

Direct Determination of Ultratrace Elements in Semiconductor Grade Hydrofluoric Acid HF (48%) using the Thermo Scientific iCAP Qs ICP-MS

Tomoko Vincent¹, Simon Chang², Vicki Wu², Fred G. Smith³

¹Thermo Fisher Scientific, Bremen, Germany, ²Joy Allied Technology, Taipei, Taiwan,

³Teledyne CETAC Technologies, Omaha, Nebraska, USA.

Key Words

Hydrofluoric Acid, Semiconductor, Cold Plasma

Goal

To directly determine ultratrace metal concentrations in undiluted 48% semiconductor grade hydrofluoric acid using the Thermo Scientific iCAP Qs ICP-MS. Demonstrate the use of cold plasma to reduce background equivalent concentrations (BEC) and improve detection limits (LoD).

Introduction

Semiconductor process chemicals with ultratrace metal content are required in order to minimize production losses due to reduced performance or defects. Hydrofluoric acid (HF), for example, is routinely used as a wet chemical etchant to remove silicon dioxide layers in semiconductor devices. Routine, automated quantification of $\text{ng}\cdot\text{L}^{-1}$ level metal contaminants in HF is therefore required in order to maintain end product quality.

HF is an extremely corrosive acid and its vapor can cause serious harm to human health. Analytical techniques that minimize exposure to manufacturing personnel are therefore preferred. Additionally, due to the extremely low target concentration levels in semiconductor process chemicals, sample handling should be minimized in order to avoid contamination.

In this application, the Thermo Scientific™ iCAP™ Qs ICP-MS, in a single cold plasma measurement mode, was used for the direct trace metal analysis of 48% semiconductor grade HF. All samples were introduced into the iCAP Qs ICP-MS using a CETAC ASX-112FR Autosampler that offers the following features:

- Compact design for close coupling, leading to reduced uptake and washout times.
- Dual flowing rinse stations with solutions supplied via a gas displacement pump to minimize contamination.
- Metal free sample area and integrated cover to prevent airborne contamination.
- Flexible configuration of (PFA) autosampler vials.
- Fully control via Thermo Scientific™ Qtegra™ Intelligent Scientific Data Solution (ISDS).

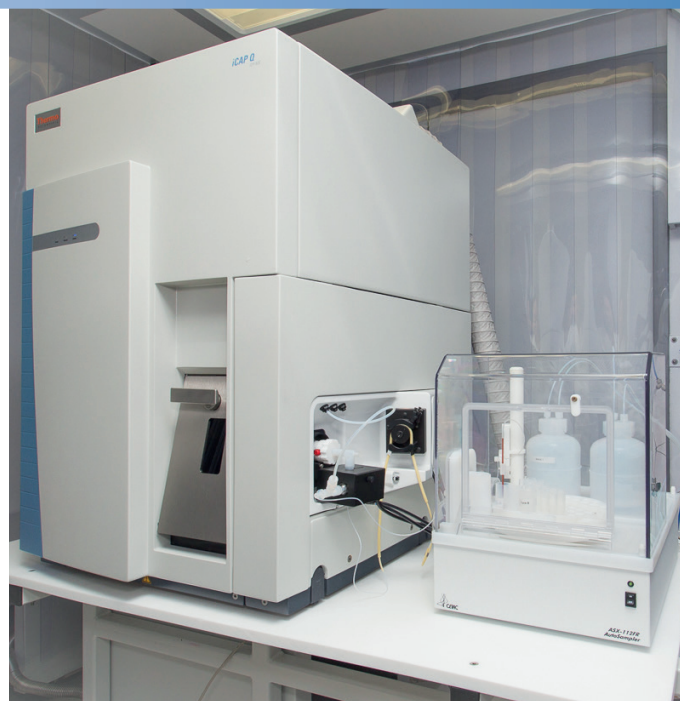


Figure 1: The Thermo Scientific iCAP Qs ICP-MS coupled with the CETAC ASX-112FR Autosampler.

Sample Preparation

Pre-cleaned PFA bottles were used for the preparation of all blanks, standards and samples. The bottles were rinsed with ultra pure water ($18.2 \text{ M}\Omega$) and left to dry in a laminar flow clean hood before use. Multi-element standards at concentrations of 10, 20, 50 and $200 \text{ ng}\cdot\text{L}^{-1}$ were prepared gravimetrically by adding the appropriate quantity of a multi-elemental SPEX Certiprep stock solution directly to the 48% HF samples. In order to assess recovery in the 48% HF, a spike recovery test at $20 \text{ ng}\cdot\text{L}^{-1}$ was performed.

Note: It is strongly recommended to pre-clean all sample vials and rinse bottles in an acid bath solution (3% HF, 2% HNO_3 , 2% H_2O_2) for 72 hours before use.

Instrument Configuration

The instrument configuration used is shown in Table 1. (Please note, the iCAP Qs ICP-MS is set up in non-clean room conditions.)

Table 1: Instrument configuration.

Parameter	Value
Spray Chamber	PFA cyclonic
Nebulizer	MicroFlow PFA-100 (self-aspirating)
Injector	2.0 mm I.D., Sapphire
Interface	Cold plasma platinum sampler and skimmer
Extraction Lens	Cold plasma lens kit

Results

Cold plasma (CP) inhibits the formation of the sample matrix induced ($^1\text{H}^1\text{F}$)₂ interference, allowing calcium to be measured at the most sensitive ^{40}Ca isotope. The improved performance that this allows is demonstrated in Figure 2 where BEC and LoD values of less than 5 ng·L⁻¹ were achieved for calcium in 48% HF.

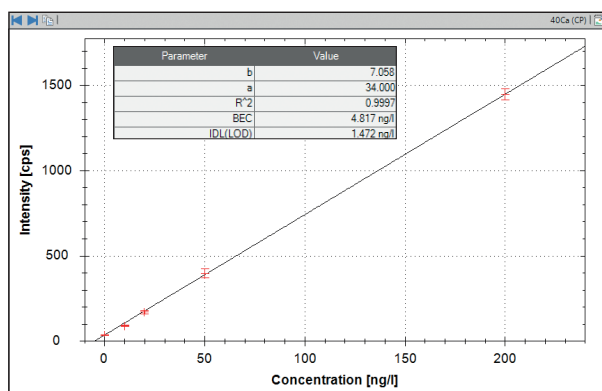


Figure 2: Cold plasma calibration curve for ^{40}Ca in 48% HF.

As can be seen in Table 2, the performance of the iCAP Qs ICP-MS for the direct analysis of 48% HF using cold plasma demonstrates:

- Excellent LoD and BEC values using a single analysis mode.
- 90% to 107% recoveries for a 20 ng·L⁻¹ spike.
- Suppression of plasma and sample matrix induced interferences for accurate analysis at ultratrace concentrations.

With a single mode analysis time of 75 seconds (including uptake and wash) high throughput analyses can be achieved.

Table 2: iCAP Qs ICP-MS performance data for 48% HF. Recovery values are shown as the percentage recovery for a 20 ng·L⁻¹ spike.

Analyte	LoD (ng·L ⁻¹)	BEC (ng·L ⁻¹)	Recovery (%)
^7Li	0.03	0.03	97
^{23}Na	0.2	1.6	100
^{24}Mg	0.5	0.6	101
^{27}Al	0.4	9.7	91
^{39}K	0.6	20.4	105
^{40}Ca	1.5	4.8	102
^{52}Cr	0.2	12.9	107
^{55}Mn	0.2	0.4	96
^{56}Fe	1.1	3.4	107
^{58}Ni	0.7	0.7	99
^{59}Co	0.5	0.7	100
^{63}Cu	1.6	4.2	91
^{66}Zn	3.5	5.0	106
^{88}Sr	0.5	0.1	99
^{107}Ag	0.6	0.7	97
^{111}Cd	0.7	0.4	90
^{115}In	0.1	0.05	96
^{205}Tl	0.1	0.03	101
^{208}Pb	0.2	0.02	93
^{209}Bi	0.5	0.1	104

Conclusion

The Thermo Scientific iCAP Qs ICP-MS has been shown to offer the high sensitivity and freedom from contamination and interference required for the direct determination of ultratrace (ng·L⁻¹) metal concentrations in semiconductor grade 48% HF. The method described is fast, simple and reliable: ideal for efficient and accurate analysis in exacting semiconductor laboratories.

www.thermofisher.com

©2016 Thermo Fisher Scientific Inc. All rights reserved. CETAC is a trademark of Teledyne Instruments Inc.; SPEX CertiPrep is a trademark of SPEX CERTIPREP, INC. 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
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