

Hydrocarbons, C₁ to C₆

Separation of light hydrocarbons on a medium polar porous polymer

Author

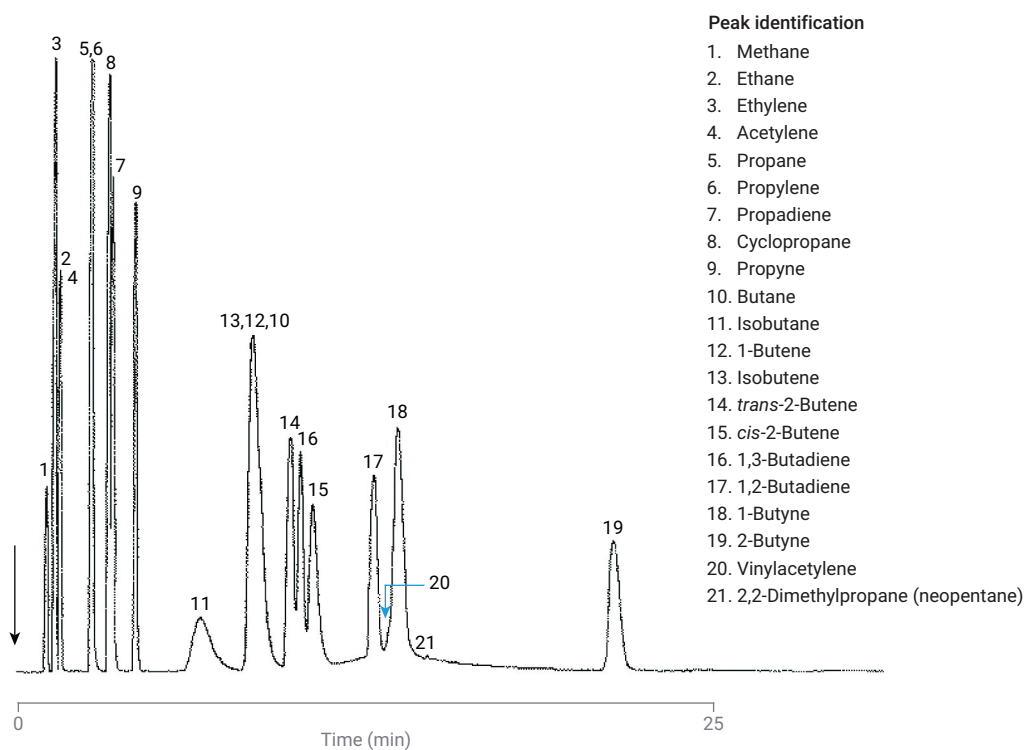
Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT U column separates 21 light hydrocarbons in 25 minutes.

Conditions

Parameter	Value
Technique	GC-wide-bore
Column	Agilent PoraPLOT U, 0.53 mm × 25 m fused silica PLOT (df = 20 µm) (p/n CP7584)
Temperature	75 °C
Carrier Gas	H ₂ , 40 cm/s
Injector	Direct, T = 100 °C
Detector	FID T = 150 °C
Sample Size	6 µL
Concentration Range	0.1% in nitrogen



www.agilent.com

DE01115181

This information is subject to change without notice.

© Agilent Technologies, Inc. 2010, 2023
Printed in the USA, January 26, 2023
A01456