

Measuring CH₄ and CO₂ with New Generation High-Precision Low-Power Low-Maintenance Closed-Path Analyzers: Lab and Field Results

K. Minish^{1*}, I. Begashaw¹, M. Johnson¹, A. Komissarov¹, W. Miller¹, D. Trutna¹, R. Walbridge¹, J. Welles¹ and G. Burba^{1,2}

¹LI-COR Biosciences, Lincoln, Nebraska USA; ²Robert B. Daugherty Water for Food Institute and School of Natural Resources, University of Nebraska, Lincoln, Nebraska, USA

Introduction

Global CO₂ and CH₄ monitoring requires instruments that must meet the stringent requirements for:

- ❑ Accuracy, precision and stability
- ❑ Low power consumption, field portability, and limited maintenance

Two new field portable analyzers were designed to address these requirements:

- ❑ LI-7810 CH₄/CO₂/H₂O
- ❑ LI-7815 CO₂/H₂O

We report on the laboratory and field performance validation of both analyzers.



LI-7810 Stability and Precision

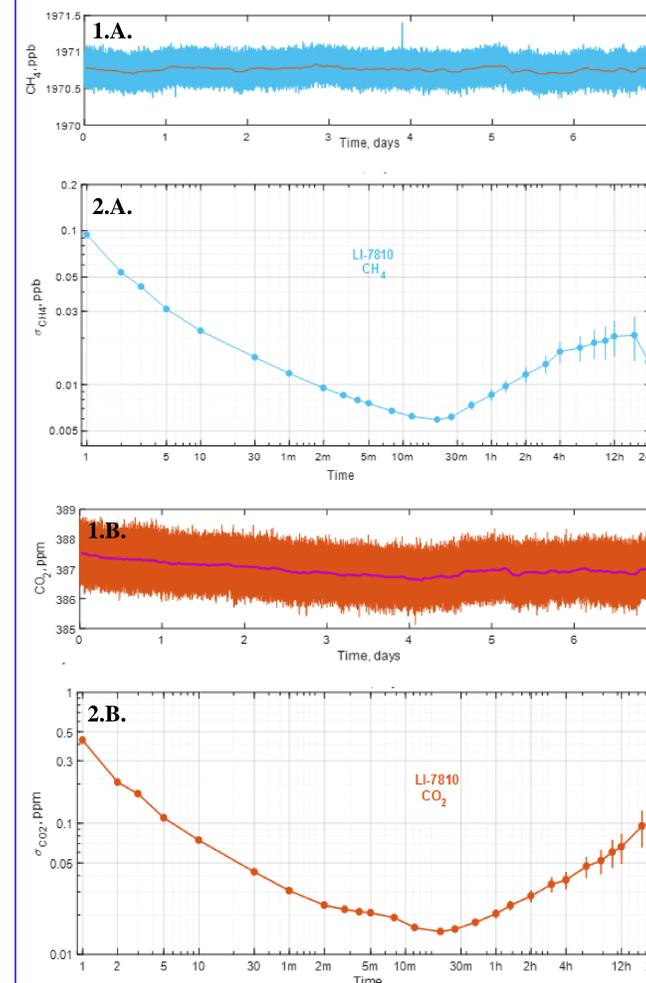


Chart 1: The instrument was operated continuously over a 7-day period with a continuous flow of tank gas. The blue line shows 1-second measurements; the orange line shows a 50 minute block average. Chart 1.A. reflects the high precision CH₄ data set, with chart 1.B. providing the CO₂ data from the same instrument. The LI-7810 provides high precision CH₄ measurements, with CO₂ data similar to other IRGAs available.

Chart 2: Allan deviation plots for the LI-7810 Analyzer. The plot shows precision at 1-second signal averaging, with improved precision as averaging time increases. Error bars represent 68% ($\pm 1\sigma$) confidence intervals based on number of averaged time series available at each τ . Chart 2.A. reflects the high precision CH₄ data set, with chart 2.B. providing the CO₂ data from the same instrument.

LI-7815 Stability and Precision

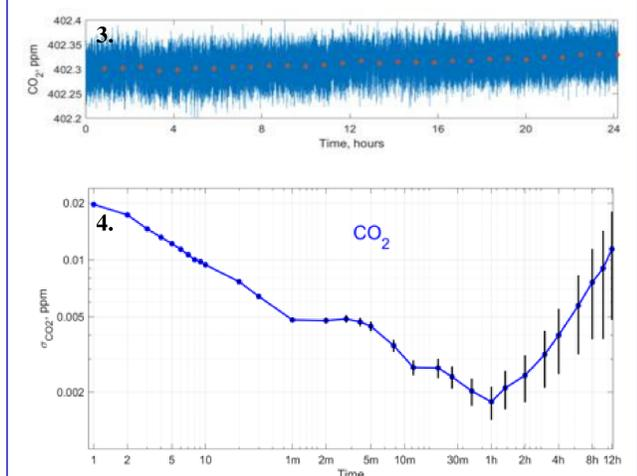
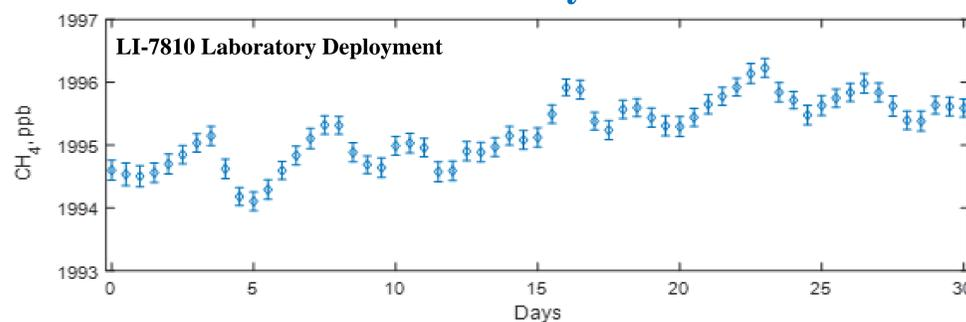


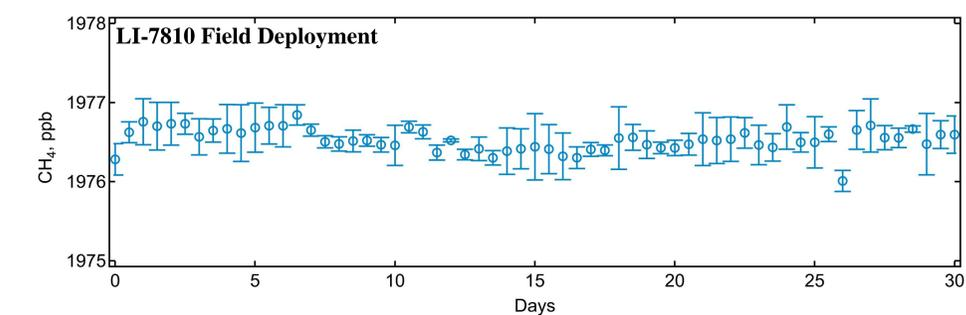
Chart 3: The instrument was operated continuously over a 24-hour period with a continuous flow of tank gas. The blue line shows 1-second measurements; the orange line shows a 50 minute block average.

Chart 4: Allan deviation plots for the LI-7815 Analyzer. The plot shows precision at 1-second signal averaging, with improved precision as averaging time increases. Error bars represent 68% ($\pm 1\sigma$) confidence intervals based on number of averaged time series available at each τ .

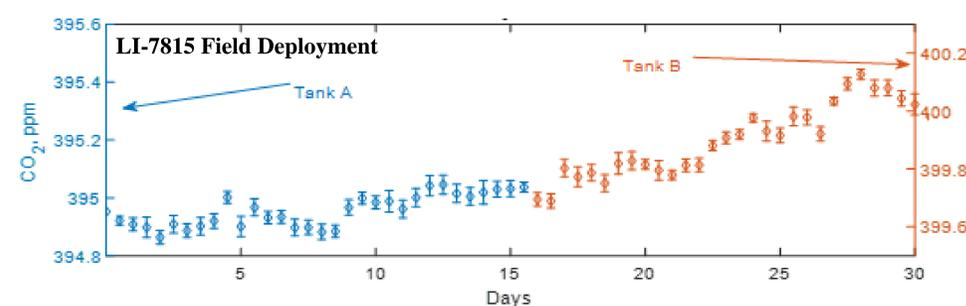
Laboratory and Field Validation



The analyzer measured ambient air between stability checks. During stability checks, the analyzer was flushed for 10 minutes with tank gas** and a time series data set was subsequently recorded for 10 minutes every 12 hours. The figure shows the mean value of the 10-minute measurements. Error bars represent a 1 σ confidence interval (68%).



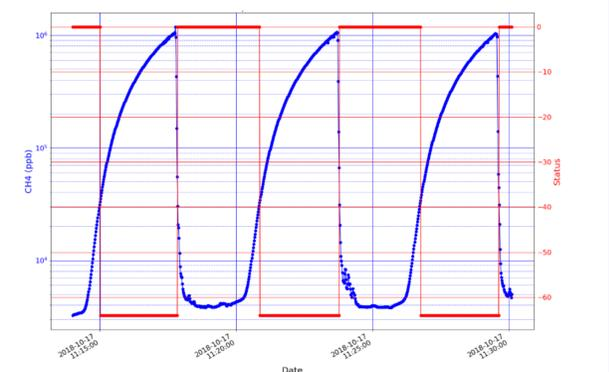
Located in an enclosed shelter, the analyzer measured ambient air between stability checks. During stability checks, the analyzer was flushed for 10 minutes with tank gas** and a time series data set was subsequently recorded for 10 minutes every 12 hours. The figure shows the mean value of the 10-minute measurements. Error bars represent a 1 σ confidence interval (68%).



Located in an enclosed shelter, the analyzer measured ambient air between stability checks. During stability checks, the analyzer was flushed for 10 minutes with tank gas** and a time series data set was subsequently recorded for 10 minutes every 12 hours. The figure shows the mean value of the 10-minute measurements. Error bars represent a 1 σ confidence interval (68%).

**Tank gas: 2 ppm CH₄, 400 ppm CO₂ mixed in air balance, 1% nominal concentration accuracy

Dynamic Range Tests



The LI-7810 was tested against higher concentrations of CH₄. The instrument responded and continued to measure, however, specifications are not guaranteed outside of the 0 – 50 ppm range for CH₄. This is reflected in the status codes given at higher concentration measurements.

Acknowledgements

The initial principles and some portions of the new technology presented herein were developed in part based on the grant from the MONITOR Program by the U.S. Department of Energy Advanced Research Projects Agency - Energy (ARPA-E), under award number DE-AR0000537.

*Contact information: kristen.minish@licor.com