

# High Sensitivity Laser Ablation MC-ICP MS

Nicholas S. Lloyd,<sup>1</sup> Claudia Bouman,<sup>1</sup> Johannes B. Schwieters,<sup>1</sup> Steve Shuttleworth<sup>2</sup> and John Roy<sup>2</sup>

<sup>1</sup>Thermo Fisher Scientific, Bremen, Germany; <sup>2</sup>Photon Machines, Redmond, WA, USA

## Keywords

NEPTUNE *Plus*, Laser Ablation, Hf Isotopes, Zircons, High Sensitivity, <sup>176</sup>Hf/<sup>177</sup>Hf, Jet Interface

## Goal

This application note demonstrates that high precision <sup>176</sup>Hf/<sup>177</sup>Hf isotope ratios can now be obtained from 25 μm diameter laser ablation spots.

## Introduction

High precision Hf isotope ratios can be obtained by simultaneous collection of Hf, Yb and Lu isotopes on Faraday cups. Typically LA spot sizes of 50 μm diameter are required to achieve targets of better than 100 ppm (epsilon unit) external precision (2RSD).

However, improved sensitivity would allow small detrital zircon grains to be analyzed, larger zircon crystals to be mapped for heterogeneity, or other accessory minerals with lower analyte concentrations to be analyzed.

The Jet Interface is an option on the Thermo Scientific<sup>™</sup> NEPTUNE<sup>™</sup> and NEPTUNE *Plus*<sup>™</sup> MC-ICP MS; it is also available on the Thermo Scientific ELEMENT 2/XR<sup>™</sup> SC-SF-ICP MS. Previously it has been shown to deliver breakthrough ICP-MS sensitivity for solution samples.<sup>1</sup>

A 25 μm diameter spot covers just 1/4 of the area of a conventional 50 μm diameter spot. In this application note, we demonstrate that precise and accurate Hf isotope ratios can be obtained from the zircon 91500 standard using a 25 μm diameter laser ablation spot size.

## Methods

### Laser Ablation

A Photon Machines Analyte.G2 laser ablation system with HelEx two-volume cell was run at 7 Hz and 6 J/cm<sup>2</sup> fluence (moderate settings) with a dwell time of 60 seconds for each spot. The excimer laser has a 193 nm wavelength and short pulse-width (ca. 4 ns). 0.9–1.0 L/min He was passed through the LA cell as a carrier gas.



Thermo Scientific NEPTUNE *Plus* MC-ICPMS and Photon Machines Analyte.G2 193 nm LA system.

## Mass Spectrometry

A NEPTUNE *Plus* MC-ICP MS was equipped with the Jet Interface option for increased sensitivity. The Jet Interface option comprises a high performance interface pump and high sensitivity cones. For desolvated solutions more than 2% of uranium ions can be detected;<sup>1</sup> accurate and precise Hf and Pb isotopes can be acquired for sub-ng solution samples.<sup>2</sup> 0.4–0.6 L/min Ar gas was mixed with the He carrier gas before the injector.

The collector configuration in Table 1 is possible with the standard NEPTUNE/NEPTUNE *Plus* MC-ICP-MS moveable Faraday collector array. By measuring <sup>173</sup>Yb/<sup>171</sup>Yb, the <sup>176</sup>Yb interference can be accurately corrected ( $\beta_{Yb} \neq \beta_{Hf}$ ).<sup>3</sup>

Table 1. Example cup-configuration for Hf isotopes on the NEPTUNE/NEPTUNE *Plus* MC-ICP MS. <sup>173</sup>Yb/<sup>171</sup>Yb is measured for accurate <sup>176</sup>Yb correction.

L4	L3	L2	L1	C	H1	H2	H3	H4
<sup>171</sup> Yb	<sup>173</sup> Yb	<sup>175</sup> Lu	<sup>176</sup> Hf	<sup>177</sup> Hf	<sup>178</sup> Hf	<sup>179</sup> Hf	<sup>180</sup> Hf	

Data were collected with 1.049 s integration times from Faraday cups equipped with 10<sup>11</sup> Ohm amplifiers (and precisely matched time constants). 1350 W RF power was used, and tuning was either for maximum LA Hf sensitivity or for lower oxide ratios (e.g. 1% UO/U).

## Nitrogen Addition

Nitrogen add gas at a rate of 7–11 mL/min was introduced by a mass flow controller installed in the NEPTUNE *Plus* MC-ICP MS. Nitrogen gas addition has previously been shown to increase LA sensitivity for most elements.<sup>4</sup>

## Data Analysis

Signal data were exported from the Thermo Scientific *multicollector* software and imported into the *Iolite* software package.<sup>5</sup> A Hf data reduction scheme was modified so that an external Yb fractionation factor could be used. The Yb fractionation factor was taken as the median value of  $^{173}\text{Yb}/^{171}\text{Yb}$  measured from all runs within each ablation session, as this is more precise than using an internal cycle-by-cycle correction from the low intensity Yb peaks (normalizing ratios<sup>6,7</sup>).

## Results

### Effect of N<sub>2</sub> Addition

N<sub>2</sub> addition increased LA Hf sensitivity by a factor of 2.4 on average. N<sub>2</sub> addition increased LA Hf sensitivity for both standard cones and for the high sensitivity cones. The optimal rate of N<sub>2</sub> addition was 7–11 mL/min, dependent on tune conditions.

### Effect of X Type Skimmer Cone

The X type skimmer cone increased LA Hf sensitivity by another factor of 1.5.

### Effect of Jet Type Sample Cone

The Jet type sample cone increased LA Hf sensitivity by another factor of 2.

### Cumulative Sensitivity Increase

The cumulative sensitivity increase from N<sub>2</sub> addition and the high sensitivity X type skimmer cone and Jet sample cone is shown in Figure 3. A sensitivity increase of 7.2 times is possible for Hf ions using this combination. The increase is slightly larger for U-Pb.

### Matrix Tolerance at High Sensitivity

Figure 1 shows a plot of measured  $^{179}\text{Hf}/^{177}\text{Hf}$  ratios for two different spot sizes. There is no difference in Hf mass fractionation between the 25 and 50  $\mu\text{m}$  diameter spot sizes, corresponding to an increase in matrix loading of 400%.

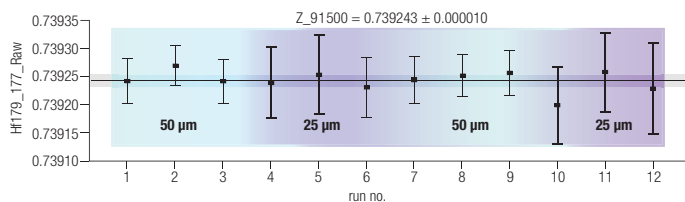


Figure 1. Raw  $^{179}\text{Hf}/^{177}\text{Hf}$  ratios measured from zircon 91500 using two different spot sizes. There is no difference in Hf mass fractionation between the 25 and 50  $\mu\text{m}$  diameter spot sizes (48 ppm 2RSD for one population).

## Effect of High Sensitivity on Accuracy and Precision

Figure 2 shows the repeatability of  $^{176}\text{Hf}/^{177}\text{Hf}$  measurements using a 25  $\mu\text{m}$  diameter spot. The results are precise and accurate, with a scatter of 86 ppm (2RSD).

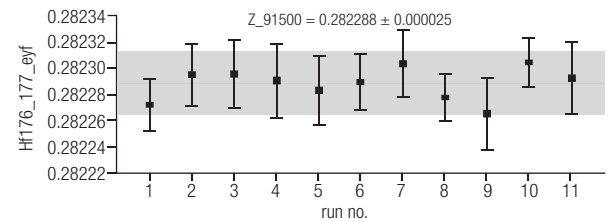


Figure 2. Plot of  $^{176}\text{Hf}/^{177}\text{Hf}$  ratios from the zircon standard 91500 using a 25  $\mu\text{m}$  diameter LA spot size. The 2RSD repeatability was 86 ppm, the results are accurate within uncertainty to accepted literature values.

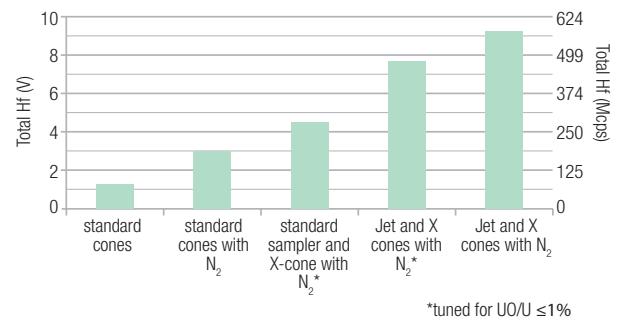


Figure 3. Plot of LA Hf sensitivity with and without N<sub>2</sub> addition and high sensitivity cones.

## Conclusions

- N<sub>2</sub> addition increases LA Hf sensitivity on the NEPTUNE *Plus* MC-ICP MS by a factor of 2.4.
- The Jet Interface option further increases sensitivity for LA Hf by a factor of 3.
- A cumulative sensitivity increase of 7.2 times is possible, whilst maintaining  $^{176}\text{Hf}/^{177}\text{Hf}$  accuracy.
- This increased sensitivity uniquely allows precise and accurate  $^{176}\text{Hf}/^{177}\text{Hf}$  ratios to be obtained from zircons using 25  $\mu\text{m}$  diameter LA spots.

## References

1. Bouman *et al.* 2009. *Geochim. Cosmochim. Acta*, 73(13), Supplement 1, A147.
2. Makishima and Nakamura 2010. *J. Anal. At. Spectrom.*, 25, 1712–1716.
3. Fisher *et al.* 2011. *Chem. Geo.*, 286, 32–47.
4. Shaheen and Fryer 2010. *J. Anal. At. Spectrom.* 25, 1006–1013.
5. Paton *et al.* 2011. *J. Anal. At. Spectrom.*, 26, 2508–2518.
6. Patchett and Tatsumoto. 1980. *Nature*, 288, 571–574.
7. Chu *et al.* 2002. *Anal. At. Spectrom.*, 17, 1567–1574.
8. Lloyd *et al.* 2013. *Min. Mag.* 6, 2029.

## Acknowledgements

The data in this application note were presented at the Goldschmidt Conference 2012.<sup>8</sup>

We would like to thank Matt Horstwood & Randy Parrish (NIGL, British Geological Survey) as well as Mohamed Shaheen & Brian Fryer (University of Windsor, CA).

## [thermoscientific.com/NeptunePlus](http://thermoscientific.com/NeptunePlus)

©2014 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. Analyte.G2 is used in trade by Photon Machines Inc and Photon-Machines is a trademark of Photon Machines Inc., Bozeman, Montana, USA. British Geological Survey is a trademark of Natural Environment Research Council, Swindon, Wiltshire, UK. The Goldschmidt Conference is a trademark of the Geochemical Society and the European Association of Geochemistry. 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.

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

AN30275\_E 05/14S

**Thermo**  
SCIENTIFIC

A Thermo Fisher Scientific Brand