



Gases

Analysis of sulfur gases

Application Note

Energy & Fuels

Authors

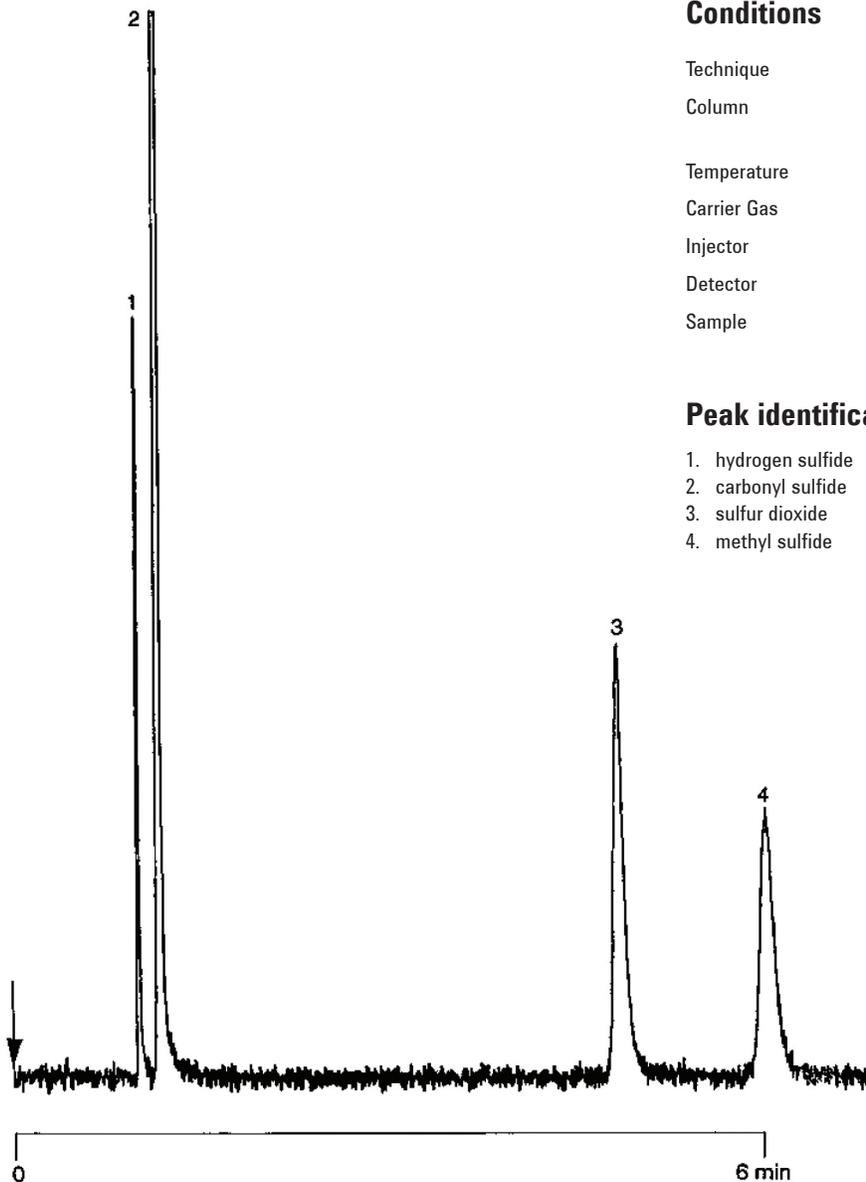
Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT U column separates four sulfur gases in six minutes.



Agilent Technologies



Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT U, 0.53 mm x 25 m (ID x L) fused silica PoraPLOT U (df = 20 μ m) (Part no. CP7584)
Temperature : 50 °C
Carrier Gas : H₂
Injector : Split, 100 mL/min
Detector : FPD
Sample : \pm 100 ppm

Peak identification

1. hydrogen sulfide
2. carbonyl sulfide
3. sulfur dioxide
4. methyl sulfide

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

A00450



Agilent Technologies