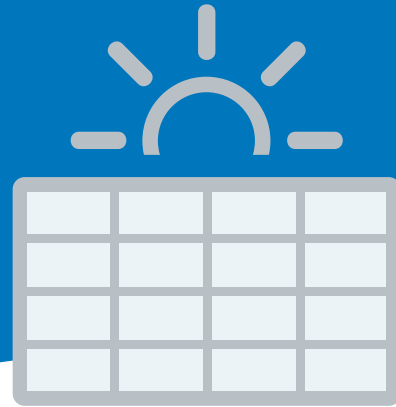


Solar Power

Safe and Reliable Power



LI-COR solar power systems are robust solutions developed to our specifications by SunWize Technologies for your eddy covariance systems. Providing up to 4 dark-days of operation, these 24-volt systems deliver power with greater efficiency than 12-volt systems. LI-COR systems are designed to ensure maximum uptime and long operation life by providing the appropriate voltage to all your instruments regardless of temperature or length of cable run.

Complete systems are available in 25, 40, 50, and 65-watt configurations for locations with 1.5 or more sun hours and are ready to deploy upon delivery. LI-COR can help determine which system is best suited for your instrument setup and location.

To learn more about the solar power options for your eddy covariance system visit licor.com/comms_power.

Key Features

- Designed to provide up to 4 dark-days of operation with LI-COR eddy covariance systems
- Robust and reliable power supply ensures maximum system uptime to avoid data loss
- Packages suitable for various eddy covariance system configurations and customized to your location
- Components provided include batteries, solar panels, and appropriately gauged wire to ensure the correct voltage reaches your instruments
- Charge controller comes preinstalled in enclosure
- Efficient 24-volt systems with available DC-to-DC converter for 24-to-12 VDC conversion

Solar Power Specifications

Typical System Includes:

- Solar Panel Array
- Lead-acid Batteries
- Vented Enclosure
- Charge Controller
- Solar Panel Power Cables
- Solar Panel Mounting Hardware
- Assembly Hardware
- Instruction Manual
- 7900-231 Power Distribution Box

25 W Solar Power Packages

Package 1: 1.5 – 1.9 sun hours

Package 2: 2.0 – 3.9 sun hours

Package 3: 4.0+ sun hours

40 W Solar Power Packages

Package 4: 1.5 – 1.9 sun hours

Package 5: 2.0 – 2.9 sun hours

Package 6: 3.0+ sun hours

50 W Solar Power Packages

Package 7: 1.5 – 2.9 sun hours

Package 8: 3.0 – 3.9 sun hours

Package 9: 4.0+ sun hours

65 W Solar Power Packages

Package 10: 2.0 – 2.9 sun hours

Package 11: 3.0 – 3.9 sun hours

Package 12: 4.0+ sun hours

Specifications subject to change without notice.