



NABT Chicago Workshop Schedule & Special Events

Join us at the Sheraton Grand Chicago Hotel at the Colorado room, for our free workshops

Friday November 15 (all at the Colorado room)

- | | |
|----------------|---|
| 8:00–9:00 AM | Are Increased Incidences of Infection the Result of Climate Change? Why does climate change matter for human health? Discover why rates of harmful infections may become more common as earth's temperature increases. |
| 10:30–11:45 AM | The Opioid Epidemic: Decoding the Genetic Associations to Opioid Abuse. As the opioid crisis surges, researchers race to decode the genetics of opioid dependence. In this hands-on workshop, use fast DNA electrophoresis to evaluate genetic links and explore personalized medicine. |
| 12:00–12:30 PM | Ready or Not, It's Coming! Biotechnology, the Science of Our Age. Are Your Students Prepared? Glowing cats? Designer babies! Empower students to be independent thinkers. Learn from a leader in biotechnology teaching how to build your lab program step-by-step with equipment, supplies, and student credentials. |
| 1:00–1:45 PM | Think Like an Engineer in Your Biology Class. Incorporate NGSS engineering practices in your biology class by challenging students to address world hunger. Students will consider constraints and design an evidence-based treatment plan (solution) for protein-energy malnutrition. |
| 2:00–3:15 PM | It's in Their DNA! Teach Personalized Medicine with Students' Own DNA. Experience a hands-on classroom activity where students work with their own genes and PCR in the context of personalized medicine, the wave of future disease treatment. |
| 3:30–4:00 PM | Precision Medicine — a Reality with CRISPR and Revolutionary Droplet Digital PCR (ddPCR) Technology! ddPCR technology is a precision medicine tool and its sensitivity makes it well-suited to "Liquid Biopsies" to detect rare cancer mutations and when combined with CRISPR technology is revolutionizing medicine. |

Friday November 15 — 12:45–1:45 PM

Proud sponsor of the **Four-Year College & University Section Luncheon** (tickets required)

Special Events

Saturday November 16 (Chicago Ballroom 8) — 11:30 AM–2:00 PM

Lab Skills: **The Escape Room! Experience an escape room like no other.** Space is limited. Get tickets at the Bio-Rad booth.

Sunday November 17 (Erie Room) — 8:30–11:30 AM

Strategies for a **More Inclusive Biology Classroom** presented by iEMBER, NABT and Bio-Rad Laboratories

Visit Us
at Booth
#508

Visit us at bio-rad.com/NABT19

Call toll free at 1-800-424-6723.

Outside the U.S. contact your local sales office.

BIO-RAD