



Anti-inflammatory/analgesic drugs

Application Note

BioPharma

Authors

Agilent Technologies, Inc.

Introduction

An Agilent Polaris C18-A HPLC column separates eight anti-inflammatory and analgesic drugs in eight minutes.



Agilent Technologies

Conditions

Column : Agilent Polaris C18-A, 4.6 x 150 mm, 5 μ m
Part No.: A2000150C046 (ChromSep)
Part No.: A2000150X046 (Conventional)

Mobile Phase : A: H₂O + 0.1% HCOOH, B: CH₃CN + 0.1% HCOOH

Gradient : 45% - 80% B in 10 min

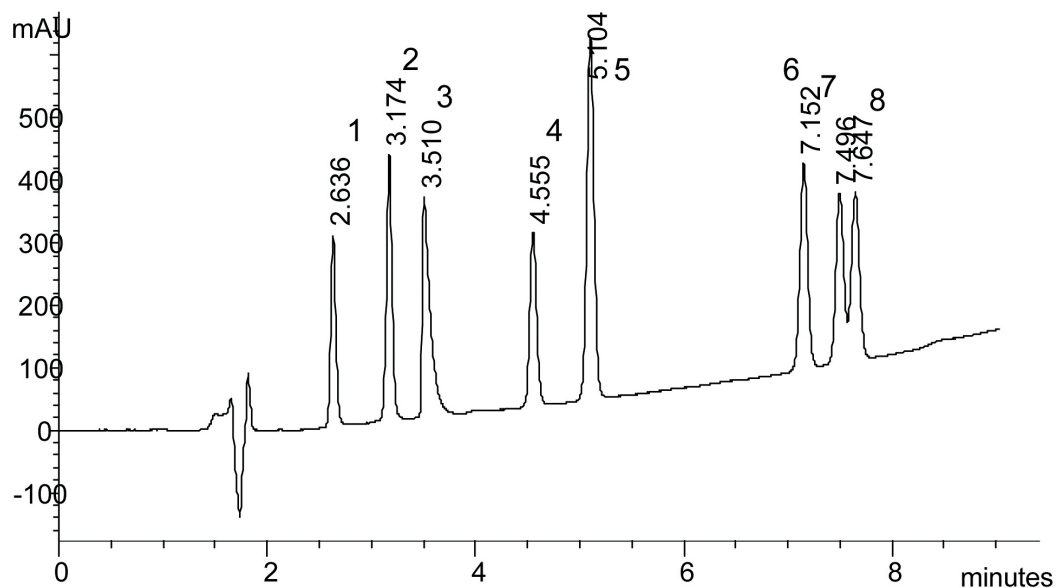
Flow Rate : 1.0 mL/min

Temperature : Ambient

Detection : 220 nm

Peak identification

1. aspirin (acetylsalicylic acid)
2. carbamazepine
3. salicylic acid
4. tolmetin
5. naproxen
6. indomethacin
7. diclofenac
8. ibuprofen



www.agilent.com/chem

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

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

A02217



Agilent Technologies