# Determination of Hexavalent Chromium in Dyes

Thunyarat Phesatcha,<sup>1</sup> Weerapong Worawirunwong,<sup>1</sup> Jeff Rohrer<sup>2</sup> <sup>1</sup>Thermo Fisher Scientific, Bangkok, Thailand; <sup>2</sup>Thermo Fisher Scientific, Sunnyvale, CA, USA

## Introduction

Hexavalent chromium [Cr(VI)] is the most toxic form of the metal chromium. In addition to regulations concerning its concentration in drinking water, many governments regulate Cr(VI) in consumer products. Leather products can be colored with dyes containing trivalent chromium, and sweat can transfer the chromium from the leather to the skin. It is possible that some of the trivalent chromium oxidizes to Cr(VI). For example, leather gloves were found to contain Cr(VI).<sup>1</sup> In this application brief, we analyze two dyes for Cr(VI) using a sensitive ion chromatography (IC) method with postcolumn reaction to produce a colored complex detected by visible absorbance.<sup>2</sup> Figure 1 shows that Cr(VI) was not found in a 1:100 dilution of the Fast Red dye while 56.1 µg/L Cr(VI) was found in a 1:100 dilution of the Metal Complex Black. Spiking both diluted dye samples with 30 µg/L Cr(VI) yielded recoveries of 99.3 and 101% for the red and black dyes, respectively, demonstrating method accuracy. This IC method delivers an easy, fast, sensitive, and accurate determination of Cr(VI) in dyes.

#### Conditions

- System Thermo Scientific<sup>™</sup> Dionex<sup>™</sup> ICS-3000 with a DP dual pump, VWD absorbance detector, and AS autosampler
- Chromatography See Figure 1.
- Sample Preparation Samples were diluted 1:100 with deionized water and then passed through a Thermo Scientific Dionex OnGuard<sup>™</sup> II P cartridge (P/N 057087) prior to analysis.



Figure 1. Chromium [Cr(VI)] in dyes.



## References

- 1. Peterson, J.; Murphy, B.; Perati, P.; Richter, B. LCGC *The Applications Notebook*, **2007**, June, 28.
- 2. Dionex (now part of Thermo Scientific) Application Update 144. *Determination of Hexavalent Chromium in Drinking Water Using Ion Chromatography*. Dionex LPN 1495, 2003, Sunnyvale, CA.

# **Acknowledgement**

Thanks to the Textile Testing Center, Thailand Textile Institute for kindly providing the samples.

www.thermoscientific.com/dionex

©2012 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 Inc. and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific Inc. 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.

Australia +61 3 9757 4486 Austria +43 1 333 50 34 0 Belgium +32 53 73 42 41 Brazil +55 11 3731 5140 China +852 2428 3282 Japan +81 6 6885 1213 Korea +82 2 3420 8600 Netherlands +31 76 579 55 55 Singapore +65 6289 1190 Sweden +46 8 473 3380



Switzerland +41 62 205 9966 Taiwan +886 2 8751 6655 UK/Ireland +44 1442 233555 USA and Canada +847 295 7500

