

**Measuring CO<sub>2</sub> Evolution in Multiple Custom Flasks?**

- Why would someone want to make measurements from flasks using the LI-8100A and multiplexer? (2:00)
- What changes need to be made in the LI-8150 in order to make these measurements? (3:03)
- How does one select the appropriate flask size for their application? (4:45)
- How would one re-plumb the LI-8150? (11:15)
- Measuring temperature in flasks? (17:52)
- What instruments are needed to make flask measurements? (22:00)