

Analysis of drinking water in compliance with the U.S. EPA Method 200.8, Revision 5.4 using the Thermo Scientific iCAP MSX ICP-MS

Authors

Sukanya Sengupta¹, Bhagyesh Surekar¹, Daniel Kutscher¹, Sabrina Antonio²

¹Thermo Fisher Scientific, Bremen, Germany

²Thermo Fisher Scientific, San Jose, CA, USA

Keywords

ICP-MS, iCAP MSX ICP-MS, U.S. EPA Method 200.8, Revision 5.4, robustness, matrix tolerance, Argon Gas Dilution, quality control, regulatory compliance

Goal

To demonstrate the performance of the Thermo Scientific[™] iCAP[™] MSX ICP-MS for the analysis of a variety of water samples, including surface waters, ground waters and drinking waters, in accordance with U.S. EPA Method 200.8, Revision 5.4

Introduction

Analytical laboratories conducting routine water analysis often experience high sample workloads that demand quick turnaround and reporting of accurate results based on requirements that adhere to local, state, and federal regulations, as applicable. Compliance with comprehensive quality control protocols and detection limits requirements ensures accurate results and data quality in such laboratories. In the U.S., the approved analytical method for drinking water compliance monitoring by inductively coupled plasma mass spectrometry (ICP-MS) is U.S. EPA Method 200.8, Revision 5.4, which includes procedures for the analysis of dissolved elements in ground waters, surface waters, and drinking water samples, as well as total recoverable elements in these types of waters and in wastewater, soils, and sludges.

thermo scientific

This application note describes an analytical workflow developed using the iCAP MSX ICP-MS that enables laboratories to perform uninterrupted analysis of water samples for extended periods. The iCAP MSX ICP-MS offers a comprehensive solution for effective handling of a variety of sample types including those with high total dissolved solids (TDS) in one analytical run using Argon Gas Dilution (AGD). This entails the automatic and reproducible online addition of argon gas, using a dedicated mass flow controller, to the sample aerosol, enhancing robustness and reliability to consistently deliver accurate and precise results. ICP-MS instrumentation must be capable of analyzing varying sample types (e.g., drinking water, seawater, wastewater, soils), which include samples with high TDS. This approach of sample dilution using argon gas reduces the deposition of matrix on the interface cones and assures consistent readouts for QC checks, as well as reliable long-term robustness with minimal downtime.

Experimental

Instrument parameters

In this application note, an iCAP MSX ICP-MS was operated in conjunction with a Thermo Scientific[™] iSC-65 Autosampler for the accurate and reliable analysis of drinking and surface

Table 1. Instrument configuration and typical operating parameters

Parameter	Value
Nebulizer	iCAP MX Series ICP-MS nebulizer
Interface cones	Ni – tipped sample and skimmer
Spray chamber	Cyclonic quartz
Injector	Quartz, 2.5 mm ID
Torch	Quartz torch
Auxiliary flow (L·min-1)	0.8
Cool gas flow (L·min-1)	14
Nebulizer flow (L·min-1)	0.447
CRC conditions	Not used as per U.S. EPA Method 200.8 v5.4 requirements
AGD setting, argon flow (L·min ⁻¹)	AGD Level 5, 0.50
RF power (W)	1,550
Sampling depth (mm)	8
Number of replicates	3
Spray chamber temp. (°C)	2.7
Dwell time	0.05 s
Sweeps	5

water samples. A standard aqueous sample introduction system was used, along with integrated argon gas dilution and typical instrument parameters, as listed in Table 1. The Thermo Scientific[™] Qtegra[™] Intelligent Scientific Data Solution[™] (ISDS) Software was used to control the ICP-MS instrument and to generate, process, and report analytical data. The method was set up as a template in the software to ensure easy generation of data files in daily analysis, where the entire workflow meets the quality control requirements described in U.S. EPA Method 200.8.

Optimized instrument conditions were obtained using the autotune routines included with the Qtegra ISDS Software to ensure that analytical robustness and detection sensitivity are achieved during analysis.

The ICP-MS was operated in standard mode, as the use of traditional correction equations (Table 2) is mandated for the mitigation of isobaric and polyatomic interferences when U.S. EPA Method 200.8 is used for drinking water compliance monitoring under the Safe Drinking Water Act and the National Primary Drinking Water Regulations (NPDWR) 40 CFR Part 141. The use of a collision/reaction cell (CRC) is not permitted. Internal standardization was employed to monitor and compensate for physical interferences, instrument drift, and signal suppression or enhancement caused by the sample matrix.

Table 2. Summary of correction equations used for the target analytes

Analytes	Correction equation	
As	3.127 [(⁷⁷ Se - 0.815 × ⁸² Se)]	
Cd	1.073 [(¹⁰⁸ Cd - 0.712 × ¹⁰⁶ Cd)]	
Pb	[²⁰⁶ Pb + ²⁰⁷ Pb + ²⁰⁸ Pb]	
Мо	[0.146 × ⁹⁹ Ru]	
V	3.127 [(⁵³ Cr - 0.113 × ⁵² Cr)]	

Calibration standards

Calibration curves were generated for 21 analytes using six calibration standards and a calibration blank. Multi-element stock standards were prepared from single element standards of each target analyte (1,000 mg·L⁻¹, SPEX[™] CertiPrep[™], Metuchen, NJ, USA). Two different stock solutions were prepared at a concentration level of 2 mg·L⁻¹ to accommodate analytes with different concentrations and chemical compatibility. The stock solutions were then diluted gravimetrically using 1% (v/v) nitric acid as a diluent to result in the concentrations specified in Table 3. An internal standard solution containing 20 µg·L⁻¹ of ⁶Li, Sc, Y, Tb, and Bi was added on-line continuously during the analysis. All solutions used for analysis consisted of 1% (v/v) HNO₃ in ultrapure water containing 100 µg·L⁻¹ Au as a stabilizer for Hg.

Table 3. List of target analytes and their concentrations ($\mu g \cdot L^{-1}$) in calibration standards

Analytes	STD 1	STD 2	STD 3	STD 4	STD 5	STD 6	STD 7
Hg	0	0.01	0.1	0.25	1	2	5
Al, Sb, As, Ba, Be, Cd, Cr, Co, Cu, Pb, Mn, Mo, Ni, Se, Ag, Tl, Th, U, V, Zn	0	1	5	25	100	200	500

Please note that the concentration of selenium used in this study was 5 times lower compared to the requirements of EPA Method 200.8, Rev. 5.4

Table 4. List of analytes and their MCLs in drinking waters allowed by NPDWR and NSDWR

0.05

0.002

MCL		SMCL		
Analyte	MCL (mg·L⁻¹)	Analyte	MCL (mg·L⁻¹)	
Sb	0.006	AI	0.05 to 0.2	
As	0.010	Ag	0.1	
Ba	2.0	Cu	1.0	
Be	0.004	Mn	0.05	
Cd	0.005	Zn	5	
Cr	0.1			
Pb	0.015			
Hg	0.002			

Sample preparation

Se Tl

In total, five different water samples, collected locally, including tap water (typical TDS level 100–300 mg·L⁻¹), ground water (typical TDS level 500–700 mg·L⁻¹), and surface water (typical TDS level up to 1,000 mg·L⁻¹) were analyzed for the determination of dissolved elements. The samples were filtered using 0.45 µm pore size membrane filters followed by acidification using nitric acid to adjust the nitric acid concentration of the samples to approximately 1%. Multiple aliquots of these samples (equally split between ground water, surface water, and tap water) were then analyzed to assess accuracy and precision, performance of duplicate measurements, and robustness over a 16-hour analysis. A certified reference material (CRM) SLRS-5 (natural river water) was included in the analysis as an additional verification of accuracy.

Results and discussion Detection limits and linearity

The detection limits of the method must be below the standards for inorganic contaminants in drinking waters set by the National Primary Drinking Water Regulations (NSDWR) and the National Secondary Drinking Water Regulations (NSDWR) (Table 4) with the maximum contaminant level (MCL) and secondary maximum contaminant level (SMCL) for each inorganic contaminant or analyte in this study. Here, the method detection limits (MDLs) for all analytes were calculated following the guidance provided in U.S. EPA Method 200.8, Revision 5.4. Seven replicate aliquots of fortified calibration blank (fortified at 2-5 times the concentration level of the expected detection limits) were analyzed, and the standard deviation of these measurements was multiplied by the Student's *t* value, which is 3.14 for seven replicate measurements. The results are summarized in Table 5. The MDLs achieved for all analytes were below the requirement of the regulations, suggesting that the developed method enables sensitive determination of all target analytes, meeting or exceeding the requirements. The correlation coefficients (R²) obtained for all analytes were greater than 0.999, which suggests excellent linear response for the established concentration range for each analyte. The analytes, with their mass-to-charge ratio (m/z) and calibration R² are also summarized in Table 5.

Table 5. List of analytes, m/z,	correlation	coefficients,	and MDLs
---------------------------------	-------------	---------------	----------

Analyte	m/z	R ²	MDL (µg·L⁻¹)
Ag	107	0.9995	0.052
Al	27	0.9992	0.228
As	75	0.9995	0.046
Ba	137	0.9998	0.025
Be	9	0.9996	0.022
Cd	111	0.9998	0.022
Со	59	0.9994	0.010
Cr	52	0.9992	0.045
Cu	63	0.9995	0.061
Hg	202	0.9993	0.005
Mn	55	0.9990	0.026
Мо	98	>0.9999	0.018
Ni	60	>0.9999	0.037
Pb	206+207+208	0.9996	0.010
Sb	123	0.9994	0.020
Se	82	0.9998	0.193
Th	232	0.9991	0.046
TI	205	0.9996	0.013
U	238	0.9993	0.024
V	51	0.9996	0.038
Zn	66	0.9995	0.122

Quality Control (QC) checks

Laboratory Reagent Blank (LRB) – LRB sample was prepared from 1% (v/v) HNO_3 in ultrapure water, which was treated similarly as real samples by exposing it to the filtration through a 0.45 µm pore diameter membrane filter. The LRB sample was analyzed every 15 samples, and resulting concentrations of all analytes were monitored. The concentrations of all analytes in subsequent LRB samples were found to be within the acceptable range as specified in U.S. EPA Method 200.8.

Laboratory Fortified Blank (LFB) – An aliquot of LRB was spiked with stock solutions to yield the final concentration of 75 μ g·L¹ for all analytes (except for Hg, which spiked at a concentration level of 0.75 μ g·L¹). The recovery of each analyte was calculated automatically within the Qtegra LabBook using a recovery test available in the comprehensive QC toolkit. The recovery values for all analytes were found to be within the range of 85–115%, which suggests that the acceptance criteria were met, and analysis could be continued. **Quality Control Sample (QCS)** – QCS solution was prepared by spiking an aliquot of the LRB using independent analyte stock solutions (different from the stock solutions used for preparation of calibration standards) to yield analyte concentrations at $50 \ \mu g \cdot L^{-1}$ level (except for Hg, which was spiked at $0.5 \ \mu g \cdot L^{-1}$ concentration level). The QCS was used to confirm the initial validity of calibration standards as well as to confirm the ongoing instrument performance during the measurement of more than 16 hours. The QCS was analyzed every 10 samples, with the percent accuracy for all analytes in each QCS found to be well within the acceptable range of $\pm 10\%$ (equivalent to 90-110%) of the true concentration.

Accuracy

Laboratory Fortified Matrix (LFM) – To assess and demonstrate method accuracy and precision, one of the water samples was spiked in duplicate at a concentration level of $50 \ \mu g \cdot L^{-1}$ for all analytes (except for Hg, which was spiked at a concentration level of $0.5 \ \mu g \cdot L^{-1}$). The percent recovery of each analyte was calculated automatically within Qtegra ISDS Software based on the measured concentration for the unspiked and spiked replicates of the LFM. The average percent recovery for all analytes was found to be within the range of 85–115%, against the acceptable range of 70–130%, with relative percent difference of less than 5% between duplicate spike measurements. Table 6 lists the analytes and their average percent recovery calculated from duplicate measurements of spiked water sample.

Analysis of Certified Reference Material (CRM) and

observed accuracy – To further demonstrate method accuracy, a certified reference material (SLRS-5, natural river water) was analyzed three times during a long-term analysis of more than 16 hours. The percent accuracy of each analyte was calculated based on the concentration data obtained during the analysis and the certified values given for each analyte in the SLRS-5 CRM. Table 7 summarizes the list of analytes, their certified concentration values, and calculated percent accuracy. The percent accuracy values given are the average calculated from three different measurements of SLRS-5 CRM.

Analyte	Concentration in unspiked water sample (µg·L ⁻¹)	Measured spike concentration (µg·L ⁻¹)	Spike recovery (%), n=2
Ag	2.4	54.1	108.2
Al	15.5	48.9	97.8
As	0.015	52.4	104.9
Ва	14.3	51.1	102.3
Be	0.036	52.5	105.1
Cd	0.032	51.0	102.1
Со	0.087	51.0	102.0
Cr	0.084	51.0	102.0
Cu	89.9	52.2	104.4
Fe	91.3	52.0	104.0
Hg	0.019	0.52	104.9
Mn	0.011	51.9	103.9
Мо	0.312	52.3	104.6
Ni	5.4	46.8	93.6
Pb	0.304	54.0	108.0
Sb	0.024	54.8	109.6
Se	0.122	49.5	99.1
Th	0.095	55.2	110.3
TI	0.158	54.0	107.9
U	0.055	55.1	110.3
V	<dl< td=""><td>51.6</td><td>103.3</td></dl<>	51.6	103.3
Zn	79.3	49.9	99.7

Table 6. Average percent recovery calculated from the duplicate measurement of spiked water sample (spike levels $0.5 \ \mu g \cdot L^{-1}$ for Hg and 50 $\ \mu g \cdot L^{-1}$ for other analytes)

Table 7. List of analytes, certified concentrations and accuracy (as percent) of the certified reference material (CRM) SLRS-5

Analyte	Certified concentration ($\mu g \cdot L^{-1}$)	Accuracy (%), n=3
Al	49.5 ± 5.0	95.2 ± 5
As	0.413 ± 0.039	94.2 ± 7
Ba	14.0 ± 0.5	112.1 ± 3
Be	0.005*	100.0 ± 3
Со	0.05*	106.8 ± 3
Cr	0.208 ± 0.023	93.8 ± 6
Cu	17.4 ± 1.3	94.4 ± 2
Fe	91.2 ± 5.8	105.1 ± 6
Mn	4.33 ± 0.18	104.2 ± 4
Мо	0.27 ± 0.04	103.4 ± 5
Ni	0.476 ± 0.064	91.6 ± 1
Pb	0.081 ± 0.006	97.3 ± 2
Sb	0.3*	86.2 ± 4
U	0.093 ± 0.006	93.0 ± 7
V	0.317 ± 0.033	101.3 ± 2
Zn	0.845 ± 0.095	97.4 ± 3

*Information values, as per CRM certificate of analysis



Figure 1. Internal standard recovery plot from a Qtegra LabBook, highlighting the consistent recovery of internal standards within 80–112% throughout the 16-hour continuous analysis

Method robustness

A robust instrument and optimized method are key requirements when analyzing varying sample matrices within an analytical run. Instrument drift and physical interferences due to the sample matrices can cause signal enhancement or suppression and must be corrected for. In addition, U.S. EPA Method 200.8 includes specific guidelines and requirements, such as internal standardization, to monitor and correct for instrument drift and physical interferences. An internal standard must be added to all standards and samples in the run sequence and must not deviate outside of the acceptable range of 60–125% compared to the response observed in the calibration blank.

As previously mentioned, an internal standard solution containing 20 µg·L⁻¹ of ⁶Li, Sc, Y, Tb, and Bi was added online during the analysis. Figure 1 shows a plot of the internal standards recovery against the calibration blank throughout the 16-hour analysis. As shown, all internal standards recoveries were within the range of 80–112%, well within the acceptance criteria of U.S. EPA Method 200.8. These results demonstrate the robustness of the method and instrument setup that included online sample dilution.

Conclusions

The quality of analytical data obtained during this study indicates that the iCAP MSX ICP-MS, equipped with a built-in AGD system for controlled and automatic dilution of the sample aerosol, is a powerful solution for the analysis of varying sample matrices within a single run sequence. Some of the important outcomes of this study are highlighted here.

- The MDLs achieved comfortably meet the requirements of the method and were below the regulations for inorganic contaminants in drinking water. This indicates that the instrument setup, with optimized sample dilution, provides the sensitivity and robustness for the analysis of a variety of water samples in a single, extended analysis.
- The accuracy and precision obtained from the analysis of the LFM and SLRS-5 CRM indicate that the developed method and instrument setup were optimized for the variety of samples analyzed. Furthermore, the ease of instrument tuning, operation, and LabBook set up through the Qtegra ISDS Software allows routine analysis of varying environmental samples with minimal effort while improving productivity.
- The accurate results obtained from the analyses of the QCS samples demonstrate on-going instrument performance and calibration over the extended analysis of varying sample matrices.
- The internal standard recoveries of 80–112% observed during the 16-hour analysis demonstrate robustness of the instrument and that the dilution applied to each sample was helpful in minimizing the effects of physical interferences.
- The QC toolkit included within the Qtegra ISDS Software enables easy set up and compliance with the comprehensive QC protocol of U.S. EPA Method 200.8. Customizable report templates are available to provide all data required for audit and regulatory purposes. The use of the flags and limits features within the software helps highlight and call out data falling outside the acceptable concentration ranges and provides easy visualization and review of data.

References

- 1. U.S. EPA Method 200.8. https://www.epa.gov/sites/default/files/2015-08/documents/ method_200-8_rev_5-4_1994.pdf
- 2. National Primary Drinking Water Regulations. https://www.epa.gov/ ground-water-and-drinking-water/national-primary-drinking-water-regulations
- 3. National Primary Drinking Water Regulations. https://www.epa.gov/sdwa/secondarydrinking-water-standards-guidance-nuisance-chemicals#what-are-secondary

General Laboratory Equipment – Not For Diagnostic Procedures. © 2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. SPEX CertiPrep is a trademark of SPEX CertiPrep, Inc. 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. AN002911 0424

thermo scientific