

Elemental Analysis: CHNS/O characterization of biomass and bio-fuels

Authors

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

Keywords

Biomass, Bio-fuels, CHNS/O,
CO₂ Emission Trade, Elemental
Analysis, Heat Values

Goal

This application note reports data on CHNS/O determination of biomass and bio-fuels for quality control purposes, performed with the FlashSmart EA.

Introduction

Biomass fuels are derived from biological material rather than being the result of geological processes. The biological material takes carbon from the atmosphere and returns it when it is burnt. Biomass derives from burning crops, manure, garbage and most commonly, wood. Burning wood, for the production of energy has been common practice both domestically and industrially for heating and cooking, or to produce steam and electricity for industrial use.

Bio-fuels are an alternative to fossil fuels and are derived mainly from plants and animals, which do not increase the carbon dioxide level in the atmosphere. These alternative fuels are used globally, most commonly for automotive transport. Many manufacturing plants and research laboratories are involved in the characterization of these bio-fuel products to establish their viability for production, carbon dioxide emission levels, impact on water resources, energy balance and efficiency. A common method to characterize these materials is by identifying their energetic value via elemental composition.

As the demand for biomass and bio-fuels testing has grown in the last years, the classical analytical methods showed to be no longer suitable, for their time-consuming sample preparation and for their use of hazardous reagents. For this reason a simple and automated technique is the requirement for modern laboratories dealing with routine analysis.

The Thermo Scientific™ FlashSmart™ Elemental Analysis (Figure 1), based on combustion method, allows the quantitative determination of the elements in a large range of concentration, without the need for sample digestion or toxic chemicals, providing advantages in terms of time and automation.

This method combines the advantages of the elemental analyzer with the sensitivity, selectivity and robustness of Flame Photometric Detector (FPD). The coupling is simple and it allows sulfur determination without any matrix effect.

The Thermo Scientific™ EagerSmart™ EA provides dedicated features for the automated calculation of the Heat Values and the CO₂ Emission Trade.

This note presents data on CHNS/O determination in biomass and bio-fuels samples to show the performance of the system and to show the reproducibility of the results obtained.

Methods

For CHNS determination the FlashSmart Elemental Analyzer operates according to the dynamic flash combustion of the sample. Samples are weighed in tin containers and introduced into the combustion reactor via the Thermo Scientific™ MAS Plus Autosampler with oxygen. After combustion the resulted gases are carried by a helium flow to a layer filled with copper, then swept through a GC column which provides the separation of the combustion gases, and finally, detected by a Thermal Conductivity Detector (TCD) (Figure 2). Total run time is less than 10 minutes.

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 and finally the sulfur is measured by the Flame Photometric Detector (FPD). Total run time 5 min. (Figure 3).

For oxygen determination samples are weighed in silver containers and introduced into the pyrolysis chamber via the MAS Plus Autosampler. The reactor contains nickel coated carbon at a temperature of 1060 °C. The oxygen in the sample, combined with the carbon, forms carbon monoxide which is 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.



Figure 1. FlashSmart Elemental Analyzer.

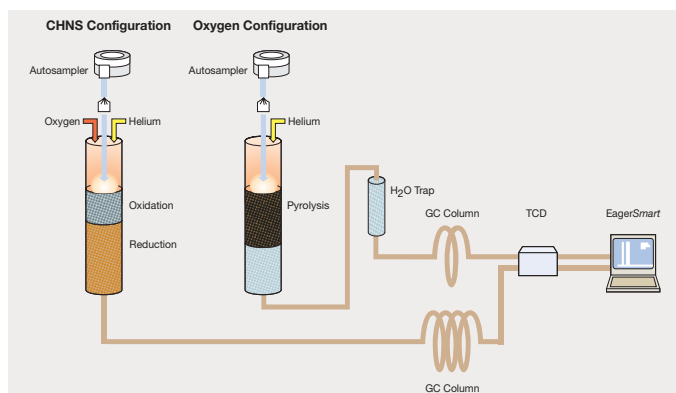


Figure 2. CHNS/O configuration.

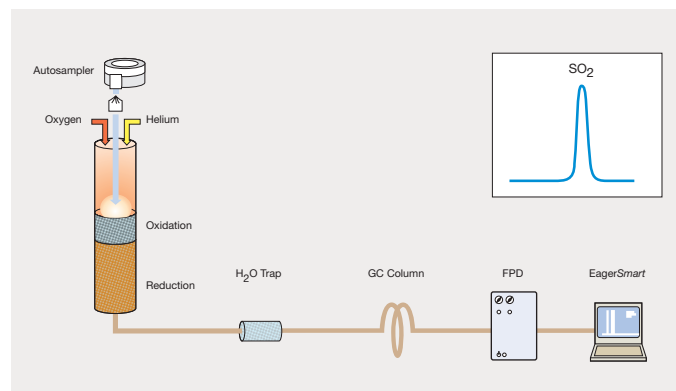


Figure 3. Sulfur configuration by FPD Detector.

Results

Different alternative fuels were chosen to show the performance obtained with the system.

The FlashSmart EA analyzer was calibrated by BBOT (2,5-Bis(5-ter-butyl-benzoxazol-2-yl)thiophene) standard for CHNS determination and by benzoic acid for oxygen determination. No matrix effect was observed when changing the nature of sample.

The EagerSmart Data Handling Software automatically calculates the GHV (Gross Heat Value in kcal/kg) and NHV (Net Heat Value in kcal/kg) using the modified Dulong equations, as well as the CO₂ Emission Trade.

Table 1 shows the CHNS data obtained from different biomass samples. Samples were homogenized by a ball mill.

Table 1. CHNS determination in biomass samples.

Sample	N%	RSD%	C%	RSD%	H%	RSD%	S%	RSD%
Biomass 1	1.382	0.977	40.091	0.025	6.421	0.326	0.087	2.012
	1.369		40.090		6.404		0.086	
	1.396		40.073		6.446		0.089	
Biomass 2	1.187	1.499	39.999	0.537	5.598	0.634	0.112	0.259
	1.181		40.193		5.574		0.111	
	1.215		39.764		5.644		0.112	
Stray Pellets	0.5282	1.452	43.278	0.075	5.695	1.008	0.149	3.631
	0.5160		43.310		5.759		0.149	
	0.5300		32.245		5.644		0.140	
Sunflower Pellets	0.712	0.498	45.531	0.158	5.716	1.049	0.125	1.791
	0.719		45.590		5.715		0.124	
	0.717		45.675		5.612		0.121	

Table 2. CHNS/O determination in vegetals as biofuels.

Sample	N%	RSD%	C%	RSD%	H%	RSD%	S%	RSD%	O%	RSD%
1	5.521	3.133	47.119	0.388	6.010	1.649	0.265	2.647	34.311	0.444
	5.790		47.276		5.821		0.277		34.045	
	5.861		47.485		5.959		0.278		34.306	
2	3.534	2.816	46.467	0.153	5.876	0.609	0.182	2.776	37.183	0.357
	3.528		46.489		5.946		0.176		37.421	
	3.706		46.600		5.926		0.186		37.199	

Table 3. Heat Values and CO₂ Emission Trade calculation in vegetals as bio-fuels.

Sample	G.H.V.	RSD%	N.H.V.	RSD%	CO ₂ e.t.	RSD%
1	4449	0.149	4145	0.167	99.883	0.157
	4461		4157		88.608	
	4450		4145		99.877	
2	4255	0.136	3952	0.139	103.056	0.145
	4245		3942		103.323	
	4255		3951		103.073	

Table 4 shows CHNS/O determinations of different wood samples while Table 5 includes the GHV (Gross Heat Value in kcal/kg), NHV (Net Heat Value in

kcal/kg) and the CO₂ emission trade. The sample data in Table 5 were obtained using the average of the element percentages.

Table 4. CHNS/O analysis of wood samples.

Sample	N%	C%	H%	S%	O%
1	0.0649	45.7746	6.1044	0.0288	47.7422
	0.0677	45.8869	6.2031	0.0274	42.7301
	0.0666	46.1076	6.1345	0.0285	42.4861
	0.0651	45.7568	6.1504	0.0275	42.5882
	0.0664	45.8578	6.0941	0.0279	42.4200
Av	0.0661	45.8767	6.1373	0.0280	42.5933
RSD%	1.7463	0.3055	0.7035	2.1913	0.3370
2	0.0667	46.6378	5.9850	0.0246	42.4833
	0.0662	47.0700	5.9247	0.0249	42.4623
	0.0659	46.8560	5.9086	0.0260	42.6421
	0.0630	46.8802	5.9816	0.0254	42.5727
	0.0660	46.9996	6.0158	0.0259	42.3615
Av	0.0656	46.8887	5.9631	0.0254	42.5044
RSD%	2.2329	0.3525	0.7520	2.6259	0.2530
3	0.0882	47.1092	6.1197	0.0312	43.2662
	0.0910	47.0881	6.1050	0.0324	43.1915
	0.0895	47.1071	6.1088	0.0315	43.1879
	0.0902	47.0995	6.1115	0.0331	43.3391
	0.0890	47.1042	6.1065	0.0319	43.2539
Av	0.0896	47.1016	6.1103	0.0320	43.2477
RSD%	1.2034	0.0178	0.0949	2.3516	0.1438

Table 5. Heat Values and CO₂ Emission Trade calculation in wood samples.

Sample	G.H.V. (kcal/kg)	N.H.V. (kcal/kg)	CO ₂ Emission Trade
1	4042	3728	108
2	4069	3763	109
3	4105	3792	109

Table 6 shows the sulfur data of wood samples obtained with the FPD Detector, where the sulfur content is at trace levels. Solid samples were weighed in tin containers with the addition of vanadium pentoxide. Instrument calibration was performed with a soil sample, of which the sulfur value (30 ppm) is the average from an International Round Robin Test (WEPAL, Wageningen Evaluating Programs for Analytical Laboratories, Wageningen University, Netherlands).

Table 6. Trace sulfur determination of wood samples by FPD Detector.

Sample	S ppm	Average S ppm	RSD%
A	36	36	3.64
	35		
	37		
	37		
	34		
B	44	42	2.60
	41		
	42		
	42		
C	52	53	2.45
	54		
	52		
	55		
	53		

Table 7 shows the CHN/O data of pyrolysis oils. Due to the low concentration of sulfur, determination was performed with the FPD Detector and calibration was

performed with a Soil Reference Material (320 ppm). The Sulfur results obtained are shown in Table 8.

Table 7. CHN/O determination in biomass pyrolysis oils.

Sample	N%	RSD%	C%	RSD%	H%	RSD%	O%	RSD%
1	0.0785	0.7788	42.1731	0.2881	7.4361	0.3150	46.9729	0.6969
	0.0774		41.9498		7.4751		47.5645	
	0.0784		42.1433		7.4330		47.0198	
2	0.0856	1.8799	46.5296	0.4357	6.9718	0.2253	43.5343	0.2671
	0.0830		46.2804		6.9984		43.6605	
	0.0859		46.6817		6.9997		43.4282	

Conclusions

For the CHNS/O determination in biomass and bio-fuels the FlashSmart Elemental Analyzer provides accurate and reproducible results.

Thanks to the dedicated features of the EagerSmart Data Handling Software, the Heat Values and the estimated CO₂ Emission Trade value are automatically provided.

The modularity of the FlashSmart Elemental Analyzer enables to perform CHNS determination in a single run, oxygen determination in a second run and TOC analysis without any modification of the analytical conditions, allowing high productivity for routine labs.

The FlashSmart Elemental Analyzer meets laboratory requirements in terms of automation and high sample throughput.

Table 8. Sulfur determination by FPD Detector in biomass pyrolysis oils.

Sample	S%	RSD%
1	0.0186	2.6961
	0.0189	
	0.0196	
2	0.0191	1.6279
	0.0185	
	0.0187	

Find out more at thermofisher.com/OEA