## thermo scientific



# Monitoring of Plant CO<sub>2</sub> Exchange Within Terrestrial Ecosystems

#### Authors

H.J. Jost,<sup>1</sup> Eric Wapelhorst,<sup>1</sup> Magda Mandic,<sup>1</sup> Nils Stoebener,<sup>1</sup> Ulf Froehlich,<sup>1</sup> Laura Gangi,<sup>2</sup> Roland Bol,<sup>2</sup> and Nicolas Brüggemann<sup>2</sup>

<sup>1</sup>Thermo Fisher Scientific, Bremen, Germany

<sup>2</sup>Forschungszentrum Jülich GmbH, Institute of Bio- and Geosciences -Agrosphere (IBG-3), Jülich, Germany

#### Keywords

Delta Ray, CO $_{_2}$  Exchange, Terrestrial Ecosystem, Monitoring,  $\delta^{_{18}}\text{O},\,\delta^{_{13}}\text{C}$ 

#### Goal

- To observe rapid changes in plant metabolism by simultaneously monitoring carbon and oxygen isotope ratios of carbon dioxide under varying ambient conditions in a plant chamber.
- To accurately estimate the CO<sub>2</sub> fluxes in terrestrial ecosystems.

#### Introduction

The Thermo Scientific<sup>™</sup> Delta Ray<sup>™</sup> IRIS (Figure 1) consists of a laser-based Isotope Ratio Infrared Spectrometer (IRIS) and a Universal Reference Interface (URI) which can perform accurate and precise, continuous, in-situ monitoring of isotopologues of trace gases at ambient concentration.





Figure 1. The Thermo Scientific Delta Ray Isotope Ratio Infrared Spectrometer (IRIS)

The Delta Ray IRIS can be integrated into a plant chamber (Figure 2) experiment to detect rapid changes in the metabolism of a plant. The objective of this experiment is to measure the impact of different plants species on the oxygen isotopic signature of atmospheric  $CO_2$  under varying environmental conditions. This is important for an accurate estimation of  $CO_2$  fluxes in terrestrial ecosystems based on  $\delta^{18}O$  in  $CO_2$ .



Figure 2. Plant chamber used in the experiment

#### Method and experimental setup

 $CO_2$  is scrubbed from ambient air to create low  $CO_2$  air, and then  $CO_2$  is added back in to create the desired  $CO_2$ concentration (Figure 3). This air is then directed through the plant chamber. The air at the input of the chamber ("chamber in") and outlet ("chamber out") is alternatingly sampled by the Delta Ray IRIS. The Delta Ray IRIS valve A was switched every 5 minutes (300 s) (Figure 3).

The signal from valve A (see Figure 3) was fed to the Delta Ray analyzer and used as a trigger to start the acquisition of the "chamber in" and "chamber out" samples. Table 1 lists the details of the sample sequence. The sequence was started with a reference measurement from the Universal Referencing Interface of 120 seconds. A flush time of 60 seconds was also included between samples, and samples and reference measurements. Measurements were taken for 230 seconds of alternating "chamber in" and "chamber out". This left a buffer of 10 seconds for Delta Ray IRIS to wait for the next trigger from valve A sent after 300 seconds.



3/2 way solenoid valve



Table 1. Labbook setup for plant chamber measurements. This30 minute sequence is repeated for 1 day, and then a new labbookstarted.

Label	Flush time [s]	Measurement time [s]	Trigger
Ref 1	60	120	
Chamber out	60	230	0
Chamber in	60	230	1
Chamber out	60	230	0
Chamber in	60	230	1
Chamber out	60	230	0
Chamber in	60	50	1

#### **Results**

Nine days of continuous measurements starting after plant watering was stopped are shown in Figure 4. Changes in isotopic composition between light on and off are clearly visible. Uptake of  $CO_2$  by photosynthesis imposes an isotopic signature on the remaining  $CO_2$  by preferentially removing the main isotopologue ( $^{12}C^{16}O^{16}O$ ). The change is >3‰ and 1.5‰ for  $\delta^{18}O$  and for  $\delta^{13}C$ , respectively. The input isotopic composition is also changing over time, due to incomplete scrubbing of  $CO_2$ from ambient air. The scientific interpretation of similar data combined with additional measurements can be found in Gangi 2015.<sup>1</sup>

#### Conclusion

The Delta Ray Isotope Ratio Infrared Spectrometer provides continuous, feature-rich data for plant research. By providing high precision, and automatically fully referenced data the Delta Ray IRIS proved to be a valuable solution for the analysis of plant metabolism. The carbon isotope ratio  $\delta^{13}$ C-CO<sub>2</sub> and oxygen isotope ratio of atmospheric carbon dioxide ( $\delta^{18}$ O-CO<sub>2</sub>) can be used to partition the gross fluxes of CO<sub>2</sub> in terrestrial ecosystems, such as plant respiration, soil respiration and plant assimilation. The characteristic  $\delta^{13}$ C value is modified by plant metabolism and photosynthesis. The  $\delta^{18}$ O is affected by the oxygen exchange between the molecules of CO<sub>2</sub> and H<sub>2</sub>O stemming from different water pools. Similar measurements could for example be used to determine the efficiency of plants in phenotyping, studying the impact of elevated CO<sub>2</sub> concentrations in a future climate or even exchanges at ecosystem level.

#### Reference

 Gangi L, Real-time quantification of oxygen isotope exchange between carbon dioxide and leaf/soil water in terrestrial ecosystems with laser-based spectroscopy. Dissertation 2015, University of Bonn, Germany L. Gangi et al., *Agricultural and Forest Meteorology 201* (2015) 128–140.



Figure 4. a) Time series of  $CO_2$  concentration of Delta Ray chamber in and chamber out; b) time series of  $\delta^{18}O$  of the Delta Ray IRIS of plant chamber in and out; c)  $\delta^{13}C$  of the Delta Ray IRIS of plant chamber in and out

### Find out more at thermofisher.com/DeltaRay

©2018 Thermo Fisher Scientific Inc. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representatives for details. **AN30291-EN 0518S** 

