Determination of Dissolved Hexavalent Chromium in Drinking Water, Groundwater and Industrial Wastewater Effluents by Ion Chromatography

Thermo Fisher Scientific Inc.

Introduction

Chromium exists in the environment in either the trivalent, Cr(III), or hexavalent, Cr(VI), form. Cr(III) is considered to be essential to mammals for the maintenance of glucose, lipid, and protein metabolism. On the other hand, Cr(VI) is known to have an adverse affect on the lungs, liver, and kidneys. The method presented here provides a sensitive and selective means of determining Cr(VI) as the chromate anion CrO_4^{2-} down to the 1 um/L level in a variety of environmental matrices.^{1,2}

This method has been documented in several standard methodologies.^{1,3,4,7} It has been validated over the range of 1–1000 µg/L in both wastewater and drinking water⁵ with a method detection limit in the range of 1 µg/L (see Figure 1). It has also been recommended for use in determining Cr(VI) in extracts of ambient air particulates⁶ as well as air particulates from waste incinerators.⁷ For a more complete discussion of this method, please refer to Dionex[™] (now part of Thermo Scientific[™]) Technical Note 26: Determination of Cr(VI) in Water, Wastewater, and Solid Waste Extracts.

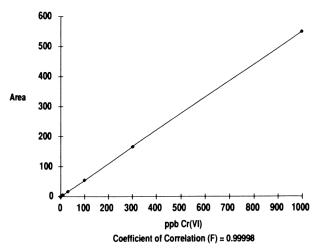


Figure 1. Area response for colorimetric chromium detection.



Equipment

Any Dionex chromatographic system* configured with:

- UV/Visible absorbance detector (VDM-2)
- Postcolumn reagent delivery system

*Equivalent or improved results can be achieved using the Thermo Scientific Dionex ICS-2100 or the Thermo Scientific Dionex ICS-5000⁺ system.

Reagents and Standards

- Ammonium sulfate
- Ammonium hydroxide
- 1,5-Diphenylcarbazide
- Methanol, HPLC grade
- Sulfuric acid



Conditions	
Columns:	Thermo Scientific™ Dionex™ IonPac™ AS7 analytical and Dionex IonPac NG1 guard
Eluent:	250 mM Ammonium sulfate, 100 mM Ammonium hydroxide
Eluent Flow Rate:	1.5 mL/min
Expected Backpressure:	900–1100 psi (6.2–7.6 kPa)
Postcolumn Reagent:	2.0 mM Diphenylcarbazide, 10% methanol, 1.0 N sulfuric acid
Postcolumn Reagent	
Flow Rate:	0.5 mL/min
Detection:	Visible, 530 nm
Sample Volume:	50 μL

Preparation of Solutions and Reagents

Eluent:

250 mM Ammonium sulfate 100 mM ammonium hydroxide

Dissolve 33 g of ammonium sulfate in 500 mL of water and add 6.5 mL of ammonium hydroxide. Dilute to 1 L with water.

Postcolumn Reagent:

2.0 mM Diphenylcarbazide10% methanol1.0 N sulfuric acid

Dissolve 0.50 g of 1,5-diphenylcarbazide in 100 mL of HPLC-grade methanol. Add 500 mL of water containing 28 mL of 98% sulfuric acid while stirring. Dilute with water to 1 L in a volumetric flask. This reagent is stable for four or five days, but should only be prepared in 1 L quantities as needed.

Standard:

1000 ppm Cr(VI)

Dissolve 0.283 g of potassium dichromate ($K_2Cr_2O_7$) that has been dried at 100 °C for one hour in water. Dilute to 100 mL in a volumetric flask.

Working standards are prepared by appropriate dilutions of the stock solutions. As an example, for a 1 ppm Cr(VI) standard, pipet 1.00 mL of the chromium stock solution into a 1 L volumetric flask. Dilute to 1 L with water.

Discussion

An aqueous sample is injected onto a high capacity anion exchange column where Cr(VI) as $CrO_4^{2^-}$ is retained and then eluted with an alkaline sulfate eluent. After this separation, a diphenylcarbazide color reagent is added to the eluent stream, which then flows through a photometric detector. The regent forms a color complex with Cr(VI), which is detected by absorbance at 530 nm. An example of hexavalent chromium in a aqueous sample is shown in Figure 2.

This method is not applicable to the determination of total chromium in environmental samples. For regulatory purposes, total chromium is traditionally defined as the sum of the free and combined chromium from both the liquid and solid portions of the sample. Accordingly, the sample must undergo a strong acid digestion to release the chromium for analysis. This step converts all chromium to the Cr(III) form. Other methods can be used to determine the total chromium in the digest solution.

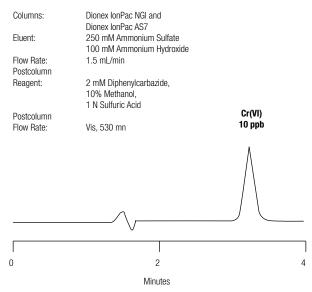


Figure 2. Determination of Cr(VI) in water, wastewater, and solid waste extracts.

References

- Arar, E.J.; Long, S.E.; Pfaff, J.D., "Method 218.6 Determination of Dissolved Hexavalent Chromium in Drinking Water, Groundwater, and Industrial Wastewater Effluents by Ion Chromatography", Nov. 1991, United States Environmental Protection Agency (U.S. EPA), Cincinnati, OH 45268.
- 2. Arar, E.J.; Pfaff, J.D. J. Chromatogr. Sci. 1991, 546, 335-340.
- 3. Proposed ASTM Method: "Dissolved Hexavalent Chromium in Water by Ion Chromatography", June 1991, American Society for Testing and Materials (ASTM), Committee D-19, Philadelphia, PA 19103.
- Proposed ASTM Method: "Collection and Analysis of Hexavalent Chromium in the Atmosphere", Nov. 1991, ASTM, Committee D-22, Philadelphia, PA 19103.

- Edgell, K.; Longbottom, J.; Joyce, R., Final Report: "Determination of Dissolved Hexavalent Chromium in Drinking Water, Groundwater, and Industrial Wastewater Effluents by Ion Chromatography: Collaborative Study", Oct. 1991, U.S. EPA, Cincinnati, OH 45268.
- 6. Pappa, R.; Castillo, N., "Determination of Ambient Levels Of Hexavalent Chromium By Ion Chromatography", California Air Resources Board, El Monte, CA. Presented at the Annual Meeting of the Air and Waste Management Association, Anaheim, CA, June 1989.
- 7. "Determination of Hexavalent Chromium Emissions from Stationary Sources", in U.S. EPA Methods Manual for Compliance With The BIF Regulations: Burning Hazardous Waste in Boilers and Industrial Furnaces, U.S. EPA/53-SW-91-010, NTIS, Springfield, VA 22161, Dec. 1990.

www.thermoscientific.com/chromatography

©2014 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. 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 Brazil +55 11 3731 5140 Canada +1 800 530 8447 China 800 810 5118 (ree call domestic) 400 650 5118 AN71358-EN 1014M 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 6 6885 1213 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 Sweden +46 8 556 468 00 Switzerland +41 61 716 77 00 Taiwan +886 2 8751 6655 UK/Ireland +44 1442 233555 USA +1 800 532 4752



A Thermo Fisher Scientific Brand