



Prostaglandins

Application Note

BioPharma

Authors

Agilent Technologies, Inc.

Introduction

GC/MS with an Agilent CP-Sil 8 CB column separates six prostaglandins in 16 minutes.



Agilent Technologies

Conditions

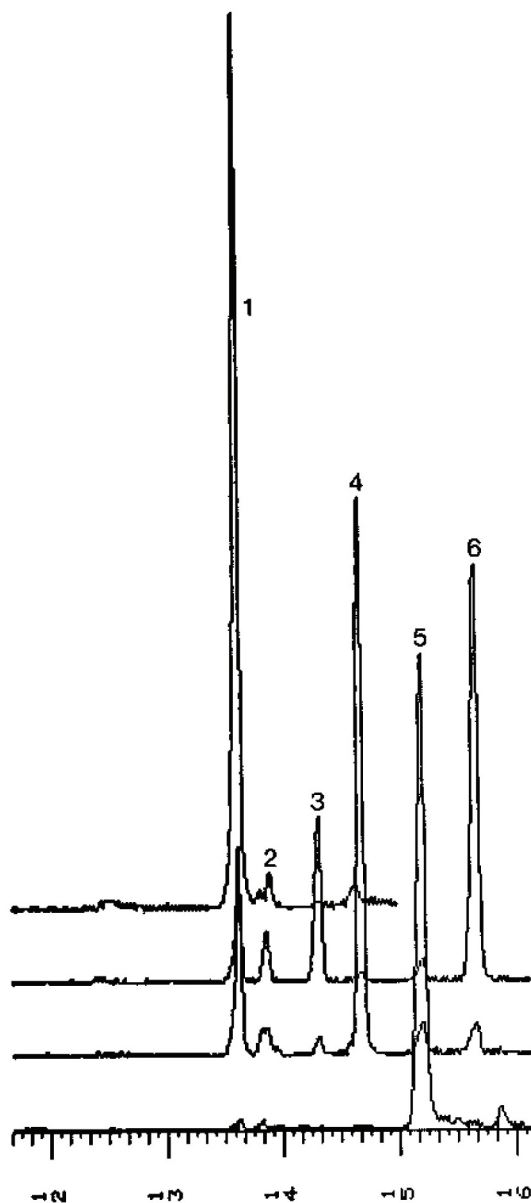
Technique : GC-capillary
Column : Agilent CP-Sil 8 CB, 0.10 mm fused silica WCOT
CP-Sil 8 CB (0.12 μ m) (Custom-made)
Temperature : 150 °C $\rightarrow$ 290 °C, 20 °C/min
Detector : selected ion monitoring MS

Courtesy : Dr. C. Chiabrando and R. Fanelli of the Marlo Negri
Institute, Milan, Italy

Peak identification

1. PGF _{1x}	Mass	3010	4610	5440	5890
2. PGE ₂ (minor isomer)	Ampl.	1	8	8	8
3. PGD ₂	Time.	20	20	20	20
4. PGE ₂ (major isomer)	Scale.	1	1	1	1
5. thromboxane B ₂	Min.	25	28		
6. 6-Keto PGF _{1x}	Max.	5808	6737	5124	10960

(pentafluorobenzyl esters, methyl oxime, trimethyl silyl ether derivatives)



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

A00204



Agilent Technologies