thermo scientific

APPLICATION BRIEF 30548

New analytical approach in monitoring of CO₂ cycle in aquatic ecosystems

Authors

Danijela Smajgl,¹ Florian Böhm,² Anton Eisenhauer,² Isabelle Taubner,² and Magda Mandic¹ ¹Thermo Fisher Scientific GmbH, Bremen, Germany ²GEOMAR Helmholtz Centre for Ocean Research Kiel, Kiel, Germany

Keywords

isotope ratio infrared analysis, carbon (δ^{13} C), oxygen (δ^{18} O), marine environment, CO₂ fluxes

Goal

To understand pathways, processes and the fate of CO_2 molecules in particular in the marine environment.

Introduction

The isotopic signatures of carbon (δ^{13} C) and oxygen (δ^{18} O) can be used as a tool to understand pathways, processes and the fate of CO₂ molecules, particularly in the marine environment (Figure 1). The reason for this can be found in the fact that δ^{13} C-values are controlled by species specific metabolic processes of respiration and photosynthesis, while the δ^{18} O-values are affected by the oxygen exchange between the molecules of CO₂ and the ambient water. Here we present a new analytical approach to determine δ^{13} C and δ^{18} O changes using a mid-infrared laser (IRIS) absorption spectrometer, Thermo Scientific[™] Delta Ray[™] IRIS with (Universal Reference Interface) URI[™] Connect.



Figure 1. Simplified view of CO₂ air-ocean flux.

Thermo Fisher

With IRIS technology, it is possible to record online changes of carbon and oxygen isotopes with time resolutions of seconds, and with this to have a better insight in CO_2 fluxes.

Delta Ray IRIS with URI Connect

The instrument is based on direct absorption spectroscopy and it uses mid-infrared laser that operates at 4.3 µm. Measurement of different carbon dioxide isotopologues and determination of stable isotope ratios from spectrum is possible due to absorption lines which are shifted relative to each other (Figures 2 and 3).



Figure 3. Thermo Scientific[™] Delta Ray[™] IRIS with URI[™] Connect and Teledyne CETAC[™] ASX-7100 Autosampler.



Figure 2. Mid-infrared spectral region at a wavelength of 4.3 μ m. The laser scans over the absorption lines at 500 Hz. The stable isotope ratios of carbon dioxide can be calculated from the respective peak areas of the different isotopologues.

Coral culturing facilities

This new approach was tested in a 600 I reef coral culturing tank at GEOMAR, Kiel.

The setup comprised three tanks with a closed recirculation system and with controlled day-night cycle (Figures 4 and 5). All water parameters are monitored to

keep the system in a stable state. Decrease of calcium concentration and carbonate alkalinity was compensated by calcium reactor, addition of bicarbonate and $CaCl_2$ solutions. Due to photosynthesis and respiration, the seawater pH ranges from 8.0 during the day and to 7.7 in the night, while temperature is kept constant at 26 ± 1°C.



Figure 4. The scheme shows day-night cycle control.



Figure 5. Coral culturing setup: 1) Main reef tank, 2) Culture tank, 3) Cleaning unit with three compartments; the first contains live rock, the second holds a protein skimmer, the third contains return pumps and sensors.

Measurements in controlled environment

IRIS technology enables us to conduct continuous measurements, with time resolution of seconds, and it also gives us opportunity to measure discrete samples.

Experimental setup: Air measurements

Equilibrator like setup was used for air measurements (Figure 6). Air was continuously pumped through the Delta Ray Analyzer with a flow of 80 mL/min. Measurement per sample was 3 minutes and with integration time of 60 seconds we achieved internal precision better than 0.05% for both δ^{13} C and δ^{18} O, per sample run.

Experimental setup: DIC measurements of discrete seawater samples

Along with continuous measurement of air, we collected water samples from the main reef tank every 1 to 3 hours during a period of 24 hours. The δ^{13} C-values of dissolved inorganic carbon (DIC) were analyzed with Delta Ray with URI Connect, using Teledyne CETACTM ASX-7100 Autosampler, following the gas evolution method with orthophosphoric acid (McCrea, 1950).





Air measurements: CO₂ (ppm), δ^{13} C and δ^{18} O

- Results present the daily average of approximately 100 days of continuous measurement.
- Due to biological production, isotopic change occur rapidly when light intensity is changed. Photosynthesis preferentially removes the light isotope ¹²C from seawater.
- Amplitude of change in CO_2 concentration during day-night cycle was approximately 740 ppm, while the average day-night amplitudes for $\delta^{13}C$ and $\delta^{18}O$ were about 2.7‰ and 0.6‰ (Figure 7, 8).







Figure 8. CO_ sources can be distinguished due to the $\delta^{\rm 13}\text{C}$ changes within a diurnal cycle.

δ^{13} C and δ^{18} O – Air and DIC measurements

- The δ¹³C-values of dissolved inorganic carbon (DIC) follow the same trend as it was observed during continuous air measurements, with an average amplitudinal change in δ¹³C of about 2.4‰. DIC in water becomes enriched in δ¹³C due to preferential uptake of ¹²C during photosynthesis.
- Measured data also showed that fractionation of CO_2 (g), relative to DIC (dissolved HCO_3^{-1}) is -7.5 ± 0.2‰, which is in a close agreement with isotope fractionation of -7.92‰ in the equilibrium system $CO_2^{-}HCO_3^{-1}$ at 26 °C (Mook et al. 1974) (Figure 9).



Figure 9. Comparison of continuous air measurements and measurements of DIC samples.

Conclusion and future work

- Delta Ray IRIS with URI Connect offers a possibility for in-situ monitoring of $\delta^{13}C$ and $\delta^{18}O$ changes in aquatic environments
- When combined with Equilibrator, Delta Ray IRIS with URI Connect offers possibility for measurements of CO₂ fluxes across air-seawater interface, and direct insight in $\delta^{13}C_{\text{DIC}}$ changes in the aquatic system
- Field experiments of this technique will show if a largescale measurement of CO₂ fluxes is possible, and whether this system may be used to monitor status of coral reef metabolism to estimate the influence of human activity in coastal areas

This project has received founding from the European Union's Horizon 2020 research and innovation programme under Marie Sklodowska-Curie grant agreement No 643084.







References

- McCrea, J. M. 1950: On the Isotopic Chemistry of Carbonates and a Paleotemperature Scale, J. Chem. Phys., vol. 18, no. 6, page 849-857 (June, 1950).
- Mook, W. G., Boommerson, J. C. and Staverman, W. H. 1974: Carbon isotope fractionation between dissolved bicarbonates and gaseous carbon dioxide. Earth and Planetary Science Letters, 22 (1974) 169-176.

Find out more at thermofisher.com/DeltaRay

©2018 Thermo Fisher Scientific Inc. All rights reserved. CETAC is a registered trademark of Teledyne Instruments Inc. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. 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. **AB30548-EN 1118S**

