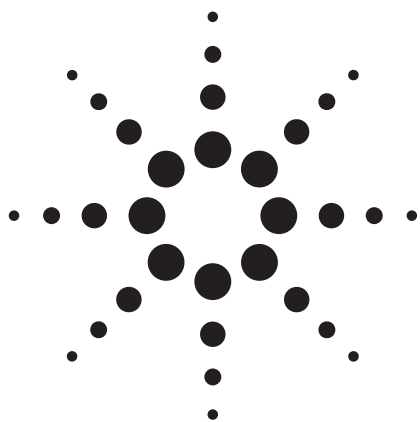




The Analysis of a Zinc-Based Alloy

Application Note

Atomic Absorption

Authors

A. W. Copley and S. Skujins

Introduction

Atomic absorption spectrophotometry has found wide application in the routine determination of both the major, as well as most of the minor, constituents of non-ferrous metal alloys.

The suitability of the technique for the analysis of non-ferrous alloys is reflected by the large number of publications [1-17].

A description is given here of a method employed at Agilent Technologies for the analysis of zinc, aluminium, copper, iron, lead, and tin in a zinc-based alloy.



Agilent Technologies

Experimental

Alloy Sample Solutions

Approximately 1 g of the alloy sample is weighed out accurately into an acid-cleaned, dry 200 mL conical flask and the alloy is reacted with 30 mL of 30% v/v hydrochloric acid.

After the reaction has almost stopped the mixture is gently heated and 5 mL of concentrated nitric acid are slowly added.

The clear sample solution is cooled and transferred to a standard volumetric flask and diluted to exactly 100 mL with distilled water.

Aliquots of this sample solution may be further diluted, if necessary, in order to bring each metal into the appropriate concentration range for measurement by atomic absorption spectrophotometry.

For example,

A 1% solution of an alloy containing approximately

will contain approximately

65% Zn	6,500 ppm Zn
30% Al	3,000 ppm Al
5% Cu	500 ppm Cu
1% Fe, Pb, Sn	100 ppm Fe, Pb, Sn

Standard Solutions

Composite standard solutions are prepared which contain all the elements that are present in the sample. The concentration range of the standards is made to cover the expected concentration of the metals in the sample solutions.

Notes on Individual Determinations

Zinc

The use of the most sensitive analytical line for zinc (0.01 ppm Zn/1% Absorption) at 2138.6 Å would require a 5,000-fold dilution of the sample solution. Such a large dilution could introduce serious errors and lead to poor accuracy in the results.

It is preferable to carry out the atomic absorption measurements at the much less sensitive (8,400X) resonance line at 3075.9 Å, using an air-C₂H₂ flame.

The recommended concentration range for the Zn standards is from 2,000-10,000 ppm.

Aluminium

For the determination of aluminium the sample is diluted approximately 30–50-fold and the atomic absorption measurements are carried out at the most sensitive Al resonance line, the 3092.7 Å, 3092.8 Å doublet, using a N₂O-C₂H₂ flame.

Aluminium standards are prepared in the range 20–150 ppm Al.

Copper

The determination of copper in an air-C₂H₂ flame can be carried out at a number of Cu resonance lines.

The use of the most sensitive line at 3247.5 Å (0.04 ppm Cu/1% Absorption) would require an approximately 100-fold dilution of the sample solution, with standards in the range 2–8 ppm Cu. If burner rotation is employed to reduce the sensitivity then such a large dilution may not be necessary.

Use may also be made of the line at 3275 Å (sensitivity 0.14 ppm Cu/1% Absorption). In this case a 40–50-fold dilution of the sample solution has to be carried out. Standards should be in the range 5–20 ppm Cu.

Other resonance lines of even lower sensitivity at 2492 Å and at 2178.9 Å may also be employed.

Iron

For an alloy containing 0.10% Fe no further dilution of the sample solution is required.

The atomic absorption measurements are made at the most sensitive line (0.08 ppm Fe/1% Absorption) at 2483.3 Å using an air-C₂H₂ flame.

The recommended range of standards is from 2–10 ppm Fe.

Lead

An air-C₂H₂ flame is used, together with the most sensitive Pb line (0.16 ppm Pb/1% Absorption) at 2170.0 Å. The use of scale expansion may be required.

The recommended range of standards is from 1–20 ppm Pb.

For samples containing less than 20 ppm Pb it is recommended that larger weights of sample be employed, for example, 2–3 g alloy/100 mL, in order to improve the accuracy of the measurements.

Tin

The best sensitivity may be obtained at the 2246.1 Å Sn resonance line, using either an air-C₂H₂ or an air-H₂ flame (0.4 ppm Sn/1% Absorption). However, in these relatively cool flames chemical interference problems are often encountered.

In order to overcome or, at least minimize chemical interferences the much hotter, although less sensitive (2.0 ppm Sn/1% Absorption), $N_2O-C_2H_2$ flame may be employed.

For samples containing less than 0.05% Sn it is recommended that large weights of sample be used, for example. 2–3 g alloy/100 mL, together with scale expansion.

The recommended range of standards is from 5–30 ppm Sn.

Some Typical Results

Approximate composition		Agilent Technologies analysis	
Zn	65%	Zn	64.6%
Al	30%	Al	29.59%
Cu	5%	Cu	4.88%
Fe	1%	Fe	0.10%
Pb		Pb	20 ppm
Sn		Sn	0.03 ppm
			99.22%.

Conclusion

The routine analysis of such non-ferrous alloys by atomic absorption spectrophotometry is relatively simple and rapid using equipment which has scale expansion and burner rotation facilities.

References

1. J. A. F. Gidley, J. T. Jones, *Analyst*, 85, 249 (1960)
2. J. A. F. Gidley, J. T. Jones, *Analyst*, 86, 271 (1961)
3. W. T. Elwell, J. A. F. Gidley, *Anal. Chim. Acta*, 24(1), 71 (1961)
4. W. Leithe, A. Hofer, *Mikrochim. Acta*, 2, 277 (1961)
5. T. R. Andrews, P. N. R. Nicholas, *Analyst*, 87, 25 (1962)
6. G. I. Nikolaev, V. B. Aleskovskii, *Zhur. Anal. Khim.*, 18(7) 817 (1963)
7. G. I. Nikolaev, *Zhur. Anal. Khim.* 19, 63 (1964)
8. K. A. Stewart, A. H. Knight, *Proc. Soc. Anal. Chem.*, 5, 116 (1965)
9. T. R. Andrews, P. N. R. Nicholas, *Analyst*, 90, 161 (1965)
10. T. Miura, *Japan Analyst*, 14, 310 (1965)
11. R. M. Dagnall, T. S. West, P. Young, *Anal. Chem.*, 38(2), 358 (1966)
12. R. C. Calkins, *Applied Spectroscopy*, 20(3), 146 (1966)
13. W. Schmidt, *Chem. Listy*, 60(4) 540 (1966)
14. R. E. Mansell, H. W. Emmel, E. L. McLaughlin, *Applied Spectroscopy*, 20, 231 (1966)
15. J. L. J. Seco, A. G. Coedeo, *Revista de Metalurgia*, 4, 621 (1968)
16. J. L. J. Seco, A. G. Coedeo, *Revista de Metalurgia*, 1, 33 (1969)
17. J. M. Hall, C. Woodward, *Spectroscopy Letters*, 2, 113 (1969)

For More Information

For more information on our products and services, visit our Web site at www.agilent.com/chem

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., 1970
Printed in the USA
November 1, 2010
A-AA05



Agilent Technologies