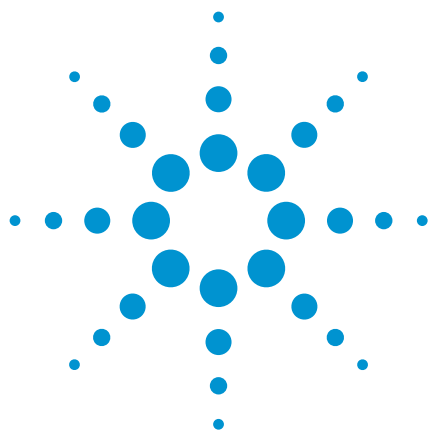




Alkaloids

Application Note

Forensic Toxicology

Authors

Agilent Technologies, Inc.

Introduction

HPLC with an Agilent Pursuit C18 5 μm packing in a ChromSep cartridge separates six alkaloids in 16 minutes.



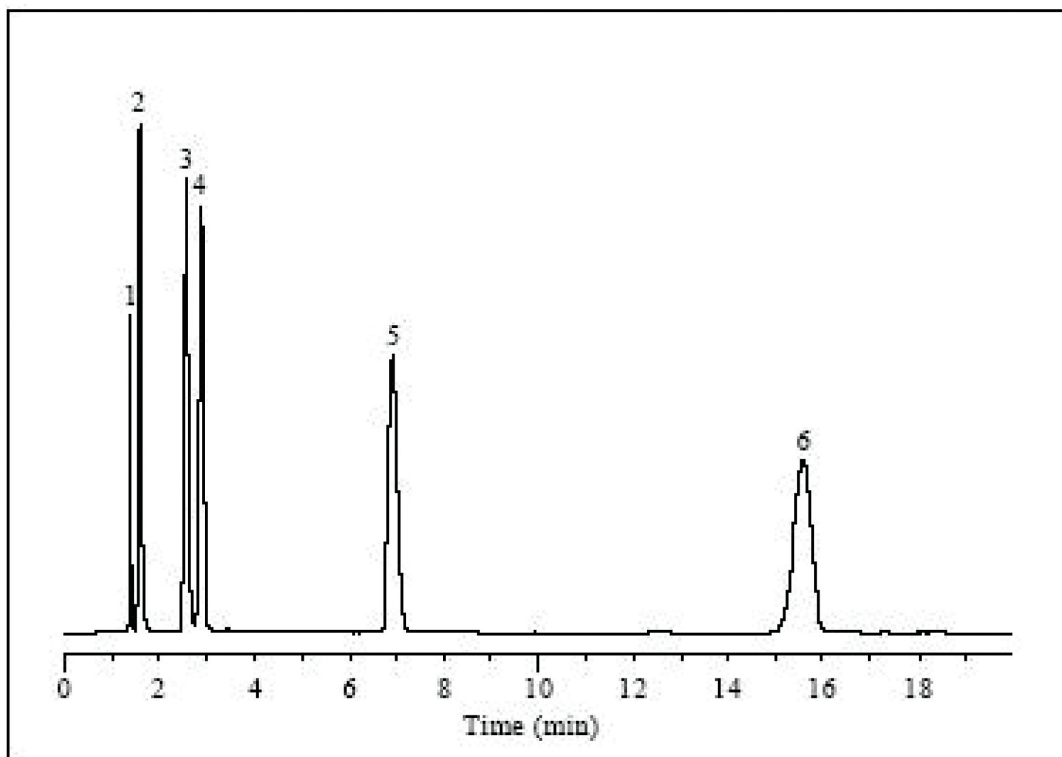
Agilent Technologies

Conditions

Technique : HPLC
Column : Agilent Pursuit C18, 4.6 x 150 mm, 5 µm
Part no. : Conventional A3000150X046
ChromSep Cartridge System A3000150C046
Temperature : Ambient
Eluent : A: MeOH
B: 25 mM KH₂PO₄, pH 6
A:B = 48:52
Flow rate : 1.5 mL/min
Detection : UV @ 214 nm

Peak identification

1. codeine
2. strychnine
3. quinidine
4. quinine
5. papaverine
6. noscapine



www.agilent.com/chem

For Forensic Use only.

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

A02308



Agilent Technologies