



# Bio-Rad Scholarship Program





# Scholarship Program Overview

The Bio-Rad Scholarship Program is designed for graduating high school students who have an interest in pursuing a career in science. The objective of the program is to provide financial assistance to those deserving students who have applied themselves in school and have decided to seek higher education to pursue their science career goals. The program is available to those students whose high school is located within a 50-mile radius of the Bio-Rad's Corporate offices in Hercules. Applications are graded on the student's financial need, their personal connection to science, as well as the originality, grammar, and sense of organization that are reflected in the applicant's answers to the questions and their essay.

Two scholarships are available:

## Bio-Rad Scholarship: \$1,000

Multiple Bio-Rad Scholarships are awarded each year.

## Ron Mardigian Scholarship: \$2,500

Only one Ron Mardigian Scholarship is awarded each year.

All applicants are automatically considered for both scholarships although only one Ron Mardigian Scholarship is awarded each year to the one applicant that stands out above the rest. The Ron Mardigian Scholarship was created in the memory of the late Ron Mardigian, who spearheaded Bio-Rad's Biotechnology Explorer™ program that continues to inspire millions of students around the world to explore their interests in science.

## Eligibility Requirements

To be eligible for this award, the applicant must:

- Be a graduating, college-bound senior from a high school within a 50-mile radius of Bio-Rad's Corporate offices in Hercules
- Have a grade point average of 2.5 or higher
- Have already been accepted at a 2- or 4-year college at the time scholarships are awarded in April of each year

## How to Apply

To apply, please complete the Scholarship Eligibility Form on the following page and return it to Bio-Rad (find address on the Scholarship Eligibility Form).

Submissions must be postmarked by March 1 each year. The applicant must also include one copy of each of the following with his or her application:

- A sealed copy of your most recent school transcript
- Two letters of recommendation from a teacher, school counselor, or community leader (the applicant may use letters of recommendation used for his or her college applications)
- The Scholarship Eligibility Form
- An essay answering one of the questions below. The essay must be at least 400 words (one and a half pages), typed double-spaced on 8.5 x 11" white bond paper in a standard typeface

## Essay Questions

Select and answer one of the following questions:

- **What can be done to encourage interest and participation in science?**
- **How can science be used to improve our world today?**





# Scholarship Eligibility Form

Please check boxes below:

I have read the eligibility requirements and I meet the criteria.

I have submitted one copy of all requested documents:

- 1. The Scholarship Eligibility Form
- 2. My high school transcripts
- 3. Two letters of recommendation
- 4. My essay

I understand that meeting the stated eligibility requirements does not guarantee that I will be selected to receive a scholarship.

By signing and submitting my application packet, I certify that the information that I am providing is complete and accurate to the best of my knowledge.

### Mail the Scholarship Eligibility Form and all required documents to:

Bio-Rad Laboratories  
Attention: Community Outreach  
1000 Alfred Nobel Drive  
Hercules, California 94547

### Deadline is March 1

Those selected to receive a Bio-Rad Scholarship will be notified no later than April 15. A Scholarship Awards ceremony is usually held the last week in April or the first week in May.

Career interests \_\_\_\_\_

Extracurricular activities \_\_\_\_\_

Please note any challenges or obstacles you face that may affect furthering your education (optional) \_\_\_\_\_

*(Note: If you do not have enough room to answer any of these questions, feel free to continue your response on a separate sheet of paper.)*

Are you applying for financial aid?

Yes  No  I am not sure yet

How did you hear about the Bio-Rad Scholarship Program? \_\_\_\_\_

Name \_\_\_\_\_

Home address \_\_\_\_\_

City \_\_\_\_\_

State \_\_\_\_\_

Zip \_\_\_\_\_

Phone (please include area code) \_\_\_\_\_

Email address \_\_\_\_\_

Full name of the high school you attend \_\_\_\_\_

If applicable, your school counselor (name and phone number)

Universities and/or colleges in which you have applied \_\_\_\_\_

Your signature \_\_\_\_\_ Date \_\_\_\_\_