

Fast Sugar Analysis for Food and Beverage Using HPAE-PAD on a Compact IC System

Hua Yang, Beibei Huang, Terri Christison, and Jeff Rohrer, Thermo Fisher Scientific, Sunnyvale, CA

OVERVIEW

Purpose: To demonstrate applications using HPAE-PAD with a High Concentration Carbohydrate Analysis kit to determine mg/L and g/L sugars in beverages, alcohol and vinegar samples.

Methods: Monosaccharide and disaccharide sugars are ionized in strong base, separated by High Performance Anion-Exchange (HPAE) chromatography, and detected by Pulsed Amperometric Detection (PAD). Analysis is facilitated by the Thermo Scientific™ Dionex™ Integri™on™ HPIC™ system.

Results: mg/L and g/L concentrations of sugars were directly determined in various diluted beverage, alcohol, and vinegar samples.³

INTRODUCTION

As a result of the Nutrition Labeling and Education Act of 1990 (NLEA), the U.S. Food and Drug Administration (FDA) requires all packaged foods and drinks to list the sugar content per serving according to Code of Federal Regulations (CFR) Title 21.¹ This information must appear on a product's nutrition facts label, along with things like calorie, fat, sodium, and fiber content. Sugars are defined as the sum of all free mono- and disaccharides (such as glucose, fructose, lactose, and sucrose). Excess sugar consumption is associated with obesity and poor health. Although labeling is not required for alcoholic beverages, consumers may need to know sugar concentrations.

As consumers' interest in health-conscious diets has grown, so has the global beverage industry – with the introduction of vitamin fortified water, energy drinks, anti-aging water, and herbal nutritional supplements. As a result, more analytical methods are needed.

High-performance anion-exchange chromatography (HPAE) coupled with pulsed amperometric detection (PAD) is a well-established technique to accurately identify and quantify carbohydrates in food and beverage samples.² By accurately determining the sugar concentrations, HPAE-PAD is used to identify contamination and adulteration, maintain product consistency, and to ensure regulatory compliance of raw ingredients (water, additives, and fruit) and the final product.

Here, we show the determination of sugars in drinks using the Thermo Scientific™ Dionex™ Integri™on™ HPIC™ system. This system allows fast determination of sugars without manual eluent preparation or sample derivatization.³

METHODS

See chromatograms

Instrument:

Dionex Integri™on™ HPIC system (Figure 1) configured for electrochemical detection, includes: eluent generation, Thermo Scientific™ Dionex™ IC PEEK Viper™ fittings (Figure 1, bottom left). The flow diagram is shown in Figure 2.

Figure 1. Dionex Integri™on™ HPIC system configured for electrochemical detection.

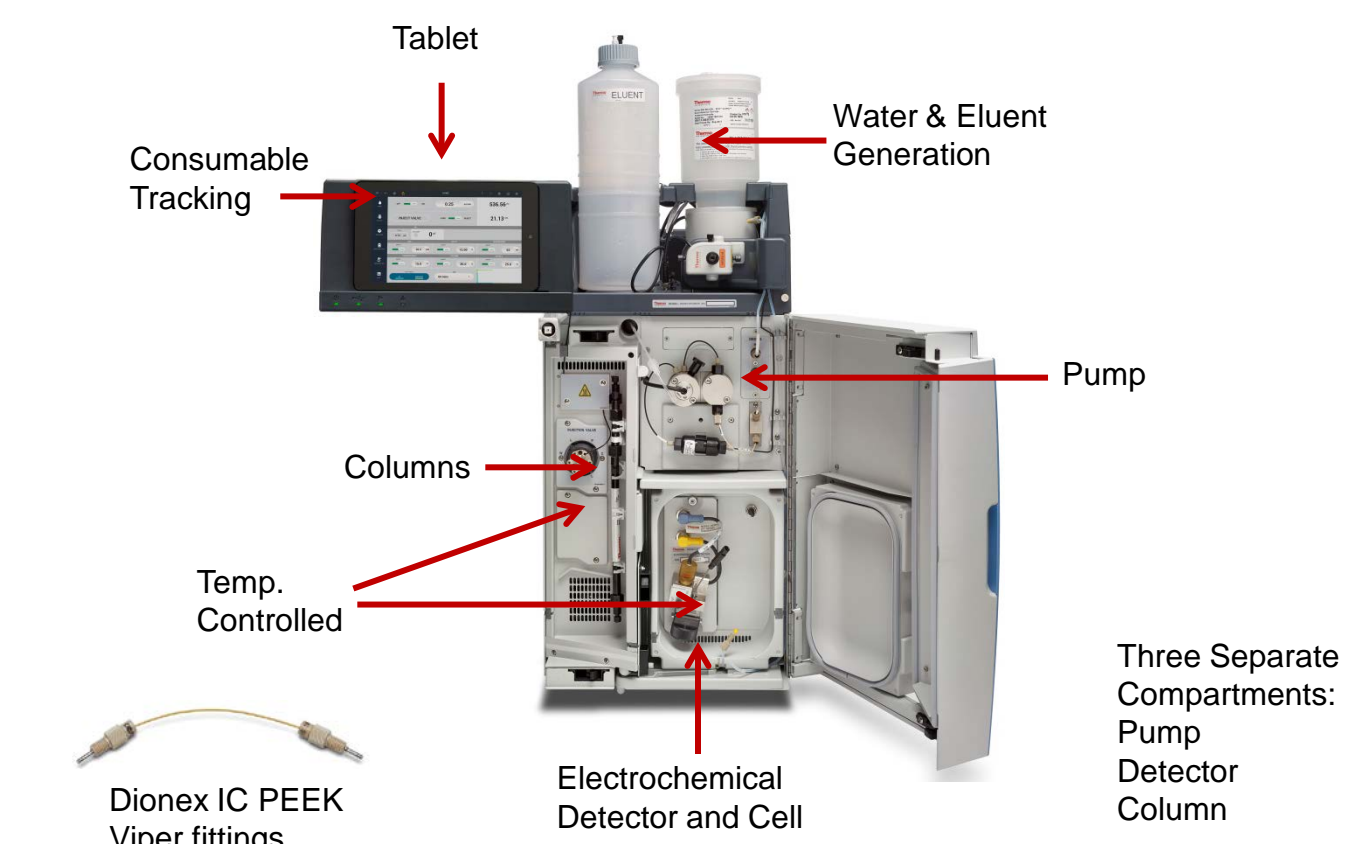
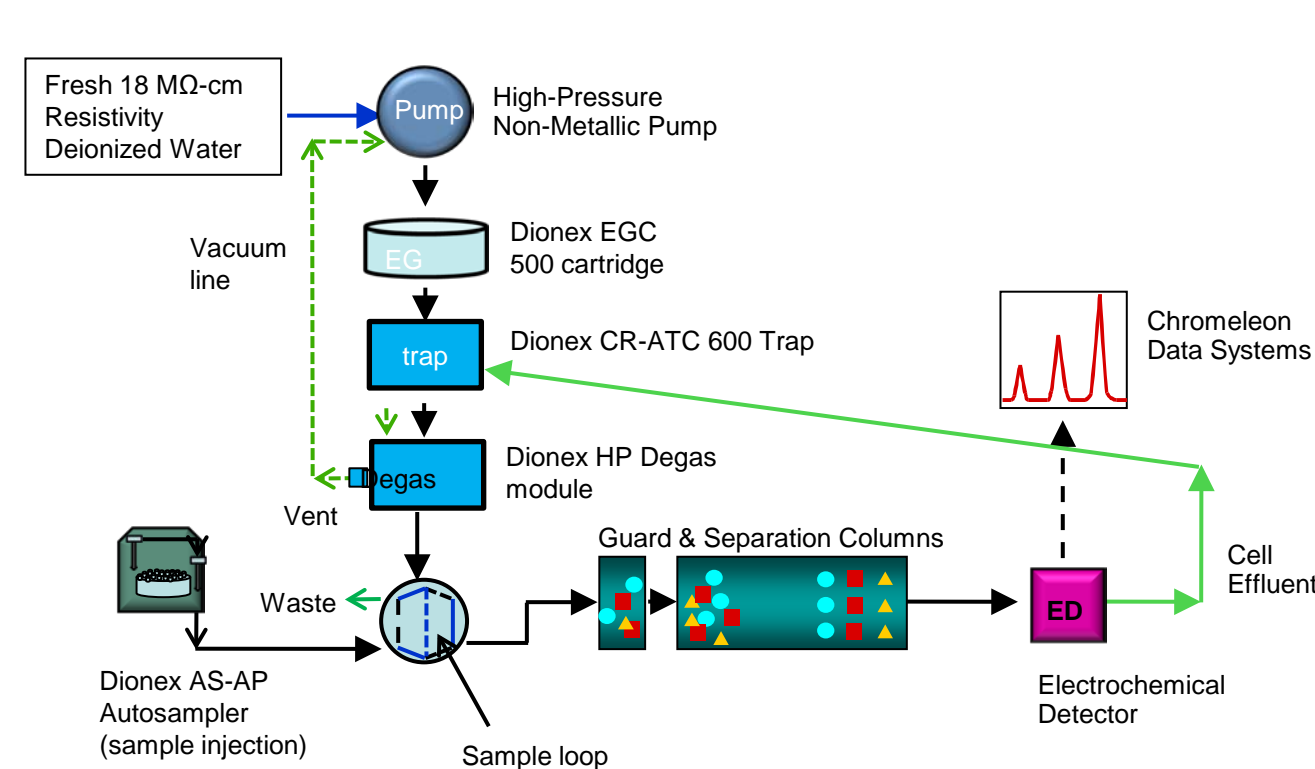
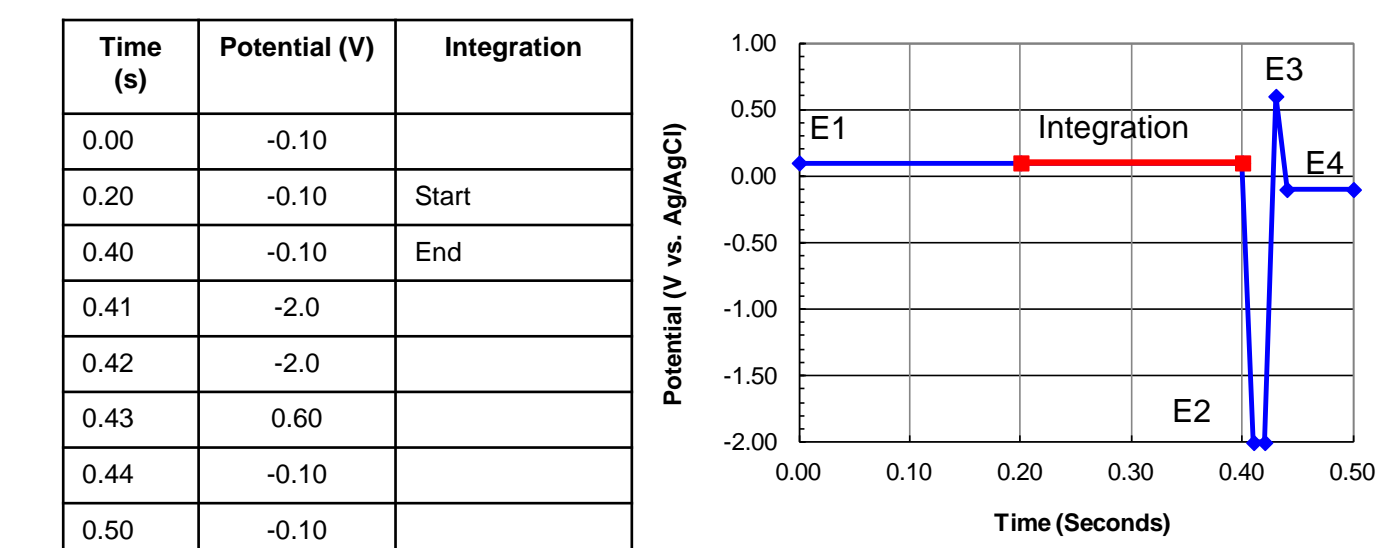


Figure 2. Flow diagram for the Dionex Integri™on™ HPIC system configured for electrochemical detection.



In PAD, using the four-potential waveform, the disposable working electrode is pulsed through the different potentials at set times, completing two cycles within one second (Figure 3). This waveform is optimized to provide a clean, stable gold-oxide layer in preparation for detection of the next eluting peak.

Figure 3. Four-potential carbohydrate waveform.



1. Detection potential (E1)
2. Duration time of E1 (t1)
3. Integration period (Integrate)
4. Reductive cleaning potential (E2)
5. Duration time of E2 (t2)
6. Oxidative cleaning potential (E3)
7. Duration time of E3 (t3)
8. Pre-detection (oxide reduction) potential (E4)
9. Duration time of E4 (t4)

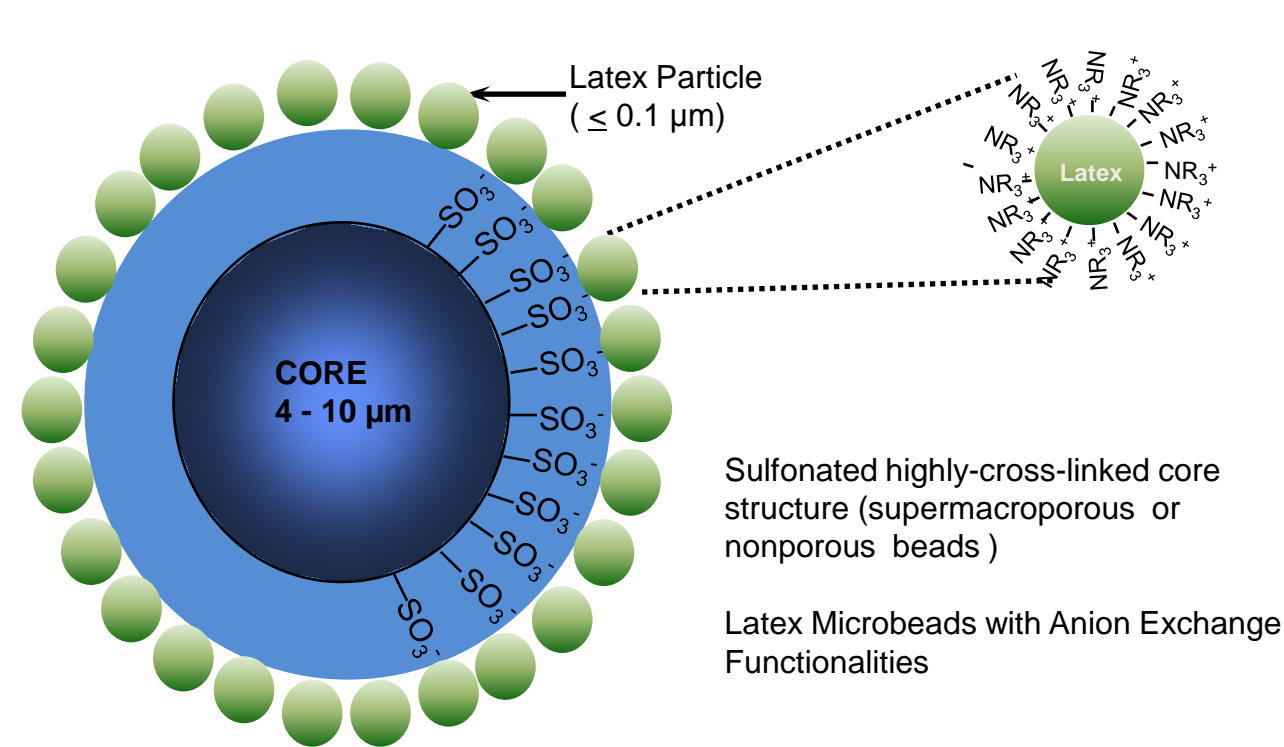
DATA ANALYSIS

Thermo Scientific™ Dionex™ Chromeleon™ Chromatography Data System (CDS) software, version 7.2 SR4.

RESULTS

The determinations of sugars in drinks using the Dionex Integri™on™ HPIC system are shown here. Carbohydrates, including sugars, are weak acids. At high pH, they are ionized, and thus can be separated by anion-exchange chromatography. Figure 4 illustrates an anion exchange bead in a Thermo Scientific™ Dionex™ CarboPac™ column. The core polymer is polystyrene divinylbenzene which has been surface sulfonated, followed by latex agglomeration. This pellicular resin structure permits excellent mass transfer, resulting in high-resolution IC chromatography. With PAD, a simple electrochemical detection, HPAE-PAD directly quantifies carbohydrates with high sensitivity and selectivity and without derivatization. With electrochemical detection, Dionex IC PEEK Viper™ fittings and Reagent-Free™ IC (RFIC), the newly introduced Dionex Integri™on™ HPIC system offers a fast and cost-effective HPAE-PAD method for sugar analysis.

Figure 4. Polymeric anion exchange particles in Dionex CarboPac columns.



Determinations of monosaccharide and disaccharide sugars were conducted in various beverages, vinegars, and flavored distilled alcohol samples. In Figures 5 and 6, sugars were determined in the same functional beverages using the Carbohydrate gold disposable working electrode. The three sugars in the 10,000-fold diluted samples were separated on the Thermo Scientific™ Dionex™ CarboPac™ PA20 column, and detected using the standard 0.003-in thick gasket (Figure 5). In contrast, and for confirmation, the 100-fold diluted samples were analyzed on the Thermo Scientific™ Dionex™ CarboPac™ SA10-4μm column using the 0.062-in thick gasket from the High Concentration Carbohydrate Analysis kit. This column was optimized for fast separations with a different selectivity as evident by the different elution order. The difference in concentration is attributed to dilution error.

Figure 5. Sugars in 10,000 × diluted functional beverage samples

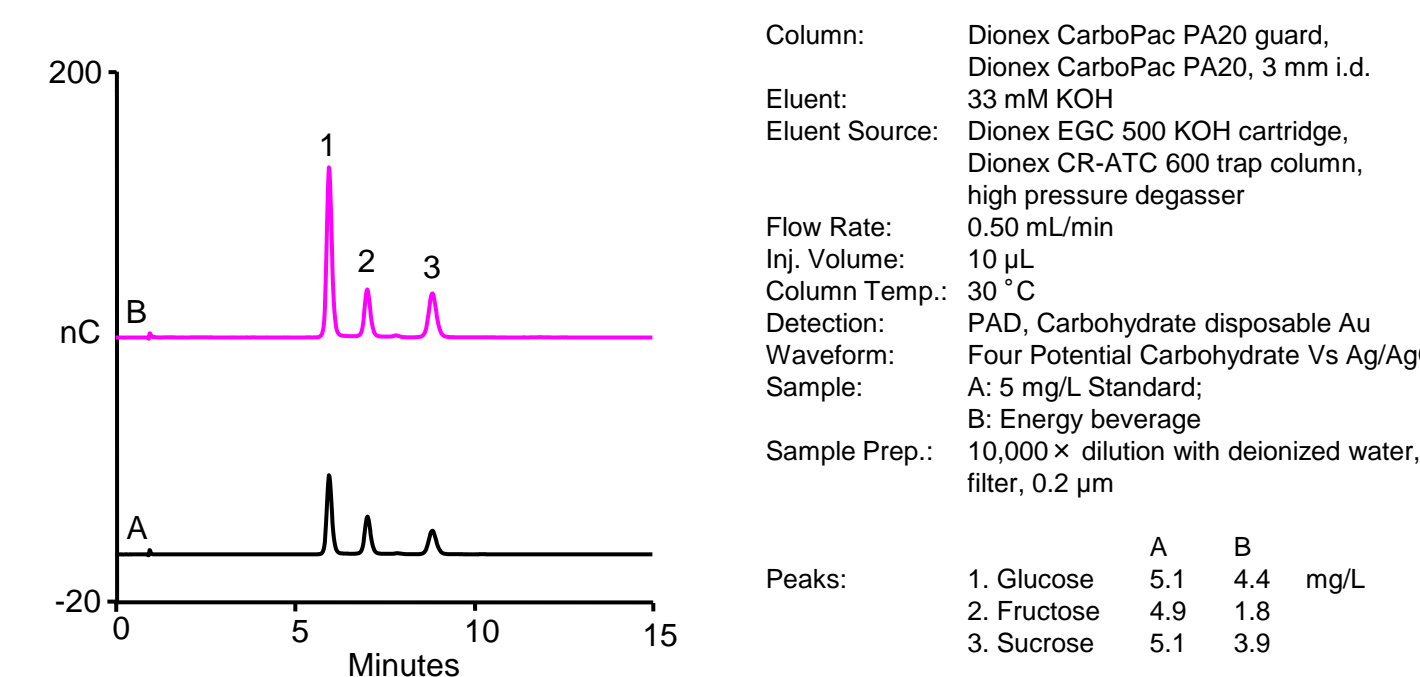
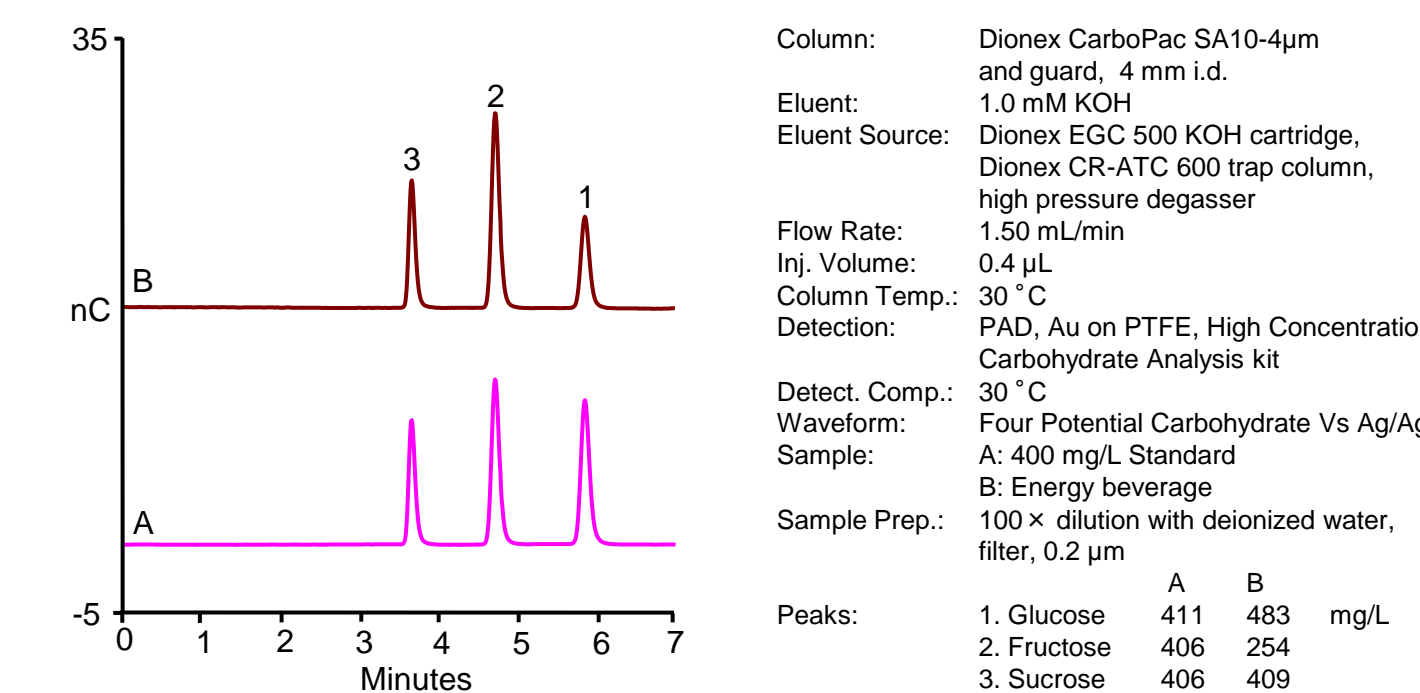


Figure 6. Sugars in 100 × diluted functional beverage samples using the Dionex CarboPac SA10-4μm column and High Concentration Carbohydrate Analysis kit.



In Figures 7 and 8, the same analysis is repeated with fruit juices and carbonated beverages on the Dionex CarboPac SA10-4μm column optimized for fast separations.

Figure 7. Sugars in 100 × diluted fruit juice samples using the Dionex CarboPac SA10-4μm column and High Concentration Carbohydrate Analysis kit.

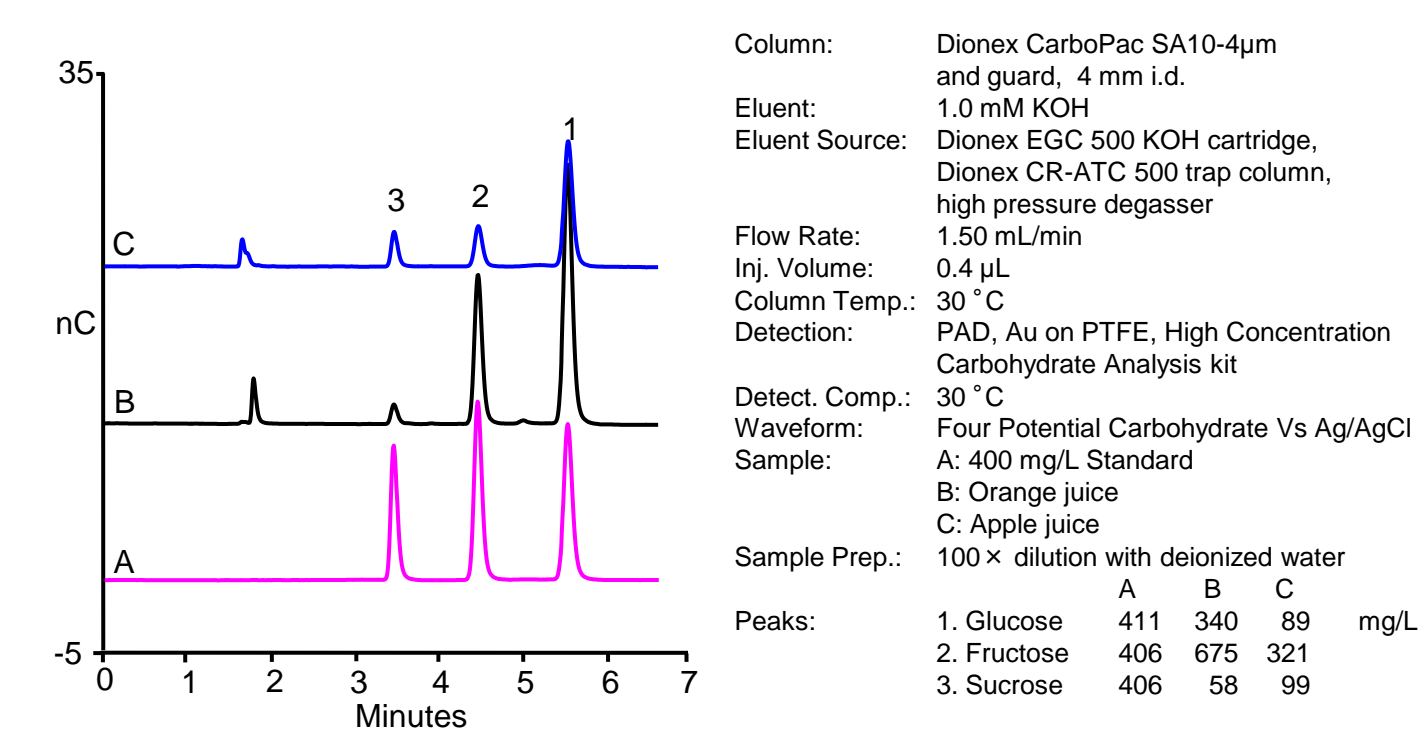
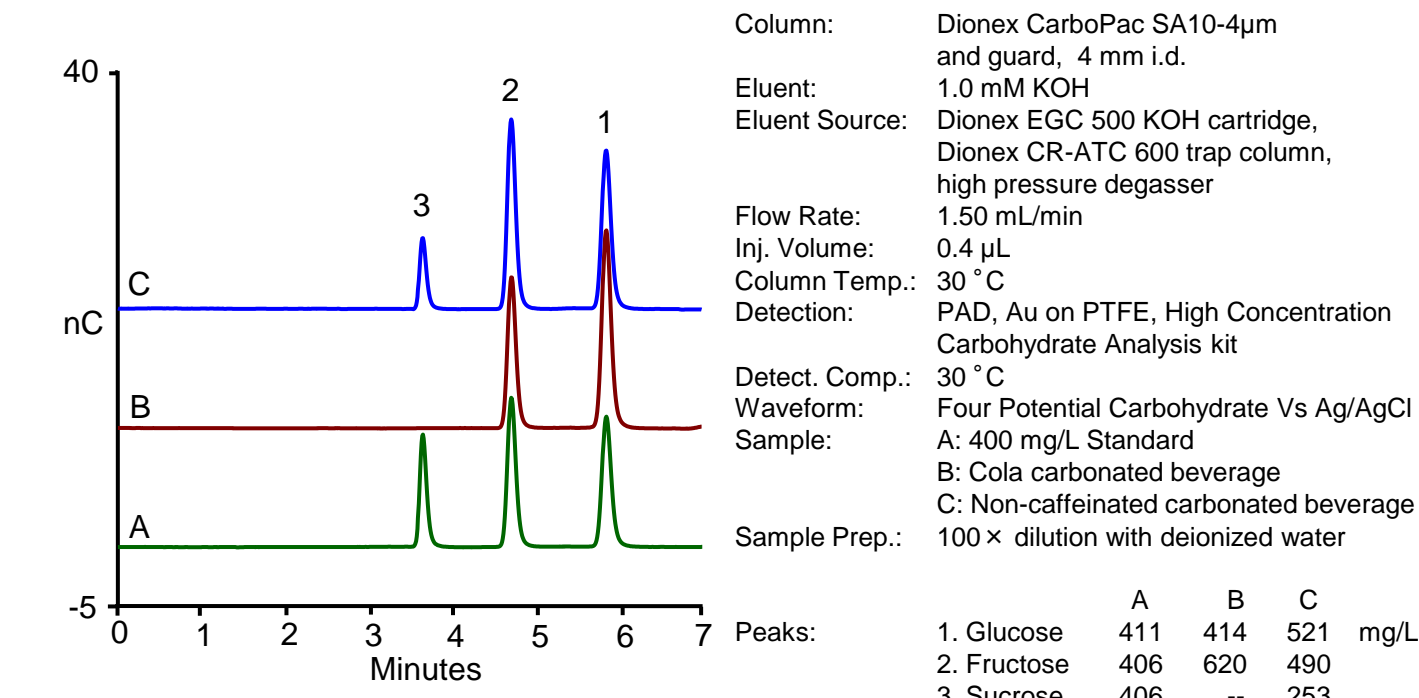


Figure 8. Sugars in 100 × diluted carbonated beverage samples using a Dionex CarboPac SA10-4μm Column and High Concentration Carbohydrate Analysis kit.



In Figures 9–11, sugars were determined in fermented and flavored distilled alcohol samples. In these examples, the Au on PTFE disposable working electrode was combined with the High Concentration Carbohydrate Analysis kit. These samples were diluted only 100-fold. For the alcoholic samples, the Dionex CarboPac PA20 column was selected because ethanol and sucrose nearly co-elute on the CarboPac SA10-4μm column. The results show good chromatography as shown by the narrow peak widths and strong response. The balsamic vinegar sample can be analyzed on either column.

Figure 9. Sugars in a balsamic vinegar sample.

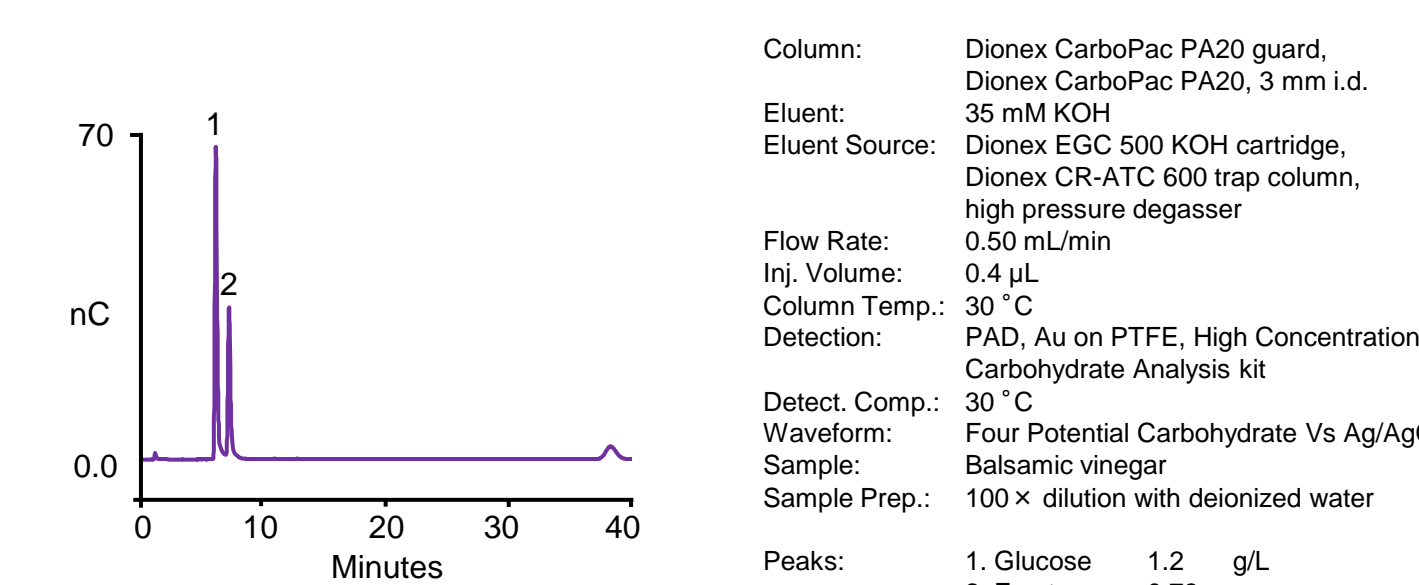


Figure 10. Sugars in a flavored rum sample.

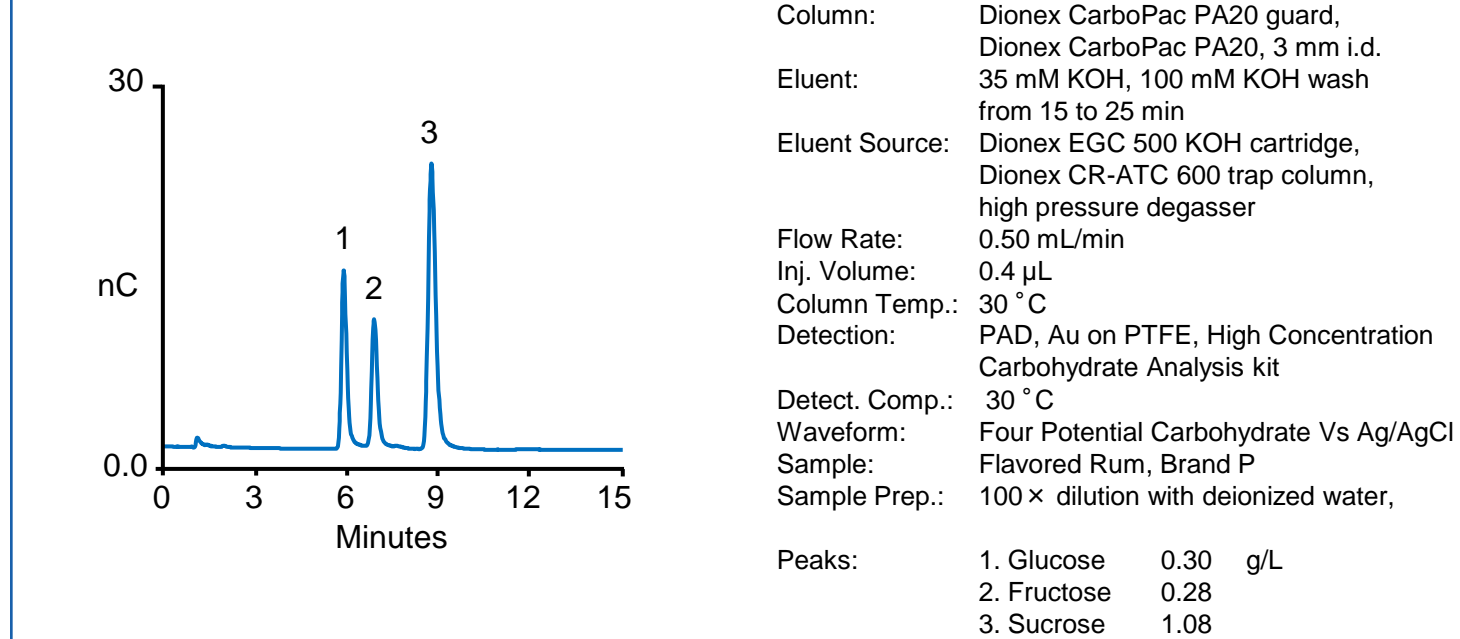
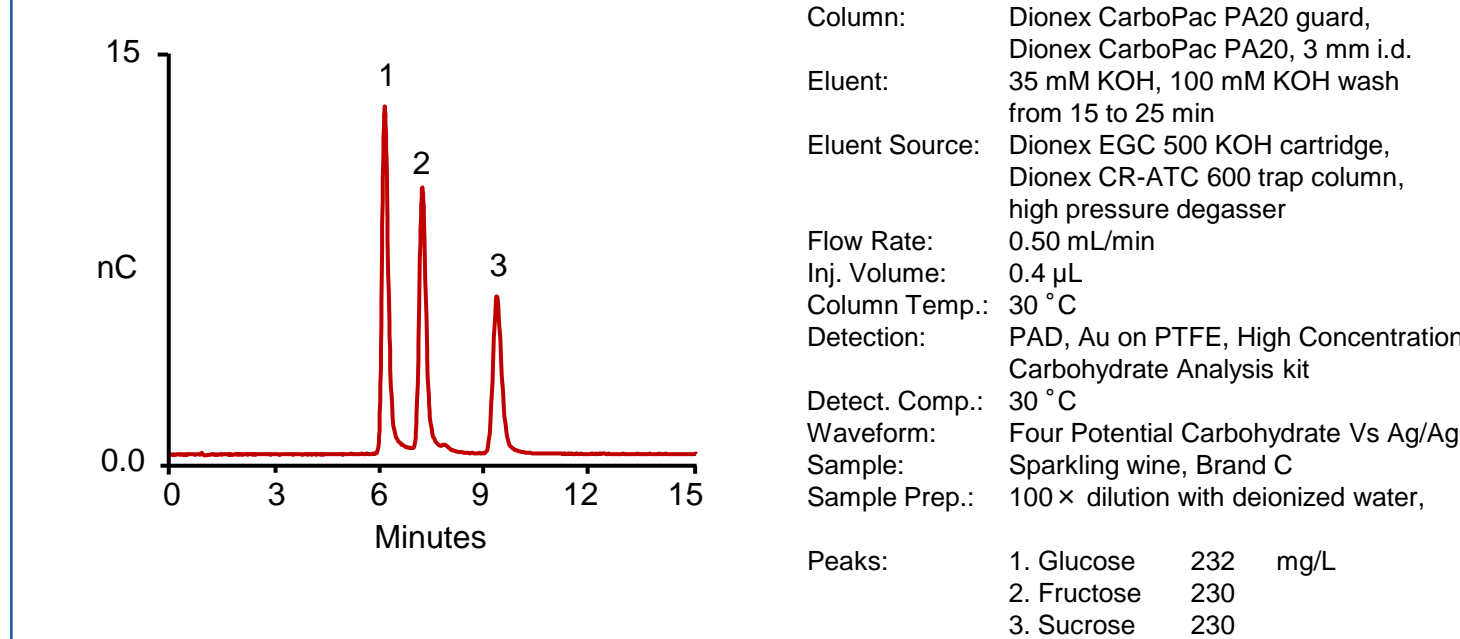


Figure 11. Sugars in a sparkling wine sample.



CONCLUSIONS

- HPAE-PAD is a sensitive method to directly determine sugars in diverse samples.
- HPAE-PAD combined with the High Concentration Carbohydrate Analysis kit extends the analytical range to mg/L and g/L concentrations with minimal dilution of samples, thus reducing dilution errors and improving reporting accuracy.
- This technique was demonstrated by the determination of mg/L to g/L sugars in sports drinks, colas, and alcoholic beverages.

REFERENCES

- 21 CFR 101.9 - NUTRITION LABELING OF FOOD
- Thermo Scientific Technical Note 20.
- Thermo Scientific™ AppsLab Library of Analytical Applications. (search word: sugars, Integri™on™)

TRADEMARKS/LICENSING

© 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.