

Elemental Analysis: CHNS/O determination in carbon

Authors

Dr. Liliana Krotz and
Dr. Guido Giazzi
Thermo Fisher Scientific,
Milan, Italy

Keywords

Coal, Coke, Carbon Black,
Graphite, CHNS/O, Heat Value

Goal

This application note reports data on CHNS/O determination on carbon samples needed for quality control purposes, performed with the FlashSmart EA.

Introduction

Carbon occurs as a variety of allotropes. There are two crystalline forms, diamond and graphite, and a number of amorphous (non-crystalline) forms, such as charcoal, coke, and carbon black. The most common use of carbon black is as a pigment and reinforcing phase in automobile tires. Coke is the solid carbonaceous material derived from destructive distillation of low-ash, low-sulfur bituminous coal. Coke is also used as a fuel and as a reducing agent in smelting iron ore in a blast furnace.

For quality control purposes, the organic elements in carbon need to be determined. For the determination of carbon, hydrogen, nitrogen, sulfur and oxygen, the combustion method is used.

The Thermo Scientific™ FlashSmart™ Elemental Analyzer (Figure 1) allows the quantitative determination of carbon, hydrogen, nitrogen and oxygen in carbon. The FlashSmart EA based on the dynamic flash combustion of the sample, provides automated and simultaneous CHNS determination in a single analysis run and oxygen determination by pyrolysis in a second run. To perform total sulfur determination at trace levels, the analyzer has been coupled with the Flame Photometric Detector (FPD).

Methods

For CHNS determination, the FlashSmart EA operates according to the dynamic flash combustion of the sample. Liquid samples are weighed in tin containers and introduced into the combustion reactor via the Thermo Scientific™ MAS Plus Autosampler. Samples can be directly injected, with a syringe via the Thermo Scientific™ AS 1310 Liquid Autosampler. In both cases a regulated amount of oxygen was used. After combustion, the resultant gases are carried by a helium flow to a layer filled with copper, then swept through a GC column that separates the combustion gases. Finally they are detected by a Thermal Conductivity Detector (TCD). Total run time is less than 10 minutes (see Figure 2). For trace sulfur determination, the gases produced by combustion are carried by a helium flow to a layer filled with copper, then swept through a water trap, a short GC column before the sulfur is measured by the Flame Photometric Detector (FPD). Total run time is 5 minutes (Figure 3).

For oxygen determination, the system operates in pyrolysis mode. Samples are weighed in silver containers and introduced into the pyrolysis chamber via the MAS Plus Autosampler or directly injected via the AI 1310 or AS 1310 Autosamplers. The reactor contains nickel coated carbon at 1060 °C. The oxygen 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). A complete report is automatically generated by the Thermo Scientific™ EagerSmart™ Data Handling Software.

Results

Different carbon samples were chosen to show the reproducibility obtained with the system. Coal, coke, graphite, lignite and carbon black samples were homogenized by a ball mill.

Tables 1 and 2 show CHNS/O and CHNS determination of different matrices. Instrument calibration was performed with 2-3 mg of BBOT (2, 5-Bis (5-ter-butyl-benzoxazol-2-yl) thiophene). No matrix effect was observed when changing the nature of sample.

Table 1 also indicates the Gross Heat Value (GHV in kcal/kg) and Net Heat Value (NHV in kcal/kg) calculated automatically by the dedicated EagerSmart Xperience software.



Figure 1. FlashSmart Elemental Analyzer coupled with Flame Photometric Detector (FPD).

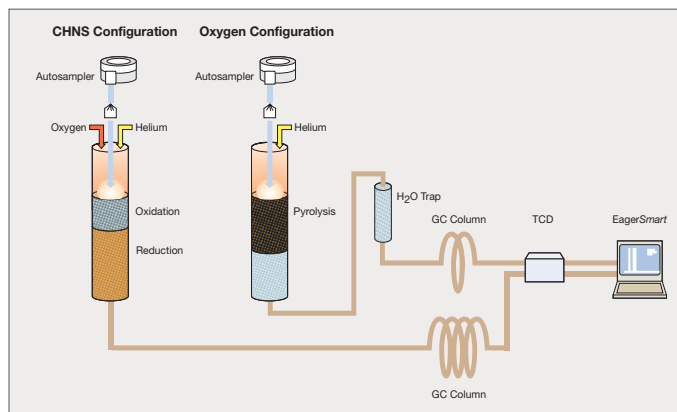


Figure 2. CHNS/O configuration with TCD Detector.

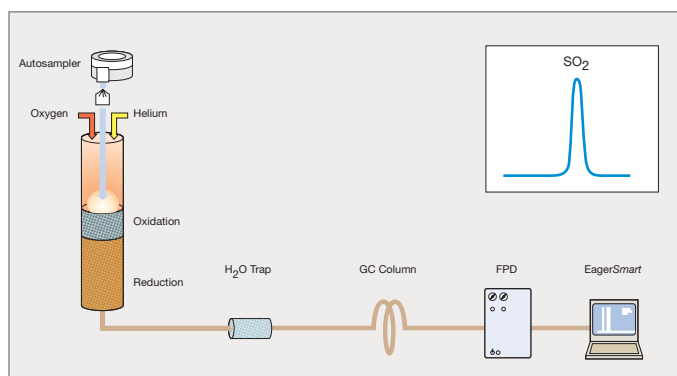


Figure 3. Sulfur configuration with FPD Detector.

Table 1. CHNS/O determination and Heat Value calculation.

Sample	N%	C%	H%	S%	O%	GHV	NHV
Coal	1.740	73.006	5.389	0.838	12.733	7316	7040
	1.732	72.950	5.358	0.857	12.731	7316	7040
	1.747	73.238	5.398	0.828	12.813	7313	7037
RSD%	0.431	0.209	0.390	1.752	0.367	0.024	0.025
Hard coal	1.287	80.137	4.621	0.488	4.903	7957	7720
	1.288	80.123	4.513	0.486	5.085	7918	7687
	1.329	80.706	4.617	0.497	5.094	8003	7766
RSD%	1.842	0.414	1.336	1.195	2.144	0.535	0.514
Brown coal	1.988	77.258	3.280	0.386	3.730	7131	7145
	2.006	77.196	3.266	0.386	3.723	7304	7136
	2.028	77.823	3.292	0.397	3.690	7364	7195
RSD%	0.998	0.446	0.397	1.630	0.575	0.442	0.444
Lignite 1	0.758	62.281	4.636	0.377	25.202	5589	5357
	0.750	62.357	4.544	0.373	25.540	5574	5343
	0.757	62.24	4.377	0.371			
RSD%	0.559	0.074	2.899	0.724	0.942	0.184	0.192

Table 2. CHNS determination.

Sample	N%	RSD%	C%	RSD%	H%	RSD%	S%	RSD%
Pet coke	1.486	2.266	96.329	0.313	0.210	8.095	0.641	0.384
	1.534		96.756		0.235		0.643	
Coke 1	1.229	0.011	86.195	0.384	4.343	0.346	0.656	4.535
	1.228		86.665		4.322		0.616	
Coke 2	0.377	1.360	98.997	0.172	---	---	0.422	1.017
	0.370		98.917				0.431	
	0.374		98.576				0.429	
	0.363		98.973				0.434	
	0.369		98.836				0.429	
Calcinated coke	1.086	0.295	96.934	0.142	0.0741	2.810	1.213	0.262
	1.086		97.161		0.0703		1.219	
	1.086		97.267		0.0716		1.213	
	1.089		96.965		0.0738		1.213	
	1.093		97.081		0.0754		1.211	
Carbon black 1	0.189	1.968	95.693	0.181	0.321	0.811	0.296	2.075
	0.186		95.629		0.323		0.288	
	0.185		95.825		0.322		0.292	
	0.181		95.835		0.319		0.280	
	0.180		96.082		0.317		0.290	
Carbon black 2	0.131	2.493	95.609	0.124	0.384	3.861	1.433	0.649
	0.137		95.542		0.361		1.452	
	0.133		95.772		0.359		1.443	
Carbon black 3	0.135	2.156	96.329	0.0465	0.420	1.496	0.711	0.505
	0.134		96.241		0.412		0.704	
	0.140		96.300		0.407		0.709	
Carbon black 4	0.172	0.331	94.591	0.0693	0.520	0.458	1.741	0.841
	0.171		94.512		0.520		1.728	
	0.172		94.642		0.524		1.757	
Carbon Black 5	0.0896	3.178	99.365	0.079	0.322	0.452	---	---
	0.0843		99.497		0.319			
	0.0884		99.506		0.321			
Graphite	---	---	99.889	0.157	---	---	---	---
			99.580					
			99.505					
			99.725					
			99.794					

Table 3. Trace sulfur determination by FPD Detector.

Sample		Sulfur		Sample		Sulfur	
Nature		ppm S	RSD%	Nature		ppm S	RSD%
Coke		398	1.507	Graphite		46	1.264
		404				46	
		392				45	

Table 3 shows trace level sulfur data obtained when using the FPD Detector. Solid samples were weighed in tin containers with the addition of Vanadium Pentoxide, a typical “oxygen donor” that allows the total conversion of sulfur.

Conclusions

For CHNS/O determination for carbon characterization samples, the FlashSmart EA performs accurate analysis. Data were obtained with good reproducibility and no matrix effect when changing the sample.

CHNS determination can be performed in a single run with the FlashSmart EA. By coupling the FlashSmart EA with the FPD Detector, trace levels of sulfur can be analyzed. Oxygen determination can also be performed. Alongside with reproducibility of the data, the Elemental Analyzer offers advantages also in terms of reduced operational costs and analysis efficiency.

Find out more at thermofisher.com/OEA