

Controlling Temperature and Humidity

- What environmental factors affect leaf gas exchange? (00:12)
- Temperature's affect on plants (00:53)
- Humidity's affect on plants (3:15)
- Controlling Relative Humidity (17:55)
- Controlling Temperature (21:58)
- What if inherent controls aren't enough? (26:00)
- LI-610 Portable Dew Point Generator (29:10)