# GC Analysis of Acylated Sugars

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

## **Key Words**

Derivatization reagent, acylation, N-methyl-bis(trifluoroacetamide) (MBTFA) glucose, fructose, sugars, TR-1701

## **Abstract**

Sugars must be derivatized to a volatile form so as to be analyzed by GC. A commonly used derivatization reagent, N-Methyl-bis (trifluoroacetamide) (MBTFA) was used for converting sugars to their volatile forms. In order to achieve separation of TFA sugar derivatives such as fructose and glucose anomers, a mid-polar 14% cyanopropylphenyl polysiloxane phase GC column was used.

### Introduction

Sugars such as glucose and fructose are very difficult to analyse by GC, as they decompose in the injector port and "crash" out on the column. The highly polar and involatile nature of the sugars reduces the efficiency of the detection of these molecules. To overcome these problems the sugars can be derivatized to remove the active hydrogens such as -OH, therefore increasing volatility and improving detectability. The most commonly used derivatization method for the analysis of sugars is an acylation reaction. The Thermo Scientific acylation reagent MBTFA is used for derivatizing sugars and it is manufactured to meet the exacting need of sensitive derivatization reactions. This involves converting the active hydrogen into trifluroesters via a carboxylic derivative. The ester in the derivatized sugar improves the volatility, which makes it easier for analysis by GC/FID MBTFA, like the majority of derivatization reagents, produces a by-product. In this case the formation of the byproduct N-methyltrifluoroacetamide does not interfere with the analysis as it elutes earlier in the chromatogram.

In order to achieve separation of fructose and glucose anomers which arise upon derivatization, a mid-polarity 14% cyanophenyl polysiloxane Thermo Scientific TRACE TR-1701 column was used.



## **Experimental Details**

# Sample Preparation

5 mg each of glucose and fructose were weighed into a Thermo Scientific Reacti-Vial containing a Reacti-Vial magnetic stirrer. To the Reacti-Vial, 0.5 mL of MBTFA was added followed by 0.5 mL of Thermo Scientific silylation grade solvent pyridine. The Reacti-Vials were then capped and placed in the Thermo Scientific Reacti-Therm Sample Incubation System and stirred for 1 hour at 65 °C. Once dissolved the reaction was complete. The final sample was then transferred to a 2 mL autosampler vial and 1  $\mu$ L was injected into the GC/FID.



Reagents		Part Number
Thermo Scientific MBTFA 10 × 1 mL ampules		TS-49700
Thermo Scientific pyridine silylation g	TS-27530	
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 clear glass reaction vials 10 mL		TS-13225
Thermo Scientific 2 mL amber vial and screw tops		60180-565
Separation Conditions		Part Number
Instrumentation:	Thermo Scientific TRACE GC Ultra	
Column:	TRACE™ TR-1701 30 m × 0.25 mm × 0.25 μm	260Q142P
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:	40 °C (1 min), 10 °C/min, 260 °C (5 min)	
Injector type:	Split/Splitless	
Injector mode:	Split	
Injector temperature:	200 °C	
Detector type:	FID	
Detector temperature:	250 °C	
Detector Hydrogen flow:	35 mL/min	
Detector Air flow:	350 mL/min	
Datastar Nitragan flaur	00 ml /min	

Detector Nitrogen flow:	30 mL/min	
Thermo Scientific TriPlus Autosampler		
Injection Volume:	1 μL	
Data Processing		
Software:	Thermo Scientific XCalibur	

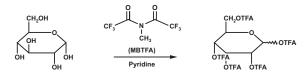


Figure 1: Acylation of glucose with MBTFA and pyridine

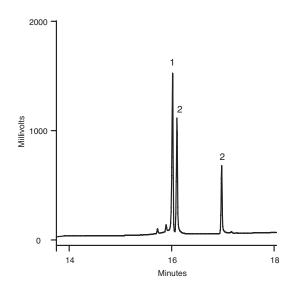


Figure 2: Chromatogram of the separation of derivatized sugars on a TR-1701 column.

Peak Number	Sugar	t <sub>R</sub> (min)
1	Fructose	16.0
2	Glucose anomers	16.1 and 17.0

## Results

The derivatized glucose produced two peaks, corresponding to the two cyclic forms of glucose existing as anomers (Figure 1) and fructose gave rise to one peak. Good baseline separation between fructose and the two anomers of derivatized glucose was observed (Figure 2) using a 14% cyanophenyl polysiloxane phase column. The stability of the sugars is improved as the acylation reagent protects the unstable groups, aiding separation on the chromatographic column.

## Conclusion

MBTFA is an ideal derivatization reagent for increasing the volatility of sugars. This enabled enhanced separation and detection of fructose and glucose anomers using a TRACE TR-1701 GC column.

## References

Thermo Scientific reagents, solvents and accessories brochure (Ref: BR20535\_E06/12). Available upon request. ACD labs software to draw chemical structures.

#### 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

**Technical Support** 



A Thermo Fisher Scientific Brand