

Bio-Rad Scholarship Program

Scholarship Program Instructions and Application Form





Bio-Rad is a global life science research and clinical diagnostics company headquartered in Hercules, CA. More than 60 years after its founding, the company remains committed to its mission of advancing research and improving lives worldwide through innovative scientific products and services.

We recognize the importance of involvement in our local community, and support the scientific education of students who may become the scientists of tomorrow.

Scholarship Overview

The Bio-Rad Scholarship Program assists local graduating high school students who wish to pursue careers in science or engineering. The program provides financial assistance to deserving students so they can continue their pursuit of science and engineering in higher education. The program is available to students attending high school within the following counties: Solano County, Alameda County, and Contra Costa County.

Applications are evaluated based on students' personal connection to science and engineering, as well as the originality, clarity, and persuasiveness of their answers to essay questions.

Three levels of scholarships are awarded:

Bio-Rad Scholarship: \$2,000

20 Bio-Rad Scholarships are awarded each year.

Ron Mardigian Scholarship: \$3,000

One Ron Mardigian Scholarship is awarded each year.

David and Alice Schwartz Scholarship: \$3,000

One David and Alice Schwartz Scholarship is awarded each year.

One Ron Mardigian Scholarship and one David and Alice Schwartz Scholarship is awarded each year by Bio-Rad's scholarship review panel to two especially deserving students. The first of these scholarships was created in the memory of the late Ron Mardigian, who spearheaded Bio-Rad's educational program in the 1990s, while the second was created to honor David and Alice Schwartz, who founded Bio-Rad in 1952 to accelerate the scientific research process.



Eligibility Requirements

To be eligible for a scholarship, the applicant must:

- Be a graduating, college-bound senior from a high school within the following counties: Solano County, Alameda County, and Contra Costa County.
- Have a grade point average of 2.5 or higher
- Have already been accepted to a 2- or 4-year college at the time scholarships are awarded.

How to Apply

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

Submissions must be postmarked by March 15. Please include one copy of each of the following with your application:

	0	,		
The Scholarship Applica	ation F	orm		
A sealed copy of your n	nost re	cent scl	hool transcri	pt
Two letters of recomme counselor, or communit			a teacher, sc	hool
An essay answering one must be at least 400 wo single sided, on 8.5 x 1	ords, ty	ped, do	ouble-space	d an

Please identify which question you are answering.

Please do not use staples.

Essay Questions

Select and answer ONE of the following questions:

- How is science <u>or</u> engineering making life better for people across the globe right now? <u>Cite specific examples.</u>
- What do you think the next big scientific or engineering advancement will be, and how will it affect society globally?
 Please be as specific as possible.
- Who are your heroes in science <u>or</u> engineering?
 <u>Please describe their contributions and</u> explain their importance.

Scholarship Awards Ceremony

The Scholarship Awards Ceremony is an exciting event held to give recognition to students who are selected as scholarship recipients for their exemplary achievements. Scholarship awards are made based on a combination of factors, demonstrated academic excellence, student life involvement, or social contributions.

Family and friends of the recipients are also invited to attend the ceremony to share in the celebration with their student. The ceremony is held at Bio-Rad Corporate Headquarters in Hercules, California.

Bio-Rad Scholarship Home address Application City State____ Please check boxes below: I have read the eligibility requirements and I meet the criteria. Phone (please include area code) ■ I have submitted one copy of all requested documents: Email address 1. The Scholarship Application Form Full name of the high school you attend _____ 2. My high school transcripts 3. Two letters of recommendation 4. My essay (remember to indentify the essay question) ■ I understand that meeting the stated eligibility requirements If applicable, your school counselor (name and phone number) does not guarantee that I will be selected to receive a scholarship. By signing and submitting my application packet, I certify that the information I am providing is complete and accurate to the Universities and/or colleges to which you have applied ____ best of my knowledge. **Mail the Scholarship Application Form** and all required documents to: Bio-Rad Laboratories Attention: Community Involvement 245 Linus Pauling Drive, Suite B Hercules, CA 94547 **Deadline is March 15** Those selected to receive a Bio-Rad Scholarship will be notified approximately 2 weeks prior to the event. A Scholarship Awards ceremony will be held 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.) Bio-Rad Scholarships are not need-based, but financial need is considered in the evaluation. How did you hear about the Bio-Rad Scholarship Program? _____

______ Date

Your signature____