



# Multi-Element Analysis of Fuel and Lubricating Oils by Simultaneous ICP-OES

## Application Note

Inductively Coupled Plasma-Optical Emission Spectrometers

### Author

Andrew Ryan

### Introduction

The determination of metal levels in both fuel oil and lubricating oils is of prime importance to the industrial machinery, transport and utility industries. In general, metal levels are determined to provide early warning of failure of components in physical contact with the oil. Oil analysis can also indicate whether oil in a system has been contaminated by another source, other than wear. For example increases in silicon or aluminum levels may indicate dirt contamination or increased levels of sodium, perhaps seawater contamination or antifreeze leakage [1]. Conversely, a decrease in the levels of additive metals (for example, Ca, Mg, Zn, Pb) may indicate a dilution of the oil by one oil not containing these additives. Many large fleet operators and mining companies rely on the analysis of used lubricating oil samples to extend the period between oil changes in their equipment and to warn of the beginning of mechanical failure so that preventive maintenance may be scheduled.

In this study various fuel oils and National Institute of Standards and Technology (N.I.S.T.), standard reference material (S.R.M.) 1084a Wear Metals in Lubricating Oil, were diluted in dekaline (decahydronaphthalene) and analyzed using an Agilent Vista simultaneous Inductively Coupled Plasma Atomic Emission Spectrometer (ICP-OES) with the axially viewed plasma.



**Agilent Technologies**

For the analysis of volatile organic solvents with an axially viewed ICP, the Auxiliary Gas Module 1 (AGM-1) oxygen accessory is required. The AGM-1 adds a small flow of oxygen to the plasma auxiliary gas, thus reducing molecular band emissions. This gives lower detection limits, and prevents the build up of carbon within the torch and on the cooled cone interface.

The use of a simultaneous ICP-OES instrument significantly improves the efficiency of analysis when a large number of elemental determinations are to be made. Typically, large laboratories specializing in used oil analysis test hundreds of samples each day. In this study 21 elements were determined at around 105 seconds per sample including sample introduction and rinse time. It should be noted that the measurement of more than this number of elements would not add further to the analysis time due to the simultaneous detection of the Vista.

## Instrumental

The Vista simultaneous ICP-OES with axially viewed plasma was used for the analysis. The Vista features a free running, air-cooled, 40 MHz RF generator and cooled cone interface. The Vista's optical system is based on an echelle polychromator with CCD detector [2]. The polychromator is thermostatted to 35 °C for stability and the unique CCD detector features 70,000 pixels (detectors) arranged to exactly match the 2 dimensional echellogram. The detector provides rapid readout and excellent detection limits. The instrument was controlled with an IBM computer with an Intel Pentium processor and Agilent's Vista software running under Microsoft Windows NT operating system.

In this work, a Vista with manual gas pressure regulator was used. Mass flow control of the nebulizer flow, which allows the nebulizer gas flow to be automatically adjusted, is available as an option.

### Instrument parameters

|                       |  |
|-----------------------|--|
| Power                 | 1.3 kW   |
| Plasma gas flow       | 15 L/min   |
| Auxiliary gas flow    | 0.75 L/min   |
| Nebulizer type        | Glass concentric type K  |
| Torch type            | Demountable with 1.8 mm quartz injector  |
| Spray chamber type    | Sturman-Masters  |
| Pump speed            | 15 rpm   |
| Pump tubing           | Inlet:–viton, black/black (0.76 mm id),<br>Outlet:–viton, white/white (1.02 mm id)   |
| Sample uptake rate    | 0.6 mL/min   |
| AGM – 1 setting       | 5 (130 mL/min)   |
| Sample delay          | 15 seconds   |
| Stabilization time    | 45 seconds   |
| Fast pump             | On   |
| Replicate time        | 5 seconds  |
| Replicates            | 3  |
| Background correction | Off peak background correction or fitted background correction. Fast Automated Curve Fitting Technique (FACT) was used for Na 588.995 nm |

## Sample Preparation

Approximately 2 g of used oil was accurately weighed into a 20 mL volumetric flask and made up to volume with dekalin.

One gram of N.I.S.T, S.R.M 1084a Wear Metals in Lubricating Oil was accurately weighed into a 50 mL volumetric flask and made up to volume with dekalin.

## Results

In this study seven fuel oil samples and one lubricating oil sample were analyzed. The lubricating oil was a standard reference material obtained from NIST, Gaithersburg, MD, USA and used to validate the method. Triplicate analyses were performed and the average result is shown.

Table 1.

|                         | Ag 328.068 nm   | Al 257.509 nm | B 249.678 nm | Ba 493.408 nm | Ca 317.933 nm | Cd 226.502 nm |
|-------------------------|-----------------|---------------|--------------|---------------|---------------|---------------|
| IC 1 (mg/kg)            | 0.011 ± 0.002   | 5.83 ± 0.03   | < 0.2        | < 0.003       | 5.35 ± 0.06   | < 0.02        |
| IC 2 (mg/kg)            | 0.0148 ± 0.0005 | 28.7 ± 0.1    | < 0.2        | < 0.003       | 26.5 ± 0.3    | < 0.02        |
| IC 3 (mg/kg)            | 0.04 ± 0.01     | 114.6 ± 1.3   | 0.270 ± 0.01 | < 0.003       | 109.0 ± 0.6   | < 0.02        |
| IC 5/4 (mg/kg)          | 0.04 ± 0.01     | 2.80 ± 0.03   | < 0.2        | 37.5 ± 0.3    | 222 ± 1       | 0.95±0.02     |
| IC 6 (mg/kg)            | 0.09 ± 0.01     | < 1           | 1.34 ± 0.1   | < 0.003       | 640 ± 4       | < 0.02        |
| IC 10 (mg/kg)           | 0.031 ± 0.002   | < 1           | 0.53 ± 0.08  | 16.3 ± 0.2    | 3210 ± 7      | 0.117 ± 0.003 |
| IC 15 (mg/kg)           | < 0.01          | < 1           | < 0.2        | < 0.003       | 0.974 ± 0.02  | < 0.02        |
| N.I.S.T S.R.M 1084a     | 100.3 ± 1.2     | 100.4 ± 0.4   | —            | —             | —             | —             |
| Certified value (mg/kg) | 101.4 ± 1.5     | (104)         | —            | —             | —             | —             |

< Values expressed as 10 times the standard deviation of background emission

Values in parenthesis are not certified but provided for information only.

— Value not certified.

Table 2:

|                         | <b>Cr 267.716 nm</b> | <b>Cu 324.754 nm</b> | <b>Fe 259.940 nm</b> | <b>Mg 279.553 nm</b> | <b>Mo 284.824 nm</b> | <b>Mn 257.610 nm</b> |
|-------------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|
| IC 1 (mg/kg)            | 6.21 ± 0.05          | 6.41 ± 0.04          | 5.95 ± 0.05          | 0.067 ± 0.007        | < 0.1                | 5.69 ± 0.03          |
| IC 2 (mg/kg)            | 29.4 ± 0.3           | 29.3 ± 0.1           | 28.5 ± 0.4           | 0.16 ± 0.03          | < 0.1                | 28.2 ± 0.3           |
| IC 3 (mg/kg)            | 113.9 ± 0.7          | 115.1 ± 0.5          | 112.8 ± 0.8          | 0.19 ± 0.04          | < 0.1                | 107.4 ± 0.5          |
| IC 5/4 (mg/kg)          | 3.05 ± 0.05          | 22.8 ± 0.3           | 50.9 ± 0.7           | 2.04 ± 0.05          | < 0.1                | 0.958 ± 0.009        |
| IC 6 (mg/kg)            | < 0.02               | 0.16 ± 0.03          | 1.11 ± 0.02          | 69.9 ± 0.9           | 21.4 ± 0.4           | 0.161 ± 0.002        |
| IC 10 (mg/kg)           | 0.131 ± 0.002        | 1.32 ± 0.02          | 6.5 ± 0.1            | 10.5 ± 0.2           | 0.46 ± 0.01          | 0.233 ± 0.003        |
| IC 15 (mg/kg)           | < 0.02               | 0.17 ± 0.01          | 0.44 ± 0.05          | 0.310 ± 0.005        | < 0.1                | 0.006 ± 0.002        |
| N.I.S.T S.R.M 1084a     | 96.8 ± 1.3           | 97.6 ± 1.1           | 98.8 ± 2.4           | 97.9 ± 1.5           | 98.8 ± 1.5           | —                    |
| Certified value (mg/kg) | 98.3 ± 0.8           | 100.0 ± 1.9          | 98.9 ± 1.4           | 99.5 ± 1.7           | 100.3 ± 1.4          | —                    |

< Values expressed as 10 times the standard deviation of background emission

— Value not certified

Table 3.

|                         | <b>Na 588.995 nm</b> | <b>Ni 216.555 nm</b> | <b>P 213.618 nm</b> | <b>Pb 220.353 nm</b> | <b>Si 251.611 nm</b> | <b>Ti 334.941 nm</b> |
|-------------------------|----------------------|----------------------|---------------------|----------------------|----------------------|----------------------|
| IC 1 (mg/kg)            | 0.436 ± 0.006        | 5.80 ± 0.03          | < 1                 | 0.98 ± 0.02          | 6.01 ± 0.08          | 5.45 ± 0.03          |
| IC 2 (mg/kg)            | 2.24 ± 0.05          | 28.3 ± 0.2           | < 1                 | 2.5 ± 0.2            | 29.4 ± 0.2           | 29.1 ± 0.2           |
| IC 3 (mg/kg)            | 2.97 ± 0.06          | 108 ± 1              | < 1                 | 48.1 ± 0.7           | 112.7 ± 0.7          | 112.7 ± 0.7          |
| IC 5/4 (mg/kg)          | 6.7 ± 0.1            | < 0.1                | 119.5 ± 0.7         | 19.1 ± 0.2           | 7.57 ± 0.09          | 0.205 ± 0.008        |
| IC 6 (mg/kg)            | 0.62 ± 0.01          | < 0.1                | 248 ± 3             | < 0.3                | 1.58 ± 0.03          | 0.011 ± 0.004        |
| IC 10 (mg/kg)           | 37.1 ± 0.8           | 3.08 ± 0.03          | 237 ± 5             | 0.91 ± 0.02          | 2.78 ± 0.03          | 0.064 ± 0.004        |
| IC 15 (mg/kg)           | 4.82 ± 0.05          | 4.58 ± 0.08          | < 1                 | < 0.3                | < 0.5                | < 0.01               |
| N.I.S.T S.R.M 1084a     | —                    | 100.2 ± 1.2          | —                   | 100.0 ± 0.9          | 96.1 ± 0.9           | 98.9 ± 1.3           |
| Certified value (mg/kg) | —                    | 99.7 ± 1.6           | —                   | 101.1 ± 1.3          | (103)                | 100.4 ± 3.8          |

< Values expressed as 10 times the standard deviation of background emission

Values in parenthesis are not certified but provided for information only

— Value not certified

Table 4:

|                         | <b>V 311.070 nm</b> | <b>Zn 213.857 nm</b> |
|-------------------------|---------------------|----------------------|
| IC 1 (mg/kg)            | 5.68 ± 0.03         | 5.76 ± 0.05          |
| IC 2 (mg/kg)            | 28.3 ± 0.3          | 28.2 ± 0.3           |
| IC 3 (mg/kg)            | 112 ± 2             | 108.1 ± 0.7          |
| IC 5/4 (mg/kg)          | < 0.1               | 32.4 ± 0.5           |
| IC 6 (mg/kg)            | < 0.1               | 250 ± 2              |
| IC 10 (mg/kg)           | 5.91 ± 0.06         | 238 ± 2              |
| IC 15 (mg/kg)           | 15.3 ± 0.2          | 0.356 ± 0.007        |
| N.I.S.T S.R.M 1084a     | 96.4 ± 1.3          | —                    |
| Certified value (mg/kg) | 95.9 ± 9.4          | —                    |

< Values expressed as 10 times the standard deviation of background emission

— Value not certified

## Discussion

For the NIST SRM 1084a sample, measured values are in good agreement with the certified values demonstrating the suitability of the axial Vista ICP-OES for the analysis of used oils.

Background structure, the result of using an organic solvent, posed some difficulties in the measurement of sodium at 588.995 nm. Using Vista's Fast Automated Curve-fitting Technique (FACT) improved the background measurement at this line and improved the levels of detection. FACT uses Gaussian curve modeling techniques to accurately characterize and correct for neighboring spectral background [3]. FACT has been proven to provide accurate background correction even when peaks are as close as 0.2 pixels and is ideal for correcting the background observed from organic solvents.

## Acknowledgment

The author wishes to thank Stewart Carter for his assistance in preparing this manuscript.

## References

T. McKenzie, "Atomic Absorption Spectrophotometry for the analysis of Wear Metals in Oil Samples", Varian Australia Pty. Ltd., Mulgrave, Victoria 3170, Australia, AA At Work No. 10, January 1981.

M. B. Knowles T. Nham, S. J. Carter, "A Charged-Coupled Device Detector for Atomic Emission Spectroscopy", American Laboratory, September, 1998.

C. Webb, A. T Zander, P. V. Wilson and G. Perlis, "A Fast Automated Spectral Curve Fitting Technique for ICP-AES". Spectroscopy, May 1999, 14(5), 58-63.

## For More Information

For more information on our products and services, visit our Web site at [www.agilent.com/chem](http://www.agilent.com/chem)

[www.agilent.com/chem](http://www.agilent.com/chem)

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc., 1999  
Printed in the USA  
November 1, 2010  
ICPES-27



**Agilent Technologies**