

# Dramatically Improved Performance of Arsenic and Selenium Determination using the Thermo Scientific iCAP Q ICP-MS with Methane Addition

Tomoko Vincent, Shona McSheehy Ducos, Julian D. Wills, Torsten Lindemann  
Thermo Fisher Scientific, Bremen, Germany

## Key Words

Arsenic, CCT, High Sensitivity, Methane Addition, Selenium

## Goal

To determine ultratrace concentration of As and Se using the Thermo Scientific iCAP Q ICP-MS and to demonstrate the use of CCT mode with CH<sub>4</sub> addition to improve analytical performance.

## Application Challenge

Arsenic and selenium ultratrace analysis is generally one of the most challenging using ICP-MS due to complicated polyatomic interferences arising from the argon plasma source and the inherent lower sensitivity of these analytes due to the high 1<sup>st</sup> ionization potential (IP).

While collision/reaction cell technology or high mass resolution enable increased performance, there is still room for improvement with regards to lowering detection limits and increasing sensitivity.

A hydrogen based reaction gas is known to improve the performance for Se analysis at  $m/z$  78. However, this method has the disadvantage of producing side reaction components such as  $^{38}\text{Ar}^{40}\text{ArH}_2^+$ , and therefore results in better BEC and LoD for  $^{78}\text{Se}$  compared to  $^{80}\text{Se}$ , despite the fact that  $^{80}\text{Se}$  is the more abundant isotope.

To further increase performance, an approach known as carbon loading has been used with the Thermo Scientific™ iCAP™ Q ICP-MS (Figure 1). Carbon loading is a technique known to improve the ionization efficiency of high IP analytes such as arsenic and selenium through what is thought to be a charge transfer process in the plasma. An approach using methane as the carbon source<sup>(1-3)</sup> has been used as it has been found in previous applications (data not shown) to be more effective than the addition of organic solvents.

Through the combination of sensitivity enhancement and interference removal, a method has been developed, which provides a significant increase in signal to noise performance enabling ultra low quantification limits for As and Se in various applications.



Figure 1. The Thermo Scientific iCAP Q ICP-MS.

## Sample Preparation

As and Se standards at concentrations of 1, 2, 3, 4 and 10 ng·L<sup>-1</sup> were prepared by diluting single stocks of SPEX CertiPrep solutions, directly in 2% HNO<sub>3</sub> prepared with Fisher Chemicals Optima grade nitric acid.

Table 1. Polyatomic interferences from an ICP source.

| Isotope          | Abundance (%) | Interference  |
|------------------|---------------|---|
| $^{75}\text{As}$ | 100           | $^{36}\text{Ar}^{38}\text{Ar}^1\text{H}^+$ , $^{40}\text{Ar}^{35}\text{Cl}$ |
| $^{74}\text{Se}$ | 0.87          | $^{36}\text{Ar}^{38}\text{Ar}$  |
| $^{76}\text{Se}$ | 9.02          | $^{40}\text{Ar}^{36}\text{Ar}$  |
| $^{77}\text{Se}$ | 7.58          | $^{36}\text{Ar}^{40}\text{Ar}^1\text{H}^+$ , $^{40}\text{Ar}^{37}\text{Cl}$ |
| $^{78}\text{Se}$ | 23.52         | $^{40}\text{Ar}^{38}\text{Ar}^+$  |
| $^{80}\text{Se}$ | 49.82         | $^{40}\text{Ar}_2^+$  |
| $^{82}\text{Se}$ | 9.19          | $^{40}\text{Ar}_2^1\text{H}_2^+$  |

## Instrument Configuration

A Thermo Scientific iCAP Qs ICP-MS was configured with a quartz, cyclonic spray chamber cooled to 2 °C and a self aspirating microflow PFA-400 nebulizer. The ICP-MS was used in CCT mode with 7% H<sub>2</sub> in He gas at 9 mL·min<sup>-1</sup>. 1% CH<sub>4</sub> in Ar gas was added to the plasma via an additional gas port on the spray chamber elbow at 120 mL·min<sup>-1</sup>. (Please note, the iCAP Qs ICP-MS was not setup in a cleanroom).

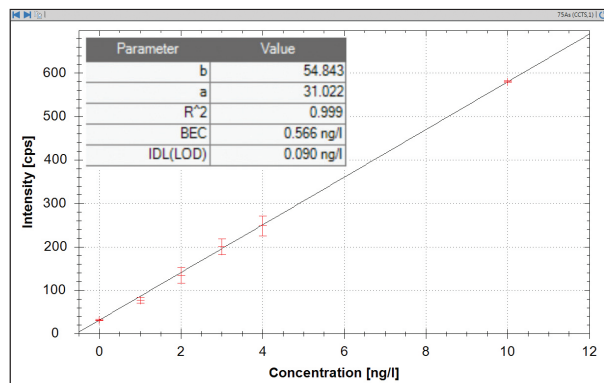


Figure 2a. Calibration curve for <sup>75</sup>As with CCT using CH<sub>4</sub> addition.

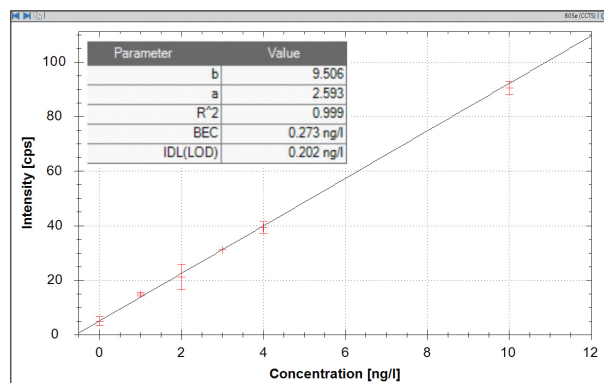


Figure 2b. Calibration curve for <sup>80</sup>Se with CCT using CH<sub>4</sub> addition.

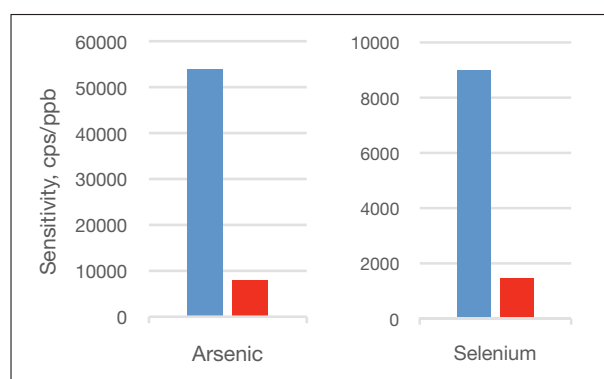


Figure 3. Sensitivity of As and Se without (red) and with (blue) the addition of methane to the plasma.

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

©2016 Thermo Fisher Scientific Inc. All rights reserved. All 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 10 3292 200  
**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

## Results

Linear calibration curves in sub ng·L<sup>-1</sup> concentration range with correlation coefficients better than 0.999 for both analytes (Figure 2a and 2b).

A dramatic improvement in sensitivity: <sup>75</sup>As 54 kcps·ppb<sup>-1</sup>, <sup>80</sup>Se 9 kcps·ppb<sup>-1</sup>, equivalent to a seven-fold increase in sensitivity compared to the same method without CH<sub>4</sub> addition (Figure 3).

Excellent LoD and BEC using a single CCT measurement mode with CH<sub>4</sub> addition.

Accurate spike recoveries from 95% to 107% for both elements at 1 ng·L<sup>-1</sup>.

Table 2. iCAP Qs ICP-MS performance data for As and Se in a 2% HNO<sub>3</sub> matrix. Recovery values are shown as the percentage recovery for a 1 ng·L<sup>-1</sup> spike.

| Analyte          | LoD (ng·L <sup>-1</sup> ) | BEC (ng·L <sup>-1</sup> ) | Recovery (%) |
|------------------|---------------------------|---------------------------|--------------|
| <sup>75</sup> As | 0.1                       | 0.6                       | 107          |
| <sup>80</sup> Se | 0.2                       | 0.3                       | 95           |

## Conclusions

The benefit of carbon loading to improve sensitivity of high IP analytes has been shown. The use of methane gas in the plasma is especially effective and has shown to give a sensitivity increase of seven fold for arsenic and selenium. Although the method has been developed with the Thermo Scientific iCAP Qs ICP-MS, the use of methane would be of equal advantage with the Thermo Scientific iCAP Qc ICP-MS and the Thermo Scientific™ ELEMENT 2 and Thermo Scientific™ ELEMENT XR. These instruments offer accurate and extremely low level quantification of arsenic and selenium with both BEC and LoD in the sub-ppt range. Methane addition is a simple technique that can be easily implemented for trace elemental analysis in various markets and matrices where ultra low detection capabilities are required.

## References

1. P. Field, M. LaVigne, K. Murphy, G. Ruiz and R. Sherrell *J. Anal. At. Spectrom.*, **22**, 1145–1151 (2007).
2. I. Rodushkin, P. Nordlund, E. Engstroem and D. Baxter *J. Anal. At. Spectrom.*, **20**, 1250–1255 (2005).
3. R. Santos *J. Anal. At. Spectrom.*, **29**, 152-161 (2014).

**Thermo**  
SCIENTIFIC

A Thermo Fisher Scientific Brand