



**Thermo Scientific**

# **Dionex CarboPac PA100**

## **Column Product Manual**

**P/N: 065545-01**

**November 2013**

## **Product Manual**

**for**

### **Dionex CarboPac PA100 Analytical Columns**

2 x 250 mm, (P/N 057182)

4 x 250 mm, (P/N 043055)

9 x 250 mm, (P/N SP2089)

22 x 250 mm, (P/N SP2667)

### **Dionex CarboPac PA100 Guard Columns**

2 x 50 mm, (P/N 057183)

4 x 50 mm, (P/N 043054)

© 2013 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.

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

Revision History:

Revision 01, November, 2013, Revised Publication;

## 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>8</b>
<b>2. System Requirements and Installation .....</b>	<b>9</b>
2.1 System Requirements.....	9
2.1.1 System Requirements for 2 mm and 4 mm Operation .....	9
2.1.2 Installation of Disposable Electrode into an ED Cell, pH-Ag/AgCl Reference Electrode or PdH Reference Electrode.....	9
2.1.3 System Void Volume .....	9
2.2 The Injection Loop.....	9
2.2.1 The 4 mm System Injection Loop, 10 - 50 $\mu$ L .....	9
2.2.2 The 2 mm System Injection Loop, 2 - 15 $\mu$ L .....	9
2.3 The Dionex CarboPac PA100 Guard Column .....	10
2.4 Installing the Dionex CR-ATC Trap Column for Use with Dionex EGC.....	10
2.5 Eluent Storage.....	10
<b>3. Purity Requirements for Chemicals.....</b>	<b>11</b>
3.1 Deionized Water .....	11
3.2 Sodium Hydroxide .....	11
3.3 Sodium Acetate .....	11
<b>4. Before You Start.....</b>	<b>12</b>
4.1 The Most Important Rules .....	12
4.2 Initial Check List.....	12
<b>5. Preparation of Eluents and Standards.....</b>	<b>13</b>
5.1 Eluent E1: Deionized Water .....	13
5.2 Eluent E2: 1 M Sodium Hydroxide.....	13
5.2.1 Sodium Hydroxide Eluent Concentration .....	14
5.3 Eluent E3: 0.1M Sodium Hydroxide / 0.1M Sodium Acetate.....	15
5.4 Eluent E4: 0.1M Sodium Hydroxide / 0.5 M Sodium Acetate.....	15
5.5 Sample Preparation .....	16
5.6 Verification of System Cleanliness.....	17
5.6.1 System Background Check .....	17
5.6.2 Verification of Column Cleanliness .....	18
5.7 Carbohydrate Sialylated N-linked Alditol Standard .....	18

<b>6. Applications .....</b>	<b>19</b>
6.1 Production Test Chromatograms .....	20
6.2 Separation of Neutral and Sialylated Oligosaccharides .....	22
6.3 Separation of Fetuin N-linked Oligosaccharide Alditols .....	23
6.4 Isocratic Separation of Food Sugars and Food Alcohols .....	25
6.5 Separation of Apple and Orange Juice Sugars Using Dionex CarboPac PA100 .....	26
6.6 Oligosaccharide Profiling During Beer Production .....	28
6.7 Analysis of Oligosaccharides in Beer .....	29
6.8 Gradient Separation of Linear Polysaccharides .....	30
6.9 Fetuin Oligosaccharide Alditol Profiling .....	31
6.10 Antibodies N-Glycan Separations.....	32
6.11 Profiling of Complex Plant Carbohydrates: Profiling of Inulins.....	33
6.12 Sialyl Lactose (N-Acetylneuraminy-Lactose).....	34
6.13 Separation of Mannose 7 D1, D3 Isomers .....	35
6.14 Separation of Neutral N-Linked Oligosaccharides: Complex Neutral Structures.....	36
6.15 High Mannose Structures.....	37
<b>7. Troubleshooting Guide.....</b>	<b>38</b>
7.1 High Back Pressure.....	38
7.1.1 Finding the Source of High System Pressure .....	38
7.1.2 Replacing Column Bed Support Assemblies for 2 mm and 4 mm columns .....	38
7.1.3 Filter Eluent.....	39
7.1.4 Filter Samples.....	39
7.2 High Background.....	39
7.2.1 Preparation of Eluents .....	40
7.2.2 CR-ATC Column .....	40
7.2.3 A Contaminated Guard or Analytical Column .....	40
7.3 Poor Resolution.....	40
7.3.1 Loss of Column Efficiency .....	41
7.3.2 Shortened Retention Times .....	41
7.3.3 Loss of Resolution for early eluting peaks .....	42
7.3.4 Spurious Peaks .....	42
7.3.5 No Peaks, Poor Peak Area Reproducibility or too Small Peak Areas .....	43
7.3.6 Large Baseline Dip in the Chromatogram.....	43
7.3.7 Unidentified Peaks Appear with Expected Analyte Peaks.....	43
7.3.8 Decreased Detection Sensitivity.....	44
7.3.9 Excessive Gradient Rise.....	44
7.4 Reconditioning or Replacement of the Gold (conventional or disposable) Electrodes or Replacement of the Reference Electrode .....	44

<b>Appendix A – CarboPac PA100 Column Care .....</b>	<b>45</b>
A.1 Recommended Operation Pressures.....	45
A.2 Column Start-Up.....	45
A.3 Column Storage .....	45
A.4 CarboPac PA100 Column Cleanup.....	45
<b>Appendix B – Quality Assurance Reports .....</b>	<b>47</b>

# 1. Introduction

The Thermo Scientific™ Dionex™ CarboPac™ family of columns is designed to address the analytical requirements of a wide range of carbohydrate chemists. Underivatized carbohydrates are separated at high pH, in an approach that is unique to Thermo Scientific Dionex, simple to reproduce, and made possible by the pH stability of the Dionex CarboPac columns. The Dionex CarboPac PA100 (2-mm and 4-mm) is a high-resolution strong anion-exchange column developed for enhanced chromatography of oligosaccharides. It uses an advanced, high-performance, pellicular anion exchange resin to offer higher efficiency than the Dionex CarboPac PA1 for the separation of neutral and anionic oligosaccharide mixtures. If desalting of collected oligosaccharides is desired, the Dionex CarboPac PA100 column is the column of choice because lower salt concentrations are required for elution compared with the Dionex CarboPac PA1.

High-pH anion-exchange chromatography coupled with pulsed amperometric detection (HPAE-PAD) has been demonstrated to be effective in the separation and detection of carbohydrates ranging from small monosaccharides to branched oligosaccharides, and large linear polysaccharides. HPAE is proven to be capable of separating monosaccharides, positional, linkage and branch isomers of branched oligosaccharides, and homopolymer oligosaccharides that differ only in length. Each class of carbohydrates has different separation characteristics requiring different column chemistries.

## **Resin Characteristics:**

Particle Size:	8.5 µm
Pore Size:	microporous (< 10 Å)
Cross-linking:	55%
Ion exchange capacity:	90 µeq per 4x250 mm column
	18 µeq per 4x50 mm column
	22.5 µeq per 2x250 mm column
	4.5 µeq per 2x50 mm column
	455 µeq per 9x250 mm column
	2722 µeq per 22x250 mm column

## **Latex Characteristics:**

Functional Group:	quaternary ammonium functionalized latex
Latex Diameter:	275 nm
Latex Cross-linking:	6%

## **Typical Operating Parameters:**

pH range:	0–14
Temperature Limit:	4–60 °C
Pressure Limit:	5000 psi
Organic Solvent Limit:	100% compatible for cleaning only
Typical eluents:	High purity water (18.2 megohm-cm), sodium hydroxide, sodium acetate



## 2. System Requirements and Installation

### 2.1 System Requirements

#### 2.1.1 System Requirements for 2 mm and 4 mm Operation

The carbohydrate separations using the Dionex CarboPac PA100 columns are optimized for use with Dionex Ion Chromatography systems equipped with electro chemical detection. It is highly recommended to ensure that the systems used for carbohydrate analysis are metal-free. Metal ions from a metal system will contaminate the CarboPac column and may contaminate the working electrode. Running a CarboPac column on a metal system voids the column warranty.

#### 2.1.2 Installation of Disposable Electrode into an ED Cell, pH-Ag/AgCl Reference Electrode or PdH Reference Electrode

The 0.002 in. thick Teflon gaskets included in each package of disposable electrodes must be used, otherwise the disposable electrode product warranty is void. In addition, the quadruple waveform must be used for carbohydrate analysis otherwise the product warranty is void. Always wear gloves when handling electrodes. Never touch the electrode surface. To install a disposable working electrode and reference electrode (pH-Ag/AgCl or PdH) refer to Product Manual for Disposable Electrodes Doc. No. 065040, ICS-5000 Ion Chromatography System Manual Doc. No. 065342 and User's Compendium for Electrochemical Detection Doc. No. 065340.

#### 2.1.3 System Void Volume

When using Dionex CarboPac PA100 columns, it is particularly important to minimize system void volume. The system void volume for 2 mm columns should be scaled down to at least 1/4 of the system volume in a standard system designed for 4 mm columns (4 mm system). For best performance, all of the tubing installed between the injection valve and detector should be 0.005" (P/N 044221) i.d. PEEK tubing for 2 mm columns and 0.010" i.d. PEEK tubing (P/N 042260) for 4 mm columns. Minimize the lengths of all connecting tubing and remove all unnecessary switching valves and couplers.

### 2.2 The Injection Loop

#### 2.2.1 The 4 mm System Injection Loop, 10 - 50 $\mu$ L

For most applications on a 4 mm analytical system, a 10 – 50  $\mu$ L injection loop is sufficient. Generally, you should not inject more than 40 nanomoles of any analyte onto the 4 mm analytical column. Injecting larger amounts of an analyte can result in overloading the column which can affect the detection linearity. For low concentrations, larger injection loops can be used to increase sensitivity.

#### 2.2.2 The 2 mm System Injection Loop, 2 - 15 $\mu$ L

For most applications on a 2 mm analytical system, a 2 – 15  $\mu$ L injection loop is sufficient. Generally, you should not inject more than 10 nanomoles of any analyte onto a 2 mm analytical column. Injecting larger amount of an analyte can result in overloading the column which can affect the detection linearity. For low concentrations, larger injection loops can be used to increase sensitivity. The Dionex CarboPac PA100 2 mm requires a microbore HPLC system configuration. Install an injection loop one-fourth or less ( $<15 \mu$ L) of the loop volume used with a 4 mm analytical system.

### 2.3 The Dionex CarboPac PA100 Guard Column

A Dionex CarboPac PA100 Guard Column is normally used with the Dionex CarboPac PA100 analytical column. Retention times will increase by approximately 20% when a guard column is placed in-line before the analytical column under isocratic conditions. A guard column is utilized to prevent sample contaminants from eluting onto the analytical column. It is easier to clean or replace a guard column than an analytical column. Replacing the Dionex CarboPac PA100 Guard Column at the first sign of peak efficiency loss or decreased retention time will prolong the life of the Dionex CarboPac PA100 analytical column.

### 2.4 Installing the Dionex CR-ATC Trap Column for Use with Dionex EGC

For Dionex CarboPac PA100 applications using the Dionex EGC KOH cartridge, a Dionex CR-ATC Continuously Regenerated Trap Column (P/N 060477) should be installed at the Dionex EGC eluent outlet to remove trace level anionic contaminants from the carrier deionized water. See the Dionex CR-TC Product Manual (Document No. 031910) for instructions.

### 2.5 Eluent Storage

Dionex CarboPac PA100 columns are designed to be used with hydroxide eluent systems. Storage under a helium atmosphere ensures contamination free operation and proper pump performance (nitrogen can be used if eluents do not contain solvents).



**NOTE**

*For assistance, contact Technical Support for Thermo Scientific Dionex Products. In the U.S., call 1-800-346-6390. Outside the U.S., call the nearest Thermo Fisher Scientific office.*

## 3. Purity Requirements for Chemicals

Obtaining reliable, reproducible and accurate results requires eluents that are free from impurities and prepared only from the chemicals recommended below. Thermo Scientific cannot guarantee proper column performance when alternate suppliers of chemicals or lower purity water are utilized.

### 3.1 Deionized Water

The deionized water used to prepare eluents should be Type I reagent grade water with a specific resistance of 18.2 megaohm-cm. The water should be free from ionized impurities, organics, microorganisms and particulate matter larger than 0.2  $\mu\text{m}$ . The availability of UV treatment as a part of the water purification unit is recommended. Follow the manufacturer's instructions regarding the replacement of ion exchange and adsorbent cartridges. All filters used for water purification must be free from electrochemically active surfactants. Expanding their period of use beyond the recommended time may lead to bacterial contamination and as a result, a laborious cleanup may be required. Use of contaminated water for eluents can lead to high background signals and gradient artifacts.

### 3.2 Sodium Hydroxide

Use 50% w/w sodium hydroxide (Certified Grade, Fischer Scientific P/N UN 1824) for preparation.

### 3.3 Sodium Acetate

Thermo Scientific highly recommends the use of Dionex Sodium Acetate Reagent (P/N 059326) for carbohydrate analysis. Thermo Scientific cannot guarantee proper detection performance when different grades or alternate suppliers of sodium acetate are utilized.

## 4. Before You Start

### 4.1 The Most Important Rules

ALWAYS	use 50% NaOH solution rather than NaOH pellets to make eluents
ALWAYS	use dedicated glassware and disposable glass or plastic ware for volume adjustments.
ALWAYS	keep your NaOH eluent blanketed with helium or nitrogen. Prepare new NaOH eluent if left unblanketed for more than 30 minutes.
ALWAYS	use EGC-KOH generated eluent when possible to avoid any eluent preparation issues
ALWAYS	pull at least 40 mL of new eluent through the lines when changing eluent or adding fresh eluent. This will ensure that your fresh eluent is primed through the lines up to the pump heads.
ALWAYS	verify the equilibration time necessary prior to injection to avoid baseline issues or artifacts or to avoid unnecessary increase in total method time.
NEVER	go to the next step of the installation if the previous step has failed.
NEVER	start an installation with any of the check list items below missing
NEVER	use ‘communal’ filtration units or filters made of unknown or unsuitable (cellulose derivatives, polysulfone) materials.
NEVER	use methanol or other organic solvents as rinse fluid in the autosampler. Use only water, replaced daily.
NEVER	run above 60°C or 5000 psi.

### 4.2 Initial Check List

These items **MUST** be available in your lab. The absence of any of these may compromise your analysis.

- ☐ Laboratory water unit delivering 18.2 megaohm-cm water at the installation site.
- ☐ Vacuum pump available for use with the vacuum filtration units.
- ☐ Sterile packed Nalgene Filtration units (pore size: 0.2 µm, filtered material: Nylon), 1 L funnel size
- ☐ Inert gas cylinder (helium or nitrogen) with a regulator valve (ca. 0–200 psi at the low pressure side) and the appropriate size adaptors plus tubing.
- ☐ Sterile-packed 10 mL and 25 mL disposable pipets and suitable pipeting bulbs or pumps.
- ☐ Disposable, plastic (PE) large-size (at least 20 mL) syringe for priming the pump.
- ☐ Plastic eluent bottles.

## 5. Preparation of Eluents and Standards

**NOTE**

*Always sanitize the entire analyzer with 2M NaOH prior to initial start-up and after idle periods.*

Obtaining reliable, consistent and accurate results requires eluents that are free from ionic and electrochemically active impurities. Chemicals and deionized water used to prepare eluents must be of the highest purity available. Maintaining low trace impurities and low particle levels in eluents also help to protect your ion exchange columns and system components. Thermo Scientific cannot guarantee proper column performance when the quality of the chemicals, solvents and water used to prepare eluents is substandard.

### 5.1 Eluent E1: Deionized Water

Vacuum degas the water by placing the eluent reservoir in a sonicator and drawing a vacuum on the filled reservoir with a vacuum pump. Vacuum degas the reservoir for 5–10 minutes while sonicating. Note: Degassing by vacuum filtration through a 0.2  $\mu\text{m}$  filter is a good alternative to degassing in a sonicator. Cap each bottle after the degassing step and minimize the length of time the bottle is opened to the atmosphere.

### 5.2 Eluent E2: 1 M Sodium Hydroxide

**NOTE**

*DO NOT prepare NaOH eluents from sodium hydroxide pellets! The pellets are coated with a layer of carbonate.*

Always store degassed NaOH eluents in plastic eluent bottles blanketed with helium or nitrogen to avoid carbon dioxide contamination from the air. Carbonate in the eluent can significantly reduce retention times for carbohydrates.

### 5.2.1 Sodium Hydroxide Eluent Concentration

#### Gravimetric Method

When formulating eluents from 50% sodium hydroxide, Thermo Scientific recommends weighing out the required amount of 50% sodium hydroxide. Use the assayed concentration value from the sodium hydroxide bottle.

Example: To make 1 L of 1 M NaOH use 80.02g of 50% sodium hydroxide:

$$\text{For 1 M: } \frac{1 \text{ mole/L} \times 40.01 \text{ g/mole}}{50\%} = 80.02 \text{ g diluted to 1 L}$$

#### Volumetric Method

Although it is more difficult to make precise carbonate-free eluents for gradient analysis volumetrically, you may choose to use the following formula to determine the correct volume of 50% sodium hydroxide to be diluted.

$$g = dvr$$

Where    g = weight of sodium hydroxide required (g)  
             d = density of concentrated solution (g/mL)  
             v = volume of the 50% sodium hydroxide required (mL)  
             r = % purity of the concentrated solution

Example: To make 1 L of 1M NaOH use 52.3 mL of 50% sodium hydroxide:

$$\text{For 1 M: } \frac{1 \text{ mole/L} \times 40.01 \text{ g/mole}}{50\% \times 1.53^* \text{ g/mL}} = 52.3 \text{ mL diluted to 1 L}$$

\* This density applies to 50% NaOH. If the concentration of the NaOH solution is significantly different from 50%, the gravimetric method should be used instead.

#### Sodium Hydroxide Eluents

Dilute the amount of 50% (w/w) NaOH Reagent specified in [Table 1](#). Prepare CarboPac Eluents with degassed, deionized water (18.2 megohm-cm) to a final volume of 1,000 mL using a volumetric flask. Avoid the introduction of carbon dioxide from the air into the aliquot of 50% (w/w) NaOH bottle or the deionized water being used to make the eluent. Do not shake the 50% (w/w) NaOH bottle or pipette the required aliquot from the top of the solution where sodium carbonate may have formed.

**Table 1** Mass or Volume of NaOH Required to Make 1 L of Common Eluents

Eluent Concentration (M)	NaOH (50%) (g)	NaOH (50%) (mL)
0.15	12.0	7.8
0.5	40.0	26.2
0.8	64.0	41.8
1.0	80.0	52.3

### 5.3 Eluent E3: 0.1M Sodium Hydroxide / 0.1M Sodium Acetate

To maintain baseline stability, it is important to keep the sodium hydroxide concentration constant during the sodium acetate gradient, because acetate has no buffering capacity at high pH. This is achieved by making the eluents as follows:

Eluent A: x mM NaOH  
 Eluent B: x mM NaOH, y mM NaOAc

To make one (1) liter of 0.1M sodium hydroxide/ 0.1M sodium acetate, dispense approximately 800 mL of DI water into a 1 L volumetric flask. Vacuum degas for approximately 5 minutes. Add a stir bar and begin stirring. Weigh out 8.2 g anhydrous, crystalline sodium acetate (Thermo Scientific Dionex Sodium Acetate Reagent, P/N 059326). Add the solid acetate steadily to the briskly stirring water to avoid the formation of clumps which are slow to dissolve. Once the salt has dissolved, remove the stir bar with a magnetic retriever. Add DI water to the flask to bring the volume to the 1 L mark.

Vacuum filter the solution through a 0.2  $\mu$ m Nylon filter. This may take a while as the filter may clog with insoluble material from the sodium acetate. Using a plastic tip volumetric pipet, measure 5.2 mL of 50% (w/w) sodium hydroxide solution. Dispense the sodium hydroxide solution into the acetate solution about 1 inch under the surface of the acetate solution. The eluent should be kept blanketed under helium or nitrogen at 34 to 55 kPa (5–8 psi) at all times, and last about 1 week.

### 5.4 Eluent E4: 0.1M Sodium Hydroxide / 0.5 M Sodium Acetate

To maintain baseline stability, it is important to keep the sodium hydroxide concentration constant during the sodium acetate gradient, because acetate has no buffering capacity at high pH. This is achieved by making the eluents as follows:

Eluent A: x mM NaOH  
 Eluent B: x mM NaOH, y mM NaOAc

To make one (1) liter of 0.1M sodium hydroxide/ 0.5M sodium acetate, dispense approximately 800 mL of DI water into a 1 L volumetric flask. Vacuum degas for approximately 5 minutes. Add a stir bar and begin stirring. Weigh out 41.0 g anhydrous, crystalline sodium acetate (Thermo Scientific Dionex Sodium Acetate Reagent, P/N 059326). Add the solid acetate steadily to the briskly stirring water to avoid the formation of clumps which are slow to dissolve. Once the salt has dissolved, remove the stir bar with a magnetic retriever. Add DI water to the flask to bring the volume to the 1 L mark.

Vacuum filter the solution through a 0.2 µm Nylon filter. This may take a while as the filter may clog with insoluble material from the sodium acetate. Using a plastic tip volumetric pipet, measure 5.2 mL of 50% (w/w) sodium hydroxide solution. Dispense the sodium hydroxide solution into the acetate solution about 1 inch under the surface of the acetate solution. The eluent should be kept blanketed under helium or nitrogen at 34 to 55 kPa (5–8 psi) at all times, and last about 1 week.

**NOTE**

*Thermo Scientific recommends the use of dedicated glassware, pipets and filtration apparatus for exclusive use in the preparation of carbohydrate eluents.*

## 5.5 Sample Preparation

The Dionex CarboPac columns are strong anion exchangers. Thus, the normal caveats applicable to ion exchange chromatography apply to these columns. High salt concentrations in the samples should be avoided wherever possible. Special care should be taken with samples containing high concentrations of anions, which are strong eluents for the Dionex CarboPac columns (e.g. chloride, carbonate, phosphate, etc.). It is best to avoid extremes of sample pH (especially extremely acidic samples). The presence of anionic detergents (e.g., SDS) in samples should be avoided entirely. Nonionic or cationic detergents may be acceptable in low concentrations.

Matrix Interferent	Effect	Possible Removal
Hydroxylated compounds (e.g. Tris buffers, alcohols)	PED-active (interferes with carbohydrate detection)	Dialysis dilution
Halides	Will bind to column, may affect retention time of analytes and interact with gold electrode.	Dialysis, dilution, or solid-phase extraction using Thermo Scientific Dionex OnGuard Ag (silver) cartridge.
Amine-containing compounds (including proteins, peptides and free amino acids).	PED active	Solid-phase extraction using Dionex OnGuard A (anion-exchange). For inline use, the Dionex AminoTrap column is used for proteins, peptides and amino acids.
Lipids	May foul column	Liquid-liquid extraction or supercritical fluid extraction.
Organic solvents	May affect analyte retention and cause diminished electrode response.	Solid-phase extraction using Dionex OnGuard RP (reverse phase).
Anionic detergents (such as SDS)	Will bind irreversibly to the column.	Solid-phase extraction using Dionex OnGuard RP.



When using pulsed electrochemical detection (PED) for detection, beware of high concentrations of electrochemically-active components (e.g., TRIS buffer, alcohols, and other hydroxylated compounds). Small amounts of organic solvents in the sample will not harm the column, although the organics may interfere with the chromatography or detection of the analytes of interest. If necessary, samples may be treated with reversed phase or ion exchange cartridges (such as the Dionex OnGuard cartridges) before analysis. However, because the Dionex CarboPac columns are extremely rugged, it is often worthwhile to analyze an aliquot of the sample directly, without any pre-column cleanup.

Sample matrices in glycoprotein analysis can be greatly simplified by performing a Western blot and selectively removing the carbohydrates from the PVDF membrane-bound proteins. Please ask for Dionex **Technical Note 30**, “Monosaccharide and Oligosaccharide Analysis of Glycoproteins Electrotransferred onto Polyvinylidene Fluoride (PVDF) membranes,” or retrieve it from our Web site at [www.thermoscientific.com](http://www.thermoscientific.com).

## 5.6 Verification of System Cleanliness

Prepare a new set of eluents as described in [Sections 5.1](#) through [5.4](#) and fill the eluent bottles. Set the eluent composition to 100% for each eluent line and draw out at least 40 mL of eluent from each eluent line.

### 5.6.1 System Background Check

This procedure is performed using the conditions of the test chromatogram. Make sure that

- A. the cell is not yet on,
- B. the pump is pumping 100 mM NaOH, 50 mM NaOAc or 200 mM KOH at 0.5 mL/min,
- C. a length of yellow tubing is installed between the injector and detector cell to generate ~1000 psi backpressure,
- D. the columns are not yet installed.

Confirm that the pH is between 12.8 and 13.4. With the pH within this range, turn on the cell using the quadruple waveform (See Table 3, Section 6.3 Disposable Electrode Manual, document number 065040) and begin monitoring the background signal from the control panel for at least 30 minutes. Confirm that the background is < 50 nC. If the background > 50 nC or the pH is out of range, see the “Troubleshooting” section at the end of this manual.

### 5.6.2 Verification of Column Cleanliness

Install the Dionex CarboPac PA100 column set only after the initial system test determines a background level within the specified range. A premature installation on a contaminated system will cause delays during the column equilibration.

The Dionex CarboPac PA100 is shipped in 100 mM NaOH + 100 mM NaOAc. Any column that is stored long-term should be stored in the same solution. Equilibrate the column set by performing two blank injections (DI water) under the test chromatogram conditions, including the column regeneration and re-equilibration steps.

Once the columns are equilibrated, inject a system suitability standard such as the column's QAR standard, to establish the performance of the column at start-up. This chromatogram can then be referred to when troubleshooting your system. Once you obtain your expected chromatography, you are ready to proceed to running your application.

Dionex recommends that the system suitability standard be run whenever you reinstall a column after long-term storage.

## 5.7 Carbohydrate Sialylated N-linked Alditol Standard

The Dionex OligoStandard, Sialylated N-Linked Alditols, P/N 043164 contains 25 nmol oligosaccharides purified from bovine fetuin. Dilute the standard prior to use, by adding a known volume of DI water (for example 1 mL). An injection of 25  $\mu$ L will then correspond to an injection of 625 pmol of the standard. Dionex recommends running this standard every time a new column is installed and subsequently anytime it becomes necessary to troubleshoot your system.

## 6. Applications

The following section provides an example of the types of applications for which the CarboPac PA100 is designed. There are several factors affecting the elution of oligosaccharides using the CarboPac PA100 column: (a) fucosylated oligosaccharides elute earlier than their non fucosylated analogs, (b) as the number of mannose residues in a high mannose oligosaccharide increases, its retention time also increases, (c) as the degree of branching increases, the retention time of the oligosaccharide increases, and (d) removal of the terminal galactose residues from a complex oligosaccharide reduces its retention time.

Separations of oligosaccharides are based on their fine structural differences such as the composition and the sequence of the oligosaccharides, linkage isomerism, degree of sialylation, and degree of branching.

## 6.1 Production Test Chromatograms

Isocratic elution of NAN-LAC on the Dionex CarboPac PA100 Analytical column has been optimized utilizing a hydroxide eluent and can be used to test the performance of the Dionex CarboPac PA100 Column. The Dionex CarboPac PA100 Analytical column should always be used with the Dionex CarboPac PA100 Guard Column; the addition of the Guard column increases elution time by ~20% when compared to the Analytical column by itself. To guarantee that all Dionex CarboPac PA100 Analytical columns meet high quality and reproducible performance specification standards, all columns undergo the following production control test. An operating temperature of 30°C is used to ensure reproducible resolution and retention.

**Figure 1**    **Dionex CarboPac PA100 Production Test Chromatograms (2mm)**

Column:                    Dionex CarboPac PA100 (2 × 250 mm)  
 Eluent:                    100mM NaOAc + 100 mM NaOH  
 Temperature:            30°C  
 Flow Rate:                0.25mL/min  
 Inj. Volume:              2.5µL  
 Detection:                Integrated Amperometry, quadruple pulse waveform  
 Working Electrode:      PTFE Gold, disposable electrode  
 Reference Electrode:    Ag/AgCl  
 Diluted Standard:        (12.0 nmol/mL / DI Water)  
     1.    Unknown  
     2.    Alpha-(2,6)-NAN-Lactose  
     3.    Alpha-(2,3)-NAN-Lactose

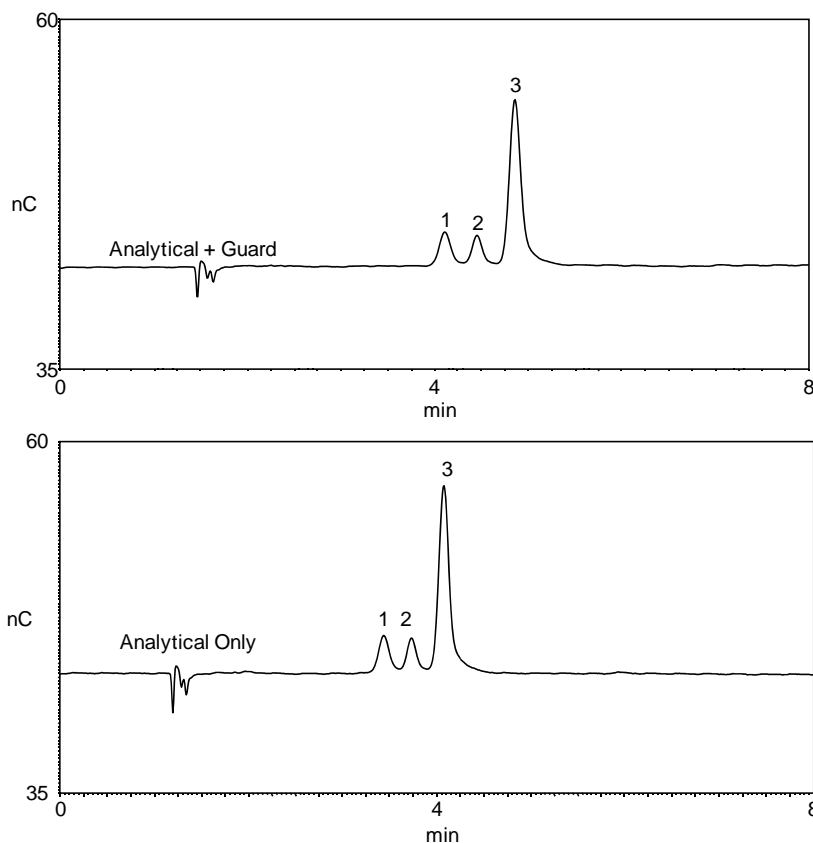
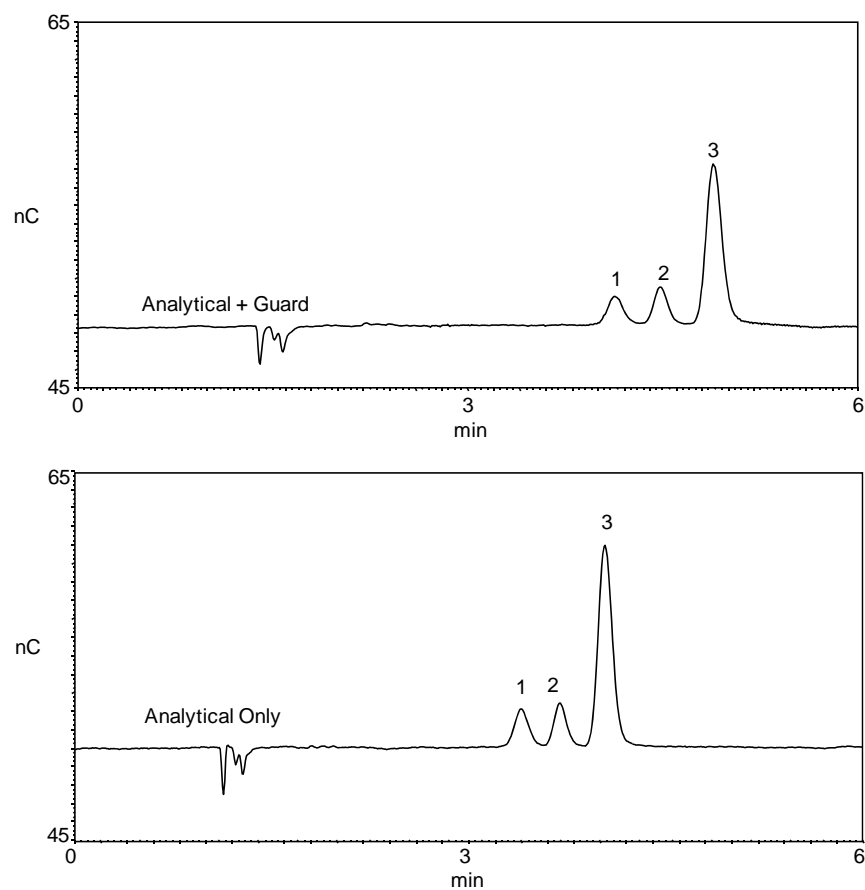


Figure 2 Dionex CarboPac PA100 Production Test Chromatograms (4mm)

Column: Dionex CarboPac PA100 (4 × 250 mm)  
Eluent: 100mM NaOAc + 100 mM NaOH  
Temperature: 30°C  
Flow Rate: 1.0 mL/min  
Inj. Volume: 10µL  
Detection: Integrated Amperometry, quadruple pulse waveform  
Working Electrode: PTFE Gold, disposable electrode  
Reference Electrode: Ag/AgCl  
Diluted Standard: (12.0 nmol/mL / DI Water)

1. Unknown
2. Alpha-(2,6)-NAN-Lactose
3. Alpha-(2,3)-NAN-Lactose



## 6.2 Separation of Neutral and Sialylated Oligosaccharides

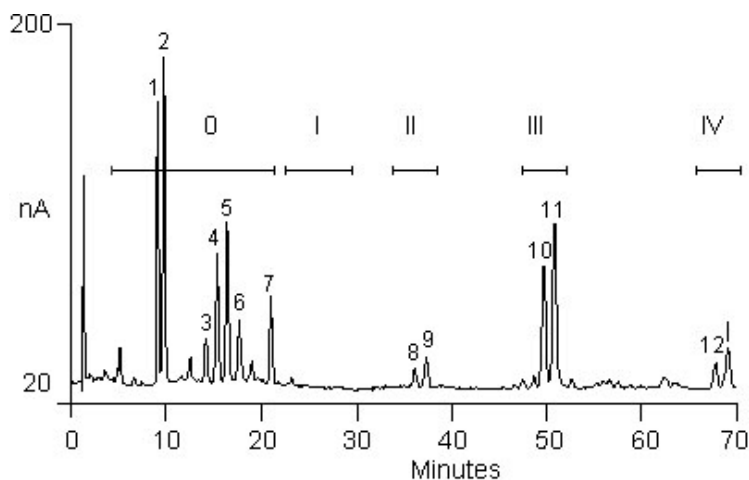
The elution of acidic sugars from the Dionex CarboPac PA100 column requires stronger eluents than those used for the elution of neutral sugars. This is usually accomplished by the addition of sodium acetate to the sodium hydroxide eluent. Sodium acetate accelerates the elution of strongly bound species without compromising selectivity or interfering with pulsed amperometric detection.

This chromatogram shows the separation of neutral and sialylated oligosaccharides in a single run, using a sodium acetate gradient in 100 mM NaOH. The neutral sugars are eluted in a group at the beginning of the profile, followed by the disialylated, the trisialylated, and finally the tetrasialylated oligosaccharides.

**Figure 3 Separation of Neutral and Sialylated Oligosaccharide Standard**

Column: Dionex CarboPac PA100 Guard and Analytical 4mm Column  
 Eluent: A – 500 mM NaOAc in 100 mM NaOH  
 B – 100 mM NaOH  
 Temperature: 30°C  
 Inj. Volume: 10µL  
 Flow Rate: 1.0 mL/min  
 Detection: Integrated Amperometry, quadruple pulse waveform  
 Working Electrode: PTFE Gold, disposable electrode  
 Reference Electrode: Ag/AgCl

Time (min)	% A	% B	Comments
0.00	0	100	Starting conditions, inject, start NaOAc gradient
110	50	50	NaOAc gradient ends
111	0	100	Back to initial conditions



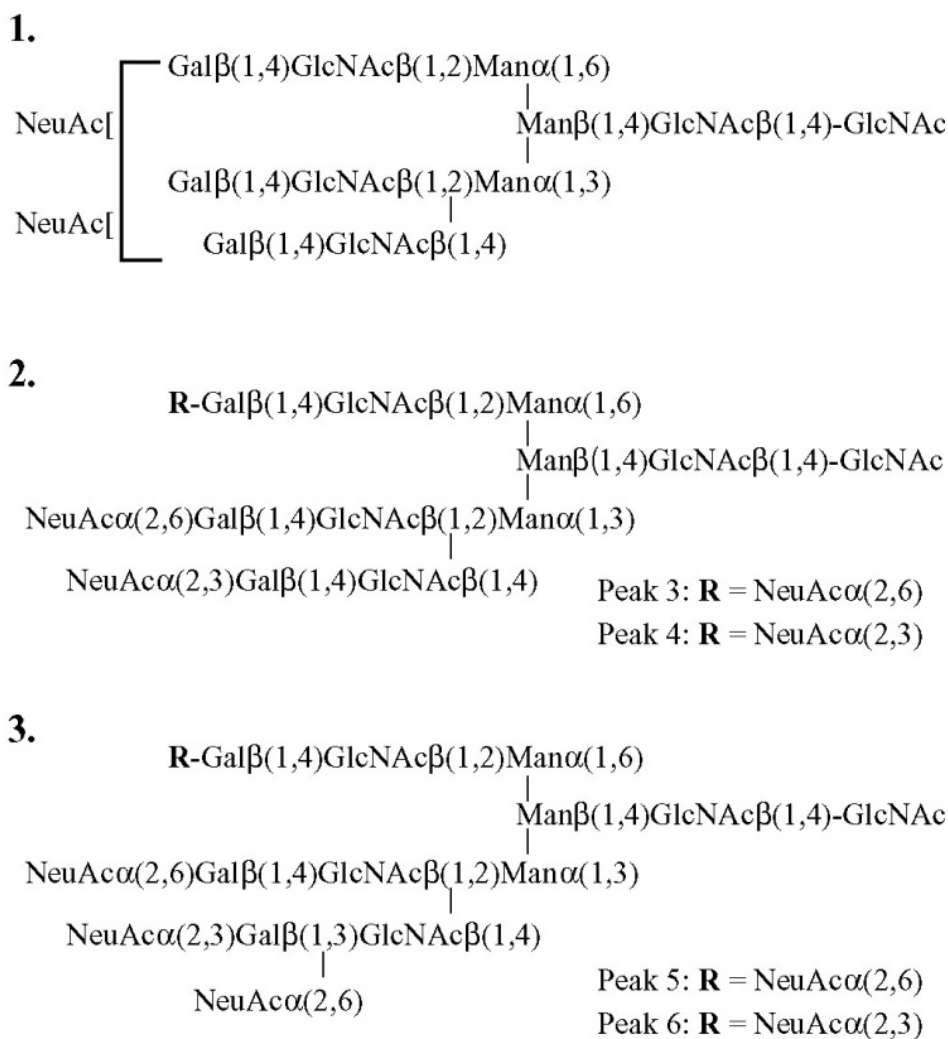
**Peaks:**

1. Fucosylated Man3GlcNAc2
2. Man3GlcNAc2
3. Asialo, agalacto bi, core fuc
4. Asialo, agalacto bi
5. Asialo bi, core fuc
6. Asialo bi
7. Man9GlcNAc2
- 8,9. Disialylated tri (reduced)
- 10,11. Trisialylated tri (reduced)
- 12,13. Tetrasialylated tri (reduced)

### 6.3 Separation of Fetuin N-linked Oligosaccharide Alditols

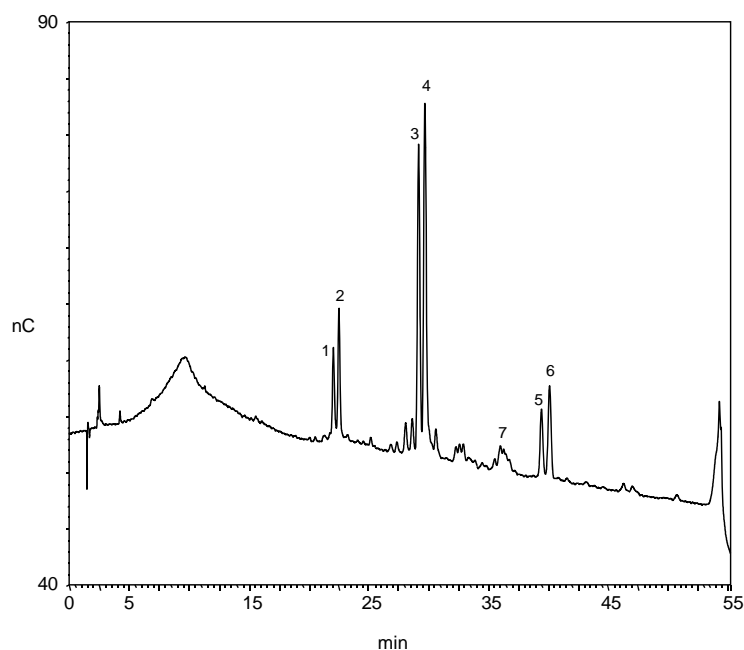
The separation of bovine fetuin oligosaccharides by HPAE-PAD is shown below. The structures for Peaks 1-6 are shown in Figure 4 below. Peak 7 is a trisialylated triantennary complex oligosaccharide. Under alkaline conditions, the technique resolves these species not only by sialic acid content, but also according to the combination of (2,3)- and (2,6)-linked sialic acids within each charge class. Oligosaccharides with the greatest proportion of (2,6)- to (2,3)-linked sialic acids are the least-retained. The neutral component of the oligosaccharides also influence separation. Of the oligosaccharides studied, those containing a Gal $\beta$  (1,3) GlcNAc sequence are retained more strongly than those with Gal $\beta$  (1,4) GlcNAc.

**Figure 4 Major Carbohydrate Structures of Bovine Fetuin**



**Figure 5 Separation of Fetuin N-linked Oligosaccharide Alditols**

Column: Dionex CarboPac PA100 Guard and Analytical 2mm Column  
 Eluent: A – 100 mM NaOH  
           B – 500mM NaOAc + 100 mM NaOH  
 Temperature: 30°C  
 Flow Rate: 0.25 mL/min  
 Inj. Volume: 2.5µL  
 Detection: Integrated Amperometry, quadruple pulse waveform  
 Working Electrode: PTFE Gold, disposable electrode  
 Reference Electrode: Ag/AgCl  
 Sample Preparation: Fetuin N-linked oligosaccharide aditol (Dionex) diluted with 500µL of DI Water

**Peaks:**

1. Disialylated, triantennary
2. Disialylated, triantennary
3. Trisialylated, triantennary
4. Trisialylated, triantennary
5. Tetrasialylated, triantennary
6. Tetrasialylated, triantennary
7. Trisialylated, triantennary

**Gradient Program**

Time (min)	%A	%B	Comments
-15.00	99	1	Equilibration
0.0	99	1	Isocratic at 5mM NaOAc
0.2	99	1	Start NaOAc gradient
10.0	90	10	Change slope of NaOAc gradient
50.0	55	45	End NaOAc gradient. Start column wash
50.1	0	100	Column wash
55.0	0	100	End column wash. Back to initial conditions
55.1	99	1	Start re-equilibration under initial conditions
70.0	99	1	End re-equilibration

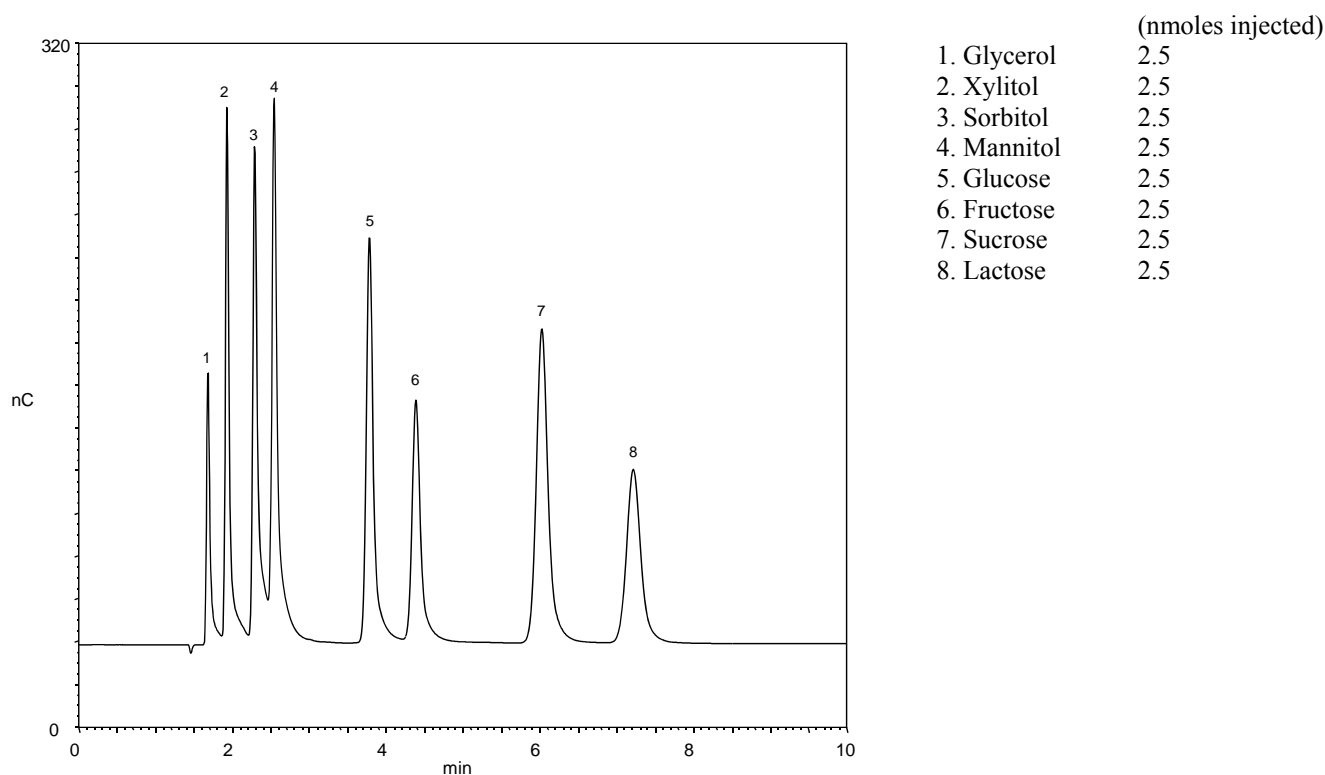


## 6.4 Isocratic Separation of Food Sugars and Food Alcohols

This application uses an isocratic elution to separate alditols, monosaccharides and disaccharides commonly found in food samples, in a single run. Alditol retention is only slightly affected by sodium hydroxide concentration between 18 mM and 50 mM. This allows for optimization of monosaccharide and disaccharide separations by increasing the hydroxide concentration while still maintaining adequate alditol retention. Use of the Dionex CarboPac BorateTrap column greatly improves the peak shape of the alditols.

**Figure 6 Separation of Food Sugars and Food Alcohols**

Column: Dionex CarboPac PA100 Guard and Analytical 2mm Column  
 Eluent: 150mM NaOH  
 Temperature: 30°C  
 Flow Rate: 0.25 mL/min  
 Inj. Volume: 2.5µL  
 Detection: Integrated Amperometry, quadruple pulse waveform  
 Working Electrode: PTFE Gold, disposable electrode  
 Reference Electrode: Ag/AgCl



## 6.5 Separation of Apple and Orange Juice Sugars Using Dionex CarboPac PA100

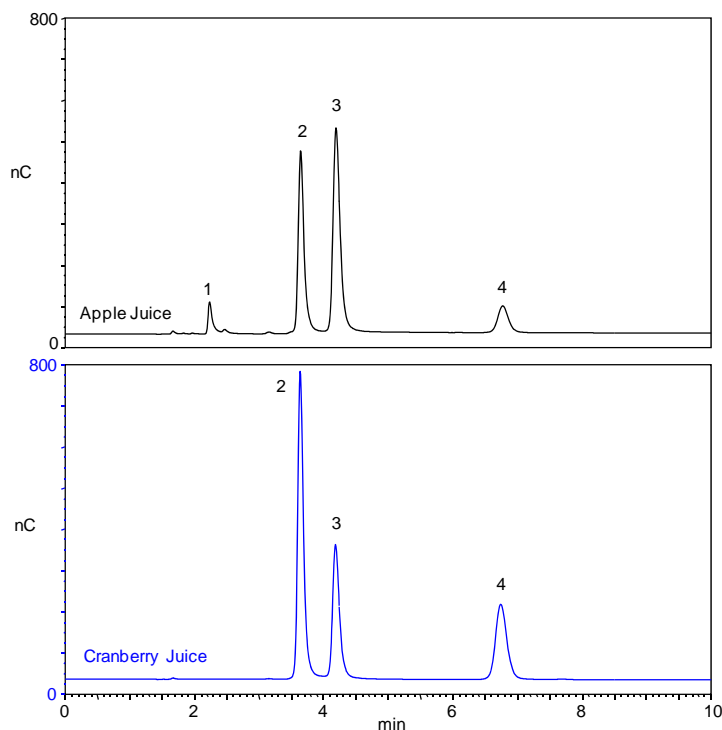
Fruit juice adulteration is an economic and regulatory problem and a matter of worldwide concern. Fruit juices are targets for adulteration because of their high cost relative to sweeteners and other ingredients. The most common forms of adulteration are dilution and blending of inexpensive and synthetically produced juices into expensive ones, and addition of pulp wash. A common adulterant used to mask the effects of dilution and addition of other adulterants is beet medium invert sugar (BMIS). BMIS is partially inverted sucrose with a glucose:fructose:sucrose ratio of 1:1:2 which closely matches the ratio found in orange juice, and it is difficult to detect.

The Figures 7 and 8 below compare the separation of juice sugars using 2mm and 4mm CPPA100 columns. Notice 2.5µL (1/4th of the 4mm) sample was injected on the 2mm column as compared to 10µL for the 4mm column, yet the peak signal is very close to the 4mm column.

The basis for this method is that invert syrups contain traces of oligosaccharides not generally found in natural juices, so the selectivity of anion exchange chromatography for oligosaccharides, and the sensitivity and specificity of PAD is uniquely suited to this analysis. The chromatograms below do not contain any peaks due to possible oligosaccharide content. The analyzed juice samples are thus unaffected by adulteration. The major peaks at the start of the chromatogram are glucose, fructose, and sucrose which are natural components of fruit juices.

**Figure 7 Separation of Juice Sugars Using Dionex CarboPac PA100 (2 mm)**

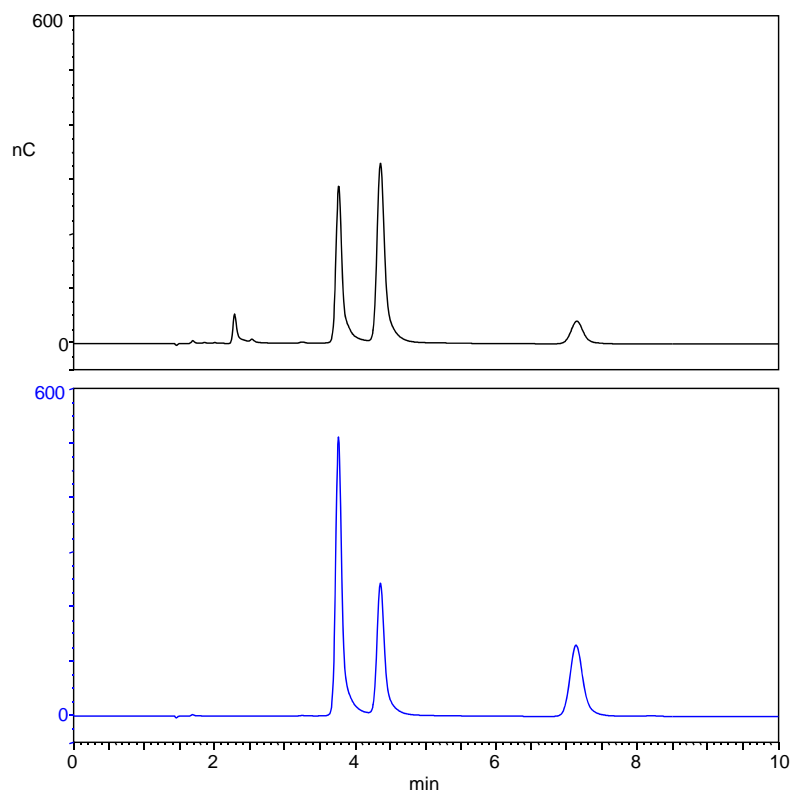
Column: Dionex CarboPac PA100 Guard and Analytical 2mm Column  
 Eluent: 150 mM NaOH  
 Temperature: 30°C  
 Flow Rate: 0.25 mL/min  
 Inj. Volume: 2.5µL  
 Detection: Integrated Amperometry, quadruple pulse waveform  
 Working Electrode: PTFE Gold, disposable electrode  
 Reference Electrode: Ag/AgCl  
 Sample: Juice samples (1:1000 dilution) / DI Water



Peaks: (nmoles injected)	Apple Cranberry	
	Apple	Cranberry
1. Sorbitol	0.05	--
2. Glucose	0.37	0.62
3. Fructose	0.72	0.48
4. Sucrose	0.12	0.41

**Figure 8** Separation of Juice Sugars Using Dionex CarboPac PA100 (4 mm)

Column: Dionex CarboPac PA100 Guard and Analytical 4mm Column  
 Eluent: 150 mM NaOH  
 Temperature: 30°C  
 Flow Rate: 1.0 mL/min  
 Inj. Volume: 10µL  
 Detection: Integrated Amperometry, quadruple pulse waveform  
 Working Electrode: PTFE Gold, disposable electrode  
 Reference Electrode: Ag/AgCl  
 Sample : Juice samples (1:1000 dilution) / DI Water



**Peaks:** (nmoles injected)

	<u>Apple</u>	<u>Cranberry</u>
1. Sorbitol	0.25	--
2. Glucose	1.56	2.68
3. Fructose	3.24	2.26
4. Sucrose	0.53	1.59

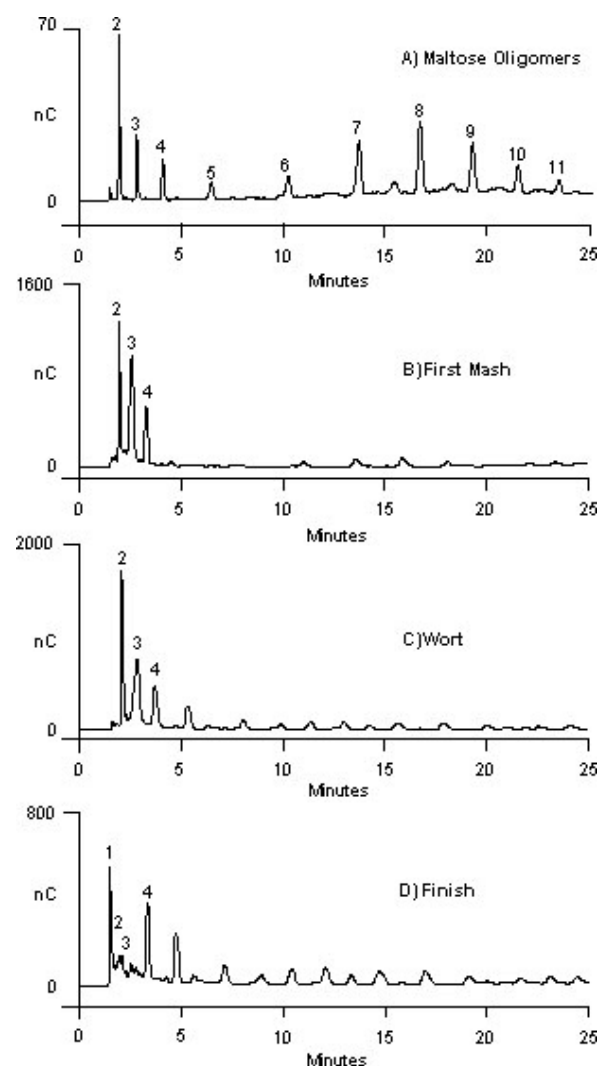
## 6.6 Oligosaccharide Profiling During Beer Production

Determining the levels of fermentable and non-fermentable sugars at every stage of beer production is important because fermentable sugars determine the final alcohol content, and non-fermentable sugars contribute to the flavor and ‘body’ of the final product. Sugars, sugar alcohols, alcohols, and glycols can be rapidly determined with high resolution at all phases of beer, wine or cider production.

A separation of maltose oligomers up to DP10 (degree of polymerization) with baseline resolution is shown in Panel A. Sugar and oligosaccharide profiles at various stages of the brewing process are shown in Panels B, C and D. All samples were diluted 1:10.

**Figure 9 Carbohydrate Profiles at Various Stages of Beer Production**

Analytical: Dionex CarboPac PA100 Guard and Analytical 4mm Column  
 Eluent A: 100 mM NaOH  
 Eluent B: 500 mM NaOAc in 100 mM NaOH  
 Flow rate: 1.0 mL/min  
 Sample Volume: 10 µL  
 Detection: Integrated Amperometry, quadruple waveform  
 Electrode: Gold



Peaks:

1. Ethanol
2. Glucose
3. Maltose
4. Maltotriose
5. Maltotetraose
6. Maltopentaose
7. Maltohexaose
8. Maltoheptaose
9. Maltooctaose
10. Maltononaose
11. Maltodecaose

Peaks:

1. Ethanol
2. Glucose
3. Maltose
4. Maltotriose

Peaks:

1. Ethanol
2. Glucose
3. Maltose
4. Maltotriose

Peaks:

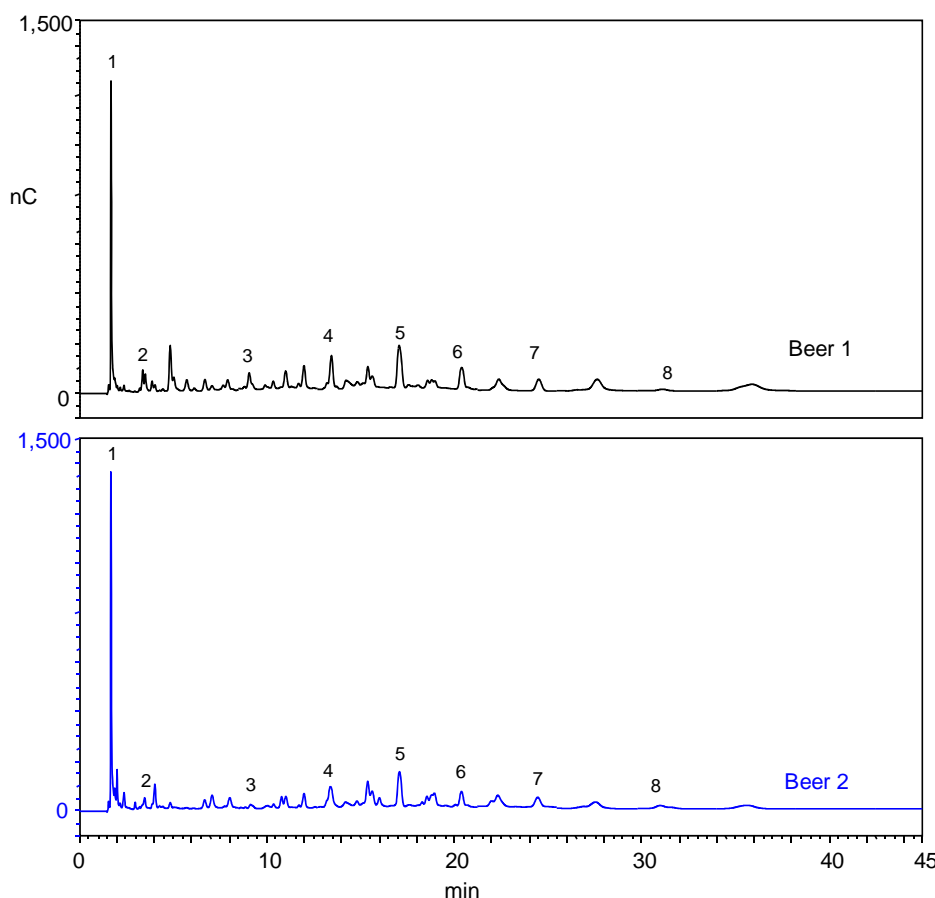
1. Ethanol
2. Glucose
3. Maltose
4. Maltotriose

## 6.7 Analysis of Oligosaccharides in Beer

The Figure 10 shows the analysis of two different beer samples using CarboPac PA100 2mm column. Notice that ethanol and glycerol co elute on the CarboPac PA100 column.

**Figure 10 Analysis of Oligosaccharides in Beer**

Column: Dionex CarboPac PA100 (2 × 250 mm) Analytical and Guard Column  
 Eluent: A – DI Water  
 B – 200 mM NaOH  
 C – 0.5 M NaOAc  
 Temperature: 30°C  
 Flow Rate: 0.25 mL/min  
 Inj. Volume: 2.5 µL  
 Detection: Integrated Amperometry, quadruple pulse waveform  
 Working Electrode: PTFE Gold, disposable electrode  
 Reference Electrode: Ag/AgCl  
 Sample: Beer samples (1:10 dilution) / DI Water



**Gradient Program**

Time (min)	%A	%B	%C
-7.00	24	75	1
0.00	24	75	1
15.00	5	75	20
45.00	5	75	20

Note Equilibration Time:  
7 min

**Peaks:**

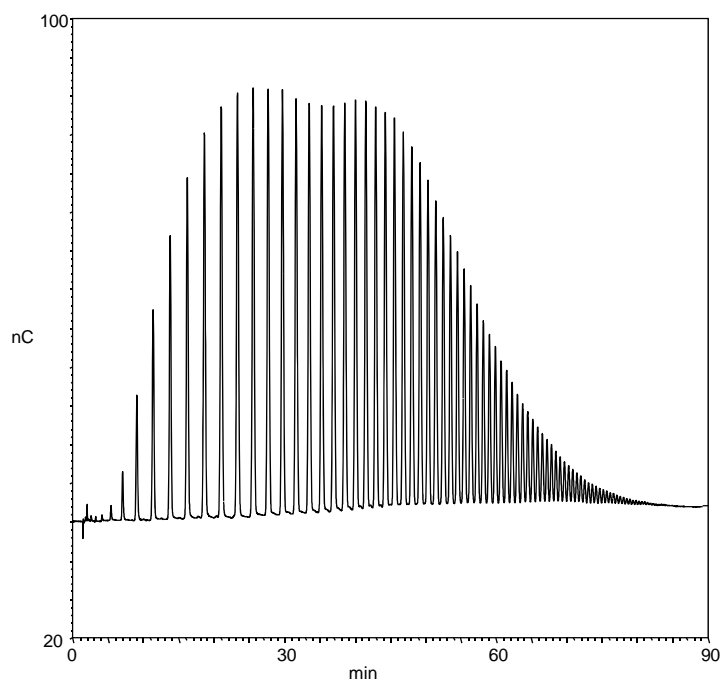
1. Ethanol+Glycerol
2. Glucose
3. Maltose
4. Maltotriose
5. Malrotetraose
6. Maltopentaose
7. Maltohexaose
8. Maltoheptaose

## 6.8 Gradient Separation of Linear Polysaccharides

Commercial inulin products have degrees of polymerization (DP) which have been tailored for a particular end use. Therefore, it is important to determine the chain length distribution during product development, production and for quality control of the end product. The following chromatogram shows the excellent resolution, which can be achieved for DP values up to 50 for inulin derived from chicory.

**Figure 11** Gradient Separation of Jerusalem Artichoke Inulin Polymers Using Dionex CarboPac PA100 Column

Column: Dionex CarboPac PA100 Guard and Analytical 2mm Column  
 Eluent: A – DI Water  
           B – 200mM NaOH  
           C – 100mM NaOH, 500mM NaOAc  
 Temperature: 30°C  
 Flow Rate: 0.25 mL/min  
 Inj. Volume: 2.5µL  
 Detection: Integrated Amperometry, quadruple pulse waveform  
 Working Electrode: PTFE Gold, disposable electrode  
 Reference Electrode: Ag/AgCl  
 Sample : Inulin from Jerusalem Artichoke (1mM dilution) / DI Water



Gradient Program

Time (min)	%A	%B	%C
-7.0	35	35	30
0.0	35	35	30
90.0	5	5	90

Note Equilibration Time:  
7 min

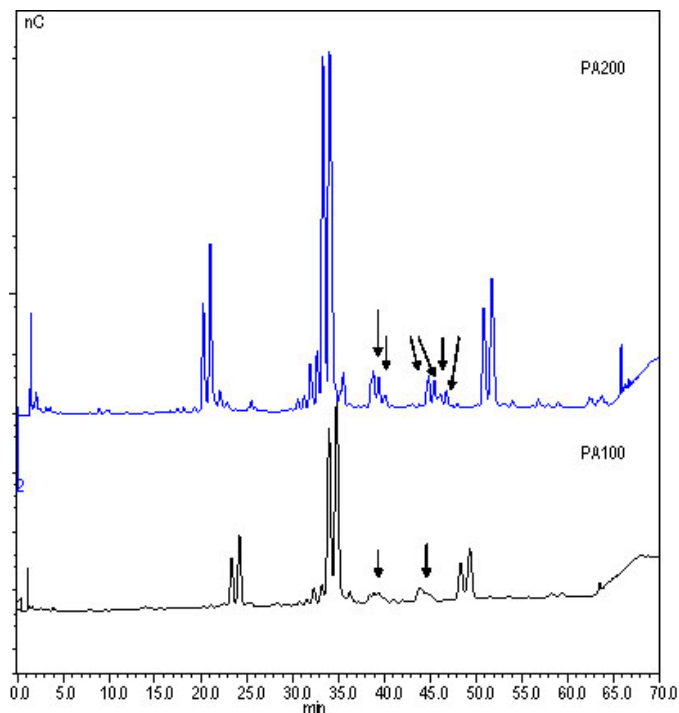
## 6.9 Fetuin Oligosaccharide Alditol Profiling

The high resolution of the Dionex CarboPac PA200 is exemplified in the following example. Dionex CarboPac columns separate mono-, oligo- and polysaccharides on the basis of fine structural differences in branching, linkage isomerism, anomericity and sialylation. In the fetuin oligosaccharide alditol standard shown below, peaks are separated according to branching, sialylation and linkage isomerism. The disialylated biantennary peaks are eluted before the trisialylated triantennary peaks which are eluted before the tetrasialylated tetraantennary peaks. In addition, within each grouping, the 2-6 isomer is eluted before, and well resolved from, the 2-3 isomer.

The Dionex OligoStandard, Sialylated N-Linked Alditols, P/N 043164 contains 25 nmol oligosaccharides purified from bovine fetuin. Dilute the standard prior to use, by adding a known volume of DI water (for example 1 mL; a 25  $\mu$ L injection then corresponds to 625 pmol of oligosaccharide). Thermo Scientific recommends running this standard every time a new column is installed and subsequently anytime it becomes necessary to troubleshoot your system.

**Figure 12** Fetuin Oligosaccharide Profiles: Dionex CarboPac PA200 vs. Dionex CarboPac PA100. The resolution of some oligosaccharides is improved on the PA200 column (see arrows)

Columns: Dionex CarboPac PA200 (3  $\times$  250 mm) and Dionex CarboPac PA100 (4  $\times$  250 mm)  
 Gradient: Dionex CarboPac PA200:  
 20–150mM NaOAc in 100 mM NaOH over 1 h  
 Dionex CarboPac PA100:  
 20–200 mM NaOAc in 100 mM NaOH over 1 h  
 Flow Rate:  
 Dionex CarboPac PA200: 0.5 mL/min  
 Dionex CarboPac PA100: 1mL/min  
 Detection: Pulsed amperometry, QP waveform, gold electrode  
 Samples: Fetuin oligosaccharide alditol standard

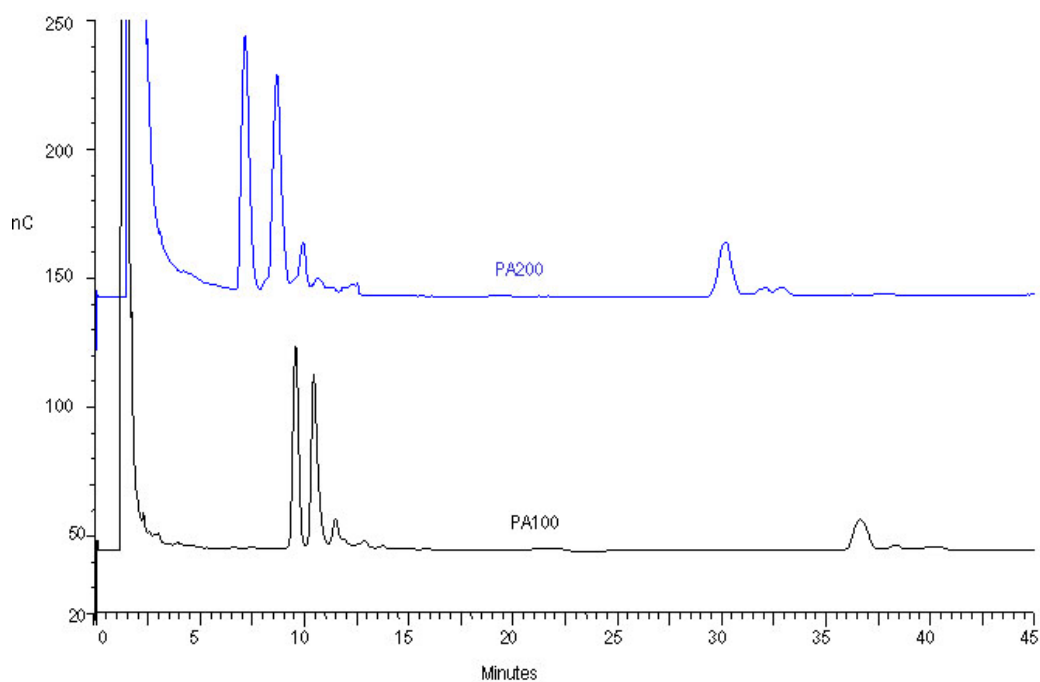


## 6.10 Antibodies N-Glycan Separations

Today, monoclonal antibodies comprise the cornerstone of numerous medical and scientific procedures and thus are produced routinely for both medical and research applications. In the medical world, they are used as diagnostics, to detect cancers or infection by certain bacteria or viruses; as vaccines, to boost the body's immune response; and as therapeutics, to target foreign bacteria, viruses or cancerous cells.

**Figure 13** Dionex CarboPac PA200 vs. Dionex CarboPac PA100 Separation of N-Glycan Released from Monoclonal Antibody N-linked Oligosaccharides

Columns: Dionex CarboPac PA200 (3 × 250 mm) and Dionex CarboPac PA100 (4 × 250 mm)  
Gradient: 10 mM or 20 mM – 150 mM NaOAc in 100 mM NaOH over 60 min  
Flow Rate: Dionex CarboPac PA200: 0.5 mL/min  
Dionex CarboPac PA100: 1 mL/min  
Detection: Pulsed amperometry, QP waveform, gold electrode  
Sample: 48 h PNGase F digest of monoclonal antibody (100 µg)



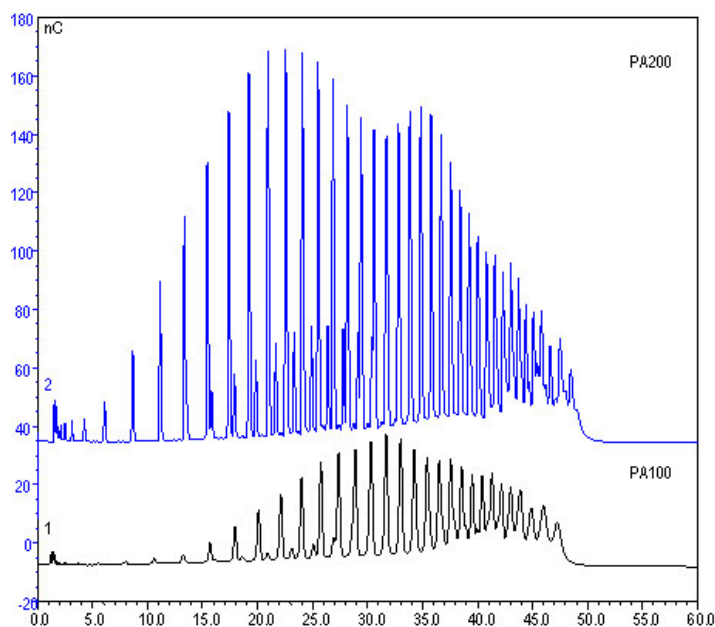


## 6.11 Profiling of Complex Plant Carbohydrates: Profiling of Inulins

Inulins and fructooligosaccharides (FOS) are increasingly being used as functional food ingredients. Chain length distribution profiles of commercial products such as those derived from inulin can be determined by using HPAE-PAD with gradient elution. By adjusting the initial gradient profile, smaller oligofructose chains can be distinguished from the inulin chains and separations up to and exceeding DP80 are possible. Quantification of individual inulin oligomers requires knowledge of the PAD response factors.

**Figure 14 Inulin Profiles: Dionex CarboPac PA200 vs. Dionex CarboPac PA100**

Columns:	Dionex CarboPac PA200 (3 × 250 mm) and Dionex CarboPac PA100 (4 × 250 mm)
Gradient:	120 – 320 mM NaOAc in 100 mM NaOH over 40 min
Flow Rate:	Dionex CarboPac PA200: 0.5 mL/min Dionex CarboPac PA100: 1 mL/min
Detection:	Pulsed amperometry, quadruple potential waveform, gold electrode
Samples:	Inulin from chicory (Sigma)

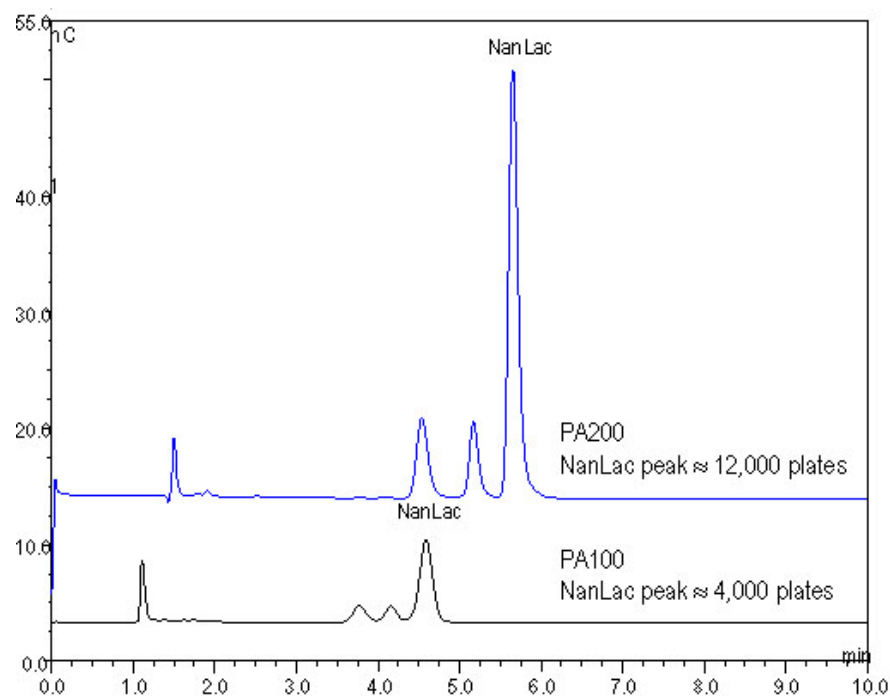


## 6.12 Sialyl Lactose (N-Acetylneuraminy-Lactose)

The following chromatography is used as the quality assurance test for the CarboPac PA200 column. Every Dionex CarboPac PA200 column is tested with respect to sialyl lactose peak efficiency. Chromatography of sialyl lactose on the Dionex CarboPac PA200 column exhibits greater than a twofold increase in peak efficiencies compared to chromatography of sialyl lactose on the Dionex CarboPac PA100 column.

**Figure 15** Chromatography of NanLac on the Dionex CarboPac PA200 exhibits > 2x the peak efficiencies compared to the Dionex CarboPac PA100

Columns:	Dionex CarboPac PA200 (3 × 250 mm) or Dionex CarboPac PA100 (4 × 250 mm)
Gradient:	
	Dionex CarboPac PA200: 100 mM NaOH, 50 mM NaOAc isocratic
	Dionex CarboPac PA100: 100 mM NaOH, 100 mM NaOAc isocratic
Flow Rate:	
	Dionex CarboPac PA200: 0.5 mL/min
	Dionex CarboPac PA100: 1 mL/min
Detection:	Pulsed amperometry, quadruple potential waveform, gold electrode
Sample:	NanLac Std

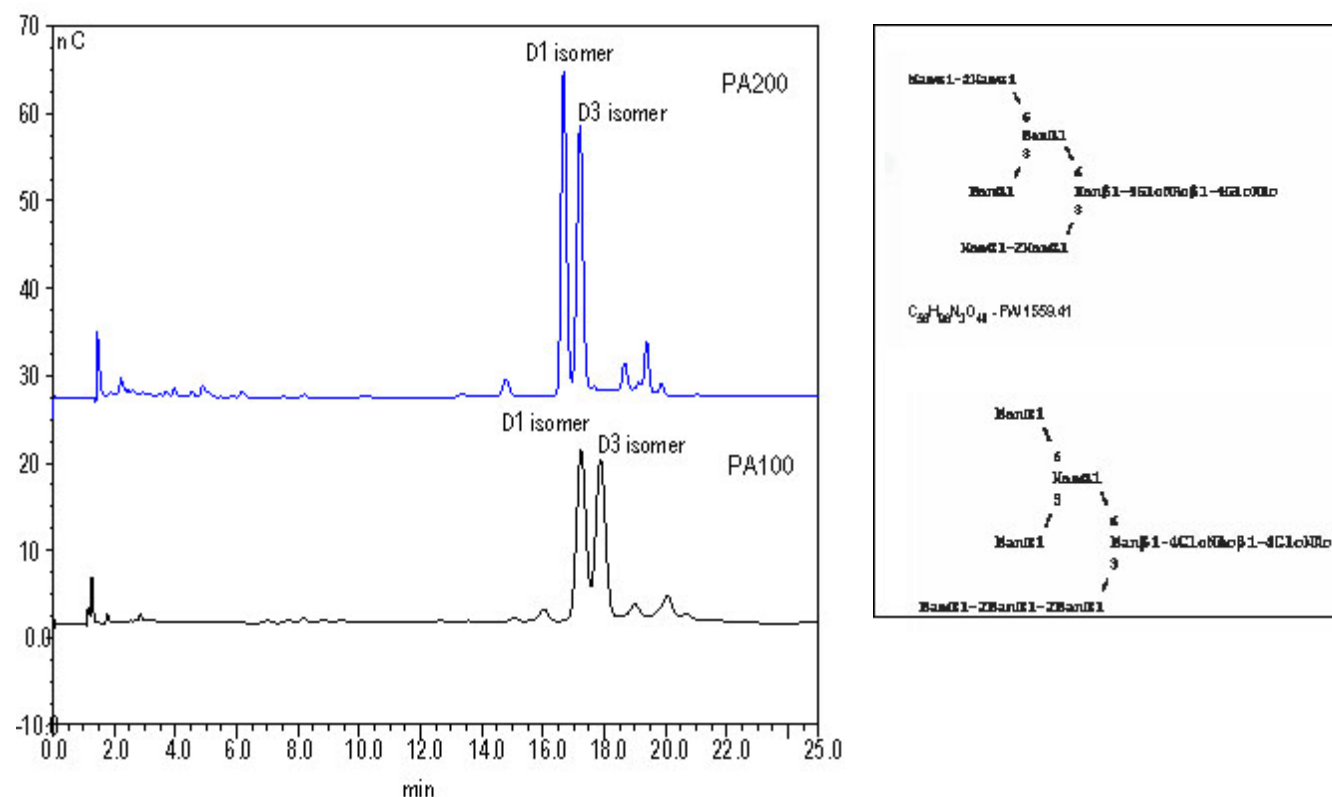


### 6.13 Separation of Mannose 7 D1, D3 Isomers

High mannose isomers that frequently would not be distinguishable by other techniques are well separated on both the CarboPac PA200 and PA100 columns. The separation of these isomers is better on the CarboPac PA200 column.

**Figure 16** Dionex CarboPac PA200 vs. Dionex CarboPac PA100 Separation of Mannose 7 D1, D3 Isomers

Columns: Dionex CarboPac PA200 (3 × 250 mm) or Dionex CarboPac PA100 (4 × 250 mm)  
 Gradient: 0–200 mM NaOAc in 100 mM NaOH over 110 minutes  
 Flow Rate: Dionex CarboPac PA200: 0.5 mL/min  
 Dionex CarboPac PA100: 1 mL/min  
 Detection: Pulsed amperometry, quadruple potential waveform, gold electrode  
 Sample: Mannose 7 isomers (Dextra Labs):



## 6.14 Separation of Neutral N-Linked Oligosaccharides: Complex Neutral Structures

N-linked glycoproteins may exhibit micro-heterogeneity with respect to the N-linked oligosaccharides that may be present. Neutral N-linked oligosaccharides released from glycoproteins could be of the complex or high mannose variety. A comparison of chromatography of neutral and high mannose oligosaccharides on the CarboPac PA200 column compared to the CarboPac PA100 column shows that the separation window is wider for the CarboPac PA200 column. Early eluting oligosaccharides on the CarboPac PA100 column tend to elute earlier on the CarboPac PA200 column.

**Figure 17** Dionex CarboPac PA200 vs. Dionex CarboPac PA100 Separation of Neutral N-linked Oligosaccharides (complex neutral structures)

Columns: Dionex CarboPac PA200 (3 × 250 mm) or Dionex CarboPac PA100 (4 × 250 mm)

Gradient: 0–200 mM NaOAc in 100 mM NaOH over 110 minutes

Flow Rate:

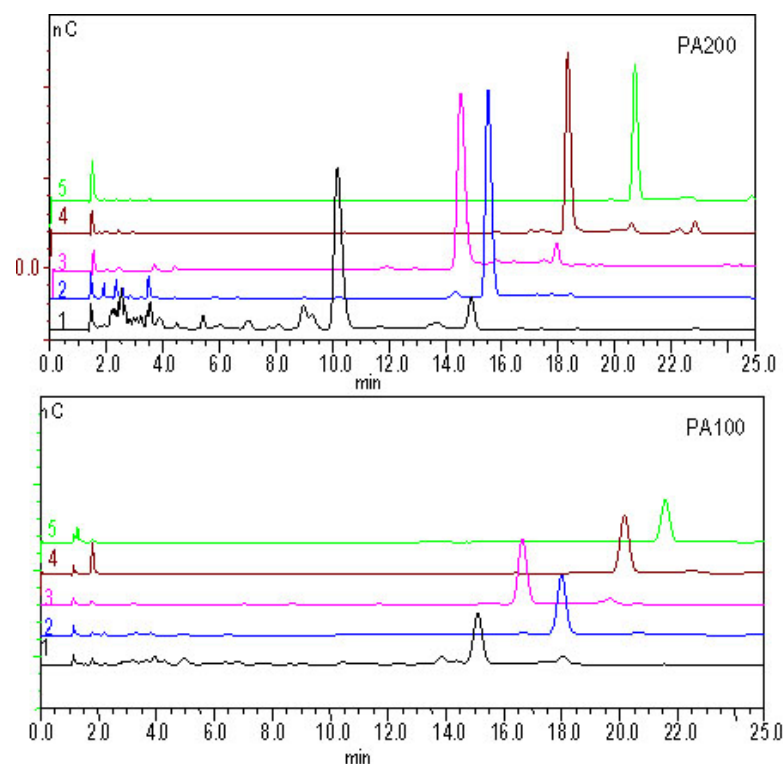
Dionex CarboPac PA200: 0.5 mL/min

Dionex CarboPac PA100: 1 mL/min

Detection: Pulsed amperometry, quadruple potential waveform, gold electrode

Samples:

1. Asialo agalacto biantennary standard
2. Asialo galacto biantennary standard
3. Asialo galacto fuco biantennary standard
4. Asialo triantennary standard
5. Asialo tetraantennary standard (All standards are from Dextra Labs)



## 6.15 High Mannose Structures

High mannose structures are well separated on both the CarboPac PA200 and PA100 columns. However, the separation and sensitivity is better on the CarboPac PA200 column.

**Figure 18** Dionex CarboPac PA200 vs. Dionex CarboPac PA100

Columns: Dionex CarboPac PA200 (3 × 250 mm) or Dionex CarboPac PA100 (4 × 250 mm)

Gradient: 0–200 mM NaOAc in 100 mM NaOH over 110 minutes

Flow Rate:

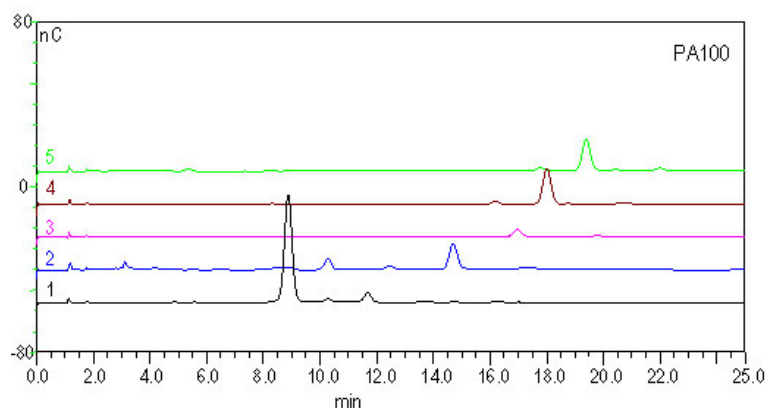
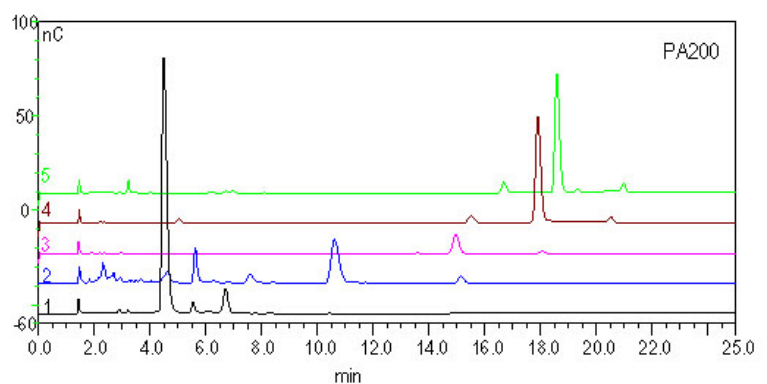
Dionex CarboPac PA200: 0.5 mL/min

Dionex CarboPac PA100: 1 mL/min

Detection: Pulsed amperometry, quadruple potential waveform, gold electrode

Samples:

1. mann 3 standard
2. mann 5 standard
3. mann 6 standard
4. mann 7 standard
5. mann 8 standard (All standards are from Dextra Labs)



## 7. Troubleshooting Guide

The purpose of the Troubleshooting Guide is to help you solve operating problems that may arise while using Dionex CarboPac columns. For more information on problems that originate with the Ion Chromatograph (IC), refer to the Troubleshooting Guide in the appropriate operator's manual. Remember that some of the problems may be related to parts of your experimental protocol (sample contamination, imprecision during sample transfer, problems during peptide or protein hydrolysis, etc.). The following text should help you to locate and eliminate problems traceable to the carbohydrate hardware and chemistries. It also provides a selection of cleanup and reconditioning procedures that have been found effective by many users.

***For assistance, contact Technical Support for Dionex Products. In the U.S., call 1-800-346-6390. Outside the U.S., call the nearest Thermo Fisher Scientific office.***

### 7.1 High Back Pressure

#### 7.1.1 Finding the Source of High System Pressure

Total system pressure for the Dionex CarboPac Guard plus the Dionex CarboPac Analytical Column when using the test chromatogram conditions should be close to the pressure listed in the QAR. If CarboPac guard and analytical column is being installed on the system, column pressure will increase by approximately 20% as compare to pressure listed in the QAR for each column type. If the total system pressure is much higher than expected, it is advisable to determine the cause of the high system pressure.

- A. Make sure that the pump is set to the correct eluent flow rate. Higher than recommended eluent flow rates will cause higher pressure. Measure the pump flow rate if necessary with an analytical balance.
- B. Determine which part of the system is causing the high pressure. High pressure could be due to a plugged tubing or tubing with collapsed or pinched walls, an injection valve with a clogged port, a column with particulates clogging the bed support, a clogged High-Pressure In-Line Filter, or the detector cell.

To determine which part of the chromatographic system is causing the problem, disconnect the pump eluent line from the injection valve and turn the pump on. Watch the pressure; it should not exceed 200 psi. (unless a backpressure coil has been installed between the pump outlet and the injection valve in which case, first disconnect the eluent line from the pump to the backpressure coil). Continue adding system components (backpressure coil (if present), injection valve, column(s), suppressor and detector) one by one, while monitoring the system pressure. The pressure should increase by the sum of the measured pressures of the individual guard and analytical columns (see product QAR) when the CarboPac Guard and Analytical columns are connected.

#### 7.1.2 Replacing Column Bed Support Assemblies for 2 mm and 4 mm columns

If the column inlet bed support is determined to be the cause of the high back pressure, it should be replaced. To change the inlet bed support assembly, refer to the following instructions, using one of the two spare inlet bed support assemblies included in the Ship Kit.

- A. Disconnect the column from the system.
- B. Carefully unscrew the inlet (top) column fitting. Use two open-end wrenches.
- C. Remove the bed support. Turn the end fitting over and tap it against a benchtop or other hard, flat surface to remove the bed support and seal assembly. If the bed support must be pried out of the end fitting, use a sharp pointed object such as a pair of tweezers, but be careful that you do not scratch the walls of the end fitting. Discard the old bed support assembly.

- D. Place a new bed support assembly (provided with each analytical column) into the end fitting. Make sure that the end of the column tube is clean and free of any particulate matter so that it will properly seal against the bed support assembly. Drop the bed support assembly into the end fitting, making sure that the bed support assembly is centered at the bottom of the end fitting. Wrap the end fitting gently on a hard surface to reorient the bed support assembly as necessary in order to properly situate the bed support assembly in the end fitting.



*If the column tube end is not clean when inserted into the end fitting, particulate matter may obstruct a proper seal between the end of the column tube and the bed support assembly. If this is the case, additional tightening may not seal the column but instead damage the column tube or the end fitting.*

- E. While holding the column in an inverted configuration, tighten the end fitting back onto the column. Tighten it finger-tight, then an additional 1/4 turn (25 in-lb). Tighten further only if leaks are observed.
- F. Reconnect the column to the system and resume operation.

### 7.1.3 Filter Eluent

Eluents containing particulate material or bacteria may clog the column inlet bed support. Filter eluents through a 0.45 µm Nylon or PTFE filter.

### 7.1.4 Filter Samples

Samples containing particulate material may clog the column inlet bed support. Filter samples through a 0.45 µm filter prior to injection.

## 7.2 High Background

While it may be possible to obtain reasonable performance even with elevated levels of detection background according to some requirements, high background frequently brings about an increased size of gradient artifacts and can be accompanied by a presence of ghost peaks. Detection sensitivity may also change suddenly when the detection background is too high. A background >50 nC with 10 mM sodium hydroxide at 0.5 mL/min and 30°C using the quadruple waveform indicates one of the following possibilities:

- A. Incorrect detection parameters.  
Verify that Ag/AgCl is specified as a reference electrode. Check all values of waveform in program against those in the Disposable Electrode Manual. If the pH reading with 10 mM NaOH or KOH is above 13.2 replace the reference electrode.
- B. Compromised working electrode surface.  
Briefly install a new working electrode and check the background as above. If the reading remains > 50 nC, remove the new electrode within 30 minutes and continue testing for column or system contamination. Otherwise continue with your work with the new electrode installed.
- C. Column contamination: Remove the column set from the system first and replace it with a length of yellow PEEK tubing, generating a pressure drop between 1000 and 2000 psi. If the background reading improves after the column is removed from the system, go to [Appendix A, “CarboPac PA100 Column Care”](#).
- D. Water contamination: Prepare eluents using a fresh ultra pure water from another source. If the background is reduced, investigate the source of contamination in the original source of water.

- E. System Contamination: If the background remains high even with fresh water and without the column, carry out the 2 M sodium hydroxide rinse. In a properly working system, the electrochemical detection (ED) background for the Dionex CarboPac PA100 QAR eluent is 50-70nC. If the background is much higher, determine the cause of high background.

### 7.2.1 Preparation of Eluents

- A. Make sure that the eluents are made correctly.
- B. Make sure that the eluents are made from chemicals with the recommended purity.
- C. Make sure that the deionized water used to prepare the reagents has a specific resistance of 18.2 megohm-cm or greater.

### 7.2.2 CR-ATC Column

- A. When using eluent generator (EGC-KOH) to generate eluent, install a Dionex CR-ATC Anion Trap Column.
- B. If the background is elevated due to contamination of the Dionex CR-ATC, please refer to Sections 5.3 and 6, in the Dionex CR-ATC Product Manual (Document No. 031910) for corrective action.

### 7.2.3 A Contaminated Guard or Analytical Column

- A. Remove the columns from the system.
- B. Install a back pressure coil that generates approximately 2000 psi and continue to pump eluent. If the background decreases, the column(s) is (are) the cause of the high background.
- C. To eliminate downtime, clean or replace the analytical column at the first sign of column performance degradation. Clean the column as instructed in, [“Appendix A, CarboPac PA100 Column Care”](#).

## 7.3 Poor Resolution

One of the unique features of Dionex CarboPac columns is the fast equilibration time in gradient applications from the last eluent (high ionic strength) to the first eluent (low ionic strength). The actual equilibration time depends on the ratio of the strongest eluent concentration to the weakest eluent concentration and application flow rate. Typically equilibration times range from 10 to 15 minutes at 1mL/min.

If increased separation is needed for early eluting peaks, reduce the initial eluent concentration.

Due to different system configurations, the gradient profile may not match the gradient shown in example applications in the product column manual. Gradient conditions can be adjusted to improve resolution or to adjust retention times either by changing the gradient timing or by changing the initial and/or final eluent concentration.

- A. Keep the eluent concentrations constant and adjust the gradient time. This is the simplest way to compensate for total system differences if resolution is the problem.
- B. Change the initial and/or final eluent concentration and adjust the gradient time. This approach requires more time to develop and more knowledge in methods development work. Its advantage is that it allows a method to be tailored for a particular application, where selectivity, resolution, and total run time are optimized. Be aware poor peak resolution can be due to any or all of the following factors.



### 7.3.1 Loss of Column Efficiency

- A. Check to see if headspace has developed in the guard or analytical column. This is usually due to improper use of the column such as exposing it to high pressures. Remove the column's inlet end fitting (see Section 7.1.2, "Replacing Column Bed Support Assemblies"). If the resin does not fill the column body all the way to the top, the column must be replaced.
- B. Extra-column effects can result in sample band dispersion, making the peaks' elution less efficient. Make sure you are using PEEK tubing with an ID of no greater than 0.010" for 4 mm systems or no greater than 0.005" for 2 mm systems to make all eluent liquid line connections between the injection valve and the detector cell inlet. Cut the tubing lengths as short as possible. Check for leaks. For capillary systems, only use precut tubing of the same type.
- C. If tubing is not connected properly from the inlet and outlet of the column, it can cause low efficiency. When installing CarboPac columns, it is recommended to turn off the pump while connecting the column inlet and the column outlet to the detector. This will avoid any slippage of the ferrule under high pressure conditions which can cause low peak efficiencies.

### 7.3.2 Shortened Retention Times



**NOTE**

*Even with adequate system and column efficiency, resolution of peaks will be compromised if analytes elute too fast.*

- A. Check the flow rate. See if the eluent flow rate is equivalent to the flow rate specified by the analytical protocol. Measure the eluent flow rate after the column using an analytical balance.
- B. Check to see if the eluent compositions and concentrations are correct. An eluent that is too concentrated will cause the peaks to elute faster. Prepare fresh eluent.



**NOTE**

*If you are using a gradient pump to proportion the eluent, components from two or three different eluent reservoirs, the resulting eluent composition may not be accurate enough for the application. Use one reservoir containing the correct eluent composition to see if this is the problem. This may be a problem when one of the proportioned eluents is less than 5%.*

- C. Column contamination can lead to a loss of column capacity. Highly retained contaminant ions will occupy a portion of the anion exchange sites limiting the number of sites available for retention of analyte ions. Refer to "[Appendix A, CarboPac PA100 Column Care](#)", for recommended column cleanup procedures.



**NOTE**

*Possible sources of column contamination are impurities in chemicals and in the deionized water used for eluents or components of the sample matrix. Be especially careful to make sure that the recommended chemicals are used. The deionized water should have a specific resistance of 18.2 megohm-cm.*

- D. Diluting the eluent will improve peak resolution, but will also increase the analytes' retention times. If a 10% dilution of the eluent is not sufficient to obtain the desired peak resolution, or if the resulting increase in retention times is unacceptable, clean the column (see [Appendix A, CarboPac PA100 Column Care](#)).

After cleaning the column, reinstall it in the system and let it equilibrate with eluent for about 30 minutes directing the column effluent to waste. Then connect the column to the electrochemical detector cell. No water wash is necessary. The column is equilibrated when consecutive injections of the standard result in reproducible retention times. The original column capacity should be restored by this treatment, since the contaminants should have been eluted from the column.

**NOTE**

*For assistance, contact Technical Support for Dionex Products. In the U.S., call 1-800-346-6390. Outside the U.S., call the nearest Thermo Fisher Scientific office.*

### 7.3.3 Loss of Resolution for early eluting peaks

If poor resolution or efficiency is observed for early eluting peaks compared to the later eluting peaks, check the following:

- A. Improper eluent concentration may be the problem if retention time is less than expected. If manually prepared eluent is used, remake the eluent as required for your application and ensure that the water and chemicals used are of the required purity. If Dionex eluent generator is used to generate the eluent, check the flow rate, as pump flow rate will affect the eluent concentration.
- B. Column overloading may be the problem. Reduce the amount of sample ions being injected onto the analytical/capillary column by either diluting the sample or injecting a smaller volume onto the column.
- C. Sluggish operation of the injection valve may be the problem due to partially plugged port faces. Refer to the valve manual for instructions.
- D. Improperly swept out volumes anywhere in the system prior to the guard and analytical/capillary columns may be the problem. Swap components, one at a time, in the system prior to the analytical/capillary column and test for early eluting peak resolution after every system change.

### 7.3.4 Spurious Peaks

- A. The columns may be contaminated. If the samples contain an appreciable level of polyvalent ions and the column is used with a weak eluent system, the retention times for the analytes will then decrease and spurious, inefficient (broad) peaks may show up at unexpected times. Clean the column as indicated in “Column Care”.

**NOTE**

*For assistance, contact Technical Support for Dionex Products. In the U.S., call 1-800-346-6390. Outside the U.S., call the nearest Thermo Fisher Scientific office.*

- B. The injection valve may need maintenance. When an injection valve is actuated, the possibility of creating a baseline disturbance exists. This baseline upset can show up as a peak of varying size and shape. This will occur when the injection valve needs to be cleaned or retorqued (see injection valve manual). Check to see that there are no restrictions in the tubing connected to the valve. Also check the valve port faces for blockage and replace them if necessary. Refer to the Valve Manual for troubleshooting and service procedures. Small baseline disturbances at the beginning or at the end of the chromatogram can be overlooked as long as they do not interfere with the quantification of the peaks of interest.

### 7.3.5 No Peaks, Poor Peak Area Reproducibility or too Small Peak Areas

- A. Check the position and filling levels of sample vials in the autosampler.
- B. Check injector needle-height setting.
- C. Check each line of the schedule for proper injector parameters. Revert to full loop and column appropriate sample loop size.
- D. Service the injection valve (check for leaks, Tefzel fragments, or sediments inside the valve).

### 7.3.6 Large Baseline Dip in the Chromatogram

A large baseline dip appearing is usually caused by oxygen in the sample injected. The ‘oxygen dip’ is normal and can be reduced in magnitude with higher NaOH concentration in the eluent.

### 7.3.7 Unidentified Peaks Appear with Expected Analyte Peaks

During the acetate or hydroxide gradient, a number of small peaks may appear. These peaks are usually due to trace contaminants in the water supply. The contaminants accumulate on the column during the isocratic section of the chromatogram and are released, frequently as irregular baseline deformations or sharp spikes, with the increasing eluent strength.

Some trace contaminants can co-elute with monosaccharides, compromising accuracy of quantitation at lower concentrations. If extraneous peaks are observed even after the water supply is excluded as a possible cause, clean the autosampler lines and sample loop. The autosampler should be cleaned using the following protocol:

- A. Disconnect the column and detector cell from the autosampler.
- B. Set the pump to 100% deionized water.
- C. Place the following solutions in the autosampler and inject in sequence. Use 25  $\mu$ L full loop injections:
  1. 1 M NaOH
  2. Deionized water
  3. IPA
  4. Deionized water
  5. 1 M HCl
  6. Deionized water

### 7.3.8 Decreased Detection Sensitivity

Always confirm the loss of response by performing at least one injection of the system suitability standard mix as described in [Section 6.1](#). This is to make sure that a decreased level of response is not being caused by system problems.

Any decrease in detection sensitivity means that the working electrode surface has been affected. The operator should install a new working electrode. Spare gold working electrodes should always be available in order to avoid unnecessary delays.

Exception:

Check the pH reading. If the value is out of range or  $>13.2$ , install a new reference electrode and then install a new gold working electrode. The system cleanup is not necessary. The decrease in sensitivity was caused by a gold-oxide-buildup on the electrode surface because the reference potential was too high. The non-disposable gold working electrode can be reconditioned by polishing.

After installing a new working electrode (with or without the complete system cleanup), confirm the normal detection sensitivity. Carry out a test with a reference standard. Should the response be too low, immediately remove the new working electrode from the system.

### 7.3.9 Excessive Gradient Rise

The magnitude of the gradient rise can be minimized by running high eluent strengths during the times when the system is not in use for sample or standard analysis. This will keep the column conditioned, free from carbonate buildup, and ready for analysis.

- A. Make sure the gradient rise is not caused by the system and/or detector cell.
- B. Increase column temperature to 40 °C and wash the guard and column with 1M NaOH or KOH for at least four hours (preferably overnight). Run a blank gradient at 30 °C and if necessary repeat the clean up with 100 mM NaOH, 950 mM sodium acetate wash at 40 °C.

## 7.4 Reconditioning or Replacement of the Gold (conventional or disposable) Electrodes or Replacement of the Reference Electrode

Refer to Product Manual for Disposable Electrodes Doc. No. 065040, Dionex ICS-5000 Ion Chromatography System Manual Doc. No. 065342 and User's Compendium for Electrochemical Detection Doc. No. 065340 for any help necessary with electrochemical detection, working and reference electrodes.

# Appendix A – CarboPac PA100 Column Care

## A.1 Recommended Operation Pressures

Operating a column above its recommended pressure limit can cause irreversible loss of column performance. The maximum recommended operating pressure for Dionex CarboPac PA100 column is 5,000 psi (34.47 MPa).

## A.2 Column Start-Up

The CarboPac columns are shipped using Sodium hydroxide (see QAR) as the storage solution. Prepare the eluent shown on the Quality Assurance Report (QAR), install the column in the chromatography module and direct the column effluent to waste for 60 minutes, and then connect to the ED cell. Test the column performance under the conditions described in the QAR. Continue making injections of the test standard until consecutive injections of the standard give reproducible retention times. Equilibration is complete when consecutive injections of the standard give reproducible retention times.

If chromatographic efficiency or resolution is poorer than the QAR, see Sections 7.3 Poor Resolution and Section 7.3.1 Loss of Column Efficiency.

### IMPORTANT

*When making any tubing connections (column installation, replacing tubing etc), it is recommended to make these connections with the pump turned off. This will avoid any slippage of the ferrule under high pressure conditions.*

## A.3 Column Storage

For short-term storage (< 1 week), use Eluent, for long-term storage (> 1 week), see QAR. Flush the column for a minimum of 10 minutes with the storage solution. Cap both ends securely, using the plugs supplied with the column.

## A.4 CarboPac PA100 Column Cleanup

The CarboPac PA100 can be readily cleaned by an approximate 50-column-volume rinse with 0.2–1 M NaOH, and/or strong sodium acetate (e.g., 1 M). More stubborn contamination problems may necessitate thoroughly cleaning the column. Use the following steps to thoroughly clean the CarboPac PA100 at a flow rate similar to application flow rate or half the application flow rate to avoid the over-pressurization of the column:

- A. Disconnect column from the ED cell and direct the column effluent to waste. If your system is configured with both a guard column and an analytical column, reverse the order of the guard and analytical column in the eluent flow path. Double check that the eluent flows in the direction designated on each of the column labels.



**CAUTION**

*When cleaning an analytical column and a guard column in series, ensure that the guard column is placed after the analytical column in the eluent flow path. Contaminants that have accumulated on the guard column can be eluted onto the analytical column and irreversibly damage it. If in doubt, clean each column separately.*

- B. Wash the CarboPac PA100 with deionized water (18.2 megohm-cm) for about 30 minutes and then clean with 1 M Methanesulfonic acid (MSA) for one to two hours at appropriate flow rate (1mL/min for 4mm and 0.25mL/min for 2mm column).
- C. Wash the column with deionized water for about 30 minutes.
- D. Then clean the CarboPac PA100 with 200 mM NaOH for at least one hour.
- E. Reconnect column to the cell and equilibrate the CarboPac PA100 to the desired initial conditions and test the performance using QAR standard and eluent.

# Appendix B – Quality Assurance Reports

Dionex CarboPac™ PA100

Analytical (2 x 250 mm)

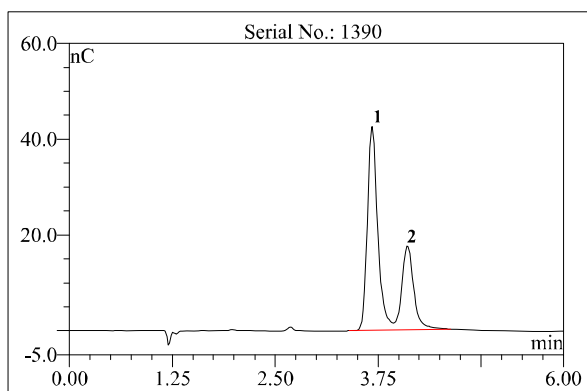
Product No. 057182

Date: 10-Dec-12 13:24

Serial No. : 001390

Lot No. : 012-26-041

**Eluent:** 100 mM NaOH/ 100mM Sodium Acetate  
**Eluent Flow Rate:** 0.25 mL/min  
**Temperature:** Ambient Temperature  
**Detection:** Electrochemical Detection  
**Injection Volume:** 2.5 µL  
**Storage Solution:** Eluent

**ED40 Operating Parameter**

Time	Potential <sup>1</sup>	Integration
0.00	0.10	
0.20	0.10	Begin
0.40	0.10	End
0.41	-2.00	
0.42	-2.00	
0.43	0.60	
0.50	-0.10	

<sup>1</sup> Reference Electrode Mode: Ag/AgCl

No.	Peak Name	Ret.Time (min)	Asymmetry (AIA)	Resolution (EP)	Efficiency (EP)	Amount Injected (nmoles)
1	alpha-(2,6)-NAN-lactosc	3.68	1.3	1.80	4470	*Unknown
2	alpha-(2,3)-NAN-lactosc	4.10	1.2	n.a.	4565	0.03

\* Exact concentration is not known

**QA Results:**

Analvte	Parameter	Specification	Results
alpha-(2,3)-NAN-lactosc	Efficiency	>=3600	Passed
alpha-(2,3)-NAN-lactosc	Retention Time	3.97-4.74	Passed
alpha-(2,3)-NAN-lactosc	Asymmetry	0.9-1.7	Passed
	Pressure	<=2420	1566

*Production Reference:*

Datasource: QAR  
 Directory: CarboPac\PA100  
 Sequence: CP\_PA100\_2X250MM  
 Sample No.: 2

6.80 SR11 Build 3161 (184582) (Demo-Installation)

Chromleon™ Thermo Fisher Scientific

066702-04 (QAR)

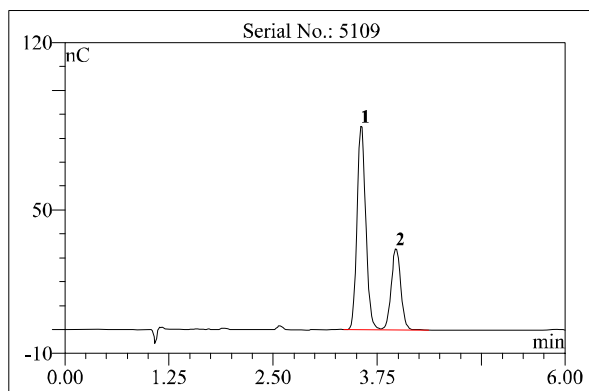
**Dionex CarboPac™ PA100**  
**Analytical (4 x 250 mm)**  
**Product No. 043055**

Date: 03-Oct-12 11:52

Serial No. : 005109

Lot No. : 011-15-088

**Eluent:** 100 mM NaOH/ 100mM Sodium Acetate  
**Eluent Flow Rate:** 1.0 mL/min  
**Temperature:** Ambient Temperature  
**Detection:** Electrochemical Detection  
**Injection Volume:** 10 µL  
**Storage Solution:** Eluent

**ED40 Operating Parameter**

Time	Potential <sup>1</sup>	Integration
0.00	0.10	
0.20	0.10	Begin
0.40	0.10	End
0.41	-2.00	
0.42	-2.00	
0.43	0.60	
0.50	-0.10	

<sup>1</sup> Reference Electrode Mode: Ag/AgCl

No.	Peak Name	Ret.Time (min)	Asymmetry (AIA)	Resolution (EP)	Efficiency (EP)	Amount Injected (nmoles)
1	alpha-(2,6)-NAN-Lactose	3.6	1.3	2.06	5326	*Unknown
2	alpha-(2,3)-NAN-Lactose	4.0	1.2	n.a.	5700	0.12

\* Exact concentration is not known

**QA Results:**

Analyte	Parameter	Specification	Results
alpha-(2,3)-NAN-Lactose	Efficiency	>=4500	Passed
alpha-(2,3)-NAN-Lactose	Retention Time	3.9-4.4	Passed
alpha-(2,3)-NAN-Lactose	Asymmetry	0.9-1.7	Passed
	Pressure	<=2420	2069

*Production Reference:*

Datasource: QAR

Directory CarboPac\PA100

Sequence: CP\_PA100\_4X250mm

Sample No.: 2

6.80 SR11 Build 3161 (184582) (Demo-Installation)

Chromleon™ Thermo Fisher Scientific

066988-04 (QAR)



**Dionex CarboPac™ PA100**  
**Semi-Prep (9 x 250 mm)**  
**Product No. SP2089**

Date: 27-Sep-12 14:31

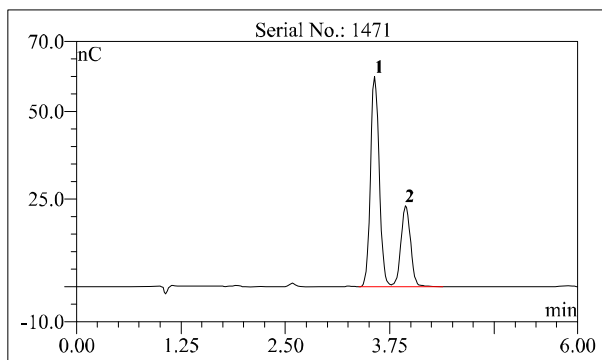
Serial No. : 001471

Lot No. : 012-26-041

**Eluent:** 100 mM NaOH/ 100mM Sodium Acetate  
**Eluent Flow Rate:** 5.0 mL/min  
 (Flow through cell approx. 1.0 mL/min with the rest flowing to the waste)  
**Temperature:** Ambient Temperature  
**Detection:** Electrochemical Detection  
**Injection Volume:** 50 µL  
**Storage Solution:** Eluent

**ED40 Operating Parameter**

Time	Potential <sup>1</sup>	Integration
0.00	0.10	
0.20	0.10	Begin
0.40	0.10	End
0.41	-2.00	
0.42	-2.00	
0.43	0.60	
0.50	-0.10	

<sup>1</sup> Reference Electrode Mode: Ag/AgCl

No.	Peak Name	Ret.Time (min)	Asymmetry (AIA)	Resolution (EP)	Efficiency (EP)	Amount Injected (nmoles)
1	alpha-(2,6)-NAN-Lactose	3.6	1.2	1.85	5686	*Unknown
2	alpha-(2,3)-NAN-Lactose	3.9	1.2	n.a.	5697	0.6

\* Exact concentration is not known

**QA Results:**

Analyte	Parameter	Specification	Results
alpha-(2,3)-NAN-Lactose	Efficiency	>=4500	Passed
alpha-(2,3)-NAN-Lactose	Retention Time	3.9-4.4	Passed
alpha-(2,3)-NAN-Lactose	Asymmetry	0.9-2.0	Passed
	Pressure	<=2420	1813

**Production Reference:**

Datasource: QAR

Directory CarboPac\PA100

Sequence: CP\_PA100\_9X250MM

Sample No.: 2

6.80 SR11 Build 3161 (184582) (Demo-Installation)

Chromeleon™ Thermo Fisher Scientific

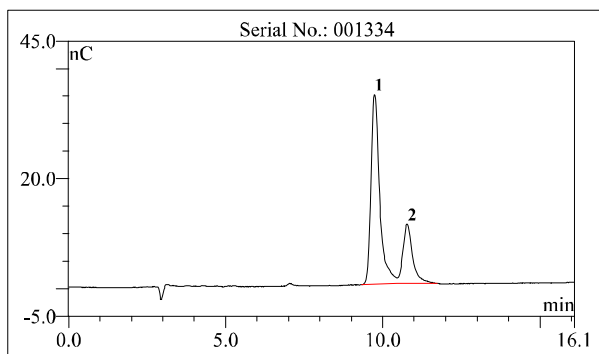
067615-04 (QAR)

**Dionex CarboPac™ PA100****Prep (22 x 250 mm)****Product No. SP2667****Date:** 13-Jan-12 16:35**Serial No. :** 001334**Lot No. :** 011-15-088

**Eluent:** 100 mM NaOH/ 100mM Sodium Acetate  
**Eluent Flow Rate:** 10.0 mL/min  
 (Flow through cell approx. 1.0 mL/min with the rest flowing to the waste)  
**Temperature:** Ambient Temperature  
**Detection:** Electrochemical Detection  
**Injection Volume:** 250 µL  
**Storage Solution:** Eluent

**ED40 Operating Parameter**

Time	Potential <sup>1</sup>	Integration
0.00	0.10	
0.20	0.10	Begin
0.40	0.10	End
0.41	-2.00	
0.42	-2.00	
0.43	0.60	
0.50	-0.10	

<sup>1</sup> Reference Electrode Mode: Ag/AgCl

No.	Peak Name	Ret.Time (min)	Asymmetry (AIA)	Resolution (EP)	Efficiency (EP)	Amount Injected (nmoles)
1	alpha-(2,6)-NAN-Lactose	9.7	1.8	2.10	6915	*Unknown
2	alpha-(2,3)-NAN-Lactose	10.8	1.4	n.a.	6548	3.0

\* Exact concentration is not known

**QA Results:**

Analyte	Parameter	Specification	Results
alpha-(2,3)-NAN-Lactose	Efficiency	>=4500	Passed
alpha-(2,3)-NAN-Lactose	Retention Time	9.7-14.1	Passed
alpha-(2,3)-NAN-Lactose	Asymmetry	0.9-2.0	Passed
	Pressure	<=2420	750

*Production Reference:*

Datasource: QAR  
 Directory: CarboPac\PA100  
 Sequence: CP\_PA100\_22x250mm  
 Sample No.: 3

6.80 SR11 Build 3161 (184582) (Demo-Installation)

Chromeleon™ Thermo Fisher Scientific

067238-04 (QAR)