

Single Mode Multi-Elemental Analysis of Traditional Chinese Medicine using the Thermo Scientific iCAP Qc ICP-MS

Jianfeng Cui, Julian D. Wills, Shona McSheehy Ducos and Lothar Rottmann,
Thermo Fisher Scientific, Bremen, Germany

Key Words

Traditional Chinese Medicine, ICP-MS, Single Mode KED

Goal

To demonstrate a simple, multi-elemental method for the routine determination of major and trace elements in traditional Chinese medicine using quadrupole ICP-MS.

Introduction

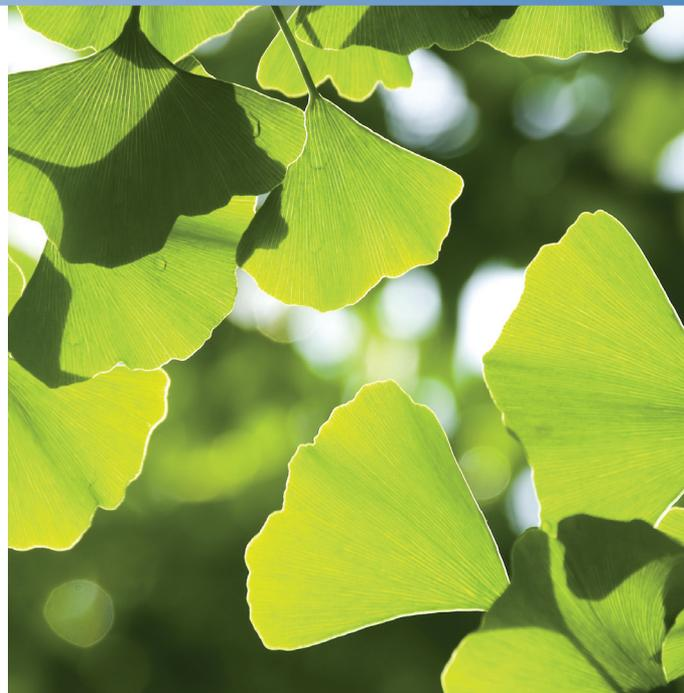
Traditional Chinese Medicine (TCM) is a group of treatments, including herbal medicines, that have been regularly employed over many hundreds of years and are still in use today. Strong tradition and culture have helped to maintain the popularity of TCM and Chinese herbal medicines but in recent years concerns have arisen that they may have been affected by mining and industrial waste. The uptake and storage of heavy metals in plant derived components of TCMs is under scrutiny as excessive exposure is linked to increased risk of metabolic disorders and other potentially degenerative effects¹.

Due to its excellent detection limits, quadrupole ICP-MS has become the standard technique for the determination of trace elements in food and plants – especially the “big four” toxic metals (arsenic, cadmium, lead and mercury). ICP-MS is therefore an important tool to assess the impact of prolonged use of TCM.

In order to demonstrate this, an authentic Ginkgo Leaf sample, a common ingredient in TCM, and two certified reference materials (CRMs) of similar matrix were prepared by microwave digestion and analyzed for their trace metal content by ICP-MS.

Method

All samples were prepared in triplicate with microwave digestion (using 0.5 g of each sample in 5 mL concentrated HNO₃ and 2 mL concentrated H₂O₂). After digestion each sample was made up to 50 mL using ultra pure water and were diluted a further 4-fold directly before analysis. Triplicate method blanks were performed for each microwave digestion batch.



All samples were quantified against an external calibration generated using multi-concentration, multi-elemental standards prepared from certified stock solutions. Final concentrations were method blank subtracted, internal standard corrected and quantified against the calibration curves.

The Thermo Scientific™ iCAP™ Qc ICP-MS was used to perform all analyses. Through its advanced low mass cut off in combination with excellent low mass sensitivity in cell modes, a single He KED measurement mode was used for the analysis of all target analytes.

Reference

- 1 Translated from the on-line source Chinese Pharmacopoeia Commission at:
<http://www.chp.org.cn/cms/business/cm/000127.html>

Table 1: Results for the NIST SRM 1515 Apple Leaves, INCT-OBTL Tobacco Leaves and an example TCM sample (Ginkgo Leaves). All concentrations are reported in mg·kg except where otherwise stated.

	LoD (ng·L ⁻¹)	MDL (µg·kg ⁻¹)	NIST SRM 1515 - Apple Leaves		INCT - OBTL Tobacco Leaves		Ginkgo Leaves
			Certified	Measured	Certified	Measured	Measured
⁷ Li	4.26	1.70		0.084 ± 0.004		24.4 ± 0.2	0.216 ± 0.019
¹¹ B	95.8	38.3	27 ± 2	27.9 ± 0.1	33.6 ± 2.2	36.4 ± 0.6	57.4 ± 4.4
²³ Na	20.3	8.12	24.4 ± 1.2	25.2 ± 0.6		213 ± 11	118 ± 8
²⁴ Mg	8.93	3.57	0.271 ± 0.008 (%)	0.273 ± 0.005 (%)	0.853 ± 0.034 (%)	0.851 ± 0.016 (%)	0.323 ± 0.018 (%)
³⁹ K	600	240	1.61 ± 0.02 (%)	1.61 ± 0.018 (%)	2.271 ± 0.076 (%)	2.385 ± 0.069 (%)	0.820 ± 0.053 (%)
⁴⁴ Ca	199	79.6	1.526 ± 0.015 (%)	1.51 ± 0.026 (%)	3.996 ± 0.142 (%)	3.902 ± 0.090 (%)	3.182 ± 0.194 (%)
⁵¹ V	0.327	0.131	0.26 ± 0.03	0.19 ± 0.04	4.12 ± 0.55	3.89 ± 0.07	0.135 ± 0.009
⁵² Cr	0.845	0.338	(0.3)	0.272 ± 0.004		4.81 ± 0.13	0.291 ± 0.029
⁵⁵ Mn	0.387	0.155	54 ± 3	53 ± 1	180 ± 6	187 ± 2	14 ± 0.7
⁵⁶ Fe	0.377	0.151	83 ± 5	77 ± 2	(0.149 %)	0.149 ± 0.003 (%)	107 ± 5
⁵⁹ Co	0.079	0.032	(0.09)	0.088 ± 0.001	0.981 ± 0.067	0.948 ± 0.016	0.022 ± 0.002
⁶⁰ Ni	0.417	0.167	0.91 ± 0.12	0.88 ± 0.04	8.50 ± 0.49	8.50 ± 0.03	0.224 ± 0.026
⁶³ Cu	0.617	0.247	5.64 ± 0.24	5.79 ± 0.15	10.1 ± 0.4	10.1 ± 0.1	4.03 ± 0.19
⁶⁶ Zn	1.52	0.608	12.5 ± 0.3	12.5 ± 0.3	52.4 ± 1.8	53.8 ± 0.4	4.58 ± 0.28
⁷⁵ As	1.17	0.47	0.038 ± 0.007	0.033 ± 0.001	0.668 ± 0.086	0.762 ± 0.055	0.087 ± 0.004
⁷⁸ Se	9.78	3.91	0.050 ± 0.009	0.052 ± 0.001		0.206 ± 0.036	0.067 ± 0.028
⁸⁵ Rb	0.224	0.090	10.2 ± 1.5	9.9 ± 0.3		16.5 ± 0.1	2.97 ± 0.08
⁸⁸ Sr	0.282	0.113	25 ± 2	25 ± 0.5	105 ± 5	102 ± 3	77 ± 2.2
⁹⁵ Mo	0.641	0.256	0.094 ± 0.013	0.092 ± 0.004	0.414 ± 0.062	0.387 ± 0.007	0.336 ± 0.024
¹⁰⁷ Ag	0.074	0.030		0.004 ± 0.001	0.053 ± 0.0105	0.055 ± 0.001	0.030 ± 0.002
¹¹¹ Cd	0.158	0.063	0.013 ± 0.002	0.013 ± 0.002	2.64 ± 0.14	2.80 ± 0.03	0.012 ± 0.001
¹²¹ Sb	0.138	0.055	(0.013)	0.012 ± 0.001	0.0755 ± 0.0125	0.070 ± 0.019	0.124 ± 0.003
¹³⁷ Ba	0.407	0.163	49 ± 2	50 ± 1	67.4 ± 3.8	67.9 ± 0.5	42.5 ± 1.3
²⁰² Hg	1.04	0.416	0.044 ± 0.004	0.051 ± 0.009	0.0209 ± 0.0013	0.0258 ± 0.0052	0.059 ± 0.006
²⁰⁵ Tl	0.024	0.009		0.012 ± 0.001	(0.0513)	0.051 ± 0.002	0.010 ± 0.001
²⁰⁸ Pb	0.035	0.014	0.470 ± 0.024	0.435 ± 0.056	2.01 ± 0.31	1.90 ± 0.03	0.899 ± 0.057
²³⁸ U	0.051	0.020	(0.006)	0.006 ± 0.001		0.077 ± 0.005	0.003 ± 0.001

Results

Figure 1 shows typical external calibration curves for Co and As in KED mode.

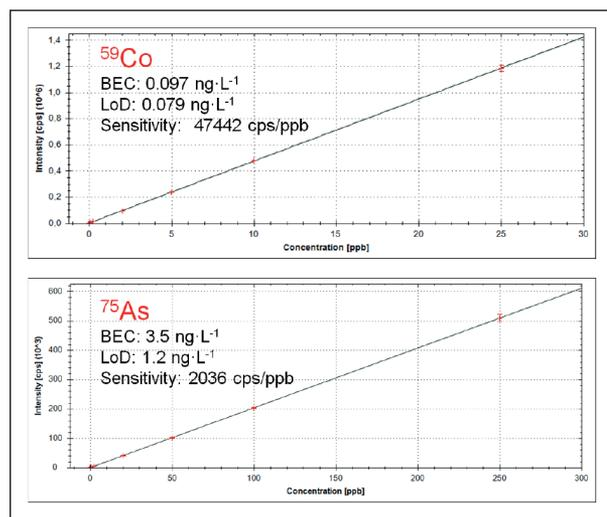


Figure 1: Fully quantitative five point calibration curves.

The results for the Apple and Tobacco Leaf CRMs and a genuine TCM (Ginkgo Leaves) are presented in Table 1. The concentrations determined for all trace metals in the two CRMs analyzed are within the certified ranges. The elemental profile of the Ginkgo Leaf sample analyzed is similar to that of the Apple Leaf CRM.

Conclusion

The iCAP Qc ICP-MS has been shown to be an excellent tool for the accurate, multi-elemental analysis of trace metals in complex samples, offering:

- A wide dynamic range (measuring µg·kg⁻¹ to % levels in a single analysis).
- High instrument sensitivity.
- Freedom from interferences using a single He KED analysis mode.
- An easy-to-use workflow enabled by the Thermo Scientific™ Qtegra™ Intelligent Scientific Data Solution (ISDS).

www.thermofisher.com

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manner 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.

Africa +43 1 333 50 34 0

Australia +61 3 9757 4300

Austria +43 810 282 206

Belgium +32 53 73 42 41

Canada +1 800 530 8447

China 800 810 5118 (free call domestic)
400 650 5118

Denmark +45 70 23 62 60

Europe-Other +43 1 333 50 34 0

Finland +358 9 3291 0200

France +33 1 60 92 48 00

Germany +49 6103 408 1014

India +91 22 6742 9494

Italy +39 02 950 591

Japan +81 45 453 9100

Latin America +1 561 688 8700

Middle East +43 1 333 50 34 0

Netherlands +31 76 579 55 55

New Zealand +64 9 980 6700

Norway +46 8 556 468 00

Russia/CIS +43 1 333 50 34 0

Singapore +65 6289 1190

Spain +34 914 845 965

Sweden +46 8 556 468 00

Switzerland +41 61 716 77 00

UK +44 1442 233555

USA +1 800 532 4752

Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand