



# Analysis of coal and coal ash using the Thermo Scientific iCAP 7400 Duo ICP-OES

## Authors

James Hannan,  
Application Specialist,  
Thermo Fisher Scientific,  
Hemel Hempstead, UK

## Keywords

Coal, Coal ash, Flue ash, Fly ash

## Goal

The goal of this application note is to discuss and demonstrate the elemental analysis of coal and coal ash samples using ICP-OES following microwave acid digestion.

## Introduction

Despite the recent improvements in, and increasing use of renewable forms of energy, approximately 40% of the world's electricity is still generated by coal fired power stations. This figure is significantly higher in some countries, for example, approximately 70% in India and over 90% in South Africa.

During coal combustion, large amounts of ash are created along with carbon dioxide and other gases. The fine particle ash that rises up with the flue gases is known as fly or flue ash while the heavier ash that does not rise is called bottom ash; collectively these are known as coal ash. Traditionally, the fly ash was released into the atmosphere, however, due to its potentially toxic effects it is now generally collected from the flue towers using electrostatic precipitators or other particle filtration equipment. It can then be disposed of or recycled into Portland cement.

The chemical makeup of fly and bottom ash varies significantly and is dependent on the source and composition of the coal being burned. This can include a wide variety of toxic substances from trace amounts to percent levels. In order to protect the environment or the quality and safety of any products it is added to, the composition of the ash product needs to be accurately analyzed before it can be recycled or disposed of.

## Instrumentation and method parameters

For this analysis the Thermo Scientific™ iCAP™ 7400 ICP-OES Duo was used, it has full control over the RF power and nebulizer gas flow rate, while the Duo view plasma allows for axial and radial viewing modes to be used for the greatest detection limits and maximum upper analysis ranges. The measurement parameters and sample introduction setup can be seen in Table 1.

**Table 1. Method parameters and sample introduction setup used.**

Parameter	Setting	
Pump Tubing (Standard Pump)	Sample Tygon® orange/white Drain Tygon® white/white Internal Standard Tygon® orange/blue	
Pump Speed	45 rpm	
Center Tube	2 mm	
Spray Chamber	Glass cyclonic	
Nebulizer	Burgener Mira Mist®	
Nebulizer Gas Flow	0.5 L·min <sup>-1</sup>	
RF Power	1350 W	
Exposure Time	<b>Axial</b>	<b>Radial</b>
	UV 15 s, Vis 5 s	UV 15 s, Vis 5 s

A 10 mg·L<sup>-1</sup> yttrium internal standard was introduced using the online internal standard kit. This allows the method to compensate for variations in sample matrices along with any changes to the sample introduction. The internal standard wavelengths used can be seen in Table 2. The Thermo Scientific™ Qtegra™ Intelligent Scientific Data Solution™ (ISDS) Software line switching feature was used for the majority of elements, allowing two or more wavelengths to be calibrated at different concentration ranges for the same element, extending the working range of the instrument.

**Table 2. Internal standard wavelengths of yttrium used.**

	Axial	Radial
UV	224.306 nm	224.306 nm
Vis	371.030 nm	371.030 nm

A range of single element solutions were prepared and analyzed in order to observe any spectral interferences that are present. The contributions of any interference present were measured, and Inter-Element Corrections (IECs) were applied. This allows the Qtegra ISDS Software to automatically compensate for these interferences on a sample by sample basis.

## Sample preparation

Two certified reference materials (CRMs) (one bituminous coal and one fly ash, both from the National Institute of Standards and Technology, USA (NIST)) were prepared in triplicate using the following method. The samples were digested by adding 6 ml of 20% hydrochloric acid (HCl) and 2 ml of 20% nitric acid (HNO<sub>3</sub>) to approximately 0.1 g of solid sample. This was then heated to 220 °C for 35 minutes using a microwave digestion system. Once the samples had cooled they were made up to 25 ml with ultrapure de-ionized water.

Calibration standards were created from 1000 mg·L<sup>-1</sup> and 10000 mg·L<sup>-1</sup> single element solutions (Fisher Scientific Chemicals, Loughborough, UK) and were matrix matched to the sample digests, i.e. 4.8% HCl and 1.6% HNO<sub>3</sub>.

## Results

The analytical results of the two CRM samples compared with the certified and reference values are shown in Table 3. An empty sample tube was subjected to the same sample digestion method as the samples, this blank was then analyzed with 10 replicates. The method detection limits were calculated by multiplying the standard deviation of these 10 replicates by a factor of 3, this was performed in triplicate and an average taken (Table 3). Table 3 shows that the analytical results are within ±5% of the certified values and within ±15% of the reference values, demonstrating the accuracy of this method for the analysis of coal and coal ash samples.

It should be noted that silicon typically constitutes a significant proportion of the composition of these sample types, particularly fly ash. Therefore if the analysis of silicon is required, then a small quantity (0.5 ml) of hydrofluoric acid (HF) will need to be added into the digestion procedure, otherwise low recovery of silicon will occur.

Table 3. Analytical results for NIST certified reference materials. Wavelengths for each element are listed in order of sensitivity.

Element	Wavelengths (nm)	Method detection limit ( $\mu\text{g}\cdot\text{L}^{-1}$ )	Units	NIST 1632d (Coal)	Certified/reference values	Recovery (%)	NIST 2690 (Fly ash)	Certified/reference values	Recovery (%)
Al	167.079 (A) 396.152 (R) 237.312 (R)	0.01	%	0.791	0.912*	86.7	12.81	12.35	103.7
As	189.042 (A)	9.4	$\text{mg}\cdot\text{kg}^{-1}$	6.026	6.1*	98.8	27.18	26*	104.5
B	249.773 (A) 249.678 (R)	6.9	$\text{mg}\cdot\text{kg}^{-1}$	58.93	62*	95.0	282.5	NA	-
Ca	422.673 (A) 184.006 (R)	0.001	%	0.133	0.144*	92.3	5.526	5.71	96.8
Cd	228.082 (A) 214.438 (R)	0.42	$\text{mg}\cdot\text{kg}^{-1}$	<0.1	0.08*	-	0.655	0.7*	93.6
Cr	283.563 (A) 267.716 (R)	4.7	$\text{mg}\cdot\text{kg}^{-1}$	12.70	13.7*	92.7	69.49	67*	103.7
Cu	324.754 (A) 224.70 (R)	3.4	$\text{mg}\cdot\text{kg}^{-1}$	5.736	5.83	98.4	64.84	NA	-
Fe	259.940 (A) 239.562 (R)	0.003	%	0.779	0.749	104.1	3.442	3.57	96.4
Hg	184.950 (A) 194.227 (A)	0.42	$\text{mg}\cdot\text{kg}^{-1}$	<0.1	0.0928	-	<0.1	<0.003*	-
K	766.490 (A+R)	0.01	%	0.106	0.1094	96.8	1.063	1.04	102.2
Mg	279.553 (A) 285.213 (R)	0.00002	%	0.034	0.039*	89.6	1.561	1.53	102.0
Mn	257.610 (A) 279.482 (R)	2.6	$\text{mg}\cdot\text{kg}^{-1}$	13.97	13.1*	106.7	313.00	300*	104.3
Na	588.995 (A) 589.592 (A)	6.4	$\text{mg}\cdot\text{kg}^{-1}$	286.5	296.9	96.5	2312.3	2400	96.3
Ni	231.604 (A) 341.476 (R)	0.23	$\text{mg}\cdot\text{kg}^{-1}$	11.99	NA	-	47.78	46*	103.9
Pb	220.353 (A) 216.999 (R)	0.96	$\text{mg}\cdot\text{kg}^{-1}$	3.750	3.845	97.5	35.75	39*	91.7
Se	196.090 (A)	2.0	$\text{mg}\cdot\text{kg}^{-1}$	1.418	1.29*	110.0	0.732	0.8*	91.5
Ti	334.941 (A) 338.376 (R)	3.8	$\text{mg}\cdot\text{kg}^{-1}$	487.7	477	102.3	4946	5200	95.1
V	282.402 (A)	0.56	$\text{mg}\cdot\text{kg}^{-1}$	22.77	23.74	95.9	148.3	NA	-
Zn	202.548 (A)	0.11	$\text{mg}\cdot\text{kg}^{-1}$	14.21	12.9*	110.2	117.9	120*	98.2

\*= non certified reference value, (A) = Axial view, (R) = Radial view

## Conclusion

The analytical results in Table 3 demonstrate that the iCAP 7400 ICP-OES Duo is capable of performing the analysis of coal and fly ash samples. The duo view combined with the features of the Qtegra ISDS Software such as line switching allow for low detection limits to be achieved whilst maintaining a large working range in a single analytical method.

Features and tools such as internal standardization and IECs combined with the adaptable, high efficiency RF generator allow for heavy matrix samples to be accurately analyzed without the need for extensive dilutions.

Find out more at [thermofisher.com/ICP-OES](https://thermofisher.com/ICP-OES)

**For Research Use Only. Not for use in diagnostic procedures.** ©2017 Thermo Fisher Scientific Inc. All rights reserved. Burgener Mira Mist is a trademark of John Burgener of Burgener Research Inc. Tygon is a trademark of Saint-Gobain Corporation. 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 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. **AN43186-EN 0717**

**ThermoFisher**  
SCIENTIFIC