



Amino acids

Application Note

BioPharma

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Chirasil Val column separates 19 amino acids (TFA/isopropylesters) in 33 minutes.



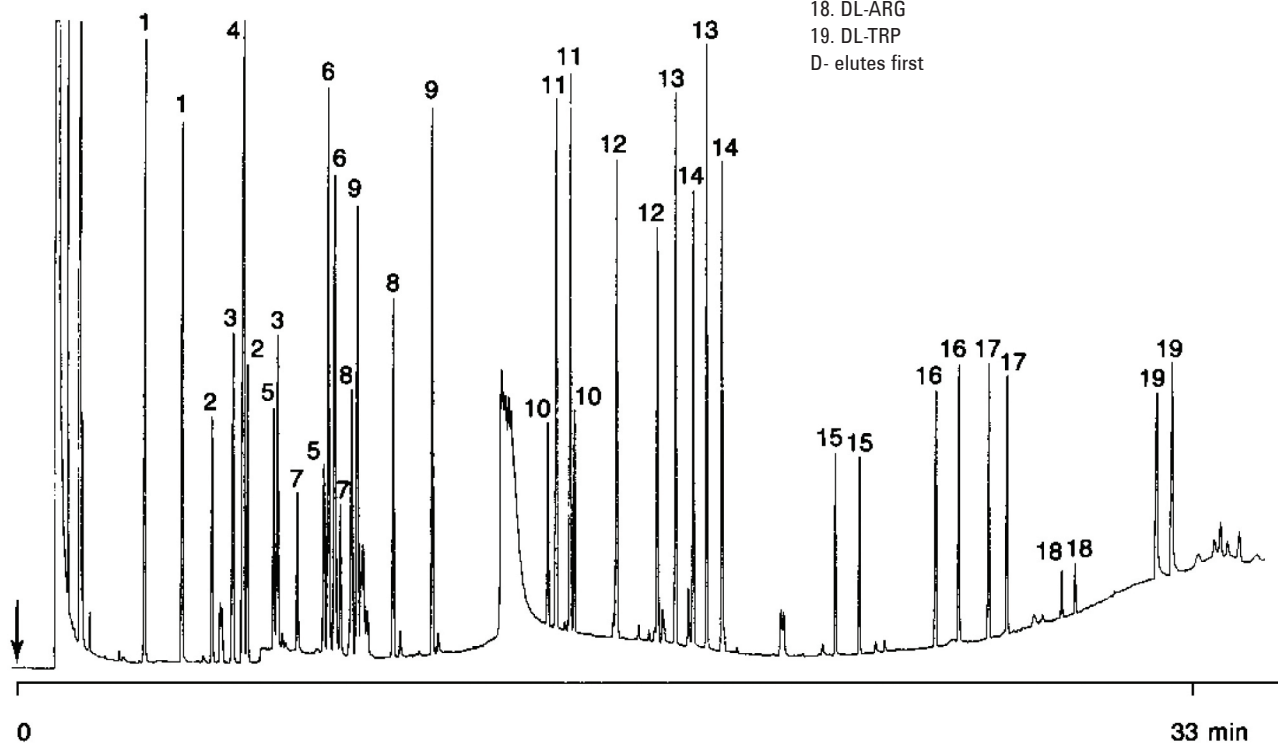
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Chirasil Val, 0.22 mm x 25 m fused silica
WCOT Chirasil Val (0.12 μ m) (Part no. CP7495)
Temperature : 75 °C (3 min) \rightarrow 195 °C, 3 °C/min
Carrier Gas : H₂, 90 kPa (0.9 bar, 13 psi), 48 cm/s
Injector : Splitter, 30 mL/min
T = 220 °C
Detector : FID, 8/4 x 10⁻¹² Afs
T = 220 °C
Sample Size : 0.05 μ L
Concentration range : 1-2 ng/component

Peak identification

1. DL-ALA
2. DL-VAL
3. DL-THR
4. GLY
5. DL- α -ILE
6. DL-PRO
7. DL-ILE
8. DL-SER
9. DL-LEU
10. DL-CYS
11. DL-ASP
12. DL-MET
13. DL-PHE
14. DL-GLU
15. DL-TYR
16. DL-ORN
17. DL-LYS
18. DL-ARG
19. DL-TRP
D- elutes first



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

A00263



Agilent Technologies