

Coming Soon

# LI-7825 CO<sub>2</sub> Isotope Analyzer

The LI-7825 CO<sub>2</sub> Isotope Analyzer measures the four most abundant CO<sub>2</sub> gas isotopologues in air and reports  $\delta^{13}\text{C}$ ,  $\delta^{17}\text{O}$ , and  $\delta^{18}\text{O}$  with high precision and accuracy.



Reveal the story of your CO<sub>2</sub> data with precise CO<sub>2</sub> isotope measurements.

## By measuring CO<sub>2</sub> isotopologues and calculating isotope ratios, you can:

- Identify the sources and sinks of atmospheric carbon
- Partition net ecosystem carbon exchange
- Gain insight into biological processes
- Evaluate carbon sequestration efforts

Backed by over thirty-five years of experience in gas analysis technology, the LI-7825 joins the LI-COR Trace Gas Analyzers trusted by researchers and networks around the world.

**LI-COR**<sup>®</sup>

# Dependable, accurate, portable, and affordable

The LI-7825 features the same rugged design, portability, accuracy, and precision as other LI-COR Trace Gas Analyzers at an affordable price. Suitable for both the bench and the field, it is easily added to an existing instrument rack. It can integrate into established systems and delivers dependable data to empower researchers to test hypotheses and produce knowledge.

## Key Specifications

### General

**Total weight:** 10.5 kg (including batteries)

**Battery life:** 8 hours typical with 2 batteries

### Preliminary Specifications

**Measurement Range:** 50 to 2,000 ppm

**Precision (1 $\sigma$ ):** <0.3 ‰ with 1 second averaging at 400 ppm CO<sub>2</sub>

**Drift:** <1 ‰ 24 hours, peak-to-peak 1 hr interval average

## Applications

- Atmospheric monitoring
- Urban emissions monitoring
- Mobile emissions monitoring
- Large area emissions monitoring
- Sensor networks
- Mud logging

**More applications information, including leaf level and soil, coming soon.**

Sign up to receive updates:  
[licor.com/CO2-isotopes](https://licor.com/CO2-isotopes)

