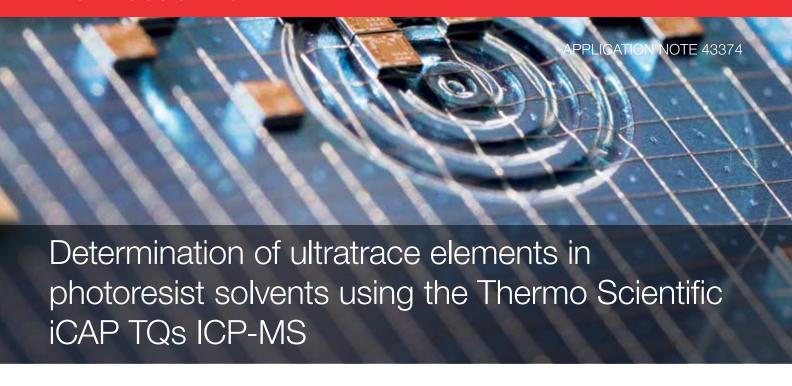
thermoscientific



Authors

Tomoko Vincent, Product Specialist, Thermo Fisher Scientific

Keywords

Cold plasma, iCAP TQs, KED, Mass shift, NMP, PGMEA, Photoresist, Triple quadrupole, Semiconductor, Ultratrace

Goal

To determine ultratrace metal concentrations in semiconductor grade NMP and PGMEA. To demonstrate the use of cold plasma and triple quadrupole technologies to reduce background equivalent concentrations (BEC) and detection limits (LOD). To demonstrate the use of the Thermo Scientific™ iCAP™ TQs ICP-MS for performing reproducible, ultratrace ng·L¹ (ppt) measurements with reliable switching between multiple analysis modes (hot/cold plasma, single/ triple quadrupole) within a single measurement.

Introduction

Propylene glycol methyl ether acetate (PGMEA) and N-methyl-2-pyrrolidone (NMP) are the base organic solvents of semiconductor photoresists used in the production of sub-µm structures in microelectronic manufacturing. As photoresists come into direct contact with wafer surfaces, their components must be controlled for trace metal purity.

Because of its high elemental sensitivity, ICP-MS is widely used in quality control analyses of materials in the semiconductor industry. A simple ICP-MS analytical method for the direct analysis of photoresist solvents would provide a useful control for ultratrace (ng·L¹) levels of analytes and avoid any contamination caused by any sample preparation.

PGMEA and NMP have historically been considered as difficult matrices to analyze directly by ICP-MS due to their high volatility (see Table 1) and high carbon content which can lead to significant polyatomic interferences.



In this application note, cold plasma, kinetic energy discrimination and triple quadrupole ICP-MS technologies are combined within a single analytical method in the Thermo Scientific iCAP TQs ICP-MS to remove all polyatomic interferences. With the use of cold plasma, the ICP ion source is run at a significantly lower forward power, effectively suppressing the ionization of argon and carbon and therefore eliminating any subsequent polyatomic species that would otherwise interfere with target analyte ions. For some analytes that are more sensitive under hot plasma conditions, a triple quadrupole mass shift analysis mode was automatically selected for accurate, interference free analysis.

Table 1. Chemical properties.

	Vapor Pressure (mm⋅Hg) at 20 °C				
H ₂ O	17.5				
PGMEA	3.7				
NMP	< 0.3				

Sample preparation

Precleaned PFA bottles were used for the preparation of all blanks, standards and samples. The bottles were rinsed with ultrapure water (18.2 MOhm cm) and left to dry in a laminar flow clean hood before use. Standards at concentrations of 200, 400 and 1000 ng·L¹ were prepared by gravimetrically adding the appropriate quantity of a multielemental stock solution (prepared from single element standards in semiconductor grade isopropyl alcohol) directly to the PGMEA and NMP samples. Semiconductor grade PGMEA was used for the rinse and blank solutions. Recovery tests were performed using 100 ng·L¹ spikes.

Instrumentation

A dedicated sample introduction system was used for the routine, direct analysis of organic solvents. The introduction system consisted of a 100 μL·min⁻¹ self-aspirating PFA micro flow concentric nebulizer (Elemental Scientific, Omaha, NE, USA) and a peltier cooled quartz spraychamber (at -10 °C). High purity oxygen (30 mL·min⁻¹) was added to the nebulizer aerosol stream via a port in the spraychamber elbow to prevent carbon matrix build up on the interface. A 1.0 mm diameter quartz injector minimized carbon loading of the plasma.

Platinum tipped sample and skimmer cones were necessary because of the oxygen addition. All samples were presented for analysis using a Teledyne CETAC Autosampler ASX-112FR System (Omaha, NE, USA). The iCAP TQs ICP-MS was operated in SQ-KED (single quadrupole He KED), SQ-CP-NH $_3$ (single quadrupole, cold plasma with NH $_3$ as a reaction gas) and TQ-O $_2$ mass shift (triple quadrupole with O $_2$ as a reaction gas) modes. All parameters are presented in Table 2. The iCAP TQs ICP-MS used in this study was not installed in a cleanroom.



Table 2. Instrument configuration and operating parameters.

Table 2. Instrument configuration and operating parameters.					
Parameter	Value				
Nebulizer	PFA concentric nebulizer 100 μL·min ⁻¹ (self-aspirating)				
Spraychamber	Quartz cyclonic spraychamber cooled at -10 °C				
Additional Gas	Pure O ₂ gas, 30 mL·min ⁻¹				
Injector	1.0 mm I.D., Quartz				
Interface	Pt sampler and Pt (high sensitivity) skimmer				
Extraction Lens	Cold plasma lens				
Measurement Mode	Sir quad	Triple quadrupole			
	SQ-KED	SQ-CP-NH ₃	TQ-O ₂		

Forward Power 1550 W 540 W 1550 W Nebulizer Gas 1.17 L·min-1 0.68 L·min-1 1.17 L·min-1 Pure O., Pure He, Pure NH_a, **CRC** Gas 4.2 mL·min-1 0.28 mL·min-1 0.3 mL·min-1 **Dwell Time** 100 to 300 ms per analyte, 5 sweeps Lens Setting Optimized using autotune

Identical instrument parameters were used for the analysis of both the NMP and PGMEA solvents for improved sample throughput without any sacrifice in performance.

Results and discussion

The Thermo Scientific iCAP TQs ICP-MS system is a powerful analytical tool for multi-element analysis in a range of routine applications. By providing the analyst with unlimited flexibility of ICP-MS technologies (cold plasma, kinetic energy discrimination and triple quadrupole analysis modes), the ultimate performance can be achieved, specifically tailored for each application. Through the implementation of Reaction Finder in the Qtegra™ Intelligent Scientific Data Solution™ (ISDS) Software, the choice of isotope and analysis mode for each target element is made automatically, making triple quadrupole ICP-MS as easy to use as single quadrupole systems.



Figure 1. iCAP TQs ICP-MS. In $TQ-O_2$ triple quadrupole mode, Q2 (collision/reaction cell) is pressurized with oxygen as a reaction gas, Q1 is set to the analyte mass (M $^{+}$) and Q3 is set to the product ion mass (MO $^{+}$) or (M $^{+}$).

For example, using the TQ- O_2 mass shift mode (schematically shown in Figure 2), the C, N, O and Ar based polyatomic interferences are efficiently removed leading to accurate analysis of As. In TQ- O_2 mass shift mode, the first quadrupole (Q1) uses intelligent mass selection (iMS) to reject unwanted ions. The second quadrupole (the Q2 collision/reaction cell) is used to selectively shift the As target analytes to the AsO+product ions using O_2 as the reaction gas, while any remaining carbon based polyatomic interference will not react with O_2 . The third quadrupole (Q3) then isolates the AsO+product ions = and removes any remaining interferences through a second stage of mass filtration allowing completely interference free quantification.

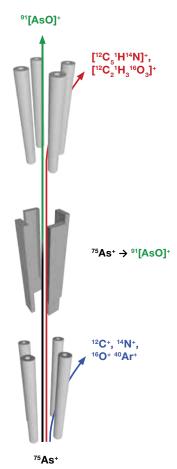


Figure 2. Schematic showing TQ-O₂ mass shift mode for the analysis of arsenic (75As as 91[AsO]*).

4	Label V9	Maniflant VS	Frauer 74	Serple Type V 4	Standard V-P
	PGMEA	10	~	ELK	-
	PGMEA =100ppx	3	8	STD	1ppb mix
1	PGMEA =200ppt	3	2	570	1ppb mix
	PGMEA -1000pet	3	2	STD	1ppb mix
	wash	3	2	UNKNOWN	
6	Recovery test PGMEA 100ppt	3	~	UNKNOWN	
1	wash	3	- 20	UNKNOWN	
*=	NAP	10		6LK	
	NMP + 100ppt	3	- 2	STD	1ppb mix
10	NMP + 200ppt	3	- 2	STD	Topt mix
#1	MMP + 1000ppt	3	· 6	STD	Toph mis
12	wash	3	₩.	UNKNOWN	*******
15	Recovery test NMP 100ppt	3	2	UNIVOYONN	

Figure 3. Sample list from Qtegra ISDS Software. NMP and PGMEA solvents were analyzed with identical instrument and method parameters within a single LabBook.

A summary of the results from the trace elemental analysis of semiconductor grade PGMEA and NMP is shown in Table 3. For each of the twenty one analytes, the analysis mode, detection limits (LOD), calculated from three times the standard deviation of ten replicate measurements of the calibration blank and spike recovery are provided.

Table 3. LOD, recovery data for the analysis of PGMEA and NMP. Please note that LOD values are dependent on the sample provided for analysis and do not necessarily demonstrate the ultimate performance possible.

Analyte	Analysis Mode	LOD (I	ng∙L ⁻¹)	Spike Recovery (100 ng·L ⁻¹)	
		PGMEA	NMP	PGMEA	NMP
⁷ Li	SQ-CP-NH ₃	0.2	0.9	97%	98%
²³ Na	SQ-CP-NH ₃	1.3	10.2	98%	97%
²⁴ Mg	SQ-CP-NH ₃	0.4	5.0	97%	91%
²⁷ AI	SQ-CP-NH ₃	1.1	3.8	97%	100%
³⁹ K	SQ-CP-NH ₃	1.3	6.4	102%	106%
⁴⁰ Ca	SQ-CP-NH ₃	1.2	8.1	96%	94%
⁵² Cr	SQ-CP-NH ₃	2.0	3.3	97%	92%
⁵⁵ Mn	SQ-CP-NH ₃	0.3	8.4	93%	99%
⁵⁶ Fe	SQ-CP-NH ₃	3.1	2.0	104%	103%
⁵⁹ Co	SQ-CP-NH ₃	0.4	0.7	95%	100%
⁶³ Cu	SQ-CP-NH ₃	1.2	3.4	98%	96%
51 V	SQ-KED	2.6	2.8	105%	96%
⁶⁶ Zn	SQ-KED	14.3	18.8	94%	100%
⁹⁸ Mo	SQ-KED	1.6	2.1	97%	111%
¹⁰⁷ Ag	SQ-KED	1.5	1.8	97%	111%
¹¹¹ Cd	SQ-KED	3.7	2.6	85%	107%
¹³⁸ Ba	SQ-KED	0.1	0.0	94%	101%
²⁰⁸ Pb	SQ-KED	0.8	1.4	97%	108%
⁴⁸ Ti at 64 <i>m/z</i>	TQ-O ₂	1.9	3.7	88%	106%
⁷⁵ As at 91 <i>m/z</i>	TQ-O ₂	2.0	2.4	99%	106%
⁸⁰ Se at 96 <i>m/z</i>	TQ-O ₂	1.9	4.6	95%	109%

Spike recoveries of between 88% and 111% were obtained for all twenty-one elements in NMP and PGMFA.

The low detection limits achieved for the more challenging analytes in NMP and PGMEA organic solvents showcase the powerful interference removal capabilities from the combination of cold plasma and triple quadrupole technologies on the iCAP TQs ICP-MS.

Calibration curves

Calibration curves for Mg, Ti, As and Ag in PGMEA are shown in Figure 4. The calibrations, performed with calibration points at 100, 200 and 1000 ng·L¹, demonstrate the excellent sensitivity and linearity achieved with different types of organic solvent using common instrument parameters.

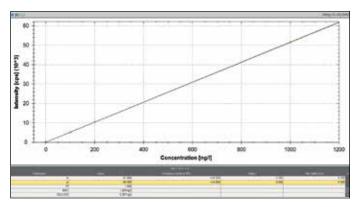


Figure 4a. Mg in PGMEA using SQ-CP-NH₃ mode.

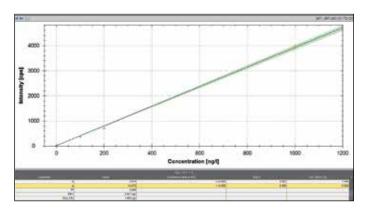


Figure 4b. Ti in PGMEA using $TQ-O_2$ mass shift mode.

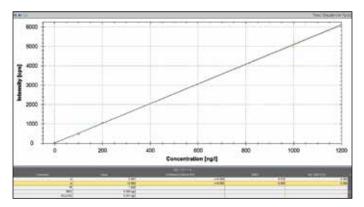


Figure 4c. As in PGMEA using TQ-O, mass shift mode.

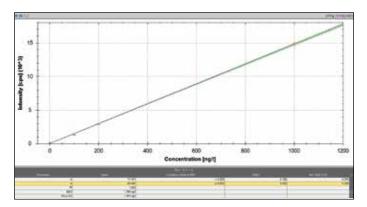


Figure 4d. Ag in PGMEA using CP-SQ-KED.

Conclusion

The Thermo Scientific iCAP TQs ICP-MS has been shown to provide sensitive and accurate multielemental analysis of NMP and PGMEA photoresist solvents at ultratrace (ng·L¹) concentration levels. The combination of single and triple quadrupole technologies with hot and cold plasma modes – even in volatile organic solvents – provides the analytical flexibility and instrumental performance to deliver the detection limits required for semiconductor level analyses.

Although NMP and PGMEA have different chemical formulae and properties, the iCAP TQs ICP-MS allows the use of a single analytical method for both solvents, simplifying the analytical procedure and improving productivity.

Find out more at thermofisher.com/TQ-ICP-MS

For Research Use Only. Not for use in diagnostic procedures. ©2018 Thermo Fisher Scientific Inc. All rights reserved. Teledyne CETAC is a trademark of Teledyne Instruments, 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 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. AN43374-EN 0218

