



Nucleotides

Application Note

BioPharma

Authors

Agilent Technologies, Inc.

Introduction

HPLC of nucleotides is straightforward with Agilent Pursuit 5 μm C8 material in a ChromSep cartridge. Six nucleotides are separated in less than five minutes



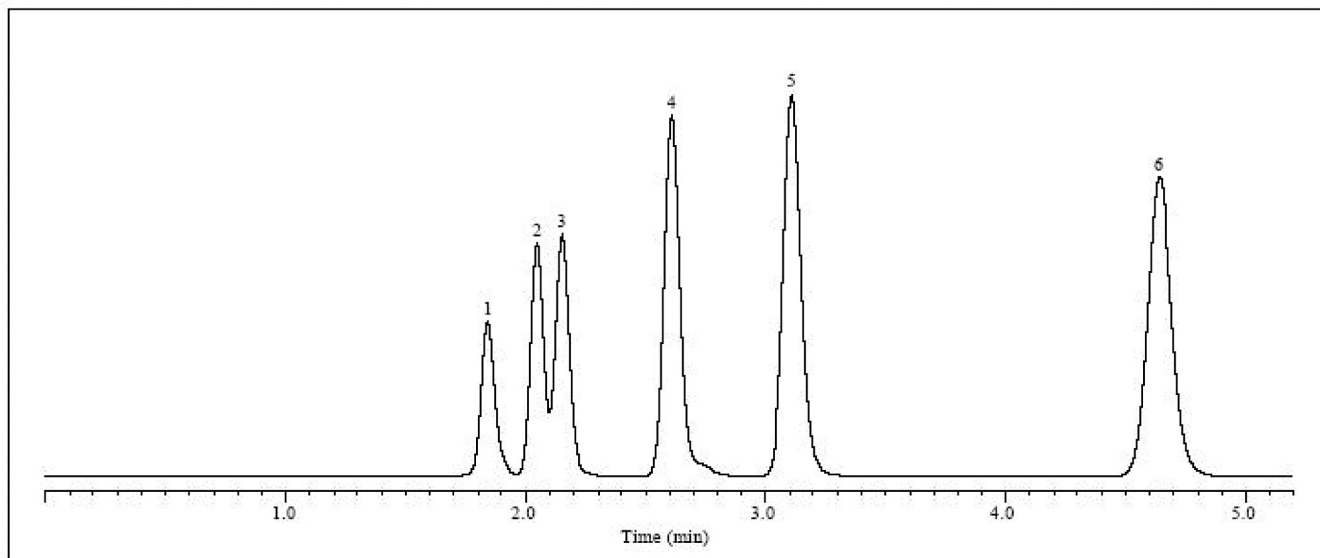
Agilent Technologies

Conditions

Technique : HPLC
Column : Agilent Pursuit C8, 4.6mm x 150 mm, 5 μ
Part no. : Conventional A3030150X046
ChromSep Cartridge System A3030150C046
Temperature : Ambient
Eluent : 50 mM K_2HPO_4 , pH 6.0
Flow rate : 1.0 mL/min
Detector : UV @ 260 nm

Peak identification

1. 5'-CTP
2. 5'-CMP
3. 5'-GDP
4. 5'-GMP
5. 5'-ADP
6. 5'-AMP



www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2011

Printed in the USA

31 October, 2011

First published prior to 11 May, 2010

A02329



Agilent Technologies