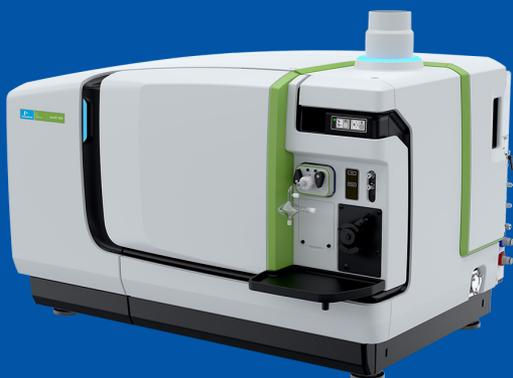


ICP - Mass Spectrometry



Novel Interference Removal Opportunities with the NexION 5000 ICP-MS

Introduction

Multi-quadrupole technology, in combination with controlled reactions in the Universal Cell

(UCT) of the NexION® 5000 ICP-MS, has opened new possibilities for spectral interference removal.¹ As was described in the technical note "Interferences in ICP-MS: Do we still have to worry about them?"², the MS/MS and Mass Shift modes, in conjunction with three cell analytical modes (Std, KED and DRC), can solve many challenging problems, not previously possible using a single quadrupole system. This work is a continuation of a discussion on spectral interference removal and presents dynamic reactions with three gases (O_2 , NH_3 and CH_4) that, in tandem with multi-quadrupole technology, are a powerful tool in interference management on the NexION 5000 ICP-MS.

Quadrupoles in the NexION 5000 ICP-MS

Each quadrupole in the NexION 5000 system plays a significant role in transmitting elements of interest to the detector. The Quadrupole Ion Deflector (QID – Q0) rejects photons and neutral species and, at the same time, transfers a selected mass range into the spectrometer, enhancing sensitivity and, in some cases, rejecting some sample matrices. The second quadrupole (Q1), placed before the Universal Cell (Q2), can work as a mass-filtering device with a nominal resolution of 0.7 amu. The mass-filtering ability of Q1 allows only ions of a specific mass to be passed into the Universal Cell, while all other ions from the matrix, solvents and plasma not residing at the mass of interest are rejected. This capability has a significant impact on the reduction of spectral interferences which may arise from reactions of lower mass matrix ions with gases in the cell. Q1 works together with the quadrupole Universal Cell where up to four reactive or non-reactive gases can be used to create desirable reactions or collisions. Q3, the fourth quadrupole in the NexION 5000 ICP-MS, works predominantly as a resolving transmission analyzer. In MS/MS mode, Q1 and Q3 are set to the same analyte mass, while in Mass Shift mode, Q1 is set to a mass of interest and Q3 is set to a higher mass where the polyatomic ion, product of a reaction of an analyte with a reaction gas, resides. In Mass Shift mode, analyte products are measured on masses free from interferences, as discussed in the following sections.

it is the opposite (Figure 3): a weak cluster at mass 254 and a strong one at mass 270. Of course, when an analyte creates multiple product ions, their sensitivity depends on the reaction gas density in the UCT. In general, the optimization of the gas flow can change the sensitivity ratio of the analyte's product ions – higher flows promote higher mass clusters.

In this next example (Figure 4), $^{98}\text{Mo}^+$ creates an interference on the Cd isotope at mass 114. However, if the UCT is pressurized with O_2 and gas density increases, MoO_2^+ at mass 130 becomes a dominant species and Cd can be determined in MS/MS mode at mass 114 free from interference.

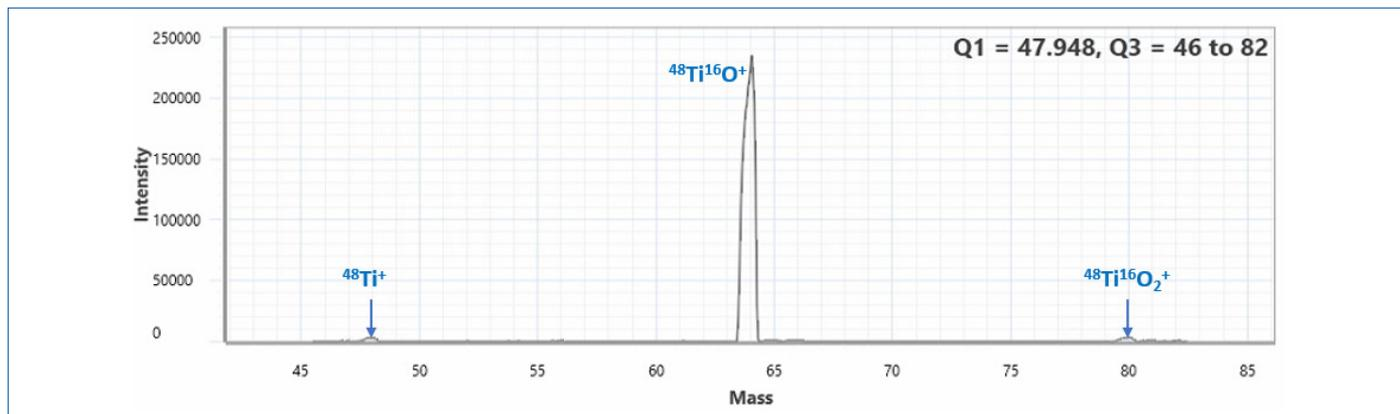


Figure 2: Ti product ion scan with O_2 .

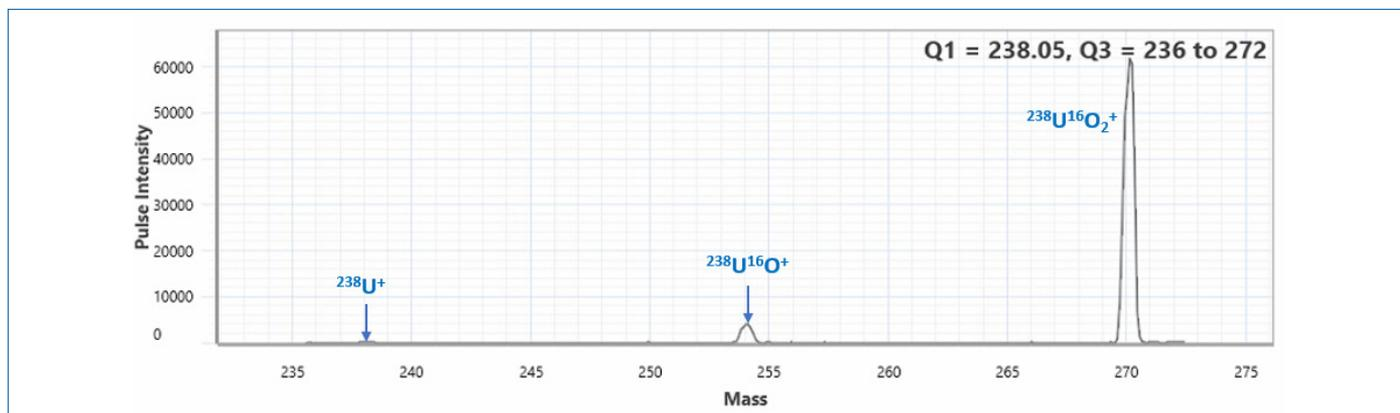


Figure 3: U product ion scan with O_2 .

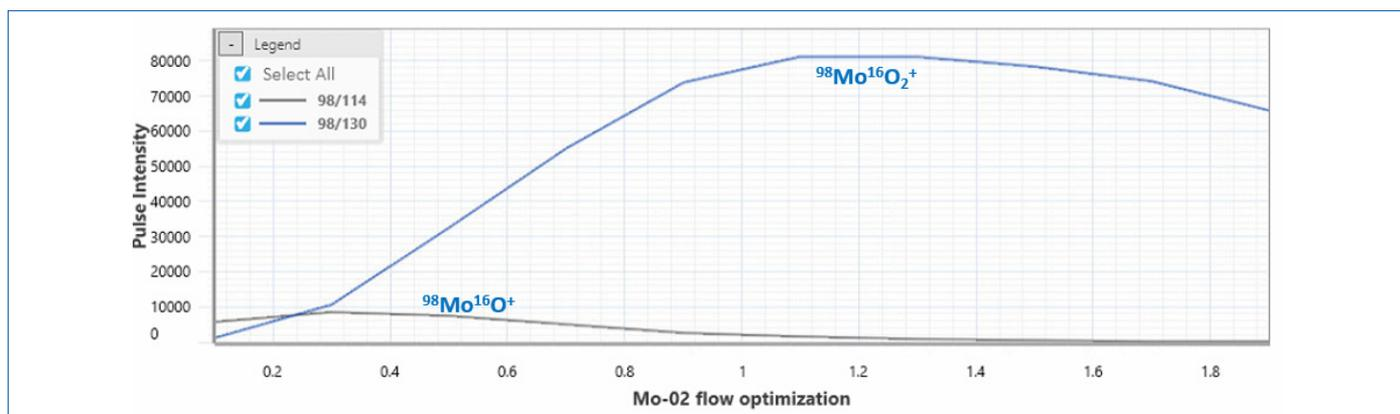


Figure 4: Optimization of O_2 flow for Mo^+ at mass 114 and MoO_2^+ at mass 130.

The dynamic bandpass, especially the RPq parameter, could promote some clustering reactions by increasing the reaction efficiency for some analytes (e.g. Cr, Mn, Fe, Co, Ni, Cu, Zn and Se) that do not react or that react weakly with O₂ under standard conditions with RPq values of 0.4-0.6.

In Figure 5, an optimization of the RPq parameter indicates that sensitivity of the CrO⁺ cluster at mass 68 is the highest at RPq=0.8. The Mass Shift approach can be an option when spectral interferences on Cr at mass 52 cannot be completely removed by MS/MS in DRC or KED modes.

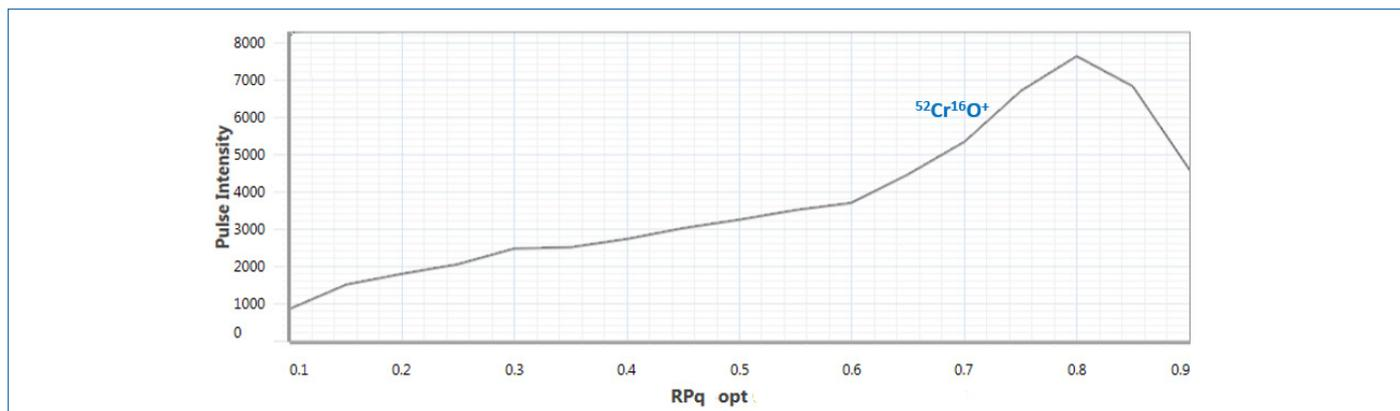


Figure 5: Optimization of Rpq parameter for CrO⁺.

Reactions with Ammonia (NH₃)

Ammonia is another reaction gas that can be very useful in moving some reactive elements to higher masses and measuring them in Mass Shift mode as clusters or products escaping spectral interferences residing on their original masses. Figure 6 shows a periodic table with elements where different shades of green indicate their reactivity with NH₃.

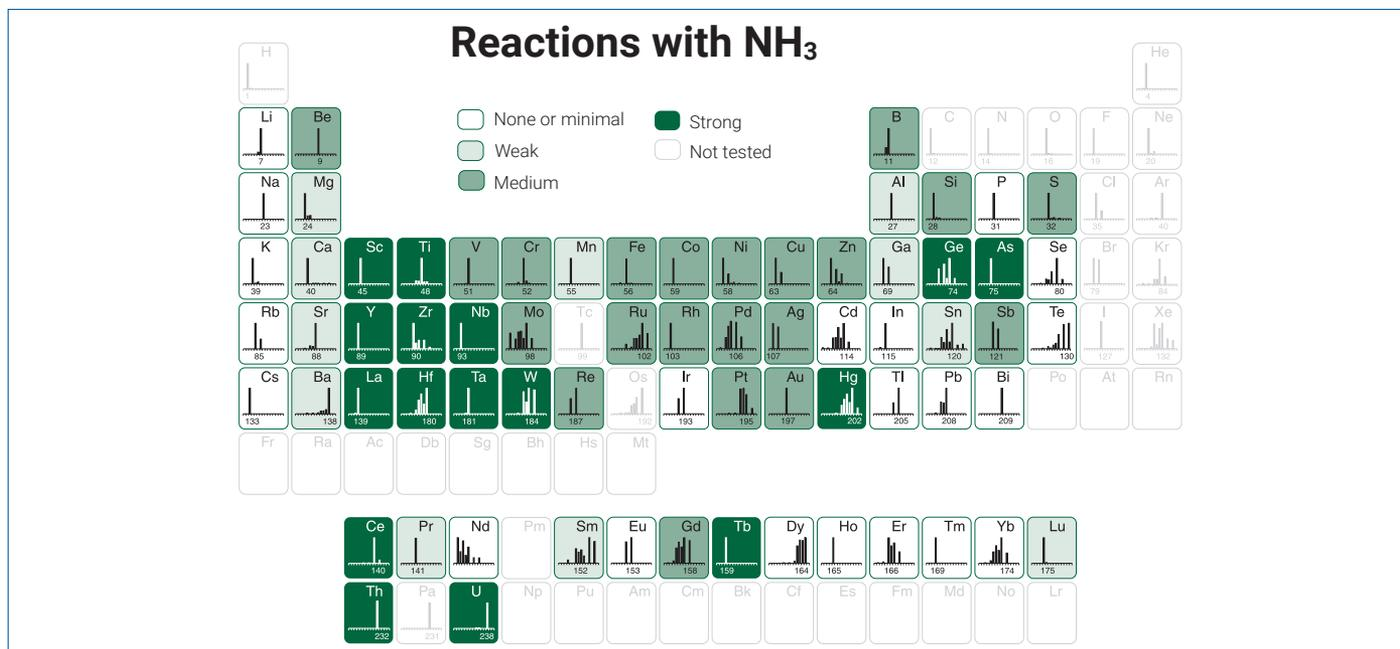


Figure 6: Elements' reactivity with NH₃.

A few elements that react strongly with NH_3 create multiple clusters with NH_3 , and the optimization of the NH_3 flow can change sensitivity of individual clusters quite dramatically. Germanium (Ge) is a good example; it creates clusters with ammonia at masses 90, 107, 124 and 141. The product ion scan run with 0.8 mL/min of NH_3 shows the highest sensitivity cluster at mass 107 (Figure 7). However, when the NH_3 flow was increased to 1.2 mL/min, the cluster at mass 124 became the most prominent (Figure 8). Such flexibility in the selection clusters can be very useful when analyzing trace elements in complex matrices.

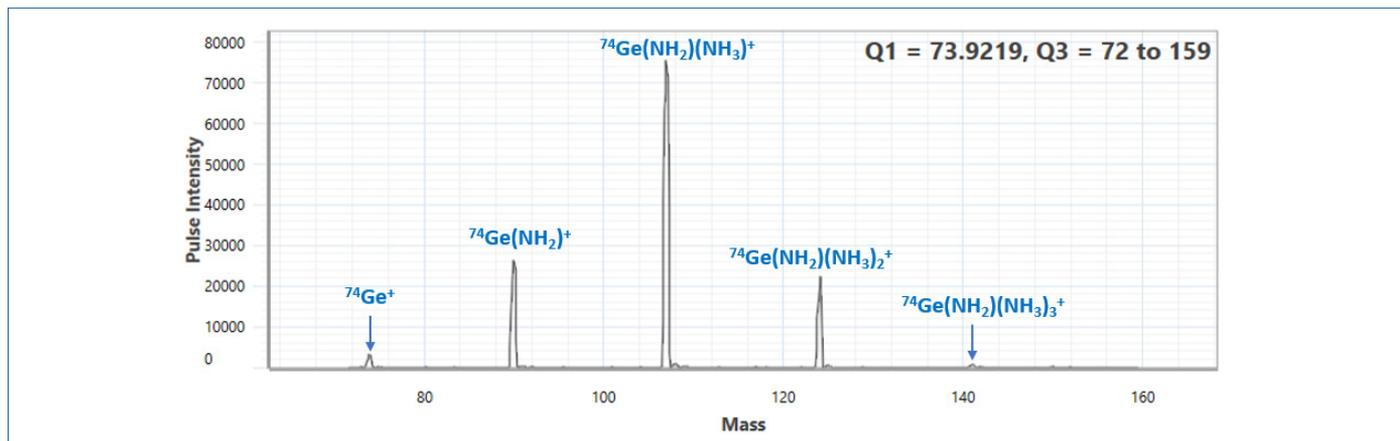


Figure 7: Ge product ion scan with 0.8 mL/min NH_3 flow.

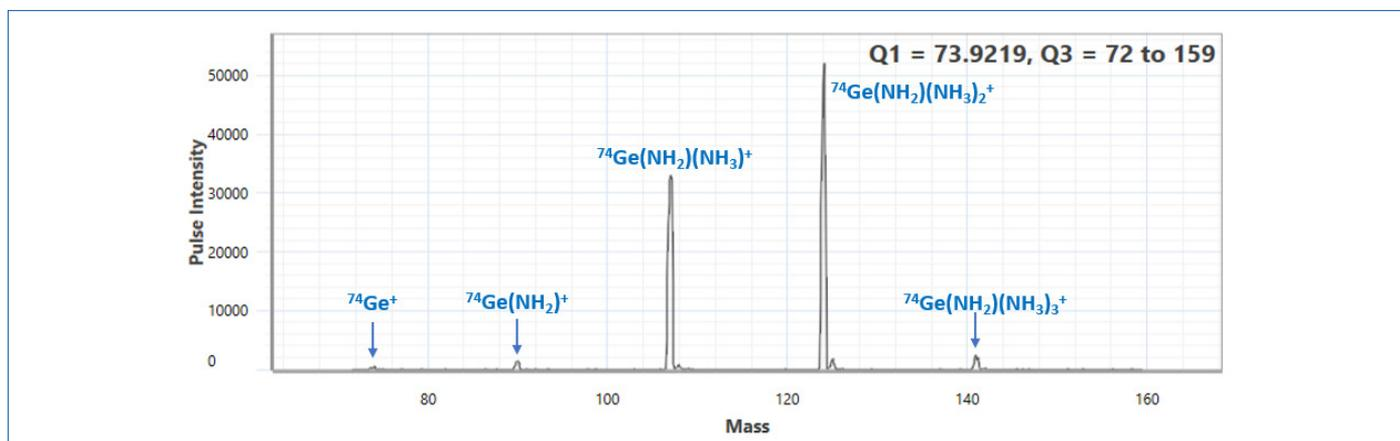


Figure 8: Ge product ion scan with 1.2 mL/min NH_3 flow.

Reactions with Methane (CH_4)

Methane is less powerful for clustering individual elements than O_2 or NH_3 , however, in certain situations, it can be useful as well. One of the CH_4 characteristics is strong reactions with some ions that create spectral interferences such as Ar^+ , ArAr^+ or ArO^+ . In such cases, instead of Mass Shift, MS/MS would be a preferable mode of analysis.

Figure 9 shows a periodic table with elements where different shades of orange indicate their reactivity with CH_4 .

A few elements, such as As, Zr, Ta, W, Ir and Pt, react relatively strongly with CH_4 and their Mass Shift results could be a confirmation of results obtained in Mass Shift mode with a different gas. For example, As is historically measured as AsO^+ at mass 91 (Figure 10), but due to a strong reaction with CH_4 , it can also be measured as AsCH_2^+ at mass 89 (Figure 11).

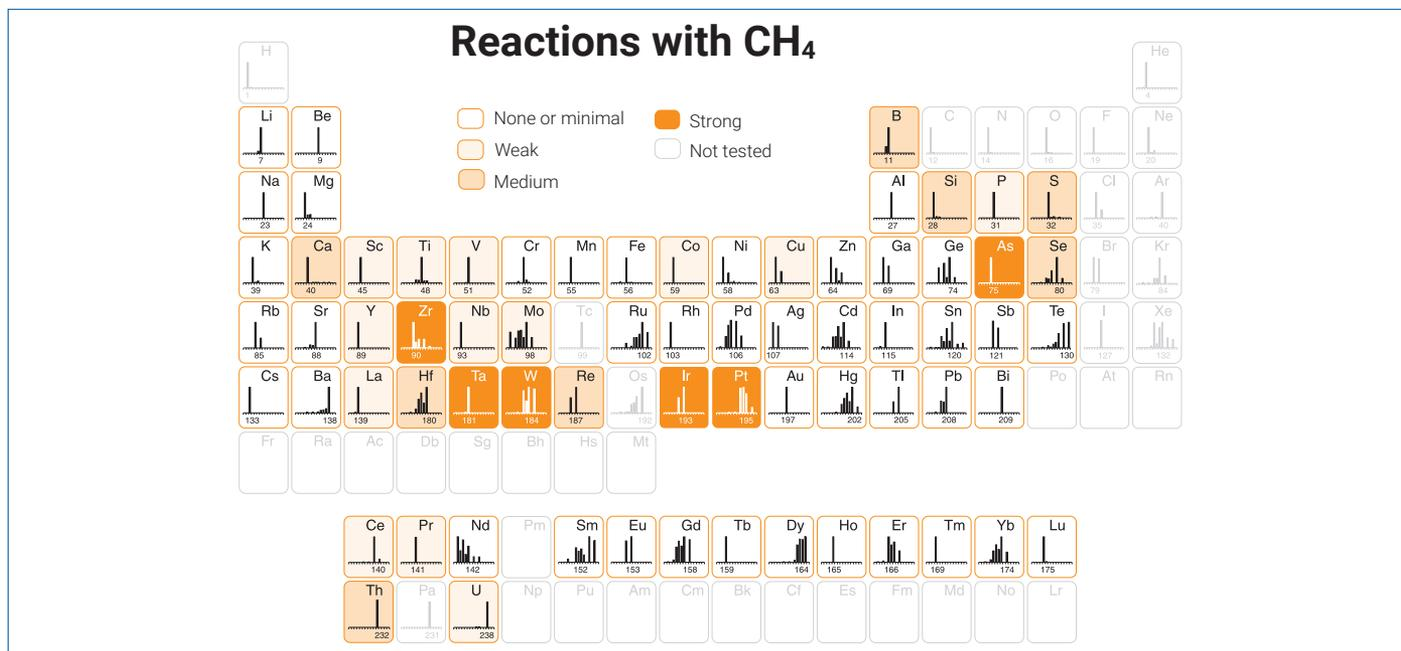


Figure 9: Elements' reactivity with CH₄.

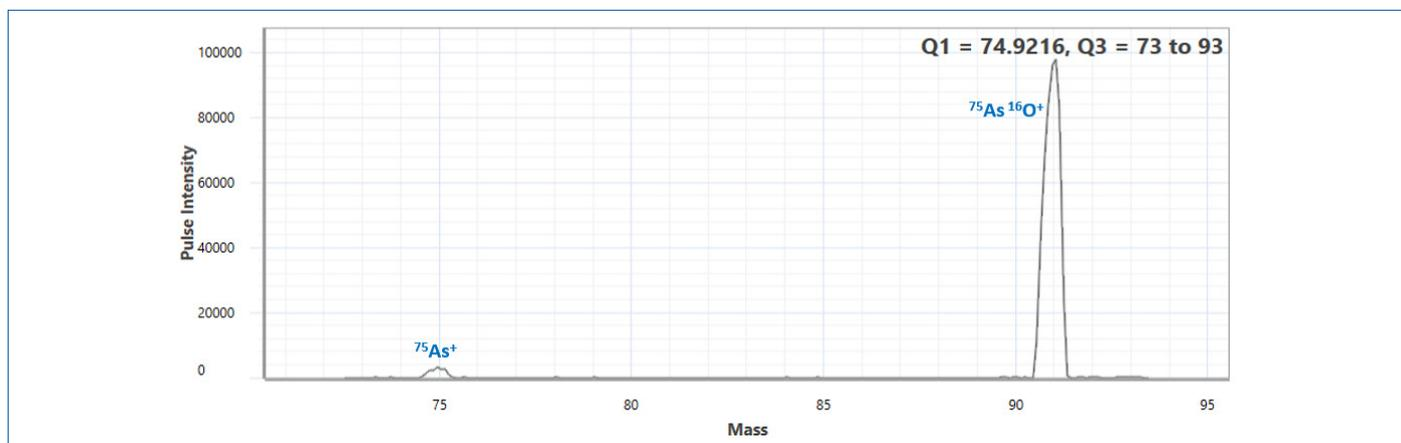


Figure 10: As product ion scan with O₂.

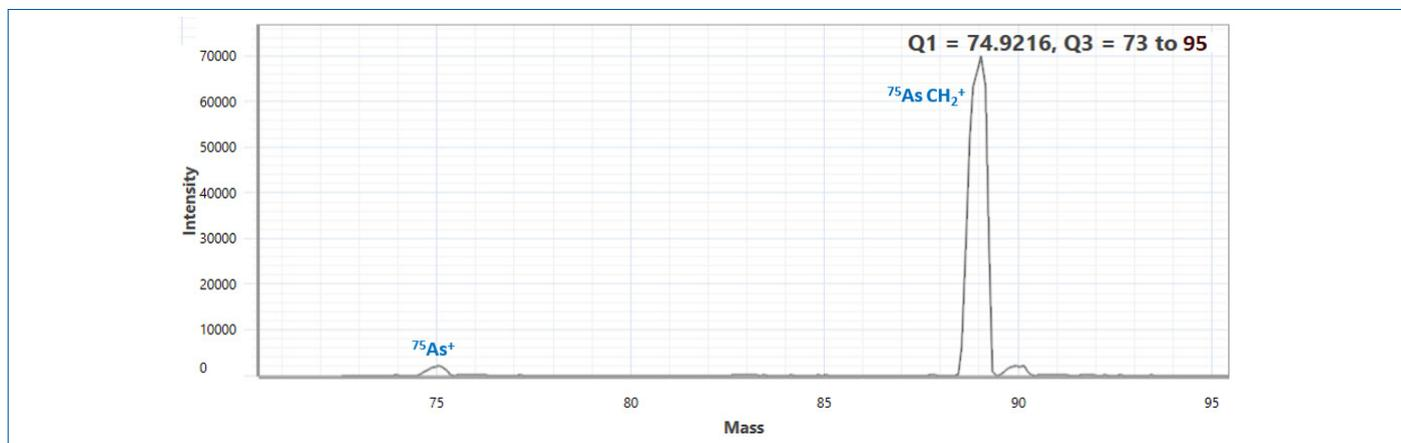


Figure 11: As product ion scan with CH₄.

Conclusion

The NexION 5000 ICP-MS with multi-quadrupole technology and Universal Cell is able to take full advantage of element reactivity with 100% pure gases by analyzing them as cluster ions at higher masses where no interferences reside, and the background is clean. The UCT, thanks to its quadrupole design in combination with dynamic bandpass tuning, provides the unique ability of controlling desirable reactions, promoting some weak reactions, rejecting interferences, and preventing side reactions from taking place.⁵ Periodic tables with color-coded elements indicating their reactivity with a specific gas are a useful tool for method development in complicated matrices, effectively predicting spectral interferences and providing a way to prevent them.

Author

Ewa Pruszkowski
PerkinElmer, Inc.
Shelton, CT USA

Reference

1. "NexION 5000 Multi-Quadrupole ICP-MS", PerkinElmer Product Note, 2020.
2. Pruszkowski E., "Interferences in ICP-MS: Do we still have to worry about them?", PerkinElmer Technical note, 2021.
3. Tanner S.D. et al., "Theory, Design, and Operation of a Dynamic Reaction Cell for ICP-MS", Atomic Spectroscopy, 20, (2), 45-52, 1999.
4. Bandura D.R. et al., "Reaction chemistry and collisional processes in multipole devices for resolving isobaric interferences in ICP-MS", Fresenius J Anal Chem, 2001.
5. Badiei H. et al., "Multimode Cells and Methods of Using Them", United States Patent 8,426,804 B2, Apr.23, 2013.
6. Anicich V.G., "An Index of the Literature for Bimolecular Gas Phase Cation-Molecule Reaction Kinetics", JPL Publication 03-19.
7. "Interactive Tables of Reaction Rates", Ion Chemistry Laboratory, York University, Toronto, Canada, http://www.yorku.ca/dkbohme/research/selection_table.html.