

Analysis of alumina powders

Authors

Joachim Hinrichs¹ and
Karol Putyera²

¹Thermo Fisher Scientific,
Bremen, Germany,

²Evans Analytical Group,
Liverpool, NY, USA

Keywords

Ceramics, GD-MS, High Purity,
Non-conductive, μ s-Pulsed

Goal

To demonstrate the capabilities of the Thermo Scientific Element GD Plus μ s-Fast-Flow Glow Discharge Mass Spectrometer for high throughput trace metal determination in high purity alumina powders with minimum sample preparation.

Introduction

Items produced from high-purity Al_2O_3 powders are found in a large variety of consumer and industrial products. With the predicted increased demand for 5N and higher purities, a fast, simple and accurate analytical technique is required to control production.

Non-conductive oxide powders in general and alumina in particular require harsh conditions for wet chemical dissolution in order to be run on ICP-MS. Direct analysis from the solid provides a cleaner sample preparation method, using a secondary electrode for analyses with DC-GD-MS.

The Thermo Scientific™ Element™ GD Plus GD-MS equipped with a pulsed power supply overcomes the analytical limitations associated with the use of a secondary electrode with high-vacuum GD sources. The μ s-pulsed fast-flow source provides state-of-the art solid sample analysis, at a sample throughput of several samples per hour.

Table 1. Instrumental parameters.

Parameter	Value
Matrix intensity	2×10^9 cps Al (MR)
Analysis time	10 min pre-sputter 10 min acquisition
Discharge voltage	1000 V
Pulse settings	~4 kHz repetition rate 50 μ s pulse duration
Anode parts	High purity graphite

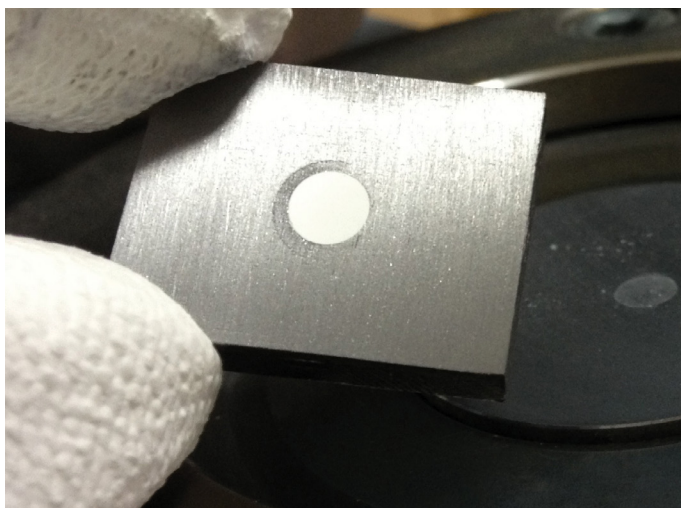


Figure 1. Sample preparation example for pressing non-conductive powder material into a secondary electrode.

Method

For sample preparation, a high purity Tantalum target was equipped with a borehole of approximately 5 mm diameter into which the sample was pressed. The target was placed on a TaW plate, and several tens of milligrams ceramic powder were filled into the borehole and pressed with a TaW pressing pin. The pressure should be adjusted to the kind of powder used. For fine-grained Al_2O_3 samples a pressure of ~ 0.4 tons, yielded stable and compact pellets, ready to be inserted without further treatment into the Element GD Plus GD-MS sample holder.

Results

- High purity alumina reference materials are reliably analyzed using Ta as a secondary electrode.
- Very good precisions are achieved (Table 2).
- The Standard RSF approach concept is shown to be valid for pulsed mode operation.
- High ionization potential Elements like boron are more efficiently ionized in pulsed mode. Therefore a dedicated RSF table should be applied.
- For the most important Elements, a matrix matched calibration can be easily established (Figure 2).

- The sample preparation method is simple, reproducible and clean.
- The Ta target used is easily resurfaced by grinding or milling for multiple use (Si contamination at low ppm level can originate from a SiC grinding step, grinding with corundum paper can serve as an alternative). Milling is therefore the preferred method for refurbishing the Ta target.
- Due to the high sensitivity of this GD-MS method ($\sim 2 \times 10^9$ cps for the matrix ion ^{27}Al , Medium Resolution), even at concentration levels as low as 0.01 ppm, good precisions are obtained (Table 1).
- Halogens are accessible for quantification at the ppm level.

Table 2. Semi-quantitative results of the high purity Al_2O_3 reference material NMIJ CRM 8007a (all concentration values in $\mu\text{g}\cdot\text{g}^{-1}$). Repeat analyses included sample preparation. Values in italics are information values.

Element	Measured conc.	Standard deviation of repeat analysis	Certified concentration
Fe	5.0	0.3	5.01 ± 0.25
Si	19.5	1.3	17.1 ± 0.4
Zr	2.5	0.6	1.80 ± 0.20
B	1.08	0.09	<i>0.21 ± 0.08</i>
Ca	2.4	1.0	<i>0.92 ± 0.14</i>
Cr	1.15	0.09	<i>0.84 ± 0.09</i>
Cu	1.25	0.06	<i>0.92 ± 0.08</i>
Mg	3.1	0.2	<i>2.8 ± 1.1</i>
Sr	0.025	0.007	<i>0.022 ± 0.009</i>
Ti	0.35	0.06	<i>0.26 ± 0.08</i>
Th	0.010	0.003	—
U	0.030	0.003	—

Conclusion

The Element GD Plus GD-MS in μs -pulsed operation mode is ideally suited for reproducible and accurate trace metal quantification of high purity alumina powders. The simple sample preparation avoids contamination and time-consuming dissolution steps, facilitating a close production control for ensuring highest quality products.

The reference material used is from the National Metrology Institute of Japan, Metrology Management Center, Reference Materials Office, 1-1-1, Umezono, Tsukuba, Ibaraki 305-8563, Japan: <http://www.nmij.jp/>

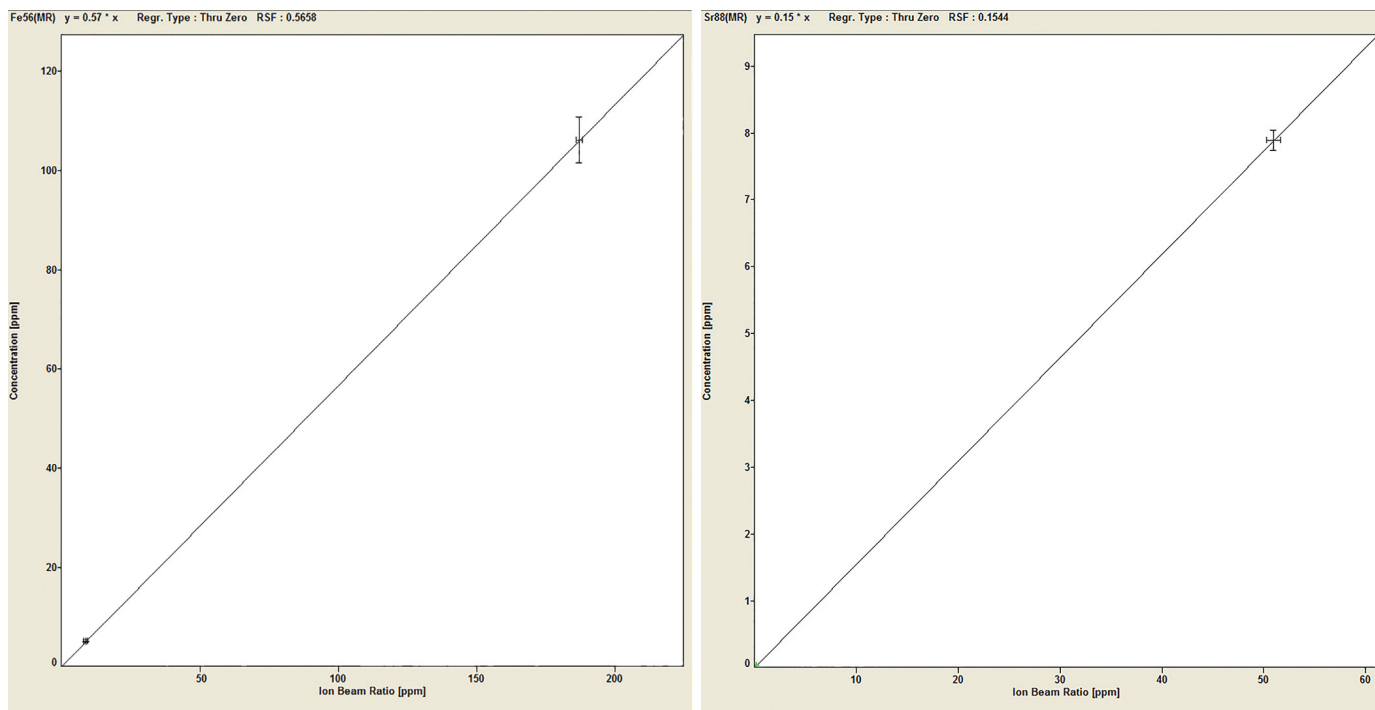


Figure 2. Calibration examples for the alumina reference materials NMIJ CRM 8006a and 8007a. Note the logarithmic scale.

Find out more at thermofisher.com/GD-MS

ThermoFisher
SCIENTIFIC

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific. 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.

AB30491-EN 0318S