

Efficient Separations Obtained by Using Smaller Particle Size Analytical Columns with a High-Pressure Ion Chromatography System

*Barbara Shao, Frank Hoefler, Maria Rey, and Linda Lopez,
Thermo Fisher Scientific, Sunnyvale, CA, USA*



Overview

Purpose: This poster demonstrates smaller-particle-size (4 μm) analytical columns combined with high-pressure ion chromatography to achieve efficient separations, fast analyses, and high throughput.

Methods: Inorganic anions in fruit juice, wine, or environmental and mineral water samples are separated using Thermo Scientific Dionex IonPac AS11-HC-4 μm , AS18-4 μm , or CS19-4 μm anion-exchange smaller particle size resins in prototype standard and microbore column formats, combined with a modified Thermo Scientific Dionex ICS-5000 HPIC system, with prototype high-pressure pumps, eluent generator cartridges, and degas modules, allowing operation in all column formats at 5000 psi system backpressures. Current eluent suppressors are compatible with high-pressure systems, and are operated using AutoSuppression™, either in recycle mode or external water mode, with detection by suppressed conductivity. The advantages of better separation of ions in complex samples and faster analysis are presented.

Results: Inorganic anions in municipal and mineral water samples are separated within 4 min with good baseline resolution using a prototype 4 μm particle size Dionex IonPac™ AS18-4 μm microbore (2 mm i.d.) format analytical column; multiple inorganic anions and organic acids are resolved in single runs in fruit juice and white wine samples, using a prototype standard (4 mm i.d.) format Dionex IonPac AS11-HC-4 μm column. Several of the separated components showed enhanced baseline resolution using the 4 μm particle-size columns relative to their 9 μm counterparts with the same stationary phase. Inorganic cations are separated within 8 min with high peak efficiency and high resolution using a prototype Dionex IonPac CS19-4 μm 4 mm i.d. column.

Introduction

Typical ion chromatography systems and analytical columns in standard 4 mm and microbore 2 mm formats operate below a maximum pressure of 3000 psi using well-established Reagent-Free Ion Chromatography (RFIC) system techniques. A smaller-particle-size analytical column, however, requires high-pressure system capability. The Dionex ICS-5000 Capillary HPIC™ system can operate continuously at up to 5000 psi using eluent generation. The modified system used here extends the high system backpressure capabilities to the entire flow rate range of the Dionex ICS-5000 system.

Running any columns at increased flow rates may cause a decrease in peak efficiencies. However, the superior chromatographic fidelity achieved using the new 4 μm columns, including the Dionex IonPac AS11-HC-4 μm , AS18-4 μm , and IonPac CS19-4 μm , minimizes such losses, to provide better separation of ions in complex matrices and can provide faster analysis times and higher throughput. Higher peak efficiencies can be achieved, even when running at faster flow rates, providing shorter analysis times and increased productivity. These columns are currently available in 0.4 mm i.d. formats, but 2 and 4 mm i.d. columns are in development, and prototypes are demonstrated here.

Methods

Sample Preparation

The samples were diluted before analysis without other pretreatment

Ion Chromatography

Equipment and Data Analysis

Dionex ICS-5000 system, modified for high-pressure operation
Dionex IonPac AS11-HC-4 μm , prototype 4 × 250 mm analytical column
Dionex IonPac AS18-4 μm , prototype 2 × 150 mm analytical column
Dionex IonPac CS19-4 μm , prototype 4 × 250 mm analytical column
Thermo Scientific Dionex Chromeleon Chromatography Data System software

Conditions

The conditions are described in the figures.

Results

Prototype 4 mm i.d. Dionex IonPac AS11-HC-4 μm Analytical Column

Small-particle-size columns such as the Dionex IonPac AS11-HC-4 μm can improve baseline resolution and separation of organic acids and inorganic anions (see Figures 1 and 2 and Table 1); the high capacity and high efficiency of this analytical column permit separation of a large number of organic acids and inorganic anions in complex sample matrices such as wine or fruit juice samples (see Figures 3 and 4). This ion-exchange column uses smaller resin particles for more efficient separations resulting in more accurate peak integration and more reliable results; therefore, a large number of inorganic anions and organic acids in fruit juice and white wine samples can be resolved in a single run using a hydroxide gradient.

FIGURE 1. Determination of organic acids and inorganic anions in standard mixture using a prototype Dionex IonPac AS11-HC-4 μ m, 4 \times 250 mm column.

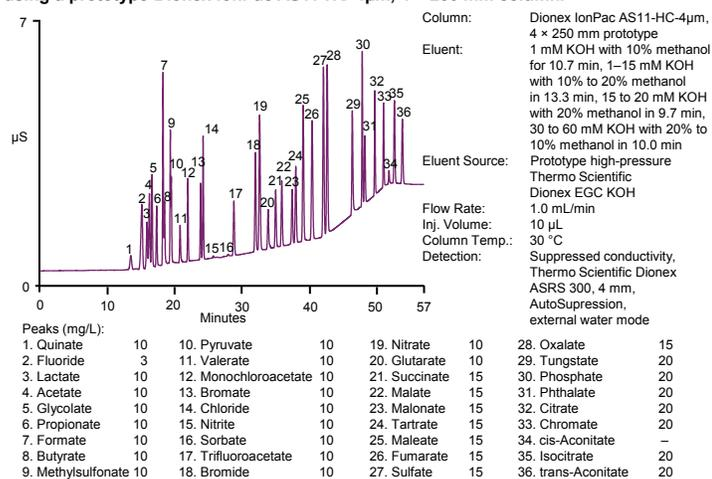


FIGURE 2. The Dionex IonPac AS11-HC smaller-particle-size (4 μ m) analytical column provides improved resolution over the AS11-HC bigger-particle-size (9 μ m) analytical column, including improved separation for monovalent carboxylic acids, such as lactate and acetate, and better baseline resolution for bromate/chloride, formate/butyrate, methylsulfonate/pyruvate, and phosphate/phthalate (see Table 1).

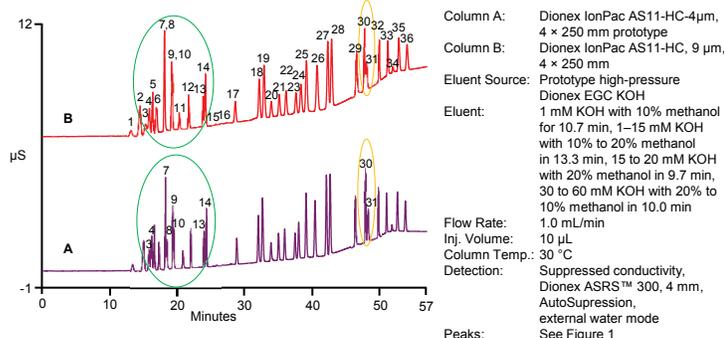


TABLE 1. Comparison of resolution obtained on both columns with different particle size

	Lactate (P3)	Bromate (P7)	Formate (P9)	Methylsulfonate (P13)	Phosphate (P30)
Resolution, 4 μ m column	1.25	1.10	0.99	1.68	1.49
Resolution, 9 μ m column	1.16	None	None	1.26	1.08

FIGURE 3. Determination of organic acids and inorganic anions in orange juice by using a prototype Dionex IonPac AS11-HC-4 μ m column with high-pressure ion chromatography (IC conditions refer to Figure 1).

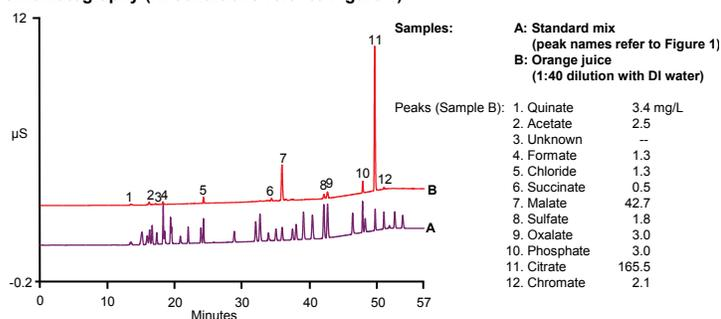
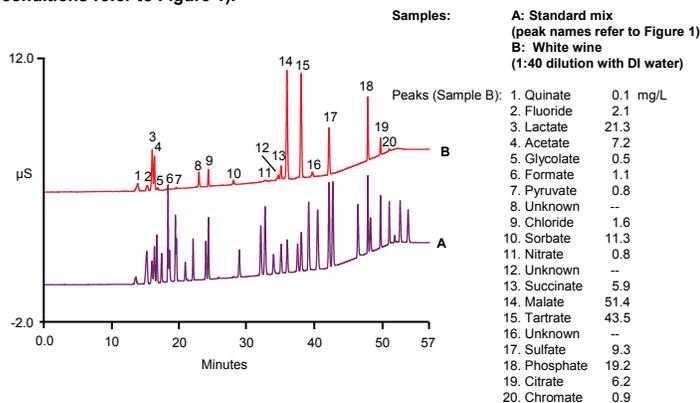


FIGURE 4. Determination of organic acids and inorganic anions in white wine using a Dionex IonPac AS11-HC-4 μ m, 4 \times 250 mm prototype anion-exchange column (for IC conditions refer to Figure 1).



Prototype 2 mm i.d. Dionex IonPac AS18-4 μ m Analytical Column

In 1993, the U.S.EPA Method 300.0 (A) defined the use of the Dionex IonPac AS4A anion-exchange column to determine inorganic anions in environmental waters using a manually prepared carbonate-based eluent and suppressed conductivity detection. Later, the development of the Dionex IonPac AS18 hydroxide-selective columns permitted use of RFIC system methods in accordance with U.S. EPA Method 300.0 (A). Reagent-Free IC systems can electrolytically generate eluent inline by just adding water, which improves selectivity and reproducibility, and eliminates errors and labor to prepare hydroxide eluent manually.

Now Reagent-Free IC methods can be employed on the Dionex IonPac AS18 chemistry using 4 μ m particle-sized columns for maximum performance. Inorganic anions that commonly exist in environmental water or drinking water samples can be analyzed using a Dionex IonPac AS18-4 μ m analytical column with an HPIC system using well-established RFIC system techniques. The modified high-pressure system combined with prototype smaller-particle-size columns permit fast run times without compromising resolution also with 2 and 4 mm columns (Figure 5).

FIGURE 5. Without sacrificing the baseline resolution and column efficiency, a shortened run time was obtained by using a prototype smaller-particle-size column (4 μ m) compared to a larger-particle-size (7.5 μ m) column for the separation of nine anions.

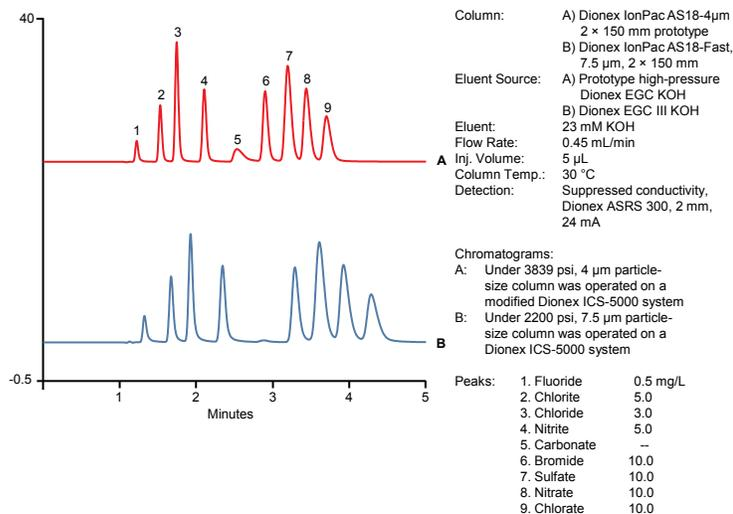


FIGURE 6. Determination of nine inorganic anions using a prototype Dionex IonPac AS18-4 μ m, 2 \times 150 mm anion-exchange column. At a high flow rate of 0.50 mL/min, a high backpressure of 4233 psi resulted with a run time of only 4 min with good baseline resolution.

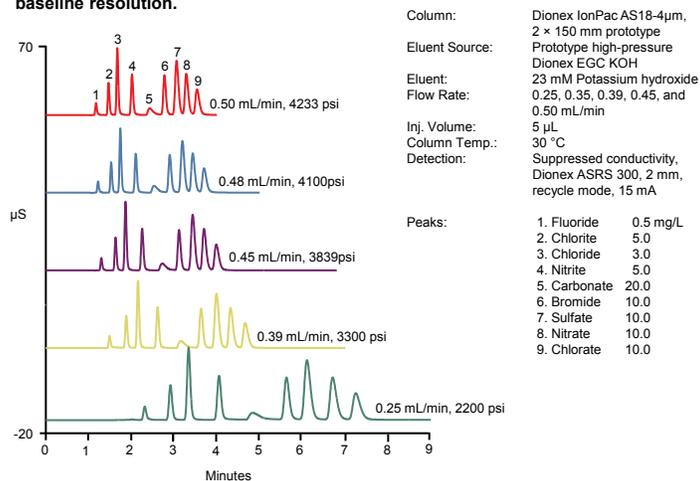
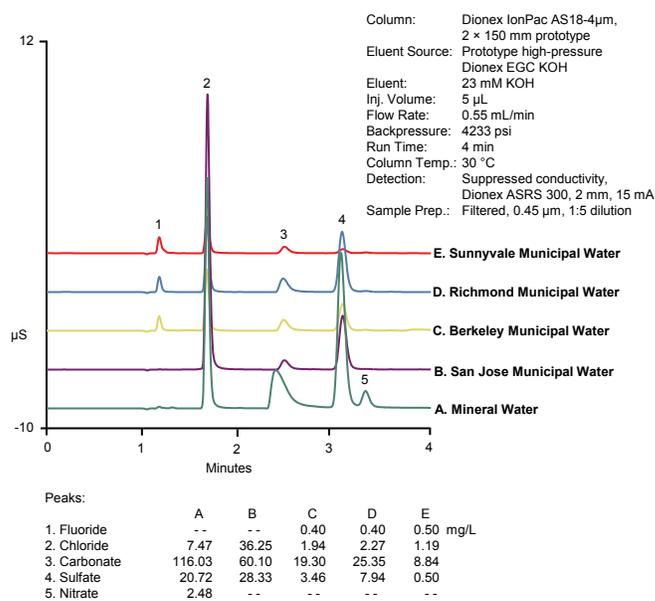
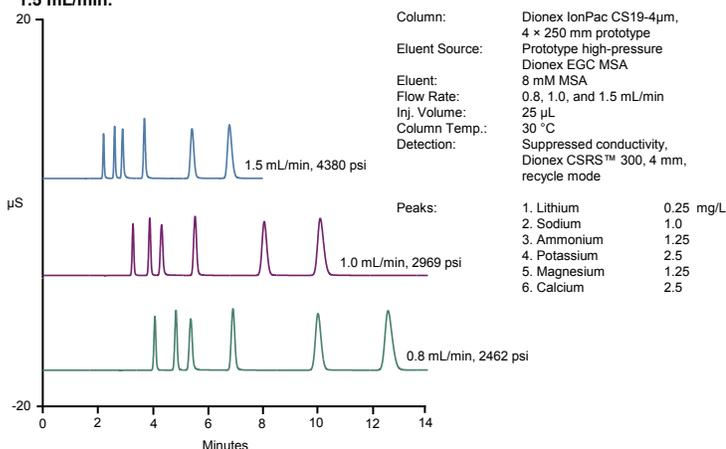


FIGURE 7. Determination of inorganic anions in water samples using a prototype Dionex IonPac AS18-4 μ m 2 \times 150 mm column. Five anions in municipal or mineral water samples were eluted within 4 min on the 4 μ m particle size column, emphasizing the fast sample runs possible with this column.



Smaller Particle Size Analytical Column for Cation Analysis

FIGURE 8. This application demonstrates resolution of the six cations within 7 min using a prototype Dionex IonPac CS19-4 μ m 4 \times 250 mm 4 μ m particle-size analytical column at high flow rates. Fast sample runs with high peak efficiency and resolution are possible with the 4 μ m particle size column; additionally, it demonstrates the requirement of high pressure (4380 psi) capability of HPIC at a flow rate of 1.5 mL/min.



Conclusion

The smaller resin particles produce a backpressure over 3000 psi at increased flow rates with the Dionex IonPac AS18-4 μ m analytical column prototype, and at standard flow rates with the Dionex IonPac AS11-HC-4 μ m and CS19-4 μ m analytical column prototypes. This requires a new, high-pressure analytical system. The modified Dionex ICS-5000 system provides the platform to operate these smaller-particle-size columns under high back pressure, yielding highly efficient separations for high resolution or fast analyses and higher throughput.

- The Dionex IonPac AS11-HC-4 μ m analytical column provides improved separation for monovalent carboxylic acids and improved resolution between formate and butyrate, methylsulfonate and pyruvate, bromate and chloride, and phosphate and phthalate, compared to the standard 9 μ m particle-size column.
- The Dionex IonPac AS18-4 μ m analytical column provides efficient anion separations at higher flow rates thereby increasing throughput for environmental and mineral water samples.
- Fast sample runs with high peak efficiency and high resolution of cations and amines are possible with the Dionex IonPac CS19-4 μ m column.

References

For more information regarding Thermo Scientific Dionex IonPac AS18, IonPac AS11-HC and IonPac CS19 analytical columns, refer to follow Thermo Fisher Scientific application notes and website.

1. Thermo Fisher Scientific. *Application Note 154, Determination of Inorganic Anions in Environmental Water Using a Hydroxide-Selective Column.*
2. Thermo Fisher Scientific. *Application Note 143, Determination of Organic Acids in Fruit Juices.*
3. <http://www.dionex.com/en-us/products/columns/ic-rfic/cation-packed/ionpac-cs19/lp-110108.html>

Acknowledgements

We would like to thank Charanjit Saini, Andy Woodruff, Cathy Tanner, Kris Bahten, Terri Christison, and Fei Pang for technical support.

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.

U.S./Canada	(847) 295 7500	Denmark	(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 3420 8600
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

Thermo
SCIENTIFIC
Part of Thermo Fisher Scientific

PN70147_E 08/16S