

# Simultaneous in situ Analysis of U-Pb Age and Hf Isotopes of Zircon by Laser Ablation Sector-Field (MC-) ICP-MS

J. B. Schwieters, C. Bouman, M. Deerberg, J. D. Wills, Thermo Fisher Scientific, Bremen, Germany  
A. Gerdes, Institut für Geowissenschaften, Johann Wolfgang Goethe Universität, Frankfurt am Main, Germany

## Introduction

The combination of laser ablation with multicollector inductively coupled plasma mass spectrometers (LA-MC-ICP-MS) has been shown to produce high quality zircon age data from U-Th-Pb isotopic studies. These results however may not correctly indicate the number of zircon forming events, especially those arising from multiple growth and alteration processes. This limitation can be overcome by combining results from Pb-U and Lu-Hf analyses made on small, well defined zircon areas.

Static measurements of Pb-U and Lu-Hf can not be achieved in a single MC-ICP-MS analysis due to the large mass range that would need to be scanned. This has previously meant sequential analyses of the Pb-U and the Lu-Hf systems at the same zircon site<sup>1</sup>, e.g. an initial Pb-U analysis by laser ablation coupled to sector-field ICP-MS (LA-ICP-SFMS) and a subsequent LA-MC-ICP-MS analysis for Lu-Hf<sup>2</sup>. The sequential analyses were made as closely as possible in the same area (growth domain) of the zircon, for example a 40 µm Lu-Hf spot was drilled on top of a 30 µm U-Pb spot<sup>2</sup>.

In order to overcome the limitations of sequential sampling of U-Th-Pb and Hf isotopes from different sites in zircons, a fast scanning single collector and a multicollector inductive coupled plasma mass spectrometer have been coupled in parallel to a New Wave UP213 213 nm UV laser ablation system. A low volume laser ablation sample chamber with fast response (< 1 s) and rapid washout time was fitted instead of the standard sample chamber in order to enable sequential sampling of heterogeneous grains during time-resolved data acquisition on both instruments. The ablated material was transported in helium, split after the ablation cell via a Y-shaped connector and introduced simultaneously into the two mass spectrometers, the Thermo Scientific ELEMENT XR for U-Th-Pb analysis and Thermo Scientific NEPTUNE for high precision Hf isotope analysis.

For accurate age dating of zircons the control of common lead contaminations is essential. This can be done by monitoring mass <sup>204</sup>Pb which is not part of the U-Pb decay scheme. Since the natural abundance of <sup>204</sup>Pb is so small every common lead contribution measured on <sup>204</sup>Pb is magnified by about a factor of 15 to 18 on <sup>206</sup>Pb and <sup>207</sup>Pb intensities. This means the small <sup>204</sup>Pb signal has to be measured precisely and therefore the highest sensitivity is required to achieve the smallest statistical error. In comparison to quadrupole ICP-MS instruments<sup>3</sup> the ELEMENT XR sector-field ICP-MS is at least one order of magnitude more sensitive and as a consequence the statistical uncertainty on <sup>204</sup>Pb is significantly improved.

## Instrumentation

	U-Th-Pb analysis	Hf isotope analysis
Instrument	ELEMENT XR	NEPTUNE
Scan mode	E-Scan	Static
Scanned masses	202, 204, 206, 207, 208, 235, 232, 238	171, 173, 175, 176, 177, 178, 179, 180
Mass resolution	300	300
Dead time	18 ns	20 ns
Dwell time/isotope	4 ms	1 s
Settling time/isotope	1 ms	–
Number of scans	1080	50
Integration time	1 s (= 15 scans)	1 s
Background	20 s	
Ablation time	50 s	50 s
Carrier gas	0.5 L/min He (+1.1 L/min Ar)	
Laser	UP213 SA (New Wave Research; Nd:YAG, 213 nm)	
Spot size	55 to 80 µm	
Laser settings	5 Hz	
Laser fluence	~ 3 J/cm <sup>2</sup>	
Drill speed	0.6 µm/s	

Table 1: Operating conditions and instrument settings.

## Key Words

- ELEMENT XR
- NEPTUNE
- Isotope Ratio
- Laser Ablation
- Zircon

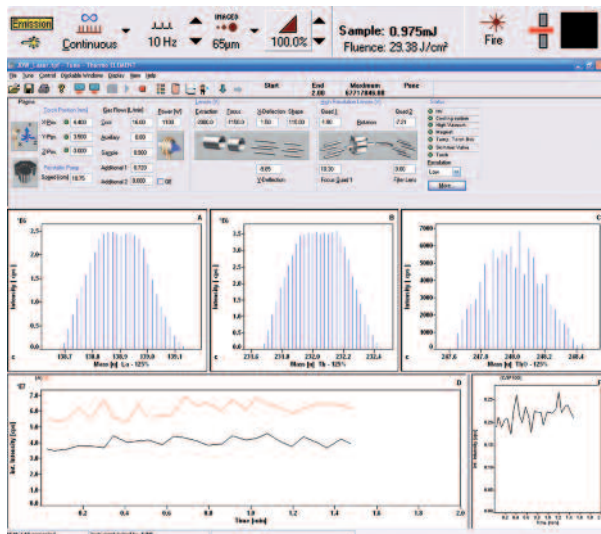


Figure 1: Screenshot from the ELEMENT software showing instrument settings and sensitivity obtained for La and Th in NIST 612 using the laser sampling parameters as shown.

The primary advantages of ICP-SFMS offers over other ICP-MS instrumentation for U-Th-Pb analyses are shown in Figure 1 and can be summarized as:

- The characteristic flat-top peaks of sector-field mass spectrometers in low mass resolution allow the reliable use of single point per peak analyses.
- The high sensitivity of ICP-SFMS allows for increased accuracy on the common lead correction that is made when measuring  $^{202}\text{Hg}$  as well as  $^{204}\text{Pb}+\text{Hg}$ . With less sensitive instrumentation small common lead contaminations detected at mass 204 would be otherwise magnified on masses  $^{206}\text{Pb}$  and  $^{207}\text{Pb}$ , leading to worsened Pb-U dating accuracy.

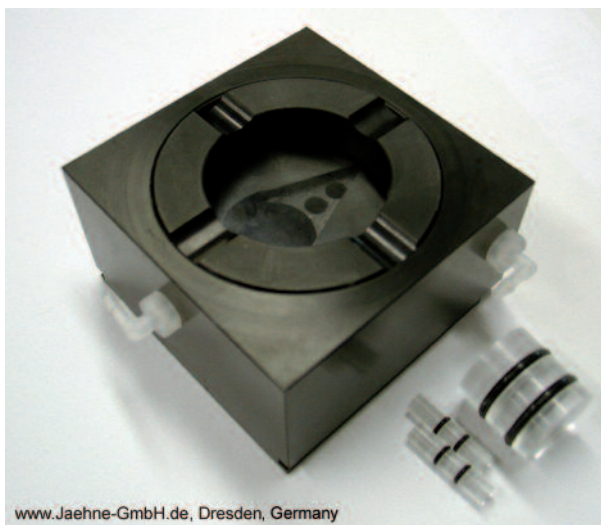
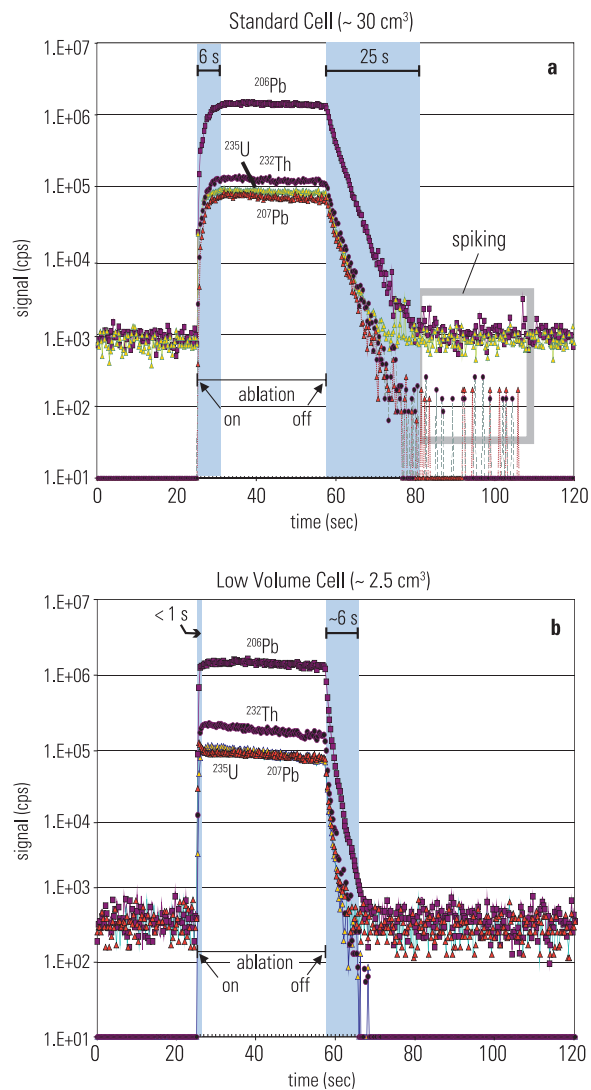


Figure 2: Low volume laser sample chamber used in all analyses.

The advantages of the low volume laser chamber<sup>4</sup> (Figure 2) for discrete laser sampling can be seen in Figure 3. Uptake delays of 1 s were achievable with the low volume cell, compared to 6 s with the standard sample cell. In the low volume cell, signal intensities after the end of the ablation dropped by 2 orders of magnitude in less than 4 s and 3 orders in about 6 s. In comparison, the signal from the standard cell needs more than 3 times longer (13 s and 25 s respectively) to drop to the same levels. Spiking is still observed in the standard cell up to 50 s after the end of ablation, while the low volume cell shows no spiking.



Frei & Gerdes, Chem. Geol. in press

Figure 3: Improvement in sample uptake and washout when using the low volume laser sample chamber.

## Results

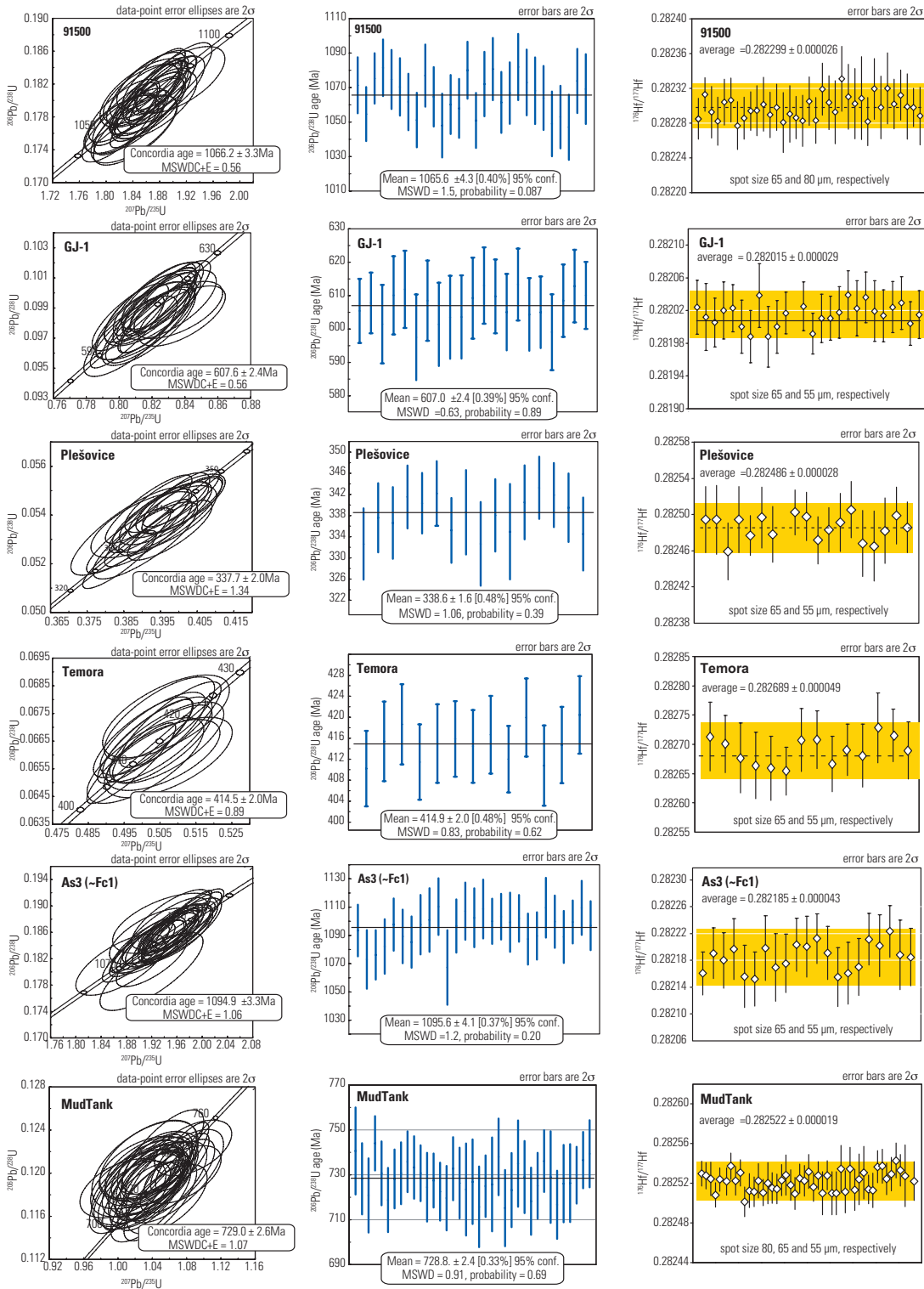


Figure 4: Simultaneous U-Pb and Hf isotope analyses in zircon.

	U-Pb age	<sup>176</sup> Hf/ <sup>177</sup> Hf
91500:	1065 Ma <sup>5</sup>	0.282306 ± 0.000008 <sup>8</sup>
GJ-1:	607 Ma <sup>2</sup>	0.282103 ± 0.000008 <sup>1</sup> (unpublished)
Plešovice:	337.1 Ma <sup>6</sup>	0.282484 ± 0.000008 <sup>6</sup>
Temora 2:	417 Ma <sup>7</sup>	0.282686 ± 0.000006 <sup>9</sup>
AS3 (-Fc1):	1099 Ma <sup>9</sup>	0.282184 ± 0.000016 <sup>7</sup>
MudTank:	732 Ma <sup>10</sup>	0.282507 ± 0.000006 <sup>9</sup>

Table 2: Reference U-Pb age and Hf isotope data for the zircons analyzed in Figure 4.

## Conclusions

Coupling the Thermo Scientific ELEMENT XR and the NEPTUNE to the laser ablation system was easy to achieve and demonstrated to be a powerful instrumental solution for the simultaneous analysis of U-Th-Pb and Hf isotope ratios in small zircons. The low volume laser cell used in this study demonstrated fast sample uptake and fast washout, resulting in low backgrounds without spiking signals. U-Pb ages and Hf isotope ratios of all zircons obtained in this study are in very good agreement with the published data.

## References

- 1 Gerdes, A., Zeh, A., 2006. Combined U-Pb and Hf isotope LA-(MC)-ICP-MS analyses of detrital zircons: comparison with SHRIMP and new constraints for the provenance and age of an Armorican metasediment in Central Germany. *Earth and Planetary Sciences Letters* 249, 47–61.
- 2 Gerdes, A., Zeh, A., 2008. Zircon formation versus zircon alteration – New insights from combined U-Pb and Lu-Hf in-situ LA-ICP-MS analyses, and consequences for the interpretation of Archean..., *Chemical Geology*, doi:10.1016/j.chemgeo.2008.03.005
- 3 Hong-Lin Yuan, Shan Gao, Meng-Ning Dai, Chun-Lei Zong, Detlef Günther, Gisela Helene Fontaine, Xia-Ming Liu, ChunRong Diwu; Simultaneous determinations of U-Pb age, Hf isotopes and trace element compositions of zircon by excimer laser-ablation quadrupole and multiple-collector ICP-MS; *Chem Geol.* 247, 110-118.
- 4 The laser ablation cell is of similar design to that described in: Davide Bleiner, Detlef Günther; Theoretical description and experimental observation of aerosol transport processes in laser ablation ICPMS, *J. Anal. At. Spectrom.*, 2001, 16, 449-456.
- 5 Wiedenbeck, M. et al, 1995. Three natural zircon standards for U-Th-Pb, Lu-Hf, trace element and REE analyses. *Geostandards Newsletter* 19, 1-23.
- 6 Slama, J. et al, 2008. Plešovice zircon – A new natural reference material for U-Pb and Hf isotopic microanalysis. *Chemical Geology* 249, 1-35
- 7 Black, L.P. et al, 2004. Improved Pb-206/U-238 microprobe geochronology by the monitoring of a traceelement related matrix effect; SHRIMP, ID-TIMS, ELA-ICP-MS and oxygen isotope documentation for a series of zircon standards. *Chemical Geology* 205, 115–140.
- 8 Woodhead, J.D., Hergt, J.M., 2005. A preliminary appraisal of seven natural zircon reference materials for in situ Hf isotope determination. *Geostandards and Geoanalytical Research* 29 (2), 183–195.
- 9 Paces, J.B., Miller, J.D., 1993. Precise U-Pb ages of Duluth Complex and related mafic intrusions, northeastern Minnesota: geochronological insights into physical, petrogenetic, paleomagnetic and tectonomagmatic processes associated with the 1.1 Ga midcontinent rift system. *Journal of Geophysical Research* 98, 13997–14013
- 10 Black, L.P., Gulson, B.L., 1978. The age of the Mud Tank carbonatite, Strangways Range, Northern Territory. *BMR Journal of Australian. Geology and Geophysics*, 3, 227–232.

*In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.*

**Africa-Other**  
+27 11 570 1840

**Australia**  
+61 2 8844 9500

**Austria**  
+43 1 333 50 34 0

**Belgium**  
+32 2 482 30 30

**Canada**  
+1 800 530 8447

**China**  
+86 10 8419 3588

**Denmark**  
+45 70 23 62 60

**Europe-Other**  
+43 1 333 50 34 0

**Finland/Norway/  
Sweden**  
+46 8 556 468 00

**France**  
+33 1 60 92 48 00

**Germany**  
+49 6103 408 1014

**India**  
+91 22 6742 9434

**Italy**  
+39 02 950 591

**Japan**  
+81 45 453 9100

**Latin America**  
+1 608 276 5659

**Middle East**  
+43 1 333 50 34 0

**Netherlands**  
+31 76 579 55 55

**South Africa**  
+27 11 570 1840

**Spain**  
+34 914 845 965

**Switzerland**  
+41 61 716 77 00

**UK**  
+44 1442 233555

**USA**  
+1 800 532 4752

**www.thermo.com**

©2009 Thermo Fisher Scientific Inc. All rights reserved. All 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 is certified  
DIN EN ISO 9001:2000

AN30170\_E 02/09C