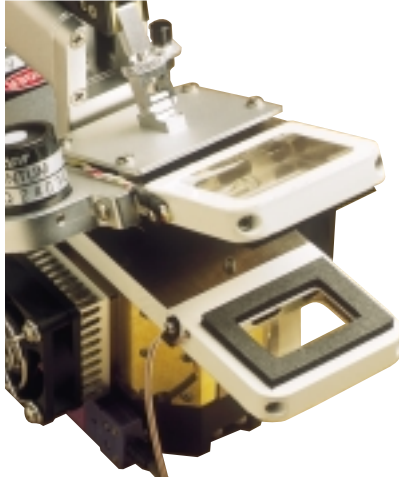


6400-08 Clear Chamber Bottom

Product Bulletin

LI-6400 Portable Photosynthesis System

Description



Illuminating the bottom of a leaf is particularly useful for survey and erect foliage measurements. The Clear Chamber Bottom has a Propafilm® window similar to the standard chamber top. Leaf temperature is computed using energy balance by measuring with a small thermocouple the temperature of air exiting the chamber.

The Clear Chamber Bottom is designed to be used with the 6400-06 PAM2000 Adapter, 6400-10 Mini-PAM Adapter or standard 2 × 3 cm chamber top.

Ordering Information

6400-08 Clear Bottom for Standard Leaf Chamber (2 x 3 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 • Offices and Distributors Worldwide