



Characterization of food and animal feed related products by the Thermo Scientific FlashSmart Elemental Analyzer

Authors

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

Keywords

Animal Feed, CHNS, Food
Quality, Food Supplements,
Labeling, QA/QC

Goal

To demonstrate the performance of the Thermo Scientific FlashSmart Elemental Analyzer for nitrogen/protein determination in food and animal feed.

Introduction

Food and animal feed is made up of chemical compounds that determine flavor, color, texture and nutritional value, and are carefully regulated by federal authorities and various international organizations to ensure that they are safe to eat and are accurately labelled.

One of the main analyses for quality control and R&D purposes is elemental characterization. The determination of nitrogen, carbon, hydrogen and sulfur, provides useful information on the characterization of these materials. It is therefore very important to have an accurate and precise technique, preferably automatic, that allows fast analysis with excellent reproducibility. The Thermo Scientific™ FlashSmart™ Elemental Analyzer (Figure 1) copes effortlessly with the wide array of laboratory requirements such as accuracy, precision and day to day reproducibility.



Figure 1. Thermo Scientific FlashSmart Elemental Analyzer.

Methods

For CHNS determinations, the elemental analyzer operates according to the dynamic flash combustion of the sample. The sample is weighed in a tin capsule and introduced into the combustion reactor via the Thermo Scientific™ MAS Plus Autosampler with oxygen. 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 being detected by a Thermal Conductivity Detector (TCD) (Figure 2). Total run time is less than 10 minutes. For NCS or for sulfur only determination, the water produced during combustion is adsorbed through a H₂O trap before entering the GC column.

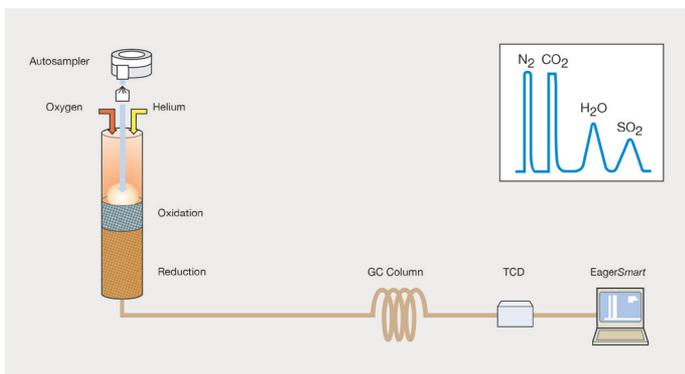


Figure 2. FlashSmart CHNS configuration.

For NC determination, after combustion, the produced gases are carried by a helium flow to a second reactor filled with copper, then swept through a H₂O trap, a GC column, before finally being detected by a thermal conductivity detector. Total run time is less than five minutes (see Figure 3).

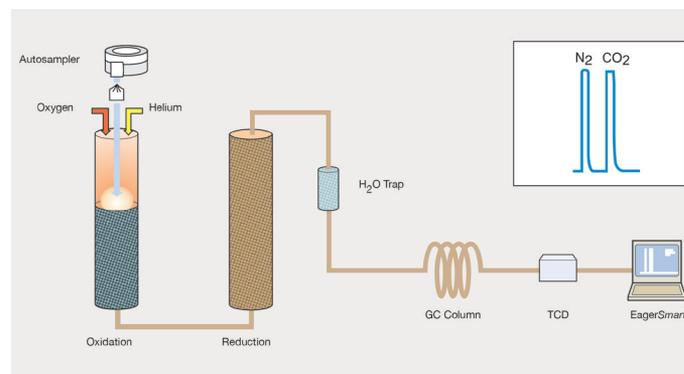


Figure 3. FlashSmart NC configuration.

A complete report is automatically generated by the Thermo Scientific™ EagerSmart™ Data Handling Software and displayed at the end of the analysis.

Results

Different food and animal feed related products with a large range of elemental concentrations were analyzed in various configurations, to show the performance of the instrument in terms of repeatability.

For the CHNS, NCS and sulfur only determinations, the addition of Vanadium Pentoxide (oxygen donor) was used for a complete conversion of the sulfur, and the instrument was calibrated with the standards BBOT* and nicotinamide. For NC determination, acetanilide and aspartic acid were used as standards to calibrate the instrument. In all cases, K factor was used as the calibration method.

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

Table 1 shows the CHNS data obtained from different sample matrices. The weight of sample was 2–3 mg for animal gelatines and food supplement A, 3–4 mg for starch and food supplement C, 8–10 mg for food supplement B.

No memory effect was observed when changing the sample nature, indicating complete combustion of all samples with quantitative determination of the elements.

Table 1. CHNS data.

Sample	N%	RSD%	C%	RSD%	H%	RSD%	S%	RSD%
Fish gelatine	16.249	0.185	43.023	0.089	6.902	2.632	0.394	2.004
	16.212		43.099		6.608		0.408	
	16.189		43.051		6.586		0.408	
Bovine gelatine	15.796	0.148	44.615	0.037	6.623	0.309	0.531	0.601
	15.835		44.647		6.658		0.536	
	15.838		44.624		6.622		0.537	
Porcine gelatine	16.088	0.226	44.460	0.096	6.631	0.585	0.531	0.970
	16.016		44.397		6.659		0.536	
	16.043		44.379		6.582		0.537	
Starch	2.530	0.329	31.008	0.204	5.396	1.004	0.399	0.902
	2.516		30.850		5.456		0.396	
	2.537		31.000		5.415		0.391	
	2.520		30.956		5.373		0.392	
	2.528		30.967		5.310		0.398	
Food supplement A	13.168	0.137	52.179	0.104	6.665	0.311		
	13.160		52.084		6.626			
	13.194		52.178		6.626			
Food supplement B	0.071	1.109	8.0197	0.036	2.207	0.254		
	0.071		8.0251		2.199			
	0.073		8.0241		2.209			
Food supplement C	0.330	0.533	40.615	0.323	6.274	0.488	0.366	1.581
	0.332		40.396		6.331		0.368	
	0.329		40.630		6.323		0.357	

Table 2 shows the NCS data obtained from different sample matrices. The weight of the sample was 3–4 mg.

No memory effect was observed when changing the sample nature, indicating complete combustion of all samples with quantitative determination of the elements.

Table 2. NCS data.

Sample	N%	RSD%	C%	RSD%	S%	RSD%
Animal feed 1	2.525	1.850	40.432	0.270	0.160	5.723
	2.523		40.342		0.143	
	2.444		40.215		0.155	
Animal feed 2	3.892	1.069	44.752	0.776	0.287	1.773
	3.845		44.893		0.282	
	3.810		44.235		0.277	
Meat 1	13.726	1.199	50.315	0.298	0.776	0.901
	14.057		50.107		0.790	
	13.931		50.026		0.784	
Meat2	12.939	0.286	51.064	0.206	0.724	0.742
	12.979		50.886		0.732	
	13.013		50.879		0.721	
Meat 3	12.255	0.094	53.381	0.118	0.764	0.798
	12.243		53.486		0.774	
	12.232		53.495		0.764	
Meat 4	12.451	0.399	52.603	0.111	0.785	0.199
	12.369		52.701		0.783	
	12.459		52.597		0.781	
Potato tuber	0.874	1.330	39.542	0.229	0.0709	1.221
	0.893		39.528		0.0698	
	0.896		39.382		0.0716	
	0.868		39.389		0.0721	
	0.883		39.348		0.0709	

Table 3 shows the sulfur data obtained from different soya and maize samples. The weight of the sample was 3–4 mg.

Table 3. Sulfur data.

Sample	S%	RSD%
Soya 1	0.356 – 0.336 – 0.338 – 0.350 – 0.343 – 0.341	2.214
Soya 2	0.351 – 0.344 – 0.343	1.635
Soya 3	0.372 – 0.373 – 0.363	1.725
Soya 4	0.366 – 0.363 – 0.364	0.388
Maize 1	0.115 – 0.114 – 0.116 – 0.113 – 0.113 – 0.106	3.137
Maize 2	0.119 – 0.111	4.919
Maize 3	0.104 – 0.104	0
Maize 4	0.112 – 0.102	6.608
Maize 5	0.102 – 0.100	1.396

Table 4 shows the NC data of blood flour, meat flour and food supplements. The weight of sample was 9–10 mg for blood flour and meat flour while for food supplements the weight was 10–20 mg.

Table 4. NC data.

Sample	N%	RSD%	C%	RSD%
Blood flour	14.536 14.307 14.427	0.797	48.791 49.238 49.052	0.458
Meat flour	7.213 7.474 7.385	1.804	32.008 33.287 32.722	1.963
Food supplement 1	5.739 5.747 5.744	0.071	14.598 14.623 14.612	0.084
Food supplement 2	11.365 11.391 11.369	0.123	45.625 45.303 45.343	0.385
Food supplement 3	5.845 5.831 5.863	0.269	17.127 17.115 17.161	0.138
Food supplement 4	3.630 3.613 3.634	0.310	37.093 36.975 36.956	0.201

Conclusions

The FlashSmart EA allows the quantitative recovery of the elements from any matrix with no memory effect observed when changing the sample. The advantage of the FlashSmart EA lies in its ability to perform NC determination or simultaneous CHNS determination in a single run. By a simple modification of the CHNS configuration, the analysis of NCS or sulfur only can be performed using the same analytical conditions.

Find out more at thermofisher.com/OEA