# Novel, Universal Approach for the Measurement of Natural Products in a Variety of Botanicals and Supplements

Ian N. Acworth<sup>1</sup>, Bruce A. Bailey<sup>1</sup>, Marc Plante<sup>1</sup>, Christopher Crafts<sup>1</sup>, David Thomas<sup>1</sup> and Mark C. Roman<sup>2</sup> <sup>1</sup>Thermo Fisher Scientific, Chelmsford, MA, USA and <sup>2</sup>Tampa Bay Analytical, Largo, FL, USA





### **Overview**

**Purpose:** To evaluate the application of HPLC with charged aerosol detection to the measurement of natural products and botanicals.

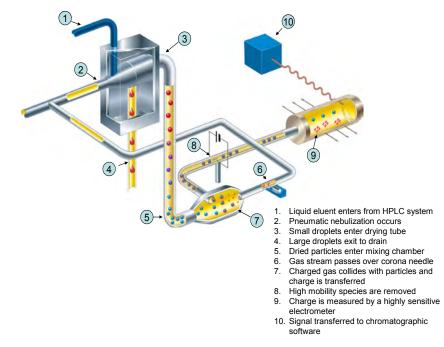
**Methods:** Several methods were developed for the analysis of black cohosh, ginkgo, ginseng, milk thistle, gotu kola and stevia.

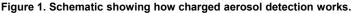
**Results:** HPLC with charged aerosol detection is sensitive (low ng levels on-column), has a wide dynamic range and minimal inter-analyte response variability. It is ideal for measuring analytes that lack a chromophore and offers superior performance over evaporative light scattering detection (ELSD) and refractive index (RI).

### Introduction

Botanicals contain a great diversity of compounds that can show extreme variation in their physicochemical properties. Analysis of potentially active components can be challenging as not all contain a chromophore or can be ionized, thereby limiting the use of UV absorbance and mass spectrometry, respectively. HPLC with charged aerosol detection is a sensitive, universal (nonselective) approach that can measure any nonvolatile and many semi-volatile compounds. A number of isocratic and gradient HPLC with the Thermo Scientific Dionex Corona Charged Aerosol Detector (CAD™) methods were developed and evaluated for the measurement of phytochemicals extracted from a variety of botanicals including: triterpene glycosides from black cohosh (Cimicifuga racemosa); ginkgolides and bilobalides from ginkgo (Ginkgo biloba); ginsenosides from ginseng (Panax ginseng); silibinins in milk thistle (Silybum marianum): ursane and oleanane triterpenes from gotu kola (Centella asiatica): and diterpene glycosides from stevia (Stevia rebaudiana). Analytes showed consistent response independent of chemical structure (typically < 10% variability between compounds corrected for gradient elution). All methods had a wide dynamic range (~four orders of magnitude), good sensitivity (typically low ng levels of detection), and excellent reproducibility (RSDs typically < 2%) even at low detection levels. Comparative data from ELS and UV detection will also be discussed.

The charged aerosol detector is a sensitive, mass-based detector, especially wellsuited for the determination of any nonvolatile analyte independent of chemical characteristics. As shown in Figure 1, the detector uses nebulization to create aerosol droplets. The mobile phase evaporates in the drying tube, leaving analyte particles, which become charged in the mixing chamber. The charge is then measured by a highly sensitive electrometer, providing reproducible, nanogram-level sensitivity. This technology has greater sensitivity and precision than ELSD and refractive index (RI), and it is simpler to operate than a mass spectrometer (MS).





### **Methods**

#### Black Cohosh

Column: Fused-Core C18 HPLC Column, 4.6 × 150 mm, 2.7  $\mu$ m Column Temp: 35 °C Flow Rate: 1.0 mL/min Mobile Phase: A – 0.1% formic acid in water; B – acetonitrile Gradient: 30% B to 40% B in 12 min; 40% B to 60% B from 12 – 36 min Injection Vol.: 10  $\mu$ L Detection: Charged aerosol detection or evaporative light scattering detection (N<sub>2</sub> pressure 2.3 bar, temperature 50 °C, gain 7) Sample Prep.: About 300 mg of sample extract was weighed into a 50 mL volumetric flask. 40 mL of methanol was added and the flask was sonicated for 15 minutes with occasional shaking. After cooling to room temperature, the flask was filled to volume with methanol and mixed well. A portion of the solution was filtered through a 0.2  $\mu$ m PTFE syringe filter into an HPLC autosampler vial.

#### Ginkgo Biloba

Column: Capcell PAK<sup>®</sup> MG 100 C18, 4.6 × 250 mm; 5 µm Column Temperature: ambient Flow Rate: 1.0 mL/min Mobile Phase A: 5% acetonitrile in 0.1% trifluoroacetic acid Mobile Phase B: 70% acetonitrile in 0.1% trifluoroacetic acid Gradient: Time, % A: 0, 100; 30, 25; 35, 25; 40, 100 Injection Vol.: 10 µL Sample Prep.: Sample (1 g powder from commercially available health supplement) was sonicated in 10 mL methanol for 30 min. The extract was centrifuged (9,000 rpm) for 5 mins and the supernatant was passed through a 0.2 µm nylon filter by centrifugation (9,000 rpm – 2 min).

#### Ginseng

Column: Fused-Core C18 HPLC Column,  $3.0 \times 100$  mm,  $2.7 \mu$ m Column Temp:  $30 \circ$ C Flow Rate: 0.67 mL/min Mobile Phase: A – water; B – acetonitrile Gradient: 15% B to 35% B in 30 minutes Injection Vol.:  $20 \mu$ L

Sample Prep.: About 400 mg of sample extract was weighed and transferred into a 100 mL volumetric flask. 15 mL of methanol was added and the flask was sonicated for 10 min with occasional shaking. About 60 mL of water was added, and the flask was sonicated for an additional 10 min. After cooling to room temperature, the flask is filled to volume with water and mixed well. A portion of the solution is filtered through a 0.2 µm PVDF syringe filter into an HPLC autosampler vial.

#### **Milk Thistle**

Column: Thermo Scientific Acclaim RSLC 120 C18 Column, 2.1 × 100 mm, 2.2  $\mu m$  Column Temp: 40 °C

Flow Rate: 0.50 mL/min

Mobile Phase A: 3 mM ammonium formate, 0.3% formic acid in water:MeOH 80:20 v:v Mobile Phase B: 3 mM ammonium formate, 0.3% formic acid in water:MeOH 20:80 v:v Gradient: 15% B to 45% B in 4.4 min.

Injection Vol.: 1 µL

Sample Prep.: About 70 mg of powdered milk thistle extract was weighed and transferred into a 250 mL amber glass bottle. 100 mL of methanol was added and the flask was sonicated for 20 minutes. A portion of the solution was transferred to a 1.5 mL polypropylene centrifuge tube and centrifuged at 5000 x g for 10 min. The supernatant was transferred into an HPLC autosampler vial.

#### Gotu Kola

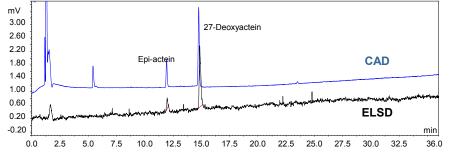
Column: Fused-Core C18 HPLC Column,  $3.0 \times 100$  mm,  $2.7 \mu$ m Column Temp:  $35 \circ$ C Flow Rate: 0.64 mL/min Mobile Phase: A – 0.1% formic acid in Water; B – acetonitrile Gradient: 18% B to 22% B in 8 min; 22% B to 45% B from 8 min to 17 min; 45% B to 80% B from 17 min to 23 min Injection Vol.: 5  $\mu$ L Sample prep.: About 250 mg of sample extract was weighed and transferred into a 100 mL volumetric flask. About 70 mL of methanol was added and the flask was sonicated for 10 min with occasional shaking. After cooling to room temperature, the flask was filled to volume with methanol and mixed well. A portion of the solution is filtered through a 0.2  $\mu$ m PTFE syringe filter into an HPLC autosampler vial.

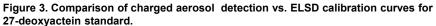
Stevia HPLC Column: Capcell PAK C18 AQ 4.6 × 250 mm, 5 µm Column Temperature: 50 °C Flow Rate: 1.0 mL/min Mobile Phase A: deionized Water (DI), acetonitrile, trifluoroacetic acid (TFA) (95:5:0.1) Mobile Phase B: DI:acetonitrile (5:95) Gradient: 0 – 90% B in 30 min. Hold for 5 min and return to starting conditions. Injection Vol.: 10 µL UHPLC Column: Acclaim<sup>™</sup> RSLC Polar Advantage II 2.1 × 250 mm, 2.2 µm Column Temperature: 40 °C Flow Rate: 0.7 mL / min Mobile Phase A: deionized Water (DI) + 0.1% formic acid Mobile Phase B: acetonitrile + 0.1% formic acid Gradient: 5% to 60% B in 9 min hold 1 min and return to 5% B Injection Vol.: 5 µL Sample Prep.: Commercially available stevia extract powder was dissolved in deionized water (0.9 mg/mL)

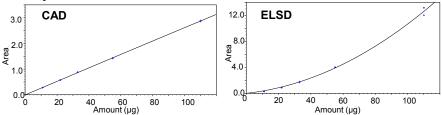
### **Results and Discussion**

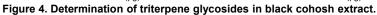
Extracts of black cohosh have been used since the 1950s to relieve symptoms of menopause. The active ingredients are believed to be related to the content of triterpene glycosides present in black cohosh, including 27-deoxyactein, actein, cimiracemoside F, and others. Many of the triterpene glycosides do not possess chromophores above 200 nm. HPLC-ELSD is the most common method for quantitation of triterpene glycosides in black cohosh products. However, ELSD suffers from poor sensitivity (Figure 2), and has a very non-linear response (Figure 3). Charged aerosol detection overcomes these issues, offers better reproducibility and can easily measure low-abundant analytes in a Black Cohosh extract (Figure 4).

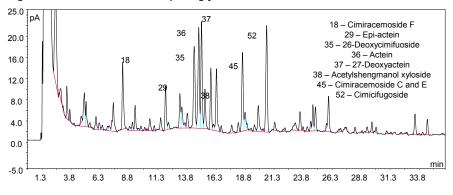
Figure 2. Comparison of charged aerosol detection vs. ELSD for determination of triterpene glycosides in black cohosh (27-deoxyactein standard solution:  $11 \mu g/mL$ ).





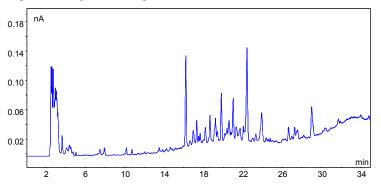






*Ginkgo biloba* is thought to possess nootropic activity, and is taken to improve memory and enhance concentration. The active ingredients are believed to be the sesquiterpenoid bilobalide and numerous diterpenoid ginkgolides. Although not labeled in Figure 5, these compounds elute between 3 and 14 minutes.





Asian Ginseng (*Panax ginseng*) has traditionally been used as a tonic to reduce the effects of stress, counteract fatigue, and increase stamina. The main bioactive ingredients found in *Panax ginseng*, and a related species *Panax quinquefolius* (American ginseng) are triterpene saponins, commonly referred to as ginsenosides. There are 7 major ginsenosides present in *Panax ginseng*: the protopanaxatriols (Rg1, Re and Rf), and protopanaxadiols (Rb1, Rc, Rb2 and Rd). HPLC with low-wavelength UV detection (203 nm) is most commonly used but suffers from poor sensitivity. HPLC with charged aerosol detection not only improves the baseline slope seen with gradient elution, but also offers improved sensitivity (Figure 6).

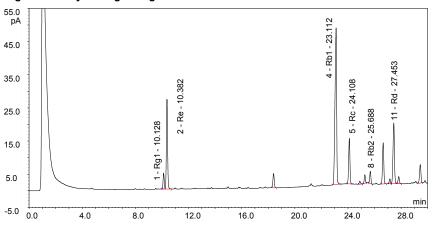


Figure 6. Analysis of ginseng extract.

Evidence suggests that milk thistle extracts may both prevent and repair damage to the liver from toxic chemicals and medications and may be of use in the treatment of mushroom poisoning. The active ingredients are thought to be numerous flavonolignans including silybinin, isosilybinin, silycristin and silydianin (Figure 7).

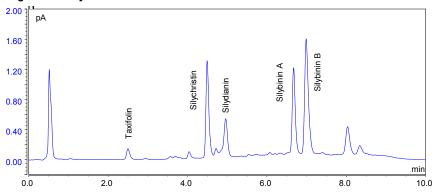
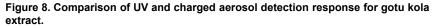
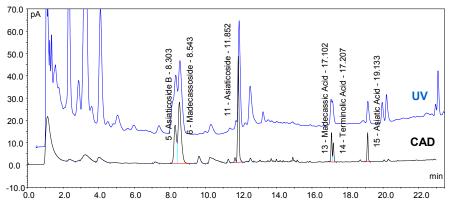


Figure 7. Analysis of milk thistle extract.

*Centella asiatica* (commonly called gotu kola) is a small herbaceous annual plant that is native to India, Sri Lanka, northern Australia, and other parts of Asia and the western Pacific. It is used as a medicinal herb in Ayurvedic medicine as well as traditional Chinese medicine for a wide variety of conditions, such as improving memory, blood flow, as a wound-healing agent, and as a topical application for skin conditions such as ulcers, wounds, and eczema. The chemical compounds of interest in gotu kola are usually considered to be the ursane- and oleanane-type triterpenes and triterpene glycosides. Although low-wavelength UV can be used to measure these compounds, it suffers from sensitivity and baseline issues – these can be readily overcome by the CAD (Figure 8). Interestingly, this figure also shows that the number of potential interferences are reduced when using the CAD, probably as the sample has a number of volatile compounds that absorb low-wavelength UV.





Stevia rebaudiana, commonly known as stevia, is currently being used as a low calorie natural sweetener. The active ingredients include the aglycone diterpene known as steviol and numerous glycosides including stevioside and the rebaudiosides. Although low-wavelength UV was initially used, it suffered from the numerous issues described above. There are a number of HPLC with charged aerosol detection methods in the literature. Here we show the improved chromatographic resolution and significantly reduced analysis time using an UHPLC method. (Figure 9).

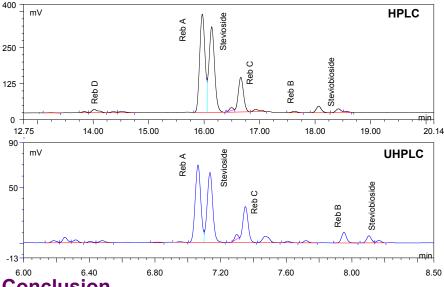


Figure 9. Comparison of HPLC and RSLC for the analysis of a stevia extract.

## Conclusion

Charged aerosol detection offers numerous analytical improvements over ELSD, RI and low-wavelength UV and is of particular use to the analysis of natural products and botanicals. The CAD:

- Measures any non-volatile and many semi-volatile species.
- Response is little affected by chemical structure.
- Allows estimation of analyte amounts even when external standards are not available.
- Has excellent sensitivity, reproducibility and a dynamic range of over 4 orders of magnitude,

#### www.thermofisher.com/dionex

Capcell PAK is a registered trademark of Shiseido. All other 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.

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



Part of Thermo Fisher Scientific

PN70021\_E 07/16S