# Measurement of Dietary Carotenoid Isomers Using HPLC-ECD

Paul Gamache and Ian Acworth Thermo Fisher Scientific, Chelmsford, MA, USA

Key Words Vitamin A, Carotenoids, HPLC-ECD, Tissue

### **Summary**

This application note describes an HPLC-ECD method using a C30 stationary phase column for the analysis of isomeric forms of carotenoids in animal and plant tissues.

### Introduction

Much of the interest in dietary carotenoids, exclusive of their pro-vitamin A activity, is related to their possible actions as preventive agents in diseases associated with oxidative stress. These electron-rich compounds can act as antioxidants *in vitro* and their possible role of protection from reactive oxygen and nitrogen species *in vivo* has received much attention.<sup>1</sup>

Dietary carotenoids can each be found as a variable mixture of geometric and positional isomers (e.g., all *trans, cis*-9-, and *cis*-13) – See Figure 1. These isomers may occur naturally or can be formed during processing<sup>2</sup> and show a variety of biological properties and chemical activities. For example,  $\beta$ -carotene isomers exhibit significant differences in pro-vitamin A activity,<sup>3</sup> tissue distribution<sup>4</sup> and detector responsivity.<sup>5</sup>

Many reversed phase HPLC methods that utilize monomeric bonded C18 stationary phases for carotenoid analysis are incapable of completely separating the various isomers and in some instances, may lead to inaccurate assessment of nutritional and other healthrelated properties.<sup>6,7,8</sup> HPLC methods that utilize a silica-based C30 stationary phase specifically designed for carotenoid separations have shown significant, and often superior, overall enhancement in shape selectivity for a variety of carotenoid isomers.<sup>5,6,7</sup> Photodiode array (PDA) detection in combination with C30 separations represents an extremely useful technique for carotenoid analysis. This approach, however, has limited sensitivity when analyzing the lower levels present in serum and animal



tissue, particularly for the less abundant isomers.<sup>7</sup> Electrochemical detection (ECD) is a highly sensitive form of HPLC detection for lipid soluble vitamins and antioxidants.<sup>9,10,11</sup> The coulometric array is a multichannel form of electrochemical detection that allows resolution of compounds based on differences in voltammetric behavior and qualitative characterization of peaks even at trace levels.<sup>12,13,14</sup> The basis of detection and resolution using coulometric array relates to differences between analytes in their ability to delocalize charge. Since the antioxidant properties of a compound are closely related to these same structural characteristics, the application of coulometric array detection to carotenoid analysis may provide some additional biochemical insight.<sup>15</sup>

This application note examines the use of coulometric array detection with HPLC using a C30 stationary phase for analysis of isomeric forms of carotenoids in animal and plant tissues. HPLC coulometric array ECD can be used both as a complementary approach to traditional methods and to provide unique information on carotenoids.







Figure 1. The structure of  $\alpha\text{-}$  and  $\beta\text{-}carotene$  and two geomrtic isomers of  $\beta\text{-}Carotene.$ 

## **Materials and Methods**

The isocratic analytical system consisted of a pump, an autosampler, a thermostatic chamber, a twelve-channel Thermo Scientific<sup>™</sup> Dionex<sup>™</sup> CoulArray<sup>™</sup> Coulometric Array Detector and a UV/vis detector placed prior to the array.



Figure 2. Analysis of a raw carrot extract.

Conditions	
LC	
Column:	C30, 5 μm, 4.6 × 250 mm
Mobile Phase:	Methanol – Methyl- <i>ter</i> t-butyl Ether (MTBE) – 1.0M Ammonium Acetate, pH 4.4 (63:35:2) (v/v/v)
Flow Rate:	1.0 mL/min
Temperature:	28 °C
Injection Volume:	10 µL
Detector	
Electrochemical Detector:	Model 5600A, CoulArray
Applied Potentials:	100, 160, 220, 280, 340, 400, 460 and 520 mV vs. Pd.
Detector Wavelength:	450 nm

#### **Standards**

Standards were prepared by dissolving ~1 mg/10 mL chloroform followed by dilution in 10 mL ethanol. Solutions were assayed spectrophotometrically and assigned a concentration value based on molar absorbtivity. Butylated hydroxyanisole was then added as a preservative. Dilutions were made in mobile phase.

# Sample Preparation

A 0.5 mL volume of serum or standard was mixed with 0.5 mL ethanol/10 mg/L BHA. After mixing for 1 min, 1.5 mL of hexane was added and after mixing for an additional 10 min. was centrifuged (4,000 rpm, 10 min). Approximately 1.0 mL of supernatant was withdrawn and the remaining sample extracted with an additional 1.5 mL of hexane. Combined hexane extracts were evaporated to dryness under a stream of nitrogen. Finally, the residue was dissolved in 0.25 mL of mobile phase.

# **Results and Discussion**

The separation of carotenoid isomers was completed within 25 minutes (Figure 2). For coulometric array ECD, the assay was linear over the range of 0.1 to 500ng on column. The limit of detection (lod) was ~20 pg (s/n 3:1) for both trans- $\alpha$ -carotene and trans- $\beta$ -carotene; a twenty fold improvement over UV detection. The within-run peak height response variability (% relative standard deviation, %R.S.D.) was 0.42 and 0.45 for 10ng of trans- $\alpha$  and trans- $\beta$ -carotene, respectively (n=8). For 200 pg, an amount well below the LOD for absorbance detection, EC response variability was 3.57 and 6.45 %R.S.D. for trans- $\alpha$  and trans- $\beta$ -carotene, respectively (n=7).

Figure 3 illustrates HPLC-coulometric array profiles of MTBE extracts of a raw (Figure 3A) and a thermally processed carrot (Figure 3B). Peak elution profiles are similar to those commonly obtained using absorbance detection where isomeric forms of  $\alpha$ - and  $\beta$ -carotene are apparent as a result of thermal processing. Figure 4 shows a representative chromatogram of a human serum extract. Levels of trans isomers, estimated using external standards, showed good agreement between EC and absorbance  $(\pm 3\%)$ . The output from 4 electrochemical channels (to illustrate voltammetric response ratios) along with absorbance at 450 nm are shown. Response ratios between adjacent channels when compared to those obtained from standards (Figure 2) are in close agreement. This is indicative of high peak purity for both *trans*- $\alpha$ and trans-\beta-carotene. The chromatogram in Figure 4 illustrates the ability to obtain useful qualitative information on low level analytes in a complex matrix. While the ability to detect and obtain spectral absorbance information is very limited, the EC response for a number of putative isomers is well above the lower limit of detection.

It is interesting to note that in Figures 3 and 4 the predominant EC response is obtained at 340mV for *trans*- $\beta$ -carotene and 400 mV for *trans*- $\alpha$ -carotene (i.e. *trans*- $\beta$ -carotene is more easily oxidized due to more effective hyper-conjugation). Furthermore, a peak eluting immediately prior to *trans*- $\alpha$ -carotene, which can be tentatively identified as 13-*cis*- $\beta$ -carotene, has dominant response at the same channel as *trans*- $\beta$ -carotene.<sup>7</sup> Other possible *cis* isomers, based on literature, may include the two unresolved peaks immediately prior to '13-*cis*- $\beta$ -carotene' and possibly the 9-*cis*- $\beta$ -carotene which elutes immediately after  $\beta$ -carotene.<sup>7</sup> This latter peak has an unique voltammetric profile with dominant response at 460 mV.

# Conclusion

In conclusion, HPLC with a silica-based C30 stationary phase in combination with coulometric array electrochemical detection allowed selective measurement of carotenoids with an estimated 20-fold improvement in sensitivity (lod ca. 20 pg) as compared to UV-Vis absorbance detection. Response was linear over a 0.1 to 500 ng range. Qualitative voltammetric data obtained across serial coulometric sensors was used to aid in peak identification and provide information on the oxidationreduction properties of analytes. The feasibility of detection and qualitative analysis of trace levels of *cis* isomeric forms of  $\alpha$  and  $\beta$ -carotene in human serum has been demonstrated. This approach has recently been applied to measurement of isomeric forms of carotenoids in tissue.<sup>16</sup>



Figure 3. Analysis of a thermally processed carrot. (\*tentative indentification).



Figure 4. Low level serum.

# References

- Krinsky, N. (1993). Actions of carotenoids in biological systems. Annu. Rev. Nutr., 13, 561-587.
- 2. Chandler, L. A., and Schwartz, S. J. (1988). HPLC separation of cis-trans carotene isomers in fresh and processed fruits and vegetables. *J. Food Sci.*, **52**, 669-672.
- Sweeney J. P., and Marsh, A. C. (1971). Effect of processing on provitamin A in vegetables. *J. Amer. Diet. Assoc.*, 59, 238-243.
- 4. Mokady, S., Avron, M., and Ben-Amotz, A (1990). Accumulation in chick livers of *9-cis* versus all-trans beta-carotene. *J. Nutr.*, **120**, 889-892.
- Bell, C. M., Sander L., C., and Wise, S. A. (1997). Temperature dependence of carotenoids on C18, C30, and C34 bonded stationary phases. *J. Chromatogr. A.*, 757, 29-39.
- Emenhiser, C., Sander, L. C., and Schwartz S. J. (1995). Capability of a polymeric C30 stationary phase to resolve cis-trans carotenoid isomers in reversed-phase liquid chromatography. *J Chromatogr. A.*, 707, 205-216.
- Sharpless, K. E., Thomas, J. B., Sander L. C., and Wise, S. A. (1996). Liquid chromatographic determination of carotenoids in human serum using an engineered C30 and a C18 stationary phase. *J. Chromatogr. B*, 678, 187-195.
- Sowell, A. L., Huff, D.L., Gunter, E. W., and Driskell, W. J. (1988). Identification of cis-carotenoids in human sera analyzed by reversed- phase high-performance liquid chromatography with diode array detection. J. Chromatogr., 431, 424-430.
- Chou, P. P., Jaynes, P. K., and Bailey, J. L. (1985). Determination of vitamin E in microsamples of serum by liquid chromatography with electrochemical detection. *Clin. Chem.*, 31, 880-882.
- Edlund, P. O. (1988). Determination of coenzyme Q10, α-tocopherol and cholesterol in biological samples by coupled-column liquid chromatography with coulometric and ultraviolet detection. *J. Chromatogr.*, 425, 87-97.
- Haroon, Y., Schubert, C. A. W., and Hauschka, P. V. (1984). Liquid chromatographic dual electrode detection system for vitamin K compounds. *J. Chromatogr. Sci.*, 22, 89-93.

- Acworth, I. N., and Gamache, P. (1996). The coulometric electrode array for use in HPLC analysis. Part 1: Theory. *American Lab.*, 5, 33-38.
- 13. Acworth, I. N., Waraska, J., and Gamache, P. (1997). The coulometric electrode array for use in HPLC analysis. Part 2: An application overview. *American Lab.*, **11**, 25-32.
- Matson, W. R., Langlais, P., Volicer, L., Gamache, P., Bird, E., and Mark, K. (1984). n-Electrode threedimensional liquid chromatography with electrochemical detection determinations of neurotransmitters. *Clin. Chem.*, 30, 1477-1488.
- Gamache, P. H., Freeto, S. M., and Acworth I.N. (1999). Coulometric array HPLC analysis of lipidsoluble vitamins and antioxidants. *Am. Clin. Lab.*, 8.
- Ferruzzi, M. G., Sander, L. C., Rock, C. L., and Schwartz, S. J. (1998). Carotenoid determination in biological microsamples using liquid chromatography with a coulometric electrochemical array. *Anal. Biochem.*, 256, 74-81.
- Epler, K. S., Zeigler, R. G., and Craft, N. E. (1993). Liquid chromatographic method for the determination of carotenoids, retinoids and tocopherols in human serum and in food. *J. Chromatogr.*, 619, 37-48.

# Acknowledgements

ESA, Inc., now part of Thermo Fisher Scientific, is grateful to Dr. S. J. Schwartz, M. G. Ferruzzi, and M. L. Nguyen Dept. Food Science and Technology, Ohio State University, for their collaboration in this study.

# **Ordering Information**

Description	Part Number
HPG-3400RS Biocompatible Binary Rapid Separation Pump with Two Solvent Selector Valves	5040.0046
WPS-3000TBRS Biocompatible Rapid Separation	5841.0020
CoulArray, Model 5600A - 8 channel	70-4325
CoulArray Organizer with Temp. Control	70-4340T
Accessory Kit, CoulArray Detector to UltiMate 3000 System	70-9191

# www.thermofisher.com/liquidchromatography

©2016 Thermo Fisher Scientific Inc. All rights reserved. All 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 (free call domestic) 400 650 5118

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