



Characterization of lubricants and oils by the Thermo Scientific FlashSmart Elemental Analyzer

Authors

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

Keywords

ASTM D5291, CHNS,
Combustion, Lubricants,
Nitrogen, Oils

Goal

This application note reports data on determination of nitrogen in lubricants and oil samples and shows the performance of the FlashSmart Elemental Analyzer and its reproducibility capabilities.

Introduction

In a typical production process of mineral oils, the nitrogen content is periodically monitored and tested for quality control. The reproducibility of data, which is measured as deviation of results from the mean value, is one of the main targets in all analytical tests for the alternative techniques accepted.

The method for nitrogen analysis in lubricants, based on combustion, is described in ASTM D5291. The ASTM method has been established for samples containing nitrogen in percentages from <0.1% to 2% (mass) with a repeatability of 0.16% (mass). The method is for the determination of carbon, hydrogen and nitrogen in laboratory samples of lubricants and petroleum products. Using the Test method D levels of 0.01 N% in lubricants can be determined.

As the demand for improved sample throughput, reduction of operational costs and minimization of human errors is steadily increasing, a simple and automated technique, which allows fast analysis with an excellent reproducibility is the key for efficient nitrogen analysis. The Thermo Scientific™ FlashSmart™ Elemental Analyzer (Figure 1), which uses the dynamic combustion of the material, requires no sample digestion or toxic chemicals, and provides important advantages in terms of time, automation and quantitative determination of nitrogen in concentrations between 0.01% (100 ppm) to higher than 5% (mass). Simultaneous determinations of nitrogen, carbon, hydrogen and sulfur are also possible with a simple upgrade.



Figure 1. Thermo Scientific FlashSmart Elemental Analyzer.

Method

The FlashSmart Elemental Analyzer uses the dynamic flash combustion of the samples, which are weighed in tin capsules and introduced into the combustion reactor via the Thermo Scientific™ MAS Plus Autosampler with oxygen.

After the combustion, for nitrogen determination only, the produced gases are carried to a second reactor filled with copper by a helium flow, then swept through CO₂ and H₂O traps, a GC column and finally detected by a Thermal Conductivity Detector (TCD). Total run time is less than 5 minutes (Figure 2).

For simultaneous CHNS, after the combustion the resulted gases are carried to a layer containing copper by a helium flow, then swept through a GC column that provides the separation of the combustion gases, and finally, detected by a TCD Detector. Total run time is less than 10 minutes (Figure 3). For NCS or sulfur determination only a water trap is installed between the reactor and the GC column.

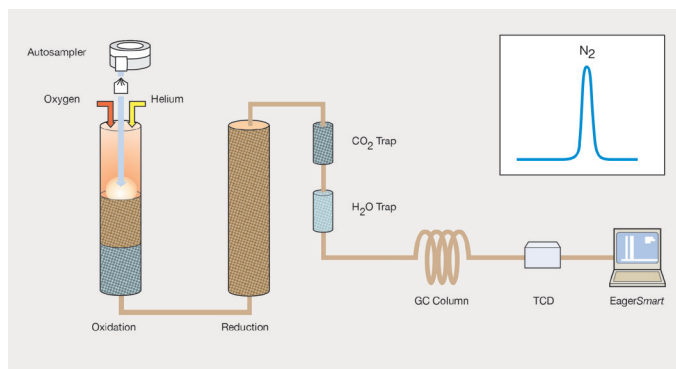


Figure 2. Nitrogen Lubricant configuration.

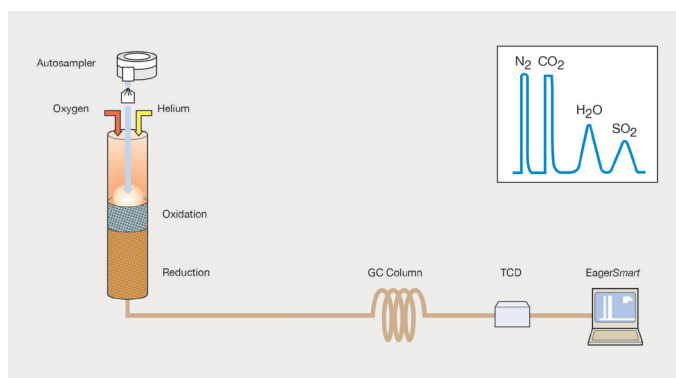


Figure 3. CHNS configuration.

A report is generated by the Thermo Scientific™ EgerSmart™ Data Handling Software.

Results

Different lubricants and oils were analyzed several times to show the reproducibility of nitrogen determination and simultaneous CHNS analysis. Samples were weighed in tin containers without pre-treatment.

In nitrogen determination only, the calibration can be performed with 4-5 mg atropine (4.84 N%) using K factor as calibration method, or by Linear Fit using a Lubricant Reference Material. Figure 4 shows a Linear Fit calibration curve. The SPX Lubricant Reference Material (1.11 N%) weighted between 0.2 mg and 10 mg. Table 1 shows the nitrogen repeatability of SPX lubricant analyzed as unknown and another lubricant at 0.08 N% analyzed in five days.

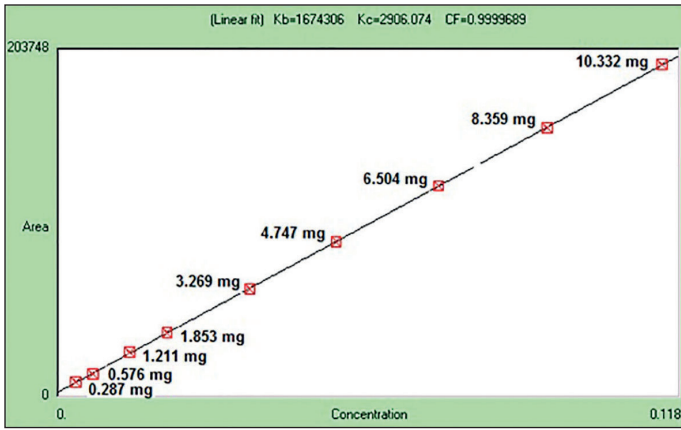


Figure 4. Nitrogen Linear Fit Calibration Curve.

Table 1. Day by day reproducibility in N determination of lubricants.

Sample	Day 1	Day 2	Day 3	Day 4	Day 5	Average N%	RSD%
SPX 1.11 N%	1.1076	1.0978	1.0994	1.0994	1.1137	1.1036	0.6
Lub 0.08 N%	0.0779	0.0815	0.0785	0.0818	0.0787	0.0797	2.2

Table 2 shows the nitrogen repeatability of lubricant samples in a wide range of concentrations analyzed in two days. No matrix effect was observed when changing sample.

Table 2. Day-by-day reproducibility in N determinations of several lubricant 3 samples.

Sample	Day 1 – N%	Day 2 – N%
1	1.130	1.130
2	1.073	1.073
3	1.596	1.594
4	1.173	1.163
5	0.874	0.866
6	1.198	1.210
7	1.123	1.117
8	0.579	0.573
9	0.878	0.888

Table 3 shows the nitrogen repeatability of three determinations of lubricant samples, containing more than 0.3% nitrogen. Table 4 shows the nitrogen repeatability of 10 determinations of lubricant samples with a nitrogen content between 300 ppm N and 0.20 N%.

Table 3. Nitrogen reproducibility of three determinations of lubricant samples.

Sample	RSD%	N%
1	0.3154	0.3047 0.3065 0.3062
2	0.6253	0.4580 0.4539 0.4594
3	0.9414	0.4700 0.4788 0.4727
4	0.4275	0.5200 0.5237 0.5197
5	0.7969	0.6400 0.6349 0.6451
6	0.7908	0.8461 0.8420 0.8331
7	0.5825	1.4617 1.4664 1.4783
8	0.3135	3.9927 3.9683 3.9760

Table 4. Nitrogen reproducibility of 10 determinations of lubricant samples.

Sample	1	2	3	4	5
N%	0.0347	0.0468	0.1051	0.1456	0.1800
	0.0323	0.0446	0.1147	0.1488	0.1819
	0.0309	0.0465	0.1158	0.1482	0.1849
	0.0298	0.0493	0.1078	0.1507	0.1860
	0.0358	0.0495	0.1147	0.1424	0.1811
	0.0323	0.0463	0.1059	0.1535	0.1849
	0.0316	0.0450	0.1084	0.1436	0.1858
	0.0304	0.0469	0.1177	0.1535	0.1894
	0.0329	0.0476	0.1089	0.1494	0.1834
	0.0292	0.0431	0.1121	0.1514	0.1886
Average N%	0.0320	0.0466	0.1101	0.1487	0.1846
RSD%	6.54	4.28	3.54	2.58	1.66

Table 5 shows the nitrogen repeatability of 10 determinations in lubricant samples. The content is to be found at trace level (less than 200 ppm N).

Table 5. Reproducibility in trace N determination of lubricant samples.

Sample	ppm N	Average ppm%	RSD%
1	193	189	5.03
	191		
	209		
	194		
	175		
	182		
	179		
	187		
	192		
	186		
2	126	129	6.63
	138		
	110		
	134		
	132		
	127		
	135		
	134		
	121		
	136		

Table 6 shows the nitrogen repeatability of three determinations of different oil samples.

Table 6. Nitrogen reproducibility of three determinations of oil samples.

Sample	N%	RSD%
Heavy Gas Oil 1	0.272	1.5851
	0.280	
	0.273	
Heavy Gas Oil 2	0.242	1.1767
	0.247	
	0.247	
Heavy Gas Oil 3	0.069	2.9182
	0.073	
	0.072	
Slurry Oil	0.213	1.2363
	0.212	
	0.217	
Residual Oil	0.276	1.9454
	0.274	
	0.266	
Vacuum Gas Oil	0.098	3.3254
	0.093	
	0.099	

To verify the performance of the FlashSmart EA results were compared with data provided by the **ASTM International Interlaboratory Program**. ASTM periodically compares test results and calculated statistical parameters with data provided by laboratories of the global petrochemical and analytical community, in order to help laboratories monitoring their performance when running **ASTM methods**. Laboratories involved in the program receive different lubricants samples and they are requested to analyze the samples according to ASTM methods. Results are first collected and then processed by ASTM and reports are sent out to each participant.

Table 7 shows the comparison of the FlashSmart EA data and the statistical results obtained by ASTM. All data obtained fall within the range indicated by the ASTM reports. The calibration of the system was performed with atropine (4.84%N) using K factor as calibration method. The samples were analyzed as received.

Table 7. Nitrogen reproducibility of ASTM lubricants.

Sample	Thermo Fisher	ASTM	ASTM Reproducibility
	N% value	Robust Mean N%	These Test Data
1	0.675	0.6685	0.0604
2	0.700	0.7088	0.0471
3	0.576	0.5790	0.0604
4	0.760	0.7916	0.0634
5	0.522	0.5272	0.0518
6	0.774	0.7937	0.0460
7	0.608	0.6516	0.0601
8	0.707	0.7166	0.0363
9	0.791	0.7923	0.0723
10	0.786	0.7843	0.0740
11	0.716	0.7182	0.0515
12	0.788	0.7738	0.0596
13	0.643	0.6466	0.0537
14	0.687	0.7079	0.0712
15	0.784	0.7875	0.0607
16	0.929	0.9431	0.0618

For the simultaneous CHNS determination, K factor was used as calibration method with BBOT* as standard. Standard and samples weighted between 2-3 mg.

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

Table 8 shows the CHNS results with crude oil samples. Table 9 shows the NCS repeatability of lubricants while Table 10 shows the repeatability of sulfur determination in oils. No memory effect was observed changing the sample nature, meaning the combustion of all samples followed by the quantitative determination of the elements.

Table 8. CHNS repeatability of crude oils.

Sample	N%	RSD%	C%	RSD%	H%	RSD%	S%	RSD%
Crude Oil 1	0.202	4.312	85.270	0.126	13.395	0.316	0.289	1.724
	0.222		85.370		13.308		0.290	
	0.200		85.359		13.319		0.283	
	0.212		85.102		13.289		0.291	
	0.214		85.283		13.355		0.297	
Crude Oil 2	0.270	2.071	85.425	0.212	11.919	0.446	2.028	0.810
	0.282		85.770		11.925		2.018	
	0.271		85.547		11.894		2.015	
	0.279		85.377		11.794		1.985	
	0.270		85.752		11.870		2.018	

Table 9. NCS repeatability of lubricants.

Sample	N%	RSD%	C%	RSD%	S%	RSD%
1	1.9692	0.4065	83.5652	0.1459	0.0334	3.9498
	1.9533		83.3596		0.0360	
	1.9623		83.3495		0.0340	
2	0.0848	1.6698	84.3010	0.3616	0.2373	1.3201
	0.0825		84.7778		0.2316	
	0.0823		84.8720		0.2324	
3	0.1205	1.0425	84.5046	0.0746	0.6492	0.4716
	0.1185		84.5403		0.6477	
	0.1208		84.4177		0.6536	

Table 10. Sulfur repeatability of oils.

Sample	S%	RSD%
Model Oil	11.461	0.540
	11.514	
	11.520	
	11.632	
	11.531	
Crude Oil 1	8.534	0.796
	8.573	
	8.499	
	8.645	
	8.471	
Crude Oil 2	4.146	0.677
	4.165	
	4.091	
	4.125	
	4.144	
Residual Oil A	2.061	0.751
	2.086	
	2.095	
	2.093	
Residual Oil B	2.937	0.918
	2.960	
	2.904	
	2.962	
Residual Oil C	4.534	0.249
	4.541	
	4.534	
	4.558	

Conclusion

The Thermo Scientific FlashSmart N Lubricant Analyzer proved to be a valuable solution for the analysis of nitrogen in lubricants in terms of accuracy, reproducibility, and sensitivity of results. Its automation, high speed of analysis and the elimination of sample preparation process allow efficient analysis and help reduce overall operational costs.

The nitrogen analysis for petroleum products and lubricants, using the Dumas method, is described in ASTM D5291. The ASTM standard has been established for samples with percentage of nitrogen from <0.1 to 2% (mass). In this demonstration the range of applicability of the technique has been extended to samples with nitrogen in trace content, up to 0.01 N% as indicated in the Test Method D.

The FlashSmart Elemental Analyzer performs the analysis without matrix effect when changing the sample and nitrogen content. With a simple upgrade the FlashSmart EA can determine CHNS, NCS or sulfur simultaneously.

Find out more at thermofisher.com/OEA

©2017 Thermo Fisher Scientific Inc. All rights reserved. ASTM is a trademark of American Society for Testing and Materials. ISO is a trademark of the International Standards Organization. All other trademarks are the property of Thermo Fisher Scientific. 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 manner 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.

AN42216-EN 0417

ThermoFisher
SCIENTIFIC