

# HPLC-UV Method for the Determination of Alkaloids Using a Synchronis aQ Column

Prakash Chander, Tushar N Mehta, Centre of Excellence for Asia Pacific Lab,  
Thermo Fisher Scientific, Ahmedabad, India

## Key Words

Nicotine, anabasine, cotinine, synchronis aQ column, stimulants in mammals, alkaloid

## Abstract

This application note demonstrates the use of the Thermo Scientific Synchronis aQ Column for the determination of alkaloids by HPLC-UV.

## Introduction

The sensitive and specific detection of nicotine, its metabolites and the tobacco alkaloid anabasine is useful in evaluating the success of smoking cessation treatments and detecting tobacco use, passive exposure and nontobacco nicotine exposure in potential transplant and elective surgical patients.

One of the key goals for the chromatographer is to achieve a consistent, reproducible separation. The selection of a highly reproducible HPLC column is essential if this goal is to be attained. The Synchronis™ column range has been engineered to provide exceptional reproducibility due to its highly pure, high surface area silica, dense bonding and double endcapping, all controlled and characterized through the use of rigorous testing.

This application note demonstrates the successful retention and separation of nicotine, anabasine and cotinine using Synchronis aQ 5 µm column. The polar endcapping used in Synchronis aQ provides a controlled mechanism for retention of these polar compounds.



## Experimental Details

Consumables	Part Number
Fisher Scientific HPLC grade water	W/0106/17
Fisher Scientific HPLC grade ammonium acetate	A/3446/50
Fisher Scientific HPLC grade acetonitrile	A/0626/17
Alkaloids purchased from Sigma Aldrich	

Sample Handling Equipment	Part Number
NSC Mass Spec Certified 2 mL clear vial with blue bonded PTFE silicone cap	MSCERT4000-34W

Separation Conditions	Part Number	
Instrumentation:	Thermo Scientific Accela UHPLC system	
Column:	Synchronis aQ 5 $\mu$ m, 150 x 2.1 mm	97305-152130
Mobile phase:	90:10 (v/v) 20 mM ammonium acetate/acetonitrile	
Flow rate:	1.0 mL/min	
Column temperature:	40 °C	
Injection details:	2 $\mu$ L partial loop	
Injection wash solvent:	90:10 (v/v) 20 mM ammonium acetate/acetonitrile	
UV detector wavelength:	260 nm	
Backpressure:	Approximately 220 bar	

## Solutions

Working standard contained 100  $\mu$ g/mL of nicotine, anabasine and cotinine in mobile phase.

## Results

The analysis was performed on a Synchronis aQ 5  $\mu$ m, 150 x 2.1 mm column. As shown in Figure 1, nicotine, anabasine and cotinine were analyzed in less than 5 minutes. Table 1 shows the results from six replicate injections.

	Anabasine	Cotinine	Nicotine
Retention time (minutes)	0.81	1.16	1.41
%RSD on retention time	0.00	0.35	0.39
Area	197049	169593	130437
%RSD on area	0.41	0.23	0.51

Table 1: Retention time and area results for anabasine, cotinine and nicotine

## Conclusion

Replicate injections of alkaloids showed that Synchronis aQ produced stable and reproducible results. This demonstrates that Synchronis aQ is an excellent choice of column for the rapid analysis of these nicotine related alkaloids.

## References

Specific detection of Anabasine, Nicotine, and Nicotine Metabolites in Urine by Liquid Chromatography-Tandem Mass Spectrometry.

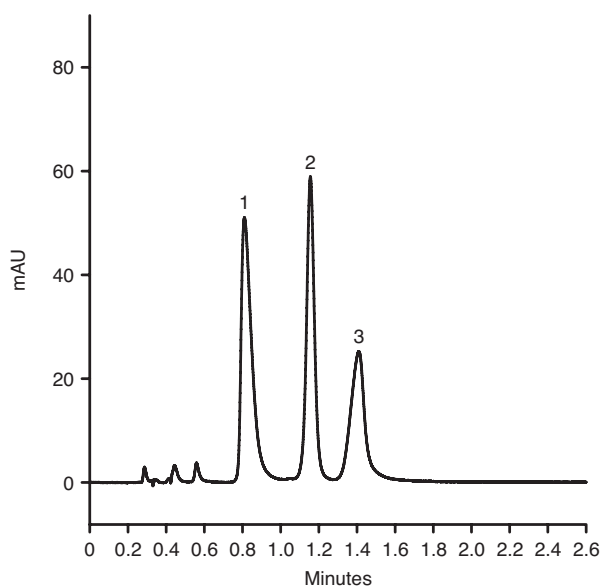


Figure 1: Chromatogram of alkaloids analyzed using a Synchronis aQ 5  $\mu$ m, 150 x 2.1 mm column 1. Anabasine 2. Cotinine 3. Nicotine

[thermoscientific.com/chromatography](http://thermoscientific.com/chromatography)

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. 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