

# 6400-11 Narrow Leaf Chamber

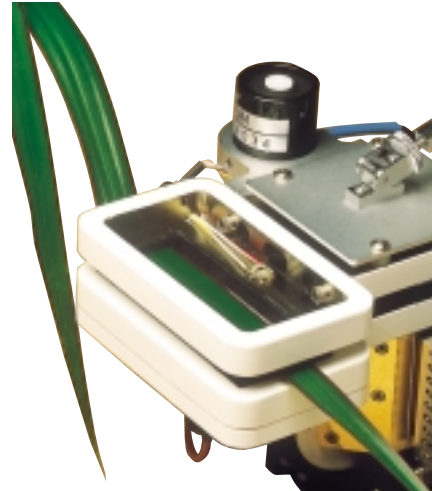
Product Bulletin

## LI-6400 Portable Photosynthesis System

### Description

The 6400-11 Narrow Leaf Chamber is ideal for foliage with narrow-bladed leaves. The 6400-11 has a 2 × 6 cm aperture like the 6400-07 Needle Chamber and uses the same chamber top. A PAR sensor for light measurement at the leaf plane is included.

The temperature of the foliage is measured with the standard LI-6400 leaf temperature thermocouple, which is inserted through the bottom of the lower half of the chamber.



### Accessories (included)

6400-32 Needle and Narrow Leaf Chamber Gaskets (10 sets)

### Ordering Information

6400-11 Narrow Leaf Chamber (2 × 6 cm)

**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](http://www.licor.com) • Offices and Distributors Worldwide