



Surfactants

Analysis of Triton X-100

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-SimDist UltiMetal column produces a fingerprint of Triton X-100 derivatized with BSTFA in 35 minutes.

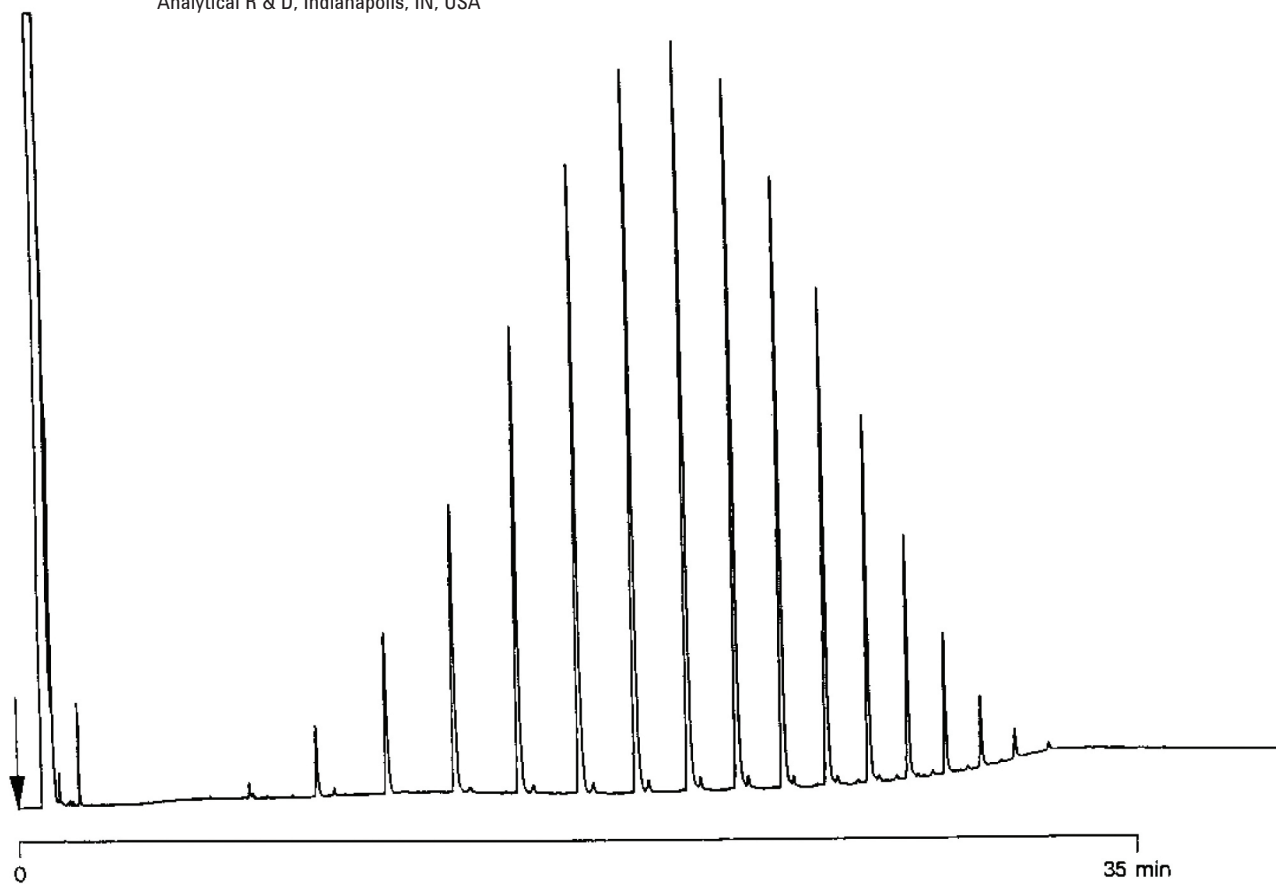


Agilent Technologies

Conditions

Technique : GC-wide-bore
Column : Agilent CP-SimDist CB UltiMetal, 0.53 mm x 15 m
WCOT CP-SimDist CB UltiMetal (df = 0.15 μ m)
This column is available under Part no. CP99921
Temperature : 80 °C \rightarrow 180 °C, 25 °C/min; 180 °C (13.5 min) \rightarrow
400 °C, 8 °C/min
Carrier Gas : He, 35 cm/s
Injector : cool on-column
Detector : FID
T = 400 °C
Solvent Sample : toluene

Courtesy : Jerry Telford, Boehringer Mannheim,
Analytical R & D, Indianapolis, IN, USA



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

A00765



Agilent Technologies