

CSTA Regional Conference 2019

# Free Teacher Workshops

Presented by Bio-Rad Laboratories

Step-by-Step, Hands-on Experience

San Jose, CA

San Jose McEnergy  
Convention Center

October 18–20

Friday | October 18

Time	Room	Title	Description
8–9 AM	220C	<b>Algae Blooms: Agriculture, Ecology, and Economy</b>	Teach photosynthesis and cellular respiration together in the context of the Gulf of Mexico dead zone. Using algae beads together with an algae bloom case study, your students engage in inquiry investigations to learn about two connected processes and their ecological and economical implications.
12:30–2 PM	220C	<b>Learn to Infuse NGSS Science/ Engineering Practices with an Engaging Activity</b>	This workshop will focus on illustrating the science and engineering practices described in the NGSS framework through the engaging pGLO Bacterial Transformation activity.

Saturday | October 19

Time	Room	Title	Description
10:00–11:30 AM	LL20A	<b>Conserving Panda Populations Through Understanding Reproductive Endocrinology (Grades 9–16)</b>	Can your students save the Giant Pandas? Explore topics such as homeostatic regulation, reproductive hormones, immunological responses, and ecosystem balance all at once as students engineer a hormone detection system that can be used for Panda conservation.
3:00–4:30 PM	LL20D	<b>Think Like an Engineer in Your Biology Class</b>	Incorporate NGSS science and engineering practices in your biology class by engaging students to define the problem of world hunger. Considering constraints, students will design a treatment plan (solution), in the form of an evidence-based argument, for protein-energy malnutrition.
4:45–6:15 PM	LL20D	<b>Investigate Photosynthesis and Cellular Respiration with Algae Beads for GenBio</b>	Use algae beads in a colorimetric assay to study photosynthesis and cellular respiration through inquiry investigations. Learn how students can observe a phenomenon, ask meaningful questions, and design and carry out their own experiments to understand how these processes are connected.

Come visit us at **Booth #223**

Can't make a session? Don't miss out — download the FREE presentations from [bio-rad.com/explorerworkshops](http://bio-rad.com/explorerworkshops)

19-2177

**BIO-RAD**

Visit us on the Web at [explorer.bio-rad.com](http://explorer.bio-rad.com)  
or call us at 1-800-4BIORAD (1-800-424-6723)

**BIO-RAD**  
**Explorer**