

# Analysis of Alumina Powders Using the Thermo Scientific ELEMENT GD PLUS GD-MS

Joachim Hinrichs,<sup>1</sup> Karol Putyera,<sup>2</sup>

<sup>1</sup>Thermo Fisher Scientific, Bremen, Germany, <sup>2</sup>Evans Analytical Group, Liverpool, NY, USA

## Key Words

Ceramics, GD-MS, High Purity, Non-conductive,  $\mu$ s-Pulsed

## Goal

To demonstrate the capabilities of the Thermo Scientific ELEMENT GD PLUS  $\mu$ 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  $\text{Al}_2\text{O}_3$  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  $\mu$ s-pulsed fast-flow source provides state-of-the-art solid sample analysis, at a sample throughput of several samples per hour.

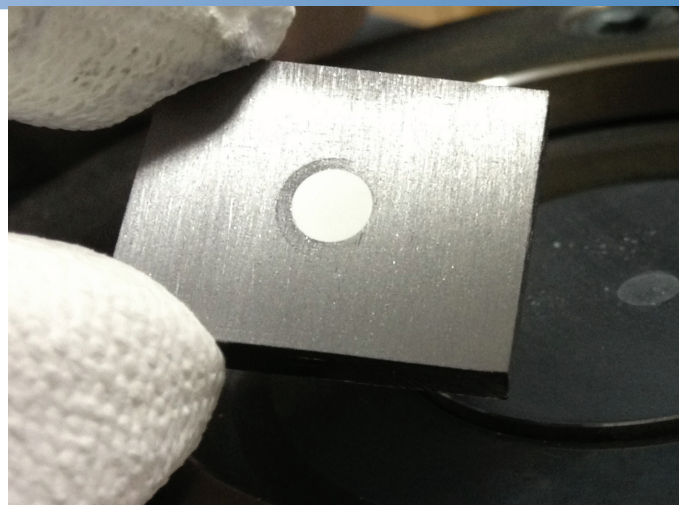


Figure 1. Sample preparation example for pressing non-conductive powder material into a secondary electrode.

## 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  $\text{Al}_2\text{O}_3$  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.

Table 1. Instrumental parameters.

Parameter	Value
<b>Matrix Intensity</b>	$2 \cdot 10^9$ cps Al (MR)
<b>Analysis Time</b>	10 min pre-sputter 10 min acquisition
<b>Discharge Voltage</b>	1000 V
<b>Pulse Settings</b>	~4 kHz repetition rate 50 $\mu$ s pulse duration
<b>Anode Parts</b>	High purity graphite

## 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 ( $\sim 2 \cdot 10^9$  cps for the matrix ion  $^{27}\text{Al}$ , 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  $\text{Al}_2\text{O}_3$  reference material NMJ 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
<b>Fe</b>	5.0	0.3	$5.01 \pm 0.25$
<b>Si</b>	19.5	1.3	$17.1 \pm 0.4$
<b>Zr</b>	2.5	0.6	$1.80 \pm 0.20$
<b>B</b>	1.08	0.09	$0.21 \pm 0.08$
<b>Ca</b>	2.4	1.0	$0.92 \pm 0.14$
<b>Cr</b>	1.15	0.09	$0.84 \pm 0.09$
<b>Cu</b>	1.25	0.06	$0.92 \pm 0.08$
<b>Mg</b>	3.1	0.2	$2.8 \pm 1.1$
<b>Sr</b>	0.025	0.007	$0.022 \pm 0.009$
<b>Ti</b>	0.35	0.06	$0.26 \pm 0.08$
<b>Th</b>	0.010	0.003	—
<b>U</b>	0.030	0.003	—

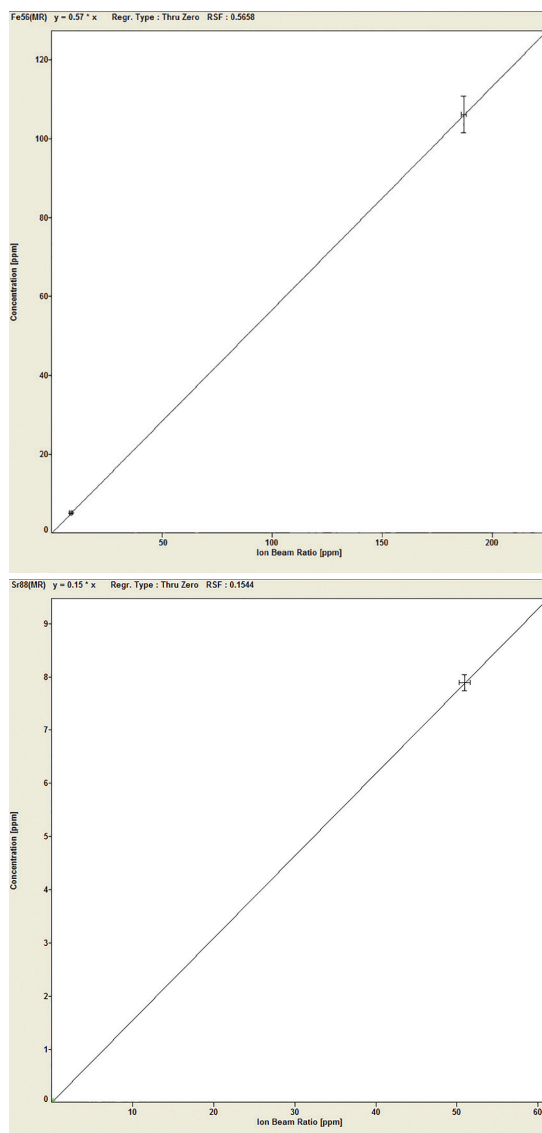


Figure 2. Calibration examples for the alumina reference materials NMJ CRM 8006a and 8007a. Note the logarithmic scale.

## Conclusion

The ELEMENT GD PLUS GD-MS in  $\mu\text{s}$ -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/>

[www.thermofisher.com/GDMS](http://www.thermofisher.com/GDMS)

©2016 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. National Metrology Institute of Japan is associated with the National Institute of Advanced Industrial Science and Technology. 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 manners 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