

Determination of LSD and Its Metabolites in Human Biological Samples by Liquid Chromatography–Tandem Mass Spectrometry

François-Ludovic Sauvage¹, Pierre Marquet^{1,2}

¹Department of Pharmacology-Toxicology, University Hospital, Limoges, France.

²Laboratory of Pharmacology, Faculty of Medicine, University of Limoges, France.

Introduction

Lysergic acid diethylamide (LSD) is a very potent hallucinogenic drug involving, particularly, behavioral disorders and is also extensively metabolized in man. Moreover, LSD and its major metabolites are present at low concentration in biological fluids, such as whole blood or urine. Identification and quantitation of such compounds for forensic use necessitate a sensitive and specific method. This study aims to describe a method using liquid chromatography/tandem mass spectrometry and permitting to quantify LSD and its metabolites at low concentrations.

Goal

The goal of this study was to identify and quantify LSD, iso-LSD, nor-LSD, nor-iso-LSD and 2-oxo-3-hydroxy-LSD in biological matrices. This report demonstrates the use of the TSQ Quantum for this application.

Experimental Conditions/Methods

Chemicals and Reagents

Lysergic acid diethylamide (LSD), d₃-LSD (internal standard), 2-oxo-3-hydroxy-LSD, iso-LSD, nor-LSD were purchased from Cerilliant (Austin, TX, USA). Ammonium formate and formic acid (>99 % pure) were purchased from Sigma. All reagents and solvents used in the extraction procedures were of analytical grade.

Sample Preparation

To 2 mL of serum, urine or whole blood content were added 100 µL of a 0.025 µg/mL aqueous solution of d₃-LSD (Internal Standard), 1 mL of a solution of pH 9.50 carbonate buffer and 8 mL of dichloromethane-isopropanol (95:5 by volume). The tubes were vortex-mixed and shaken on an oscillatory mixer. After centrifugation at 3,400 g for 5 min, the organic phase was poured in a conical glass tube and evaporated under a stream of nitrogen at 37°C. The dried extracts were reconstituted in 25 µL of acetonitrile : pH 3.0, 2 mmol/L ammonium formate (30:70 by volume) and 10 µL were injected into the chromatographic system.

Instrumentation Methods

HPLC Conditions

The chromatographic system consisted of a Shimadzu 10ADvp micro-flow rate, high-pressure gradient pumping system with a Rheodyne® Model 7725 injection valve equipped with a 5 µL internal loop. A C18, 5 µm (50×2.1 mm) column, maintained at 25°C, was used with a linear gradient of mobile phase A (pH 3.0, 2 mmol/L ammonium formate) and mobile phase B (acetonitrile:pH 3.0, 2 mmol/L ammonium formate [90:10; v/v]), flow rate of 200 µL/min, programmed as follows: 0-1.5 min, 5% B; 1.5-9 min, 5 to 50% B; 9-10 min, 50 to 90% B; 10-10.5 min, decrease from 90 to 5% B; 10.5-13 min, equilibration with 5% B.

MS Conditions

Mass Spectrometer: Thermo Scientific TSQ Quantum
Source: ESI mode
Ion Polarity: Positive
Spray Voltage: 4000 V
Sheath/Auxiliary gas: Nitrogen
Sheath gas pressure: 25 (arbitrary units)
Auxiliary gas pressure: 15 (arbitrary units)
Ion transfer tube temperature: 250°C
Scan type: SRM
Collision gas: Argon
Collision gas pressure: 1.5 mTorr

SRM Conditions

Settings were optimized by infusing at 5 µL/min a 1 µg/L solution containing the studied compound in acetonitrile: pH 3.0, 2 mmol/L ammonium formate (30:70, by volume). The structure of these compounds is shown in Figure 1.

Key Words

- TSQ Quantum
- Drugs of Abuse
- Forensic analysis
- LC-MS/MS
- LSD (Lysergic acid diethylamide)
- LSD metabolites
- Toxicology

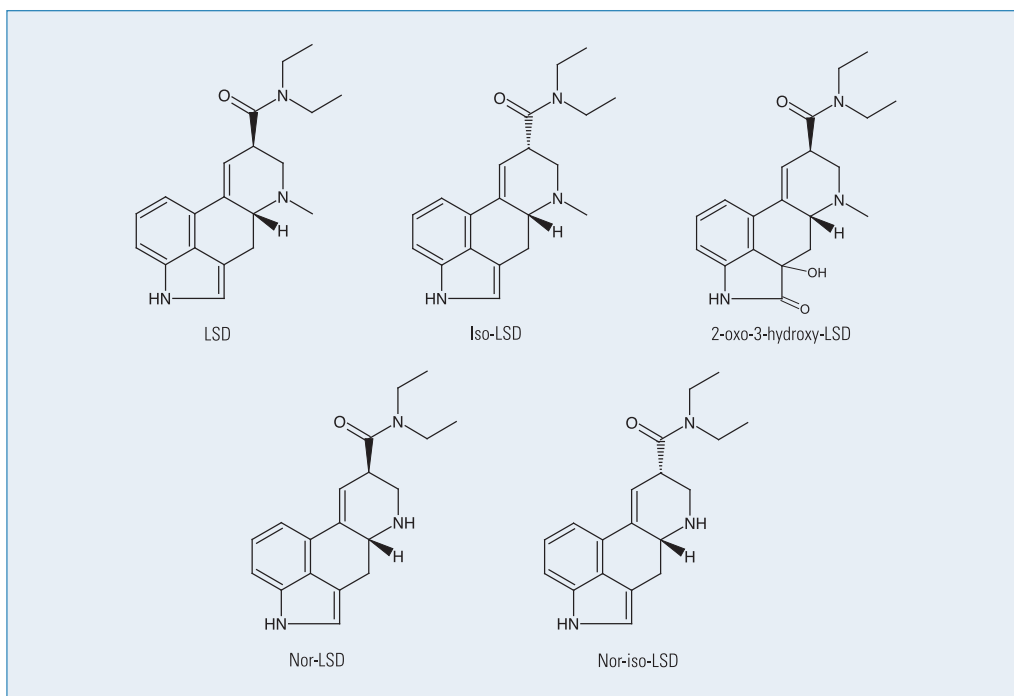


Figure 1: Structures of investigated compounds

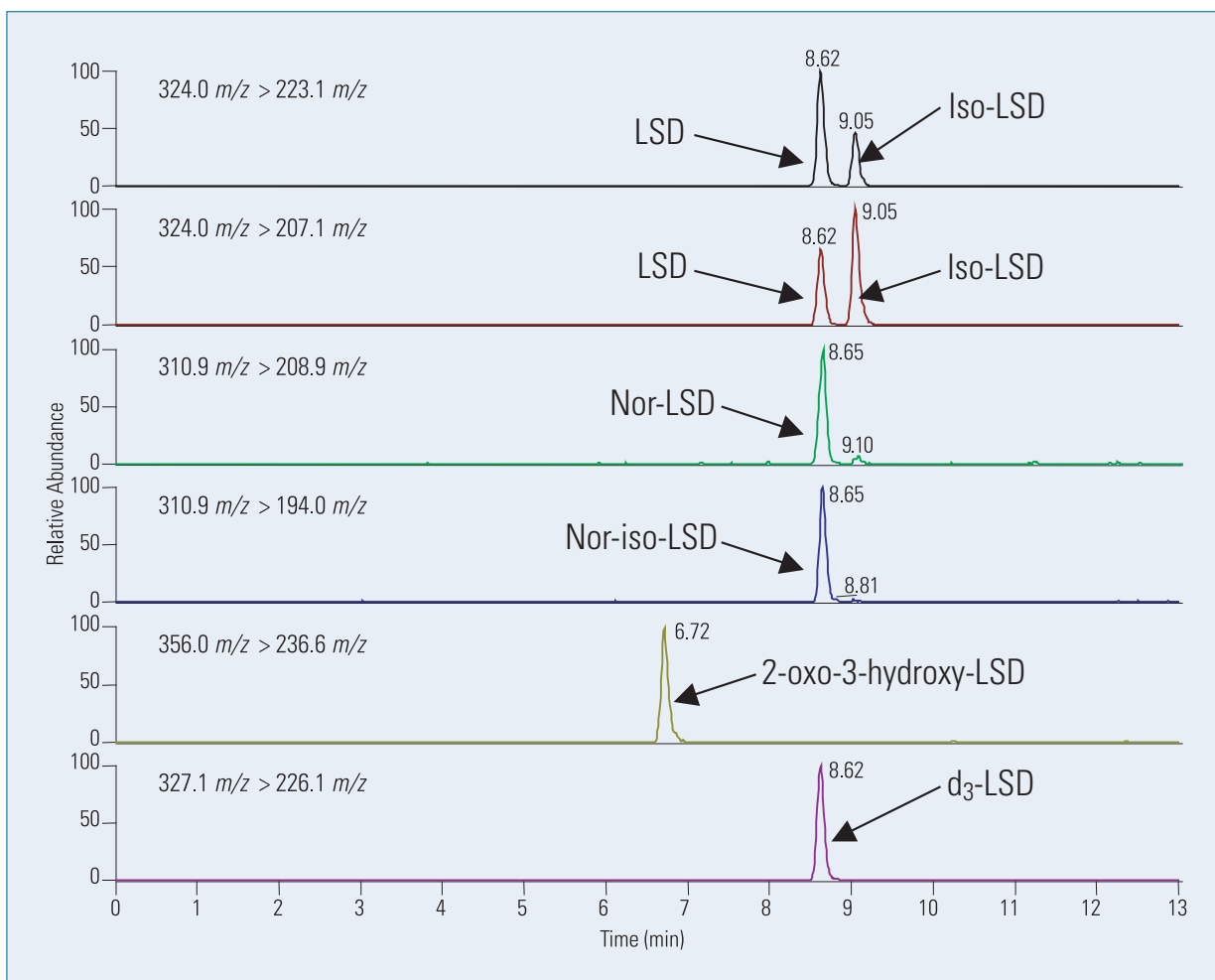


Figure 2: Chromatogram of a urine spiked at 0.5 ng/mL

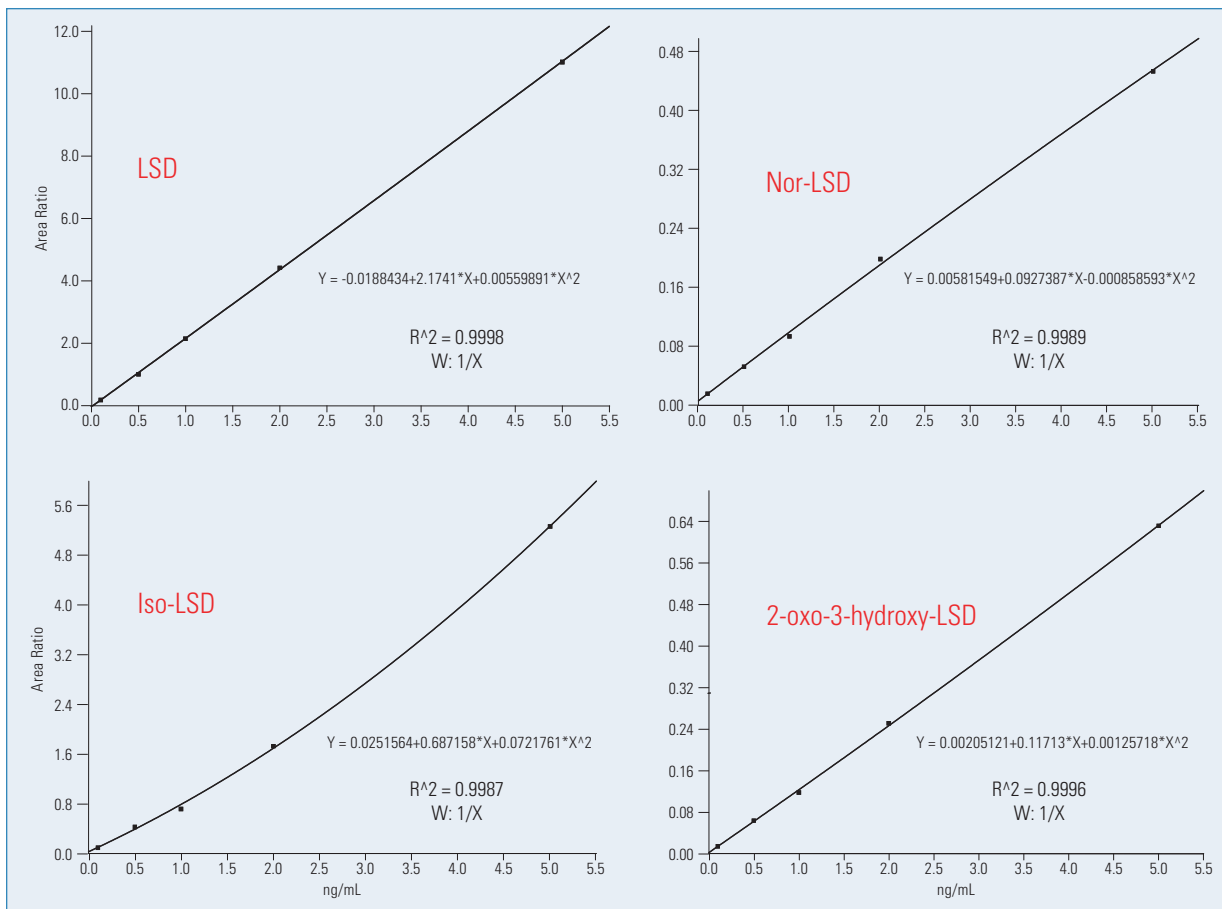


Figure 3: Representative calibration curves from standards spiked in urine

Compounds	Quantification transition	Collision energy	Confirmation transition	Tube lens voltage
LSD	324.0/223.1	30	324.01/207.1	50
Iso-LSD	324.0/223.1	30	324.01/207.1	50
Nor LSD	310.9/208.9	28	310.91/194.0	54
Nor-iso-LSD	310.9/208.9	28	310.91/194.0	54
2-oxo-3-hydroxy-LSD	356.0/236.6	30	356.01/222.0	36
d3-LSD	327.1/210.1	50	327.11/226.2	30

Results and Discussion

The LC-ESI/SRM chromatograms obtained for a blank urine spiked at 0.5 ng/mL are shown in Figure 2. As presented, LSD and iso-LSD are separated using the chromatographic conditions described previously. Identification of LSD is performed using two characteristic transitions and the retention time given by its deuterated internal standard.

Linearity

Calibration curves obtained for each compound spiked in urine samples are presented in Figure 3. Concentration ranges were comprised between 0.1 ng/mL and 5 ng/mL.

Conclusion

This application note described a sensitive, specific method developed for the quantitation of lysergide and metabolites in various biological matrices for forensic use.

Laboratory Solutions Backed by Worldwide Service and Support

Tap our expertise throughout the life of your instrument. Thermo Scientific Services extends its support throughout our worldwide network of highly trained and certified engineers who are experts in laboratory technologies and applications. Put our team of experts to work for you in a range of disciplines – from system installation, training and technical support, to complete asset management and regulatory compliance consulting. Improve your productivity and lower the cost of instrument ownership through our product support services. Maximize uptime while eliminating the uncontrollable cost of unplanned maintenance and repairs. When it's time to enhance your system, we also offer certified parts and a range of accessories and consumables suited to your application.

To learn more about our products and comprehensive service offerings, visit us at www.thermo.com.

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 2 8844 9500

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
+1 608 276 5659

Middle East
+43 1 333 50 34 0

Netherlands
+31 76 579 55 55

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

www.thermo.com

©2007-2009 Thermo Fisher Scientific Inc. All rights reserved. Rheodyne is a registered trademark of IDEX Health & Science group. All other 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.

View additional Thermo Scientific LC/MS application notes at: www.thermo.com/appnotes



Thermo Fisher Scientific,
San Jose, CA USA is ISO Certified.

AN62232_E 12/09S