

6400-07 Needle Chamber

Product Bulletin

LI-6400 Portable Photosynthesis System

Description

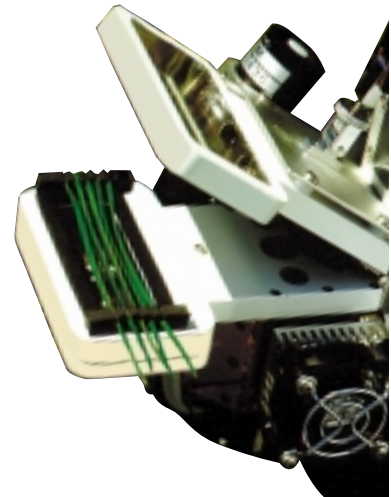
Photosynthesis rates of long-needled conifers are easily measured using the 6400-07 Needle Chamber. The 6400-07 has a 2 × 6 cm aperture and specially designed foam mounting pads that hold up to five needles in the chamber. The 6400-07 allows each needle to be pressed into place in the chamber, where it is held securely in the channels of the foam mounting pad. (The foam mounting pads can be ordered separately for measuring shorter needles with the standard 2 × 3 cm leaf chamber.)

The chamber top has a clear Propafilm® window for excellent light penetration. A PAR sensor in the chamber provides light measurement at the plane of the needles.

The chamber bottom is clear, allowing light to penetrate both the top and bottom of the chamber. The chamber bottom has a Propafilm® window similar to the chamber top. Leaf temperature is computed using energy balance by measuring with a small thermocouple the temperature of air exiting the chamber

Accessories: (included with 6400-07)

- | | |
|---------|--|
| 6400-32 | Needle and Narrow Leaf Chamber Gaskets (10 sets) |
| 6400-34 | Needle Holder Kit (5 sets) |



Ordering Information

- | | |
|---------|----------------|
| 6400-07 | Needle 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