

Advantages of Automated Soil CO₂ Flux Measurements

- What do we mean by automated soil CO₂ flux measurements? (2:59)
- Why is understanding the soil flux important in an ecosystem? (3:47)
- Why is measuring soil flux difficult? (5:53)
- Why do we need to get the right answer? (6:55)
- What are some measurement methods used for measuring soil CO₂ flux? (8:15)
 - Survey Measurements (8:20)
 - Long-term Measurements (9:25)
 - Multiplexed Long-term Measurements (9:55)
 - What can be learned from Multiplexed systems? (10:44)
- Summary of main points in the webinar (19:37)