

**In partnership with InnovATEBIO**, the Bio-Rad Explorer program developed this Life Science Careers poster to help you recruit students into your science and biotechnology courses and programs. We hope that by sharing this poster, you can bring awareness to the breadth and depth of career opportunities that are possible for students with a life science or biotechnology background.

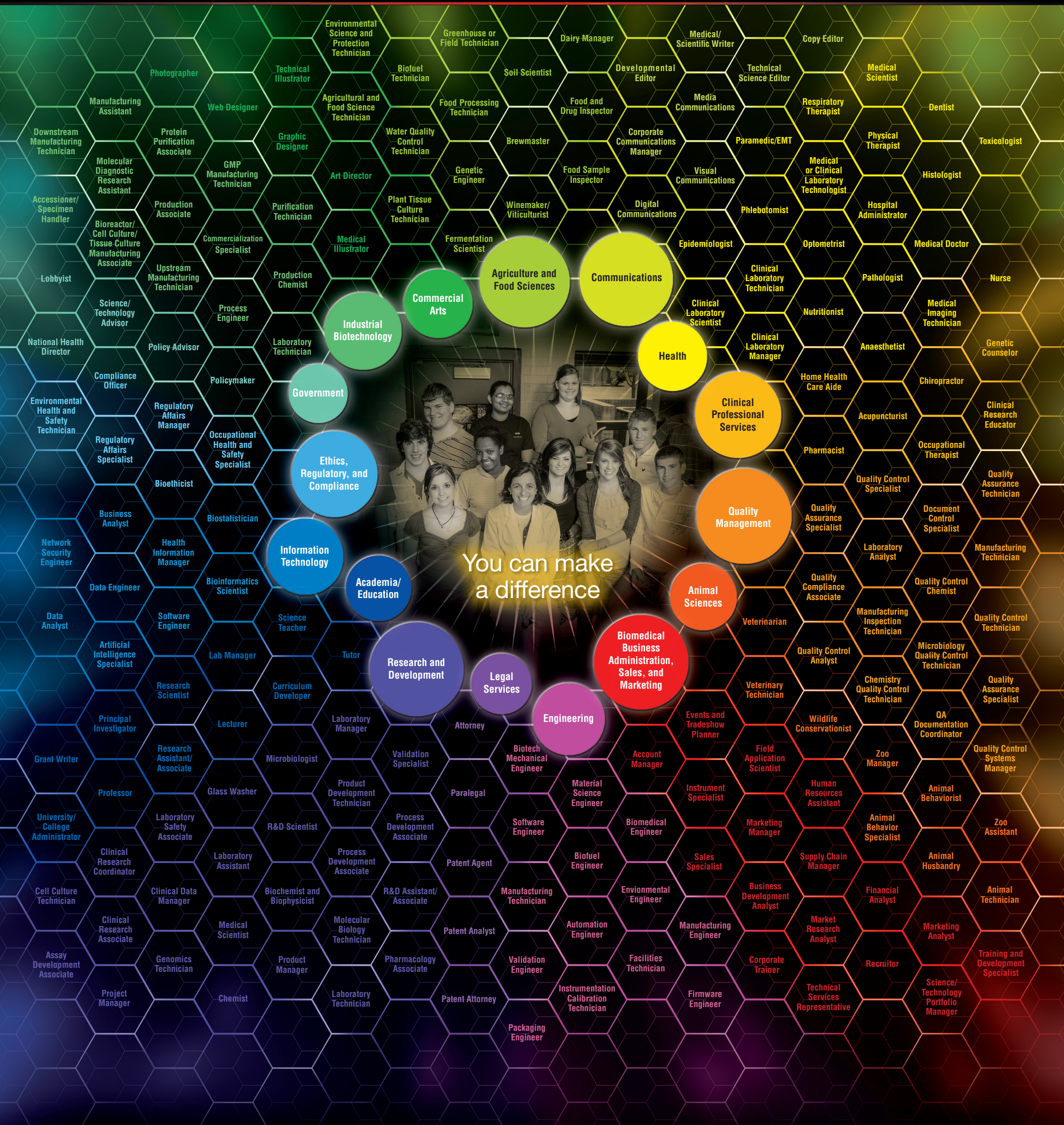
**Using Acrobat Reader to customize this poster for printing**, add your own logo, branding, and/or contact information into the editable boxes at the bottom of the PDF. (Poster optimum size: 11x17")





BE PART OF THE SOLUTION WITH A

# Life Science Career



You can make a difference



[bio-rad.com/explorer](http://bio-rad.com/explorer)



Made in partnership with InnovATEBIO and California Community Colleges. Visit [bio-rad.com/innovatebiocareers](http://bio-rad.com/innovatebiocareers) to learn more about these and other career options.

The Bio-Rad Explorer Program brings you the next generation of inquiry-based biotechnology curricula, kits, equipment, and supplies. 1-800-4BIORAD