



LI-6400XT Training CD

LI-COR's strong commitment to outstanding technical support for the LI-6400XT Portable Photosynthesis System is expanding with the introduction of LI-6400XT Training CD.

WHAT IS IT?

The Training CD (part number 6400-512) is intended to serve as an introduction and overview of the operation of the LI-6400XT Portable Photosynthesis System. It contains information on both the hardware and software operation of the system. The CD contains more than three hours of instruction and information ranging from assembly of the system to daily maintenance, configuration for different experimental protocols, software overview, making measurements, and more.

WHO BENEFITS?

The first-time user of the LI-6400XT, new graduate and/or undergraduate students, technicians or researchers who want to learn the proper use and maintenance of the LI-6400XT, as well as anyone who wants a refresher course about the instrument.

CD CONTENT

The CD is organized chronologically into individual information modules, starting with what to expect when you first unpack the instrument, and concluding with how to efficiently manage the data collected.



A summary of the CD contents:

Introduction

- Initial Setup
- Installing Accessories

Software Introduction

- Open Screen
- Welcome Screen
- Shortcuts
- Utility Screen
- Configuration Screen

Getting Ready

- Calibration
- Daily Preparation Checklist

Making a Measurement

- Measurements
- New Measurements Screen
- Making Measurements Screen

Data Handling

- File Exchange Screen
- Recomputing data

<u>Part Number</u>	<u>Description</u>
6400-512	LI-6400XT Training CD
6400-513	LI-6400XT Training Set (5-CD Package)

For more information or to order the training CD, please contact LI-COR.



4647 Superior Street • P.O. Box 4425 • Lincoln, Nebraska 68504 USA
North America: 800-447-3576 • International: 402-467-3576 • Fax: 402-467-2819
envsales@licor.com • envsupport@licor.com • www.licor.com

In Germany and Norway - LI-COR GmbH: +49 (0) 6172 17 17 771
envsales-gmbh@licor.com • envsupport-gmbh@licor.com
In United Kingdom and Ireland - LI-COR UK Ltd: +44 (0) 1223 422102
envsales-UK@licor.com • envsupport-UK@licor.com