

**Quality Assurance and Quality Control of Soil CO<sub>2</sub> Flux Datasets**

- What do QA and QC mean? (00:58)
- Five Elements of Quality Assurance(1:25)
- Five Elements of Quality Control (4:54)
  - Instrument diagnostics (8:49)
  - Flux variables (9:46)
  - Computed flux values (11:48)
- Quality Control Example (13:18)