



Engelhardt test

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

The Engelhardt test uses eight compounds to assess the hydrophobicity and silanophilic activity of Agilent Pursuit C18 5 μm reversed-phase material in a ChromSep cartridge.



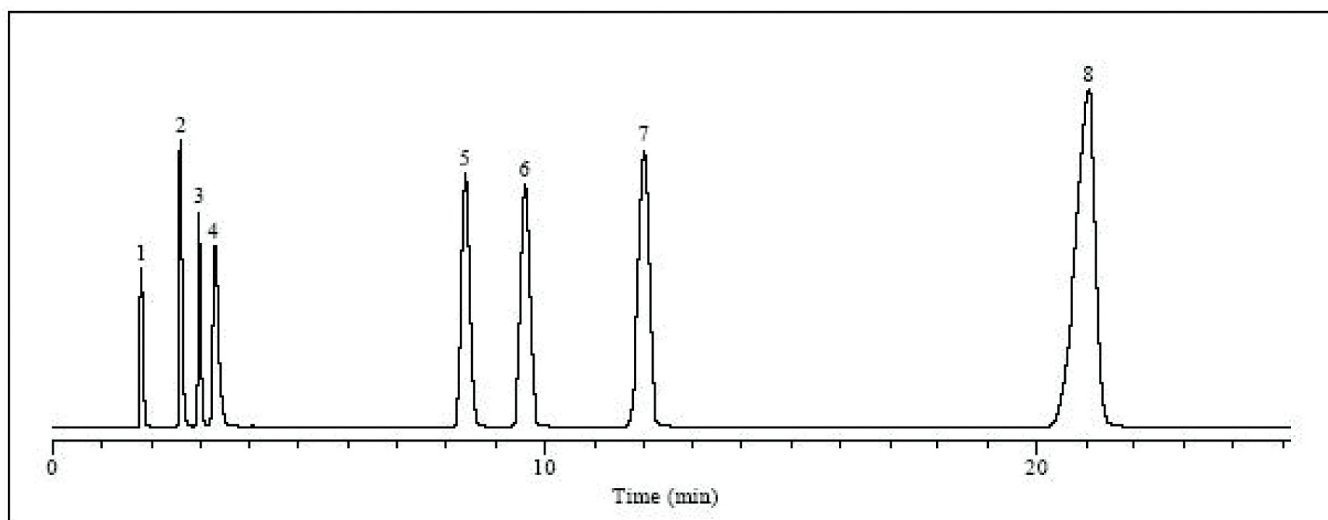
Agilent Technologies

Conditions

Technique : HPLC
Column : Pursuit C18, 4.6 x 150 mm, 5 μ m
Part no. : Conventional A3000150X046
ChromSep Cartridge System A3000150C046
Temperature : Ambient
Eluent : A: MeOH
B: Water
A:B = 55:45
Flow rate : 1.0 mL/min
Detection : UV @ 254 nm

Peak identification

1. thiourea
2. aniline
3. phenol
4. toluidine (o-,m-,p-)
5. n,n-dimethylaniline
6. ethylbenzoate
7. toluene
8. ethylbenzene



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

A02316



Agilent Technologies