

Quantifying Co in Doping Control Urine Samples – A Pilot Study

Daniel Kutscher,¹ Oliver Krug,² Mario Thevis,² Shona McSheehy Ducos¹

¹Thermo Fisher Scientific, Bremen, Germany;

²German Sports University, Cologne, Germany



Overview

Purpose: To demonstrate the power of ICP-MS for the analysis of different elements in biological matrices and demonstrate the use of trace elemental analysis in clinical applications.

Methods: Total elemental determinations of Co and other trace elements in urine were performed using the Thermo Scientific™ iCAP™ Qc ICP-MS in a single, optimized operation mode. The excretion of Co in urine was measured in an elimination study after intake of Co²⁺ and vitamin B12 respectively.

Results: Good agreement between the results determined using a certified reference material was achieved. The elimination study showed that the determination of Co in urine can be a potential tool to screen for its misuse as a performance enhancing substance.

Introduction

The role of Inductively Coupled Plasma Mass Spectrometry (ICP-MS) in clinical research is relatively unknown in the mass spectrometry community. However, it is a powerful tool for trace elemental analysis thanks to high sensitivity and robustness when faced with high matrix samples. ICP-MS provides a fully quantitative technique for almost all elements in the periodic table, offers isotopic information and isotope ratio determinations and can also provide species specific information when coupled with a separation device such as Ion Chromatography (IC) or Liquid Chromatography (LC).

The analysis of trace metals in urine is a way to address potentially harmful exposure to contaminants. One of the advantages for urine as a sample matrix is that the specimen is easily derived. Apart from screening for potential exposure to toxic elements like As, Cd, Hg and Pb, the determination of Co in urine has gained interest, as Co²⁺ could potentially be used as a performance enhancing substance in endurance sports¹.

Methods

Sample Preparation

Urine samples were diluted 10-fold in 2% HNO₃, and 50 µl of an internal standard solution (5 ng·mL⁻¹ Sc, Ge, 2.5 ng·mL⁻¹ Rh and Ir) was added to each sample.

Mass Spectrometry

The iCAP Qc ICP-MS (Figure 1) was used for acquisition of all data.

An SC-4Q autosampler (Elemental Scientific, Omaha, NE, USA) was used for all measurements. The iCAP Qc ICP-MS was operated in single He KED mode for all analyte ions.

All instrumental parameters can be found in table 1.



FIGURE 1. Thermo Scientific iCAP Qc ICP-MS.

TABLE 1. Instrumental Conditions.

| Parameter | Value |
|--------------------|--------------------------------------|
| Nebulizer | PFA-ST |
| Nebulizer Gas Flow | 1.06 L·min ⁻¹ |
| RF Power | 1550 W |
| Interface Set-Up | Ni Cones, High Matrix Skimmer Insert |
| QCell conditions | |
| Cell Gas Flow | 4.8 mL·min ⁻¹ 100% He |
| KED Voltage | 3 V |

Data Analysis

Thermo Scientific™ Qtegra™ Intelligent Scientific Data Solution™ software was used for quantitative assessment of the data.

Results

The underlying mechanism in which Co^{2+} may lead to improved performance of especially endurance athletes is displayed in figure 2. In 2014, Co^{2+} was not listed on the World Anti-Doping Agency's list of prohibited substances², so that dishonest sportsmen could potentially use this option to increase their oxygen transport capacity. In contrast, testing for elevated levels of Co is already done in e.g. horse racing³.

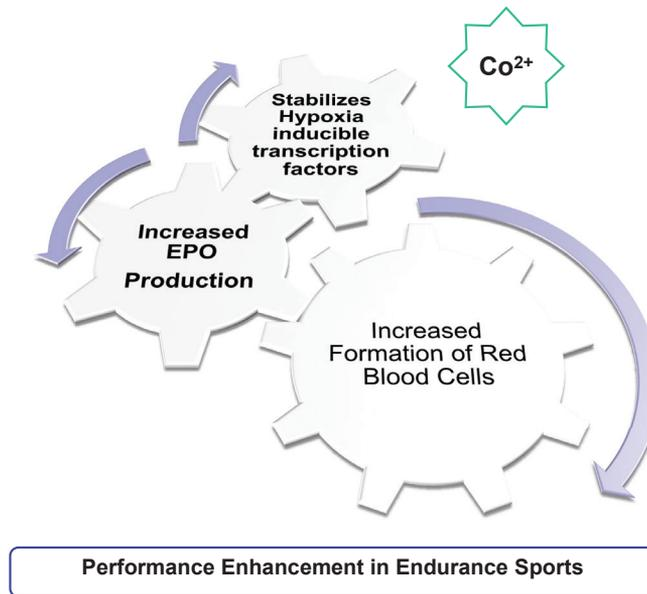


FIGURE 2. Proposed mechanism for Co^{2+} induced performance enhancement in enduring sports.

- The effect of 5 μmol Co is equivalent to stimulation of Erythropoiesis as 1 IU rhEPO.
- Typically, 6-25 $\text{IU}\cdot\text{L}^{-1}$ of EPO are circulating in humans.
- 150 mg CoCl_2 leads to an increase of RBC from 0.5 to 1.19 $\text{Mio cells}\cdot\text{mm}^{-2}$ in 7-22 days.

Analysis of Trace Elements in Urine

In a pilot study, approx. 200 samples were analyzed for Co and various other trace elements in order to verify baseline levels of Co in urine.



Verify baseline Co levels in urine

- 100 healthy non-elite athletes
- 96 endurance sports athletes

- Co concentrations in urine were found to be in the typically observed range [0.1 to $2 \text{ ng}\cdot\text{mL}^{-1}$]

All calibration standards were prepared in diluted nitric acid and a six point calibration curve was generated between 0.025 to 25 ng·L⁻¹ (verification of reference levels) and 0.25 to 100 ng·mL⁻¹ (elimination study). The attainable instrumental limit of detection for ⁵⁹Co was found to be 0.6 ng L⁻¹. UTAK® Urine Control Samples (Low and High Range, UTAK Laboratories, Valencia, CA, USA) were regularly interspersed in the sample list (every 10 samples) to verify the accuracy of the method. The obtained results for the repeated analysis of both control samples within a batch (containing 70 samples) are displayed in table 2.

TABLE 2. Results for the repeated analysis of urine controls as a quality control within one sample batch, all values given in µg·L⁻¹.

| | As | Cd | Ca | Cr | Co | Cu | Fe | Pb | Mn | Mo | Ni | Se | Zn |
|------------------|-------|--------|-----|-------|---------------|--------|-------|--------|-------|--------|-------|------|-------|
| Reference | 16.5 | 0.4 | 96 | 1.3 | 1.9 | 32 | 30 | 0.6 | 3.5 | 52 | 2.6 | 40 | 269 |
| Minimum | 14 | 0.34 | 82 | 1.1 | 1.6 | 27 | 26 | 0.5 | 3.0 | 44 | 2.2 | 34 | 229 |
| Maximum | 19 | 0.46 | 110 | 1.5 | 2.2 | 37 | 35 | 0.7 | 4.0 | 60 | 3.0 | 46 | 309 |
| Recovery | 94% | 95% | 98% | 95% | 95% | 103% | 97% | 122% | 95% | 94% | 117% | 90% | 117% |
| RSD | 1.8 | 6.6 | 1.6 | 4.6 | 3.3 | 2.9 | 4.9 | 4.7 | 2.9 | 2.2 | 3.0 | 2.7 | 2.5 |
| Reference | 118 | 4.6 | 490 | 6.2 | 7.6 | 81 | 565 | 137 | 4.2 | 71 | 27 | 55 | 859 |
| Minimum | 100 | 3.9 | 417 | 5.3 | 6.5 | 69 | 480 | 116 | 3.6 | 60 | 23 | 47 | 730 |
| Maximum | 136 | 5.3 | 564 | 7.1 | 8.7 | 93 | 650 | 158 | 4.8 | 82 | 31 | 63 | 988 |
| Recovery | 94% | 96% | 91% | 105% | 98% | 95% | 87% | 89% | 100% | 95% | 107% | 91% | 100% |
| RSD | 3.6 | 2.9 | 1.7 | 1.9 | 1.6 | 3.2 | 3.1 | 1.0 | 4.1 | 3.8 | 1.7 | 5.2 | 2.8 |
| IDL | 0.007 | 0.0006 | 0.3 | 0.001 | 0.0006 | 0.0008 | 0.002 | 0.0003 | 0.002 | 0.0007 | 0.002 | 0.02 | 0.006 |

Utak 12111 Normal Range

Utak 12110 High Range

Good agreement between the certified and the obtained values was achieved. The low relative standard deviation between the different replicates highlights the robustness of the method.

The obtained results for both groups (mean values 0.28 ng·mL⁻¹ for non elite athletes and 0.36 ng·mL⁻¹ for listed endurance athletes) were found to be within the expected range, but showing a modest but statistically significant increase for professional endurance athletes (Wilcoxon rank sum test, p<0.01). However, strongly increased urinary Co concentrations were not observed.

Feasibility of Co Urinary Concentration as a Tool for Anti Doping Testing

In a second step, an elimination study was performed to verify that intake of Co^{2+} leads to an increase of the urinary Co concentration. The results obtained for both groups are displayed in figure 3.

Investigate excretion after intake of Co^{2+}

- 3 participants. single dose; 500 μg
- 3 participants. multiple doses of 500 μg

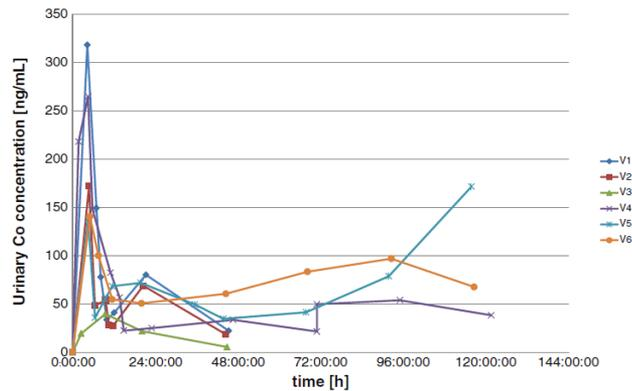


FIGURE 3. Proposed mechanism for Co^{2+} induced performance enhancement in enduring sports.

The results show that the Co concentration in urine is increased in comparison to the reference values found between $0.1\text{-}2.0\text{ ng}\cdot\text{mL}^{-1}$. A similar picture was found after normalization of the determined Co concentration in order to compensate for the specific gravity of the samples. There are two conclusions to be drawn from the results:

- A strong increase of the urinary Co concentration is observed within 6 h (between 40 and $318\text{ ng}\cdot\text{mL}^{-1}$) after intake of Co^{2+} .
- Multiple doses of Co^{2+} lead to elevated levels of Co^{2+} observed even several days after the intake was stopped.

Furthermore, the excretion of Co in urine was monitored upon regular (daily) intake of Vitamin B12 ($500\text{ }\mu\text{g}\cdot\text{day}^{-1}$, accounting for $22\text{ }\mu\text{g}$ Co). However, Vitamin B12 does not seem to affect the urinary Co level.

Conclusion

- The iCAP Q ICP-MS is a perfect tool to determine trace elements in a challenging matrix such as urine. The combination of a unique robust interface design and a single analysis mode for all analytes greatly simplifies method development and helps to improve throughput when it comes to routine analysis of a large number of samples.
- The detection of Co in urine has been shown to be a potential tool to investigate its abuse in sports doping.
- Further investigation is required to elucidate the role of additionally influencing factors and to establish tests for complementary sample matrices like erythrocytes.

References

1. O. Krug et al., Quantifying cobalt in doping control urine samples – a pilot study, *Drug Testing Analysis* **6** (2014), 1186-1190
2. The Prohibited List, published by the World Anti Doping Agency, <https://www.wada-ama.org/en/resources/science-medicine/prohibited-list>
3. P. Bartley: in *The Sidney Morning Herald*, February 4th 2014.

www.thermofisher.com

©2016 Thermo Fisher Scientific Inc. All rights reserved. Utak is the property of Utak Laboratories, Inc. 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 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.

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
Korea +82 2 3420 8600
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