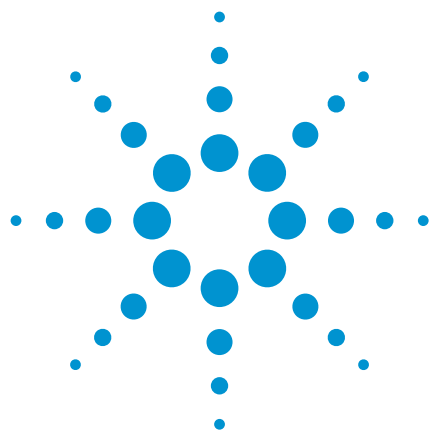




Phthalate esters

Analysis of phthalate esters to EPA 606

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

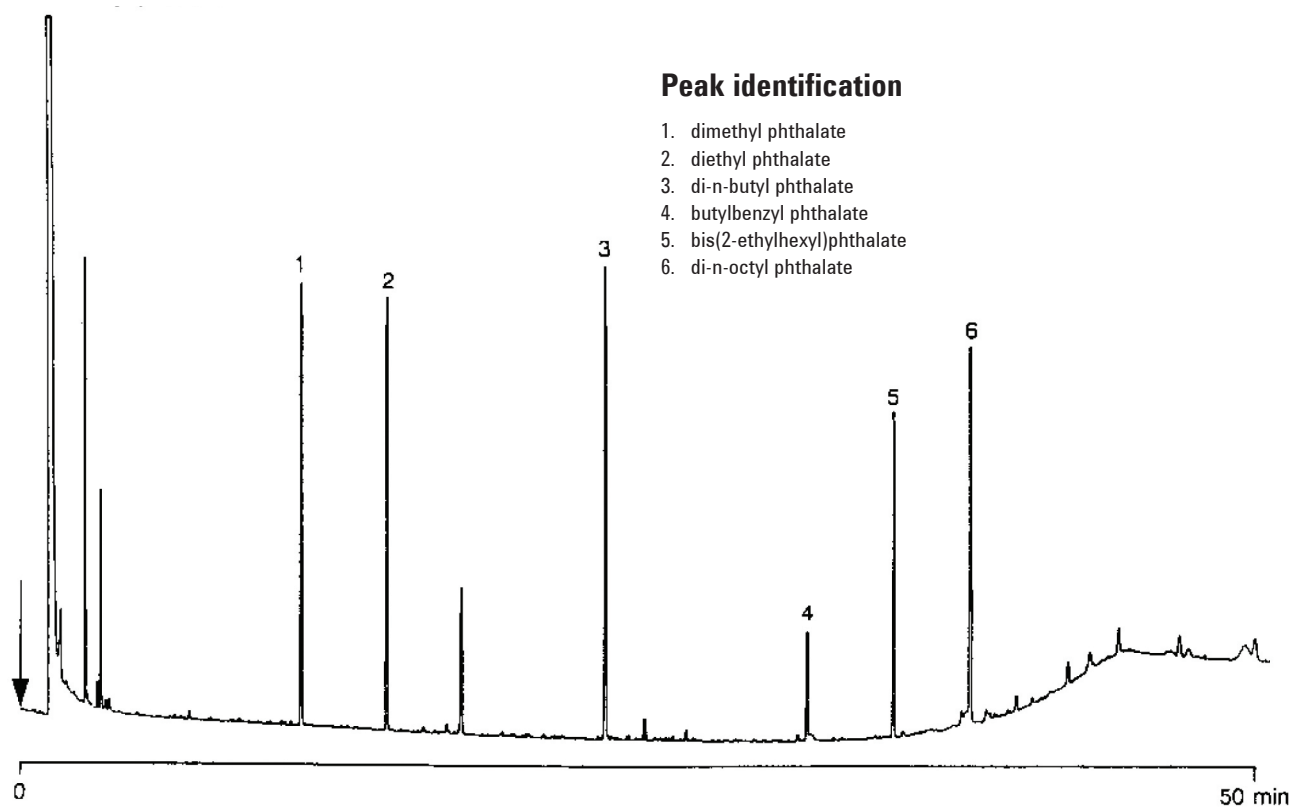
Gas chromatography with an Agilent CP-Sil 8 CB column separates six phthalate esters according to EPA 606 in 50 minutes.



Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 8 CB, 0.32 mm x 50 m fused silica
WCOT CP-Sil 8 CB (df = 0.4 μ m) (Part no. CP7789)
Temperature : 100 °C $\rightarrow$ 300 °C, 10 °C/min
Carrier Gas : H₂, 150 kPa (1.5 bar, 21 psi)
Injector : Split, 100 mL/min
T = 265 °C
Detector : FID
T = 300 °C



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

A00784



Agilent Technologies