

Robust single method determination of major and trace elements in foodstuffs using the Thermo Scientific iCAP PRO X Duo ICP-OES

Author

Nora Bartsch,
Application Specialist,
Thermo Fisher Scientific,
Bremen, Germany

Keywords

Foodstuffs, Micronutrients,
Microwave digestion,
Toxic elements

Goal

This application note demonstrates the ability of the Thermo Scientific™ iCAP™ PRO Series ICP-OES to determine trace elements and major components in foodstuffs to comply with regulations.

Introduction

Food protection from potentially hazardous contaminants has become a major topic of public interest. As well as the standard regulatory testing, it is necessary to account for contaminants, which may enter the food chain via many means, such as industrial pollution or environmental contamination, i.e., polluted rainfall on crops. Once toxic elements are in the food chain, they can pose significant health risks. With this in mind and the increasing number of micronutrients requiring determination, it is critical that the method of testing is a rigorous and reliable one. The organization charged with the development of the Codex standards (worldwide food and consumer regulations) is the Codex Alimentarius Commission, which is an intergovernmental body jointly sponsored by the Food and Agriculture Organization and the World Health Organization.

Instrumentation

The Thermo Scientific iCAP PRO Series ICP-OES employs a high-resolution Echelle spectrometer with a charge injection device (CID) detector. Advancements in CID technology allow this detector to feature lower noise and better separation of spectral orders than any of its predecessors. With the new optical design, the full spectrum from 167 nm up to 852 nm can be measured with a single exposure. A Thermo Scientific™ iCAP™ PRO X Duo ICP-OES was chosen for this analysis as this enables maximum sensitivity using axial view whilst maintaining excellent matrix tolerance in radial viewing mode. The instrument parameters used are listed in Table 1.

Table 1. Instrument parameters.

Parameters	Setting	
Pump tubing	Sample Tygon® orange/white	Drain Tygon® white/white
Spray chamber	Glass cyclonic	
Nebulizer	Glass concentric	
Center tube	2.0 mm	
Pump speed	45 rpm	
Nebulizer gas flow	0.6 L·min ⁻¹	
Auxiliary gas flow	0.5 L·min ⁻¹	
Coolant gas flow	12.5 L·min ⁻¹	
RF power	1150 W	
Exposure time	Axial	Radial
	15 s	15 s

Method development and analysis

Initially, more than one wavelength was selected for each element (using both axial and radial view). The subarrays for each wavelength were then examined and the most appropriate wavelength for the application was chosen (Table 2) based on the presence of interferences, calibration curve quality, QCs, and CRMs and the required linearity for the element. The subarray plots for each element can be easily manipulated by the analyst, allowing the optimum peak integration and background correction points to be selected. The Thermo Scientific™ Qtegra™ Intelligent Scientific Data Solution™ (ISDS) Software was used for data acquisition and provides easy options for post-analysis data manipulation. Qtegra ISDS Software has integrated Quality Control (QC) checks that allows the user to define automatically controlled actions to ensure data integrity, as well as a Flags and Limits function which flags samples according to user specifications.

Table 2. Element, wavelength and plasma view used.

Element	Wavelength (nm)	Plasma view
Ca	317.933	Radial
Cu	324.754	Axial
Fe	259.940	Axial
Mg	285.213	Radial
Mn	257.610	Axial
Ni	231.604	Axial
P	178.284	Radial
Pb	220.353	Axial
Zn	206.200	Axial

Sample preparation

Samples were prepared using certified reference material (CRM) standards. 0.5 g aliquots of rice flour (NBS1568b) and bovine liver (NBS1577c) were digested in 9 mL TraceMetal™ grade nitric acid (Fisher Chemicals, Loughborough, UK) using a standard food method program in a high pressure microwave system. The final digests were made up to 50 mL with deionized water before analysis.

Standard preparation

High purity standards (1000 mg·kg⁻¹ single element standards) were used to prepare the calibration standards for this method. All samples and standards were acid matched. Table 3 indicates the concentration of each of the standards, selected to cover the concentration range of the samples.

Table 3. Standard concentrations in mg·kg⁻¹.

Element	STD 1	STD 2	STD 3	STD 4
Pb	0.01	0.05	0.2	0.5
Mn, Ni, Zn, Cu	0.1	0.5	2	5
Ca, P, Fe, Mg	1	5	20	50

Results

The instrument was calibrated using a blank and at least two standards for each element. After inspection, a linear fit was applied to all elements. The calibration standards and samples were analyzed in a single sequence with an acid rinse between samples. The sample data was measured by interpolation and results are shown in Table 4. Suitable dilutions were made to over-range elements to ensure they fell within the calibration range. Method detection limits (MDLs) were also established by analyzing the acid matched calibration blank using a 10 replicate analysis and multiplying the standard deviation of the analysis by 3. This was repeated three times and the average values for method detection limits are also summarized in Table 4.

Conclusion

The results table shows that major and trace elements were measured with equal success. Precise, accurate results for digested foodstuff samples are easily attained on the Thermo Scientific iCAP PRO X Duo ICP-OES. The full wavelength coverage of the unique CID detector allowed the optimum wavelengths to be selected and the whole spectrum to be measured with one exposure, while the sensitivity of the axial view provided the low detection limits for this application.

Table 4. Results of measurement, CRM and MDLs in mg·kg⁻¹ and calculated recoveries in %.

Element	SRM1577c			SRM1568b			MDL µg·kg ⁻¹
	CRM	Measured	Recovery (%)	CRM	Measured	Recovery (%)	
Ca	131	124	95	118.4	107	90	1.90
Cu	275.2	272.3	99	2.35	2.13	91	0.4
Fe	197.94	205.97	104	7.42	7.78	105	1.5
Mg	580	579	100	549	478	87	1.23
Mn	10.46	10.12	97	19.2	18.1	94	0.04
Ni	0.0353	0.0325	92	-	-	-	0.52
P	-	-	-	1490	1302	87	12
Pb	0.0628	0.0623	99	< MDL	-	-	2.9
Zn	179.1	163.3	91	19.16	15.59	81	0.28

Find out more at thermofisher.com/icp-oes

ThermoFisher
SCIENTIFIC

©2020 Thermo Fisher Scientific Inc. All rights reserved. TraceMetal is a trademark of Fisher Scientific. Tygon is a trademark of Saint-Gobain Corporation. All other 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.

AN44423-EN 0320C