GC Analysis of Derivatized Amino Acids

Anila I Khan, Thermo Fisher Scientific, Runcorn, Cheshire, UK

Key Words

Derivatization reagent, silylation, N-methyl-N-(trimethylsilyl) trifluoroacetamide (MSTFA), L-alanine, L-leucine, L-lysine, amino acid

Abstract

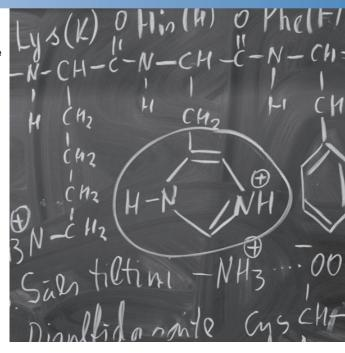
The derivatization of amino acids was achieved using the MSTFA silylation reagent. The TFA derivatized compounds were then analysed on a 5% phenyl methylpolysiloxane column using FID repetition of detection.

Introduction

Amino acids such as L-alanine, L-leucine and L-lysine are very difficult to analyse by GC as they decompose in the injector port and crash out on the column. The highly polar amino acid moiety, aids their solubility in water, but not in organic solvents which reduces the efficiency of their detection. To overcome these problems the amino acids can be derivatized to remove the active hydrogens and improve detectability. One of the most commonly used derivatization methods for the analysis of amino acids is a silylation reaction.

The Thermo Scientific Silylation reagent, N-methyl- N –(trimethylsilyl)trifluoroacetamide (MSTFA) is commonly used for derivatizing amino acids. The main advantage of using MSTFA is the formation of very volatile by-products such as TMS-amide, whilst the reagent itself elutes with the solvent front. MSTFA is manufactured to meet the exacting need for sensitive derivatization reactions. This reagent protects the functional groups present in the molecule aiding separation on the chromatography column. MSTFA involves converting the active hydrogen into a stable and volatile derivative which makes it easier for detection by FID.

In order to achieve separation and highly symmetrical peak shapes of these derivatized amino acids, a Thermo Scientific TRACE TR-5 column was used.



Experimental Details

Sample Preparation

2.5 mg each of L-alanine, L-leucine and L-lysine were weighed into a Thermo Scientific Reacti-Vial containing a Reacti-Vial magnetic stirrer. To the Reacti-Vial, 0.5 mL of MSTFA was added followed by 1 mL of acetonitrile. The Reacti-Vials were capped and placed in the Reacti-Therm Sample Incubation System and stirred for 20 minutes at 70 °C. Once dissolved the reaction was complete. The final sample was then transferred to a 2 mL autosampler vial and 1 μ L was injected into the GC/FID.



Reagents		Part Number
Thermo Scientific MSTFA 10 × 1 mL ampules		TS-48910
Thermo Scientific acetonitrile silylation grade solvent		TS-20062
Sample Handling Equipment		Part Number
Thermo Scientific Reacti-Therm III Heating/Stirring Module		TS-18823
Thermo Scientific Reacti-Vap III Evaporator		TS-18826
Thermo Scientific Reacti-Block Q-1 (Holds 8 × 10 mL Reacti-Vials)		TS-18814
Thermo Scientific Reacti-Vial reaction cl	ear glass vials 10 mL	TS-13225
Thermo Scientific 2 mL amber vial and s	screw tops	60180-565
Separation Conditions		Part Number
Instrumentation:	Thermo Scientific TRACE GC Ultra	
Column:	TRACE TR-5 30 m \times 0.25 mm \times 0.25 μ m	260E142P
Thermo Scientific BTO 17 mm septa		31303211
5 mm ID focus split liner, 105 mm long		453T1905
Graphite liner seal		29033406
10 μL, 50 mm needle length gauge 25 syringe		36500525
Graphite ferrules to fit 0.1-0.25 mm ID columns		29053488
Carrier gas:	Helium	
Split flow:	60 mL/min	
Column flow:	1.2 mL/min, Constant flow	
Split ratio:	50:1	
Oven temperature:	100 °C, 15 °C/min, 300 °C	
Injector type:	Split/Splitless	
Injector mode:	Split	
Injector temperature:	240 °C	
Detector type:	FID	
Detector temperature:	280 °C	
Detector Hydrogen flow:	35 mL/min	
Detector Air flow:	350 mL/min	

Thermo Scientific TriPlus Autosampler	
Injection Volume:	1 μL

30 mL/min

Detector Nitrogen flow:

Data Processing	
Software:	Thermo Scientific XCalibur

Results

Separation of derivatized amino acids was achieved using a 5% phenyl methylpolysiloxane (TRACE TR-5) column (Figure 1). The stability of the amino acids and enhanced detection is observed following derivatization with MSTFA.

Conclusion

MSTFA increases the volatility and stability of amino acids, allowing for enhanced separation and detection using the TRACE TR-5 GC column.

References

Thermo Scientific reagents, solvents and accessories brochure (Ref: BR20535_E 06/12S). Available upon request.

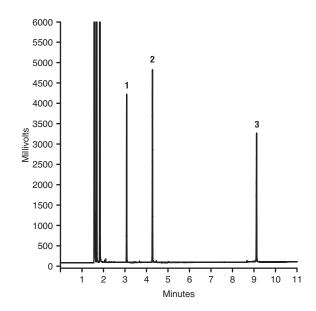


Figure 1: Chromatogram of the separation of the derivatized amino acids

Peak Number	Derivatised Amino acid	t _R (min)
1	L-alanine	3.1
2	L-leucine	4.3
3	L-lysine	9.1

thermoscientific.com/chromatography

©2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

USA and Canada +1 800 332 3331 Australia 1300 735 292 (free call domestic) China 800 810 5118 (free call domestic) 400 650 5118

France +33 (0)1 60 92 48 34 Germany +49 (0) 2423 9431 20 or 21 India +91 22 6742 9494 +91 27 1766 2352
 Japan
 0120 753 670 (free call domestic)

 0120 753 671 (fax)

 Korea
 +82 2 3420 8600

 United Kingdom
 +44 (0) 1928 534 110

United Kingdom +44 (0) 1928 534 110 New Zealand 0800 933 966 (free call domestic) Singapore +65 6289 1190

All Other Enquiries +44 (0) 1928 534 050



For advice and support, please visit our website: www.thermoscientific.com/chromexpert



A Thermo Fisher Scientific Brand