



6400-05 Conifer Chamber

Product Bulletin

LI-6400 Portable Photosynthesis System

Description

- ❑ Teflon lined to minimize water sorption.
- ❑ Accommodates needles up to 3.5 cm long and stems up to 1 cm in diameter.
- ❑ A 5 cm length shoot can be enclosed in the chamber.
- ❑ Light measurement via external LI-COR quantum sensor.

The 6400-05 Conifer Chamber is used to measure short needle conifers in a chamber environment where temperature, humidity and CO₂ are controlled by the LI-6400.

The conifer chamber is constructed of acrylic for durability and lined with teflon to minimize water sorption. Shoots with needles up to 3.5 cm long can be accommodated in the 7.5 cm diameter chamber. The 6400-05 mounts directly to the sample cell of the LI-6400's CO₂ and H₂O IRGAs in order to minimize response time.



Leaf temperature is computed by the LI-6400 using energy balance based on the exit air temperature measured by a thermocouple suspended in the air exhaust port. Modifications to the LI-6400's software for operation with the conifer chamber are stored in memory for easy setup.

A bracket is provided on the exterior of the chamber to mount a LI-COR quantum sensor to measure photosynthetically active radiation (PAR).

Accessories:

- | | |
|----------|--|
| 6400-31 | Conifer Chamber Gaskets (5 sets, included) |
| 9901-013 | External Quantum Sensor (not included) |

Ordering Information

6400-05 Conifer Chamber

LI-COR

Biosciences

4421 Superior Street • P.O. Box 4425 • Lincoln, Nebraska 68504 USA
US & Canada: 800-447-3576 • Int'l.: 402-467-3576 • FAX: 402-467-2819
www.licor.com • Offices and Distributors Worldwide

PPS 232 08/01