



# Give Students a Competitive Edge and Job-Ready Skills

## Biotechnology: Growing Employment Opportunities

Biotechnology applies biological molecules and processes for the improvement of human lives. It is at the core of innovation in many industries, including medicine, food processing, textiles, agriculture, energy, and environmental conservation. Biotechnology touches almost every aspect of our lives.

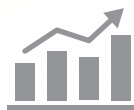
These life science and biotechnology industries are growing and continue to hire at a rapid pace, creating an ongoing need for talent. The time to educate your students and prepare them for this exciting future is now.

## Prepare Students by Partnering with Bio-Rad

With over 25 years of experience in science education and almost 75 years as a global biotechnology leader, Bio-Rad Laboratories can help you develop a 21st-century biotechnology course or program.

## Think Biotechnology. Think Bio-Rad.

Partner with us to create a robust and scalable curriculum that can adapt to your changing needs. Visit [bio-rad.com/explorer](http://bio-rad.com/explorer) for more information about the Bio-Rad Explorer program.



### Sustainable job growth

Opportunities expected to grow by at least 9% through 2029



### Competitive salaries

Almost twice the private sector national average



### College degree not required

Multiple entry points and opportunities for career advancement

## Industry-Quality Products That Work

Success creates confidence and ignites passion. This is why Bio-Rad partners with master teachers and researchers to create relevant and robust educational products that work. All products undergo rigorous quality control to guarantee reliability and longevity.



### Biotechnology: A Laboratory Textbook

Aligned to the Biotechnology Assistant Credentialing Exam (BACE), our textbook can guide development of a new course or supplement existing programs.



### Innovative Hands-On Lab Activities

From DNA extraction to protein purification and real CRISPR gene editing, our hands-on lab activities complement the textbook and are guaranteed to work — set your students up for success!



### Real, Industry Research Equipment

Students learn on the same instruments used in research centers and laboratories.



### Professional Development

Online or in-person, we offer training to match each teacher's background and training needs.

## Credentialing and CTSO Alignments

The Bio-Rad Explorer program partners with industry leaders to offer the alignments you need for a robust CTE program.



### Biotechnician Assistant Credentialing Exam (BACE)

We can help you develop a program that prepares students for this industry-recognized credential and gives them a competitive advantage when seeking employment — in industrial or in academic labs!



### HOSA: A Career and Technical Student Organization (CTSO)

The HOSA Competitive Event in Biotechnology uses the Bio-Rad textbook as a primary resource, allowing you to build a CTE program that aligns with this popular and impactful CTSO.

Visit [info.bio-rad.com/ExplorerContactUs](http://info.bio-rad.com/ExplorerContactUs) to connect with a curriculum training specialist near you.

## Program Solutions and Support

Whether you are looking to add just a taste of biotechnology or to start a full course or program, we work with you to tailor your curriculum to your needs.

**Flexible** — Use a modular approach to fit all school schedules and timeframes.

**Adaptable** — Customize your program to fit your vision for school, student, and community needs.

**Scalable** — Choose from simple integrations to multi-course certification pathways.

**Economical** — With generous discounts and no subscription costs, you can grow your program sustainably and within budget.

© 2023 Bio-Rad Laboratories, Inc. BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Website** [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23  
**France** 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050  
**Italy** 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 080 007 7373 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670  
**The Netherlands** 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23  
**Russian Federation** 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23  
**Switzerland** 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

