

Routine determination of trace rare earth elements in high purity Nd₂O₃ using the Agilent 8800 ICP-QQQ

Application note

Geochemistry, mining and materials science

Authors

Juane Song¹, Xiang-Cheng Zeng¹,
Dong Yan¹ and Wei-ming Wu²

¹ Agilent Technologies, China

² Jiangxi University of Science and
Technology, Jiangxi, China



Introduction

Advanced technology products containing rare earth elements (REEs) continue to evolve at a rapid pace. Consequently, the use of REEs has extended from well-established applications such as glass polishing, to high-performance magnets, high-tech catalysts, electronics, glass, ceramics, and alloys. An alloy of the second most abundant REE, neodymium (Nd), with iron and boron (NIB) is an important material in the super permanent magnets that are used in vehicle components, computer data storage devices, MRI scanners, and loudspeakers. Nd is also used to make glass and safety glasses (didymium) used by glassblowers and welders. As REEs are increasingly used in high-tech products, control of any impurities is important. For example, the presence of other REEs as contaminants in a purified single-element REE material may affect the functionality of the final product, so impurities in the REE oxide raw material must be carefully controlled.



Agilent Technologies

Inductively Coupled Plasma Mass Spectrometry (ICP-MS) is the most commonly used atomic spectrometry technique for the measurement of trace REEs. This is partly because the mass spectrum of the REEs is relatively simple, particularly when compared to the complex REE optical emission spectrum produced by techniques such as ICP-OES. The measurement of mid- and high-mass REEs in a low-mass REE matrix is, however, very challenging for ICP-MS because REEs have high metal-oxide (M-O) bond strengths, and the oxide ions of the low mass REEs overlap the preferred isotopes of the mid-mass and high-mass REEs. For example, in the analysis of trace REEs in high-purity Nd₂O₃, ¹⁴⁵Nd¹⁶OH₂⁺ and ¹⁴⁶Nd¹⁶OH⁺ overlap the preferred isotope of dysprosium (¹⁶³Dy⁺), ¹⁴³Nd¹⁶O⁺ overlaps the only isotope of terbium (¹⁵⁹Tb⁺) and ¹⁴⁸Nd¹⁶OH⁺ overlaps the sole isotope of holmium (¹⁶⁵Ho⁺). While separation of the trace REEs from the REE matrix can be performed using a chelating resin, this technique is time-consuming and needs to be customized to the particular analyte and matrix under investigation. Clearly there is a requirement for a method capable of the direct analysis of trace REEs in a variety of high purity REE matrices.

In this study, an Agilent 8800 Triple Quadrupole ICP-MS (ICP-QQQ) was used for the direct analysis of trace REEs in a high-purity Nd₂O₃ sample. In MS/MS reaction mode, the Nd-based interferences were all removed effectively, enabling the 8800 ICP-QQQ to determine all 13 REE impurities at trace levels in the high purity REE-matrix material.

Experimental

A standard configuration Agilent 8800 ICP-QQQ (G3663A #100) with nickel interface cones was used. The sample introduction system consisted 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. The 8800 ICP-QQQ features High Matrix Introduction (HMI) aerosol dilution technology, which reduces the total sample load introduced to the plasma by accurately and reproducibly diluting the sample aerosol using an additional argon gas flow. A preset plasma condition (HMI-Low) was selected in the MassHunter software

to achieve a robust plasma to permit the routine (long-term) analysis of samples containing a 500 ppm Nd₂O₃ matrix. To compare the efficiency of removing Nd-based polyatomic ions, five different collision/reaction cell (CRC) modes (no gas, He, O₂ mass-shift, NH₃ on-mass and NH₃ mass-shift) were tested. Table 1 summarizes the tuning conditions used in the study.

Table 1. Agilent 8800 ICP-QQQ tuning parameters (NH₃ as 10% NH₃ in He)

| Parameter | | No gas | He | O ₂ mass shift | NH ₃ on mass | NH ₃ mass shift |
|-----------------|--------|--------|-----|---------------------------|-------------------------|----------------------------|
| Scan mode | | SQ | | MS/MS | | |
| RF power | W | 1550 | | | | |
| Sampling depth | mm | 8 | | | | |
| Carrier gas | L/min | 0.6 | | | | |
| Dilution gas | L/min | 0.5 | | | | |
| Octopole bias | V | -8 | -18 | -4 | -18 | |
| Octopole RF | V | 190 | | 150 | 190 | |
| Cell entrance | V | -30 | -40 | -50 | -120 | |
| Cell exit | V | -50 | -60 | -70 | -120 | |
| Deflect | V | 10 | 0 | 4 | -5 | |
| Plate bias | V | -50 | -40 | -70 | -120 | |
| KED | V | 5 | 3 | -7 | -15 | |
| He | mL/min | - | 5 | - | 1 | |
| O ₂ | mL/min | - | - | 0.3 | - | |
| NH ₃ | mL/min | - | - | - | 8 | 3 |

Five calibration standards [0, 0.1, 0.5, 2.0 and 5.0 µg/kg (ppb)] were prepared from an Agilent REE mixed standard (Part no. 8500-6944). Rhodium (Rh) and rhenium (Re) were used as internal standards (ISTD). High purity (99.999%) Nd₂O₃ (Baotou Research Institute of Rare Earths, China) was dissolved gently in semiconductor grade HNO₃, diluted to 500 ppm as Nd₂O₃ and analyzed for 13 REE impurities by ICP-QQQ.

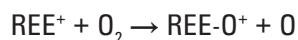
Results and discussion

All thirteen trace REEs were measured in the Nd₂O₃ sample using the no gas, He and O₂ cell modes. The results are summarized in Table 2. Pr and Sm were evidently present as impurities but, as expected, the Background Equivalent Concentration (BEC) of the other low- and mid-mass REEs, La, Ce, Pr, Sm, Eu and Gd, were comparable in all modes, as these elements are free from interferences due to Nd. In contrast, the BECs for the high-mass REEs were lower in He mode than in

no gas mode, suggesting that the high-mass REEs suffer interferences from Nd-derived polyatomic ions.

O₂ mass-shift mode

All 13 trace REEs react with O₂ efficiently to form REE-oxide ions, as shown below:



This reaction pathway, combined with the MS/MS capability of the 8800 ICP-QQQ, allows spectral interferences on the REEs to be avoided by using “mass-shift” mode. In O₂ mass-shift mode, all 13 REEs can be detected as their REE-O⁺ reaction product ions at 16 amu higher than the original elemental mass (M+16). This mode can be used in the CRC of a conventional quadrupole ICP-MS (ICP-QMS), but in practice the reaction chemistry is not consistent enough for routine analysis of real samples. In ICP-QMS, all of the ions produced in the plasma are able to enter the reaction cell, so other analyte ions or cell-formed reaction product ions may overlap the analyte ion of interest. For example, in the case of Dy measured as ¹⁶³Dy¹⁶O⁺ using O₂ mass-shift mode, ¹⁴⁶Nd⁺ also reacts with the O₂ cell gas to form ¹⁴⁶Nd¹⁶O₂H⁺ that would overlap the ¹⁶³Dy¹⁶O⁺ product ion at *m/z* 179. Furthermore, hafnium has a natural isotope at *m/z* 179 which could also overlap the target product ion. To avoid the formation of the overlapping NdOH⁺ product ion, the precursor ion ¹⁴⁶Nd⁺ (and any native ions such as Hf⁺ that would appear at the analyte product ion mass) must be removed before they enter the reaction cell; this can only be achieved with MS/MS mode on the ICP-QQQ. In MS/MS mode, the first quadrupole (Q1) operates as a single amu mass filter so that only ions with the targeted *m/z* enter the reaction cell and can react with the cell gas. All other masses are rejected by Q1, so they are not able to pass through the cell or undergo the reactions that would lead to interfering product ions.

In addition to the significant improvement in the analysis of the elements Tb, Dy and Ho, that suffer intense interference from NdO⁺, O₂ mass-shift mode also delivers improved BECs for the other high-mass REEs Er, Tm, Yb and Lu, indicating that these elements were also affected by interferences from Nd-based polyatomic ions in no gas mode. Likely overlaps are ¹⁵⁰NdOH₃⁺ on ¹⁶⁹Tm⁺; ¹⁴²NdN₂⁺, ¹⁴²NdCO⁺ and ¹⁴⁴NdCN⁺

on ¹⁷⁰Er⁺; ¹⁴²NdO₂⁺ on ¹⁷⁴Yb⁺; ¹⁴³NdO₂⁺, ¹⁴⁴NdONH⁺ and ¹⁵⁰NdC₂H⁺ on ¹⁷⁵Lu⁺. The Nd-based interferences on Er, Tm, Yb and Lu are not very significant (less than 0.2 µg/kg in the Nd₂O₃ matrix), but nevertheless O₂ mass-shift mode was shown to be an effective approach for the removal of all the polyatomic ion interferences, typically providing around a 5-10x lower BEC compared to no gas mode and He mode.

Table 2. BECs of 13 REEs in 500 ppm Nd₂O₃. All units are ppb.

| Element | Isotope | No gas | He | O ₂ mass shift | NH ₃ on mass | NH ₃ mass shift |
|---------|---------|--------|-------|---------------------------|-------------------------|----------------------------|
| La | 139 | 0.143 | 0.127 | 0.143 | - | - |
| Ce | 140 | 0.018 | 0.012 | 0.011 | - | - |
| Pr | 141 | 1.376 | 1.202 | 1.056 | - | - |
| Sm | 152 | 1.061 | 0.950 | 0.999 | - | - |
| Eu | 153 | 0.032 | 0.026 | 0.028 | - | - |
| Gd | 155 | 0.035 | 0.046 | 0.033 | - | - |
| Tb | 159 | 442.6 | 74.6 | 1.258 | - | 0.022 |
| Dy | 163 | 250.3 | 196 | 1.161 | 0.040 | - |
| Ho | 165 | 20.43 | 16.2 | 0.101 | 0.004 | - |
| Er | 170 | 0.065 | 0.020 | 0.013 | - | - |
| Tm | 169 | 0.084 | 0.031 | 0.003 | - | - |
| Yb | 174 | 0.251 | 0.120 | 0.058 | - | - |
| Lu | 175 | 0.014 | 0.006 | 0.004 | - | - |

NH₃ on-mass mode for Dy and Ho

A previous study [1] showed that NH₃ cell gas reacts with many of the polyatomic ions that interfere with the REEs. However, NH₃ also reacts quickly with some of the REE ions, leading to reduced sensitivity of < 1 cps/ppt for La, Ce, Nd, Sm, Gd, Tb and Lu. For the remaining REEs that react less quickly (or not at all) with NH₃, i.e. Pr, Eu, Dy, Ho, Er, Tm and Yb, on-mass measurement with NH₃ cell gas could therefore be a valuable approach for avoiding the Nd-based overlaps. The results in Table 2 show that NH₃ on-mass mode gave excellent results for Dy and Ho in the Nd₂O₃ matrix, with an improvement in BECs of 20x compared to O₂ mass-shift mode.

NH₃ mass-shift mode for Tb

For the REEs that react efficiently with NH₃ (La, Ce, Nd, Sm, Gd, Tb and Lu), NH₃ cell gas can be used in combination with mass-shift mode, where the target REE element is measured as its reaction product ion.

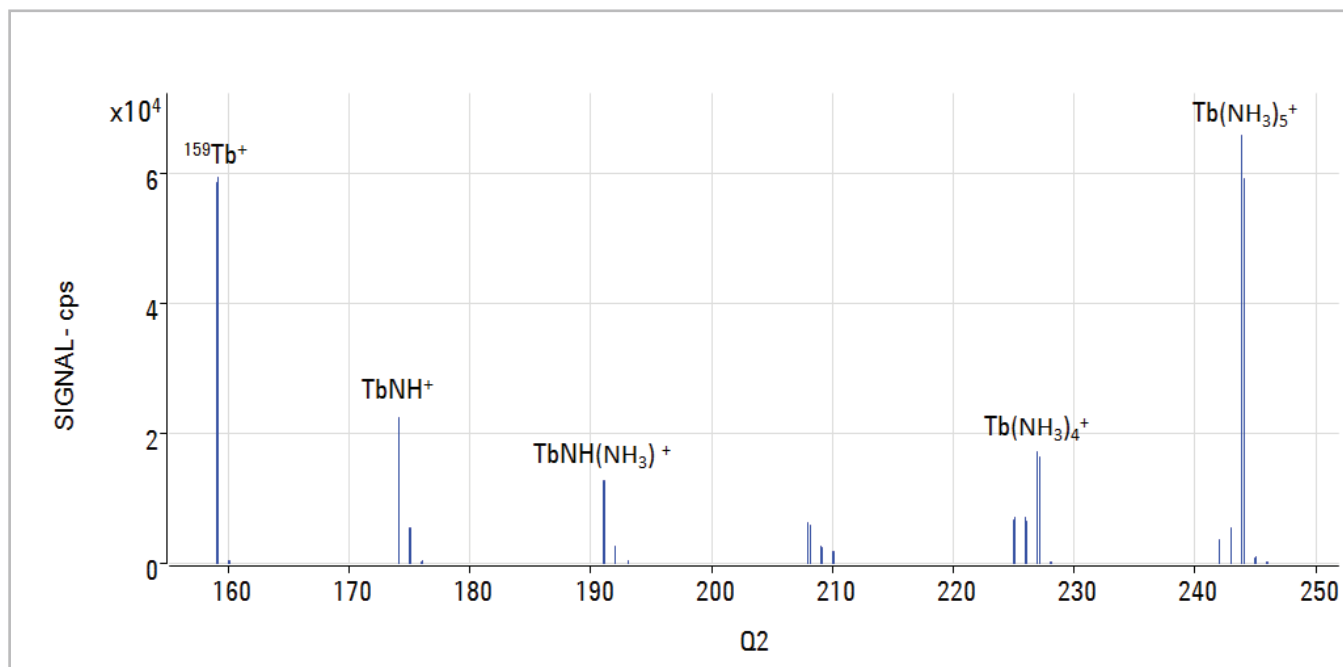


Figure 1. Product ion scan for ^{159}Tb in NH_3 mode

In this study, NH_3 mass-shift mode was investigated for the determination of Tb. In order to find the most appropriate ammonia-cluster product ion for Tb, a 10 ppb Tb solution was introduced and a product ion scan was performed. Q1 was set at m/z 159, allowing only ions at the mass of the target precursor ion (^{159}Tb) to enter the cell. Q2 was scanned over a selected mass range to measure all of the product ions formed in the cell via NH_3 reaction with ^{159}Tb (Figure 1). The scan revealed the four most abundant cluster ions: TbNH^+ (at $Q1 + 15$ amu), $\text{TbNH}(\text{NH}_3)^+$ (at $Q1 + 32$ amu), $\text{Tb}(\text{NH}_3)_4^+$ (at $Q1 + 68$ amu) and $\text{Tb}(\text{NH}_3)_5^+$ (at $Q1 + 85$ amu). For each of the four candidate cluster ions, the BEC of Tb in the Nd matrix was studied and TbNH^+ (m/z 174) was found to give the lowest BEC of 22 ppt for Tb in a 500 ppm Nd_2O_3 solution. This BEC is 50x lower than the result achieved in O_2 mass-shift mode, indicating the effective avoidance of the NdO^+ overlap.

Spike recoveries and long term stability

To validate the method, a spike-recovery test was carried out by analyzing a spiked solution containing 0.5 ppb mixed REE standard in a 500 ppm Nd_2O_3 sample. NH_3 on-mass mode was used for the determination of Dy and Ho, NH_3 mass-shift mode was used for Tb, and O_2 mass-shift mode for the remaining REEs. The results, which are shown in Figure 2, indicate that the

recoveries for all elements were good, demonstrating the sensitivity and effectiveness of the method.

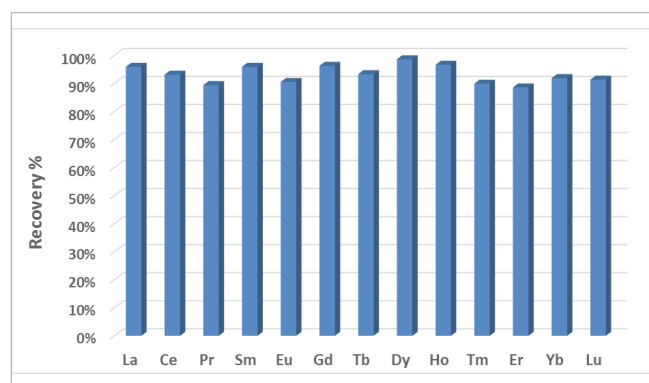


Figure 2. Recovery of 0.5 ppb REE spike added to 500 ppm Nd_2O_3 sample solution.

Long term stability over 2 hours was also measured using the same spiked sample. The stability of each REE signal is illustrated in Figure 3 and the average signal and RSD of each signal over the test period is summarized in Table 3. The data shows excellent stability (<5% RSD), demonstrating the applicability of the method for the routine analysis of high purity REE materials using the Agilent 8800 ICP-QQQ.

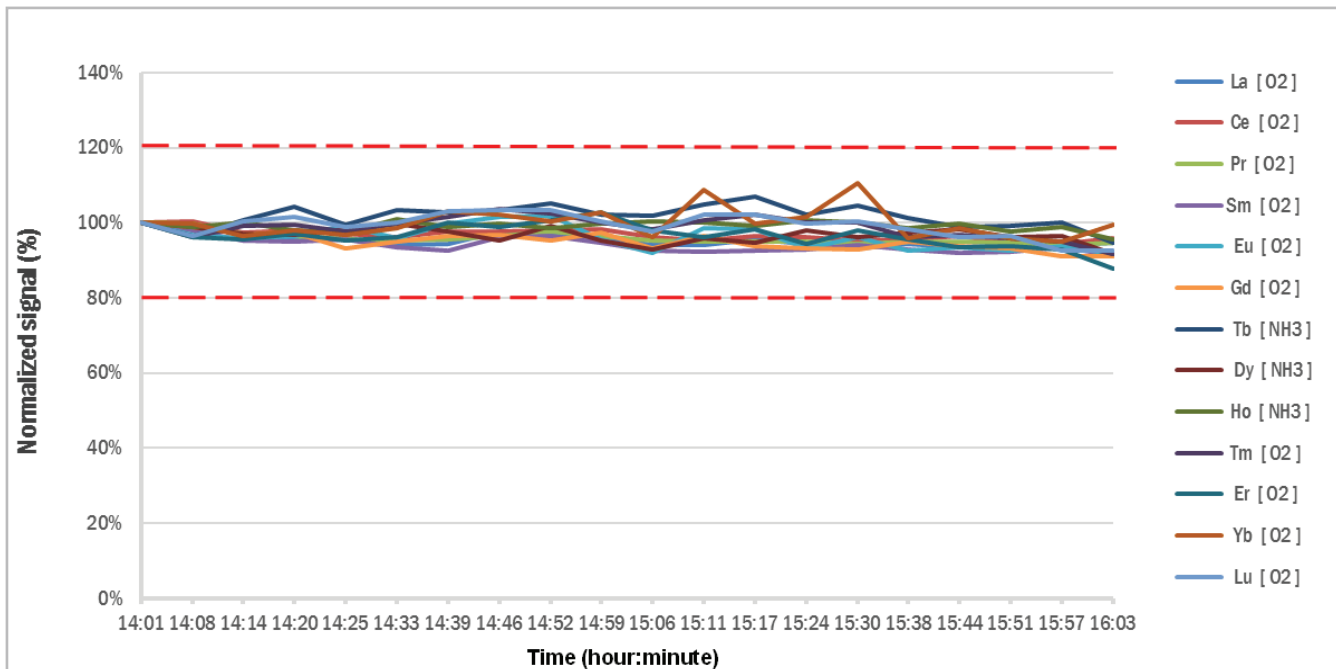


Figure 3. Long-term (2 hour) stability of 13 trace REEs spiked at 0.5 ppb in 500 ppm Nd₂O₃ solution

Table 3. Long-term (2 hour) %RSD for 13 trace REEs spiked at 0.5 ppb in 500 ppm Nd₂O₃ solution

| Element | Mass pair | Reaction gas | CPS | | Concentration (ppb) | |
|---------|-----------|-----------------|-------|------|---------------------|------|
| | | | Mean | RSD% | Mean | RSD% |
| La | 139->155 | O ₂ | 33443 | 2.0 | 0.625 | 1.1 |
| Ce | 140->156 | O ₂ | 24217 | 1.7 | 0.480 | 1.4 |
| Pr | 141->157 | O ₂ | 93453 | 1.7 | 1.512 | 1.2 |
| Sm | 152->168 | O ₂ | 24081 | 2.4 | 1.483 | 1.2 |
| Eu | 153->169 | O ₂ | 2171 | 3.3 | 0.485 | 3.2 |
| Gd | 155->171 | O ₂ | 4268 | 2.6 | 0.517 | 1.9 |
| Tb | 159->174 | NH ₃ | 1549 | 3.0 | 0.495 | 2.4 |
| Dy | 163->163 | NH ₃ | 939 | 2.3 | 0.530 | 3.1 |
| Ho | 165->165 | NH ₃ | 3508 | 1.3 | 0.484 | 2.3 |
| Er | 170->186 | O ₂ | 3735 | 3.4 | 0.463 | 3.7 |
| Tm | 169->185 | O ₂ | 18649 | 2.9 | 0.452 | 3.8 |
| Yb | 174->190 | O ₂ | 548 | 4.1 | 0.529 | 5.2 |
| Lu | 175->191 | O ₂ | 20623 | 3.2 | 0.470 | 4.0 |

Conclusion

The Agilent 8800 ICP-QQQ with MS/MS capability was used to successfully measure 13 REE impurities in a high-purity Nd₂O₃ sample solution. Oxygen mass-shift mode was effective at removing all of the polyatomic interferences from the Nd₂O₃ matrix. However, NH₃ on-mass mode further improved the BECs for Dy and Ho, and NH₃ mass-shift mode improved the BEC of Tb. Using NH₃ reaction mode, overlaps from NdO⁺, NdOH⁺, NdOH₂⁺ on Dy, Tb and Ho respectively were reduced by around two orders of magnitude compared to O₂ mass shift mode, allowing the trace (ppt) measurement of Dy, Tb and Ho in a 500 ppm Nd₂O₃ matrix.

With the combination of Agilent's HMI and MS/MS reaction cell mode, the 8800 ICP-QQQ provided good long term stability over 2 hours for all trace REEs present in the 500 ppm Nd₂O₃ sample, demonstrating the validity of the analytical method for the direct analysis of high purity Nd₂O₃.

Reference

1. Naoki Sugiyama and Glenn Woods, Direct measurement of trace rare earth elements (REEs) in high-purity REE oxide using the Agilent 8800 Triple Quadrupole ICP-MS with MS/MS mode, Agilent publication 5991-0892EN (2012).

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. 2015
Published January 20, 2015
Publication number: 5991-5400EN