

**Monitoring Instrumentation for Carbon Capture and Storage Projects**

- Why surface monitor? (2:00)
- What instruments from LI-COR can be used to measure CO<sub>2</sub> at the surface? (3:15)
- How can the LI-8100A System help detect leaks? (6:00)
- How do soil CO<sub>2</sub> fluxes vary in time and space? (19:00)
- Where are LI-COR instruments being used for CCS? (21:40)