

High Precision Nickel Alloy Analysis by the Thermo Scientific ELEMENT GD PLUS GD-MS

Joachim Hinrichs
Thermo Fisher Scientific, Bremen, Germany

Key Words

GD-MS, Nickel Alloys, Pulsed

Goal

To achieve superior performance of the Thermo Scientific ELEMENT GD PLUS GD-MS pulsed for the analysis of Ni alloys, compared to the bulk analysis of conductive metals with continuous DC GD-MS.

Introduction

The Thermo Scientific™ ELEMENT™ GD GD-MS is used and well accepted for trace metal determination in Ni alloys controlled by aerospace regulations. With the upgrade to a pulsed supply of the GD source, the precision achievable on the Thermo Scientific™ ELEMENT™ GD PLUS GD-MS needs to be verified. The accuracy is dominated by the calibration materials used and is equivalent to continuous DC operation.

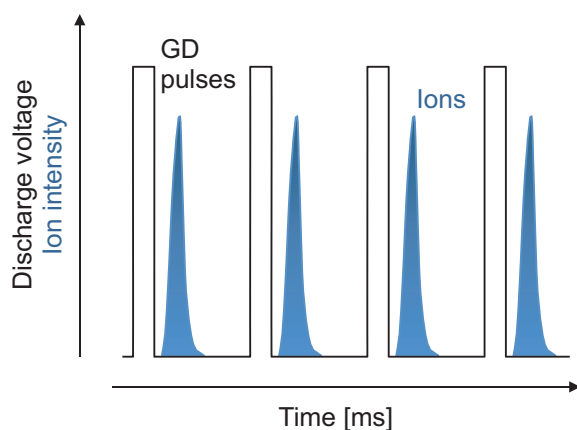


Figure 1. Schematics of pulsed discharge operation.



Method

The Certified Reference Material (CRM) BAS346A is a nickel superalloy (IN100) with a well characterized set of trace metals important for materials used in the aerospace industry. Therefore this CRM has been used exemplarily to investigate the precision achievable in routine GD-MS quality control tasks.

Repeat analyses were performed including sample preparation by wet grinding the sample on SiC paper grit 80, followed by a washing with de-ionized water. As a finish, the sample surface was soaked with iso-propanole and blown dry in a nitrogen stream.

Table 1. Instrumental parameters.

Parameter	Value
Discharge Voltage	800V
Pulse Frequency	4 kHz
Pulse Duration	40 μ s
Average Discharge Current	~15 mA
Discharge Gas	500 mL argon min ⁻¹
Matrix Intensity	~1·10 ¹⁰ cps (MR)
Preputter Time	6 min
Acquisition Time	6 min
Anode Consumables	Stainless steel

Table 2. Precision data of repeat analysis of CRM BAS346A.

Element	Isotope (Resolution)	Certified value [$\mu\text{g}\cdot\text{g}^{-1}$]	RSD of all 16 spots	RSD of 6 spots with flow tube and cap exchange	RSD of 4 spots with cap exchange	RSD of 6 spots without any part exchanged
B	11 (MR)	~200	4.2%	4.0%	1.3%	1.2%
C	12 (MR)	~1500	3.2%	4.3%	2.9%	0.8%
Mg	24 (MR)	130 \pm 5%	1.2%	1.1%	0.7%	1.2%
Ca	44 (MR)	~20	8%	11%	6%	6%
Zn	66 (MR)	28.8 \pm 5%	2.5%	2.3%	1.1%	1.1%
Ga	69 (MR)	49.6 \pm 4%	1.5%	1.8%	0.7%	0.9%
As	75 (MR)	50.4 \pm 5%	1.2%	0.6%	1.0%	1.1%
Se	82 (MR)	5.7 \pm 14%	10%	13%	8%	9%
Ag	107 (MR)	42.5 \pm 2%	2.4%	2.0%	1.6%	0.7%
Cd	111 (MR)	0.37 \pm 11%	9%	9%	9%	11%
In	115 (MR)	~20	1.6%	1.6%	1.1%	1.1%
Sn	119 (MR)	93 \pm 9%	0.9%	1.0%	1.0%	0.9%
Sb	121 (MR)	45 \pm 9%	1.2%	0.9%	1.1%	1.2%
Te	130 (MR)	9.3 \pm 9%	7.7%	10.6%	5.5%	3.8%
Tl	205 (MR)	1.9 \pm 16%	3.0%	3.4%	0.7%	3.3%
Pb	208 (MR)	22.2 \pm 5%	1.4%	1.7%	0.6%	1.1%
Bi	209 (MR)	10.3 \pm 7%	2.4%	3.0%	1.1%	2.4%

Results

- The results in Table 2 indicate excellent reproducibility for 16 repeat runs including sample preparation.
- Presputter and acquisition times are comparable to continuous (non-pulsed) operation.
- Best precisions are observed when no source parts are exchanged. This allows for extended analysis (>2 h sputter time, equivalent to approximately 10 sample runs) of similar sample types.
- The typical procedure of changing the anode cap does not significantly impact the precision.
- The exchange of the flow tube shows a visible impact, but still gives precisions similar or better than the reference values itself.
- A long-term test run on one spot without sample change showed that even after 4 hours the measured concentrations are within a 10% limit. About half of this variation originates from the deep crater eroded, which is never achieved in routine work with ~10-15 minutes sputter time per sample.
- The methodology shown here is subject to ongoing and further optimization, depending on individual lab requirements.

Conclusion

The ELEMENT GD PLUS GD-MS offers a solution for highly precise trace metal determination in nickel super alloys. Its much lower sputter rate at a similar level of intensity compared to the non-pulsed mode yields an excellent spot-to-spot as well as per-spot precision.

www.thermofisher.com/GDMS

©2016 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards 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 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.

Africa +43 1 333 50 34 0
Australia +61 3 9757 4300
Austria +43 810 282 206
Belgium +32 53 73 42 41
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 9 3291 0200
France +33 1 60 92 48 00
Germany +49 6103 408 1014
India +91 22 6742 9494
Italy +39 02 950 591

Japan +81 45 453 9100
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
Spain +34 914 845 965
Sweden +46 8 556 468 00
Switzerland +41 61 716 77 00
UK +44 1442 233555
USA +1 800 532 4752

Thermo
 SCIENTIFIC

A Thermo Fisher Scientific Brand