





Figure 3: Thermo Scientific ELEMENT 2 HR-ICP-MS calibration curve for Sr.

## Results and Discussion

The calibration curves for the determination of Ca and Sr concentrations in the coral sample are shown in Figures 1 and 3.

The precision for each calibration point was < 1% RSD throughout the analysis sequence.

The correlation coefficients for all calibration curves were > 0.99996 (Figure 3).

The results of the Sr/Ca and Mg/Ca elemental ratio calculations are summarized in Tables 2 and 3.

ISOTOPE	CONCENTRATION [NG/ML]				
	RUN1	RUN 2	RUN 3	RUN 4	RUN 5
<sup>48</sup> Ca (Low Resolution)	1665.2	1652.2	1647.3	1643.2	1640.5
<sup>88</sup> Sr (Low Resolution)	32.64	32.34	32.28	32.21	32.17
Sr/Ca0.019601	0.019576	0.019594	0.019602	0.019612	

  

RATIO	AVER	STD. DEV.	RSD[%]
Sr/Ca	0.019597	0.0000134	0.07

Table 2: Sr/Ca elemental ratio calculations.

ISOTOPE	CONCENTRATION [NG/ML]				
	RUN1	RUN 2	RUN 3	RUN 4	RUN 5
<sup>48</sup> Ca (Low Resolution)	878.7	877.1	879.3	879.9	879.7
<sup>24</sup> Mg (Low Resolution)	2.793	2.790	2.793	2.800	2.798
Mg/Ca	0.003178	0.003181	0.003176	0.003182	0.003181

  

RATIO	AVER	STD. DEV.	RSD [%]
Mg/Ca	0.003180	0.0000023	0.07

Table 3: Mg/Ca elemental ratio calculations.

Under routine analytical conditions, an external precision of 0.07% RSD has been achieved for the measurement of both Sr/Ca and Mg/Ca elemental ratios. These values are comparable to previously reported values using a Thermo Scientific ELEMENT<sup>TM</sup> and superior to those obtained using quadrupole ICP-MS<sup>[1]</sup>. The improved precision obtained with the ELEMENT 2 shown in this work can be directly attributed to the characteristic 'flat top' peak shape in low resolution rather than the round shouldered peak of ICP-QMS. An additional factor that influences the precision of ratio measurements by ICP based techniques, is the instability of the plasma (so-called 'plasma flicker'). By using a stabilized sample introduction system the droplet size distribution of particles introduced into the plasma is minimized, significantly reducing 'plasma flicker'.

The comparatively simple analysis procedure presented in this report does not require time consuming sample preparation steps and is therefore suitable especially for a high sample throughput: 5 samples an hour (5 replicate analysis per sample). Because of the high sensitivity of the ELEMENT 2 (> 1 Mcps/ppb Sr) only a small sample amount (~ 2 µg) is required per analysis.

## Conclusions

The Thermo Scientific ELEMENT 2 is shown to be a powerful tool for the high precision measurement of Sr/Ca and Mg/Ca elemental ratios in corals. Using solution analysis, Sr/Ca and Mg/Ca elemental ratios in corals were determined with an external precision of 0.07% RSD (1σ). An autosampler enabled fully automatic unattended analysis. The complete measurement procedure is highly efficient and is suitable for high sample throughput.

## References

- [1] F. Le Cornec, T. Corrège, J. Anal. Atom. Spectrosc. 12 (1997), 969-973
- [2] J.W. Beck, R.L. Edwards, E. Ito, F.W. Taylor, J. Recy, F. Rougerie, P. Joannot, C. Henin, Science, 257 (1992), 644-647
- [3] D.W. Lea, P.A. Martin, Geochimica et Cosmochimica Acta, 60 (1996), 3143-3149
- [4] D.P. Schrag, in press
- [5] D.J. Sinclair, L.P.J. Kinsley, M.T. McCulloch, 62 (1998), 1889-1901
- [6] S.R. Hart, A.L. Cohen, Geochimica et Cosmochimica Acta, 60 (1996), 3075-3084
- [7] Y. Rosenthal, M.P. Field, R.M. Sherrell, Winter Conference on Plasma Spectrochemistry, 1999, Pau, France
- [8] T.M. Quinn, D.E. Sampson, Fall AGU, 1998, San Francisco

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

**Australia**  
+61 2 8844 9500

**Austria**  
+43 1 333 50340

**Belgium**  
+32 2 482 30 30

**Canada**  
+1 800 532 4752

**China**  
+86 10 5850 3588

**Denmark**  
+45 70 23 62 60

**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

**Netherlands**  
+31 76 587 98 88

**South Africa**  
+27 11 570 1840

**Spain**  
+34 91 657 4930

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

**Switzerland**  
+41 61 48784 00

**UK**  
+44 1442 233555

**USA**  
+1 800 532 4752

[www.thermo.com](http://www.thermo.com)

## Legal Notices

©2003, 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific Inc. products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

AN30001\_E 08/16C