

Nitrogen and Carbon Determination of Soils and Plants by the Thermo Scientific FLASH 2000 Analyzer Using a Single Reactor

Dominique Chevalier¹, Francesco Leone², Liliana Krotz,² and Guido Giazzi²

¹Thermo Fisher Scientific, Courtaboeuf, France, ²Thermo Fisher Scientific, Milan, Italy

Overview

Purpose: Demonstrate the nitrogen and carbon determination by Organic Elemental Analysis (OEA) using a single combustion/reduction reactor.

Methods: Soils and plants reference materials were analyzed through an elemental analyzer with automated autosampler using a single reactor.

Results: Data collected of nitrogen and carbon from different soils and plants reference material are discussed to assess the performance of the OEA analyzer using a single reactor without matrix effect.

Introduction

Nitrogen and carbon determination by combustion analysis is very commonly used for soils, plants, leaves, sediments, deposit on filters, animal tissues analysis and marine sciences analysis because these elements provides important information for their characterization in agricultural and environmental research.

The importance of soil and plant testing has increased in the last years, as many of the traditional methods are no longer suitable for routine analysis, due to their time consuming preparation and the required use of environmentally hazardous reagents. For this reason the need for an efficient analytical technique has become critical. As the demand for improved sample throughput, reduction of operational costs and minimization of human errors has increased dramatically, a simple and automated technique which allows fast analysis with an excellent reproducibility is the key for efficient nitrogen and carbon determination.

The Thermo Scientific™ FLASH™ 2000 Analyzer (Figure 1), operating with the dynamic flash combustion of the sample, meets the laboratory requirements, such as accuracy, reproducibility and high sample throughput. The standard configuration is based on a double reactors system: first reactor for combustion and catalytic oxidation of the combustion gases, the second is used to reduce nitrous oxides as N₂. The superior performance of the Thermo Scientific FLASH 2000 Analyzer allows reducing the amount of oxidation catalyst needed for NC analysis using a single reaction tube. The reactor filled with less amount of oxidation catalyst and copper ensures the complete conversion of gases produced by the combustion and offers advantages as large numbers of analysis before the maintenance, shorter analytical circuit allowing faster analysis, and the possibility to install two analytical circuits which are used alternatively (for example NC in the left furnace and CHNS/NCS in the right furnace, or NC in the left furnace and Sulfur by FPD detector in the right furnace) in a single analyzer. Using two autosamplers, the switching time from one analytical circuit to the second is only of few seconds, without the need for tools or mechanical modification.



FIGURE 1. Thermo Scientific FLASH 2000 NC Analyzer.

Methods

The FLASH 2000 Analyzer operates according to the dynamic flash combustion method. Samples are weighed in tin containers and introduced into the combustion reactor via the Thermo Scientific™ MAS 200R Autosampler together with oxygen. For NC *single reactor configuration*, after combustion, the resultant gases are carried by a helium flow to oxidation catalyst and to reduced copper, and finally pass through a halogen and sulfur absorber located inside the reaction tube.

After the reaction tube, water is trapped; a GC column separate gases and finally gases are detected by a sensitive and stable thermal conductivity detector (TCD) (Figure 2). A complete report is automatically generated by the dedicated Thermo Scientific™ Eager Xperience Data Handling Software and displayed at the end of the analysis.

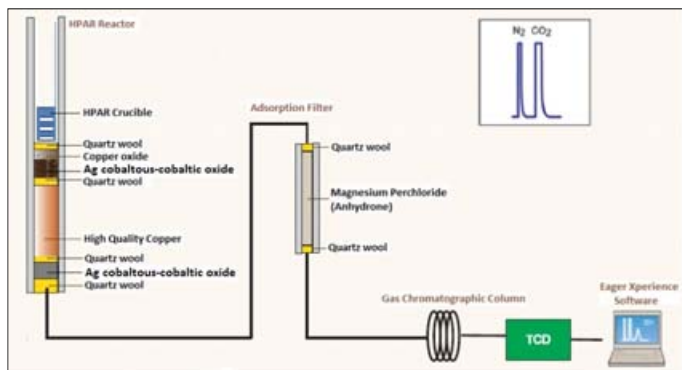


FIGURE 2. FLASH 2000 NC configuration.

To allow a large range of amount of nitrogen and carbon detected with a high quality of separation of these two elements, a three-meter GC column installed in the oven is used. Figure 3 shows a typical NC chromatogram.

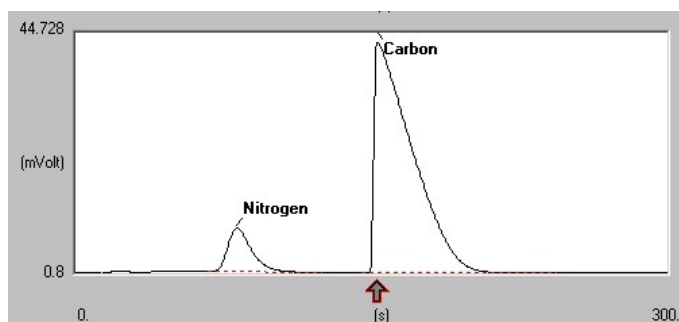


FIGURE 3. Typical NC chromatogram.

Results

Soils and plants reference materials with different nitrogen and carbon concentrations were analyzed in order to evaluate the performance of the single reactor. Table 1 shows the certified N % and C % and the relative uncertainty.

Reference Materials Description	Specification			
	N %	Uncertainty (±)	C %	Uncertainty (±)
Thermo Scientific Soil	0.21	0.01	2.29	0.07
Low Organic Content Soil	0.133	0.023	1.61	0.09
Medium Organic Content Soil	0.27	0.02	3.19	0.07
Loamy Soil	0.27	0.02	2.75	0.12
Chalky Soil	0.35	0.02	5.39	0.09
Sandy Soil	0.07	0.01	0.83	0.05
Birch Leaves	2.12	0.06	48.09	0.51
Orchard Leaves	2.28	0.04	50.40	0.40
Alfalfa	3.01	0.20	na	-

TABLE 1. Certified Nitrogen and Carbon of Reference Materials.

Table 2 shows the calibration sequence for soil analyses while Table 3 shows the weight of Soils Reference Materials analyzed and the experimental results obtained using Quadratic Fit or Linear Fit as calibration method. All results are obtained with a good repeatability and the averages values are inside the specification of the reference materials certificates.

Run	Standard	Type	Weight	Theoretical Value	
			mg	N %	C %
1	Benzoic acid	STD	5.974	-	68.85
2	Benzoic acid	STD	6.455	-	68.85
3	Aspartic acid	STD	6.548	10.52	36.09
4	Aspartic acid	STD	5.799	10.52	36.09
5	Aspartic acid	STD	2.528	10.52	36.09
6	Aspartic acid	STD	2.092	10.52	36.09
7	TF Soil ref. mat	STD	6.288	0.21	2.29

TABLE 2. Calibration sequence for soils analysis.

Sample Information	Quadratic Fit Calibration Method					Linear Fit Calibration Method				
	Reference Material	W (mg)	N%	RSD%	C%	RSD %	N%	RSD %	C%	RSD%
Thermo Scientific Soil	54.879	0.21			2.26		0.20		2.27	
	62.212	0.20	2.84		2.23	0.68	0.20		2.25	
	59.210	0.20			2.25		0.20	0	2.27	0.51
Low Organic Content Soil	52.310	0.12			1.57		0.12		1.58	
	59.495	0.12	4.68		1.56	0.97	0.12	4.68	1.56	0.97
	56.449	0.13			1.59		0.13		1.59	
Medium Organic Content Soil	57.837	0.27			3.15		0.27		3.17	
	66.428	0.27	0		3.15	0.18	0.27	2.17	3.16	0.32
	70.375	0.27			3.16		0.26		3.18	
Loamy Soil	66.068	0.26			2.67		0.25		2.68	
	56.356	0.26	0		2.69	0.37	0.25	2.28	2.71	0.57
	65.322	0.26			2.68		0.26		2.70	
Chalky Soil	53.043	0.37			5.32		0.36		5.34	
	59.841	0.37	0		5.34	0.47	0.36	1.59	5.35	0.28
	67.876	0.37			5.37		0.37		5.37	
Sandy Soil	240.058	0.065			0.813		0.064		0.818	
	255.114	0.066	0.88		0.817	0.32	0.065	0.89	0.821	0.31
	229.087	0.066			0.818		0.065		0.823	

TABLE 3. Experimental Nitrogen and Carbon data of Soils Reference Materials.

Table 4 shows the calibration sequence for plants analyses while Table 5 shows the weight of Plant Reference Materials and the experimental results obtained using Quadratic Fit or Linear Fit as calibration method. All results are obtained with a good repeatability and the values of the reference materials are inside the specification.

Run	Standard	Type	Weight	Theoretical Values	
			mg	N %	C %
1	Benzoic acid	STD	6.442	-	68.85
2	Benzoic acid	STD	6.749	-	68.85
3	Aspartic acid	STD	7.796	10.52	36.09
4	Aspartic acid	STD	7.03	10.52	36.09
5	Aspartic acid	STD	2.342	10.52	36.09
6	Aspartic acid	STD	2.21	10.52	36.09
7	Acetanilide	STD	11.756	10.36	71.09
8	Acetanilide	STD	12.772	10.36	71.09
9	Acetanilide	STD	5.233	10.36	71.09

TABLE 4. Calibration sequence for plant analysis.

Sample Information			Quadratic Fit Calibration Method			Linear Fit Calibration Method			
Reference Material	W (mg)	N%	RSD%	C%	RSD%	N%	RSD%	C%	RSD%
Birch Leaves	16.185	2.17	0.27	48.45	0.08	2.19	0.26	48.47	0.08
	15.952	2.16		48.37		2.19		48.39	
	16.746	2.16		48.41		2.20		48.44	
Orchard Leaves	16.777	2.28	0.25	50.40	0.22	2.30	0.43	50.43	0.22
	15.214	2.29		50.34		2.32		50.36	
	15.878	2.29		50.19		2.31		50.21	
Alfalfa	18.079	3.03	0.19	43.06	0.12	3.01	0.19	43.07	0.11
	20.441	3.03		43.06		3.01		43.06	
	20.585	3.02		43.15		3.00		43.15	

TABLE 5. Experimental Nitrogen and Carbon data of Plants Reference Materials.

Table 6 shows the NC data of other matrices analyzed to evaluate the repeatability. No memory effect was observed when changing the sample matrix indicating that the combustion is complete at this higher weight of sample.

Sample	W (mg)	N%	RSD%	C%	RSD%
France Maize	19.393	1.21	0	42.02	0.18
	19.908	1.21		41.96	
	20.273	1.21		41.87	
Russian Maize	19.580	1.55	0	43.14	0.14
	18.960	1.55		43.12	
	19.001	1.55		43.03	
Soja	19.626	7.93	0.15	41.82	0.77
	19.036	7.95		41.93	
	19.226	7.95		41.43	
Alfalfa	19.858	2.53	0.68	44.00	0.59
	19.767	2.56		43.48	
	19.742	2.56		43.74	
Sunflower	19.222	5.81	0.46	43.68	0.10
	19.765	5.77		43.61	
	20.313	5.82		43.69	
Oatmeal	20.354	1.87	1.43	42.56	0.22
	20.067	1.83		42.54	
	20.152	1.82		42.39	
Soja Bean	20.333	7.59	0.33	43.20	0
	20.656	7.56		43.20	
	20.179	7.54		43.20	
Barley flour	20.188	1.83	1.09	41.56	0.06
	20.349	1.81		41.55	
	20.119	1.85		41.60	
Wheat Flour	20.441	1.37	0.42	40.00	0.07
	20.513	1.38		40.03	
	19.914	1.37		40.06	
Rice Flour	20.162	1.35	1.26	40.74	0.13
	20.267	1.33		40.70	
	20.201	1.38		40.75	

TABLE 6. Experimental Nitrogen and Carbon data of other matrices.

Conclusion

Using one large reaction tube coupled with the 3 m length GC column, instead of the classical two reactors configuration, extended the analytical capabilities of the Thermo Scientific FLASH 2000 Elemental Analyzer:

- High quality of results
- Faster NC analysis
- Larger sample weight for vegetal
- Easier maintenance
- High productivity: possibility to use a single reactor in the left furnace and another reactor in the right furnace for NC/NC or NC/Sulfur determinations and others.



www.thermoscientific.com

©2016 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. All other 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 representative for details.

COMPANY WITH QUALITY MANAGEMENT
SYSTEM CERTIFIED BY DNV
= ISO 9001:2008 =

Thermo Fisher Scientific S.p.A. Milan, Italy
is ISO 9001:2008 Certified.

Africa +43 1 333 50 34 0
Australia +61 3 9757 4300
Austria +43 810 282 206
Belgium +32 53 73 42 41
Brazil +55 11 2730 3006
Canada +1 800 530 8447
China 800 810 5118 (free call domestic)
400 650 5118

Denmark +45 70 23 62 60
Europe-Other +43 1 333 50 34 0
Finland +358 10 3292 200
France +33 1 60 92 48 00
Germany +49 6103 408 1014
India +91 22 6742 9494
Italy +39 02 950 591

Japan +81 6 6885 1213
Korea +82 2 3420 8600
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
Norway +46 8 556 468 00

Russia/CIS +43 1 333 50 34 0
Singapore +65 6289 1190
Sweden +46 8 556 468 00
Switzerland +41 61 716 77 00
Taiwan +886 2 8751 6655
UK/Ireland +44 1442 233555
USA +1 800 532 4752
PN42248-EN 0216S

Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand