High-Throughput Analysis of Dextromethorphan and Dextrorphan

Kimberly Phipps, Thermo Fisher Scientific, Runcorn, Cheshire, UK

Key Words

Accucore C18, dextromethorphan, dextrorphan

Abstract

A liquid chromatography method for the analysis of dextromethorphan and dextrorphan was developed on a Thermo Scientific™ Accucore™ C18 2.6 µm, 50 x 2.1 mm HPLC column for a fast separation with a cycle time of 2 minutes and excellent peak shape and resolution.

Introduction

Dextromethorphan (Figure 1) is used as a cough suppressant and can be found in many over-the-counter cough and cold remedies. Dextromethorphan is quickly metabolized to dextrorphan (Figure 2) in the liver. This process can be used to assess the speed of metabolism. The retention, peak shape, and resolution of dextromethorphan and dextrorphan are demonstrated in this application.

Thermo Scientific Accucore HPLC columns use Core Enhanced Technology TM to facilitate fast and highly efficient separations. The 2.6 μ m 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 carbon loading of the Accucore C18 HPLC column provides high retention of non-polar analytes via a predominantly hydrophobic interaction mechanism. The tightly controlled 2.6 μ m diameter of Accucore particles results in much lower backpressures than typically seen with sub-2 μ m materials.

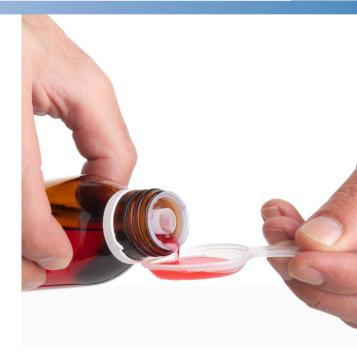


Figure 1. Dextromethorphan

Figure 2. Dextrorphan



Experimental Details

Sample Handling	Part Number
Fisher Scientific™ HPLC grade water	W/0106/17
Fisher Scientific HPLC grade acetonitrile	A/0626/17
Fisher Scientific HPLC grade ortho phosphoric acid (OPA)	J/5655/PB08
Thermo Scientific™ National™ Mass Spec Target DP Certified 2 mL clear vial with ID patch and blue bonded cap with PTFE/Silicone septum	MSCERT4000-34W

Sample Preparation

A primary standard of dextromethorphan and dextrorphan was prepared in methanol at 1 mg/mL

The working standard contained 1 µg/mL dextromethorphan and dextrorphan in 80:20 (v/v) water / acetonitrile

Separation Conditions			Part Number
Instrumentation:	Thermo Scientific Dionex™ UltiMate™ 3000 HPLC system		
Column:	Accucore C18 2.6 μm, 50 x 2.1 mm		17126-052130
Mobile phase A:	Water + 0.1% OPA		
Mobile phase B:	Acetonitrile + 0.1% OPA		
Gradient:	Time (min)	%B	
	0	5	
	1	95	
	1.01	5	
	2	5	
Flow rate:	1.4 mL/min		
Column temperature:	40 °C		
Pressure:	360 Bar		
Injection details:	1 μL		
Injection wash solvent:	Water / acetonitrile (60:40 v/v)		
Wavelength:	194 nm		

Results

This application was run at a high flow rate to provide a fast separation suitable for high-throughput laboratories. The Thermo Scientific Accucore C18 HPLC column gave excellent peak shape and resolution, as demonstrated in Figure 3. The calculated time of T_0 for this method is 7.43 seconds or 0.12 minutes. This can be clearly seen on the chromatogram and provides dextrorphan with a k' of 6.8, proving the peaks are well retained even with a short runtime.

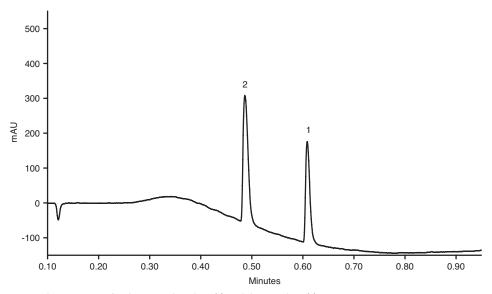


Figure 3: Chromatogram for dextromethorphan (1) and dextrorphan (2)

Replicate injections of the standard containing dextromethorphan and dextrorphan showed that the Accucore C18 HPLC column produced stable and reproducible results (Table 1).

	Dextromethorphan	Dextrorphan
Retention time (minutes)	0.60	0.48
% RSD on retention time	1.05	1.57
Asymmetry	1.43	1.46
Resolution	-	7.58

Table 1: Results obtained from an Accucore C18 column, based upon 6 replicate injections

Conclusion

- Analysis of dextromethorphan and dextrorphan can be achieved using an Accucore C18 HPLC column.
- The Accucore C18 HPLC column provides good resolution for dextromethorphan and dextrorphan in less than 2 min.
- The Accucore C18 HPLC column gives excellent peak shape for dextromethorphan and dextrorphan.
- The Accucore C18 HPLC column performs reproducibly for the analysis of dextromethorphan and dextrorphan.

For Research Use Only.

thermoscientific.com/accucore

© 2013 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

