

irm-GC/MS: $^{15}\text{N}/^{14}\text{N}$ Isotope Ratio Monitoring Chlorophyll Derivatives

Andreas Hilkert, Thermo Fisher Scientific, Bremen, Germany
First published as Finnigan MAT Application Flash Report No. G13 (1995)

Introduction

Tetrapyrrole ring structures (porphyrines) are the backbone of important biomolecules like haemoglobin, chlorophyll, and vitamin B₁₂. Because of the inherent stability of porphyrines, they can hold information on biological and environmental processes expressed in their isotope ratios for long periods of time.

Irm-GC/MS allows to read the ^{13}C and ^{15}N information in derivatized porphyrines in the low nmol to pmol range either for natural abundances or by conducting tracer experiments.¹ In irm-GC/MS, all compounds eluting from a GC column are converted on-line into CO₂ and/or N₂ in the combustion interface and transferred on-line into an Isotope Ratio Mass Spectrometry (IRMS). ^{15}N determination of compounds with high molecular weight (e.g. M* = 822 g/mol for the chlorophyll derivative described in this report) is a rigorous test of all components of an irm-GC/MS system. The GC/C interface and IRMS meet the requirements for highest sensitivity, linearity and stability while maintaining chromatographic integrity.

The ^{15}N determination of porphyrines is made difficult by the relatively low nitrogen content of the compounds and their derivatives. In the current case, the chlorophyll derivative contains 6.8% N, corresponding to 3.4% N₂. Ion statistics and ionization efficiency then ask for 43 times more sample for ^{15}N determination than for ^{13}C measurement with the same precision. In addition, the GC separation of such relatively high molecular weight compounds requires high temperature and high carrier gas flow.

This, together with the limited injection capacity of the GC column, reduces the amount of N₂ transferred to the IRMS and increases the risk of nitrogen background contribution formed by micro leaks in the GC connections.

Experimental

IRMS	Finnigan MAT 252
GC	Hewlett Packard 5890 series II
Injector	On column
Column	HT 5, 12 m, 0.32 mm i.d., film thickness 0.10 μm
GC program	1 min at 40°C; 35°C/min to 80°C/min; 20°C/min to 275°C/min; 6°C/min to 315°C/min; 10 min at 315°C; constant flow
GC/C interface	Standard GC/C II interface

Results

The chlorophyll derivative was prepared in duplicate (A₁, A₂) to demonstrate the external reproducibility of the entire ^{15}N determination procedure.¹ The injected amounts ranged from 750 pmol to 1.3 nmol corresponding to 120 - 220 pmol N₂ into the IRMS. Although the N₂ background is extremely low and stable, it forms a permanent (rectangular) contribution of about 30% to the integrated porphyrine peak (see Figure 1). Optimized and fully automated background correction algorithms during acquisition

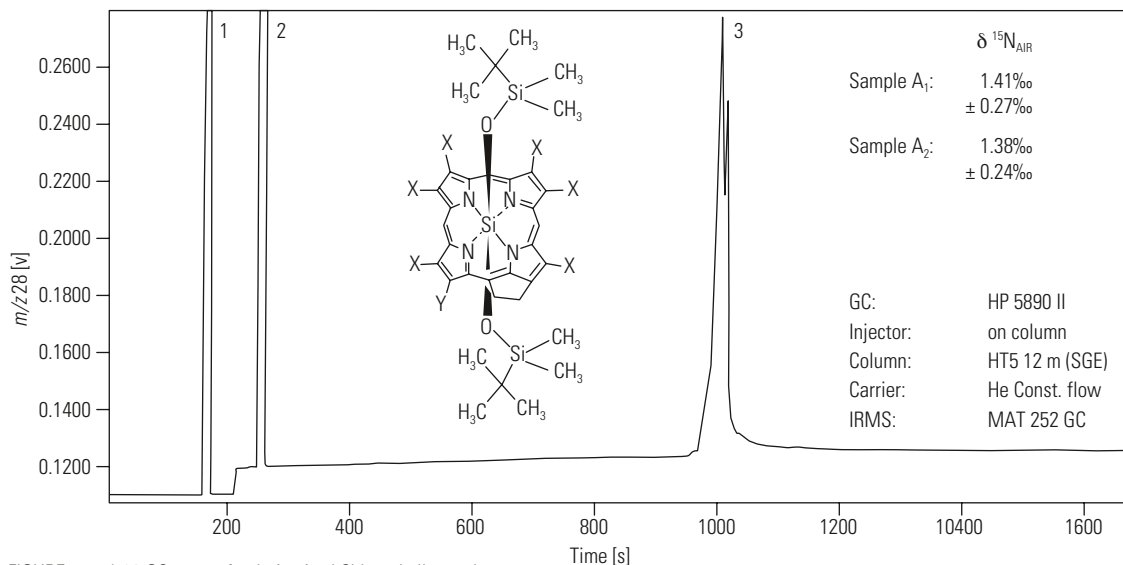


FIGURE 1. m/z 28 GC trace of a derivatized Chlorophyll sample.

Key Words

- Biomarkers
- GC Combustion
- Isotope Ratio MS
- Polyphyrines

tion and in post-acquisition data review of the ISODAT data system ensure a correct isotope ratio determination. The $\delta^{15}\text{N}$ value versus atm. air (international standard) of sample A¹ was $1.41\text{‰} \pm 0.27\text{‰}$, of sample A₂ $1.38\text{‰} \pm 0.24\text{‰}$. The corresponding standard deviation (n = 6) in atom percent was ± 0.00009 at %. The absolute values were within analytical error of the off-line determination. The results prove the high sensitivity in combination with linearity and stability of the MAT 252 IRMS coupled to the GC/C combustion interface.

Similar results can be expected from a DELTA V IRMS or a MAT 253 IRMS coupled with the GC Isolink unit via the ConFlo IV interface.

References

- ¹ J.P. Sachs, D.J. Repeta, A. Hilkert, Measurement and Assessment of the Carbon and Nitrogen Isotopic Composition of Chlorophyll Derivatives, presented at the *17th International Meeting on Organic Geochemistry*, San Sebastian, Spain, 1995

In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.

Africa-Other

+27 11 570 1840

Australia

+61 3 9757 4300

Austria

+43 1 333 50 34 0

Belgium

+32 53 73 42 41

Canada

+1 800 530 8447

China

+86 10 8419 3588

Denmark

+45 70 23 62 60

Europe-Other

+43 1 333 50 34 0

Finland/Norway/ Sweden

+46 8 556 468 00

France

+33 1 60 92 48 00

Germany

+49 6103 408 1014

India

+91 22 6742 9434

Italy

+39 02 950 591

Japan

+81 45 453 9100

Latin America

+1 561 688 8700

Middle East

+43 1 333 50 34 0

Netherlands

+31 76 579 55 55

New Zealand

+64 9 980 6700

Russia/CIS

+43 1 333 50 34 0

South Africa

+27 11 570 1840

Spain

+34 914 845 965

Switzerland

+41 61 716 77 00

UK

+44 1442 233555

USA

+1 800 532 4752



www.thermoscientific.com/irms

© 2012 Thermo Fisher Scientific Inc. All rights reserved. Comparable Hewlett Packard hardware is now sold under the Agilent brand. Agilent is a registered trademark of Agilent. HT5 is a trademark of SGE Analytical Science Pty Ltd.. ISO is a trademark of the International Standards Organisation. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

Thermo Fisher Scientific (Bremen)
GmbH Management System
Registered to ISO 9001:2008

AN30087_E 02/12G