

# A New Hydroxide Selective Anion Exchange Phase for Ion Chromatography

Charanjit Saini, Chris Pohl, Yan Liu  
Thermo Fisher Scientific, 1228 Titan Way, Sunnyvale, CA 94085

## Overview

**Purpose:** Development and evaluation of a new anion exchange phase for Ion Chromatography

**Methods:** Ion Chromatography

**Results:** Dionex IonPac® AS28-4 $\mu$ m offers superior performance in terms of selectivity, peak efficiency and sensitivity.

## Introduction

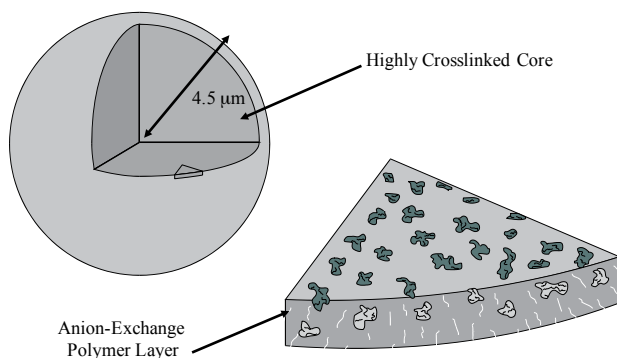
In the area of ion chromatography of anions, hydroxide based eluent systems provide the highest detection sensitivity as well as the lowest noise, resulting in substantially lower detection limits. Hydroxide eluent is ideally suited to the application of gradients in anion exchange ion chromatography in terms of detection properties. As a result, we have developed a series of new hydroxide selective phases over the past 15 years aimed at producing materials well suited to the chemical and chromatographic properties of hydroxide eluents. In this presentation, we will demonstrate recent advances in hydroxide selective stationary phases.

## Experimental Conditions

- All separations were performed using an ICS-5000 system (Dionex, Sunnyvale, CA) equipped with dual pump, eluent generator, conductivity detectors and AERS in recycle mode
- Columns used were Dionex IonPac® AS15 and AS28-4 $\mu$ m
- Dionex Chromeleon 6.8 chromatography management software was used for controlling system and data processing

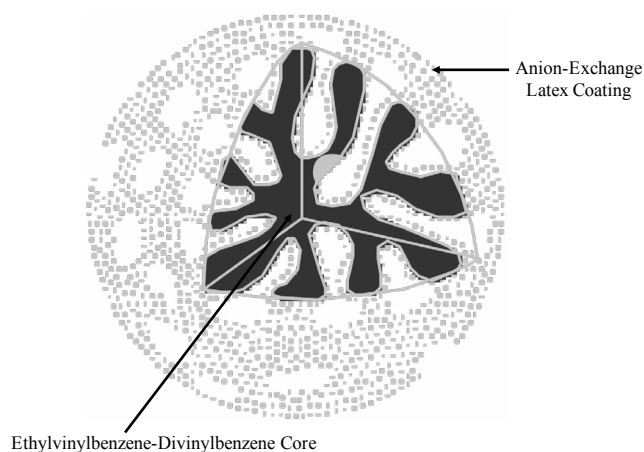
## Results

FIGURE 1. Resin Bead Structure for the Dionex IonPac® AS15



Dionex IonPac® AS15 packing is composed of a highly crosslinked core and an anion-exchange layer grafted to the surface, as shown in Figure 1. The AS15 column substrate is composed of a 9 $\mu$ m macroporous resin bead consisting of ethylvinylbenzene crosslinked with 55% divinylbenzene. The anion-exchange layer is functionalized with alkanol quaternary ammonium groups.

FIGURE 2. Resin Bead Structure for the Dionex IonPac® AS28-4 $\mu$ m



Dionex IonPac® AS28-4 $\mu$ m packing is a unique structure composed of a highly crosslinked core and a MicroBead™ anion-exchange layer attached to the surface, as shown in Figure 2. The AS28-4 $\mu$ m column substrate is a 4 $\mu$ m supermacroporous resin bead consisting of ethylvinylbenzene crosslinked with 55% divinylbenzene. The anion-exchange layer is functionalized with alkanol quaternary ammonium groups and has a controlled thickness, which results in excellent mass transfer characteristics and consequently, highly efficient peaks.

**FIGURE 3. Dionex IonPac® AS28-4µm 4x150mm Versus AS15 4x250mm: Comparison of Selectivity Using Isocratic Eluent**

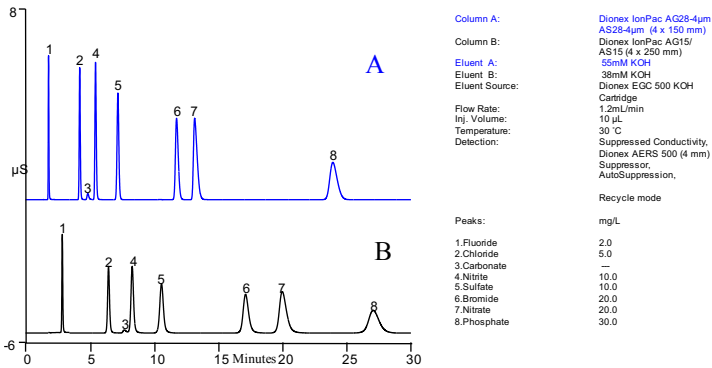


Figure 3A and B illustrates the common anion selectivity for the IonPac®AS15 and AS28-4µm column using recommended eluent at the same linear velocity. Notice the excellent peak shape for the AS28-4µm due to controlled thickness of MicroBead™ and 4µm supermacroporous resin beads.

**FIGURE 4. Dionex IonPac® AS28-4µm 4x150mm Versus AS15 4x250mm Column: Comparison of Selectivity and Sensitivity Using Gradient Chromatography**

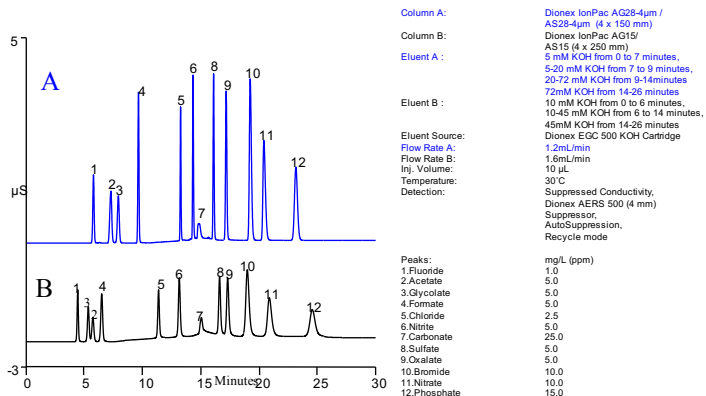


Figure 4A and B demonstrates common inorganic anion and organic acid selectivity for the IonPac®AS15 and AS28-4µm column. The IonPac®AS15 and AS28-4µm are high-capacity, hydroxide-selective anion-exchange columns designed for the determination of trace-level concentrations of inorganic anions and low molecular weight organic acids in high-purity water matrices encountered in the semiconductor and power generation industries. Both columns offer excellent retention of fluoride from the water dip and use a simple hydroxide gradient easily supplied by the Eluent Generator (EG). However, notice the excellent peak shape and better sensitivity offered by the IonPac AS28-4µm column.

**FIGURE 5. Effect of Increasing Flow Rate on Analysis Speed Using Dionex IonPac AS28-4µm and Isocratic Eluent**

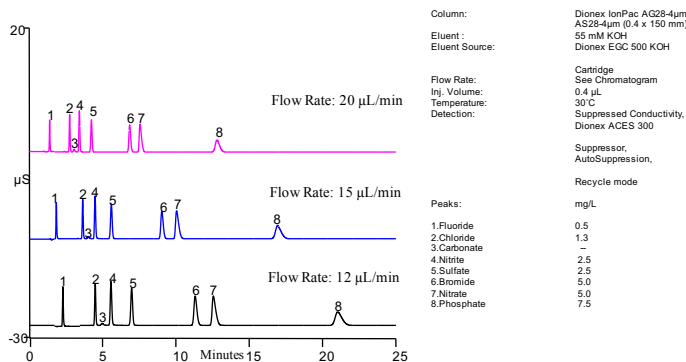


Figure 5 shows how analysis time can be reduced by increasing the flow rate without changing the common inorganic anion selectivity.

**FIGURE 6. Optimization of a Hydroxide Gradient and Flow Rate to Reduce the Analysis Time**

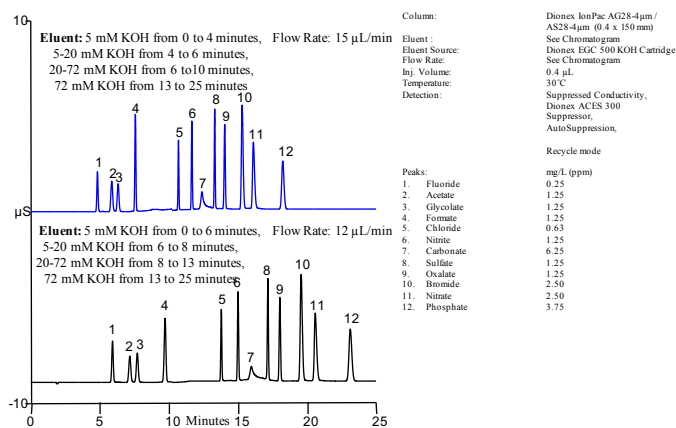


Figure 6 demonstrates optimization of a hydroxide gradient and flow rate to reduce analysis time for the common inorganic anions and organic acids encountered in the semiconductor and power generation industries.

**FIGURE 7. Analysis of Trace Organic Acids and Inorganic Anions Using a Hydroxide Gradient, Large Loop Injection, and Dionex IonPac® AS28-4µm 2 mm Column**

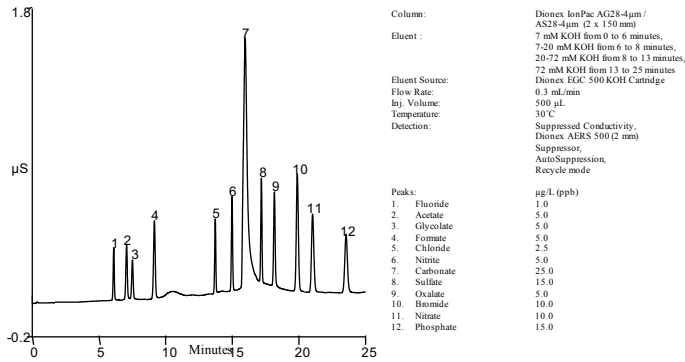


Figure 7 illustrates the separation of inorganic anions and low molecular weight organic acids in a high-purity water matrix using large loop injection. Separation was achieved using a potassium hydroxide gradient with suppressed conductivity detection at a temperature of 30 °C. Low µg/L (ppb) levels of these analytes can easily be quantified using a 500µL injection on an IonPac® AS28-4µm (2x 150 mm) column. The conductivity of the hydroxide eluent can be suppressed to a very low background, facilitating trace-level analysis.

**FIGURE 8. Analysis of Toothpaste Samples Using a Hydroxide Gradient and Dionex IonPac® AS28-4µm 4 mm Column**

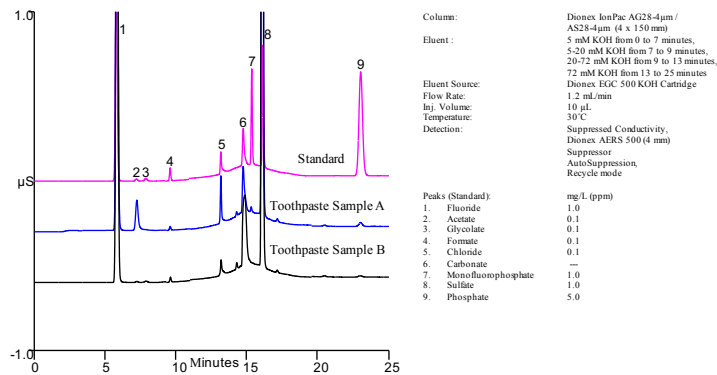


Figure 8 illustrates that the IonPac® AS28-4µm column can also be used for analysis of tooth paste samples.

**FIGURE 9. Separation of Various Environmental Anions Using a Hydroxide Gradient and Dionex IonPac® AS28-4µm 0.4 mm Column**

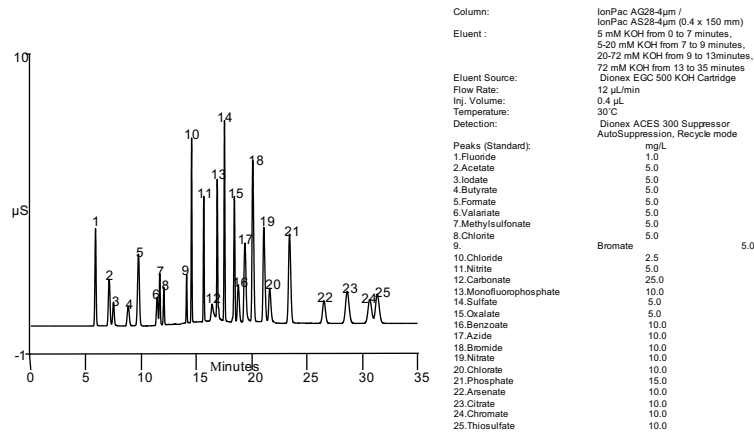


Figure 9 illustrates the separation of various environmental anions using hydroxide gradient.

## Conclusion

Dionex IonPac® AS28-4µm offers the following advantages relative to IonPac® AS15 column:

- Improved peak shape, efficiency and resolution
- Superior column selectivity with improved sensitivity
- Superior resolution of early-eluting anions (fluoride, glycolate, acetate, and formate)
- Use large-loop injection for ppb-level analysis
- Recommended column for high-purity water matrices encountered in the semiconductor and power generation industries

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

©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 2730 3006  
**Canada** +1 800 530 8447  
**China** 800 810 5118 (free call domestic)  
 400 650 5118

**Denmark** +45 70 23 62 60  
**Europe-Other** +43 1 333 50 34 0  
**Finland** +358 10 3292 200  
**France** +33 1 60 92 48 00  
**Germany** +49 6103 408 1014  
**India** +91 22 6742 9494  
**Italy** +39 02 950 591

**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

**Thermo**  
 SCIENTIFIC

A Thermo Fisher Scientific Brand

PN71981-EN 08/16S