

A Fully Automated LC-MS Screening System using Automated Online Sample Preparation for Forensic Toxicology

Mueller, D.M.¹; Duretz, B.²; Espourteille, F.A.³; Rentsch, K.M.¹

¹Institute for Clinical Chemistry, University Hospital, Zurich, Switzerland; ²Thermo Fisher Scientific, Les Ulis, France;

³Thermo Fisher Scientific, Franklin, USA

Introduction

Liquid chromatography-mass spectrometry (LC-MS) is a powerful tool widely used for forensic targeted drug screening. However, the quality of the results is highly affected by the sample preparation. Offline solid phase extraction (SPE) and liquid-liquid extraction (LLE) are widely used, but these methods are often time-consuming and costly. To provide a fast and sensitive approach, an automated online sample preparation method using Thermo Scientific Transcend TLX-1 system powered by TurboFlow™ technology for the forensic toxicological screening of more than 400 acidic, neutral, and basic drugs in urine with LC/MSⁿ has been developed.

Goal

To evaluate the performance of an automated online sample preparation method for an LC/MSⁿ screening approach.

Experimental

Sample preparation was performed by an online sample extraction method utilizing Thermo Scientific TurboFlow technology. Two TurboFlow columns (Cyclone, C18XL) were connected in series and used for sample extraction. Urine samples were run both natively and after enzymatic hydrolysis. The eluent was then transferred to the LC column (Thermo Scientific Betasil Phenyl-Hexyl, 100 x 3 mm, 3 μm) for separation.

A 30-minute gradient from 1% to 98% organic was employed for separation of the analyte with flow rates of 300 μL/min. All samples were then analyzed on a Thermo Scientific LXQ linear ion trap mass spectrometer with the atmospheric pressure chemical ionization (APCI) source. A data-dependent polarity switching method was used for data acquisition. MS² and MS³ spectra were acquired. Since polarity switching was used, a single injection of a sample containing unknown compounds was sufficient to detect both substances ionizing in negative and positive mode. The data was automatically processed, post-acquisition, by Thermo Scientific ToxID automated screening software.

Results and Discussion

The method using online extraction has been fully validated. A minor matrix effect (suppression < 5%) was observed for over 98% of the compounds. A recovery of more than 90% was seen in 90% of the substances. The limit of identification (LOI) was below 10 ng/mL for 60% of the substances and 90% could be identified at a concentration of 100 ng/mL. The 400-compound library contains both MS² and MS³ spectra. MS³ spectra bring an additional level of specificity, although in most cases, the analytes can be easily identified by using only the MS² spectra. However, some analytes may have the same molecular weight, very similar MS² spectra, and a very close retention time. For these reasons, MS³ data have to be used for the identification. One example is the isobaric compounds O-desmethylvenlafaxine and tramadol. The two analytes have the same molecular weight, very close retention times (see details in Table 1), and the same MS² spectra (Figure 1). Therefore, by running only MS² experiments, it is impossible to properly differentiate the two analytes. When MS³ spectra are recorded, tramadol does not fragment ions while O-desmethylvenlafaxine gives a specific spectrum (Figure 1). Therefore, the analytes can be properly identified. Total run time of the analysis is 30 minutes. An example of a chromatogram obtained from a sample is presented in Figure 2.

Table 1. Tramadol and O-desmethylvenlafaxine information

	O-Desmethylvenlafaxine	Tramadol
Precursor mass	264.3	264.3
MS ² Fragment	246.3	246.3
Retention Time	10.6 min	10.3 min

Key Words

- Transcend TLX-1 system
- TurboFlow Technology
- LXQ Linear Ion Trap
- Illicit Drugs
- Forensic Toxicology

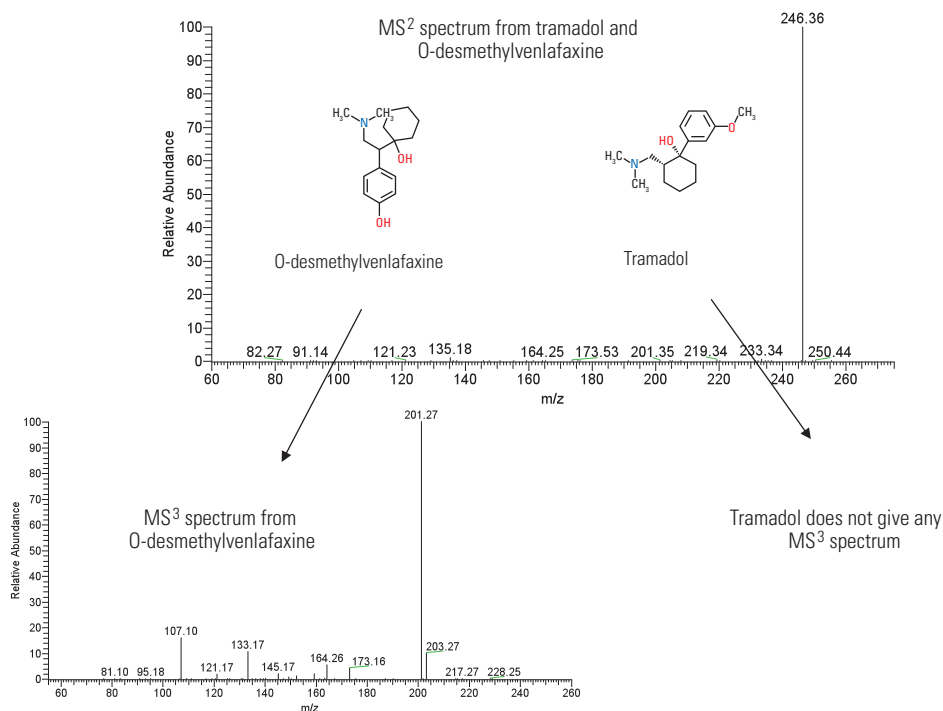


Figure 1. Fragmentation of tramadol and O-desmethylenlafaxine in MS² and MS³

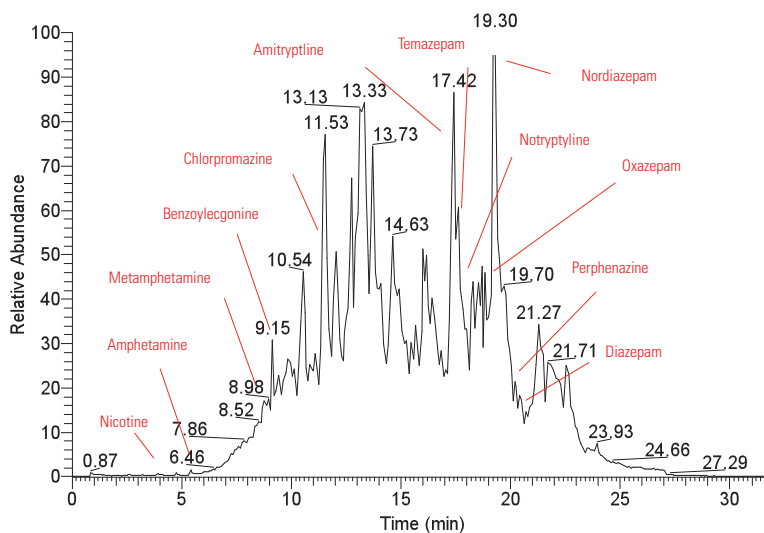


Figure 2. Full scan MS chromatogram of a sample containing 12 different analytes

Conclusion

The automated online TurboFlow method with the LXQ™ linear ion trap mass spectrometer allows a fast and specific approach for the identification of a broad range of compounds in positive and negative mode in a single run. The sample preparation time is 15 minutes

with this method as compared to 2 hours with an offline approach. The LOIs are below 100 ng/mL for more than 90% of the analytes. MS³ spectra acquisition brings an additional level of specificity for forensic toxicology laboratories.

www.thermofisher.com

Legal Notices: ©2016 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.

In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.

Africa-Other
+27 11 570 1840

Australia
+61 3 9757 4300

Austria
+43 1 333 50 34 0

Belgium
+32 53 73 42 41

Canada
+1 800 530 8447

China
+86 10 8419 3588

Denmark
+45 70 23 62 60

Europe-Other
+43 1 333 50 34 0

Finland/Norway/Sweden
+46 8 556 468 00

France
+33 1 60 92 48 00

Germany
+49 6103 408 1014

India
+91 22 6742 9434

Italy
+39 02 950 591

Japan
+81 45 453 9100

Latin America
+1 561 688 8700

Middle East
+43 1 333 50 34 0

Netherlands
+31 76 579 55 55

New Zealand
+64 9 980 6700

Russia/CIS
+43 1 333 50 34 0

South Africa
+27 11 570 1840

Spain
+34 914 845 965

Switzerland
+41 61 716 77 00

UK
+44 1442 233555

USA
+1 800 532 4752

AN63402_E_06/16S