

# The Spectro-Electro Array: A Novel Platform for the Measurement of Secondary Metabolites in Botanicals, Supplements, Foods and Beverages - Part 4: Beer Polyphenols, Proanthocyanidins and Bitter Acids

*Paul A. Ullucci, Ian N. Acworth, Marc Plante, Bruce A. Bailey, and Christopher Crafts  
Thermo Fisher Scientific, Chelmsford, MA, USA*

## Overview

**Purpose:** To develop gradient HPLC methods using a spectro-electro array platform for use to either measure specific analytes in beer samples or in a metabolomic approach to distinguish between different beer samples, as well as study beer stability.

**Methods:** Gradient HPLC with diode-array detection and electrochemical array detection were used.

**Results:** A wide variety of polyphenols and bitter acids could easily be identified and quantified and their stability measured using the bitter acid targeted method. The metabolomic approach could readily distinguish between different beer types when using EC data, but less so when using UV data.

## Introduction

Beer is the most widely consumed alcoholic beverage in the world and the third most popular drink after water and tea. Beer is typically brewed from four basic ingredients: water, a starch source (e.g., malted barley), brewer's yeast, and a flavoring agent such as hops. Many varieties of beer result from differences in these ingredients, the additives used and the brewing process. Beer contains a complex mixture of phenolic compounds extracted from starch and hops. A recent trend from microbreweries in the USA is the development of a wide range of extremely bitter beers created by the addition of extra hops during the brewing process. The hop-derived xanthohumol and the iso-alpha acids formed are primarily responsible for the perceived bitterness. Many of these secondary metabolites are not only purported to offer health benefits but are essential to the flavor and stability of the beer itself. Conversely, some secondary metabolites contribute to the degradation of beer during storage with the formation of haze (e.g., catechins and their polymers the proanthocyanidins).

Presented here are gradient HPLC Spectro-Electro Array methods for studying beer composition. Two targeted assays were developed for the measurement of either polyphenols (including catechins and proanthocyanidins) or xanthohumols and iso-alpha acids. The bitter acid method was used to study beer stability. Also discussed is a metabolomics approach where patterns of both known and unknown analytes can be used to study differences between beer samples—an approach that is relevant to quality control.

## Methods

### Liquid Chromatography Method 1: Polyphenols

Pump: Thermo Scientific Dionex LPG-3400BM with SR-3000 Solvent Rack  
Autosampler: Thermo Scientific Dionex WPS-3000TBSL  
UV Detector: Thermo Scientific Dionex DAD-3000RS diode-array detector  
Channel 1: 218 nm Channel 2: 240 nm  
Channel 3: 254 nm Channel 4: 275 nm  
EC Detector: Thermo Scientific Dionex CoulArray Detector with Thermal Organizer  
EC Parameters: 16 channel array from 0 to +900 mV in +60 mV increments  
Mobile Phase A: 20 mM monobasic sodium phosphate, 3% acetonitrile, 0.2% tetrahydrofuran, pH 3.35  
Mobile Phase B: 20 mM monobasic sodium phosphate, 50% acetonitrile, 10% tetrahydrofuran, pH 3.45  
Mobile Phase C: 90% methanol  
Gradient: 0-2 min: 2%B/3%C. 30 min: 97%B/3%C, Concave. 45 min 97%B/3%C  
Column: Thermo Scientific Acclaim 120, C18, 3  $\mu$ m, 3.0  $\times$  150 mm  
Flow Rate: 0.65 mL/min  
Injection: 10 or 20  $\mu$ L

### Liquid Chromatography Method 2: Bitter Acids

Column: Acclaim™ 120, C18, 3  $\mu$ m, 3.0  $\times$  150 mm,  
Flow Rate: 0.65 mL/min  
Injection: 20  $\mu$ L  
Mobile Phase A: 25 mM sodium perchlorate, 50% acetonitrile, 2.5 mM perchloric acid  
Mobile Phase B: 25 mM sodium perchlorate, 90% acetonitrile, 2.5 mM perchloric acid  
Mobile Phase C: 90% methanol  
Gradient: 0-3 min: 0%B/3%C. 30 min: 40%B/3%C  
40 min: 97%B/3%C. 45 min: 97%B/3%C  
EC Parameters: Thermo Scientific Dionex model 5011A Dual Channel Cell  
E1: +550 mV E2: +850 mV

### Data Analysis

Data were analyzed using Thermo Fisher Dionex Chromeleon Chromatography Data System 6.8 (SR 9) and CoulArray™ software 3.1. EC-array data were transferred to Pirouette® software for chemometric analysis using a CoulArray version 2.0 Software Utility (Pattern Recognition Setup Wizard). UV data were tabularized prior to transfer to Pirouette.

### Standard Preparation

For the polyphenol method, standards, depending upon solubility, were prepared in ethanol, methanol, or methanol/water solutions at 1000 or 100 µg/mL. Working standards were prepared at 0.20, 0.50, and 1.0 µg/mL in 10% methanol containing 0.2% ascorbic acid/0.02% EDTA.

For the bitter acids method, standard mixes were obtained from the American Society of Brewing Chemists: International Calibration Standards for the HPLC Analyses of Isomerized alpha-acids including: DCHA-Iso, ICS-12; DCHA-Rho, ICS-R2; Tetra, ICS-T2 and DEHA-Hexa, ICS-H1. All acids were obtained as their dicyclohexylamine salt. Stock standard solutions were prepared by dissolving 100 mg of the standard material with 100 mL of acidified methanol (0.2% phosphoric acid in methanol) and sonicating for 15 mins. Calibration Standards were prepared in 50% acetonitrile containing 0.2% phosphoric acid in the range of 0.10– 2.0 µg/mL.

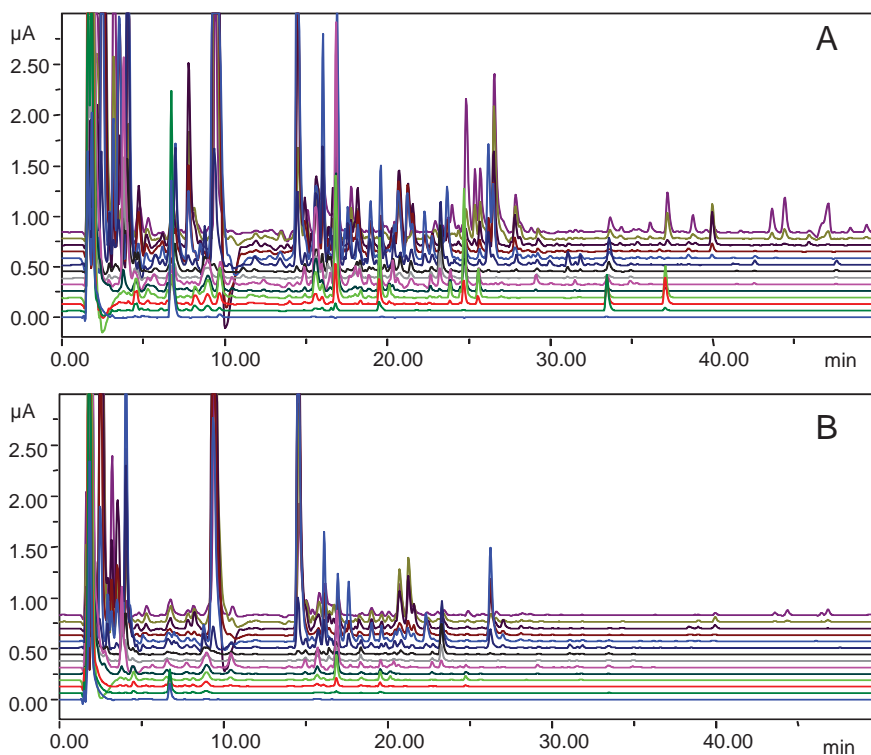
### Sample Preparation

A random variety of beer samples were obtained from the local liquor store and included domestic beer and a light equivalent, numerous domestic microbrews, European examples (Bavaria and Belgium), and a domestic extremely bitter (highly-hopped) beer. Beer samples were treated as follows: 0.50 mL of beer + 0.50 mL of acidified acetonitrile (0.2% phosphoric acid in acetonitrile) were mixed, centrifuged, and the clear supernatant analyzed. For the stability study, beer samples were transferred to a sealed container and kept at 4 °C in the dark, and processed as needed.

## Results and Discussion

The Spectro-Electro Array makes use of both spectrophotometric and electrochemical data.<sup>1</sup> While UV data provides identification and quantitation of the major components in a sample, EC array detection provides additional information. First, the EC array is incredibly sensitive with low pg limits of detection (LOD) and it is capable of measuring compounds missed by UV. Second, it voltammetrically resolves compounds that co-elute chromatographically. Third, the EC array is fully gradient compatible, thereby extending the number of analytes that can be measured in a sample. Fourth, the redox behavior of a compound reacting across the array provides qualitative information and can be used for analyte identification/authentication.

**FIGURE 1. A. Polyphenol Method Chromatogram of a High-hop Beer and B. Regular Domestic Beer.**



### Polyphenol Method – Targeted Analysis

The analytical figures of merit for this assay were described previously.<sup>1</sup> Briefly, the limits of detection were typically 10–50 pg on column by ECD and 100–500 pg by UV. The limits of quantification were 200–1000 pg on column by ECD and 500–5000 pg by UV. Response range was over seven orders of magnitude by ECD and five by UV. Typical R<sup>2</sup> values were ~0.99 or better for all compounds. Average intra-day retention time precision for all analytes averaged 0.55% RSD over a 10 day period, with a range of 0.30–1.22%. Multi-channel EC array chromatograms for two different beer samples are presented in Figure 1A with high-hops beer and 1B with regular domestic beer. As can be seen, the high-hops beer contains a great abundance of analytes, as confirmed in Table 1. Analytes are in agreement with previously published data.

**TABLE 1. Polyphenol Data Presented in mg/L. Beer 1 – High-hop; Beer 3 – Regular.**

Compound	Beer 1	Beer 3	Compound	Beer 1	Beer 3
4-Hydroxybenzyl alcohol	2.32	ND	Ferulic acid	1.42	2.54
2-Hydroxybenzyl alcohol	0.04	ND	Gentisic acid	0.24	0.06
3,4-Dihydroxybenzoic acid	2.10	ND	Hesperidin	0.40	0.04
4-Hydroxybenzaldehyde	0.16	ND	Isorhamnetin	0.94	ND
4-Hydroxybenzoic acid	0.38	0.36	Isoxanthohumol	1.16	0.14
4-Hydroxyphenyl acetic acid	0.14	0.24	Kaempferol	1.04	ND
4-Hydroxycoumarin	11	0.46	Myricetin	0.5	0.2
Apigenin	0.2	0.02	Naringin	2.08	5.1
Caffeic acid	0.14	0.1	p-Coumaric acid	2.02	2.8
Carnosol	0.38	ND	Quercetin	1.56	ND
Catechin	4.48	1.8	Salicylic acid	0.18	0.04
Cavarcrol	1.2	0.25	Sinapic acid	0.68	0.58
Chlorogenic acid	0.52	0.04	Syringaldehyde	1.39	ND
Chrysin	0.12	0.26	Syringic acid	0.08	0.1
Epicatechin	1.76	0.40	Thymol	0.78	0.09
Epicatechingallate	1.54	0.26	Umbelliferone	0.34	0.64
Epigallocatechin	0.14	0.16	Vanillic acid	0.22	0.15
Ellagic acid	1.38	0.64	Vanillin	1.06	0.26
Ethyl vanillin	0.2	0.02	Xanthohumol	0.26	0.04

### Polyphenol Method – Metabolomic Study

A simple metabolomics experiment was conducted to evaluate whether the Spectro-Electro Array platform could be used to differentiate between different beer types including: matched domestic normal and light beers, a variety of domestic microbrews, two European beers (from Bavaria and Belgium), an Irish stout, and a domestic Ultra (extra hopped) IPA. Metabolomic profiles containing several hundred analytes, including both known (e.g., see Table 1) and unknown compounds, were measured in each sample. Principle component analysis (PCA) was then used to differentiate samples for both EC data (Figure 2A) and UV data (Figure 2B). The EC data gave the best differentiation between samples and could distinguish between the light and normal beer, Irish stout, Belgian beer and the extra hopped IPA. UV data were less effective with no ability to distinguish between light and normal beers.

### Bitter Acid Method – Targeted Analysis

This method showed similar figures of merit to the polyphenol method above. Rather than using an EC array approach, a simple two channel cell was used to detect xanthohumol and the alpha and beta bitter acids on the first lower potential channel, and isoxanthohumol, prenylnaringenin, and the cis- and trans-iso bitter acids on the second higher potential channel. Example chromatograms are shown in Figure 3A (standards) and Figure 3B (extra hopped IPA). As expected, the Ultra IPA beers were abundant in the bitter acids and related compounds, more so than matched normal and light beers (Table 2).

FIGURE 2. Principal Component Plots for A) ECD and B) UV Data.

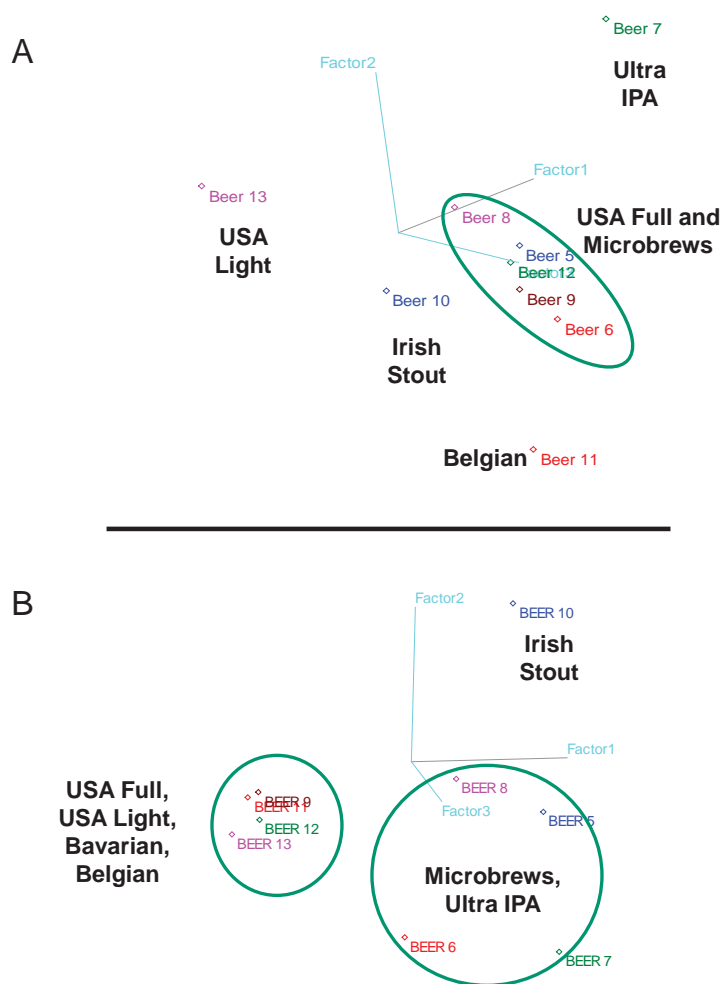
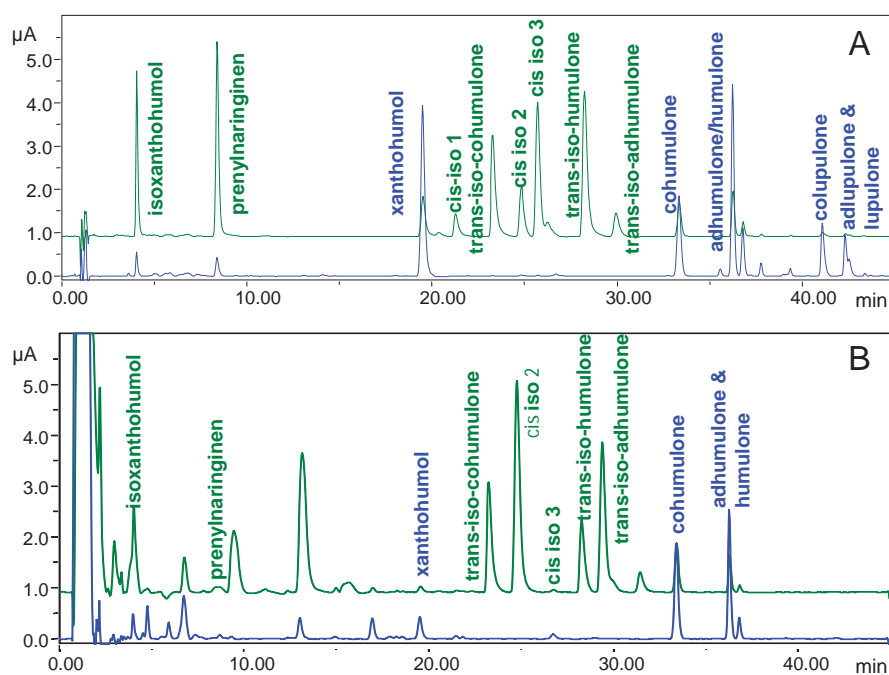


FIGURE 3. Chromatograms of A. Bitter Acids Standard Mixture; B. Ultra IPA Beer Sample. Blue Trace at 550 mV, Green Trace at 850 mV.



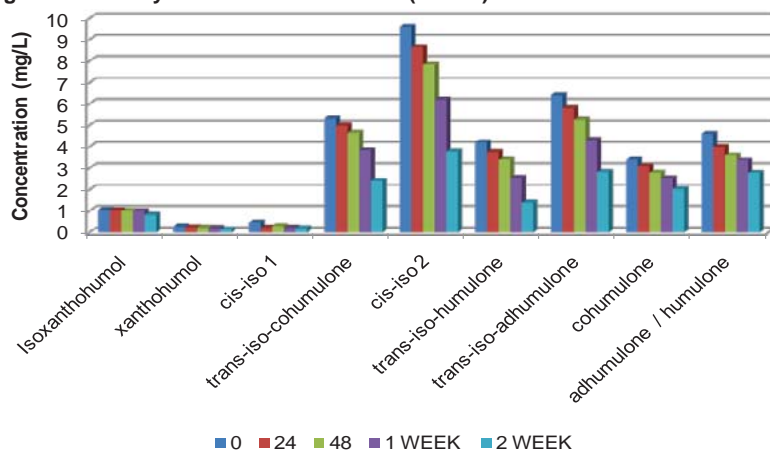
**TABLE 2. Hops Bitter Acids Data Presented in mg/L.**

Compound	Beer 1 Ultra IPA	Beer 2 Ultra IPA	Beer 3 Regular	Beer 4 Light Beer
Isoxanthohumol	2.10	1.3	0.38	0.28
Xanthohumol	0.52	0.48	ND	ND
Cis-iso-acid 1	0.90	0.4	ND	ND
Trans-iso-cohumulone	10.6	7.0	ND	ND
Cis-iso-acid 2	19.1	12.2	3.8	1.6
Trans-iso-humulone	8.4	7.6	0.20	ND
Trans-iso-adhumulone	12.8	11.0	3.2	2.6
Cohumulone	6.8	6.4	0.03	0.02
Adhumulone\Humulone	9.2	9.0	ND	ND

### Beer Stability Study

Auto-oxidation, including decomposition of iso-alpha-acids, plays an important role in the deterioration of taste and flavor qualities of beer during aging. It is believed that degradation of iso-alpha-acids is the cause of gradual decrease of beer bitterness. The current bitter acid method was used to evaluate the stability of a selection of the beers. Data for the stability of a variety of analytes in an Ultra IPA sample tested over a two week period are presented in Figure 4. Marked decreases in many analytes was seen but particularly for cis-iso 2 (-61%), trans-iso-humulone (-67%) , trans-iso-adhumulone (-56%), and trans-iso-cohumulone (-55%). This data is consistent with previous studies.<sup>2</sup>

**Figure 4. Stability Data for Ultra IPA Beer (Beer 1).**



## Conclusions

- The polyphenol method can be used in a targeted approach to accurately and sensitively measure numerous phenols, phenolic acids and polyphenols in beer and other samples not possible by methods employing UV detection alone.
- Metabolomic approaches using the patterns of numerous known and unknown analytes can be used to differentiate between different samples. Such an approach can be used to study fermentation, product stability, and authenticity.
- The bitter acid method enabled the sensitive targeted measurement of isoxanthohumol, xanthohumol, prenylnaringenin, the trans- and cis-iso-alpha bitter acids and the alpha and beta bitter acids in a single run. Use of EC detection eliminated the need for solid phase extraction procedures for sample pre-concentration commonly used in UV detection methods. This approach was used to measure beer stability over a two week period.

## References

1. Ullucci, P.A.; Bailey, B.; Acworth, I.N.; Crafts, C.; Plante, M. The Spectro-Electro Array: A Novel Platform for the Measurement of Secondary Metabolites in Botanicals, Supplements, Food, and Beverages. Part 1. Theory and Concepts, Pittsburgh Conference, **2012**, Poster 1770-1 P.
2. De Cooman, L.; Aerts, G.; Overmeire, H. Alteration of the Profiles of Iso-alpha-Acids During Beer Ageing, Marked Instability of Trans-Iso-alpha-Acids and Implications for Beer Bitterness, Consistency in Relation to Tetrahydroiso-alpha-acids , *Journal of the Institute of Brewing*. **2000**, 106: 169-178.

[www.thermofisher.com/dionex](http://www.thermofisher.com/dionex)

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

---

<b>U.S./Canada</b>	(847) 295 7500	<b>Denmark</b>	(45) 36 36 90 90	<b>Sweden</b>	(46) 8 473 3380	<b>India</b>	(91) 22 2764 2735
<b>Brazil</b>	(55) 11 3731 5140	<b>France</b>	(33) 1 39 30 01 10	<b>Switzerland</b>	(41) 62 205 9966	<b>Japan</b>	(81) 6 6885 1213
<b>Austria</b>	(43) 1 616 51 25	<b>Germany</b>	(49) 6126 991 0	<b>United Kingdom</b>	(44) 1276 691722	<b>Korea</b>	(82) 2 2653 2580
<b>Benelux</b>	(31) 20 683 9768	<b>Ireland</b>	(353) 1 644 0064	<b>Australia</b>	(61) 2 9420 5233	<b>Singapore</b>	(65) 6289 1190
	(32) 3 353 42 94	<b>Italy</b>	(39) 02 51 62 1267	<b>China</b>	(852) 2428 3282	<b>Taiwan</b>	(886) 2 8751 6655

**Thermo**  
S C I E N T I F I C

Part of Thermo Fisher Scientific

PN70045\_E 07/16S