB/amp/ara plate.
2. The total amount of pGLO plasmid DNA used for bacterial transformation that was spread on the LB/amp/ara plate.

#### 1. Determining the Total Number of Transformed Green Fluorescent Colonies

Place the LB/amp/ara plate near a UV light source. Count the number of green fluorescent colonies that glow under UV light on the plate.

Enter that number here →

#### 2. Determining the Amount of pGLO DNA in the Cells Spread on the LB/Amp/Ara Plate

You need two pieces of information to determine the amount of pGLO DNA in the bacterial cells spread on the LB/amp/ara plate in this experiment: (a) the total amount of DNA we began the experiment with, and (b) the fraction of the DNA (in the bacteria) that was spread onto the LB/amp/ara plates.

Once you calculate these data, you multiply the total amount of pGLO DNA used in this experiment by the fraction of DNA you spread on the LB/amp/ara plate. This will tell you the amount of pGLO DNA in the bacterial cells that were spread on the LB/amp/ara plate.

##### A. Determining the total amount of pGLO plasmid DNA

The total amount of DNA we began with is equal to the product of the concentration and the total volume used, or

$$(\text{DNA in } \mu\text{g}) = (\text{concentration of DNA in } \mu\text{g}/\mu\text{l}) \times (\text{volume of DNA in } \mu\text{l})$$

In this experiment you used 10  $\mu\text{l}$  of pGLO at a concentration of 0.08  $\mu\text{g}/\mu\text{l}$ . This means that each microliter of solution contained 0.08  $\mu\text{g}$  of pGLO DNA. Calculate the **total amount of DNA** used in this experiment

Enter that number here →

**How will you use this piece of information?**

**B. Determining the fraction of pGLO plasmid DNA (in the bacteria)**

Not all the DNA you added to the bacterial cells was transferred to the agar plate; therefore, you must determine the fraction of the DNA actually spread onto the LB/amp/ara plate. To do this, divide the volume of DNA you spread on the LB/amp/ara plate by the total volume of liquid in the test tube containing the DNA:

$$\text{Fraction of DNA used} = \frac{\text{Sample volume spread on LB/amp/ara plate, } \mu\text{l}}{\text{Total sample volume in microcentrifuge tube, } \mu\text{l}}$$

You spread 100  $\mu\text{l}$  of cells containing DNA from a test tube containing a total volume of 510  $\mu\text{l}$  of solution. Do you remember why there is 510  $\mu\text{l}$  total solution? Look in the laboratory procedure and locate all the steps where you added liquid to the reaction tube. Add the volumes.

Use the above formula to calculate the **fraction of pGLO plasmid DNA** you spread on the LB/amp/ara plate.

Enter that number here →

**How will you use this piece of information?**

**C. So how many micrograms of pGLO DNA did you spread on the LB/amp/ara plate?**

To answer this question, you will need to multiply the total amount of pGLO DNA used in this experiment by the fraction of pGLO DNA you spread on the LB/amp/ara plate.

$$\text{pGLO DNA spread } (\mu\text{g}) = \text{amount of DNA used } (\mu\text{g}) \times \text{fraction of DNA}$$

Enter that number here →

**What will this number tell you?**

Look at all your calculations above. Decide which of the numbers you calculated belong in the table below. Fill in the following table.

Number of colonies on LB/amp/ara plate = _____
pGLO DNA spread on the plate, $\mu\text{g}$ = _____



Now use the data in the table to calculate the efficiency of the pGLO transformation:

$$\text{Transformation efficiency} = \frac{\text{Total number of colonies growing on the agar plate}}{\text{Amount of DNA spread on the agar plate (in } \mu\text{g)}}$$

Enter that number here → Transformation efficiency = \_\_\_\_\_ transformants/ $\mu$ g

### **The Analysis**

Transformation efficiency calculations result in very large numbers. Scientists often use a mathematical shorthand, referred to as scientific notation, to express large numbers. For example, if the calculated transformation efficiency is 1,000 bacteria/ $\mu$ g of DNA, they often report this number as:

$10^3$  transformants/ $\mu$ g      ( $10^3$  is another way of saying  $10 \times 10 \times 10$  or 1,000)

**1. How would scientists report 10,000 transformants/ $\mu$ g in scientific notation?**

Carrying this idea a little further, suppose scientists calculated an efficiency of 5,000 bacteria/ $\mu$ g of DNA. This would be reported as:

$5 \times 10^3$  transformants/ $\mu$ g      ( $5 \times 1,000$ )

**2. How would scientists report 40,000 transformants/ $\mu$ g in scientific notation?**

One final example: If 2,600 transformants/ $\mu$ g were calculated, then the scientific notation for this number would be:

$2.6 \times 10^3$  transformants/ $\mu$ g      ( $2.6 \times 1,000$ )

Similarly:

$5,600 = 5.6 \times 10^3$      $271,000 = 2.71 \times 10^5$      $2,420,000 = 2.42 \times 10^6$

**3. How would scientists report 960,000 transformants/ $\mu$ g in scientific notation?**

4. Report your calculated transformation efficiency in scientific notation.

5. Use a sentence or two to explain what your calculation of transformation efficiency means.

Biotechnologists generally agree that the transformation protocol you just completed usually has a transformation efficiency of between  $8.0 \times 10^2$  and  $7.0 \times 10^3$  transformants per microgram of DNA.

6. How does your transformation efficiency compare with the above?

7. In the table below, report the transformation efficiency of several of the teams in the class.

<i>Team</i>	<i>Efficiency</i>

How does your transformation efficiency compare to theirs?

**IMPORTANT:** Save the **transformed pGLO plates** for the inquiry labs that follow.

### III. Investigations #2–5: Guided and Open Inquiry

Below is a general list of materials that could be used for the following inquiry investigations. It will be up to you to determine what materials and supplies you'll need to design and develop your inquiry investigations. Check with your teacher to make sure it is okay to use all of the materials you think you need.

#### **Guided and Open Inquiry Materials Checklist**

- E. coli* starter plate
- pGLO transformed plate(s) from the pGLO Bacterial Transformation lab
- pGLO transformed plate(s) with satellite colonies from the pGLO Bacterial Transformation lab
- Poured agar plates
  - LB nutrient agar plates
  - LB/amp agar plates
  - LB/amp/ara agar plates
- Transformation solution
- LB nutrient broth
- Inoculation loops
- Disposable plastic transfer pipets
- Foam microcentrifuge tube holder/float
- Container (such as foam cup) full of crushed ice (not cubed ice)
- Marking pen
- Microcentrifuge tubes
- pGLO plasmid ( $0.8 \times 10^{-1}$  µg/µl stock solution)
- Ampicillin solution (100 mg/ml stock solution)
- Arabinose solution (300 mg/ml stock solution)
- Sterile water

#### **Guided and Open Inquiry Investigation Accessories**

- 42°C water bath and thermometer
- UV pen light
- 37°C incubator
- Clock or timer for counting seconds
- Forceps
- 2–20 µl adjustable volume micropipet (optional)
- 2–20 µl micropipet tips (optional)

## Investigation #2: Transformation Efficiency (Guided Inquiry)

### Overview

Earlier, you determined how well your bacterial transformation worked. You did this by calculating the transformation efficiency: You calculated how many colonies of transformed bacteria you could get for every microgram of DNA you used.

$$\text{Transformation efficiency} = \frac{\text{Total number of colonies growing on the agar plate}}{\text{Amount of DNA spread on the agar plate (in } \mu\text{g)}}$$

Knowing transformation efficiency has many useful applications. For example, bacterial transformation is used as a method to produce more plasmid (cloning) or to manufacture gene products (enzymes and other proteins, such as insulin). In these situations, optimizing your transformation efficiency allows you to produce the most plasmid or protein possible. Similarly, if you are working with a new strain of bacteria, you may need to adjust parameters within the transformation protocol to help maximize efficiency.

In this guided set of inquiry experiments using the pGLO Bacterial Transformation protocol and supplies, you will examine which variables in the protocol can be adjusted (and by how much) to maximize transformation efficiency.

### Designing the Procedure

In this section, you will develop a procedure to increase transformation efficiency by changing the Bio-Rad protocols you used previously. Due to the amount of material provided in the kit and the number of variables you could change to increase transformation efficiency, you will need to work within the following parameters:

- You will have 2 agar plates: 2 LB/amp –OR– 2 LB/amp/ara
- One (1) of your agar plates must be used for the control condition
- All the agar plates will have been prepared in a standard manner
- You will be provided with an *E. coli* starter plate
- You will use the materials provided by your teacher, unless otherwise noted

Based on the inquiry investigation materials checklist on page 16 and the parameters given above, list the materials and equipment that you will need to conduct your experiment.

**Materials needed:**

**Equipment needed:**

## ThINQ! Exercises

**Collaborate and use outside resources to answer the following questions:**

What are some specific examples where bacterial transformation is commonly used to manufacture gene products?

What is a control condition and why it is important?

### **Steps for Designing a Scientific Investigation**

1. Make observations about the natural world.

Before you can begin to design your experiment or develop your hypothesis, you need to determine the variables that affect bacterial transformation. In the pGLO bacterial transformation lab, which step(s) of the protocol do you think had the greatest effect on the efficiency of bacterial transformation?

<b>Critical Protocol Steps</b>	<b>Variable(s)</b>

2. Ask questions about those observations.

Develop a question you would like to explore, given the observations and variable(s) you noted above, that could affect transformation efficiency. It will be impossible to test all these conditions at once, so focus on a single protocol step and explore one of its variables. Write your question below:

3. Formulate a reasonable hypothesis to explain your observations.

Formulate a reasonable hypothesis that you can test, given the supplies for inquiry investigations. The simplest form of a hypothesis is an if/then format. Think about what you are trying to determine with this experiment and how you will know whether your hypothesis is supported. Write your hypothesis below:

4. Create and execute experiments to test the hypothesis and generate data.

Design your experiment. What variables will you manipulate and what variables should you keep constant to test your hypothesis? (Hint: Don't forget to describe what controls you have designed for the experiment.)

Consider your hypothesis and your experiment and predict what you think you will observe.

5. Analyze your data, compare to the hypothesis, and communicate your findings.

To be able to illustrate results graphically and make comparisons, you will need to interpret your results quantitatively (numerically) as well as qualitatively (with words). Please explain how you will quantify your results (that is, measure them).

### Writing the Procedure

To help keep track of your experiment, fill in the table below as a quick guide. We have provided some of the critical steps for transformation. For the best comparison, vary only the experimental variable(s) chosen while keeping all other steps the same as the control.

<b>Critical Protocol Step</b>	<b>Control Condition</b>	<b>Experimental Conditions</b>	
Transformation solution (CaCl <sub>2</sub> )	50 mM CaCl <sub>2</sub>	(CaCl <sub>2</sub> ) (mM) =	
Amount pGLO DNA	0.8 µg (1 loopful)	Amount (µg) =	
Pre-incubation of DNA	10 min on ice	Time(s) =	Temperature(s) =
Heat shock	50 sec @ 42°C	Time(s) =	Temperature(s) =
Heat shock recovery	2 min on ice	Time(s) =	Temperature(s) =
Growth recovery	10 min @ room temperature	Time(s) =	Temperature(s) =

When you develop a protocol, it is helpful to draw pictures for each step you carry out. Think about the materials you have and the question you are trying to answer. You can use outside resources to help generate ideas and guide you before you begin to write your protocols.

Write or illustrate your proposed protocols in a lab notebook or on a sheet of lab notebook paper. Ask your teacher to review the experiment before you begin.

**Results**

Use this page to restate your hypothesis and summarize your results.

**Hypothesis****Results**

What did you do to test the hypothesis?

What were the results? Calculate and compare the transformation efficiency for each of the conditions you tested.

Do the data support your hypothesis?

List any observations relevant to your experiment.

List any ideas you have for refining your hypothesis and testing your experiment.



### Investigation #3: Effect of Ampicillin on Bacterial Growth (Open Inquiry)

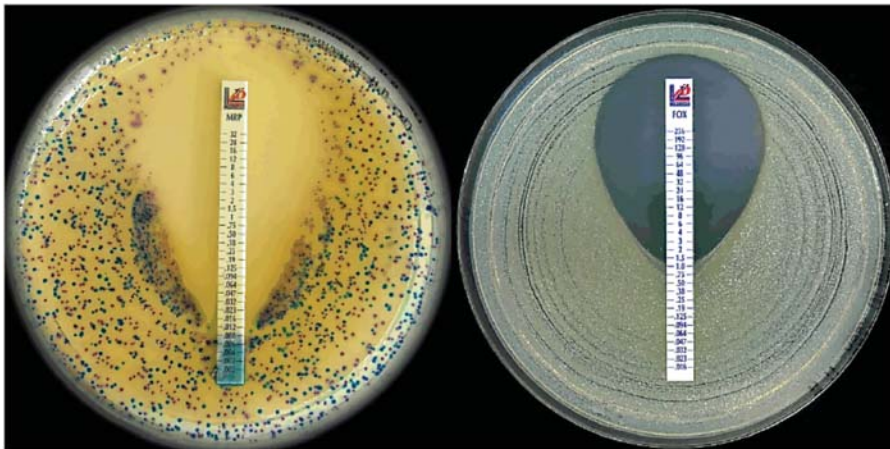
#### Overview

Imagine you are planning to work with a new strain of bacteria. How would you know which antibiotic to use, and how much of it, for selection in your transformation experiments? In this set of experiments, you will use the supplies in the pGLO Transformation and Inquiry Kit and some extra information to design an experiment that tests the effects of ampicillin on bacterial growth.

Ampicillin is an antibiotic used to treat a number of bacterial infections, such as bronchitis, sinus and ear infections. Ampicillin reduces the growth of bacteria by acting as an irreversible inhibitor of the enzyme transpeptidase, which is needed by bacteria to make cell walls. Though ampicillin is often a first line of treatment for these common infections, some strains of bacteria have developed resistance to it. Antibiotics like ampicillin are also used routinely in bacterial transformation experiments to help select for transformants.

Kirby-Bauer (KB) and Epsilonometer (E) testing (diffusion antibiotic sensitivity testing) provide simple but effective ways of quickly assessing whether a particular bacterium is affected by an antibiotic. In the KB test, a filter disk is impregnated with an antibiotic and placed on an agar plate with bacteria and the plate is left to incubate. If the antibiotic stops growth of (or kills) the bacteria, there will be a clear area around the filter disk where the bacteria have not grown (the zone of inhibition). The zone of inhibition includes the disk and is measured to the nearest millimeter using a ruler across the diameter of the circular zone of inhibition. Bacterial growth touching or around the edge of the disk is reported as a zone of 0 mm. The size of this zone depends on how effective the antibiotic is at stopping the growth of the bacterium.

The E test is a bit more quantitative. A rectangular strip with an exponential scale printed on it is impregnated with the drug to be studied. The strip is laid on a prepared plate and, after a 24-hour incubation period, an elliptical zone of inhibition is produced. The point at which the ellipse meets the strip indicates the minimum inhibitory concentration (MIC) of the drug. The information from these tests can be used to choose appropriate antibiotics and concentrations of those antibiotics for selection in transformation experiments.



Mixed culture on Chromatic MH  
MIC Test Strip Meropenem 0.002-32 µg/ml.

Staphylococcus ATCC® 29213  
MIC Test Strip Cefoxitin 0.016-256 µg/ml.  
M.I.C. = 1.5 µg/ml.

Examples of Epsilonometer (E) tests from a Liofilchem MIC Test Strip.

## ThINQ! Exercises

Collaborate and use outside resources to answer the following questions:

Define the following terms:

Transpeptidase:

Transformant:

In your own words, describe how the Kirby-Bauer test relates to antibiotic resistance:

### ***Designing the Procedure***

In this section, you will formulate a hypothesis and develop a procedure to test the effect of ampicillin on bacterial growth. Due to the amount of material provided in the kit, you will need to work within the following parameters:

- You will have 1 LB agar plate
- You will be provided with an *E. coli* starter plate
- You will need to use the materials provided by your teacher, unless otherwise noted

Based on the inquiry investigation material checklist on page 16 and the parameters given above, list the materials and equipment that you will need to conduct your experiment.

**Materials needed:**

**Equipment needed:**

### ***Steps for Designing a Scientific Investigation***

1. Make observations about the natural world.

Think about what you have read about antibiotic resistance, the Kirby-Bauer test, and the E test. Make some observations about the natural world and the pGLO experiment that may relate to antibiotic dosage and resistance.

2. Ask questions about those observations.

Develop a question you would like to explore, given the observations you have made and components of the pGLO bacterial transformation kit. Write your question below:

3. Formulate a reasonable hypothesis to explain your observations.

Formulate a reasonable hypothesis that you could test, given the supplies for inquiry investigations. The simplest form of a hypothesis is an if/then format. Think about what you are trying to determine with this experiment and how you will know whether your hypothesis is supported. Write your hypothesis below:

4. Create and execute experiments to test the hypothesis and generate data.

Design your experiment. What variables will you manipulate and what variables should you keep constant to test your hypothesis?

5. Analyze your data, compare to the hypothesis, and communicate your findings.

Consider your hypothesis and your experiment and predict what you think you will observe.

To illustrate results graphically and make comparisons, you will need to interpret your results quantitatively (numerically) as well as qualitatively (with words). Please explain how you will quantify your results (that is, measure them).

**Writing the Procedure**

When you develop a protocol, it is helpful to draw pictures for each step you carry out. Think about the materials you have and the question you are trying to answer. You can use outside resources to help generate ideas and guide you before you begin to write your protocols.

Write or illustrate your proposed protocols in a lab notebook or on a sheet of lab notebook paper. Ask your teacher to review the experiment before you begin.

**Results**

Use this page to restate your hypothesis and summarize your results.

**Hypothesis****Results**

What did you do to test the hypothesis?

What were the results?

Do the data support your hypothesis?

List any observations relevant to your experiment.

List any ideas you have for further refining your hypothesis and testing your experiment.

## Investigation #4: Effect of Arabinose on GFP Expression (Open Inquiry)

### Overview

#### Why Is Gene Expression Regulated?

Gene expression is carefully regulated in part to allow organisms to adapt to different environmental conditions. Genes can be turned on when particular proteins are needed and off when they are no longer necessary, preventing wasted effort and overproduction of proteins.

For bacteria, the genes involved in the breakdown of different food sources are often highly regulated. The simple plant sugar arabinose, for example, is a source of both energy and carbon for bacteria. The bacterial genes that encode digestive enzymes that break down arabinose for food are not expressed when arabinose is not in the environment. When arabinose is present, however, these genes are turned on, and when the arabinose runs out, the genes are turned off again. Arabinose initiates transcription of these genes by promoting the binding of RNA polymerase.

In the genetically engineered pGLO plasmid, some of the genes involved in arabinose breakdown have been replaced by the jellyfish gene that encodes GFP. Therefore, the expression of GFP is induced by arabinose. When bacteria transformed with the pGLO plasmid are grown in the presence of arabinose, expression of the GFP gene is turned on, and the bacteria glow brilliant green when exposed to UV light.

#### Why Regulate Expression of a Gene We Transform into Bacteria?

In the wild, regulation of gene expression is clearly beneficial. But why would it be necessary in the lab, when we are trying to get our bacteria to glow green?

One reason is that certain proteins, when overexpressed, may be toxic to the bacteria. If production of these proteins is not regulated, the toxic proteins would quickly accumulate and kill the bacteria that are producing them.

Another reason is that bacteria expend significant amounts of energy and environmental resources to grow and divide. Consumption of these energy and resource pools by constant production of a large amount of a foreign protein may interfere with the bacteria's ability to grow and divide. In this situation, bacteria would overexpress the foreign protein at the expense of their ability to make other proteins needed for growth and population expansion, resulting in a smaller bacterial population.

## ThINQ! Exercises

**Collaborate and use outside resources to answer the following questions:**

Define the following terms:

Transcription:

RNA polymerase:

What are clinical trials?

What is the FDA?

What is the FDA's role in clinical trials?

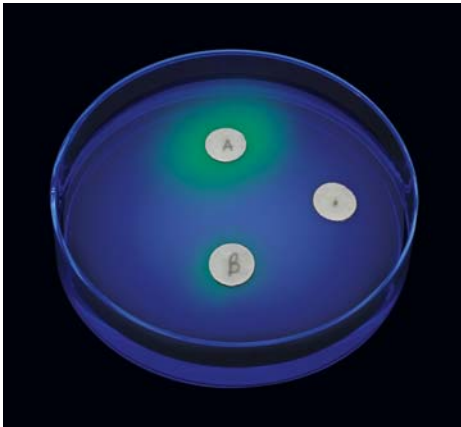
Why are FDA-approved clinical trials important?

### **Why Does It Matter if Bacterial Growth or Survival Is Negatively Affected by Expression of a Foreign Protein?**

Imagine a scenario in which you have transformed bacteria with plasmid DNA that encodes a new and potentially powerful drug (for example, insulin). You will need to make large amounts of this peptide for FDA-regulated clinical trials as a potential new treatment.

If it is difficult to isolate large amounts of this peptide and costly to synthesize large amounts of it chemically, you may wish to express large quantities of it in bacteria. However, if constant expression of this peptide either is toxic to the bacteria or prevents the bacteria from growing and dividing, you would not be able to grow enough bacteria to make enough of the peptide to carry out your FDA-regulated clinical trials.

But what if peptide production could be regulated? You could grow very large numbers of bacteria, and then activate gene expression for the plasmid-encoded peptide, which would allow you to obtain massive amounts of the peptide without worrying about its negative effects on bacteria growth or survival.



*Examples of GFP regulation with varied arabinose dosage.*

### ***Designing the Procedure***

In this section, you will develop a procedure to test your question and hypothesis about GFP gene regulation and arabinose dosage. Due to the amount of material provided in the kit, you will need to work within the following parameters:

- You will have 1 LB/amp agar plate
- You will be provided with a plate containing pGLO transformed *E. coli*
- You will need to use the materials provided by your teacher, unless otherwise noted

Based on the inquiry investigation material checklist on page 16 and the parameters given above, list the materials and equipment that you will need to conduct your experiment.

**Materials needed:**

**Equipment needed:**



### **Steps for Designing a Scientific Investigation**

1. Make observations about the natural world.

Think about what you have read about gene expression regulation and similar experiments that test antibiotic concentration dosage. Make observations about the natural world and the pGLO experiment that may relate to gene expression regulation.

2. Ask questions about those observations.

Develop a question you would like to explore given the observations you have made and materials available in the pGLO Transformation and Inquiry Kit. Write your question below:

3. Formulate a reasonable hypothesis to explain your observations.

Formulate a reasonable hypothesis that you could test given the supplies for inquiry investigations. The simplest form of a hypothesis is an if/then format. Think about what you are trying to determine with this experiment and how you will know whether your hypothesis is supported. Write your hypothesis below:

4. Create and execute experiments testing the hypothesis and generating data.

Design your experiment. What variables will you manipulate and what variables should you keep constant to test your hypothesis?

5. Analyze your data, compare to the hypothesis, and communicate your findings.

Consider your hypothesis and your experiment and predict what you think you will observe.

In order to illustrate results graphically and make comparisons, you will need to interpret your results quantitatively (numerically) as well as qualitatively (with words). Please explain how you will quantify your results (that is, measure them).

**Writing the Procedure**

When you develop a protocol, it is helpful to draw pictures for each step you carry out. Think about the materials you have and the question you are trying to answer. You can use outside resources to help generate ideas and guide you before you begin to write your protocols.

Write or illustrate your proposed protocols in a lab notebook or on a sheet of lab notebook paper. Ask your teacher to review the experiment before you begin.

**Results**

Use this page to restate your hypothesis and summarize your results.

**Hypothesis****Results**

What did you do to test the hypothesis?

What were the results?

Do the data support your hypothesis?

List any observations relevant to your experiment.

List any ideas you have for further refining your hypothesis and testing your experiment.

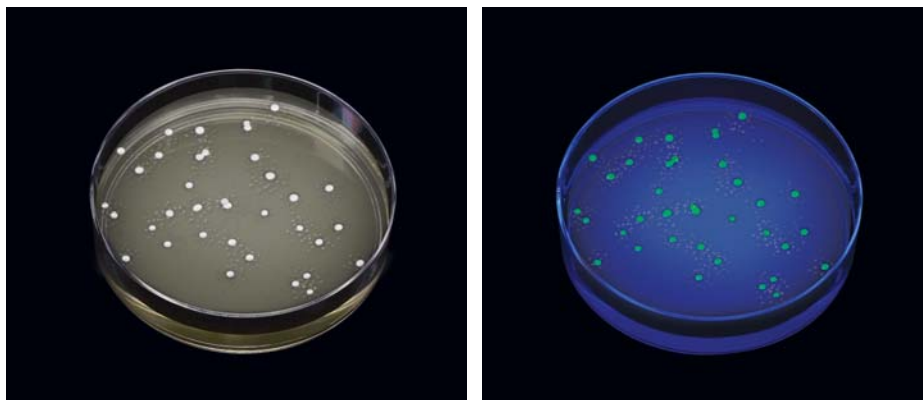
## Investigation #5: Satellite Colonies (Open Inquiry)

Recall that the pGLO plasmid DNA contains *bla*, a gene that confers resistance to the antibiotic ampicillin. Ampicillin was included in the growth medium following transformation so that only transformed bacteria would grow. This inclusion of antibiotic applies selective pressure to the bacterial population; it makes the environment on the plate inhospitable to any bacteria that cannot express genes that encode antibiotic resistance. Without this selection strategy, all bacteria would grow on the plate, and it would be difficult to determine which bacteria were transformed.

Over time, the amount of antibiotic present in the growth medium may become depleted for a number of reasons. When this happens, the selective pressure on the bacterial population is removed, and bacteria that are not antibiotic resistant may grow on the plate.

You may have noticed that your LB/amp/ara plates contain different types of colonies: the primary (larger) colonies of transformed bacteria and smaller “satellite” colonies surrounding them. This occurs especially if the plates are kept at room temperature for several days. Why might that be?

Just as a scientist in a lab producing a powerful drug like insulin would, you need to investigate what the properties of these satellite colonies are. Develop a question and design an experiment to understand what these foreign satellite colonies are and why they grew on your transformed pGLO colony plates.



**Plate containing pGLO transformants and satellite colonies.** In normal light (left), satellite colonies appear as noticeably smaller colonies surrounding larger colonies. Under UV light (right), the larger colonies glow bright green, whereas the satellite colonies do not.

## ThiNQ! Exercises

**Collaborate and use outside resources to answer the following questions:**

Can you see different types of colonies on your LB/amp/ara plates? Compare and contrast the physical characteristics of these different types of colonies.

What is selective pressure?

What are some reasons the antibiotic in the growth medium might become depleted over time?

How does the pGLO plasmid give rise to ampicillin resistance? Refer back to the plasmid map and the introductory information in the pGLO Bacterial Transformation exercise.

Why would you expect to see satellite colonies if your growth plate contains a large number of transformed colonies on a plate containing ampicillin?

### ***Designing the Procedure***

In this section, you will develop your own hypothesis about satellite colony formation. You can ask questions about whether satellite colonies contain a plasmid, or if they are resistant to an antibiotic. Try to think outside the box, but keep in mind you will need to work within the following parameters:

- You will have 2 agar plates: 1 LB/amp/ara and 1 LB
- You will be using an already prepared plate that contains satellite colonies
- You will need to use the materials provided by your teacher, unless otherwise noted

Based on the inquiry investigation material checklist on page 16 and the parameters given above, list the materials and equipment that you will need to conduct your experiment.

**Materials needed:**

**Equipment needed:**

### **Steps for Designing a Scientific Investigation**

1. Make observations about the natural world.

Think about what you have read and researched about satellite colonies and what you might need to explore about these unexpected colonies. Make observations about the natural world and the pGLO experiment that may relate to satellite colonies.

2. Ask questions about those observations.

Develop a question you would like to explore given the observations you have made and materials available in the pGLO Transformation and Inquiry Kit. Write your question below:

3. Formulate a reasonable hypothesis to explain your observations.

Formulate a reasonable hypothesis that you could test given the supplies for inquiry investigations. The simplest form of a hypothesis is an if/then format. Think about what you are trying to determine with this experiment and how you will know whether your hypothesis is supported. Write your hypothesis below:

4. Create and execute experiments testing the hypothesis and generating data.

Design your experiment. What variables will you manipulate and what variables should you keep constant to test your hypothesis?

5. Analyze your data, compare to the hypothesis, and communicate your findings.

Consider your hypothesis and your experiment and predict what you think you will observe.

In order to illustrate results graphically and make comparisons, you will need to interpret your results quantitatively (numerically) as well as qualitatively (with words). Please explain how you will quantify your results (that is, measure them).

**Writing the Procedure**

When you develop a protocol, it is helpful to draw pictures for each step you carry out. Think about the materials you have and the question you are trying to answer. You can use outside resources to help generate ideas and guide you before you begin to write your protocols.

Write or illustrate your proposed protocols in a lab notebook or on a sheet of lab notebook paper. Ask your teacher to review the experiment before you begin.

**Results**

Use this page to restate your hypothesis and summarize your results.

**Hypothesis****Results**

What did you do to test the hypothesis?

What were the results?

Do the data support your hypothesis?

List any observations relevant to your experiment.

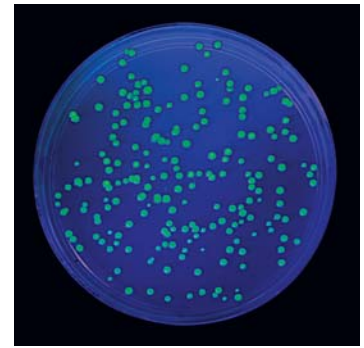
List any ideas you have for further refining your hypothesis and testing your experiment:



## Post-Lab Assessment

### Question 1

Calculate transformation efficiency from the results of the plate pictured. One loopful (10  $\mu$ l) of plasmid DNA with a concentration of 10 ng/ $\mu$ l was used. The cells were incubated in a total of 500  $\mu$ l of solution, and 100  $\mu$ l of this solution was used to inoculate the plate.



If the transformation efficiency was supposed to be  $10^8$  colonies/ $\mu$ g, explain whether the transformation was successful.

### Question 2

A biologist is interested in purifying a large amount of a protein she is studying. The protein is a fascinating new type of enzyme from a rare bacterium that grows only in deep sea hydrothermal vents. She clones the gene and expresses it in *E. coli*, hoping to use the *E. coli* as a factory for that protein. The plasmid she uses is similar to the pGLO plasmid — it includes the *bla* selectable marker (encoding ampicillin resistance) and it uses the arabinose-dependent inducer of GFP expression (*araC*) to regulate expression of the gene.

Her initial transformation efficiency is high and she continues growing transformants on LB/amp/ara plates. After several weeks, however, she notices that the transformants are less viable — they are growing more slowly, and then not at all. Through a quick test, she sees the bacteria still appear to have the plasmid.

What might be going wrong? Can you suggest a solution that might help?

**Question 3**

Olivia Hamilton, a clinical lab technician, is testing a panel of antibiotics against a new strain of bacteria that is causing infections in the local area. The goal is to help find the best treatment strategy. She performs a Kirby-Bauer test by growing lawns of the bacterial strain on agar plates with four paper disks impregnated with the same concentration of three antibiotics (A, B, and C) and a control with sterile water. A clear zone forms around the paper disks if the antibiotic inhibits bacterial growth.



Why would it be useful to have more than one antibiotic to treat these infections?

Rank the bacterial resistance to these antibiotics from highest to lowest resistance.

**Question 4**

Of the following potential applications for genetic engineering, which one is not yet done regularly and why do you think that is?

- Production of tomatoes in which ripening is delayed, making them easier to ship
- Production of hormones for treating diabetes
- Genetic testing for harmful alleles in adults (such as alleles associated with diseases like breast cancer)
- Genetic testing for harmful alleles in unborn infants (prenatal genetic testing)
- Transformation of engineered genes into human gametes (germ-line gene therapy)

## IV. Class Discussion

### Science Case Study: Can Bacterial Transformation Stop the Spread of Malaria?

#### I. The Global Impact of Malaria

Lerato sits in the clinic, watching her son Baruti, age 4, writhe with fever. He is sleeping for the first time in days, and Lerato is anxious for signs that the medication the doctor had given her son is working. She worries she may have waited too long to bring Baruti to the clinic.

Baruti had fallen ill a week before with fever, chills, and body aches, all vague flu-like symptoms that Lerato had assumed would clear in a few days, as so many illnesses had in the past. Lerato and her family live an hour's walk away from the medical clinic in Seronga, a remote village in Okavango, Botswana. Making that long trip with a sick child is difficult, but it was a trip she had to make when her son's symptoms grew more severe. He was now in a clinic bed, suffering from extreme anemia secondary to (brought on by) malaria.

Malaria is spread by mosquitoes, and like so many others in the Okavango region, Lerato's family had been issued mosquito nets for their beds. The nets are covered with insecticidal chemicals and are an effective and relatively inexpensive method for controlling mosquitoes. The nets, though, do not allow much air circulation and so are very hot to sleep under, and little Baruti tends to kick them away while he sleeps, exposing his limbs to the bites of mosquitoes.

#### Questions

1. Malaria is the third leading cause of infectious disease death in the world, after tuberculosis and AIDS. According to the World Health Organization, 3.4 billion people — nearly half the global population — are currently at risk for malaria. Most prevalent in African or tropical Asian countries, malaria is often considered a “disease of the developing world.” Though vaccines are not yet available, it can be cured if diagnosed and treated promptly.

Given this information, what might the biggest hurdles be in fighting malaria? Consider the regions the disease affects and the challenges faced by the people living there.

2. Malaria can be spread only through the bite of a mosquito, and it was nearly eliminated in the U.S. back in the 1950s. Despite this, as many as 1,500–2,000 new cases of malaria are reported in the U.S. annually. How can this be? How might malaria be coming into the country?

3. Considering that malaria can be spread only from infected blood and through mosquito bites, how might malaria eventually be eradicated in a particular region, such as the U.S.?

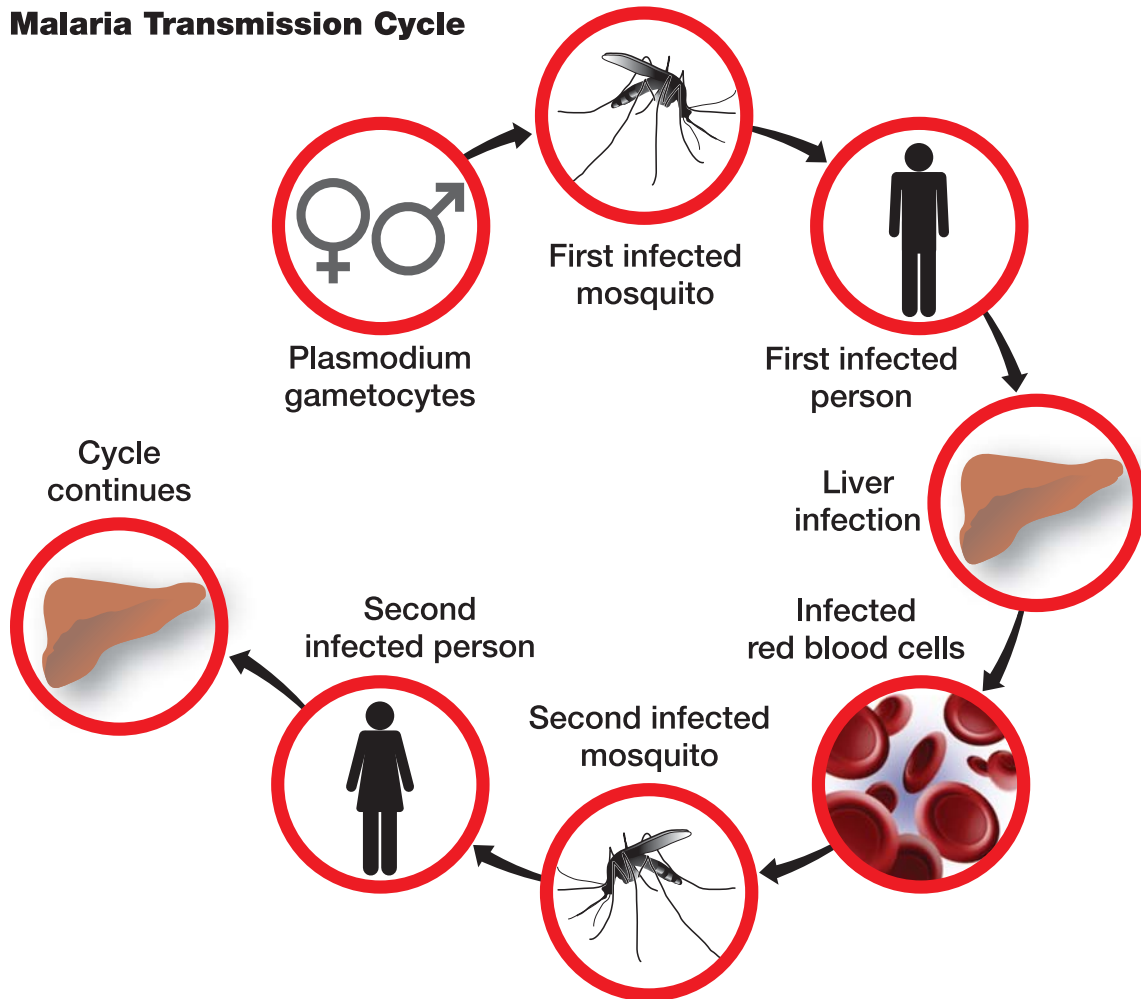
## II. Mosquitoes — Flying Factories of Malaria

In order to expand the discussion of malaria and possible methods for its treatment, control, and eradication, it is important to understand the biology behind the disease.

### *Plasmodium* — the parasitic protist behind malaria

Malaria is a parasitic infection caused by single-celled protists in the genus *Plasmodium*. Of the more than 100 species of *Plasmodium*, only four infect humans and cause disease. *Plasmodium* is a member of the phylum *Apicomplexa*, a fascinating group of protists believed to have evolved from photosynthetic dinoflagellates (forms of plankton). It has a complex life cycle that will not be discussed here except to say that to complete that life cycle, *Plasmodium* requires two hosts: (1) humans and (2) the female *Anopheles* mosquito.

### Malaria Transmission Cycle



**The *Plasmodium* life cycle depends on two hosts: humans and mosquitoes.**

In humans, the *Plasmodium* parasite invades the cells of the liver, lymph nodes, and red blood cells (erythrocytes) in the bloodstream, where it replicates and eventually causes cells to rupture. The human immune system also responds to the invasion, producing the high fevers, nausea, diarrhea, and other flu-like symptoms characteristic of malaria. In mosquitoes, *Plasmodium* lives in the gut and salivary glands; it has no known negative effects on the mosquito host.

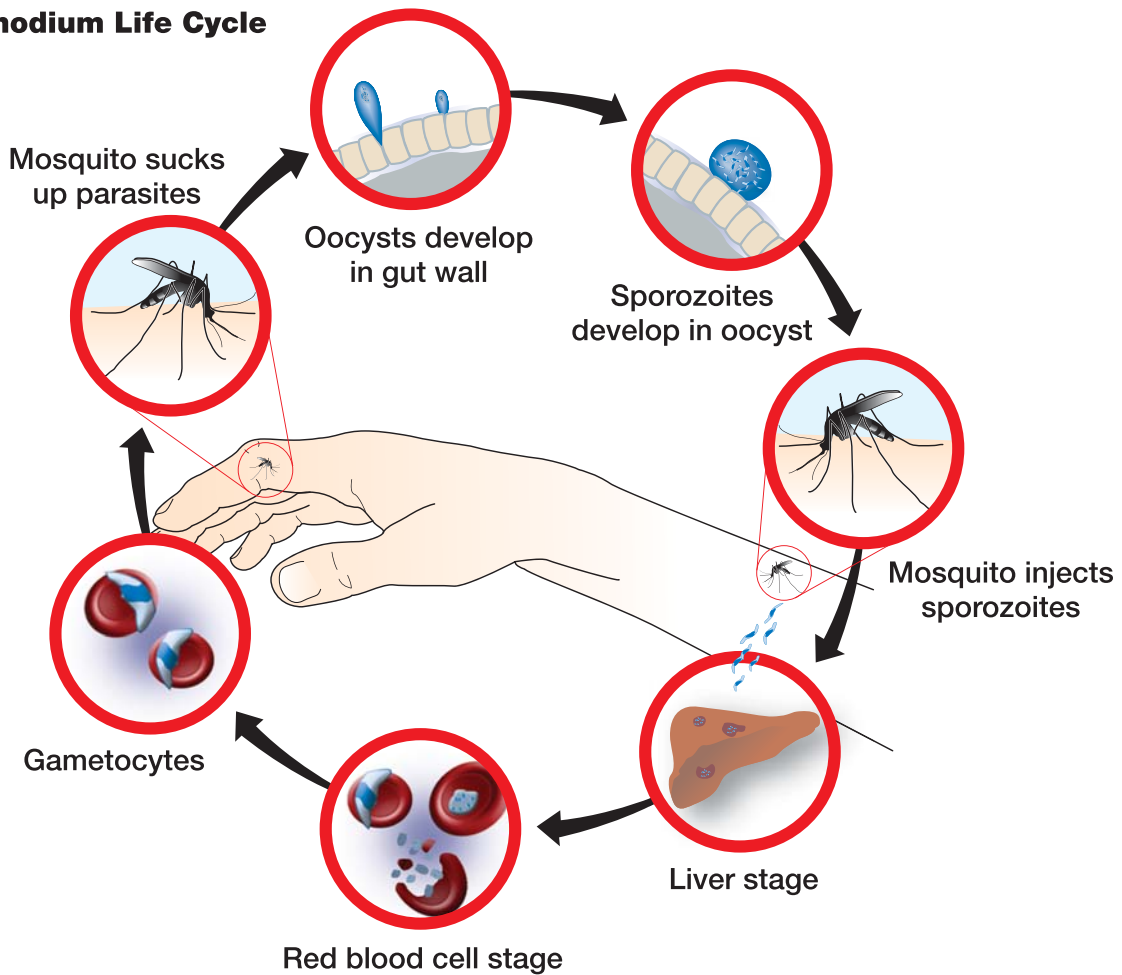
As mentioned, *Plasmodium* is mainly transmitted between infected humans by mosquitoes. Specifically, it is spread only by female mosquitoes of the genus *Anopheles*. This is because only female mosquitoes bite humans. They ingest human blood to obtain the proteins necessary for egg development. Most female *Anopheles* mosquitoes are nocturnal feeders (they bite only at night).

When a female mosquito bites and takes blood from a person infected with *Plasmodium*, the microscopic parasite moves along with the human's red blood cells to the mosquito's gut, where it continues through its life cycle. It then moves to the mosquito's salivary gland; when the mosquito takes a bite from another human, *Plasmodium* is injected along with the mosquito's saliva. (It is the proteins in mosquito saliva that trigger an immune response from your body, causing bites to itch.) The parasites can then be ingested by another mosquito, completing the life cycle and transmitting the disease from human to human to human.



**Blood smear showing the presence of the *Plasmodium* parasite (crescent shapes).** In the absence of more sophisticated tests, microscopic analysis of blood samples is a common diagnostic approach for malaria in clinics.

### Plasmodium Life Cycle



*Malaria is passed from person to person through mosquito bites.*

**Questions**

1. Malaria is most often transmitted by mosquitoes, but considering *Plasmodium* lives in erythrocytes, what are other ways in which the disease might spread from human to human?

2. Understanding the biology of malaria, it is not surprising that the most common methods of reducing outbreaks involve mosquito control. Relatively inexpensive and simple to perform, indoor spraying with insecticides kills mosquitoes for 3–6 months, and insecticidal bed nets provide additional protection from bites, with the insecticidal qualities of the nets lasting 3–5 years. What are the main benefits and drawbacks of both these approaches?

**Indoor spraying**

Benefits:

Drawbacks:

**Bed nets**

Benefits:

Drawbacks:

3. Malaria is only one of many parasite-mediated, mosquito-borne illnesses affecting the world's population. For this reason, some researchers advocate mosquito elimination — completely killing off mosquito populations in affected regions. The benefits of elimination are obvious (many of us would love to see mosquito-free lakes, streams, and ponds). However, what might some of the concerns be in eliminating a species from a habitat? Do we know what roles mosquitoes play in the environment? Does it make a difference if we know or don't know?

4. Many regions in Africa are reporting increasing populations of *Anopheles* mosquito species that show resistance to insecticides. Even more troubling are findings that climate change is expanding the habitat of the mosquitoes into regions where malaria had not been a health concern. How does your answer to the previous question change if the mosquito in question is an invasive species (for example, in the spread of dengue fever, the mosquito vector in question is often an invasive species new to habitats in North America)?

### **III. A Novel Approach Involves Bacterial Transformation**

Biotechnology and genetic engineering methods are also being investigated as mechanisms for eliminating malaria. Theoretically, any of the species involved — the *Plasmodium* parasite or the human or mosquito host — can be the target of genetic modifications that disrupt either the life cycle or transmission of the parasite. Practically, however, these systems have their drawbacks in terms of their ability to be either cultured or manipulated genetically. Therefore, researchers have turned their attention to a more familiar subject: bacteria.

#### **How can DNA “transform” bacteria? A background to bacterial transformation**

In the 1920s, scientists demonstrated how to turn a harmless strain of bacteria into a virulent strain, just by mixing the two strains together (Griffith 1928). What is truly incredible about this experiment is that the virulent strain had been killed prior to mixing, so something in the dead bacteria could “transform” the harmless bacteria, making them virulent.

It wasn't until the 1940s that scientists understood the chemical basis for this transformation. A team of scientists led by Oswald Avery at the Rockefeller Institute found that an extract of the bacteria was unaffected by treatment with protein-digesting enzymes, but was destroyed by a DNA-digesting enzyme. This showed that the agent that transformed the harmless bacteria was DNA (Avery et al. 1944).

Today, we understand that genes within DNA encode proteins that give rise to certain traits. We also know how to exploit the fact that many bacteria can acquire new genes by taking up DNA molecules encoding those genes (for instance, a plasmid) from their surroundings. The process is optimized by adding salts to the transformation medium and using a heat shock step, steps we use deliberately to transform bacteria and other microorganisms. The ability to transform the bacterium *E. coli*, for example, has made possible the cloning of genes, the cornerstone of many modern advances in sciences and of the biotechnology industry.

So how does bacterial transformation relate to our battle against malaria?

#### **Mosquitoes also have gut microbiota?**

It is surprising to many that, like humans, mosquitoes harbor a number of symbiotic bacteria within their gut. These symbiotic bacteria can be engineered, using procedures like those you used to transform *E. coli* bacteria, to produce proteins. However, in this case the symbiotic bacteria can be engineered to produce and secrete proteins that interfere with the life cycle of the *Plasmodium* parasite.

In one experiment (Wang et al. 2012), researchers used a bacterium called *Pantoea agglomerans*, which grows abundantly inside *Anopheles* mosquitoes. *P. agglomerans* can be grown and transformed using the same culturing and transformation techniques used with other more common bacteria, like *E. coli*. Researchers used these techniques to engineer *P. agglomerans* to express the genes of the hemolysin (hly) A system of *E. coli* bacteria, three proteins that cause red blood cells to lyse. The transformed bacteria were fed to mosquitoes through sugar solutions. The idea was that when the transformed bacteria colonized the mosquito gut, they would produce the toxic proteins. If that host mosquito then fed upon a human infected with malaria, the toxins produced by the transformed bacteria would cause the red blood cells (from the human blood) to burst. This would halt the life cycle of the *Plasmodium* parasite and stop the spread of malaria.

However, the researchers also hypothesized that transformation and expression of foreign genes might affect the ability of the transformed *P. agglomerans* bacteria to grow or colonize the mosquito gut (in other words, their fitness for that environment might be reduced). This could jeopardize the effectiveness of this strategy for fighting malaria in the wild. So they carried out another experiment: they transformed the bacteria with a plasmid that contains a green fluorescent protein (GFP), derived from the jellyfish *Aequorea victoria*, and fed the transformed bacteria to mosquitoes. The researchers then monitored how much fluorescence came from the mosquito gut. They found that after the host mosquitoes were given a blood meal, the GFP fluorescence in their guts increased, indicating the number of transformed bacteria there had rapidly increased. This demonstrated that transformed *P. agglomerans* could grow in the mosquito gut and, more importantly, replicate quickly when the mosquito ingested a blood meal. This meant the bacteria would also likely produce more hly A proteins when the host mosquito ingested potentially infected blood cells. In terms of the efficacy of the transformed bacteria against the *Plasmodium* parasite, when mosquitoes with the transformed bacteria were fed a blood meal containing the *Plasmodium* parasite, the development of the parasite was inhibited by nearly 98% (Wang et al. 2012).



## Questions

1. Scientists have also demonstrated that it is possible to genetically modify the *Anopheles* mosquito to produce a substance in their gut that kills off *Plasmodium*. Why might symbiotic bacteria be a more suitable subject for transformation than the *Anopheles* mosquito?
2. As in the human gut, many different bacterial species inhabit the *Anopheles* gut. If you were the researcher, how would you pick the best species for use in this transformation experiment? What factors should you consider?
3. Why was GFP used in the bacterial transformation experiment?
4. *E. coli* was also used in this experiment for plasmid production (to grow more copies of the plasmids), and the plasmid also contained genes for antibiotic resistance. After transformation, the bacteria were grown on plates with antibiotic in them. Why do you think this is a common step in bacterial transformation?
5. The experiment demonstrated that in the lab the transformed bacteria could survive and proliferate within the mosquito gut after transformation and that they could inhibit *Plasmodium* growth and development by nearly 98%. What other experiments might be needed to demonstrate this is a viable option for malaria elimination in the wild? Consider the life cycles and roles of all the key players in the spread of disease.
6. The hly A system used in the experiment described causes lysis of red blood cells. Considering this, would you have concerns about releasing these transformed bacteria into the environment? Under what conditions might these concerns be alleviated?
7. As a final thought, what do you think the greatest hurdles will be to successfully implement this bacteria-based approach in the wild? What are the technical, ethical, or regulatory challenges, and how might they be handled?

#### **IV: Prognosis**

Baruti opens his eyes and sees his mother sitting at his side. She takes his hand and tells him she loves him. She has waited for this moment for three days. She is exhausted from the sleepless nights and constant worry.

The doctors come into the room to check on his progress. They are cautiously optimistic that the mixture of antimalarial drugs is working, and they tell Lerato to be patient. They tell her that her boy was very ill and that treatment takes time, but they are encouraged by the progress he is making. They expect him to recover.

#### **Questions**

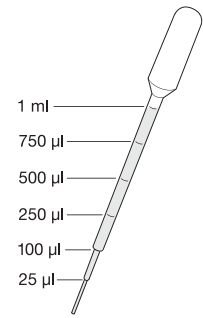
1. Most antimalarial drugs target the red blood cell (erythrocytic) stage of malaria infection, which is the phase of infection that causes symptomatic illness. Why might it be important to research other medications targeting other stages (for example, the liver stage) of the life cycle? Refer to the figure describing the life cycle of *Plasmodium*.
2. When *Plasmodium* becomes resistant to antimalarial drugs, this results in a delayed or incomplete clearance of the parasite from the patient's blood. How might an organism develop resistance to a chemical that can otherwise kill it?
3. The problem of antimalarial drug resistance can be compounded by cross-resistance, in which resistance to one drug confers resistance to other drugs that belong to the same chemical family or have similar modes of action. How and why do you think this might happen?
4. Current practice in treating cases of malaria is based on combination therapy, in which several different classes of drugs are combined. What might some advantages of this approach be?
5. Many people take antibiotics to treat bacterially mediated illnesses like strep throat or sinus infections. When you take antibiotics, you are told you must take the entire course of the medication in order to reduce the risk of developing antibiotic resistance. Why is a full course needed?

# Appendix A: Serial Dilution and Dilution Instructions

## How to Perform Serial Dilutions

A serial dilution is a series or chain of dilutions of a substance in a solution, typically distilled or sterile water. Serial dilutions are used to create accurate, highly diluted solutions intended to result in a concentration change. Many of the serial dilutions that are needed in the pGLO Transformation and Inquiry Kit will start from an original stock solution and will be successively diluted to create the desired concentration(s).

Serial dilutions require a degree of precision that is difficult to obtain using a DPTP, but not impossible. The dilution series depicted here can be done with a DPTP using the following gradient patterns.



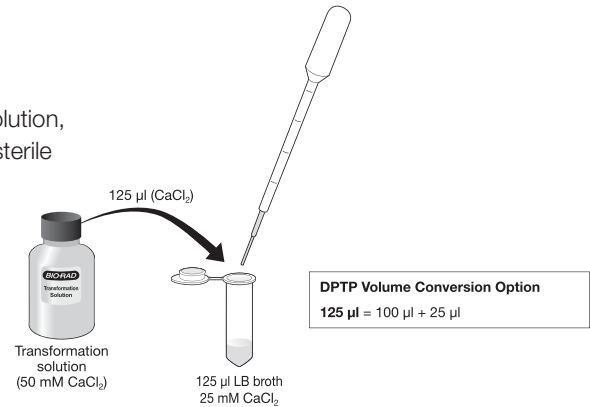
## Inquiry Investigation #1: Transformation Efficiency

### Transformation solution (CaCl<sub>2</sub>) concentration

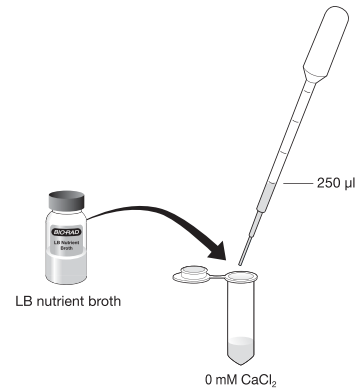
If you have chosen to vary the concentration of CaCl<sub>2</sub> transformation solution, you will need to make either a dilution of transformation solution using sterile water or a 0 mM solution using LB nutrient broth.

To make experimental dilutions of CaCl<sub>2</sub> you will use transformation solution and LB nutrient broth.

- a. For a 25 mM dilution of CaCl<sub>2</sub> use a sterile DPTP to transfer 125 µl of 50 mM CaCl<sub>2</sub> into a tube of 125 µl of LB broth.



- b. For a 0 mM dilution of CaCl<sub>2</sub> use a sterile DPTP to transfer 250 µl of LB broth into a tube.

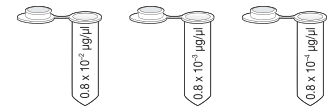


### Amount of pGLO plasmid DNA

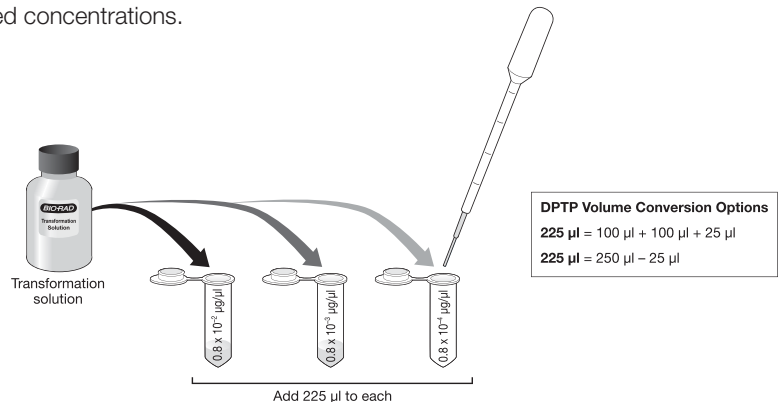
#### Transforming with less pGLO plasmid DNA

To make experimental dilutions of the pGLO plasmid DNA use the pGLO stock (0.8 × 10<sup>-1</sup> µg/µl) and transformation solution. A sample dilution series is illustrated.

- a. Use the rehydrated stock pGLO plasmid DNA (0.8 × 10<sup>-1</sup> µg/µl)
- b. Label blank microcentrifuge tubes with the expected concentrations.



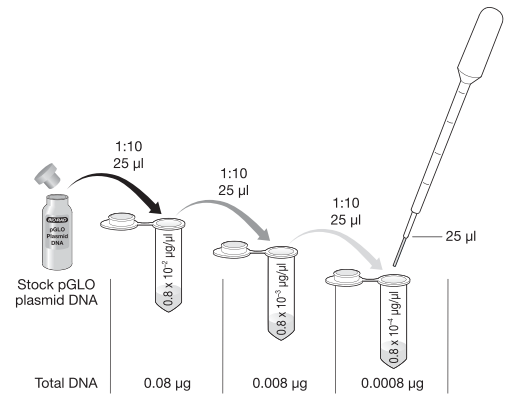
- c. Using a sterile DPTP, transfer 225 µl of transformation solution (the diluent for pGLO plasmid DNA) to the tubes.



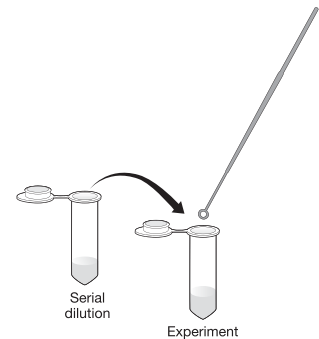
- d. Transfer 25  $\mu\text{l}$  of the stock pGLO plasmid DNA ( $0.8 \times 10^{-1} \mu\text{g}/\mu\text{l}$ ) to the first dilution, flick the tube to mix.

Then transfer 25  $\mu\text{l}$  of the first dilution ( $0.8 \times 10^{-2} \mu\text{g}/\mu\text{l}$ ) to the second dilution, flick the tube to mix.

Continue making serial dilutions until the series is complete. Use a new sterile DPTP or sterile pipet tip every time.



- e. Choose which serial dilution will be used as the experiment variable and use an inoculation loop to transfer 10  $\mu\text{l}$  into the experiment tube.



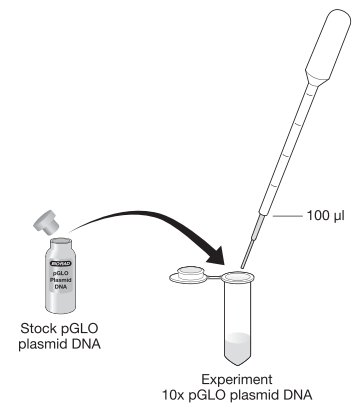
**Transforming with more pGLO plasmid DNA (micropipet is required)**

If you want to investigate the effect of transforming bacteria with more pGLO plasmid DNA (rather than less, in the form of a dilution), you must use a larger volume of pGLO plasmid DNA stock solution relative to the total volume in your experimental tube. This means you will need to reduce the amount of transformation solution and increase the amount of pGLO plasmid DNA stock solution. Following is an example of how to make a pGLO plasmid DNA stock that is ten times more concentrated than the standard (1x) pGLO plasmid DNA solution.

- a. Use a micropipet to transfer 160  $\mu\text{l}$  of transformation solution into the experiment tube. Use a 20–200  $\mu\text{l}$  micropipet to perform this accurately.



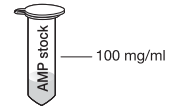
- b. Transfer 100  $\mu\text{l}$  of the pGLO plasmid DNA stock ( $0.8 \times 10^{-1} \mu\text{g}/\mu\text{l}$ ) into the experiment tube to make 10x pGLO plasmid DNA solution.



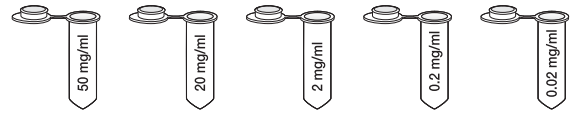
## Inquiry Investigation #2: Effect of Ampicillin on Bacterial Growth

In order to test different concentrations of the antibiotic ampicillin, you will need to make serial dilutions from a 100 mg/ml ampicillin (amp) stock solution. Set up dilutions of amp as described:

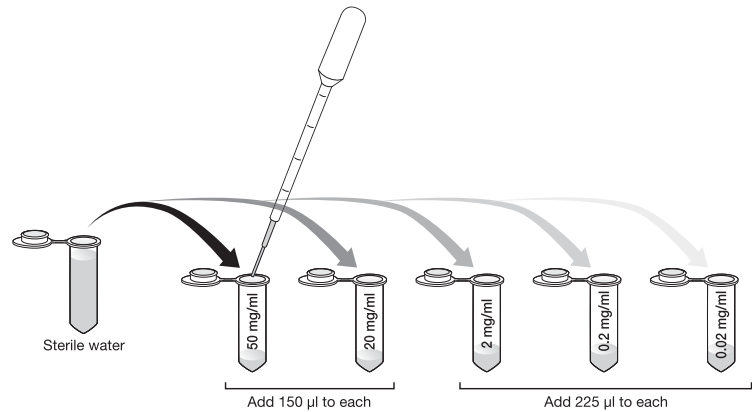
a. You should have a 100 mg/ml stock solution of ampicillin.



b. Label blank microcentrifuge tubes with the appropriate concentrations (for instance, 50 mg/ml, etc.).



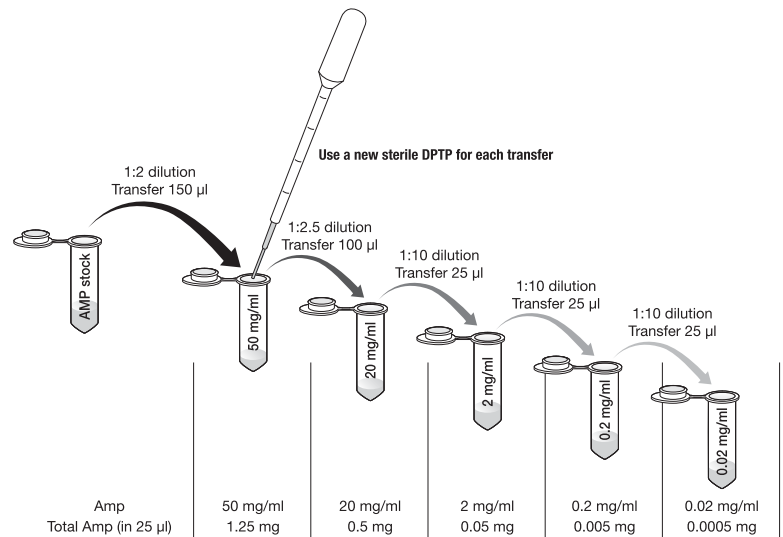
c. Using a sterile DPTP, transfer 150 or 225  $\mu$ l of sterile water (the diluent for ampicillin) to your tubes as shown.



d. Transfer 150  $\mu$ l of ampicillin from the stock tube to the first dilution as shown. Flick the tube to mix.

e. Transfer 100  $\mu$ l of ampicillin from the first dilution to the second dilution. Flick the tube to mix.

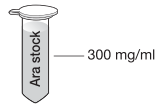
f. Continue making serial dilutions until your series is complete. Use a new sterile DPTP every time. Label each DPTP with the appropriate concentration for use later.



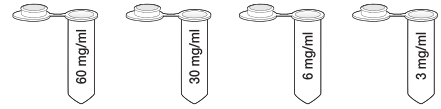
### Inquiry Investigation #3: Effect of Arabinose on GFP Expression

In order to test different concentrations of the inducer arabinose you will need to make serial dilutions from a 300 mg/ml arabinose (ara) stock solution. Set up dilutions of ara as described:

a. You should have a 300 mg/ml stock solution of arabinose.

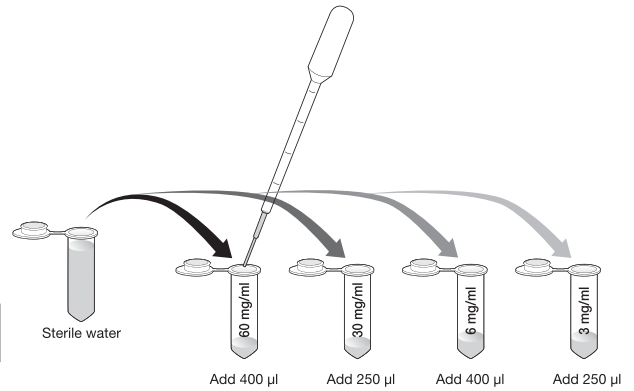


b. Label blank microcentrifuge tubes with the appropriate concentrations (for instance, 60 mg/ml, etc).



c. Using a sterile DPTP transfer 400 or 250 µl of sterile water (the diluent for arabinose) to your tubes.

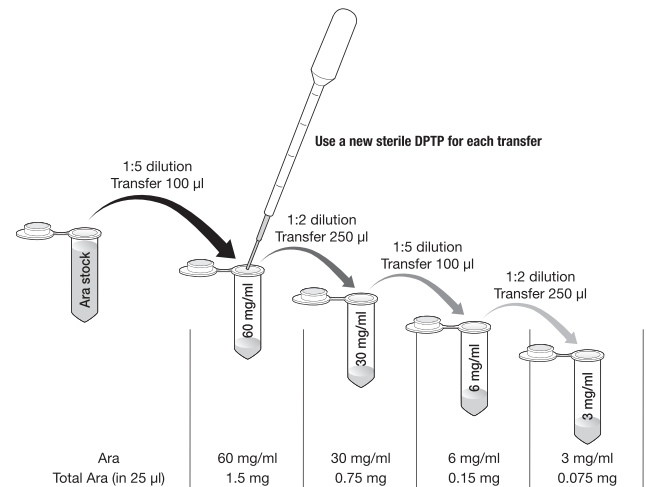
**DPTP Volume Conversion Option**  
 $400\ \mu\text{l} = 100\ \mu\text{l} + 100\ \mu\text{l} + 100\ \mu\text{l} + 100\ \mu\text{l}$



d. Transfer 100 µl of arabinose from the stock tube to the first dilution as shown. Flick the tube to mix.

e. Transfer 100 µl of arabinose from the first dilution to the second dilution. Flick the tube to mix.

f. Continue making serial dilutions until your series is complete. Use a new sterile DPTP or sterile pipet tip every time. Label each DPTP with the appropriate concentration for use later.



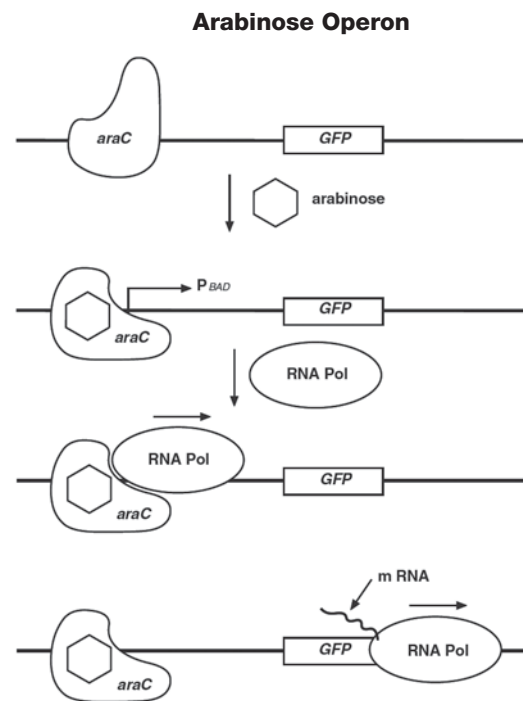
## Appendix B: Gene Regulation

Our bodies contain thousands of different proteins, which perform many different jobs. Digestive enzymes are proteins; some of the hormone signals that run through our bodies and the antibodies protecting us from disease are proteins. The information for assembling a protein is carried in our DNA. The section of DNA that contains the code for making a protein is called a gene. There are 30,000–100,000 genes in the human genome. Each gene codes for a unique protein. The gene that codes for a digestive enzyme in your mouth is different from one that codes for an antibody or the pigment that colors your eyes.

Organisms regulate expression of their genes and ultimately the amounts and kinds of proteins present within their cells for a myriad of reasons, including developmental changes, cellular specialization, and adaptation to the environment. Gene regulation not only allows for adaptation to differing conditions, but also prevents wasteful overproduction of unneeded proteins, which would put the organism at a competitive disadvantage. The genes involved in the transport and breakdown (catabolism) of food are good examples of highly regulated genes. For example, the sugar arabinose is both a source of energy and a source of carbon. *E. coli* bacteria produce three enzymes (proteins) needed to digest arabinose as a food source. The genes that code for these enzymes are not expressed when arabinose is absent, but they are expressed when arabinose is present in their environment. How is this so?

Regulation of the expression of proteins often occurs at the level of transcription from DNA into RNA. This regulation takes place at a very specific location on the DNA template, called a promoter, where RNA polymerase sits down on the DNA and begins transcription of the gene. In bacteria, groups of related genes are often clustered together and transcribed into RNA from one promoter. These clusters of genes controlled by a single promoter are called operons.

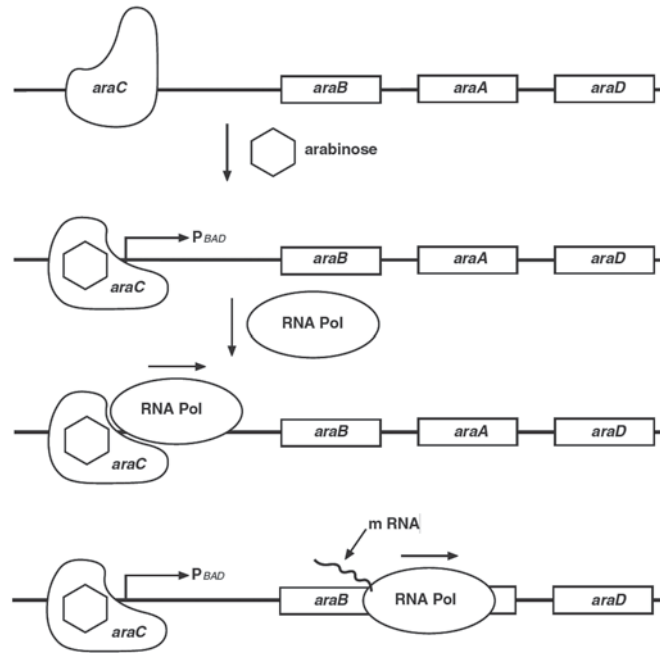
The three genes (*araB*, *araA*, and *araD*) that code for three digestive enzymes involved in the breakdown of arabinose are clustered together in what is known as the arabinose operon. These three proteins depend on initiation of transcription from a single promoter,  $P_{BAD}$ . Transcription of these three genes requires the simultaneous presence of the DNA template (promoter and operon), RNA polymerase, a DNA binding protein called *araC*, and arabinose. The *araC* binds to the DNA at the binding site for the RNA polymerase (the beginning of the arabinose operon). When arabinose is present in the environment, bacteria take it up. Once inside, the arabinose interacts directly with *araC*, which is bound to the DNA. The interaction causes *araC* to change its shape, which in turn promotes (actually helps) the binding of RNA polymerase, and the three genes *araB*, *araA*, and *araD* are transcribed. Three enzymes are produced, they break down arabinose, and eventually the arabinose runs out. In the absence of arabinose, the *araC* returns to its original shape and transcription is shut off.



The DNA code of the pGLO plasmid has been engineered to incorporate aspects of the arabinose operon. Both the promoter ( $P_{BAD}$ ) and the *araC* gene are present. However, the genes that code for arabinose catabolism, *araB*, *araA*, and *araD*, have been replaced by the single gene that codes for GFP. Therefore, in the presence of arabinose, *araC* protein promotes the binding of RNA polymerase and GFP is produced. Cells fluoresce brilliant green as they produce more and more GFP. In the absence of arabinose, *araC* no longer facilitates the binding of RNA polymerase and the GFP gene is not transcribed. When GFP is not made, bacteria colonies will appear to have a wild-type (natural) phenotype of white colonies with no fluorescence.

This is an excellent example of the central dogma of biology in action: **DNA > RNA > PROTEIN > TRAIT**

### Expression of Green Fluorescent Protein





### References

Avery OT et al. (1944). Studies on the chemical nature of the substance inducing transformation of pneumococcal types: Induction of transformation by a desoxyribonucleic acid fraction isolated from *Pneumococcus* Type III. *J Exp Med* 79, 137–158.

Griffith F (1928). The significance of pneumococcal types. *J Hyg* 27, 113–159.

Wang S et al. (2012). Fighting malaria with engineered symbiotic bacteria from vector mosquitoes. *Proc Natl Acad Sci* 109, 12734–12739.

### Suggested Reading, Listening

Kill 'Em All (NPR Radiolab podcast) <http://www.radiolab.org/story/kill-em-all/>

What if We Don't Kill 'Em All (NPR Radiolab podcast) <http://www.radiolab.org/story/what-if-we-dont-kill-em-all/>

### Legal Notice

Liofilchem is a trademark of Liofilchem S.R.L., LLC.



**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 424 6723 Australia 61 2 9914 2800 Austria 43 1 877 89 01 Belgium 03 710 53 00 Brazil 55 11 3065 7550  
Canada 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00  
France 01 47 95 69 65 Germany 49 89 31 884 0 Greece 30 210 9532 220 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300  
Israel 03 963 6050 Italy 39 02 216091 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666  
New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04  
Singapore 65 6415 3188 South Africa 27 (0) 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 026 674 55 05  
Taiwan 886 2 2578 7189 Thailand 1800 88 22 88 United Kingdom 020 8328 2000