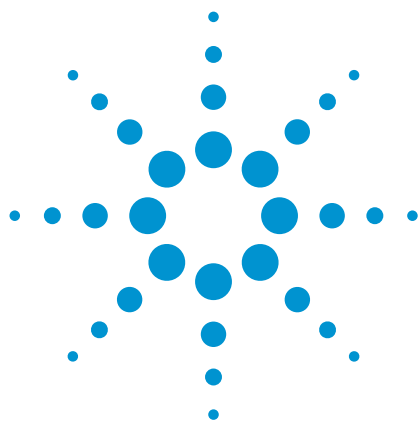




Analysis of Nucleotides Using an Agilent Polaris C18-A Column

Application Note

Pharmaceutical

Author

Anita Der Martirosian
Agilent Technologies, Inc.

Introduction

An HPLC separation of ten nucleotides was achieved in 24 minutes using an Agilent Polaris C18-A column.



Agilent Technologies

Conditions

Column Agilent Polaris C18-A, 4.6 × 150 mm, 5 µm (p/n A2000150X046)
Sample Nucleotide mix, 10
Mobile phase A: H₂O + 0.01 M TBAH + 0.01 M KH₂PO₄
B: CH₃OH + 0.01 M TBAH
Gradient 5% - 50% B in 25 min, hold for 2 min,
ramp to 95% B in 0.1 min, hold for 3 min,
back to 5% B in 0.1 min, hold for 2 min
Flow rate 1.5 mL/min
Temperature Ambient
Detector 254 nm

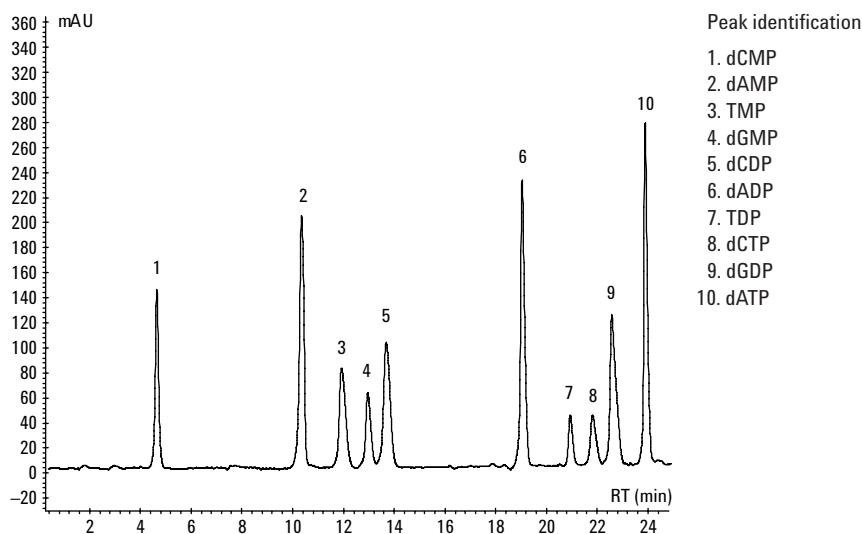


Figure 1. Analysis of nucleotides using an Agilent Polaris C18-A column.

For More Information

These data represent typical results. For more information on our products and services, visit our Web site at www.agilent.com/chem.

www.agilent.com/chem

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice. This publication was originally published in 2008.

© Agilent Technologies, Inc., 2011
Published in USA, July 8, 2011
SI-01261



Agilent Technologies