

Webinar Outline: An Introduction to Soil CO₂ Flux

- Why is CO₂ soil flux important? (1:40)
- Where does soil CO₂ come from? (3:59)
- What affects soil CO₂ production? (4:44)
- What affects soil CO₂ movement? (7:20)
- What are some challenges in taking soil CO₂ flux measurements? (11:00)
- What is a closed system approach? And what are the advantages and disadvantages of that approach? (12:45)
- How does the pressure vent on the LI-8100A decrease pressure fluctuations from the wind? (15:23)