

Nitrogen Determination in Lubricants by Flash Combustion using Argon as Carrier Gas

Liliana Krotz, Francesco Leone, and Guido Giuzzi
Thermo Fisher Scientific, Milano, Italy

Overview

Purpose: Demonstrate nitrogen determination in lubricants by Organic Elemental Analysis

Methods: Lubricant samples were analyzed using an elemental analyzer with automated autosampler using helium and argon as carrier gas.

Results: Nitrogen data from different lubricant samples are presented to demonstrate the performance of the Elementar Analyzer.

Introduction

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

The method for Nitrogen analysis in lubricants, based on combustion, is described in ASTM D5291. The method covers the instrumental determination of carbon, hydrogen and nitrogen in laboratory samples of petroleum products and lubricants. 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 increasing, it is very important to have a simple and automated technique which allows fast analysis with excellent reproducibility.

The Thermo Scientific™ FLASH™ 2000 Elemental Analyzer (Figure 1), based on the dynamic combustion of the material (Dumas method), requires no sample digestion or toxic chemicals, while providing important advantages in terms of time, automation and quantitative determination of nitrogen in a large range of concentration. However, as there is a possible worldwide shortage and an increase in costs for helium, the FLASH 2000 Analyzer can work with an alternative gas, argon which is readily available.



FIGURE 1. The Thermo Scientific FLASH 2000 Nitrogen Lubricant Analyzer.

Methods

The FLASH 2000 Organic Elemental Analyzer operates according to the dynamic flash combustion of the sample. Samples are weighed into tin capsules and introduced into the combustion reactor via the Thermo Scientific™ MAS 200R Autosampler together with oxygen.

After combustion, the produced gases are carried by a argon flow to a second reactor filled with copper, then swept through CO₂ and H₂O traps, a GC column and finally detected by a thermal conductivity detector. (Figure 2).

A report is automatically generated by the Thermo Scientific™ Eager Xperience Data Handling Software and displayed at the end of the analysis. The dedicated software allows the automatic calculation of the empirical formula.

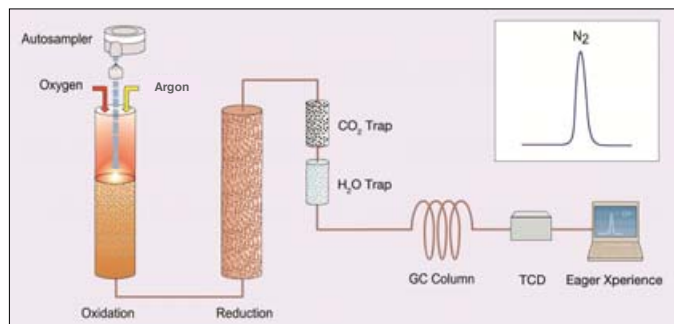


FIGURE 2. Nitrogen/Lubricant configuration.

Table 1 shows the analytical conditions using helium and argon as carrier gas.

| Parameter | Helium Gas | Argon Gas |
|--------------------------------|------------|------------|
| Combustion Reactor Temperature | 950°C | 950°C |
| Reduction Reactor Temperature | 840°C | 840°C |
| Oven Temperature | 50°C | 50°C |
| Carrier Gas Flow | 140 ml/min | 60 ml/min |
| Reference Gas Flow | 100 ml/min | 60 ml/min |
| Sample Delay | 10 sec | 12 sec |
| Oxygen Flow | 300 ml/min | 300 ml/min |
| Oxygen Injection Time | 8 sec | 20 sec |

TABLE 1. Analytical Conditions.

Note: The Eager Xperience Data Handling Software provides a new option AGO (Argon Gas Option) which allows the user to manage the flow of argon gas during the run.

Results

To evaluate the repeatability of the system using helium and argon as carrier gases, the Thermo Scientific Lubricant Reference Material (1.12 N%, accepted range 1.02-1.22 N%) was analyzed several times. Using helium as carrier gas, the instrument calibration was performed with about 4-5 mg atropine (4.84 N%) with K factor as calibration method, the Lubricant Reference Material analyzed as unknown was weighed between 8 and 10 mg. While using argon as carrier gas, the calibration was performed with 19-21 mg of atropine with K factor as calibration method, the Lubricant Reference Material analyzed as unknown was weighed between 30- 35 mg. Table 2 shows the Nitrogen data obtained with both gases.

| Gas | Helium Carrier Gas | Argon Carrier Gas |
|------------|--------------------|-------------------|
| N % | 1.054 | 1.082 |
| | 1.057 | 1.089 |
| | 1.061 | 1.091 |
| | 1.054 | 1.069 |
| | 1.045 | 1.066 |
| | 1.052 | 1.087 |
| | 1.054 | 1.093 |
| | 1.066 | 1.071 |
| | 1.058 | 1.093 |
| | 1.071 | 1.091 |
| Average N% | 1.057 | 1.083 |

TABLE 2. Nitrogen repeatability of Thermo Scientific Lubricant Reference Material with helium and argon gases.

The performance of the FLASH 2000 Analyzer was evaluated through the analysis of lubricant samples coming from an ASTM International Interlaboratory Program. The participating laboratories receive different lubricants samples by year and they are requested to analyze the samples according to their own procedures. The results are collected and processed by ASTM and a complete report is sent to each participants.

Table 3 shows the comparison of the calculated statistical results obtained by ASTM while Table 4 shows the FLASH 2000 data using helium and argon as carrier gases. All data obtained are inside in the range indicated in the ASTM reports. The samples were analyzed as received. The Sample Code include the name ALA follow by the year and the month in which the sample was analyzed and evaluated by ASTM.

Using helium as carrier gas, the instrument calibration was performed with about 4-5 mg atropine (4.84 N%) with K factor as calibration method, the ASTM Lubricants were weighed between 8 and 10 mg.

While using argon as carrier gas, the calibration was performed with atropine (13-14 mg, 4.84 N%), ALA 0902 (9-35 mg, 0.561 N%) and Thermo Scientific Lubricant Reference Material (36-37 mg) with Linear Fit as calibration method; the ASTM Lubricants were weighed between 30-35 mg.

Table 5 shows the nitrogen data of other lubricant samples obtained using helium and argon as carrier gas.

| Sample Name –Year-Month | Number of laboratories participants | ASTM Robust Mean N% | ASTM Range Accepted N% |
|-------------------------|-------------------------------------|---------------------|------------------------|
| ALA 0702 | 26 | 0.708 | 0.669 – 0.769 |
| ALA 0706 | 31 | 0.788 | 0.748 – 0.848 |
| ALA 0802 | 25 | 0.701 | 0.650 – 0.747 |
| ALA 0902 | 31 | 0.561 | 0.500 – 0.606 |
| ALA 1102 | 32 | 0.931 | 0.854 – 0.974 |

TABLE 3. ASTM information.

| Sample | Helium Carrier Gas | | | Argon Carrier Gas | | |
|----------|--------------------|------------|-------|-------------------|------------|-------|
| | N% | Average N% | RSD % | N% | Average N% | RSD % |
| ALA 0702 | 0.729 | 0.726 | 0.59 | 0.710 | 0.719 | 1.20 |
| | 0.727 | | | 0.721 | | |
| | 0.721 | | | 0.727 | | |
| ALA 0706 | 0.785 | 0.781 | 0.41 | 0.784 | 0.781 | 0.63 |
| | 0.780 | | | 0.783 | | |
| | 0.779 | | | 0.775 | | |
| ALA 0802 | 0.716 | 0.723 | 0.91 | 0.718 | 0.712 | 1.01 |
| | 0.729 | | | 0.708 | | |
| | 0.724 | | | 0.714 | | |
| ALA 0902 | 0.565 | 0.564 | 0.14 | 0.574 | 0.578 | 0.61 |
| | 0.563 | | | 0.571 | | |
| | 0.564 | | | 0.568 | | |
| ALA 1102 | 0.941 | 0.939 | 0.49 | 0.947 | 0.946 | 0.54 |
| | 0.933 | | | 0.950 | | |
| | 0.941 | | | 0.940 | | |

TABLE 4. Nitrogen reproducibility of ASTM lubricants with helium and argon gases.

| Sample Number | Helium Carrier Gas | | | Argon Carrier Gas | | |
|---------------|--------------------|------------|-------|-------------------|------------|-------|
| | N% | Average N% | RSD % | N% | Average N% | RSD % |
| 1 | 0.436 | 0.437 | 0.52 | 0.454 | 0.455 | 0.34 |
| | 0.440 | | | 0.455 | | |
| | 0.436 | | | 0.457 | | |
| 2 | 0.470 | 0.474 | 0.95 | 0.495 | 0.497 | 1.74 |
| | 0.479 | | | 0.489 | | |
| | 0.473 | | | 0.506 | | |
| 3 | 0.520 | 0.521 | 0.43 | 0.529 | 0.526 | 0.67 |
| | 0.524 | | | 0.522 | | |
| | 0.520 | | | 0.526 | | |
| 4 | 0.186 | 0.188 | 0.92 | 0.191 | 0.194 | 1.19 |
| | 0.189 | | | 0.195 | | |
| | 0.189 | | | 0.195 | | |
| 5 | 0.147 | 0.146 | 1.19 | 0.148 | 0.152 | 2.18 |
| | 0.147 | | | 0.148 | | |
| | 0.144 | | | 0.153 | | |
| 6 | 0.640 | 0.640 | 0.80 | 0.645 | 0.65 | 1.05 |
| | 0.635 | | | 0.648 | | |
| | 0.645 | | | 0.658 | | |

TABLE 5. Nitrogen reproducibility of lubricants with helium and argon gases.

The stability of the FLASH 2000 Analyzer using argon as carrier gas was verified though the analysis of 19-21 mg of atropine standard (4.84 N%, accepted range 4.79-4.91 N%) and 30-35 mg of Thermo Scientific Lubricant Reference Material (accepted range 1.184-1.224 N%) analyzed as unknown for 12 days. Calibration was performed with 19-21 mg of atropine using K factor as calibration method. During these 12 days maintenance was performed as follows:

- on day 3: the argon cylinder was replaced.
- on day 8: oxygen cylinder was replaced
- on day 11: the crucible was cleaned and the CO₂ and H₂O traps filling were replaced

Every evening and every weekend the instrument was set in Stand-By mode to reduce the consumption of argon gas. Table 6 shows the data obtained which falls within the accepted range of the technical specification of the analyzer, meaning excellent stability of the FLASH 2000 Elemental Analyzer.

| Run Day | February Date | Sample | N% | Av. N% | RSD % | Sample | N% | Av. N% | RSD % |
|---------|---------------|----------|-------|--------|-------|-----------|-------|--------|-------|
| 1 | 10 | Atropine | 4.868 | 4.87 | 0.560 | Lubricant | 1.178 | 1.186 | 0.999 |
| 2 | 11 | | 4.894 | | | | 1.177 | | |
| 3 | 12 | | 4.888 | | | | 1.178 | | |
| 4 | 13 | | 4.885 | | | | 1.188 | | |
| 5 | 14 | | 4.895 | | | | 1.182 | | |
| 6 | 17 | | 4.886 | | | | 1.184 | | |
| 7 | 18 | | 4.900 | | | | 1.198 | | |
| 8 | 19 | | 4.836 | | | | 1.187 | | |
| 9 | 20 | | 4.893 | | | | 1.204 | | |
| 10 | 21 | | 4.816 | | | | 1.188 | | |
| 11 | 24 | | 4.895 | | | | 1.182 | | |
| 12 | 25 | | 4.789 | | | | 1.195 | | |

TABLE 6. Day by day reproducibility in N determination using argon as carrier gas.

Conclusion

The FLASH 2000 N Lubricant Analyzer is an excellent solution for the analysis of nitrogen in lubricant in terms of:

- accuracy,
- reproducibility,
- sensitivity,
- automation,
- speed of analysis,
- cost per analysis.

All data showed were obtained with an acceptable repeatability and no matrix effect was observed when changing the sample. Good repeatability was obtained with the FLASH 2000 Analyzer using argon as carrier gas. The data are comparable with those obtained when using helium.

By using the FLASH 2000 Analyzer it is possible to characterize the different lubricant samples according to the nitrogen content coming from the additive. Knowledge of its concentration can be used to predict performance.



www.thermoscientific.com

©2016 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. AOAC is a trademark of The Association of Official Analytical Chemists. AOCS is a trademark of The American Oil Chemists' Society. AACC is a trademark of The American Association of Cereal Chemists. ASBC is a trademark of The American Society of Brewing Chemists. IFFO is a trademark of The International Fishmeal and Fish Oil 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.

COMPANY WITH QUALITY MANAGEMENT
SYSTEM CERTIFIED BY DNV
= ISO 9001:2008 =

Thermo Fisher Scientific S.p.A. Milan, Italy
is ISO 9001:2008 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 2730 3006
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
PN42247-EN 0216S

Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand