

Fast Analysis of Coffee Bean Extracts Using a Solid Core HPLC Column

William Mullen¹, Joanna Freeke², Valeria Barattini²,

¹University of Glasgow, Glasgow, UK; ²Thermo Fisher Scientific, Runcorn, Cheshire, UK

Key Words

Accucore RP-MS, fast method, coffee, phenols, polyphenols, chlorogenic acids, caffeic acid, p-coumaric acid, ferulic acid

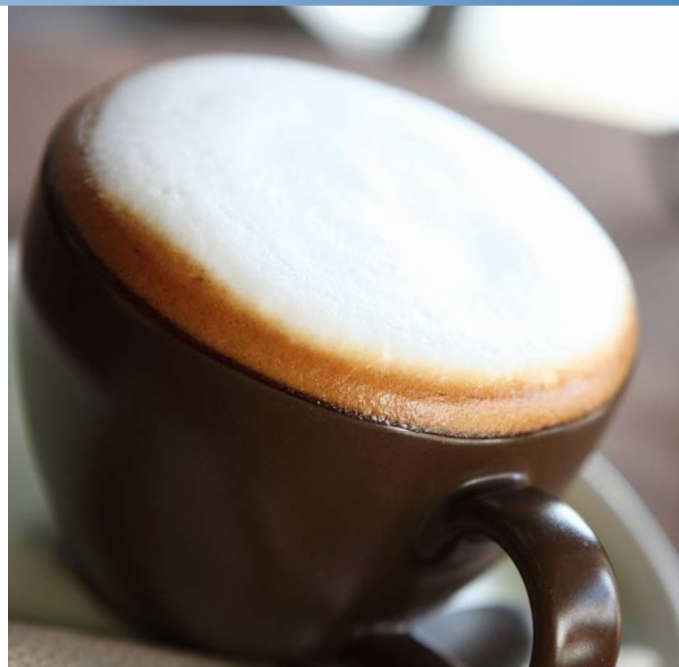
Abstract

This application note demonstrates the benefits of using Thermo Scientific™ Accucore™ RP-MS HPLC columns for the analysis of coffee extracts. By transferring the established method to the Accucore phase a 75% reduction in analysis time was achieved while maintaining full resolution of all components.

Introduction

It is believed that dietary phenols and polyphenols have the potential to protect the consumer's health and as such there has been a great deal of research into plant products that contain these compounds. One rich source of these antioxidant compounds is coffee, where a range of chlorogenic acids is found. For people who consume coffee beverages regularly, whether prepared from soluble powders or freshly brewed, daily intakes of 500 mg or even more of chlorogenic acids are easily achieved [1]. Chlorogenic acids are a family of esters formed between quinic acid and certain trans-cinnamic acids, most commonly caffeic, p-coumaric and ferulic. As many as 45 of these chlorogenic acids have been reported in coffee [2]. Thus quantification of individual compounds by absorbance detection requires good chromatographic separation.

Accucore HPLC columns use Core Enhanced Technology™ to facilitate fast and highly efficient separations. The 2.6 µm diameter particles are not totally porous, but instead have a solid core and a porous outer layer. The optimized phase bonding creates a series of high-coverage, robust phases. The tightly controlled 2.6 µm diameter of the Accucore particles results in much lower backpressures than typically seen with sub-2 µm materials. In this application an Accucore HPLC column was used to speed up the separation of a range of chlorogenic acids from a coffee extract.



Experimental Details

Consumables	Part Number
Fisher Scientific™ LC-MS grade water	W/0112/17
Fisher Scientific Pierce LC-MS grade acetonitrile	TS-511001
Fisher Scientific Pierce LC-MS grade formic acid	PI-28905
9 mm screw thread vial, 300 µL, fused insert	03-FISV
9 mm open top short screw cap, 6 mm hole, blue silicone/PTFE septum	9-SC(B)-ST101

Separation Conditions		Part Number
Instrumentation:	Thermo Scientific Accela HPLC system	
Column 1:	Fully porous C18, 4 µm, 250 x 2.0 mm	
Column 2:	Accucore RP-MS, 2.6 µm, 150 x 3.0 mm	17626-153030
Mobile phase A:	0.1% formic acid in water	
Mobile phase B:	0.1% formic acid in acetonitrile	
Gradient (column 1)	Time (min)	% B
	0	5
	20	10
	60	35
Flow rate:	200 µL/min	
Gradient (column 2)	Time (min)	% B
	0	2
	8	7
	15	50
Flow rate:	700 µL/min	
Backpressure at starting conditions for both columns:	Approx. 200 bar	
Column temperature:	40 °C	
Injection details:	5 µL	
Detection details:	UV absorbance detection at 325 nm	

Sample Preparation

Soluble coffee extract (Futureceuticals Inc. USA) in water/methanol (90:10 v/v)

Data Processing

Software: Thermo Scientific Xcalibur Software v1.3

Results

Current methods for the measurement of chlorogenic acid content of commercial coffee bean extracts [3] typically require a gradient method with an analysis time of 60 minutes. Such methods use a 250 x 2.0 mm C18 column packed with conventional 4 µm particles (Figure 1) to achieve adequate separation. Here the first analysis run of a similar coffee bean extract using a 150 x 3.0 mm Accucore RP-MS column is reported. The method was modified to take advantage of the very high performance offered by Accucore columns and reduce the run time. An equivalent separation at a similar backpressure was achieved with an analysis time of only 15 minutes. These excellent results, achieved from the first analysis, meant that no further method development was undertaken (Figure 2).

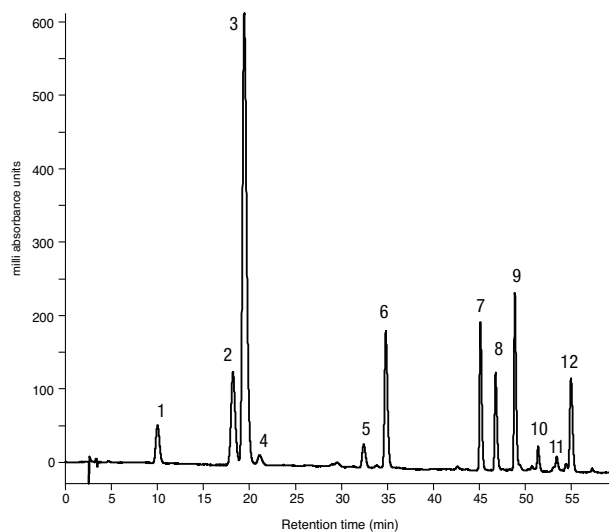


Figure 1: Analysis of a coffee bean extract using a 60 minute gradient on a fully porous C18 HPLC column (4 μm particle size, 250 x 2 mm)

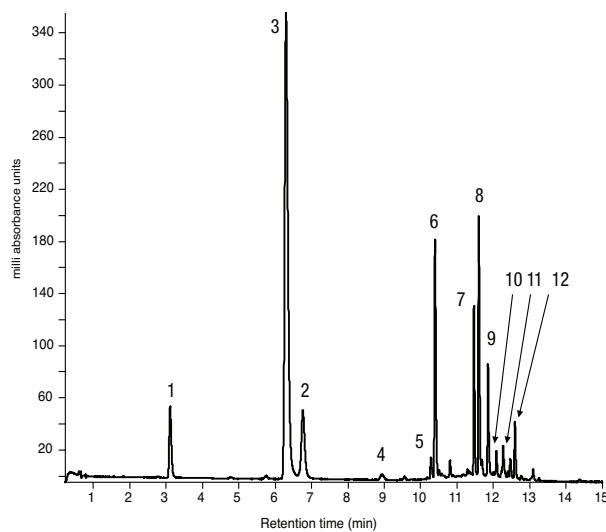


Figure 2: Analysis of a coffee bean extract using a 15 minute gradient on a solid core Accucore RP-MS HPLC column (2.6 μm particle size, 150 x 3.0 mm)

Along with an improvement in the resolution between peaks 2 and 3, the Accucore column also exhibits a change in selectivity, and compound 3 elutes before compound 2.

Peak number	Compound	Peak number	Compound
1	3-O-Caffeoylquinic acid	7	3,4-O-Dicaffeoylquinic acid
2	4-O-Caffeoylquinic acid	8	3,5-O-Dicaffeoylquinic acid
3	5-O-Caffeoylquinic acid	9	4,5-O-Dicaffeoylquinic acid
4	3-O-Feruloylquinic acid	10	3-O-Feruloyl-4-O-caffeoylquinic acid
5	4-O-Feruloylquinic acid	11	3-O-Caffeoyl-5-O-feruloylquinic acid
6	5-O-Feruloylquinic acid	12	4-O-Caffeoyl-5-O-feruloylquinic acid

Table 1: Peak numbers and identities of compounds in coffee extracts

Conclusion

- The Accucore RP-MS HPLC column is suitable for analyzing coffee extracts and quantifying the amount of different chlorogenic acids.
- Transferring the established analysis method to an Accucore RP-MS HPLC column enabled a 75% reduction in analysis time (60 minutes down to 15 minutes).
- Good resolution of all components was observed without additional method development.
- Backpressure was suitable for use on a conventional HPLC systems.

References

1. Clifford, M. N., Chlorogenic acids and other cinnamates - nature, occurrence, dietary burden, absorption and metabolism. *J. Sci. Food Agric.* 2000, 80 (7), 1033-1043.
2. Clifford, M. N.; Johnston, K. L.; Knight, S.; Kuhnert, N., Hierarchical scheme for LC-MSn identification of chlorogenic acids. *J. Agric. Food Chem.* 2003, 5(10), 2900-11.
3. Mullen, W.; Nemzer, B.; Ou, B.; Stalmach, A.; Hunter, J.; Clifford, M. N.; Combet, E., The antioxidant and chlorogenic acid profiles of whole coffee fruits are influenced by the extraction procedures. *J. Agric. Food Chem.* 2011, 59(8), 3754-62.

thermoscientific.com/chromatography

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific Inc. 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
France +33 (0)1 60 92 48 34
Germany +49 (0) 2423 9431 20 or 21
United Kingdom +44 (0)1928 534110
Japan +81 3 5826 1615

China +86 21 68654588 +86 10 84193588
 +86 20 83145199 800 810 5118
India +91 22 6742 9494 +91 27 1766 2352
Australia 1 300 735 292 (free call domestic)
New Zealand 0800 933 966 (free call domestic)
All Other Enquiries +44 (0) 1928 534 050

Technical Support
North America +1 800 332 3331
Outside North America +44 (0) 1928 534 440

Thermo
 SCIENTIFIC

Part of Thermo Fisher Scientific