

LI-7200RS

Enclosed CO₂/H₂O Gas Analyzer

Water vapor and carbon dioxide are the two most influential greenhouse gases affecting global climate change. The LI-7200RS is the next generation high-performance, enclosed CO₂/H₂O gas analyzer chosen by leading flux networks around the world including NEON and ICOS.



The LI-7200RS is designed to deliver the high-speed CO₂/H₂O measurements of a closed path analyzer with the ruggedness and low power demands of an open path analyzer. Temperature-controlled optics ensure stability – even in very harsh environments.

To learn more about the LI-7200RS visit [licor.com/LI-7200RS](https://www.licor.com/LI-7200RS).

Key Features

- Minimize data gaps with continuous measurements through rain, snow, and fog
- An insulated or heated intake tube ensures high frequency response of water vapor measurements, even in high humidity
- SmartFlux[®] System processes real-time fluxes on-site using EddyPro[®] Software
- SmartFlux syncs instrument clocks using GPS and PTP for precise data synchronization within and across sites
- Improved, temperature-controlled optics deliver stable measurements, even in extreme conditions
- Compatible with nearly all fast response, three-dimensional sonic anemometers to meet your unique research needs

LI-7200RS Specifications

CO₂ Measurements

Calibration range:

0 to 3000 $\mu\text{mol mol}^{-1}$

Accuracy: Within 1% of reading**Zero drift (per °C):**

- ± 0.1 ppm typical
- ± 0.3 ppm maximum

RMS noise (typical @ 370 ppm CO₂ and 10 mmol mol⁻¹ H₂O):

- @5 Hz: 0.08 ppm
- @10 Hz: 0.11 ppm
- @20 Hz: 0.16 ppm

Gain drift (% of reading per °C @ 370 ppm):

- $\pm 0.02\%$ typical
- $\pm 0.1\%$ maximum

Direct sensitivity to H₂O (mol CO₂/mol H₂O):

- $\pm 2.00\text{E-}05$ typical
- $\pm 4.00\text{E-}05$ maximum

H₂O Measurements

Calibration range: 0 to 60 mmol mol⁻¹**Accuracy:** Within 2% of reading**Zero drift (per °C):**

- ± 0.03 mmol mol⁻¹ typical
- ± 0.05 mmol mol⁻¹ maximum

RMS noise (typical @ 370 ppm CO₂ and 10 mmol mol⁻¹ H₂O):

- @5 Hz: 0.0034 mmol mol⁻¹
- @10 Hz: 0.0047 mmol mol⁻¹
- @20 Hz: 0.0067 mmol mol⁻¹

Gain drift (% of reading per °C @ 20 mmol mol⁻¹):

- $\pm 0.15\%$ typical
- $\pm 0.30\%$ maximum

Direct sensitivity to CO₂ (mol H₂O/mol CO₂):

- ± 0.02 typical
- ± 0.05 maximum

General

Analysis Type: Absolute, non-dispersive infrared spectroscopy

Data Storage: 16 GB removable industrial grade USB storage device

Data Communication: Ethernet, Synchronous Devices for Measurement (SDM; >50 Hz), RS-232 (115,200 baud; 20 records per second max), 6 DACs (0-5 V; 300 Hz)

Inputs: Four analog input channels (differential; bi-polar; ± 5 V; 300 Hz)

Operating Temperature Range: -25 to 50 °C (-40 to 50 °C verification test available on request)

Power Requirements: 10.5 to 30 volts DC

Power Consumption: 12 W nominal (up to 30 W during startup)

Detector: Thermoelectrically cooled lead selenide

Bandwidth: 5, 10, or 20 Hz, user-selectable

User Interface: Windows® based

Cable length:

5 meters (all cables); 5-meter head cable extension available

Analyzer Head

Size: 7.5 cm (3") diameter, 31 cm (12.2") length

Weight: 1.8 kg (3.95 lbs.)

LI-7550 Analyzer Interface Unit

Size: 35 cm × 30 cm × 15 cm (13.8" × 12" × 6")

Weight: 4.4 kg (10 lbs)

7200-101 Flow Module

Operating Temperature Range: -25 to 50 °C

Power Requirements: 10.5 to 30 VDC

Power Consumption: < 16 W nominal

Flow Rate: 15 LPM nominal

Size: 35 cm × 30 cm × 15 cm (13.8" × 12" × 6")

Weight: 6.15 kg (13.55 lbs)

Insulated Intake Tube

Length: Up to 1 meter

Inside Diameter: 5.33 mm

Outside Diameter: 6.35 mm

Heated Intake Tube

Length: 71.1 cm

Inside Diameter: 5.33 mm

Outside Diameter: 6.35 mm

Operating Temperature Range: -40 to 50 °C

Output Wattage: 0.1 to 6 W

Heat Density Ratio: (short tube to long tube) 2:1

Weight: 0.54 kg (1.2 lbs)

7550-101 Auxiliary Sensor Interface

Size: 11.5 cm × 6.5 cm × 4.2 cm (4.5" × 2.6" × 1.7")

Weight: 0.39 kg (0.85 lbs) including mounting bracket

Specifications subject to change without notice.