

Highly-Sensitive Uranium Isotopic Analysis for Nuclear Safeguards, Using the NEPTUNE *Plus* with Jet Interface Option

Nicholas S. Lloyd, Claudia Bouman, and Johannes B. Schwieters
Thermo Fisher Scientific, Bremen, Germany

Keywords

NEPTUNE *Plus*, MC-ICP-MS, Uranium Isotope, Nuclear Safeguards

Introduction

This application note describes the analytical performance of the Thermo Scientific™ NEPTUNE™ *Plus* multicollector ICP-MS (MC-ICP-MS) for measuring uranium isotopes from 2 ng uranium samples.

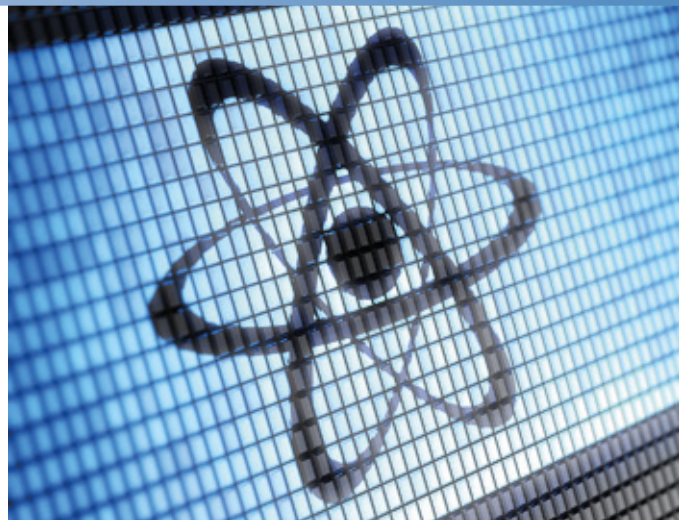
Instrumentation

Jet Interface for Ultimate Sensitivity

The NEPTUNE *Plus* MC-ICP-MS was equipped with the Jet Interface option for highest sensitivity (see also Application Note: 30187). The Jet Interface option comprises a high performance interface pump with Jet sample and X skimmer cones. Solutions were introduced by an ESI μ Flow PFA-100 self-aspirating nebulizer and a CETAC Aridus II desolvating nebulizer system. The sequence was automated using a CETAC ASX-112 FR autosampler.

The Thermo Scientific NEPTUNE *Plus* MC-ICP-MS was equipped with a multi-ion-counting array configured for nuclear applications. This array uses a combination of compact discrete dynode (CDD) type ion counters (see also Application Note: 30192) and full-sized SEM ion counters. The CDD type ion counters offer similar performance characteristics to the full-sized SEM.

The beams ^{238}U and ^{235}U were measured on Faraday cups. ^{234}U and ^{236}U were measured simultaneously in SEM ion counters with RPQ ion energy filters for improved abundance sensitivity (dual RPQ option). The abundance sensitivity from ^{238}U on m/z 236 (2 a.m.u.) was estimated to be ca. 1×10^{-7} .



Outline of Methodology

The NBS standard reference material U-030 was run as a standard, bracketing samples of U-010 (currently available as CRM U-010, New Brunswick Laboratory, Argonne, IL, USA). The uranium concentration of the U-010 'sample' solution was 2.3 ng/g (ppb) in 3 wt. % HNO_3 .

A certified 1 $\mu\text{g/L}$ (ppb) uranium concentration solution gave a ^{238}U signal of 1.22 V on a standard $10^{11} \Omega$ amplifier (1220 V/ppm at 87 $\mu\text{L/min}$). The sample utilization (ions detected/atoms in solution consumed) is estimated to be 2.1%.

Results

Combination of Highest Sensitivity and Stability

The data for each of the standards and isotopes are plotted in Figure 1, and statistics are summarized in Table 1. It can be seen that for each isotope the scatter of measured data is small and within the range of uncertainty of the certified reference materials. More recent and precise reference values from Richter and Goldberg (2003) are plotted for comparison against the certified values. Accuracy of the analytical data was not compromised by using the high-sensitivity Jet Interface option.

These data demonstrate that the analytical performance of the NEPTUNE Plus MC-ICP-MS for ^{235}U abundance analysis is well within the International Target Values for Measurement Uncertainties in Safeguarding Nuclear Materials (IAEA 2010).

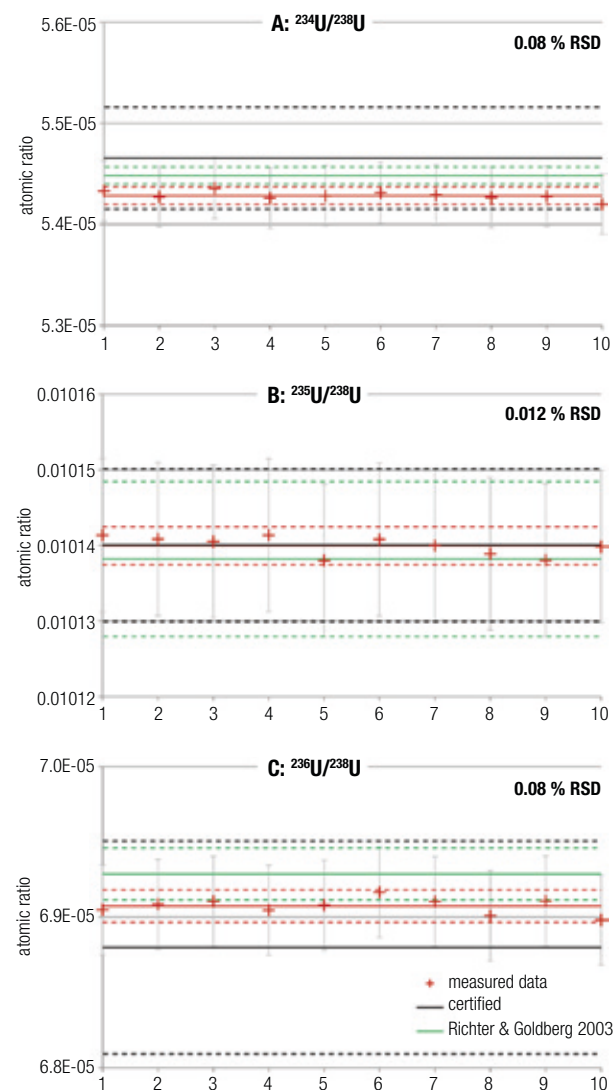


Figure 1. The scatter of data for the measurements is small compared to the uncertainty of the certified reference values (dashed lines show uncertainty at 95%). The 95% C.I. error bars for the measured data include the uncertainty propagated from the NBS U-030 standard's certified reference values.

Table 1. Mean and external reproducibility for 10 repeat analyses of NBS U-010 (externally corrected using NBS U-030).

2 ng NBS U-010, n=10	$^{234}\text{U}/^{238}\text{U}$	$^{235}\text{U}/^{238}\text{U}$	$^{236}\text{U}/^{238}\text{U}$
mean	5.43E-05	0.01014	6.91E-05
1RSD	0.08%	0.012%	0.08%

Conclusions

The data show that the NEPTUNE Plus MC-ICP-MS with Jet Interface option is a MC-ICP-MS that is capable of measuring high-precision $^{235}\text{U}/^{238}\text{U}$ ratios from small uranium samples.

The NEPTUNE Plus MC-ICP-MS is supported by a global network of dedicated service engineers, vital for continuous industry critical operations.



Figure 2. Thermo Scientific NEPTUNE Plus MC-ICP-MS.

Related Products

The Thermo Scientific TRITON Plus thermal ionization mass spectrometer (TIMS) is well established within the nuclear industry. It can optionally be equipped with the same ion counting configurations that are available on the NEPTUNE Plus MC-ICP-MS. The TRITON Plus TIMS offers new capabilities for the 'modified total evaporation' methodology, for improved minor isotope analysis. (Richter et al. 2011).



Figure 3. Thermo Scientific TRITON Plus TIMS.

The Thermo Scientific URANUS™ multi-collector gas source instrument is dedicated to the high-precision analysis of uranium isotopes directly from UF_6 samples. This system can be integrated into uranium enrichment production processes for online isotope ratio analysis without sample preparation.



Figure 4. Thermo Scientific URANUS multi collector mass spectrometer for uranium hexafluoride analysis.

Three high-precision analytical instruments cater for the needs of uranium isotope analysis within the nuclear industry, all of which are supported by a global network of dedicated service engineers.

References

- Application Note: 30136. *Improvements in TIMS High Precision Isotope Ratio Measurements for Small Sample Sizes*. Thermo Fisher Scientific, Bremen, Germany. Thermo Fisher Scientific, Bremen, Germany.
- Application Note: 30187. *NEPTUNE and NEPTUNE Plus: Breakthrough in Sensitivity using a Large Interface Pump and New Sample Cone*. Thermo Fisher Scientific, Bremen, Germany.
- Application Note: 30192. *New Compact Discrete Dynode Multipliers Integrated into the Thermo Scientific TRITON Variable Multicollector Array*. Thermo Fisher Scientific, Bremen, Germany.
- Application Note: 30238. *Uranium Isotopic Analysis for the Nuclear Industry Using the Thermo Scientific NEPTUNE Plus MC-ICP-MS*. Thermo Fisher Scientific, Bremen, Germany.
- IAEA. 2010. *International Target Values 2010 for Measurement Uncertainties in Safeguarding Nuclear Materials*. STR-368. International Atomic Energy Agency.
- Richter, S. & Goldberg, S.A. 2003. *Int. J. Mass Spectrom.* 229, 181–197.
- Richter, S. Kühn, H. Aregbe, Y. et al. 2011. *J. Anal. At. Spectrom.*, 26, 550–564.

www.thermoscientific.com/IRMS

©2014 Thermo Fisher Scientific Inc. All rights reserved. Teledyne Cetac Technologies formerly Cetac are trademarks of Teledyne Technologies Incorporated; ESI is a trademark of Elemental Scientific Inc.; ISO is a trademark of the International Standards Organization. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

Thermo Fisher Scientific (Bremen) GmbH
Management System Registered to
ISO 9001:2008

Africa +43 1 333 50 34 0	Denmark +45 70 23 62 60	Japan +81 45 453 9100	Singapore +65 6289 1190
Australia +61 3 9757 4300	Europe-Other +43 1 333 50 34 0	Latin America +1 561 688 8700	Spain +34 914 845 965
Austria +43 810 282 206	Finland +358 9 3291 0200	Middle East +43 1 333 50 34 0	Sweden +46 8 556 468 00
Belgium +32 53 73 42 41	France +33 1 60 92 48 00	Netherlands +31 76 579 55 55	Switzerland +41 61 716 77 00
Canada +1 800 530 8447	Germany +49 6103 408 1014	New Zealand +64 9 980 6700	UK +44 1442 233555
China 800 810 5118 (free call domestic) 400 650 5118	India +91 22 6742 9494	Norway +46 8 556 468 00	USA +1 800 532 4752
	Italy +39 02 950 591	Russia/CIS +43 1 333 50 34 0	

AN30269_E 02/14S

Thermo
SCIENTIFIC

Part of Thermo Fisher Scientific