



# Bio-Rad Science Teacher Grants Program

For science that goes beyond the textbook!





# Science Teacher Grants Program Overview

*As part of Bio-Rad's Community Outreach effort, the Science Teacher Grants Program is dedicated to enhancing science education in public schools in our local community at the K-12 level. The program was created to support innovative and creative approaches to science education and provide students with the opportunity to learn about and experience research techniques in a hands-on way.*



## Funding Priorities

The Science Teacher Grants Program encourages sustainable projects that improve the way science is taught at the K–12 level. Priority is given to requests involving Bio-Rad products and technologies provided by the company's Biotechnology Explorer™ program\* and/or involve Bio-Rad employees. Priority is also given to inquiry-based approaches to science education.

*\* Bio-Rad's Biotechnology Explorer program offers research-grade equipment and inquiry-based educational kits with curricula at a discount to educators. For more information please visit or call 1 800-4BIORAD (1 800-424-6723). Catalogs are available upon request.*

## Grant Timeline and Deadlines

- Teachers may apply for grants of up to \$5,000; this amount may be cash, Bio-Rad products, or combination of the two (cash and Bio-Rad products).
- The deadline for applications is April 15 (applications must be postmarked on or before April 15).
- Applicants who have been selected to receive a grant will be notified by May 31.
- Funds are distributed in the fall of the same year the application was submitted.\*\*

*\*\*Due to unforeseen budget considerations for the company during the year applications are submitted, the applications may be held over until the following year.*

## Eligibility of the Bio-Rad Science Teacher Grants

- Only public schools (and organizations) are eligible for the grants; the school must be located within a 50 mile radius of Bio-Rad's headquarters in Hercules, California.
- Only one grant will be awarded per school, per year.
- The grants may *not* be used toward general operations of a school, school maintenance, capital projects, endowments, conferences, independent study, fundraising events (or similar activities), salaries, textbooks, field trips, research, or for any type of religious or political causes.
- Organizations that discriminate on the basis of sex, race, age, disability, or religion are not eligible for the grants.



## Instructions For Preparing Your Application

Please prepare your grant application according to the instructions provided below. Teachers must be involved in the planning of projects. There are 4 sections of the application. Please submit your application according to the following format:

### Section 1: Application Cover Page

Complete form and attach evidence that the school (or organization) is public.

### Section 2: Project Summary

Briefly summarize your proposal and list its two most important learning objectives.

### Section 3: Endorsement

Have your proposal endorsed by a senior school official at the level of assistant principal or above.

### Section 4: Proposal

Create your proposal according to the format provided below. Your answers to sections (a-f) should be clear and succinct. Your proposal should not be longer than 3 pages, typed single space, using a 12 point font size. When you have completed your proposal, attach it to your Application Cover Page.

#### a. Project Description

Include a title for your proposal and a brief description of how you plan to use the grant. Indicate if other teachers or individuals are involved in your project; if so, please provide their names, positions, and roles in the project.

#### b. Learning Objectives

Describe the learning objectives for the students that will be met by your proposal. The objectives should be realistic, directly related to the project, and attainable within the period of time of the project.

#### c. Teaching Strategies

Describe how you propose to meet the learning objectives stated above. Please also provide proposed teaching strategies, the duration of the project, and the classroom time and resources that will be dedicated to the project. How will your project differ from current teaching strategies?

#### d. Experience/Additional Resources

Do you (or any of the other individuals involved in your project) have any additional resources, experience, or training that will ensure the success of the project? For example: professional development, resources already procured, funding commitments from other sources, etc.

#### e. How You Will Determine Success

Upon completion of the project, describe how you will determine if your project met its objectives and was a success.

#### f. Appendices

Please feel free to include any additional information that you feel is relevant to your proposal in this section. For example, you may wish to include letters of recommendation or support for your proposal. Please keep this supplementary material to a minimum.

### Section 5: Proposal Budget

Use this section to explain specifically how the funds will be used for your project.

### Section 6: Project Success Report

This section is to be completed only by those who receive grants. Please provide this Project Success Report within 30 days after the completion of your project.

## Grant Application Submission

Please submit your application along with the required Sections; be sure to obtain necessary endorsements (signatures) requested in Section 3 and mail to:

#### Bio-Rad Laboratories

Attn: Community Outreach  
1000 Alfred Nobel Drive  
Hercules, CA 94547

If you have any questions about this application or the submission process, please feel free to contact us at [community\\_outreach@bio-rad.com](mailto:community_outreach@bio-rad.com).



# Section 1: Application Cover Page

(Note: it may be easier to complete the Application Cover Page once you have completed the other sections of the application.)

<b>PROJECT TITLE (MAXIMUM 140 CHARACTERS INCLUDING SPACES)</b>		
<b>PROJECT MANAGER (NAME/TITLE)</b>		
<b>SCHOOL OR ORGANIZATION NAME</b>	<b>SCHOOL DISTRICT (IF APPLICABLE)</b>	
<b>ADDRESS</b>		
<b>CITY</b>	<b>STATE</b>	<b>ZIP</b>
<b>HAVE YOU RECEIVED A GRANT/SPONSORSHIP FROM BIO-RAD IN THE PAST? IF SO, PLEASE EXPLAIN.</b>		
<b>THE NUMBER OF MILES FROM YOUR SCHOOL/ORGANIZATION TO BIO-RAD'S HEADQUARTERS IN HERCULES, CALIFORNIA*</b>		
<b>PHONE (WORK)</b>	<b>PHONE (AFTER SCHOOL OR WHEN SCHOOL IS NOT IN SESSION)</b>	
<b>EMAIL ADDRESS</b>	<b>ALTERNATE EMAIL (WHEN SCHOOL IS NOT IN SESSION)</b>	
<b>SCHOOL/ORGANIZATION ADMINISTRATOR (NAME/TITLE)</b>	<b>ADMINISTRATOR'S PHONE</b>	
<b>GRANT AMOUNT REQUESTED</b>	<b>NUMBER OF STUDENTS WHO WILL BENEFIT FROM YOUR PROJECT</b>	
<b>GRADE LEVEL</b>	<b>PROJECT BEGINNING &amp; END DATES</b>	

Only public schools and organizations are eligible to apply for Bio-Rad's Science Teacher Grants. Applicants must provide evidence that their school is public as recognized by an appropriate local or state government agency. For example, a copy of the state tax exempt certificate could be used to satisfy the eligibility requirement.

We confirm we are a public school or organization and have attached proof with this application.

\* Your school must be located within a 50 mile radius of Bio-Rad's headquarters in Hercules, California.



# Section 2: Project Summary

Summarize your project idea in a short paragraph and list the two most important learning objectives (maximum 200 words):

# Section 3: Endorsement

The project described must be acknowledged and endorsed by a school official at the level of assistant principal or above.

I endorse this grant proposal, including the proposed budget.

<b>PROJECT MANAGER</b>	<b>DATE</b>
<b>SCHOOL OFFICIAL</b>	<b>DATE</b>
<b>SCHOOL OFFICIAL'S TITLE</b>	



## Section 4: Proposal

Use this section to write your proposal according to instructions provided on page 2 “Section 4: Proposal” (your answers to questions a-f). Your answers should be clear and succinct. Your proposal should be not be longer than 3 pages, single space text, using a 12 point font size. When you have completed your proposal, attach it to your Application Cover Page.

## Section 5: Proposal Budget

The Bio-Rad Science Teacher Grants Program is dedicated to promoting sustainable science projects for the classroom that may be incorporated into the school’s science education curriculum for years to come. We understand that to begin a project some training and consumables may be required. However, we request that the majority of the grant funds be used toward equipment that has sustainability and may be used for science education on a continuing basis. Grant applicants that have sustainability built into their projects are highly encouraged. Please base your cost estimates (requested below) on current market value of equipment (obtained from vendors, catalogs, the internet, etc.); be sure to ask suppliers if they offer special pricing for teachers. If you would like to receive a Bio-Rad quote for the company’s products that you are requesting to help with your budget, please call 1-800-4BIORAD (1-800-424-6723) for pricing.

<b>PROJECT TITLE</b>
<b>PROJECT MANAGER (NAME/TITLE)</b>

Provide a cost estimate for each area in which you seek funds (attach additional pages if necessary); indirect charges are not allowed).

### Costs of your Proposal

ITEM:	REQUESTED AMOUNT:	DISTRICT LOCAL MATCH (CASH OR IN-KIND) IF AVAILABLE:	PRIVATE, PUBLIC, NON-PROFIT SECTORS MATCH (CASH OR IN-KIND) IF AVAILABLE:	ITEM TOTAL:
EQUIPMENT				
CONSUMABLES (MAXIMUM 25% OF TOTAL GRANT AMOUNT)				
PROFESSIONAL DEVELOPMENT (MAXIMUM 25% OF TOTAL GRANT AMOUNT)				
<b>TOTAL</b>				



## Budget Justification

*Please provide an explanation of your budget request, addressing proposed expenditures in each category. Specify any Bio-Rad products by catalog number (maximum 500 words)..*



# Section 6: Project Success Report

In this section, please describe the outcome of your project. Please send your response to Bio-Rad within 30 days after your project is complete. If you need more room, feel free to attach additional sheets as necessary.

<b>PROJECT TITLE</b>
<b>PROJECT MANAGER (NAME/TITLE)</b>

<b>1) BRIEFLY SUMMARIZE THE OUTCOME OF YOUR PROJECT</b>
<b>2) WOULD YOU CONSIDER YOUR PROJECT A SUCCESS? WHY OR WHY NOT?</b>
<b>3) WILL THE PROJECT CONTINUE? IF NO, WHY NOT? IF YES, PLEASE DESCRIBE</b>
<b>4) WHAT WAS THE HIGHLIGHT OF YOUR PROJECT EXPERIENCE?</b>
<b>5) IS THERE ANY STUDENT FEEDBACK THAT YOU WOULD LIKE TO SHARE?</b>

Please submit your Project Success Report to:

**Bio-Rad Laboratories**  
Attn: Community Outreach  
1000 Alfred Nobel Drive  
Hercules, CA 94547