

Separation of Large Double-Stranded DNA (dsDNA) Fragments Using a Wide-Pore Reversed-Phase Chromatography Column

Julia Baek, Jim Thayer, Shanhua Lin, Xiaodong Liu, Thermo Fisher Scientific, Sunnyvale, CA

Key Words

Ion-pair reversed-phase chromatography, double-stranded DNA (dsDNA), DNAPac RP, HPLC

Goal

To demonstrate the separation of dsDNA fragments and a PCR product using a wide-pore, polymer-based DNAPac RP column.

Introduction

Restriction nucleases are enzymes that generate dsDNA fragments of various sizes, and help researchers construct plasmids for expression of modified or novel proteins. In addition, DNAs that contain sequences for alternative gene editing systems (such as CRISPR-Cas9) may also generate plasmids.¹ Restriction enzymes may also be used to identify single nucleotide polymorphisms (SNPs) where the base change alters the specific restriction nuclease site.² For all these applications, restriction nuclease DNA fragments are typically purified prior to subsequent use. Agarose and polyacrylamide gels are most often used for separating and recovering dsDNA fragments. Gel electrophoresis is followed by manual excision of the target dsDNA size from the gel and extraction of the DNA from the excised gel. The method is laborious and time consuming and produces relatively low yields. It also requires handling of toxic materials such as acrylamide or ethidium bromide. HPLC provides reliably higher yields, allows direct DNA quantitation, and can be automated using fraction collection. Here we describe the use of a new reversed-phase column for the separation and recovery of dsDNA molecules.

The Thermo Scientific™ DNAPac™ RP column is based on a 4 μm , wide-pore polymer resin, which is stable at high pH and high temperatures. The wide-pore resin (1000–2000 Å) separates large dsDNA molecules as well as single-stranded oligonucleotides. This application note demonstrates the separation of DNA fragments generated from restriction enzymes and DNA ladder components, with automated purification and collection of the PCR product using ion-paired reversed-phase chromatography on the DNAPac RP column.



Experimental

Chemicals and Reagents

- Deionized (DI) water, 18.2 M Ω -cm resistivity
- Acetonitrile (Fisher Scientific™ P/N A955-4)
- Triethylammonium acetate (TEAA) 2.0 M (Applied Biosystems P/N 400613)

Samples

- pBR322-BsuRI digest (Thermo Scientific P/N SM0252)
- Φ X174-BsuRI digest (Thermo Scientific P/N SM0271)
- Thermo Scientific™ FastRuler™ Low Range DNA Ladder, ready-to-use (P/N SM1103)
- Thermo Scientific FastRuler Middle Range DNA Ladder, ready-to-use (P/N SM1113)
- Thermo Scientific FastRuler High Range DNA Ladder, ready-to-use (P/N SM1123)
- WPRE PCR product (Generous gift from Sanjay Vasu)

Sample Handling and Electrophoresis Equipment

Vial and closures: 300 μ L PP 8-425 Screw Thread Vial (P/N C4013-11), Pre-Assembled Black PP 8-425 Cap with Pre-Slit PTFE/Silicone Septa (P/N C4013-77A)

Agarose Gels: Thermo Scientific™ Invitrogen™ 1.2% E-Gel (P/N G501801) with Invitrogen E-Gel Base

Power supply: FisherBiotech™ FB200Q Electrophoresis Power Supply or equivalent

Visualization: UV/Vis transilluminator

Sample Preparation

pBR322-BsuRI and Φ X174-BsuRI digests were diluted five-fold with DI water to make 100 μ g/mL solutions.

FastRuler DNA Ladders and the PCR product were used as delivered.

Separation Conditions

Instrumentation	1. Thermo Scientific™ Vanquish™ Flex UHPLC system consisting of:
	System Base (P/N VF-S01-A)
	Quaternary Pump Flex (P/N VF-P20-A)
	Split-Loop Sampler FT (P/N VF-A10-A)
	Column Compartment H (P/N VH-C10-A)
	Active Pre-heater (P/N 6732.0110)
	Diode Array Detector HL (P/N VH-D10-A)
	Thermo Scientific™ LightPipe™ Flow Cell, Standard, 10 mm (P/N 6083.0100)
	2. Thermo Scientific™ Dionex™ UltiMate™ 3000 BioRSLC system equipped with:
	SR-3000 Solvent Rack (without degasser) (P/N 5035.9200)
	LPG-3400RS Biocompatible Quaternary Rapid Separation Pump with degasser (P/N 5040.0036)
	WPS-3000TBFC Thermostatted Biocompatible Pulled-Loop Well Plate Autosampler with Integrated Fraction Collection (P/N 5825.0020)
	TCC-3000RS Rapid Separation Thermostatted Column Compartment (P/N 5730.0000) equipped with 11 μ L pre-column heater (P/N 6723.0252)
	VWD-3400RS Rapid Separation Variable Wavelength Detector (P/N 5074.0010) equipped with a micro flow cell (PN 6074.0300)
Columns	DNAPac RP, 2.1 \times 100 mm (P/N 088923) DNAPac RP, 2.1 \times 50 mm (P/N 088924)
Mobile Phase A	0.1 M TEAA, pH 7.0
Mobile Phase B	0.1 M TEAA in water/acetonitrile (75:25 v/v)
Gradient	As specified in Figures
Flow Rate	As specified in Figures
Column Temperature	As specified in Figures
UV Detector Wavelength	260 nm

Data Processing

The Thermo Scientific™ Dionex™ Chromeleon™ 7.2 Chromatography Data System was used for data analysis.

Results and Discussion

Ion-pair reversed-phase chromatography on the DNAPac RP column was used to separate restriction digest components (Figure 1). At 55 °C, all of the dsDNA digestion fragments from pBR322 plasmid (Figure 1a) and Φ X174 viral DNA (Figure 1b) with BsuRI enzyme were separated. The DNAPac RP column was able to separate fragments differing by as little as a single base

pair (bp) up to 100 bp. In addition, the retention of the fragments correlated to the length of the DNA. Retention versus log (number of base pairs) of both pBR322 and Φ X174 digests was plotted in Figure 1c. The data fit well to a second-order polynomial equation, indicating good correlation between the retention time and the length of the fragments.

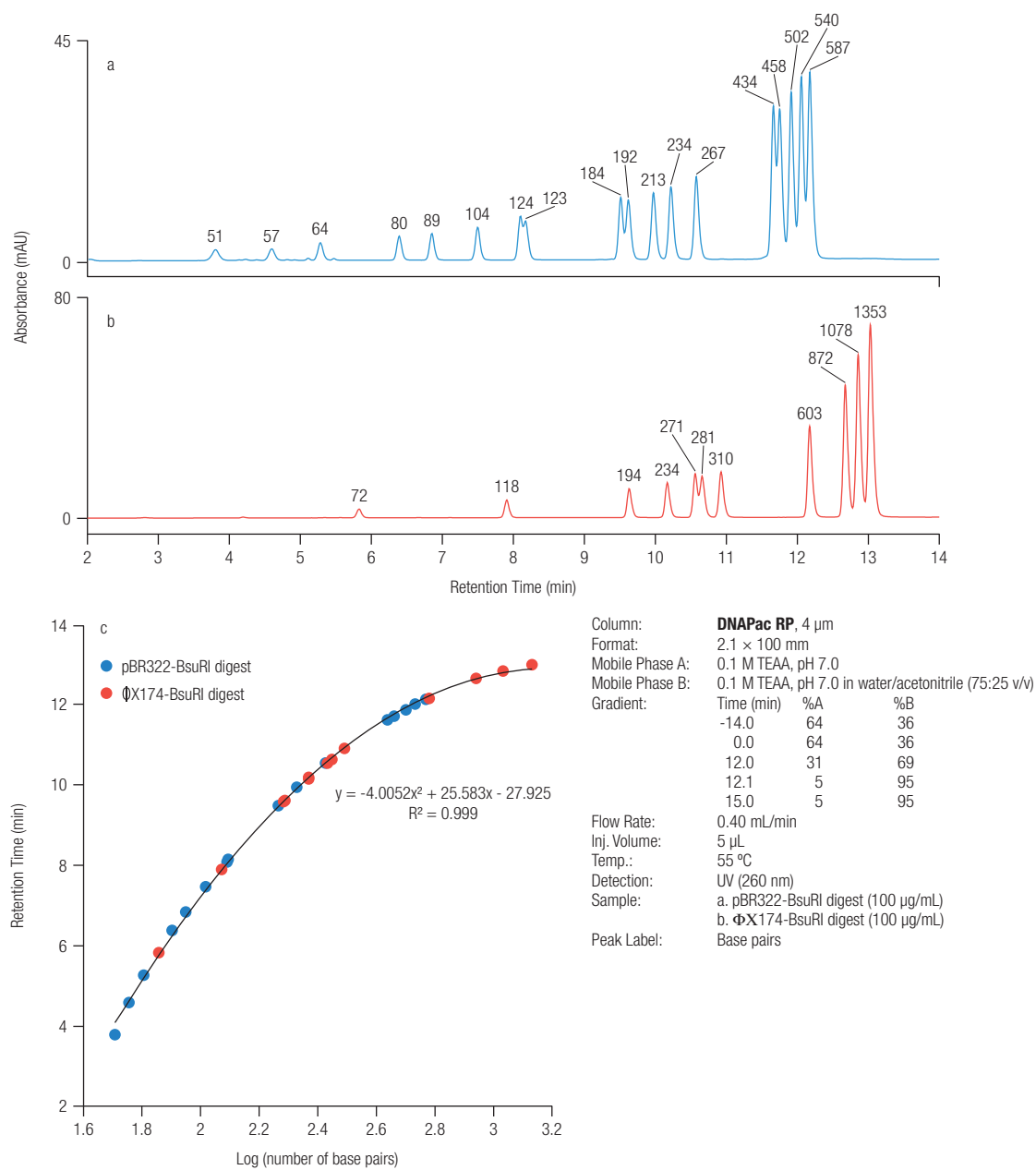
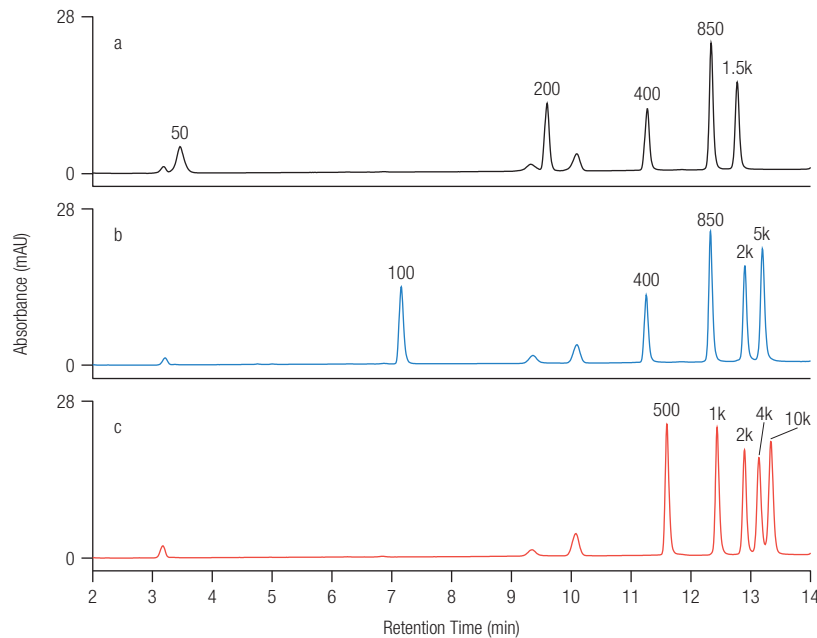


Figure 1. Separation of restriction enzyme digests.

To further investigate the DNAPac RP column resolution of large dsDNA fragments, DNA ladders with components up to 10,000 base pairs were examined (Figure 2). All of the DNA ladder components were resolved in less than 14 minutes with the resolution of

components differing by approximately 10% for fragments up to around 10,000 bp. The DNA ladder retention time can be compared to the retention time of a DNA fragment to estimate its size.



Column:	DNAPac RP , 4 μ m		
Format:	2.1 \times 100 mm		
Mobile Phase A:	0.1 M TEAA, pH 7.0		
Mobile Phase B:	0.1 M TEAA, pH 7.0 in water/acetonitrile (75:25 v/v)		
Gradient:	Time (min)	%A	%B
	-14.0	65	35
	0.0	65	35
	12.0	30	70
	12.1	10	90
	15.0	10	90
Flow Rate:	0.40 mL/min		
Inj. Volume:	10 μ L		
Temp.:	55 $^{\circ}$ C		
Detection:	UV (260 nm)		
Sample:	DNA ladders		
	a. Low range		
	b. Middle range		
	c. High range		
Peak label:	Base pairs		

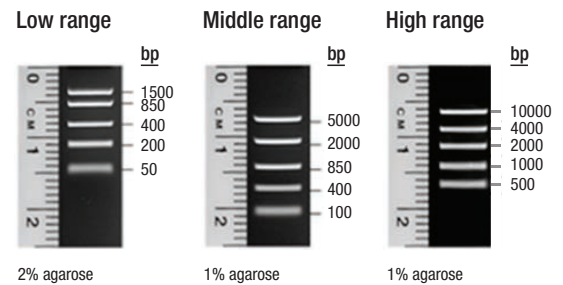


Figure 2. Separation of FastRuler DNA ladders.

Separation of Φ X174-BsuRI digest, which contains components from 72 to 1353 bp, on a DNAPac RP column was compared to competitive columns of similar format (all equivalent 2.1 mm columns, except for brand C, which has a 2.0 mm ID). The same mobile phase and linear velocity were used. The larger pore size

DNAPac RP column gave significantly better separation of the larger dsDNA fragments. Differences in separation can be attributed to larger pore size; however, selectivity, particle size, and particle chemistry may have also contributed to the difference in the separation of dsDNA fragments (Figure 3).

The Φ X174-BsuRI digest was used to study the temperature effect on the dsDNA retention and resolution (Figure 4). Resolution and the retention of these dsDNA fragments increased with temperature up to 50 °C without changing the peak intensity pattern. However, at 60 °C, the retention pattern changed and the resolution of the peaks appeared to be higher. Since the intensities of several peaks change at 60 °C, partial denaturation of some dsDNA fragments is likely induced (hyperchromic shift). Two peaks also eluted earlier at 60 °C than at

50 °C, which is consistent with exposure of ssDNA sections within the fragment. At 70 °C, denaturation is obvious as the number of peaks increased from 11 to 18, and many peaks eluted much earlier than at lower temperatures, suggesting full denaturation of some dsDNA into single stranded DNA (ssDNA) forms. At even higher temperatures (80–90 °C), more peaks were observed and all eluted much earlier, indicating full denaturation of all of the dsDNA fragments.

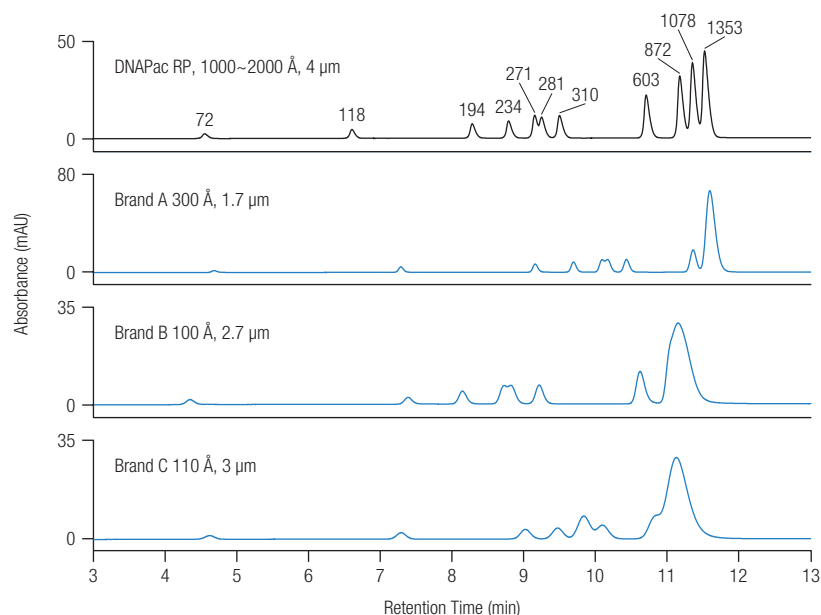


Figure 3. Separation of Φ X174-BsuRI digest on columns with different pore sizes.

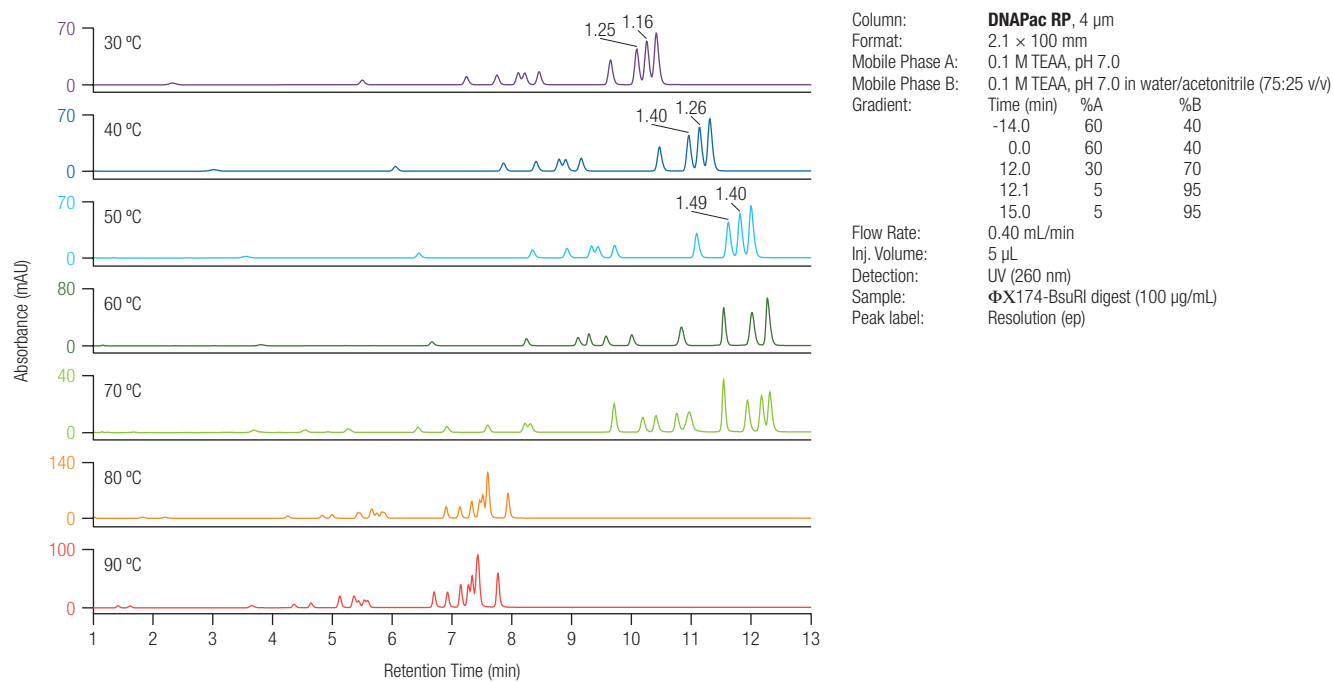
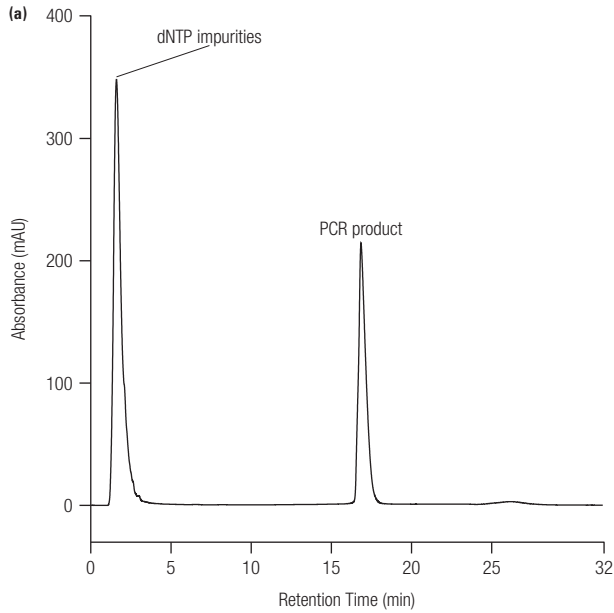


Figure 4. Separation of Φ X174-BsuRI digest at different temperatures.

PCR is routinely used in molecular biology labs for amplification of DNA and sequencing. PCR products are typically purified from the dNTPs and other impurities in the reaction prior to use. Using an UltiMate 3000 system with an UltiMate WPS-3000TBFC (Well Plate Autosampler configured for Fraction Collection), a PCR product was separated from its impurities and

automatically collected into a well plate (Figure 5).^{4,5} The collected sample was then analyzed with a 1.2% agarose E-gel to confirm its purity (Figure 5b). This example demonstrates automated HPLC with fraction collection of dsDNA using the DNAPac RP column, showing facile purification from PCR-related components in 20 minutes.



Column:	DNAPac RP, 4 μm																					
Format:	2.1 \times 100 mm																					
Mobile Phase A:	0.1 M TEAA, pH 7.0																					
Mobile Phase B:	0.1 M TEAA in water/acetonitrile (75:25 v/v)																					
Gradient:	<table border="0"> <thead> <tr> <th>Time (min)</th> <th>%A</th> <th>%B</th> </tr> </thead> <tbody> <tr> <td>0.0</td> <td>56</td> <td>44</td> </tr> <tr> <td>20.0</td> <td>32</td> <td>68</td> </tr> <tr> <td>20.1</td> <td>10</td> <td>90</td> </tr> <tr> <td>22.0</td> <td>10</td> <td>90</td> </tr> <tr> <td>22.1</td> <td>56</td> <td>44</td> </tr> <tr> <td>32.0</td> <td>56</td> <td>44</td> </tr> </tbody> </table>	Time (min)	%A	%B	0.0	56	44	20.0	32	68	20.1	10	90	22.0	10	90	22.1	56	44	32.0	56	44
Time (min)	%A	%B																				
0.0	56	44																				
20.0	32	68																				
20.1	10	90																				
22.0	10	90																				
22.1	56	44																				
32.0	56	44																				
Flow Rate:	0.25																					
Inj. Volume:	50 μ L																					
Temp.:	50 $^{\circ}$ C																					
Detection:	UV (260 nm)																					
Sample:	PCR product																					

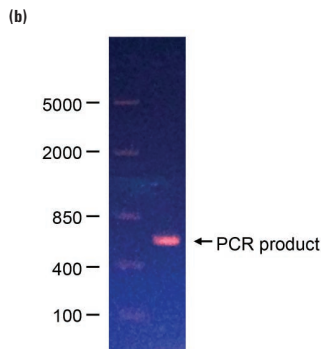


Figure 5. Purification of a PCR product.

Conclusion

- The combination of the DNAPac RP column and Vanquish Flex UHPLC system delivers rapid, high-resolution dsDNA fragment separation up to 10,000 bp.
- Elution of the dsDNA fragments is essentially proportional to fragment length (in bp).
- The UltiMate 3000 system equipped with the WPS-3000 TBFC Well Plate Autosampler configured for Fraction Collection facilitates purification and recovery of dsDNA fragments.

Reference

1. Sander, J.D.; Joung, J.K. CRISPR-Cas systems for editing, regulating and targeting genomes, *Nature Biotechnology*, 2014, 32, 347.
2. Wolff, J.N.; Gemmell, N.J. Combining allele-specific fluorescent probes and restriction assay in real-time PCR to achieve SNP scoring beyond allele ratios of 1:1000, *BioTechniques*, 2008, 44, 193.
3. Baek, J.; Thayer, J.; Lin, S.; Liu, X. Separation of Mixed-Base Oligonucleotides Using a High-Resolution, Reversed-Phase Chromatography Column, Thermo Fisher Scientific AN 21476 (2016).
4. Thayer, J.R., Puri, N., Burnett, C., Hail, M. and S. Rao. Identification of linkage isomers by anion exchange purification with electrospray ionization mass spectrometry of automatically desalted phosphodiesterase-II digests. *Analytical Biochemistry*, 2010, 399, 110.
5. Thayer, J.R., Flook, K.J., Woodruff, A., Rao, S., Pohl, C. New monolith technology for automated anion-exchange purification of nucleic acids. *Journal of Chromatography B*, 2010, 878, 933.

Useful Links

AppsLab Library

The eWorkflow and the Chromeleon Backup (cmbx) file can be downloaded at AppsLab Library:

www.thermofisher.com/appslab

For Research Use Only. Not for use in diagnostic procedures

www.thermofisher.com/BioLC

© 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. Specifications, terms and pricing are subject to change.
Not all products are available in all countries. Please consult your local sales representative for details.

Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand