

# Analysis of elemental impurities in electronic grade copper sulfate using the Thermo Scientific iCAP 7400 ICP-OES Duo

## Authors

James Hannan,  
Application Specialist,  
Thermo Fisher Scientific,  
Hemel Hempstead, UK  
Thomas Huang,  
Application Engineer,  
Joy Allied Technology Inc., Taiwan  
Vicki Wu, Application Manager,  
Joy Allied Technology Inc., Taiwan

## Keywords

Electronic grade copper sulfate,  
 $\text{CuSO}_4$ , Raw materials testing

## Goal

This application note focuses on the analysis of electronic grade copper sulfate demonstrating the superior performance and high optical resolution of the Thermo Scientific iCAP 7400 ICP-OES Duo for the analysis of this complex matrix sample.

## Introduction

As technology advances, increasingly sophisticated manufacturing techniques are required for electronic goods. The main electrical conductor of these technologies is high purity copper. Components such as High Density Interconnect (HDI) printed circuit boards, flip chips used in bumping connections and wafer components, are manufactured using high purity copper sulfate. Electro-plating is used to bind the positive copper ions, from the copper sulfate, onto the required surface. As a result of this electroplating technique, any other positive ions present in the copper sulfate, such as iron, calcium or chromium, are also plated to the surface, producing impurities in the copper and reducing its conductive efficiency. Therefore any impurities in the copper sulfate must be quantified, prior to use, in order to maintain the high quality required.

Analysis of impurities in copper sulfate by Inductively Coupled Plasma-Optical Emission Spectrometry (ICP-OES) is an ideal choice of technique as it enables accurate, high throughput, multi-element analysis while requiring minimal sample preparation.

## Instrumentation

For this analysis the Thermo Scientific™ iCAP™ 7400 ICP-OES Duo was used as it provides the most cost effective solution and speed of analysis, while its axial plasma view provides the highest sensitivity required for trace analysis.

## Method development and analysis

The solid electronic grade copper sulfate ( $\text{CuSO}_4$ ) samples were dissolved in ultra-pure deionized (DI) water (resistivity  $>18 \text{ M}\Omega/\text{cm}$ ). 5 g of solid sample was dissolved in 50 ml of DI water and made up to a final volume of 250 ml; these samples were then ready for analysis. Multi-element calibration standards were made at concentrations of 0, 20, 50 and  $100 \mu\text{g}\cdot\text{L}^{-1}$  (ppb) and matrix matched to the samples (2%  $\text{CuSO}_4$ ).

A LabBook was created using the Thermo Scientific™ Qtegra™ Intelligent Scientific Data Solution™ (ISDS) Software to analyze the elements of interest, the elements and analytical wavelengths selected can be seen in Table 2 and the measurement modes, acquisition parameters and sample introduction accessories used are shown in Table 1. Instrument calibration plots were created using the calibration standards (0, 20, 50 and  $100 \mu\text{g}\cdot\text{L}^{-1}$ ).

**Table 1. Instrument parameters.**

Parameter	Setting
Pump Tubing (Standard Pump)	Sample Tygon® orange/white Drain Tygon® white/white
Spray Chamber	Glass cyclonic
Nebulizer	AeroSalt
Center Tube	2.0 mm
Pump Speed	50 rpm
Nebulizer Gas Flow	$0.45 \text{ L}\cdot\text{min}^{-1}$
Auxiliary Gas Flow	$0.5 \text{ L}\cdot\text{min}^{-1}$
Coolant Gas Flow	$12 \text{ L}\cdot\text{min}^{-1}$
RF Power	1150 W
Exposure Time	UV 20 s, Vis 5 s

## Results

The method detection limits for this method (MDL) were calculated by analyzing a matrix matched blank with ten repeats, then multiplying the standard deviation (SD) by a factor of 3. The calculated detection limits can be seen in Table 2.

The suitability of the analytical method was demonstrated by analyzing a series of matrix spiked samples. Five electronic grade copper sulfate samples were spiked with  $50 \mu\text{g}\cdot\text{L}^{-1}$  of all the target elements. Both the spiked and unspiked samples were analyzed and their measured concentrations compared. The results of this analysis along with the spike recoveries are shown in Table 3. The spike recoveries of all elements, for all samples, are between 92% and 105%.

In order to demonstrate the method and instrument stability, a sample of electronic grade copper sulfate was spiked with all of the target elements at  $50 \mu\text{g}\cdot\text{L}^{-1}$  and then analyzed repeatedly over a period of 60 minutes, the results of which can be seen in Figure 1. The average results, standard deviation (SD) and relative standard deviation (RSD) can be seen in Table 2. The relative standard deviations for all elements were less than 2%, with the exception of lead which was below 4%.

**Table 2. Average, SD and RSD of spiked  $\text{CuSO}_4$  solution and detection limits (DL) achieved.**

Element	Wavelength (nm)	Plasma view	Average ( $\mu\text{g}\cdot\text{L}^{-1}$ )	SD ( $\mu\text{g}\cdot\text{L}^{-1}$ )	RSD (%)	DL ( $\mu\text{g}\cdot\text{L}^{-1}$ )
Ag	328.068	Axial	49.95	0.69	1.38	1
As	189.042	Axial	52.15	0.57	1.09	2.3
Ca	393.366	Axial	52.9	0.56	1.07	0.04
Cd	226.502	Axial	50.21	0.24	0.47	0.15
Co	228.616	Axial	53.35	0.34	0.63	0.33
Cr	267.716	Axial	51.71	0.49	0.94	0.47
Fe	259.940	Axial	62.5	0.55	0.88	0.82
In	230.606	Axial	53.47	1.02	1.9	4.2
K	766.490	Axial	48.16	0.86	1.79	0.44
Mg	279.553	Axial	51.76	0.47	0.91	0.03
Mn	257.610	Axial	50.94	0.43	0.84	0.17
Na	589.592	Axial	51.99	1	1.92	0.2
Ni	231.604	Axial	52.72	0.31	0.59	0.39
Pb	168.215	Axial	50.69	1.58	3.13	4.1
Sn	189.989	Axial	53.13	0.33	0.62	0.67
Ti	336.121	Axial	50.97	0.47	0.92	0.16
Tl	190.856	Axial	48.63	0.69	1.43	2.2
Zn	206.200	Axial	51.37	0.25	0.48	0.23

Table 3. Spiked/unspiked sample concentrations in  $\mu\text{g}\cdot\text{L}^{-1}$  and spike recoveries in %.

Element	346 50X	346 50X spiked	Spike recovery (%)	348 50X	348 50X spiked	Spike recovery (%)	484 50X	484 50X spiked	Spike recovery (%)	486 50X	486 50X spiked	Spike recovery (%)
Ag	<DL	48.26	96.52	<DL	49.3	98.6	<DL	48.95	97.9	<DL	48.66	97.32
As	<DL	50.51	101.03	<DL	50.24	100.48	<DL	51.77	103.54	<DL	52.08	104.16
Ca	0.62	50.96	100.69	1.16	51.47	100.62	0.56	52.04	102.96	1.77	52.31	101.07
Cd	<DL	49.08	98.16	<DL	49.09	98.18	<DL	50.38	100.76	<DL	49.97	99.94
Co	<DL	50.31	100.62	<DL	50.38	100.76	<DL	51.73	103.46	<DL	51.2	102.4
Cr	0.49	49.56	98.14	<DL	49.55	99.1	<DL	50.65	101.3	0.46	50.94	100.96
Fe	8.38	57.57	98.38	9.38	58.01	97.25	7.35	57.58	100.47	12.63	62.57	99.88
In	<DL	47.21	94.42	<DL	47.7	95.4	<DL	48.19	96.38	<DL	48.32	96.64
K	<DL	49.19	94.38	<DL	46.24	92.48	<DL	47.33	94.66	<DL	46.7	93.4
Mg	0.18	49.73	99.09	0.25	49.43	98.37	0.2	50.57	100.74	0.36	50.35	99.98
Mn	<DL	49	98	<DL	48.88	97.76	<DL	49.91	99.82	<DL	49.74	99.48
Na	0.82	51.71	101.77	1.13	50.33	98.41	1.62	51.83	100.42	1.09	50.94	99.69
Ni	<DL	50.53	101.06	<DL	50.79	101.58	<DL	51.98	103.96	ND	51.74	103.48
Pb	<DL	47.27	94.54	<DL	46.05	92.1	<DL	50.72	101.44	<DL	47.65	95.3
Sn	<DL	48.16	96.32	<DL	48.33	96.66	<DL	49.8	99.6	<DL	49.45	98.9
Ti	<DL	48.83	97.66	<DL	48.65	97.3	<DL	49.78	99.56	<DL	49.89	99.78
Tl	<DL	46.18	92.36	<DL	46.18	92.36	<DL	47.05	94.1	<DL	46.78	93.56
Zn	<DL	49.71	99.42	<DL	51.25	102.5	<DL	51.13	102.26	<DL	50.3	100.6

\*<DL: measured concentration below method detection limit.

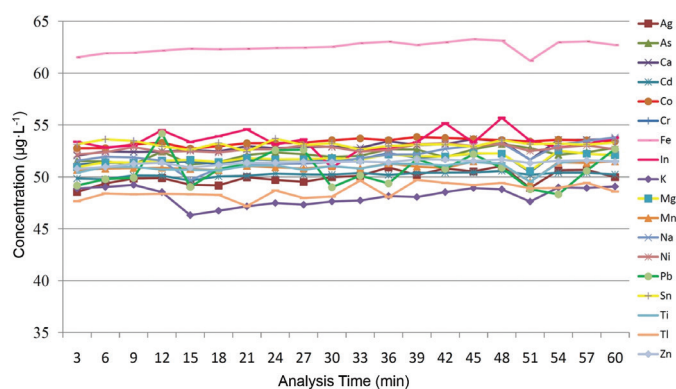


Figure 1. Stability test of spiked  $\text{CuSO}_4$  solution.

## Conclusion

The low relative standard deviations derived from the stability study, combined with the % recovery values of the sample spikes, show the high level of accuracy and precision of analytical results obtained with the iCAP 7400 ICP-OES Duo. The good method detection limits achieved also show that the iCAP 7400 ICP-OES Duo can easily perform the analysis of electronic grade copper sulfate required for electronic factories, semiconductor plants, precious metal technology sites and other electronic grade raw materials testing.

Find out more at [thermofisher.com/ICP-OES](https://thermofisher.com/ICP-OES)

**For Research Use Only. Not for use in diagnostic procedures.** ©2017 Thermo Fisher Scientific Inc. All rights reserved. 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. **AN43178-EN 0717**

**ThermoFisher**  
SCIENTIFIC