



Promoting Biotechnology Education Through Partnership

Bio-Rad and the InnovATEBIO National Biotechnology Education Center have partnered to help meet the growing need for a skilled technical workforce in the biotechnology industry.

Together, we aim to redefine biotechnology education by offering:

- Professional development (teacher workshops, webinars, and on-site technical training)
- Teaching resources (curricula, video tutorials, on-site presentations, and other instructional materials)
- Sponsorships and other types of support

Together, we leverage our unique expertise and strengths to:

- Identify and address emerging workforce and industry trends
- Ensure students have access to and learn to work with industry-based, research-quality equipment and supplies
- Build a broad community of educators, mentors, employers, and others throughout the biotechnology community

BIO-RAD Explorer

Bio-Rad + InnovATEBIO: Helping to meet the growing need for a skilled technical workforce in the biotechnology industry

The InnovATEBIO Center

Funded by the National Science Foundation in 2019,¹ InnovATEBIO helps educate highly skilled technicians for the nation's biotechnology workforce. By supporting educational initiatives and strengthening the connections between high school and community college biotechnology programs, InnovATEBIO strives to address the changing needs of the biotechnology community by:

- Developing a collaborative infrastructure that supports innovation
- Monitoring and addressing emerging trends in the biotechnology industry
- Coordinating and leveraging outputs from ATE-funded biotechnology projects
- Nurturing partnerships and collaborations that accelerate innovation in biotechnology education
- Engaging underrepresented populations in biotechnology labs and related emerging technologies

To learn more about InnovATEBIO, visit innovatebio.org

Bio-Rad Explorer Program

Bio-Rad Laboratories, a global leader in the development and manufacture of products for life science research and clinical diagnostics, understands its responsibility to support the education of a science-literate population. To meet this responsibility, the company established the Bio-Rad Explorer Program.

For more than 20 years, the mission of the Bio-Rad Explorer Program has been to make science accessible and to keep students at the center of instruction. By leveraging the expertise within Bio-Rad and working in partnership with educators, researchers, and industry leaders, the program brings state-of-the-art science and technology into the classroom. It offers:

- Hands-on laboratory activities that prepare students with foundational skills and expand their experience with cutting-edge techniques — all in a format that is appropriate for classroom laboratories
- Engaging lessons that connect techniques with scientific concepts and provide the context of real-world scenarios
- Industry-quality instruments and reagents — the same tools scientists use in their research
- Professional development opportunities and teaching resources that support teachers in their quest to spark interest in the life sciences
- Industry insights and decades of experience that can help shape biotechnology programs

To learn more about the Bio-Rad Explorer program, visit explorer.bio-rad.com.

¹ DUE 1901984

© 2020 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.



(L to R) **Dr. Molly Beth Malcolm**, Executive Vice President of Campus Operations and Public Affairs at Austin Community College; **Dr. Linnea Fletcher**, Department Chair of Biology at Austin Community College and NBEC Founder; **Andre Williams**, AP Biology Teacher; and **Dr. Ingrid Hermanson-Miller**, Bio-Rad Explorer Marketing Manager; meet at Anderson High School in Austin, TX.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website 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 2 3473 4460 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

