



# Dionex Guardcap Products

065705 Revision 04 • March 2018

**thermo**scientific

*For Research Use Only. Not for use in diagnostic procedures.*

# **Product Manual**

**for**

## **Dionex Guardcap H**

(Package of 125, Item # 302504)

(Package of 12 with sample transfer line, Item # 302765)

## **Dionex Guardcap Na**

(Package of 125, Item # 302797)

(Package of 12 with sample transfer line, Item # 302796)

## **Dionex Guardcap HRP**

(Package of 125, Item # 302799)

(Package of 12 with sample transfer line, Item # 302798)

© 2018 Thermo Fisher Scientific Inc. All rights reserved.

All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

Thermo Fisher Scientific Inc. provides this document to its customers with a product purchase to use in the product operation. This document is copyright protected and any reproduction of the whole or any part of this document is strictly prohibited, except with the written authorization of Thermo Fisher Scientific Inc.

The contents of this document are subject to change without notice. All technical information in this document is for reference purposes only. System configurations and specifications in this document supersede all previous information received by the purchaser.

Thermo Fisher Scientific Inc. makes no representations that this document is complete, accurate or error free and assumes no responsibility and will not be liable for any errors, omissions, damage or loss that might result from any use of this document, even if the information in the document is followed properly.

This document is not part of any sales contract between Thermo Fisher Scientific Inc. and a purchaser. This document shall in no way govern or modify any Terms and Conditions of Sale, which Terms and Conditions of Sale shall govern all conflicting information between the two documents.

Revision History:

Revision 01, December, 2016, Original Publication;

Revision 02, March, 2017, Added the item number for 12 pack (# 302765).

Revision 03, June, 2017, Added Dionex Guardcap Na products.

Revision 04, March, 2018, Added Dionex Guardcap HRP products.

## Safety and Special Notices

Make sure you follow the precautionary statements presented in this guide. The safety and other special notices appear in boxes.

Safety and special notices include the following:



**SAFETY**

*Indicates a potentially hazardous situation which, if not avoided, could result in death or serious injury.*



**WARNING**

*Indicates a potentially hazardous situation which, if not avoided, could result in damage to equipment.*



**CAUTION**

*Indicates a potentially hazardous situation which, if not avoided, may result in minor or moderate injury. Also used to identify a situation or practice that may seriously damage the instrument, but will not cause injury.*



**NOTE**

*Indicates information of general interest.*

**IMPORTANT**

*Highlights information necessary to prevent damage to software, loss of data, or invalid test results; or might contain information that is critical for optimal performance of the system.*

**Tip**

*Highlights helpful information that can make a task easier.*

# Contents

|  |           |
|--|-----------|
| <b>1. Introduction .....</b>   | <b>6</b>  |
| 1.1 Overview of Automatic Sample Preparation using Dionex Guardcap™<br>Products and the Dionex AS-DV Autosampler .....                 | 6         |
| 1.2 Deionized Water Requirements for Ion Chromatography Methods .....  | 6         |
| 1.3 Setup .....  | 7         |
| 1.3.1 Vials and Caps .....   | 7         |
| 1.3.2 Setup of AS-DV .....   | 7         |
| 1.3.3 Preparing Sample Vials and Dionex Guardcap vial caps .....   | 10        |
| 1.3.4 Filling the PolyVial and Installing the Dionex Guardcap vial cap .....   | 10        |
| 1.4 In-line Sample Preparation using Dionex Guardcap vial caps<br>with PolyVials in the AS-DV Autosampler.....                         | 11        |
| 1.4.1 Dionex Guardcap H – Removal of Metals and Other Cations.....   | 11        |
| 1.4.2 Dionex Guardcap H – Neutralization of Base .....   | 12        |
| 1.4.3 Dionex Guardcap Na – Determination of pH-sensitive Analytes with Removal of Metals<br>and Other Cationic Matrix Components ..... | 14        |
| 1.4.4 Dionex Guardcap Na – Determination of Anions in Matrices Containing Components that<br>Precipitate at low pH, Humic Acids .....  | 14        |
| 1.4.5 Dionex Guardcap HRP – Removal of Surfactants and Other Hydrophobic Matrix Components .....                                       | 15        |
| 1.5 Filtering of Particulates .....  | 15        |
| <b>2. Troubleshooting .....</b>  | <b>16</b> |
| <b>3. Reordering Information.....</b>  | <b>17</b> |
| 3.1 Dionex Guardcap Introductory Packs (each pack includes a 70 cm<br>reduced volume sample transfer line).....                        | 17        |
| 3.2 Dionex Guardcap Vial Caps .....  | 17        |
| 3.3 Accessories .....  | 17        |
| <b>Appendix A – Installing the Sample Transfer Line .....</b>  | <b>18</b> |

# 1. Introduction

## 1.1 Overview of Automatic Sample Preparation using Dionex Guardcap™ Products and the Dionex AS-DV Autosampler

Thermo Scientific™ Dionex™ Guardcap™ vial caps are designed to work with 5-mL Thermo Scientific Dionex PolyVials™ in Thermo Scientific Dionex AS-DV autosamplers to achieve inline removal of cations and neutralization of bases with and without sample acidification.

The Dionex AS-DV holds 50 vials that can be sampled in any order and multiple samples can be taken from each vial. A full description of operating features and instructions for use can be found in the Dionex AS-DV Autosampler Operator's Manual, Document No. 065259. The unique PolyVial sample vial incorporates a cap that acts as a simple piston to force sample out of the vial. This eliminates the need for an external sampling pump. This vial/cap system can be used for inline sample preparation and filtration by using Dionex Guardcap vial caps.

Dionex Guardcap vial caps fit 5-mL PolyVials and contain either H<sup>+</sup>-form cation exchange resin, Na<sup>+</sup>-form cation exchange resin, or a hydrophilic reversed phase resin in a high density polyethylene filtration matrix. The Dionex Guardcap H and Guardcap Na filters are brown in color, as compared to the white Dionex Guardcap HRP and Dionex Filtercap filters. Note that if filtration-only is needed, the standard filter cap contains a 20 µm filter to remove solid materials from the sample before injection. By default, the Dionex AS-DV comes with a sample transfer line that is 142 cm long. To maximize the ion exchange capacity in Dionex Guardcap products, the Dionex AS-DV autosampler should be configured with an onboard 6-port high pressure injection valve (Item # 068920) and 35 cm reduced volume sample transfer line (Item # 22120-60051) in order to minimize sample volumes required for adequate flushing between samples. If an onboard injection valve is not available, direct connection to injection valves mounted in IC systems should be made with a 70 cm reduced volume sample transfer line (Item # 22120-60055).

## 1.2 Deionized Water Requirements for Ion Chromatography Methods

For all procedures requiring deionized water, use ASTM Type 1 (18 megohm-cm) filtered and deionized water that meets the following specifications.

**Table 1 Specifications for ASTM Type I water (with addition of a limit on metals)**

|                        |                  |
|------------------------|------------------|
| Resistivity            | > 18.0 megohm-cm |
| Organics-TOC           | <10 ppb          |
| Iron-Transition metals | <1 ppb           |
| Pyrogens               | <0.03 Eu/mL      |
| Particulates >0.2 µm   | < 1 units/mL     |
| Colloids-silica        | <10 ppb          |
| Bacteria               | 1 cfu/mL         |

## 1.3 Setup



NOTE

*The standard 100  $\mu$ L sample transfer line in the AS-DV autosampler should be replaced with a shorter and narrower, reduced volume sample transfer line in order to make optimum use of the capacity of Dionex Guardcap products. Please see the Setup of AS-DV and Appendix A for instructions to change the sample transfer line.*

### 1.3.1 Vials and Caps

Dionex Guardcap vial caps are available for 5-mL PolyVials. Use of an insertion tool (Item # 037987), included in the Dionex AS-DV Ship Kit (Item # 068887), simplifies insertion of the caps. Section 3, Reordering Information, contains part numbers for Dionex Guardcap vial caps, PolyVials, and other accessories.

Dionex Guardcap H and Guardcap Na filters are brown in color, as compared to the white Dionex Guardcap HRP and Filtercap filters. The brown color may appear darker for wet Dionex Guardcap vial caps than for dry vial caps, however color does not affect performance.

Dionex Guardcap products are shipped slightly damp. While Dionex Guardcap H and Guardcap Na perform equally well from a dry state, Dionex Guardcap HRP must remain damp and instructions are provided for re-hydrating Dionex Guardcap HRP vial caps that have dried out in Section 1.4.5.

The product packaging is an opaque, resealable bag and it is recommended to re-seal the bag when not in use to avoid airborne contamination (such as from amines) and to maintain moisture.

### 1.3.2 Setup of AS-DV

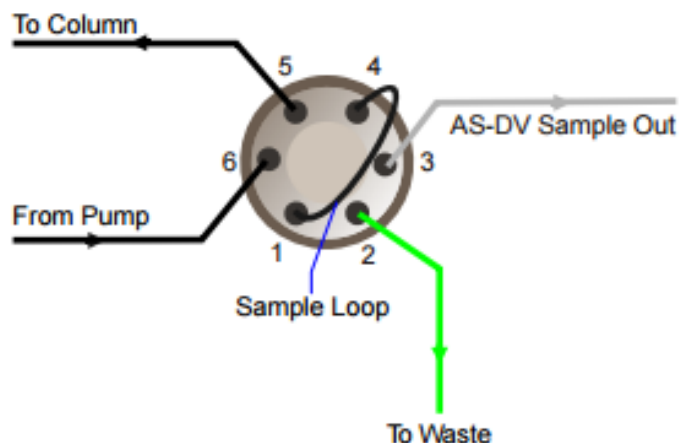
#### 1.3.2.1 Injection valve installed in AS-DV

For best performance and the highest use of the capacity of Dionex Guardcap products, the AS-DV autosampler should be setup with the on-board injection valve and a reduced volume sample transfer line assembly (Item # 22120-60051). Other configurations with direct connection to injection valves located in IC instruments are discussed in section 1.3.2.2. The standard AS-DV sample line will reduce the capacity of Dionex Guardcap products by up to 80%. Large flush volumes and extra tubing volume will exhaust the capacity, depending on the ionic strength of the sample matrix.

Dionex Guardcap products can be used at flow rates up to 3 mL/min but optimum performance is achieved at 1 mL/min. The following instructions describe setup of the hardware and method parameters for accomplishing sample preparation using Dionex Guardcap products.

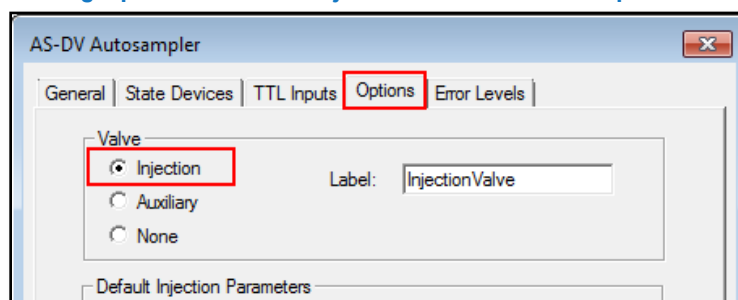
1. Install a 6-port injection valve in the AS-DV autosampler (see Figure 1). Note that the valve installs upside down.

Figure 1 AS-DV 6-port Injection Valve Plumbing



- a. Install a new, reduced volume AS-DV sample transfer line assembly (Item # 22120-60051). This is a narrower Teflon line shortened to 35 cm. See Appendix A or the AS-DV Operators manual for detailed installation instructions.
2. Add an AS-DV to the Chromeleon Instrument Configuration. Confirm the onboard valve is identified as an Injection Valve in the AS-DV options (Figure 2):

Figure 2 Confirming Optional Onboard Injection Valve in AS-DV Options



3. The sample delivery volume in the AS-DV will be controlled by the Chromeleon Instrument Method/Program using the command line below (Figure 3). The injection volume in the Sequence is not used. The actual injection volume is the volume of the sample loop installed on the injection valve.

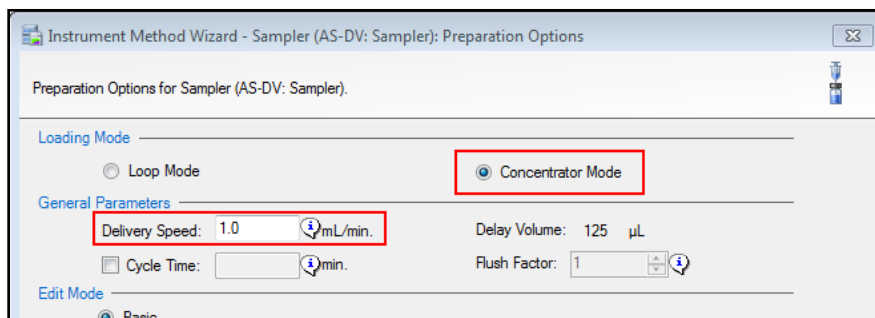
Figure 3 Setting Sample Delivery Volume to 175 µL in the Instrument Method

| Command               | Value           |
|-----------------------|-----------------|
| Sampler.FlushFactor   | 1               |
| Sampler.DeliverSpeed  | 1.0 [ml/min]    |
| Pump_ECD.Flow         | 1 [ml/min]      |
| Sampler.LoadPosition  |                 |
| Sampler.DeliverSample | Volume=175 [µl] |
| Sampler.EndSamplePrep |                 |



4. The loading flow rate (Delivery Speed) can be specified either in the Chromeleon Instrument Method Wizard (Figure 4) or directly in the Instrument Method/Program (Figure 5).
  - a. Set Loading Mode = **Concentrator Mode**
  - b. Set Delivery Speed = **1.0 mL/min**

**Figure 4 Confirming Delivery Speed in Instrument Method Wizard**



**Figure 5 Confirming Delivery Speed in the Instrument Method/Program**

| Command               | Value           |
|-----------------------|-----------------|
| Sampler.FlushFactor   | 1               |
| Sampler.DeliverSpeed  | 1.0 [ml/min]    |
| Pump_ECD.Flow         | 1 [ml/min]      |
| Sampler.LoadPosition  |                 |
| Sampler.DeliverSample | Volume=175 [µl] |
| Sampler.EndSamplePrep |                 |

This configuration allows about 10x flushing per sample while maximizing the capacity available in Dionex Guardcap products.

The actual injection loop volumes can be in the range of 2–25 µL for 2-mm or 4-mm operation.

### 1.3.2.2 Injection valve installed in IC system

1. Install a new, reduced volume AS-DV sample transfer line assembly (Item # 22120-60055). This is a narrower Teflon line shortened to 70 cm. Note: the AS-DV must be placed very close to the IC in order to use this sample line.
2. Setup is the same as steps 3 and 4 of section 1.3.2.1 except the sample delivery volume is defined as 350 µL in the Chromeleon Instrument Method/Program (Figure 6).

Figure 6 Setting Sample Delivery Volume to 350  $\mu\text{L}$  in the Instrument Method

| Command               | Value                        |
|-----------------------|------------------------------|
| Sampler.FlushFactor   | 1                            |
| Sampler.DeliverSpeed  | 1.0 [ml/min]                 |
| Pump_ECD.Flow         | 1 [ml/min]                   |
| Sampler.LoadPosition  |                              |
| Sampler.DeliverSample | Volume=350 [ $\mu\text{L}$ ] |
| Sampler.EndSamplePrep |                              |

### 1.3.3 Preparing Sample Vials and Dionex Guardcap vial caps

Thermo Scientific recommends thoroughly rinsing all PolyVials and caps with deionized water before filling to remove any traces of dust. Thermo Scientific does not recommend reusing vials or caps and does not guarantee their cleanliness or proper operation if reused.

To prevent the introduction of contaminants, extreme care must be used in cleaning. Use only ASTM Type I (18 megohm-cm) filtered and deionized water that meets the specifications listed in Table 1 for the cleaning procedures.

Dionex Guardcap vial caps are provided in a clean state (see Figure 7). However, for trace analysis it may be necessary to pre-rinse the vial caps. Dionex Guardcap vial caps can be flushed with deionized water using the same procedure used for filter caps in section 3.1 of the AS-DV Operator's manual. Using a clean syringe, force 10-50 mL of deionized water through the Dionex Guardcap from the top of the cap (Luer inlet). Be careful not to dislodge the Dionex Guardcap filter. Rinsed vial caps can be stored in a clean polyethylene container, filled with deionized water and capped.

Dionex PolyVials should be rinsed 10 times with a jet of deionized water. For trace analysis, rinsed vials should be allowed to soak in deionized water for at least 24 hours. When ready to use, empty the vial and rinse two times with sample.

### 1.3.4 Filling the PolyVial and Installing the Dionex Guardcap vial cap



**NOTE**

*To prevent contamination when doing trace-level analysis, use forceps to handle the vials and caps. Also, avoid touching any surface that will be wetted with sample.*

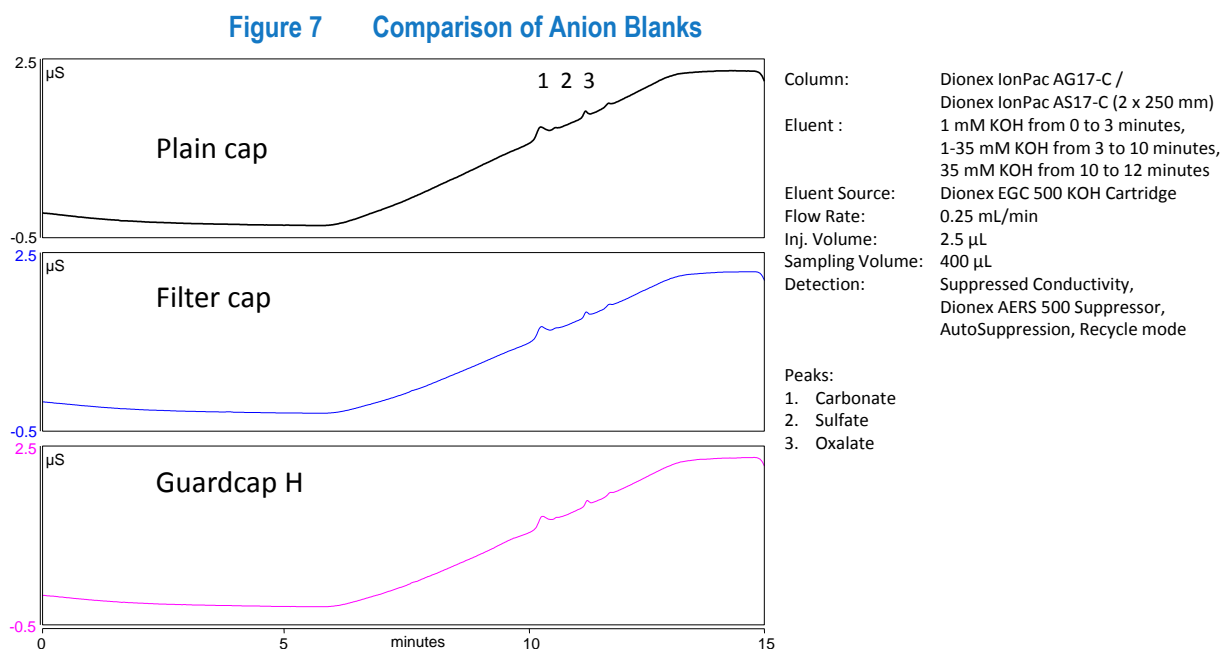
It is best to minimize the initial air bubble under the Dionex Guardcap vial cap when installed in the 5-mL PolyVial. The procedure below assures that adequate sample is placed in the vial, but does not overfill the vial.

1. Weigh 6.1-6.4 g sample/vial and install the Dionex Guardcap vial cap using the flat end of the installation tool so the cap is flush with the top of the vial. Note that this amount of sample is above the top of the vial tray. The top of the Dionex Guardcap vial cap must be flush with the top of the PolyVial in order for the vial to be properly sensed by the

autosampler. The backpressure applied by the user to install Dionex Guardcap products is higher than when installing a filter or plain cap. This is normal since Dionex Guardcap products contain ion exchange resin in the pores of the filter.

2. After installation of the Dionex Guardcap vial cap, invert the vial to confirm that no air bubble is freely floating under the cap.
3. Before loading the vial into the carousel, shake off any liquid that has been forced into the cap socket. Do not use laboratory wipes to blot liquid from the cap sockets as wipes can leave fibers that can clog the liquid flow path and cause backpressure.

Figure 7 shows carbonate, sulfate, and oxalate from a system blank. All peaks are less than 50 ppb. The chromatograms in the figure show a comparison using a plain cap, filter cap, and Dionex Guardcap H.



## 1.4 In-line Sample Preparation using Dionex Guardcap vial caps with PolyVials in the AS-DV Autosampler

The following application examples illustrate several types of methods for sample preparation.

### 1.4.1 Dionex Guardcap H – Removal of Metals and Other Cations

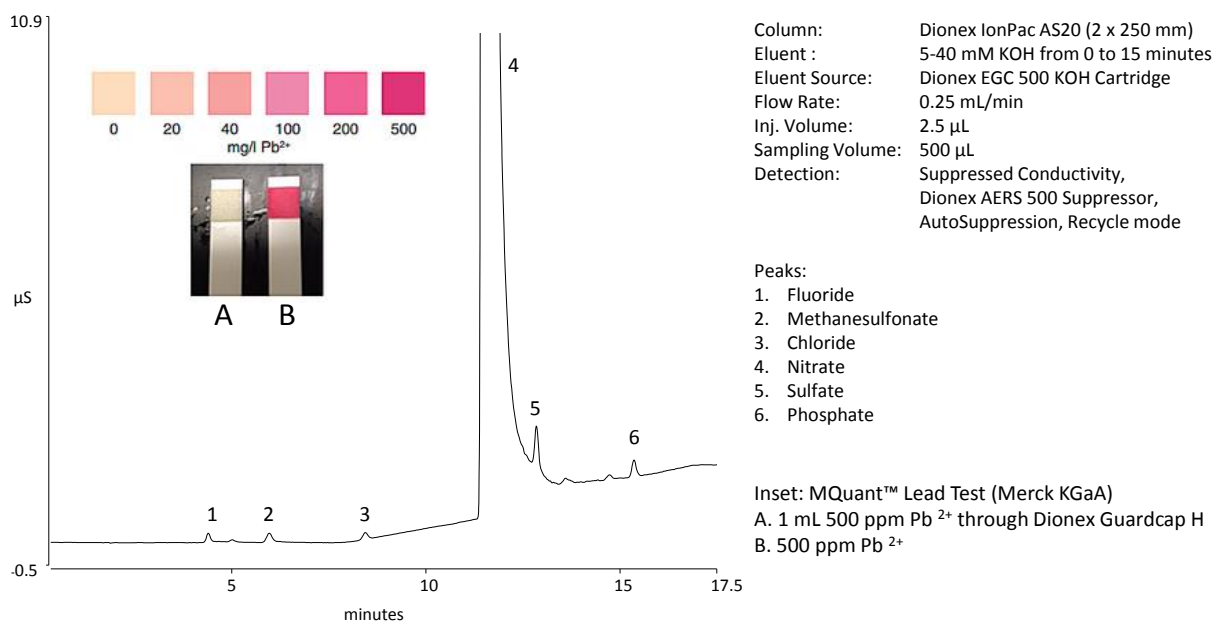
Dionex Guardcap H vial caps contain strong cation exchange resin in the hydronium ion form. One Dionex Guardcap H can neutralize up to 500 µL of 50 mM sodium hydroxide or remove 1000 mg/L divalent cation, i.e. calcium or transition metals, from 800 µL of sample. The following application examples illustrate several types of methods for sample preparation.

The Dionex Guardcap H resin is provided in the H<sup>+</sup>-form, meaning that alkali, alkaline earth, and transition metals in the cation form, and protonated weak bases including amines are exchanged for hydronium ion. One Dionex Guardcap H vial cap contains about 50 µEq of cation exchange capacity, available for removing cations from samples. As an example, one Dionex Guardcap H vial cap can remove 1000 ppm calcium from 800 µL of sample, or 500 ppm of a divalent cation from 1.6 mL of sample. It is important to note that the necessary capacity needed for an application must be calculated based on the sample volume passing through the Dionex Guardcap H vial cap rather than the volume of the injection loop.

The sample can become acidic in this mode as cations are exchanged for the hydronium ion. This application is useful to remove cations that can foul the anion suppressor, which is a cation exchanger, or cations that can build up by precipitation as insoluble salts on system components. This method of use is exactly analogous to using Dionex OnGuard™ II H Cartridges for offline sample preparation removal of metals and other higher valency cations.

Figure 8 below shows the determination of methanesulfonate in a lead nitrate matrix using Dionex Guardcap H sample preparation. The inset in Figure 8 shows the results of passing 1 mL of a 500 ppm Pb<sup>2+</sup> standard through a Dionex Guardcap H vial cap. Dionex Guardcap H effectively removes all of the Pb<sup>2+</sup> from the standard.

**Figure 8** Determination of Methanesulfonate in a Lead Nitrate Matrix Using Dionex Guardcap H Sample Preparation



#### 1.4.2 Dionex Guardcap H – Neutralization of Base

The Dionex Guardcap H resin is provided in the H<sup>+</sup>-form. When a cation such as sodium is removed from the sample, it is replaced by the hydronium ion and can neutralize bases that are present in the sample. This is especially useful for lowering the pH of high pH samples that will be injected into a gradient anion system.

Figure 9 below shows the band broadening effects of a high pH matrix injected onto a Dionex™ IonPac™ AS17-C column and the improvement using Dionex Guardcap H. The pH of the sample

is lowered to a pH that more closely approximates the initial conditions of the gradient, thus eliminating disturbances in the ion exchange equilibria. This is also useful for large loop injections, which for 2-mm systems, are injections volumes in the range of 25  $\mu\text{L}$ . When using standard loop injection, the most common effect of high hydroxide concentration in the sample is peak broadening of early-eluting peaks, most commonly fluoride.

**Figure 9 Neutralization of Base to Improve Peak Efficiency**

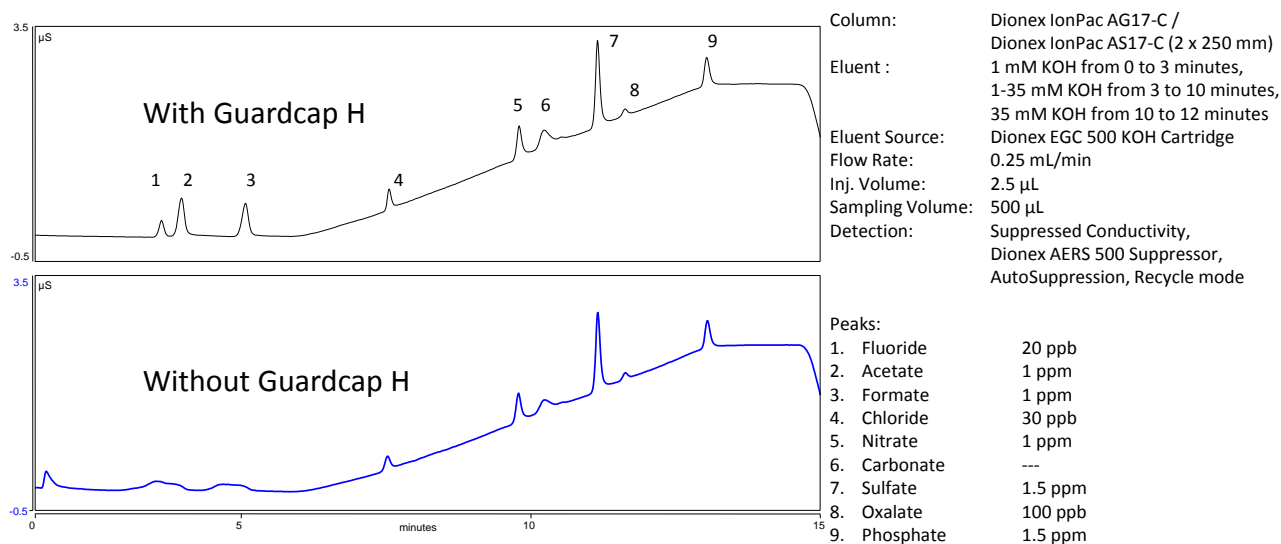
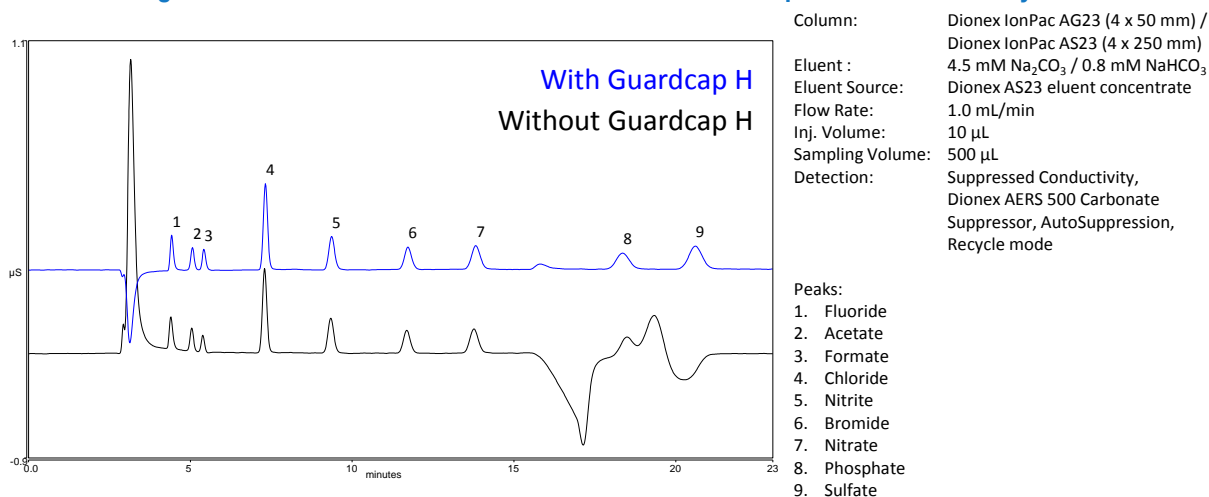


Figure 10 below shows the baseline disturbance caused by a high pH matrix injected onto a Dionex™ IonPac™ AS23 column and the improvement using Dionex Guardcap H.

**Figure 10 Neutralization of a 50 mM NaOH Matrix to Improve Peak Efficiency**



The capacity of Dionex Guardcap H for neutralization using the AS-DV flow rate of 1 mL/min is about 500  $\mu\text{L}$  of 50 mM NaOH. If the concentration of base in the sample is higher than about 50 mM, the sample should be diluted prior to treatment using Dionex Guardcap H.

### 1.4.3 Dionex Guardcap Na – Determination of pH-sensitive Analytes with Removal of Metals and Other Cationic Matrix Components

Dionex Guardcap Na vial caps contain the same strong cation exchange resin found in Dionex Guardcap H but in the sodium ion form. One Dionex Guardcap Na vial cap can remove 1000 mg/L divalent cation, i.e. calcium or transition metals, from 800  $\mu$ L of sample without acidification of the sample. The following application examples illustrate several types of methods for sample preparation.

The Dionex Guardcap Na resin is provided in the Na<sup>+</sup>-form, meaning that higher alkali, alkaline earth, and transition metals in the cation form, and protonated weak bases including amines are exchanged for sodium ion. The sodium content of the sample increases after treatment with Dionex Guardcap Na. One Dionex Guardcap Na vial cap contains about 50  $\mu$ Eq of cation exchange capacity, available for removing more highly retained cations from samples. As an example, one Dionex Guardcap Na vial cap can remove 1000 ppm calcium from 800  $\mu$ L of sample, or 500 ppm of a divalent cation from 1.6 mL of sample. It is important to note that the necessary capacity needed for an application must be calculated based on the sample volume passing through the Dionex Guardcap Na vial cap rather than the volume of the injection loop.

Some analytes are affected by the pH change through H<sup>+</sup>-form resin, most notably nitrite. Removal of metals from a sample requiring determination of nitrite must be accomplished through sodium-form resin, using Dionex Guardcap Na. Table 2 below shows a comparison of recoveries for nitrite using Dionex Guardcap H and Dionex Guardcap Na sample preparation. These are the same results found using Dionex OnGuard H and Dionex OnGuard Na products.

**Table 2 Comparison of Nitrite Recovery using Dionex Guardcap H and Dionex Guardcap Na**

| Sample   | Dionex Guardcap H |             |                            | Dionex Guardcap Na |             |                            |
|--|-------------------|-------------|----------------------------|--------------------|-------------|----------------------------|
|  | %Recovery*        | %RSD<br>N=7 | pH after<br>sample<br>prep | %Recovery*         | %RSD<br>N=7 | pH after<br>sample<br>prep |
| 1 ppm nitrite<br>in 500 ppm<br>calcium<br>(CaSO <sub>4</sub> ) | 66                | 9.3         | 2                          | 101                | 2.7         | 6                          |

*\*compared to 1 ppm nitrite standard in deionized water matrix*

### 1.4.4 Dionex Guardcap Na – Determination of Anions in Matrices Containing Components that Precipitate at low pH, Humic Acids

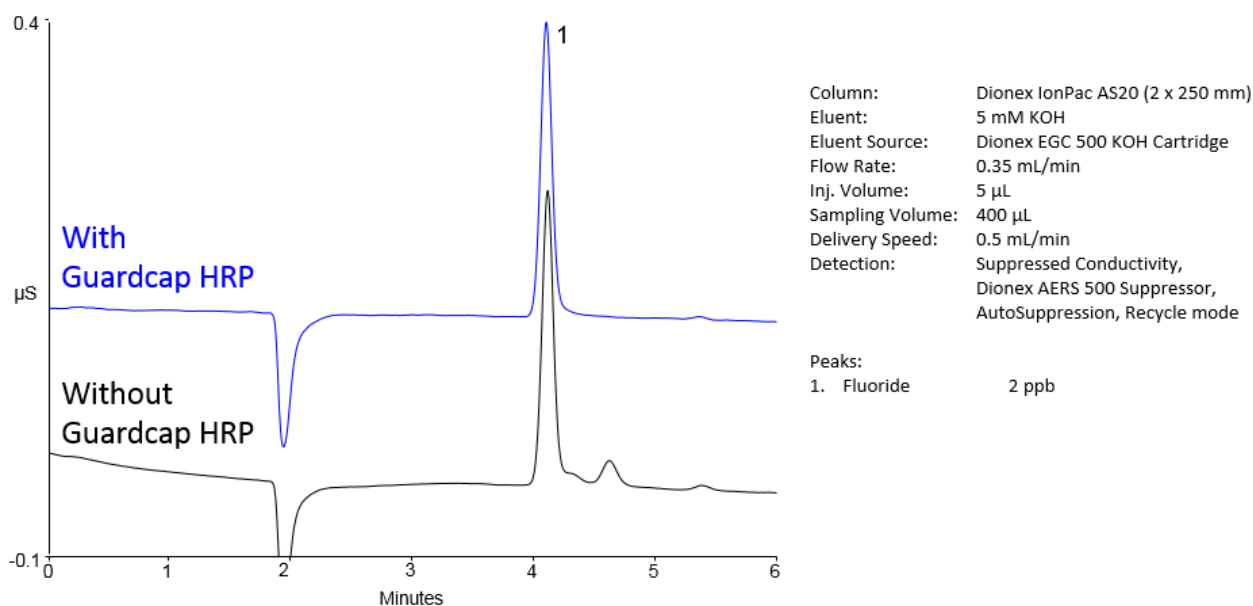
In another type of application, Dionex Guardcap Na allows removal of metals from samples containing other matrix components that can precipitate at low pH. An example is a natural water sample containing humic acids. Many humic acids precipitate below about pH 4, depending on their molecular weight and carboxylate:phenolate ratios.

### 1.4.5 Dionex Guardcap HRP – Removal of Surfactants and Other Hydrophobic Matrix Components

Dionex Guardcap HRP vial caps contain a proprietary styrene-divinylbenzene resin grafted with a hydrophilic monomer that helps maintain the pre-wetted state required for maximum retention of hydrophobic compounds. Dionex Guardcap HRP has the same selectivity as Dionex OnGuard RP. This resin is embedded in a high density polyethylene filtration matrix. Dionex Guardcap HRP filters are white. Dionex Guardcap HRP vial caps are pre-washed and treated so that they are pre-wetted and ready for use when opened. These products are shipped in re-sealable moisture-barrier bags so it is recommended to immediately reseal the bag after use. However, if they are open to air for more than 24 hours they can dry out and should be re-wet before use. Freshly opened Dionex Guardcap HRP vial caps can be removed from the bag, rinsed and stored in a sealed container such as a plastic tub with cap, as in Section 1.3.3 if desired. If Dionex Guardcap HRP vial caps have been open to the air for more than overnight, it is recommended to re-wet them with 1 mL 50% v/v methanol/deionized water then 1 mL deionized water using a Luer-slip syringe and then store them in a closed container.

Dionex Guardcap HRP can remove 0.01% Triton x-405 surfactant from 500  $\mu$ L of sample. Optimum matrix removal using Dionex Guardcap HRP is achieved using a reduced sampling flow rate of 0.25 – 0.5 mL/min. Use of higher flow rates will reduce the concentration of hydrophobic compounds that can be removed from the sample. Figure 11 shows the separation and detection of fluoride in a synthetic wastewater sample containing 0.005% Triton x-405.

**Figure 11** Determination of Fluoride in 0.005% Triton x-405 Using Dionex Guardcap HRP Sample Preparation



## 1.5 Filtering of Particulates

Similar to filter caps, all Dionex Guardcap products will remove better than 80% of particulate matter greater than 20  $\mu$ m in particle diameter.

## 2. Troubleshooting

| Symptom  | Possible Issue  | Action  |
|--|---|---|
| AS-DV chatters, grinds, or stalls during sample loading      | Sample loading flow rate (delivery speed) is set too high.  | Reduce loading flow rate to 1 mL/min.   |
|  | There is a clog in the sample loading path.   | Work through each connection to identify the source of the clog. Replace lines or components as necessary, taking care not to overtighten fittings.                                       |
|  |   | Clean or replace injection valve components (rotor and stator face assembly).   |
| Carryover  | Insufficient sample volume to adequately flush lines between samples.   | Ensure that a reduced volume sample transfer line was installed and that the instrument method/program specifies the appropriate sample delivery volume (see Section 1.3.2.1 or 1.3.2.2). |
|  |   | Increase sample delivery volume by 100 $\mu$ L in the instrument method/program.  |
| High blank   | Sample flow path may be contaminated.   | Flush the AS-DV with 5 mL deionized water using a clean PolyVial and plain vial cap.  |
|  | Dionex Guardcap vial caps may have become contaminated.   | Pre-rinse Dionex Guardcap vial caps with clean deionized water as noted in Section 1.3.3.   |
|  | Dionex Guardcap vial caps may be beyond the “Best If Used By” date.   | Check the “Best If Used By” date printed on the box label.  |
|  | Shared AS-DV used for both anions and cations can become contaminated with insoluble salt deposits that contribute to background peaks. | Replace sample loop, sample transfer line, and connecting PEEK tubing from the injection valve to the column.   |
|  | Large carbonate peak in blank.  | Use a Carbonate Removal Device (CRD) and/or degas samples and deionized water used for eluent.  |
| Hydrophobic compounds are not removed by Dionex Guardcap HRP | Sample loading flow rate (delivery speed) is set too high.  | Reduce the sample loading flow rate in the instrument method/program to 0.25 – 0.5 mL/min, or by half.  |



## 3. Reordering Information

### 3.1 Dionex Guardcap Introductory Packs

(each pack includes a 70 cm reduced volume sample transfer line)

| Part Number | Description   | Quantity |
|-------------|---|----------|
| 302765      | 5 mL Dionex Guardcap H vial caps with transfer line   | 12       |
| 302796      | 5 mL Dionex Guardcap Na vial caps with transfer line  | 12       |
| 302798      | 5 mL Dionex Guardcap HRP vial caps with transfer line | 12       |

### 3.2 Dionex Guardcap Vial Caps

| Part Number | Description                        | Quantity |
|-------------|------------------------------------|----------|
| 302504      | 5 mL Dionex Guardcap H vial caps   | 125      |
| 302797      | 5 mL Dionex Guardcap Na vial caps  | 125      |
| 302799      | 5 mL Dionex Guardcap HRP vial caps | 125      |

### 3.3 Accessories

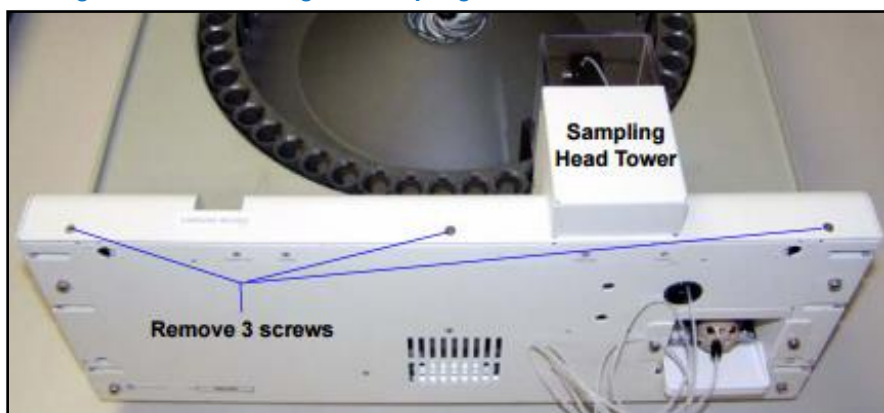
| Part Number | Description  | Quantity |
|-------------|--|----------|
| 22120-60051 | 35 cm Sample Transfer Line Assembly<br>For use with injection valves mounted in AS-DV        | 1        |
| 22120-60055 | 70 cm Sample Transfer Line Assembly<br>For use with injection valves mounted in an IC system | 1        |
| 038008      | 5 mL Dionex PolyVials  | 250      |
| 037987      | Tool, vial cap insertion   | 1        |
| 068920      | 6-port High Pressure Valve Kit   | 1        |
| 068947      | Vial holder tray for 5 mL vials  | 1        |

## Appendix A – Installing the Sample Transfer Line

The following tools are needed to complete this procedure: #1 Phillips screwdriver and 3-mm Allen wrench.

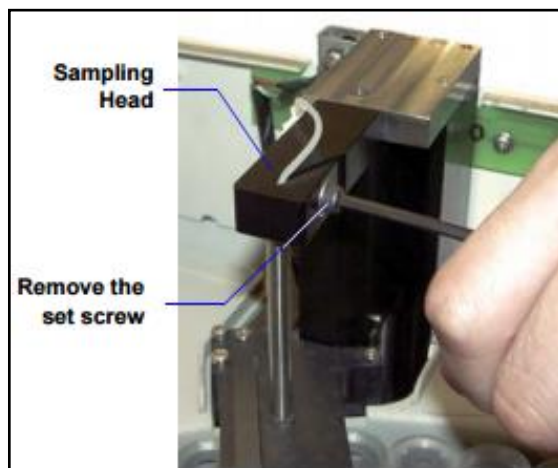
1. Disconnect the sample out line from the injection valve. Remove the fitting.
2. Press the Carousel Release button to retract the sampling tip to its fully up position.
3. Turn off the Dionex AS-DV power and disconnect the power cord. Raise the top cover.
4. Remove the Dionex AS-DV front cover by gripping the indentations at each side and pulling the cover straight off toward you.
5. Use a #1 Phillips screwdriver to remove the three flathead screws from the sampling head tower cover (see Figure 11). Lift up the cover to remove it.

**Figure 11** Removing the Sampling Head Tower Cover



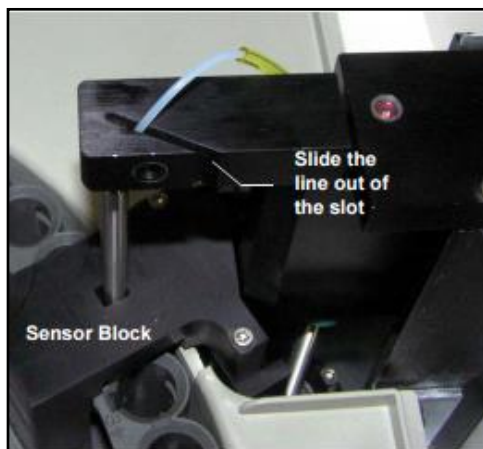
6. Use a 3-mm Allen wrench to remove the set screw on the sampling head (Figure 12).

**Figure 12.** Removing the Set Screw



7. Pull slightly on the sample line to gain a small amount of slack in the line. Then, slide the tubing out of the slot in the sampling head (Figure 13).

Figure 13 Sliding the Sample Transfer Line out of the Sampling Head Slot



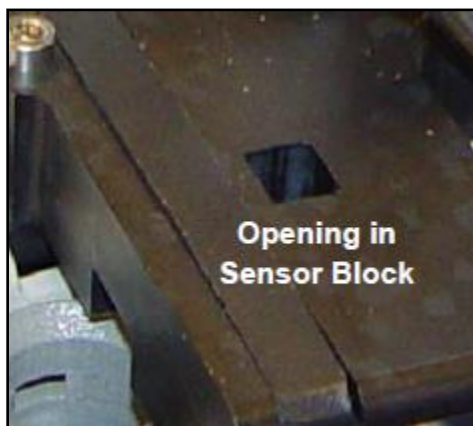
8. Push down on the sampling head to move the needle into the carousel a small amount. Twist the needle to allow the sampling tip to exit through the opening in the sensor block.
9. Unscrew the sampling tip from the needle and pull out and discard the tip and tubing.
10. Locate the new sample transfer line assembly. This assembly consists of a length of tubing with a new sampling tip installed on one end.
11. Thread the free end of the tubing up through the needle and screw the tip into place on needle (Figure 14).

Figure 14 Sampling Tip



12. Continue threading the free end of the tubing up through the sampling head and out the aperture on the front of the Dionex AS-DV.
13. Slide the needle into place and slide the sampling needle tubing back into the slot in the sampling head.
14. Align the flat sides of the sampling tip so that they fit through the opening in the sensor block (Figure 15).

Figure 15 Sensor Block Opening



15. Make sure that the needle is pushed as far as it will go up into the sampling head and then tighten the set screw firmly.

***NOTE: The set screw must be very tight to secure the sampling needle. This will require a firm turn of the set screw. Test the line to make sure it is tightly held.***

16. Reinstall the sampling head tower cover.
17. Reinstall the front cover and close the top cover.
18. Reconnect the sample out line to the valve.
19. Reconnect the power cord and turn on the power.