

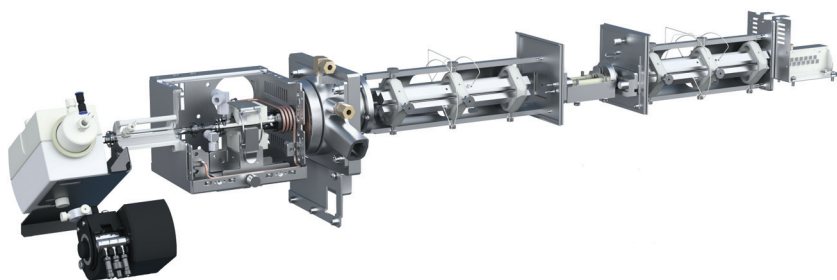
Reaction data for 70 elements using O₂, NH₃ and H₂ gases with the Agilent 8800 Triple Quadrupole ICP-MS

Technical note

Authors

Naoki Sugiyama and
Kazumi Nakano

Agilent Technologies, Japan



Introduction

Currently, methods based on Collision/Reaction Cell (CRC) technology are the most commonly used approach to resolve spectral interferences in ICP-MS. Collision mode uses a chemically inert gas, usually helium (He), to discriminate an analyte ion from any interfering polyatomic ions on the basis of Kinetic Energy Discrimination (KED). KED is effective for all polyatomic ion interferences and rarely requires optimization or customization for particular interferences or sample types, so collision mode using He is the most widely applicable approach for resolving interferences in elemental analysis by ICP-MS. The alternative approach is reaction mode, which relies on a chemical reaction taking place between the interfering ion (or the analyte) and the reactive cell gas in the CRC to resolve a specific interference.



Agilent Technologies

Various reactive gases are used in reaction cell mode including hydrogen (H_2), oxygen (O_2), ammonia (NH_3), methane (CH_4), nitrous oxide (N_2O), carbon dioxide (CO_2) etc, with the choice of gas depending on the individual interference that needs to be addressed. As reactions occur very quickly in the cell (often proceeding with the first collision between the ion and reaction gas molecule), a reaction gas may be more effective than a collision gas for the removal of a given interference. As a result, reaction mode is often used to remove the most intense plasma-based interferences, such as $^{40}Ar^+$ on $^{40}Ca^+$, $^{40}Ar^{16}O^+$ on $^{56}Fe^+$, and $^{40}Ar_2^+$ on $^{80}Se^+$, when ultra-trace level analysis is required, for example in high purity semiconductor reagents. However, conventional quadrupole ICP-MS (ICP-QMS) has been unable to utilize the full potential of reaction mode for more complex or variable samples, because the reaction gas will also react with matrix elements, other analytes, and polyatomic species, resulting in the formation of new and unwanted interferences, which often leads to inaccurate and inconsistent results. For this reason, the application of reactive gases with ICP-QMS has been limited to the determination of a small number of analytes with predictable and consistent interferences in simple, well characterized matrices.

To meet the growing demand from industry for more reliable determination of a wider range of elements at lower concentrations and in a greater variety of matrices, a new type of ICP-MS, a triple quadrupole ICP-MS (ICP-QQQ) has been developed. The Agilent 8800 ICP-QQQ realizes the full potential of reaction chemistry in ICP-MS, by using MS/MS mode to control the reaction chemistry, eliminating the errors and variability of reaction cell operation in ICP-QMS. This enables the Agilent 8800 ICP-QQQ to remove interferences far more effectively, offering significantly improved detection limits and much more consistent reaction mode performance than conventional quadrupole ICP-MS. In MS/MS mode, both quadrupoles are operated as mass filters (unit mass resolution), with the first quadrupole (Q1, positioned in front of the cell) controlling which ions enter the CRC and take part in the reactions. This ensures that the chemical reactions

in the cell are consistent and predictable even when the sample composition is variable, allowing the operator to use reaction chemistry to achieve much more reliable results and lower detection limits than is possible when the same reaction chemistry is used with ICP-QMS.

The MS/MS configuration of ICP-QQQ also simplifies reaction mode method development by eliminating the interference- or analyte/isotope-specific customization of cell conditions that is often required when using reactive cell gases with ICP-QMS.

Information available in the literature relating to reaction cell technology [1, 2, 3] and thermochemical reaction data [4, 5] can be used as a reference to aid development of reaction cell methods. In this technical note, specific guidance on the important tuning parameters and some fundamental reaction data is provided to guide method development when using an 8800 ICP-QQQ in reaction mode.

Reaction cell methods

On-mass mode and mass-shift mode

There are two distinct reaction cell modes of operation for the 8800 ICP-QQQ: on-mass mode and mass-shift mode.

On-mass mode detects the analyte ion at its true isotope mass e.g., $^{51}V^+$ would be detected at m/z 51. This method is used when the analyte ion is relatively unreactive with the chosen cell gas, while the interfering ion reacts efficiently with the cell gas. The interfering ion is therefore removed from the mass of the analyte ion, either by being converted to a neutral species, or by forming a reaction product ion at a new mass away from the analyte ion.

By contrast, mass-shift mode detects the analyte ion as a reaction product ion at the product ion mass. For example, using O_2 as the cell gas, $^{51}V^+$ would be detected as VO^+ at m/z 67. This method is used when the analyte ion reacts efficiently with the cell gas to form a new reaction product ion, while the interfering ion reacts slowly or not at all with the cell gas, so does not contribute significantly to the signal at the

new mass of the analyte product ion. Both methods are illustrated schematically in Figure 1 using the determination of vanadium in a chlorine matrix as an example.

Evaluation of different reaction gases

The three reaction gases that are most commonly used with the Agilent 8800 ICP-QQQ are O_2 , NH_3 (used as a mixture of 10% NH_3 in He), and H_2 . As the characteristics of the three gases differ, the reaction of 70 elements with these gases was investigated.

H_2 reacts with argide ions (ArM^+) at a relatively fast rate compared to its reaction rate with several interfered elements. It is therefore an effective gas to remove argide-interferences such as $^{40}Ar^+$ on ^{40}Ca , $^{38}ArH^+$ on $^{39}K^+$, $^{40}Ar^{12}C^+$ on $^{52}Cr^+$, $^{40}Ar^{40}Ar^+$ on $^{80}Se^+$, allowing measurement of the analyte ions at their original mass using on-mass mode. In addition, some elements that form MH_n^+ product ions can be measured with H_2 cell gas using mass-shift mode. For example phosphorus (P) can be detected as PH_4^+ to avoid the interference from $^{30}SiH^+$ [6] at m/z 31. Likewise, $^{35}Cl^+$ can be measured as $^{35}ClH_2^+$ to avoid interference by $^{16}O^{18}OH^+$ at m/z 35 [7].

O_2 is often used in mass-shift methods to “move” the target analyte from its elemental ion mass to its oxide product ion (MO^+) mass, by setting Q2 to 16 amu higher

than Q1. For example, sulfur (S) analyte ions react with O_2 cell gas, so O_2 mass shift mode can be used to move $^{32}S^+$ to its product ion $^{32}S^{16}O^+$ at m/z 48. Since O_2^+ doesn't react with O_2 cell gas to form O_3^+ , this mass shift mode can be used to avoid the intense interference from O_2^+ on ^{32}S at its original elemental ion mass of m/z 32.

NH_3 (Agilent 8800 ICP-QQQ uses a mix of NH_3 in He) is highly reactive due to its lone electron pair. It is commonly used as a cell gas in both on-mass mode and mass-shift mode depending on the interference to be resolved. Titanium (Ti) suffers a severe interference from S-based polyatomic ions. To avoid these interferences, NH_3 is used to convert Ti^+ to an ammonium cluster product ion, usually $Ti(NH_3)_6^+$. Using this approach, Ti can be measured at low levels in samples with a matrix containing a large amount of S and Ca [8].

Experimental

An Agilent 8800 Triple Quadrupole ICP-MS (option #100) including an Octopole Reaction System (ORS³) cell was used. The standard 8800 ICP-QQQ configuration features Ni interface cones, x-lens and a sample introduction system consisting of a MicroMist glass concentric nebulizer, a Peltier-cooled quartz double-pass Scott-type spray chamber, and a quartz torch with 2.5 mm injector.

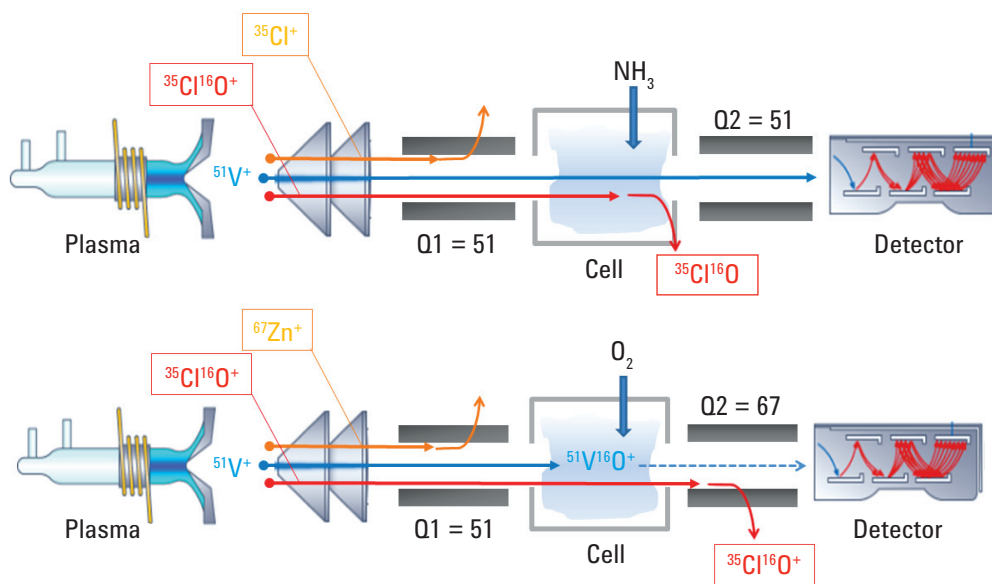


Figure 1. On-mass and mass-shift methods for vanadium measurement in a chlorine matrix: $^{35}Cl^{16}O^+$ interferes with $^{51}V^+$. In on-mass mode (upper), the interfering ion ClO^+ reacts with NH_3 to form neutral ClO (no-charge). ClO is removed while the V^+ analyte ion passes to the detector. In mass-shift mode (lower), the analyte ion V^+ reacts with O_2 to form the product ion VO^+ which is detected at m/z 67. The interfering ion ClO^+ remains at m/z 51 and is rejected by Q2.

The performance of MS/MS mode was investigated for 3 different reaction gases, O₂, NH₃ and H₂. Soft extraction lens tune (Ex1 = 0 V, Ex2 = -180 V, Omega Bias = -80 V and Omega Lens = +10 V) was applied. The preset plasma condition “low matrix” was used throughout the study, providing robust plasma conditions with a CeO⁺/Ce⁺ ratio of < 1 %.

Among the various operational settings of the ICP-QQQ, two key tuning parameters affect reaction efficiency: “octopole bias” and “cell gas flow rate”. Octopole bias determines an ion’s kinetic energy before any reaction occurs in the cell or, if the ion is unreactive with the cell gas, its ion kinetic energy after undergoing a series of collisions in the cell. Cell gas flow rate affects cell gas density, i.e., the probability of an ion experiencing a (reactive) collision with the cell gas, or the number of collisions for an unreactive ion. In this study, these settings were optimized in each cell gas mode as follows: no gas tune was optimized for maximum counts of Co⁺; H₂ tune was optimized for maximum counts of PH⁺, O₂ tune was optimized for maximum counts of SeO⁺, and NH₃ tune was optimized for maximum counts of Ti(NH₂)⁺.

The 8800 ICP-QQQ includes a deflection lens positioned after the ORS³. An ion’s trajectory within the deflection lens is determined by the kinetic energy of the ion and

Table 1. 8800 ICP-QQQ tuning parameters for each gas mode

	No gas mode	H ₂ mode	O ₂ mode	NH ₃ mode
Cell exit (V)	-90			
Deflect (V)	20	-10	-4	3
Plate bias (V)	-100			
Cell gas	N/A	H ₂	O ₂	NH ₃ /He
Cell gas flow rate (mL/min)	N/A	5.0	0.5	1.5
Octopole bias (V)	-8	-25	-18	-8
Octopole RF (V)	180			
KED (V)	5	-8		

the voltage applied to the lens. Changes in octopole bias and/or cell gas flow rate alter the kinetic energy of the ion at the deflection lens, so the “deflect” voltage was tuned to maximize ion transmission for each set of cell conditions. Tuning parameters for each of the three gas modes are summarized in Table 1.

Plasma conditions were tuned for robustness. Theoretically, the plasma conditions (which determine the plasma temperature) could affect reaction efficiency since each ionic species may exist in multiple electronic states (ground state and excited states) in the plasma, and ions in different electronic states are known to have different reactivity [9, 10]. However, it is unlikely that small changes in plasma conditions would have a significant effect on reaction efficiency as long as the plasma is operated consistently in “hot plasma conditions” as defined by Ce⁺/CeO⁺ of < a few %.

Sample preparation

Six multi-element solutions were prepared from multi-element standards purchased from SPEX Certiprep (New Jersey, USA) and single standards bought from Kanto Kagaku (Saitama, Japan) as detailed below.

- Solution 1: 1 ppm of each of Li, Be, Na, Mg, Al, K, Ca, V, Cr, Mn, Fe, Co, Ni, Cu, Zn, Ga, As, Se, Rb, Sr, Ag, Cd, Cs, Tl, Pb, Bi, Th and U in 1% HNO₃
- Solution 2: 1 ppm of each of Sc, Y, La, Ce, Pr, Nd, Sm, Eu, Gd, Tb, Dy, Ho, Er, Tm, Yb and Lu in 1% HNO₃
- Solution 3: 1 ppm of each of B, Si, P, S, Ti, Ge, Zr, Nb, Mo, Ta, W and Re in 0.1% HF
- Solution 4: 1 ppm of each of Ru, Rh, Pd, Sn, Sb, Te, Hf, Ir, Pt and Au in 1% HNO₃/0.5% HCl
- Solution 5: 1 ppm of each of Cl, Br and I in 1% TMAH
- Solution 6: 100 ppb Hg in 1% HNO₃/0.5% HCl

Blank solutions matching the acid content of each standard were also prepared for blank subtraction to determine the net sensitivity in each mode. For each element, several mass-pairs (combinations of Q1 set mass and Q2 set mass) were surveyed in order to determine the net sensitivity of the element and its various product ions. So for each analyte mass-pair, Q1 was set to the mass of the natural isotope (the precursor ion), and Q2 was set to either the same mass

(Q1 = Q2) or the mass-shifted m/z of the product ion (Q2 = Q1+ Δm). The common mass-shift settings (the mass difference between Q1 and Q2) associated with the most common reaction product ion transitions (e.g., +16 amu for O₂ reaction mode) are predefined in the ICP-MS MassHunter software, to aid method development.

Reaction data

The sensitivity of 70 elements and their respective product ions was studied in no gas, O₂, NH₃ and H₂ modes and the results are summarized in Table 2A (absolute sensitivity) and Table 2B (sensitivity as % relative to no gas mode).

It should be emphasized that the sensitivity data given here has not been optimized for individual elements; all results were obtained using the tuning parameters given in Table 1. Tuning cell parameters for a target analyte ion would lead to an increase in sensitivity of that ion.

The data given in the tables is designed to provide an insight on the reactivity and product ion formation of a wide range of elements with O₂, NH₃ and H₂. For example, we can see from the data that some elements form not only oxides but also dioxides and trioxides. In certain applications, it may be possible to achieve lower detection limits by measuring higher cluster ions.

Conclusions

This note provides fundamental reaction data that was obtained experimentally using an Agilent 8800 Triple Quadrupole ICP-MS (ICP-QQQ) with ORS³ collision reaction cell. The tandem mass spectrometer layout allows the 8800 ICP-QQQ to use MS/MS mode, where both quadrupoles (Q1 and Q2) are operated as unit mass filters. This allows precise control of reaction chemistry in either on-mass or mass-shift mode, to eliminate problematic spectral interferences that cannot be removed with conventional ICP-QMS. With MS/MS, mass selection is performed by the first quadrupole (Q1), allowing the cell to operate much more efficiently and selectively, and opening up the potential of reaction

cell chemistry in ICP-MS to a wide range of elements. The sensitivity data for 70 elements obtained in different reaction gas modes (oxygen, ammonia and hydrogen) has been investigated in order to provide valuable insight into reaction cell methodology and to aid method development when using the 8800 ICP-QQQ in reaction mode. By referring to the data in the tables, you can see at a glance which elements form potentially useful product ions.

References

1. D.R. Bandura, V.I. Baranov and S.D. Tanner, *Fresenius J. Anal. Chem.*, 2001, 370, 454-470
2. John W. Olesik and Deanna R. Jones, *J. Anal. At. Spectrom.*, 2006, 21, 141-159
3. *Agilent 8800 ICP-QQQ application handbook*, 5991-2802EN
4. V.G. Anicich, *The Astrophysical Journal Supplement Series*, 1993, 84, 215-315
5. York University web site, http://www.chem.yorku.ca/profs/bohme/research/selection_table.html
6. Agilent application note, Improvement of ICP-MS detectability of phosphorus and titanium in high purity silicon samples using the Agilent 8800 Triple Quadrupole ICP-MS, 5991-2466EN
7. Agilent application note, Trace level analysis of sulfur, phosphorus, silicon and chlorine in NMP using the Agilent 8800 Triple Quadrupole ICP-MS, 5991-2303EN
8. Lieve Balcaen and et al, *Anal. Chim. Acta*, 2014, 809, 27, 1-8
9. M. Hamdan and et al., *J. Phys. B: At. Mol. Phys.*, 1984, 17, 849-856
10. P. B. Armentrout, *J. Anal. At. Spectrom.*, 2004, 19, 571-580

Table 2A. Sensitivity of elemental ions and their reaction product ions in no gas mode and three reaction cell modes. The green circle represents sensitivity > 1000 cps/ppb and the yellow circle represents sensitivity > 100 cps/ppb

Element	Atomic Num	Mass Num	No gas (cps/ppb)	H ₂ (cps/ppb)				O ₂ (cps/ppb)			
				M ⁺	MH ⁺	MH ₂ ⁺	MH ₃ ⁺	M ⁺	MO ⁺	MO ₂ ⁺	MO ₃ ⁺
				ΔM	0	1	2	3	0	16	32
Li	3	7	● 41759	● 4521	0	0	0	● 232	0	0	0
Be	4	9	● 10193	● 3898	● 526	1	1	● 136	18	6	0
B	5	11	● 5266	● 2193	5	39	0	● 317	6	1	0
Na	11	23	● 57839	● 39584	0	0	0	● 16182	0	0	0
Mg	12	24	● 39815	● 24945	9	0	0	● 13593	● 1031	8	2
Al	13	27	● 56763	● 34045	1	0	0	● 22139	● 1398	3	22
Si	14	28	● 16421	● 10884	● 5137	22	1	● 3386	● 2482	● 187	7
P	15	31	● 2486	● 1285	● 485	95	6	21	● 902	3	0
S	16	32	● 6181	● 4137	● 631	97	4	● 131	● 2038	2	0
Cl	17	35	● 226	2	6	● 102	0	0	6	0	0
K	19	39	● 43539	● 33850	0	0	0	● 20183	0	0	0
Ca	20	44	● 2142	● 1249	0	0	0	● 987	80	22	0
Sc	21	45	● 95839	● 48343	● 807	0	0	● 1186	● 45969	● 524	16
Ti	22	47	● 5107	● 2940	21	0	0	77	● 2502	49	17
V	23	51	● 74095	● 43074	● 132	0	0	● 2118	● 46960	● 1605	● 170
Cr	24	52	● 64552	● 34482	7	0	0	● 28691	● 11279	● 254	52
Mn	25	55	● 81996	● 33599	40	0	0	● 37512	● 13080	22	24
Fe	26	56	● 66340	● 32190	● 143	0	0	● 29356	● 15154	66	14
Co	27	59	● 69707	● 45629	57	1	0	● 29359	● 16187	59	31
Ni	28	60	● 14472	● 12522	2	0	0	● 7848	● 1877	22	2
Cu	29	63	● 34480	● 27423	0	0	0	● 20842	● 294	35	0
Zn	30	66	● 11313	● 3930	0	0	0	● 6120	● 303	2	1
Ga	31	71	● 38632	● 14743	0	0	0	● 22568	0	0	0
Ge	32	72	● 17998	● 7688	92	0	0	● 7062	● 1719	● 161	1
As	33	75	● 10603	● 6438	● 131	1	0	● 334	● 6013	6	0
Se	34	82*	● 732	● 282	9	0	0	● 212	● 204	0	0
Br	35	81	● 1847	● 799	● 260	8	0	● 116	● 173	1	0
Rb	37	85	● 76018	● 32869	0	0	0	● 43163	0	0	0
Sr	38	88	● 95671	● 40573	15	0	0	● 52149	● 5582	● 1471	0
Y	39	89	● 150007	● 81923	98	0	0	● 978	● 77354	● 1877	11
Zr	40	90	● 58386	● 31671	35	0	0	● 188	● 26699	● 5106	● 1432
Nb	41	93	● 102753	● 76003	57	0	0	● 557	● 4744	● 48482	14
Mo	42	95	● 17028	● 10677	3	0	0	● 752	● 956	● 6125	6
Ru	44	101	● 16832	● 10171	1	0	0	● 4789	● 3427	● 121	46
Rh	45	103	● 85372	● 47902	5	0	0	● 32214	● 7102	40	6
Pd	46	105	● 15176	● 11137	1	0	0	● 8697	● 161	7	0
Ag	47	107	● 42300	● 24631	0	0	0	● 29437	1	1	0
Cd	48	111	● 10075	● 3724	0	0	0	● 6238	55	0	0
Sn	50	118	● 28708	● 8585	2	0	0	● 10789	● 1496	88	0
Sb	51	121	● 34183	● 12507	4	0	0	● 8256	● 8954	31	0
Te	52	125	● 1413	● 554	0	0	0	● 394	● 324	0	0
I	53	127	● 20843	● 10059	5	0	0	● 9338	● 2819	2	0

*⁸²Se was used in order to allow measurement of the no gas mode sensitivity

Table 2A. continued showing H₂ and O₂

Element	Atomic Num	Mass Num	No gas (cps/ppb)		H ₂ (cps/ppb)				O ₂ (cps/ppb)			
					M ⁺	MH ⁺	MH ₂ ⁺	MH ₃ ⁺	M ⁺	MO ⁺	MO ₂ ⁺	MO ₃ ⁺
					ΔM	0	1	2	3	0	16	32
Cs	55	133	● 100507	● 45313	0	0	0	● 72832	0	0	0	
La	57	139	● 132424	● 98130	24	0	0	● 320	● 87357	● 2007	0	
Ce	58	140	● 130831	● 53284	18	0	0	● 115	● 80926	● 1311	72	
Pr	59	141	● 181014	● 47241	18	0	0	● 139	● 88626	● 767	0	
Nd	60	146	● 26895	● 9859	3	0	0	30	● 16631	● 175	0	
Sm	62	147	● 21771	● 8404	2	0	0	● 315	● 14053	● 305	0	
Eu	63	153	● 85415	● 28133	7	0	0	● 24774	● 19372	● 1969	0	
Gd	64	157	● 28661	● 9333	4	0	0	● 1282	● 14006	● 362	1	
Tb	65	159	● 164457	● 56971	17	0	0	● 511	● 95468	● 1918	5	
Dy	66	163	● 36080	● 14809	6	0	0	● 175	● 28535	● 420	1	
Ho	67	165	● 181582	● 56523	27	0	0	● 680	● 99092	● 2435	11	
Er	68	166	● 50501	● 18679	9	0	0	● 296	● 34575	● 982	7	
Tm	69	169	● 172859	● 51606	73	0	0	● 8973	● 100996	● 2411	17	
Yb	70	172	● 34696	● 11596	30	0	0	● 10624	● 9081	● 415	1	
Lu	71	175	● 153754	● 55401	11	0	0	● 1994	● 104484	● 3932	32	
Hf	72	178	● 36060	● 13109	6	0	0	15	● 13849	● 7833	0	
Ta	73	181	● 150282	● 63258	48	0	0	63	● 5121	● 75315	● 235	
W	74	182	● 36381	● 17171	20	0	0	60	● 2035	● 17187	● 266	
Re	75	185	● 48114	● 19353	36	0	0	● 3173	● 4413	● 10111	● 5323	
Ir	77	193	● 50400	● 41464	57	3	0	● 17592	● 12528	● 664	● 105	
Pt	78	195	● 18826	● 17165	12	3	0	● 10722	● 1409	11	1	
Au	79	197	● 32902	● 31730	4	3	0	● 18775	● 114	4	0	
Hg	80	202	● 8205	● 3859	0	0	0	● 4710	4	0	0	
Tl	81	205	● 67878	● 25667	0	0	0	● 42998	0	0	0	
Pb	82	208	● 46191	● 18386	0	0	0	● 26069	● 589	0	0	
Bi	83	209	● 77596	● 29212	1	0	0	● 41364	● 3776	0	0	
Th	90	232	● 87948	● 31544	22	2	0	2	● 16035	NA	NA	
U	92	238	● 90868	● 32366	14	0	0	2	● 559	NA	NA	

Table 2A. continued showing NH₃

Element	Atomic Num	Mass Num	NH ₃ (cps/ppb)																			
			M+	M(NH) ⁺	M(NH ₂) ⁺	M(NH ₃) ⁺	MNH(NH ₃) ⁺	MNH ₂ (NH ₃) ⁺	M(NH ₃) ₂ ⁺	MNH(NH ₃) ₂ ⁺	MNH ₂ (NH ₃) ₂ ⁺	M(NH ₃) ₃ ⁺										
			0	15	16	17	32	33	34	49	50	51										
Li	3	7	●	1543	0	0	2	0	0	1	0	0	1									
Be	4	9	●	366	0	38	3	0	●	116	16	0	●	190	●	249						
B	5	11	●	131	0	12	2	51	1	0	●	258	3	0								
Na	11	23	●	40985	0	0	80	0	0	15	0	0	0									
Mg	12	24	●	28157	0	0	●	178	0	0	55	1	0	36								
Al	13	27	●	35328	0	0	80	0	2	3	1	0	0									
Si	14	28	●	1511	0	●	438	3	1	0	0	1	0	0								
P	15	31		9	1	2	0	0	0	0	0	0	0	0								
S	16	32		0	0	0	0	0	0	0	0	0	0	0								
Cl	17	35		16	0	0	0	0	0	0	0	0	0	0								
K	19	39	●	36532	0	0	21	0	0	1	0	0	0	0								
Ca	20	44	●	1562	0	1	5	1	0	1	0	0	0	0								
Sc	21	45	●	18843	●	7197	●	2139	64	●	2167	●	372	47	●	1256	●	406	●	115		
Ti	22	47	●	1148	●	469	78	1	100	13	0	84	17	2								
V	23	51	●	39430	●	831	●	4927	●	185	●	301	●	362	124	●	338	●	668	●	130	
Cr	24	52	●	48910	8	9	●	182	3	1	●	405	4	3	14							
Mn	25	55	●	64783	0	4	●	278	0	1	●	187	0	1	26							
Fe	26	56	●	50691	2	27	●	280	1	8	●	2474	2	7	●	254						
Co	27	59	●	48441	3	20	●	493	2	5	●	3499	2	4	●	800						
Ni	28	60	●	10516	0	1	96	0	0	●	317	0	0	●	566							
Cu	29	63	●	25533	0	1	●	210	0	0	●	1856	0	0	17							
Zn	30	66	●	9000	0	0	●	119	0	0	61	0	0	12								
Ga	31	71	●	30730	0	0	68	0	0	11	0	0	0	0								
Ge	32	72	●	6211	0	●	2571	25	0	●	1418	73	0	●	185	●	588					
As	33	75	●	2569	8	●	1599	59	1	●	230	32	1	16	4							
Se	34	82*		347	0	0	0	0	0	0	0	0	0	0	0							
Br	35	81		1	0	0	0	0	0	0	0	0	0	0	0							
Rb	37	85	●	54979	1	0	0	0	0	0	0	0	0	0	0							
Sr	38	88	●	79654	0	48	●	147	15	2	16	3	1	11								
Y	39	89	●	47349	●	22301	●	5740	●	587	●	5828	●	974	●	154	●	1903	●	672	●	187
Zr	40	90	●	11579	●	10060	●	790	14	●	2172	●	155	6	●	1433	●	255	29			
Nb	41	93	●	19857	●	6968	●	206	5	28	1	0	22	1	0							
Mo	42	95	●	10344	23	12	31	0	0	●	143	0	0	1								
Ru	44	101	●	11426	1	2	47	0	0	●	137	0	0	10								
Rh	45	103	●	55102	0	3	●	226	0	1	●	548	0	0	27							
Pd	46	105	●	11381	0	0	49	0	0	●	127	0	0	130								
Ag	47	107	●	37325	0	0	81	0	0	●	247	0	0	3								
Cd	48	111	●	8290	0	0	21	0	0	6	0	0	0	1								
Sn	50	118	●	16684	0	5	76	0	1	69	0	0	0	3								
Sb	51	121	●	24016	3	13	●	1147	1	3	●	179	1	1	4							
Te	52	125	●	1096	0	0	0	0	0	0	0	0	0	0								
I	53	127		5	0	0	0	0	0	0	0	0	0	0								

⁸²Se was used in order to allow measurement of the no gas mode sensitivity

Table 2A. continued showing NH₃

Element	Atomic Num	Mass Num	NH ₃ (cps/ppb)										
			M+	M(NH) ⁺	M(NH ₂) ⁺	M(NH ₃) ⁺	MNH(NH ₃) ⁺	MNH ₂ (NH ₃) ⁺	M(NH ₃) ₂ ⁺	MNH(NH ₃) ₂ ⁺	MNH ₂ (NH ₃) ₂ ⁺	M(NH ₃) ₃ ⁺	
			0	15	16	17	32	33	34	49	50	51	
Cs	55	133	● 80994	0	0	3	0	0	0	0	0	0	0
La	57	139	● 51900	● 31021	● 12190	● 868	● 1145	● 284	96	● 245	● 123	76	
Ce	58	140	● 47439	● 22882	● 8447	● 521	● 915	● 328	68	● 238	● 130	71	
Pr	59	141	● 96491	● 3154	● 7649	● 505	● 107	● 158	47	26	47	39	
Nd	60	146	● 18452	● 190	● 1053	62	10	22	8	2	8	7	
Sm	62	147	● 14593	1	● 513	32	1	16	7	0	7	7	
Eu	63	153	● 59006	0	● 150	79	4	3	8	1	0	7	
Gd	64	157	● 11875	● 4066	● 840	56	● 589	77	13	● 160	44	12	
Tb	65	159	● 78912	● 13589	● 4297	● 160	● 3202	● 498	67	● 941	● 318	77	
Dy	66	163	● 28299	● 121	● 635	33	11	42	8	3	18	7	
Ho	67	165	● 112235	● 826	● 2120	● 133	93	262	35	20	98	31	
Er	68	166	● 32776	● 367	● 730	38	45	89	10	9	32	10	
Tm	69	169	● 113722	27	● 749	● 103	4	64	20	1	16	17	
Yb	70	172	● 22193	0	24	22	0	1	3	0	0	3	
Lu	71	175	● 115121	● 1507	● 3205	● 259	● 359	● 449	56	85	● 153	47	
Hf	72	178	● 6021	● 13915	● 703	15	● 2018	99	2	● 470	54	4	
Ta	73	181	● 23148	● 29292	● 2691	32	● 230	16	1	● 117	5	0	
W	74	182	● 9277	● 6208	● 315	3	13	1	0	1	0	0	
Re	75	185	● 39126	● 289	19	2	1	0	0	0	0	0	
Ir	77	193	● 25337	● 1931	20	● 143	89	22	31	4	5	1	
Pt	78	195	● 15021	7	4	82	2	2	● 200	1	2	● 121	
Au	79	197	● 23788	5	1	● 112	2	0	● 595	1	0	7	
Hg	80	202	2	0	0	0	0	0	0	0	0	0	
Tl	81	205	● 52670	0	0	2	0	0	0	0	0	0	
Pb	82	208	● 35788	0	1	5	0	0	1	0	0	0	
Bi	83	209	● 60593	2	4	11	0	0	1	0	0	0	
Th	90	232	● 15603	● 32219	● 6181	● 229	NA	NA	NA	NA	NA	NA	
U	92	238	● 29649	● 24271	● 4917	● 248	NA	NA	NA	NA	NA	NA	

Table 2B. Sensitivity (relative to no gas mode sensitivity) of elemental ions and their respective product ions in three reaction cell modes. The green circle represents sensitivity > 2 % and the yellow circle represents sensitivity > 0.5%.

Element	Atomic Num	Mass Num	No gas (cps/ppb)	H ₂ (%)				O ₂ (%)				
				M ⁺	MH ⁺	MH ₂ ⁺	MH ₃ ⁺	M ⁺	MO ⁺	MO ₂ ⁺	MO ₃ ⁺	
				ΔM	0	1	2	3	0	16	32	48
Li	3	7	41759	● 10.8%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 0.6%	● 0.0%	● 0.0%	● 0.0%
Be	4	9	10193	● 38.2%	● 5.2%	● 0.0%	● 0.0%	● 0.0%	● 1.3%	● 0.2%	● 0.1%	● 0.0%
B	5	11	5266	● 41.6%	● 0.1%	● 0.7%	● 0.0%	● 0.0%	● 6.0%	● 0.1%	● 0.0%	● 0.0%
Na	11	23	57839	● 68.4%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 28.0%	● 0.0%	● 0.0%	● 0.0%
Mg	12	24	39815	● 62.7%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 34.1%	● 2.6%	● 0.0%	● 0.0%
Al	13	27	56763	● 60.0%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 39.0%	● 2.5%	● 0.0%	● 0.0%
Si	14	28	16421	● 66.3%	● 31.3%	● 0.1%	● 0.0%	● 0.0%	● 20.6%	● 15.1%	● 1.1%	● 0.0%
P	15	31	2486	● 51.7%	● 19.5%	● 3.8%	● 0.2%	● 0.0%	● 0.8%	● 36.3%	● 0.1%	● 0.0%
S	16	32	6181	● 66.9%	● 10.2%	● 1.6%	● 0.1%	● 0.0%	● 2.1%	● 33.0%	● 0.0%	● 0.0%
Cl	17	35	226	● 1.0%	● 2.8%	● 45.3%	● 0.0%	● 0.0%	● 0.1%	● 2.6%	● 0.0%	● 0.0%
K	19	39	43539	● 77.7%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 46.4%	● 0.0%	● 0.0%	● 0.0%
Ca	20	44	2142	● 58.3%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 46.1%	● 3.7%	● 1.0%	● 0.0%
Sc	21	45	95839	● 50.4%	● 0.8%	● 0.0%	● 0.0%	● 0.0%	● 1.2%	● 48.0%	● 0.5%	● 0.0%
Ti	22	47	5107	● 57.6%	● 0.4%	● 0.0%	● 0.0%	● 0.0%	● 1.5%	● 49.0%	● 1.0%	● 0.3%
V	23	51	74095	● 58.1%	● 0.2%	● 0.0%	● 0.0%	● 0.0%	● 2.9%	● 63.4%	● 2.2%	● 0.2%
Cr	24	52	64552	● 53.4%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 44.4%	● 17.5%	● 0.4%	● 0.1%
Mn	25	55	81996	● 41.0%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 45.7%	● 16.0%	● 0.0%	● 0.0%
Fe	26	56	66340	● 48.5%	● 0.2%	● 0.0%	● 0.0%	● 0.0%	● 44.3%	● 22.8%	● 0.1%	● 0.0%
Co	27	59	69707	● 65.5%	● 0.1%	● 0.0%	● 0.0%	● 0.0%	● 42.1%	● 23.2%	● 0.1%	● 0.0%
Ni	28	60	14472	● 86.5%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 54.2%	● 13.0%	● 0.1%	● 0.0%
Cu	29	63	34480	● 79.5%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 60.4%	● 0.9%	● 0.1%	● 0.0%
Zn	30	66	11313	● 34.7%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 54.1%	● 2.7%	● 0.0%	● 0.0%
Ga	31	71	38632	● 38.2%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 58.4%	● 0.0%	● 0.0%	● 0.0%
Ge	32	72	17998	● 42.7%	● 0.5%	● 0.0%	● 0.0%	● 0.0%	● 39.2%	● 9.5%	● 0.9%	● 0.0%
As	33	75	10603	● 60.7%	● 1.2%	● 0.0%	● 0.0%	● 0.0%	● 3.2%	● 56.7%	● 0.1%	● 0.0%
Se	34	82*	732	● 38.5%	● 1.3%	● 0.0%	● 0.0%	● 0.0%	● 29.0%	● 27.9%	● 0.1%	● 0.0%
Br	35	81	1847	● 43.3%	● 14.1%	● 0.4%	● 0.0%	● 0.0%	● 6.3%	● 9.4%	● 0.1%	● 0.0%
Rb	37	85	76018	● 43.2%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 56.8%	● 0.0%	● 0.0%	● 0.0%
Sr	38	88	95671	● 42.4%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 54.5%	● 5.8%	● 1.5%	● 0.0%
Y	39	89	150007	● 54.6%	● 0.1%	● 0.0%	● 0.0%	● 0.0%	● 0.7%	● 51.6%	● 1.3%	● 0.0%
Zr	40	90	58386	● 54.2%	● 0.1%	● 0.0%	● 0.0%	● 0.0%	● 0.3%	● 45.7%	● 8.7%	● 2.5%
Nb	41	93	102753	● 74.0%	● 0.1%	● 0.0%	● 0.0%	● 0.0%	● 0.5%	● 4.6%	● 47.2%	● 0.0%
Mo	42	95	17028	● 62.7%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 4.4%	● 5.6%	● 36.0%	● 0.0%
Ru	44	101	16832	● 60.4%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 28.4%	● 20.4%	● 0.7%	● 0.3%
Rh	45	103	85372	● 56.1%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 37.7%	● 8.3%	● 0.0%	● 0.0%
Pd	46	105	15176	● 73.4%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 57.3%	● 1.1%	● 0.0%	● 0.0%
Ag	47	107	42300	● 58.2%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 69.6%	● 0.0%	● 0.0%	● 0.0%
Cd	48	111	10075	● 37.0%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 61.9%	● 0.5%	● 0.0%	● 0.0%
Sn	50	118	28708	● 29.9%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 37.6%	● 5.2%	● 0.3%	● 0.0%
Sb	51	121	34183	● 36.6%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 24.2%	● 26.2%	● 0.1%	● 0.0%
Te	52	125	1413	● 39.2%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 27.9%	● 22.9%	● 0.0%	● 0.0%
I	53	127	20843	● 48.3%	● 0.0%	● 0.0%	● 0.0%	● 0.0%	● 44.8%	● 13.5%	● 0.0%	● 0.0%

*⁸²Se was used in order to allow measurement of the no gas mode sensitivity

Table 2B. continued showing H₂ and O₂

Element	Atomic Num	Mass Num	No gas (cps/ppb)	H ₂ (%)				O ₂ (%)			
				M ⁺	MH ⁺	MH ₂ ⁺	MH ₃ ⁺	M ⁺	MO ⁺	MO ₂ ⁺	MO ₃ ⁺
				ΔM	0	1	2	3	0	16	32
Cs	55	133	100507	● 45.1%	0.0%	0.0%	0.0%	● 72.5%	0.0%	0.0%	0.0%
La	57	139	132424	● 74.1%	0.0%	0.0%	0.0%	0.2%	● 66.0%	● 1.5%	0.0%
Ce	58	140	130831	● 40.7%	0.0%	0.0%	0.0%	0.1%	● 61.9%	● 1.0%	0.1%
Pr	59	141	181014	● 26.1%	0.0%	0.0%	0.0%	0.1%	● 49.0%	0.4%	0.0%
Nd	60	146	26895	● 36.7%	0.0%	0.0%	0.0%	0.1%	● 61.8%	● 0.6%	0.0%
Sm	62	147	21771	● 38.6%	0.0%	0.0%	0.0%	● 1.4%	● 64.5%	● 1.4%	0.0%
Eu	63	153	85415	● 32.9%	0.0%	0.0%	0.0%	● 29.0%	● 22.7%	● 2.3%	0.0%
Gd	64	157	28661	● 32.6%	0.0%	0.0%	0.0%	● 4.5%	● 48.9%	● 1.3%	0.0%
Tb	65	159	164457	● 34.6%	0.0%	0.0%	0.0%	0.3%	● 58.1%	● 1.2%	0.0%
Dy	66	163	36080	● 41.0%	0.0%	0.0%	0.0%	● 0.5%	● 79.1%	● 1.2%	0.0%
Ho	67	165	181582	● 31.1%	0.0%	0.0%	0.0%	0.4%	● 54.6%	● 1.3%	0.0%
Er	68	166	50501	● 37.0%	0.0%	0.0%	0.0%	● 0.6%	● 68.5%	● 1.9%	0.0%
Tm	69	169	172859	● 29.9%	0.0%	0.0%	0.0%	● 5.2%	● 58.4%	● 1.4%	0.0%
Yb	70	172	34696	● 33.4%	0.1%	0.0%	0.0%	● 30.6%	● 26.2%	● 1.2%	0.0%
Lu	71	175	153754	● 36.0%	0.0%	0.0%	0.0%	● 1.3%	● 68.0%	● 2.6%	0.0%
Hf	72	178	36060	● 36.4%	0.0%	0.0%	0.0%	0.0%	● 38.4%	● 21.7%	0.0%
Ta	73	181	150282	● 42.1%	0.0%	0.0%	0.0%	0.0%	● 3.4%	● 50.1%	0.2%
W	74	182	36381	● 47.2%	0.1%	0.0%	0.0%	0.2%	● 5.6%	● 47.2%	● 0.7%
Re	75	185	48114	● 40.2%	0.1%	0.0%	0.0%	● 6.6%	● 9.2%	● 21.0%	● 11.1%
Ir	77	193	50400	● 82.3%	0.1%	0.0%	0.0%	● 34.9%	● 24.9%	● 1.3%	0.2%
Pt	78	195	18826	● 91.2%	0.1%	0.0%	0.0%	● 57.0%	● 7.5%	0.1%	0.0%
Au	79	197	32902	● 96.4%	0.0%	0.0%	0.0%	● 57.1%	0.3%	0.0%	0.0%
Hg	80	202	8205	● 47.0%	0.0%	0.0%	0.0%	● 57.4%	0.0%	0.0%	0.0%
Tl	81	205	67878	● 37.8%	0.0%	0.0%	0.0%	● 63.3%	0.0%	0.0%	0.0%
Pb	82	208	46191	● 39.8%	0.0%	0.0%	0.0%	● 56.4%	● 1.3%	0.0%	0.0%
Bi	83	209	77596	● 37.6%	0.0%	0.0%	0.0%	● 53.3%	● 4.9%	0.0%	0.0%
Th	90	232	87948	● 35.9%	0.0%	0.0%	0.0%	0.0%	● 18.2%	NA	NA
U	92	238	90868	● 35.6%	0.0%	0.0%	0.0%	0.0%	● 0.6%	NA	NA

Table 2B. continued showing NH₃

Element	Atomic Num	Mass Num	NH ₃ (%)															
			M+	M(NH) ⁺	M(NH ₂) ⁺	M(NH ₃) ⁺	MNH(NH ₃) ⁺	MNH ₂ (NH ₃) ⁺	M(NH ₃) ₂ ⁺	MNH(NH ₃) ₂ ⁺	MNH ₂ (NH ₃) ₂ ⁺	M(NH ₃) ₃ ⁺						
			0	15	16	17	32	33	34	49	50	51						
Li	3	7	●	3.7%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%				
Be	4	9	●	3.6%	0.0%	0.4%	0.0%	0.0%	●	1.1%	0.2%	0.0%	●	1.9%	●	2.4%		
B	5	11	●	2.5%	0.0%	0.2%	0.0%	●	1.0%	0.0%	0.0%	●	4.9%	0.1%	0.0%			
Na	11	23	●	70.9%	0.0%	0.0%	0.1%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%			
Mg	12	24	●	70.7%	0.0%	0.0%	0.4%	0.0%	0.0%	0.0%	0.1%	0.0%	0.0%	0.0%	0.1%			
Al	13	27	●	62.2%	0.0%	0.0%	0.1%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%			
Si	14	28	●	9.2%	0.0%	●	2.7%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%			
P	15	31		0.3%	0.1%	0.1%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%			
S	16	32		0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%			
Cl	17	35	●	7.3%	0.0%	0.0%	0.1%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%			
K	19	39	●	83.9%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%			
Ca	20	44	●	72.9%	0.0%	0.1%	0.2%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%			
Sc	21	45	●	19.7%	●	7.5%	●	2.2%	0.1%	●	2.3%	0.4%	0.0%	●	1.3%	0.4%	0.1%	
Ti	22	47	●	22.5%	●	9.2%	●	1.5%	0.0%	●	2.0%	0.2%	0.0%	●	1.7%	0.3%	0.0%	
V	23	51	●	53.2%	●	1.1%	●	6.7%	0.3%	0.4%	●	0.5%	0.2%	0.5%	●	0.9%	0.2%	
Cr	24	52	●	75.8%	0.0%	0.0%	0.3%	0.0%	0.0%	0.0%	●	0.6%	0.0%	0.0%	0.0%	0.0%		
Mn	25	55	●	79.0%	0.0%	0.0%	0.3%	0.0%	0.0%	0.0%	0.2%	0.0%	0.0%	0.0%	0.0%	0.0%		
Fe	26	56	●	76.4%	0.0%	0.0%	0.4%	0.0%	0.0%	0.0%	●	3.7%	0.0%	0.0%	0.0%	0.4%		
Co	27	59	●	69.5%	0.0%	0.0%	●	0.7%	0.0%	0.0%	●	5.0%	0.0%	0.0%	●	1.1%		
Ni	28	60	●	72.7%	0.0%	0.0%	●	0.7%	0.0%	0.0%	●	2.2%	0.0%	0.0%	●	3.9%		
Cu	29	63	●	74.1%	0.0%	0.0%	●	0.6%	0.0%	0.0%	●	5.4%	0.0%	0.0%	0.0%	0.0%		
Zn	30	66	●	79.6%	0.0%	0.0%	●	1.1%	0.0%	0.0%	●	0.5%	0.0%	0.0%	0.0%	0.1%		
Ga	31	71	●	79.5%	0.0%	0.0%	0.2%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%		
Ge	32	72	●	34.5%	0.0%	●	14.3%	0.1%	0.0%	●	7.9%	0.4%	0.0%	●	1.0%	●	3.3%	
As	33	75	●	24.2%	0.1%	●	15.1%	●	0.6%	0.0%	●	2.2%	0.3%	0.0%	0.2%	0.0%		
Se	34	82*	●	47.4%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%		
Br	35	81		0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%		
Rb	37	85	●	72.3%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%		
Sr	38	88	●	83.3%	0.0%	0.0%	0.2%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%		
Y	39	89	●	31.6%	●	14.9%	●	3.8%	0.4%	●	3.9%	●	0.6%	0.1%	●	1.3%	0.4%	0.1%
Zr	40	90	●	19.8%	●	17.2%	●	1.4%	0.0%	●	3.7%	0.3%	0.0%	●	2.5%	0.4%	0.0%	
Nb	41	93	●	19.3%	●	6.8%	0.2%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	
Mo	42	95	●	60.7%	0.1%	0.1%	0.2%	0.0%	0.0%	0.0%	●	0.8%	0.0%	0.0%	0.0%	0.0%		
Ru	44	101	●	67.9%	0.0%	0.0%	0.3%	0.0%	0.0%	0.0%	●	0.8%	0.0%	0.0%	0.0%	0.1%		
Rh	45	103	●	64.5%	0.0%	0.0%	0.3%	0.0%	0.0%	0.0%	●	0.6%	0.0%	0.0%	0.0%	0.0%		
Pd	46	105	●	75.0%	0.0%	0.0%	0.3%	0.0%	0.0%	0.0%	●	0.8%	0.0%	0.0%	●	0.9%		
Ag	47	107	●	88.2%	0.0%	0.0%	0.2%	0.0%	0.0%	0.0%	●	0.6%	0.0%	0.0%	0.0%	0.0%		
Cd	48	111	●	82.3%	0.0%	0.0%	0.2%	0.0%	0.0%	0.0%	0.1%	0.0%	0.0%	0.0%	0.0%	0.0%		
Sn	50	118	●	58.1%	0.0%	0.0%	0.3%	0.0%	0.0%	0.0%	0.2%	0.0%	0.0%	0.0%	0.0%	0.0%		
Sb	51	121	●	70.3%	0.0%	0.0%	●	3.4%	0.0%	0.0%	●	0.5%	0.0%	0.0%	0.0%	0.0%		
Te	52	125	●	77.6%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%		
I	53	127		0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%		

*⁸²Se was used in order to allow measurement of the no gas mode sensitivity

Table 2B. continued showing NH₃

Element	Atomic Num	Mass Num	NH ₃ (%)												
			M+	M(NH) ⁺	M(NH ₂) ⁺	M(NH ₃) ⁺	MNH(NH ₃) ⁺	MNH ₂ (NH ₃) ⁺	M(NH ₃) ₂ ⁺	MNH(NH ₃) ₂ ⁺	MNH ₂ (NH ₃) ₂ ⁺	M(NH ₃) ₃ ⁺			
			0	15	16	17	32	33	34	49	50	51			
Cs	55	133	●	80.6%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	0.0%	
La	57	139	●	39.2%	●	23.4%	●	9.2%	●	0.7%	●	0.9%	0.2%	0.1%	0.1%
Ce	58	140	●	36.3%	●	17.5%	●	6.5%	●	0.4%	●	0.7%	0.3%	0.1%	0.1%
Pr	59	141	●	53.3%	●	1.7%	●	4.2%	●	0.3%	●	0.1%	0.1%	0.0%	0.0%
Nd	60	146	●	68.6%	●	0.7%	●	3.9%	●	0.2%	●	0.0%	0.1%	0.0%	0.0%
Sm	62	147	●	67.0%	●	0.0%	●	2.4%	●	0.1%	●	0.0%	0.1%	0.0%	0.0%
Eu	63	153	●	69.1%	●	0.0%	●	0.2%	●	0.1%	●	0.0%	0.0%	0.0%	0.0%
Gd	64	157	●	41.4%	●	14.2%	●	2.9%	●	0.2%	●	2.1%	0.3%	0.0%	0.0%
Tb	65	159	●	48.0%	●	8.3%	●	2.6%	●	0.1%	●	1.9%	0.3%	0.0%	0.0%
Dy	66	163	●	78.4%	●	0.3%	●	1.8%	●	0.1%	●	0.0%	0.1%	0.0%	0.0%
Ho	67	165	●	61.8%	●	0.5%	●	1.2%	●	0.1%	●	0.1%	0.1%	0.0%	0.0%
Er	68	166	●	64.9%	●	0.7%	●	1.4%	●	0.1%	●	0.1%	0.2%	0.0%	0.0%
Tm	69	169	●	65.8%	●	0.0%	●	0.4%	●	0.1%	●	0.0%	0.0%	0.0%	0.0%
Yb	70	172	●	64.0%	●	0.0%	●	0.1%	●	0.1%	●	0.0%	0.0%	0.0%	0.0%
Lu	71	175	●	74.9%	●	1.0%	●	2.1%	●	0.2%	●	0.2%	0.3%	0.0%	0.0%
Hf	72	178	●	16.7%	●	38.6%	●	2.0%	●	0.0%	●	5.6%	0.3%	0.0%	0.0%
Ta	73	181	●	15.4%	●	19.5%	●	1.8%	●	0.0%	●	0.2%	0.0%	0.0%	0.0%
W	74	182	●	25.5%	●	17.1%	●	0.9%	●	0.0%	●	0.0%	0.0%	0.0%	0.0%
Re	75	185	●	81.3%	●	0.6%	●	0.0%	●	0.0%	●	0.0%	0.0%	0.0%	0.0%
Ir	77	193	●	50.3%	●	3.8%	●	0.0%	●	0.3%	●	0.2%	0.0%	0.1%	0.0%
Pt	78	195	●	79.8%	●	0.0%	●	0.0%	●	0.4%	●	0.0%	0.0%	0.0%	0.0%
Au	79	197	●	72.3%	●	0.0%	●	0.0%	●	0.3%	●	0.0%	0.0%	0.0%	0.0%
Hg	80	202	●	0.0%	●	0.0%	●	0.0%	●	0.0%	●	0.0%	0.0%	0.0%	0.0%
Tl	81	205	●	77.6%	●	0.0%	●	0.0%	●	0.0%	●	0.0%	0.0%	0.0%	0.0%
Pb	82	208	●	77.5%	●	0.0%	●	0.0%	●	0.0%	●	0.0%	0.0%	0.0%	0.0%
Bi	83	209	●	78.1%	●	0.0%	●	0.0%	●	0.0%	●	0.0%	0.0%	0.0%	0.0%
Th	90	232	●	17.7%	●	36.6%	●	7.0%	●	0.3%	●	NA	NA	NA	NA
U	92	238	●	32.6%	●	26.7%	●	5.4%	●	0.3%	●	NA	NA	NA	NA

www.agilent.com

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc. 2014

Published May 30, 2014

Publication number: 5991-4585EN



Agilent Technologies