

### **How Low Can You Go?**

- Why control chamber CO<sub>2</sub> at low concentrations?(00:01)
- How do C4 plants differ from C3 plants? (00:23)
- Controlling LI-6400/XT chamber concentrations (1:24)
- Needed time, supplies, and tools to modify the flow path (4:08)
- Techniques and Warnings to know before modifying the flow path(5:06)
- System Disassembly (7:41)
- Installing Modified Plumbing (9:24)
- Software Edits and Calibration (14:15)
- Making measurements with modified flow path (19:05)
  - Flow Control (19:43)
  - Water Control (20:40)
- Diffusion at Low CO<sub>2</sub> concentrations (22:09)
- Modification Summary (23:40)