

# Characterization of Carbon Black with the Thermo Scientific FLASH 2000 Organic Elemental Analyzer

Liliana Krotz and Guido Giazzi  
Thermo Fisher Scientific, Rodano, Milan, Italy

## Overview

**Purpose:** Characterization of carbon black by Organic Elemental Analysis.

**Methods:** Samples were analyzed using an elemental analyzer with automated autosampler.

**Results:** CHNS/O data are shown to assess the performance of the FLASH 2000 OEA.

## Introduction

Carbon black is a material produced by the incomplete combustion of heavy petroleum products such as coal tar, ethylene cracking tar, and a small amount from vegetable oil. It is mainly used as a reinforcing filler in tires and other rubber products while in plastics, paints, and inks carbon black is used as a color pigment.

In a typical production process of carbon black the elemental content is periodically monitored and tested for quality control, in particular carbon quantification. The total carbon content is required for the calculation of carbon dioxide emissions and it can be also used in calculations to estimate yield of the process.

The Thermo Scientific™ FLASH™ 2000 CHNS/O Analyzer (Figure 1), which operates with dynamic combustion of the sample (Dumas method), provides automated and simultaneous CHNS determination in a single analysis run and it provides oxygen determination by pyrolysis in a consequent run. The dedicated Thermo Scientific™ Eager Xperience Data Handling Software calculates automatically the Gross Heat and Net Heat Values (GHV and NHV), both expressed in kcal/kg, and the CO<sub>2</sub> Emission Trade data.

## Methods

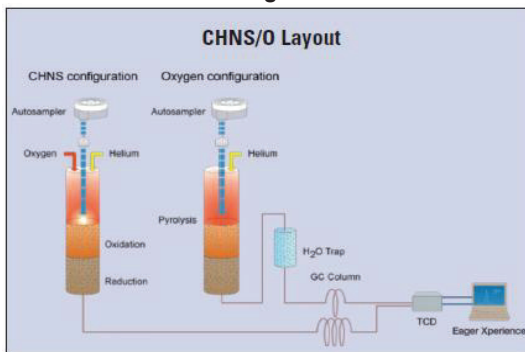
For CHNS determination, samples are weighed in tin capsules and introduced into the combustion reactor via the Thermo Scientific MAS 200R Autosampler. After combustion the resulted gases are carried by a helium flow to a layer filled with copper, then swept through a GC column that provides the separation of the combustion gases, and finally, detected by a thermal conductivity detector (TCD) (Figure 2).

For oxygen determination, samples are weighed in silver capsules and introduced into the pyrolysis chamber via the MAS 200R Autosampler. The reactor contains nickel coated carbon at 1060°C. The oxygen present in the sample, combined with the carbon, forms carbon monoxide which is then chromatographically separated from other products and detected by the TCD detector (Figure 2).

FIGURE 1. FLASH 2000 OEA CHNS/O Analyzer.



FIGURE 2. CHNS/O configuration.



## Results

Different types of carbon black samples were analyzed according to the different element content, in order to demonstrate the performance of the system and show the repeatability (without any matrix effect).

To verify the complete combustion of high containing carbon samples, carbon determination was evaluated in a Carbon Mesoporous Reference Material (99.95 C%, Aldrich No. 699624). The system was calibrated with BBOT\* as standard, 2,5-Bis (5-tert-butyl-benzoxazol-2-yl) thiophene.

Table 1 shows the data obtained with good repeatability. The average is comparable with the expected value, indicating the complete oxidation of the sample.

\*BBOT: 2,5-Bis (5-tert-butyl-benzoxazol-2-yl) thiophene

TABLE 1. Carbon determination.

| C %   | Av. C % | RSD % |
|-------|---------|-------|
| 99.91 | 99.71   | 0.143 |
| 99.58 |         |       |
| 99.86 |         |       |
| 99.71 |         |       |
| 99.68 |         |       |
| 99.69 |         |       |
| 99.82 |         |       |
| 99.72 |         |       |
| 99.69 |         |       |
| 99.41 |         |       |

Table 2 shows the repeatability of CHNS determination of carbon black samples while Table 3 shows the oxygen data and the GHV and NHV values (both expressed in kcal/kg), and the CO<sub>2</sub> Emission Trade data calculated automatically by the Eager Xperience Data Handling Software. For CHNS, carbon black samples were run in two series of analyses. In each series samples were analyzed in duplicate, the statistical data showed in the table is the average of the four determinations.

TABLE 2. CHNS data of carbon black samples.

| Sample   | N %   | RSD % | C %    | RSD % | H %   | RSD % | S %   | RSD % |
|----------|-------|-------|--------|-------|-------|-------|-------|-------|
| <b>A</b> | 0.188 | 0.676 | 96.559 | 0.111 | 0.300 | 1.843 | 0.829 | 2.193 |
|          | 0.185 |       | 96.734 |       | 0.295 |       | 0.848 |       |
|          | 0.186 |       | 96.817 |       | 0.304 |       | 0.868 |       |
|          | 0.186 |       | 96.706 |       | 0.308 |       | 0.868 |       |
| <b>B</b> | 0.273 | 0.810 | 96.178 | 0.159 | 0.267 | 2.086 | 0.823 | 0.493 |
|          | 0.277 |       | 95.899 |       | 0.280 |       | 0.814 |       |
|          | 0.272 |       | 95.351 |       | 0.272 |       | 0.815 |       |
|          | 0.273 |       | 96.079 |       | 0.277 |       | 0.817 |       |
| <b>C</b> | 0.239 | 2.155 | 95.768 | 0.196 | 0.343 | 2.697 | 1.016 | 0.311 |
|          | 0.229 |       | 96.126 |       | 0.350 |       | 1.020 |       |
|          | 0.239 |       | 95.826 |       | 0.341 |       | 1.019 |       |
|          | 0.232 |       | 95.700 |       | 0.328 |       | 1.013 |       |
| <b>D</b> | 0.148 | 1.681 | 97.073 | 0.076 | 0.282 | 1.465 | 1.273 | 0.583 |
|          | 0.146 |       | 97.066 |       | 0.285 |       | 1.267 |       |
|          | 0.149 |       | 96.966 |       | 0.292 |       | 1.256 |       |
|          | 0.152 |       | 97.146 |       | 0.286 |       | 1.261 |       |

The performance of the FLASH 2000 CHNS/O Analyzer was evaluated through the comparison of the repeatability of the CHN data obtained in comparison to the ASTM D5373 requirements showed in Table 4.

The precision of the FLASH 2000 OEA for CHN determination was evaluated from the results showed in Table 2. Table 5 shows the difference (Diff.) calculated between both data of each series of analyses. All differences are acceptable and less than the Repeatability Limit indicates in the Official Method indicating the good homogeneity and the complete combustion of the samples.

**TABLE 3. Oxygen data, Heat Values and CO<sub>2</sub> value of carbon black samples.**

| Sample   | O %   | RSD % | GHV (kcal/kg) | RSD % | NHV (kcal/kg) | RSD % | CO <sub>2</sub> E.T. | RSD % |
|----------|-------|-------|---------------|-------|---------------|-------|----------------------|-------|
| <b>A</b> | 1.149 | 0.397 | 8005          | 0.003 | 7990          | 0.003 | 105.96               | 0.000 |
|          | 1.158 |       | 8005          |       | 7989          |       | 105.96               |       |
|          | 1.152 |       | 8005          |       | 7990          |       | 105.96               |       |
| <b>B</b> | 1.661 | 0.623 | 7915          | 0.006 | 7901          | 0.006 | 106.37               | 0.005 |
|          | 1.679 |       | 7916          |       | 7901          |       | 106.38               |       |
|          | 1.661 |       | 7915          |       | 7901          |       | 106.37               |       |
| <b>C</b> | 1.597 | 0.181 | 7933          | 0.002 | 7916          | 0.002 | 106.00               | 0.005 |
|          | 1.597 |       | 7934          |       | 7916          |       | 106.00               |       |
|          | 1.602 |       | 7934          |       | 7916          |       | 106.01               |       |
| <b>D</b> | 0.624 | 0.990 | 8062          | 0.003 | 8048          | 0.003 | 105.59               | 0.006 |
|          | 0.612 |       | 8063          |       | 8048          |       | 105.58               |       |
|          | 0.616 |       | 8063          |       | 8048          |       | 105.58               |       |

**TABLE 4. Concentration Range and Limit of Repeatability accepted by ASTM D5373.**

| Element  | Concentration range (%) | Repeatability Limit r | Repeatability Limit (r): the value below which the absolute difference between two test results calculated to a dry basis of separate and consecutive test determinations, carried out on the same sample, in the same laboratory, by the same operator, using the same apparatus. |
|----------|-------------------------|-----------------------|--|
| Carbon   | 48.6 to 90.6            | 0.64                  |  |
| Hydrogen | 0.14 to 5.16            | 0.16                  |  |
| Nitrogen | 0.69 to 1.57            | 0.11                  |  |

**TABLE 5. CHN data of carbon black samples according to ASTM D5373 requirements.**

| Sample   | Serie    | N %   | Diff. | C %    | Diff. | H %   | Diff. |
|----------|----------|-------|-------|--------|-------|-------|-------|
| <b>A</b> | <b>1</b> | 0.188 | 0.003 | 96.559 | 0.175 | 0.300 | 0.005 |
|          |          | 0.185 |       | 96.734 |       | 0.295 |       |
|          | <b>2</b> | 0.186 | 0.000 | 96.817 | 0.111 | 0.304 | 0.004 |
|          |          | 0.186 |       | 96.706 |       | 0.308 |       |
| <b>B</b> | <b>1</b> | 0.273 | 0.004 | 96.178 | 0.279 | 0.267 | 0.013 |
|          |          | 0.277 |       | 95.889 |       | 0.280 |       |
|          | <b>2</b> | 0.272 | 0.001 | 95.851 | 0.228 | 0.272 | 0.005 |
|          |          | 0.273 |       | 96.079 |       | 0.277 |       |
|          |          |       |       |        |       |       |       |

The Thermo Scientific FLASH 2000 Elemental Analyzer proved to be a valuable solution for the elemental analysis of carbon black in terms of:

- Accuracy, repeatability and sensitivity of results.
  - Its automation and high speed of analysis allow efficient analysis and help reduce overall operational costs.
  - All data were obtained according to the ASTM D5373 method.
- No matrix effect was observed when changing the sample, indicating complete combustion.

[www.thermoscientific.com](http://www.thermoscientific.com)

©2015 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.



Thermo Fisher Scientific,  
Sunnyvale, CA USA is  
ISO 9001 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 3731 5140  
**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  
PN42237-EN 1015S

**Thermo**  
SCIENTIFIC  
A Thermo Fisher Scientific Brand