



6400-70 AC Module

Product Bulletin

Description

This optional accessory allows the LI-6400/6400XT to be powered from mains power. It consists of a transformer, and a battery-shaped box that can fit in the console (Figure 1).



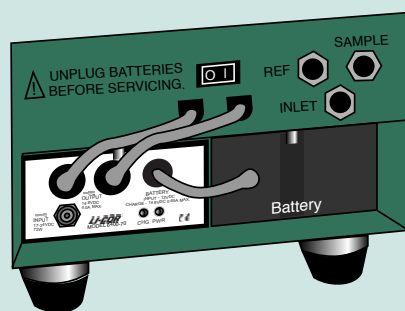
Figure 1. The 6400-70 AC Module fits in a battery compartment in the console. If you install a battery in the other compartment, and plug it into the module (as shown at left), that battery will trickle charge, but more importantly, provide continued operation for 1-2 hours in the event of a mains power failure.

Notes on using the 6400-70 AC Module

- **A battery is not required**
The module will power the LI-6400/6400XT by itself. However, plugging a battery into the module will provide 1 or 2 hours backup in case of mains power failure.
- **Use both plugs - most of the time**
The AC module has two plugs that go into the console's battery connectors. For normal operations, plug both of them in.
- **Hot swapping**
You can switch from AC operation to battery operation by unplugging one of the module's plugs from the console, and replacing it with a fresh battery. Then disconnect the other plug, and replace it with second battery if you wish. Reverse the process to switch from battery to AC operation. Either way, observe the caution in Figure 2 below.

Safe for extended operations

Battery safely trickle charges, and is available to run the instrument if mains power fails.



Don't do this for more than a minute

Battery is being charged in an uncontrolled manner. Damage could result to the battery and/or the transformer.

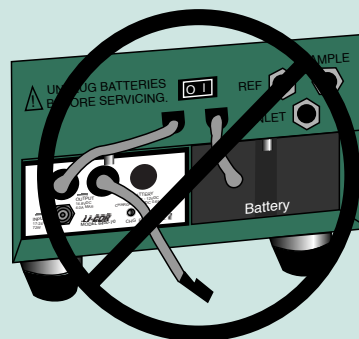


Figure 2. Avoid extended (more than a minute) operations with a battery and one of the AC module's plugs connected to the console. In this configuration, you will be providing uncontrolled charging of the battery, which could damage the battery and/or transformer.

LI-6400/6400XT Portable
Photosynthesis System

LI-COR®

Biosciences

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 Biosciences GmbH: +49 (0) 6172 17 17 771 • envsales-gmbh@licor.com • envsupport-gmbh@licor.com

In UK and Ireland – LI-COR Biosciences UK Ltd.: +44 (0) 1223 422102 • envsales-UK@licor.com • envsupport-UK@licor.com