

Preparing your LI-6400/XT for Field Season

- Checking the Power Supply (1:18)
- Other parts to check on the instrument (7:38)
- Scrub Tubes (10:50)
- Checking the filter in the CO₂ source (21:30)
- Checking parts on the sensor head (25:00)
 - Chamber Gaskets (25:25)
 - Leaf Temperature Thermocouple (29:13)
- Troubleshooting the match valve (38:40)
- AGC (automatic gain control) voltages(47:00)
- Cleaning the analyzers (51:20)