

LI-6800 Portable Photosynthesis System

The LI-6800 Portable Photosynthesis System measures gas exchange from leaves and other samples. When equipped with the fluorometer, it also measures chlorophyll *a* fluorescence. This document provides performance specifications for a typical LI-6800 and accessories.

Distinguishing specifications

CO₂ gas analyzers

CO₂ Analyzer Precision: Within 0.1 $\mu\text{mol mol}^{-1}$ RMS with 4-second averaging at 400 $\mu\text{mol mol}^{-1}$

H₂O gas analyzers

H₂O Analyzer Precision: Within 0.01 mmol mol^{-1} RMS with 4-second averaging at 10 mmol mol^{-1}

System

Bulk Flow Rate Range: 680 – 1700 $\mu\text{mol s}^{-1}$ at SATP¹

Light Source Uniformity: $<\pm 10\%$ variation over 90% of aperture

Chamber Temperature Control Range: 10 °C above or below ambient

Fluorometer

Modulation Frequency: 1 Hz – 250 kHz

Measuring Light Peak Wavelength: 625 nm

Red Actinic and Saturating Flash Peak

Wavelength: 625 nm

Blue Actinic Peak wavelength: 475 nm

Far-red Peak Wavelength: 735 nm

Actinic Light Output:

0 – 3000 $\mu\text{mol m}^{-2} \text{s}^{-1}$ total at 25 °C

0 – 1000 $\mu\text{mol m}^{-2} \text{s}^{-1}$ blue at 25 °C

0 – 2000 $\mu\text{mol m}^{-2} \text{s}^{-1}$ red at 25 °C

Saturation Light: 0 – 16,000 $\mu\text{mol m}^{-2} \text{s}^{-1}$ at 25 °C

Far-red Light: 0 – 20 $\mu\text{mol m}^{-2} \text{s}^{-1}$ at 25 °C

Full specifications

CO₂ gas analyzer

Type: Absolute non-dispersive infrared gas analyzer

Measurement range: 0 – 3100 $\mu\text{mol mol}^{-1}$

Precision (signal noise) at 400 $\mu\text{mol mol}^{-1}$

RMS with 4-second signal averaging: $\leq 0.1 \mu\text{mol mol}^{-1}$

Accuracy: Within 1% of reading at 200 $\mu\text{mol mol}^{-1}$ or above, $\pm 2 \mu\text{mol mol}^{-1}$ at $< 200 \mu\text{mol mol}^{-1}$

Orientation sensitivity: $\leq \pm 1 \mu\text{mol mol}^{-1}$ variation at 400 $\mu\text{mol mol}^{-1}$ from any orientation

H₂O gas analyzer

Type: Absolute non-dispersive infrared gas analyzer

Measurement range: 0 – 75 mmol mol^{-1}

Precision (signal noise) at 10 mmol mol^{-1}

RMS with 4-second signal averaging: $\leq 0.01 \text{mmol mol}^{-1}$

Accuracy: within 1.5% of reading at $> 5 \text{mmol mol}^{-1}$; $\pm 0.08 \text{mmol mol}^{-1}$ at $< 5 \text{mmol mol}^{-1}$

Temperatures

Operating temperature range: 0 – 50 °C

Storage temperature range: -20 °C – 60 °C

Temperature control range

Leaf temperature: $\pm 10 \text{ }^\circ\text{C}$ from ambient with 3 cm \times 3 cm chamber

Setpoint resolution: 0.1 °C

Chamber exhaust air temperature and temperature control block

Type: Thermistor

Range: -10 – 60 °C

Accuracy: ±0.15 °C

Leaf temperature sensor

Type: Type E fine-wire thermocouple

Sensitivity range: -10 – 60 °C

Accuracy: <±0.5 °C total; ±0.2 °C cold junction reference; ±0.3 °C thermocouple when within ±10 °C of cold junction temperature

Air flow rates

Bulk flow rate range: 680 – 1700 $\mu\text{mol s}^{-1}$ at SATP¹

Leaf chamber flow rate range: 0 – 1400 $\mu\text{mol s}^{-1}$ at SATP

Pressure

Console pressure sensor

Operating range: 50 – 110 kPa

Accuracy: ±0.4 kPa

Resolution: 1.5 Pa typical

Signal noise: ≤ 0.004 kPa peak-to-peak with 4-second signal averaging

Chamber pressure sensor

Range: -2 – 2 kPa

Resolution: < 1 Pa typical

Signal noise: 1 Pa peak-to-peak with 4-second signal averaging

Setpoint resolution: 1.0 Pa

Control range: 0 – 0.1 kPa (dependent on flow rate through the chamber)

Communication

RJ-45 Ethernet; IP/TCP for networks and computers: 1

Head connections: 2

Accessory connections: 2

Batteries

Weight: 0.435 kg

Capacity: 6800 mAh

Type: Lithium ion

Storage: -20 – 60 °C; ≤ 80% RH

CO₂ control

CO₂ control range: 0 – > 2000 $\mu\text{mol mol}^{-1}$ (with pump set to low; dependent on bulk flow rate)

Sources:

8 gram CO₂ cartridge

External tank with optional adapter kit

CO₂ cartridge lifetime: > 8 hours after puncture (dependent on setpoint)

CO₂ scrubber: Soda lime

H₂O control

H₂O control range: 0 – 90% RH (noncondensing)

Humidifier substrate: Pall Stuttgarter Masse ceramic substrate (Pall Corporation)

Desiccant: Sorbead[®] Orange CHAMELEON[®] silica gel beads

Light measurement

Chamber and light source PAR sensors:

Detector: Gallium arsenide

Sensitivity range: 0 – 3000 $\mu\text{mol m}^{-2} \text{s}^{-1}$

Resolution: < 1 $\mu\text{mol mol}^{-1}$

Calibration accuracy: ±5% of reading; traceable to the U.S. National Institute of Technology (NIST)

External LI-190R PAR Sensor:

Detector: Silicon photodiode

Sensitivity: 5 – 10 μA per 1000 $\mu\text{mol s}^{-1} \text{m}^{-2}$

Calibration accuracy: ±5% of reading; traceable to NIST

Console

Processor: Arm[®] Cortex^{®2} A9 Quad Core running at 1 GHz.

Memory: 2 GB RAM; 8 GB Flash memory

Display: Sunlight-readable TFT LCD with capacitive touch screen

Resolution: 1024 × 600 pixels

Dimensions: 26 cm diagonally

Size: 18.5 cm × 27.5 cm × 21 cm; (D × W × H)

Weight: 6.1 kg

Power requirements: 12 to 18 VDC or 24 VDC

Head

Size with 3x3 cm chamber: 37 cm × 11.5 cm × 21.6 cm (L × W × H)

Weight: 2.15 kg without chamber

Display resolution: 128 × 128 pixels

Display dimensions: 3.15 cm diagonally

Sensor head inputs:

Leaf temperature thermocouple: 2

LI-190R light sensor: 1

Sensor head light source connections: 1

Leaf Chambers, Light Sources, and Accessories

The following components may be ordered separately or as a package with the LI-6800.

Fluorometer chamber

Modulated light: Software controlled, selectable frequencies of 1 Hz – 250 kHz

Measuring light peak wavelength: 625 nm

Blue actinic peak wavelength: 475 nm

Red actinic and saturating flash peak wavelength: 625 nm

Far-red peak wavelength: 735 nm

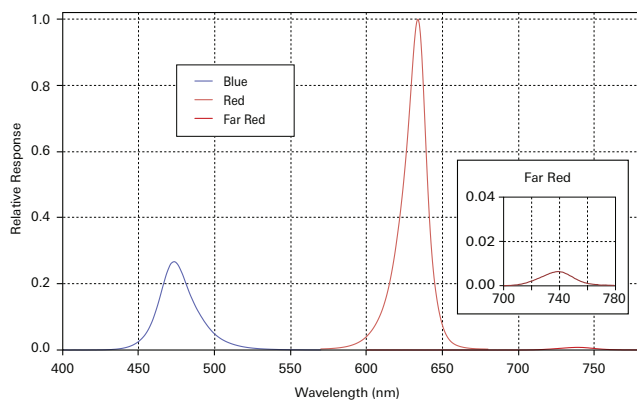


Figure 1. Typical spectral output of the blue, red, and far red LEDs in the fluorometer.

Actinic light output:

0 – 3000 $\mu\text{mol m}^{-2} \text{s}^{-1}$ total at 25 °C

0 – 1000 $\mu\text{mol m}^{-2} \text{s}^{-1}$ blue at 25 °C

0 – 2000 $\mu\text{mol m}^{-2} \text{s}^{-1}$ red at 25 °C

Saturation light: Software controlled intensity; 0 – 16,000 $\mu\text{mol m}^{-2} \text{s}^{-1}$ at 25 °C

Far-red light: Software controlled intensity; 0 – 20 $\mu\text{mol m}^{-2} \text{s}^{-1}$ at 25 °C

Fluorescence signal temperature dependence: - 0.25% per °C

Uniformity:

<±10% over 92% of the aperture with white top gasket

<±10% over 90% of the aperture with black top gasket

Power consumption:

< 18 W at 25 °C with 3000 $\mu\text{mol m}^{-2} \text{s}^{-1}$ actinic light

< 60 W at 25 °C with 16,000 $\mu\text{mol m}^{-2} \text{s}^{-1}$ saturating flash

Leaf area: 6 cm² or 2 cm²; Round apertures

Dimensions: 16.6 cm × 11.5 cm × 13.6 cm (L × W × H)

Weight: 0.86 kg

Small leaf chamber (3 × 3 cm)

Aperture areas: 9 cm² (3 cm × 3 cm), 6 cm² (2 cm × 3 cm), 3 cm² (1 cm × 3 cm)

Size: 15.4 cm × 11.5 cm × 5.9 cm (L × W × H)

Weight: 0.3 kg

Large leaf and needle chamber (6 × 6 cm)

Maximum leaf area: 36 cm²

External dimensions: 11.5 cm wide (spring to spring) × 16.8 cm long (front of chamber to back of chamber link) × 5.9 cm tall (bottom of chamber manifold to top of chamber)

Weight: 0.35 kg

Large leaf and needle chamber with conifer sprig blocks

Chamber Volume: 420.8 cm³ (internal)

Maximum planar leaf area: 36 cm²

Internal chamber height: 6.7 cm

External dimensions: 11.5 cm wide (spring to spring) × 16.8 cm long (front of chamber to back of chamber link) × 7.2 cm tall (bottom to top of chamber)

Small plant chamber

Chamber volume: 193.2 cm³ (internal)

Internal dimensions: 7 cm diameter; 4.46 cm deep

External dimensions: 8.4 cm wide (chamber block) × 12.7 cm long (chamber manifold interface to tip of chamber) × 6.47 cm (bottom of chamber manifold to top of chamber)

Weight: 0.60 kg

Bryophyte chamber

Chamber volume: 193.2 cm³ (internal)

Internal dimensions: 7 cm diameter; 4.45 cm deep

External dimensions: 8.4 cm wide (chamber block) × 12.7 cm long (chamber manifold interface to tip of chamber) × 6.47 cm (bottom of chamber manifold to top of chamber)

Weight: 0.60 kg

Insect respiration chamber

Chamber volume: 49.9 cm³ (not including tubing)

Exterior dimensions: 11.25 cm length; 3 cm diameter

Weight: 0.07 kg

Small light source (3 × 3 cm)

The small light source is compatible with the small leaf chamber.

Total output range: 0 – >2000 $\mu\text{mol m}^{-2} \text{s}^{-1}$ at 25 °C

Blue output range: 0 – > 400 $\mu\text{mol m}^{-2} \text{s}^{-1}$ at 25 °C

Red output range: 0 – > 1600 $\mu\text{mol m}^{-2} \text{s}^{-1}$ at 25 °C

Red peak wavelength: 660 nm

Blue peak wavelength: 453 nm

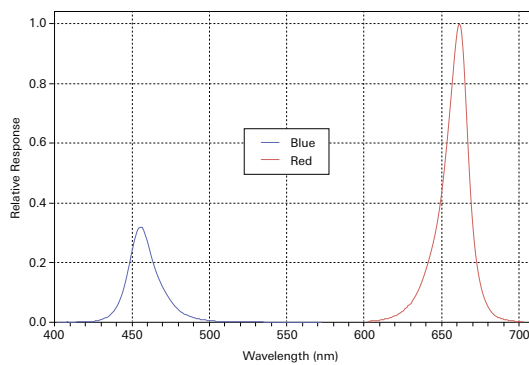


Figure 2. Typical spectral output of the LEDs in the 3x3 cm light source.

Uniformity:

±10% over 90% of the aperture with white top gasket, typically

±10% over 77% of the aperture with black gasket, typically

Power consumption at 2000 $\mu\text{mol m}^{-2} \text{s}^{-1}$: < 5 W

Operating temperature range: 0 – 50 °C

Size: 6.6 cm × 5.9 cm × 5.8 cm (L × W × H)

Weight: 0.21 kg

Large light source (6 × 6 cm)

The large light source is compatible with the large leaf chamber, small plant chamber, sprig adapter kit, and the bryophyte chamber.

Total output range: 0 – >2500 $\mu\text{mol m}^{-2} \text{s}^{-1}$ at 25 °C

Blue output range: >2000 $\mu\text{mol m}^{-2} \text{s}^{-1}$ at 25 °C

Green output range: >1000 $\mu\text{mol m}^{-2} \text{s}^{-1}$ at 25 °C

Red output range: >2400 $\mu\text{mol m}^{-2} \text{s}^{-1}$ at 25 °C

White output range: >1500 $\mu\text{mol m}^{-2} \text{s}^{-1}$ at 25 °C

Blue peak wavelength: 453 nm

Green peak wavelength: 523 nm

Red peak wavelength: 660 nm

White color temperature: 4000 K

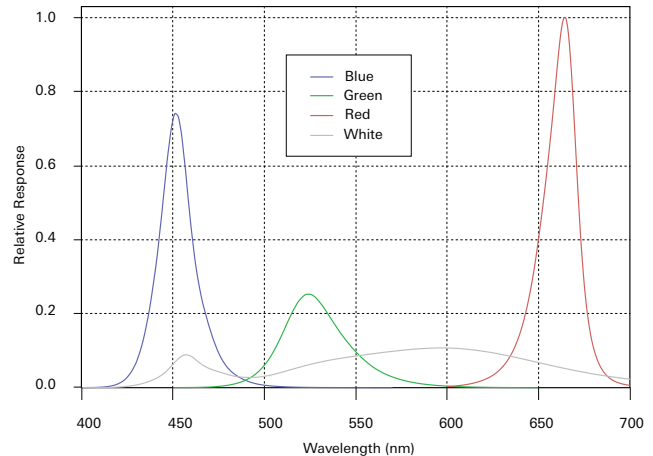


Figure 3. Typical spectral output of the blue, green, red, and white LEDs in the 6x6 cm large light source.

Uniformity: ±10% over 90% of the aperture

Power consumption: 15 W at 2000 $\mu\text{mol m}^{-2} \text{s}^{-1}$

equal parts red, green, blue, and white

Operating temperature range: 0 – 50 °C

Operating relative humidity range: 0 – 85%

Size: 11.7 cm × 11 cm × 13 cm (L × W × H)

Weight: 0.54 kg

Aquatic chamber

Sample volume: >0 to 20 mL; 15 mL recommended

Wetted materials: 316 stainless, float glass, Viton, PTFE, silicone, acetal

Operating temperature range: Non-freezing to 50 °C with no solar load

Sample temperature control: User supplied water bath

Storage temperature range: -20 to 60 °C with clean and dry chamber

Sample salinity range: 0 to 35 ‰

pH probe port

pH probe opening 12 mm diameter O-ring sealed port

Compatible probe type: Passive glass-electrode based pH probe

BNC Connector: Nominal -59mV/pH slope, integrated amplifier with user-settable calibration

Septum port

Material: Silicone-PTFE

Size: 3/8" diameter

Soil chamber

System

Chamber volume: 4244 cm³

IRGA volume: 57 cm³

Soil area: 317.8 cm²

Air temperature thermistor:

Operating range: -20 to 45 °C

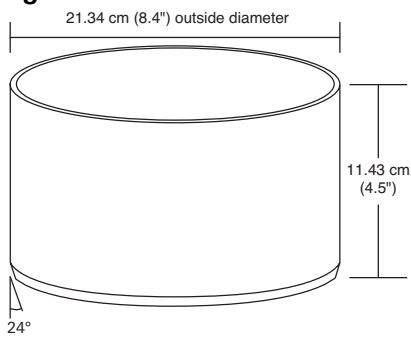
Accuracy: ±0.5 °C over 0 °C to 70 °C

Soil Collars

Inside diameter: ~20 cm

Outside diameter: ~21 cm

Height: 11.43 cm



¹ SATP is defined as Standard Ambient Temperature (25 °C) and Pressure (100 kPa).

² Arm and Cortex are registered trademarks of Arm Limited (or its subsidiaries) in the US and/or elsewhere.



LI-COR Biosciences

4647 Superior Street
Lincoln, Nebraska 68504
Phone: +1-402-467-3576
Toll free: 800-447-3576 (U.S. and
Canada)
envsales@licor.com

LI-COR Distributor Network:
www.licor.com/envdistributors

Regional Offices

LI-COR Biosciences GmbH
Siemensstraße 25A
61352 Bad Homburg
Germany
Phone: +49 (0) 6172 17 17 771
envsales-gmbh@licor.com

LI-COR Biosciences UK Ltd.
St. John's Innovation Centre
Cowley Road
Cambridge
CB4 0WS
United Kingdom
Phone: +44 (0) 1223 422102
envsales-UK@licor.com

Copyright © 2021 LI-COR, Inc.
All rights reserved.
982-16875 10/2021

Specifications subject to change. LI-COR is an ISO 9001 registered company. LI-COR is a registered trademark of LI-COR, Inc. in the United States and other countries. All other trademarks belong to their respective owners. For trademark information, visit https://www.licor.com/corp/copyrights_trademarks. For patent information, visit www.licor.com/patents.