

THC-COOH Quantification in Urine Using Dilute and Shoot LC-MS/MS Method for Forensic Toxicology

Samuele Scurati, Thermo Fisher Scientific, Rodano, Italy

Maura Brambilla, Ludovico D'Amato, Paolo Brambilla, University Department of Laboratory Medicine, Hospital of Desio, Desio, Italy

Introduction

Cannabis sativa is a widely used drug of abuse. Tetrahydrocannabinol (THC) is the major psychoactive chemical compound in the cannabis plant. After smoke inhalation, THC is absorbed and distributed in blood. Subsequently, it is rapidly metabolized to THC-COOH, conjugated with glucuronic acid, and excreted through urine. Liquid chromatography-tandem mass spectrometry (LC-MS/MS) is considered a useful tool to establish the consumption of cannabis by the assessment of THC-COOH in urine for forensic toxicology purposes.

Goal

To develop a reliable and fast analytical method for the quantitative determination of THC-COOH in urine using a Thermo Scientific TSQ Quantum Access MAX triple stage quadrupole mass spectrometer.

Experimental

Sample Preparation

A urine sample was hydrolyzed with 10M NaOH and heated at 60 °C for 15 minutes. The pH was restored with Fisher Chemical acetic acid. Hydrolyzed samples as well as calibrators were diluted 1:10 in Fisher Chemical water/acetonitrile (1:1). Then, 10 µL were directly injected. Quantitative analysis was performed on the basis of calibration curves prepared in urine, ranging from 7.8 to 1000 ng/mL. Calibrators were injected in duplicate.

UHPLC conditions

Liquid chromatography separation was performed using a Thermo Scientific Accela autosampler and pump. The sample was injected directly on a Thermo Scientific Hyperasil GOLD column (50 × 2.1 mm, 1.9 µm). A gradient LC method used mobile phases A (0.1% aqueous formic acid) and B (Fisher Chemical Optima LC/MS acetonitrile) at a flow rate of 300 µL/min. The run time was 6 minutes.

Mass Spectrometry

MS analysis was carried out on a TSQ Quantum Access MAX™ triple stage quadrupole mass spectrometer equipped with a Thermo Scientific Ion Max source with a heated electrospray ionization (HESI) probe. The MS conditions were as follows:

Scan type:	SRM
Divert valve:	2 - 4 min to source
Selected ions for quantification:	m/z 343 → 299 + 245 for THC-COOH in negative mode

Results and Discussion

Figures 1 and 2 show the ion chromatograms of the lowest and highest calibration points. Excellent linearity ($r^2 = 0.99$) fits for the calibration curve were observed over the range of 7.8-1000 ng/mL urine, with a Coefficient of Variation (%CV) at the lower end of 6.5%. The limit of quantitation (LOQ) was established as 7.8 ng/mL in urine.

Figure 4 reports an ion chromatogram of a real urine sample positive for cannabinoids (225 ng/mL urine), analyzed as described.

To examine the difference between hydrolyzed and non-hydrolyzed urine, we analyzed the same urine sample without the hydrolysis step. When urines were not hydrolyzed, the portion excreted as free THC-COOH was detected at 3.06 minutes, while THC-COOH-glucuronide was detected at 2.58 minutes (Figure 5). The precursor ion m/z 343 was generated as result of an in-source fragmentation and a consequent loss of glucuronic acid.

Because THC-COOH is mainly excreted as glucuronic acid conjugate, it is always necessary to perform urine hydrolysis before the LC-MS analysis to obtain an accurate quantification of THC-COOH.

Key Words

- TSQ Quantum Access MAX
- Accela Pump
- Cannabinoids
- Forensic Toxicology

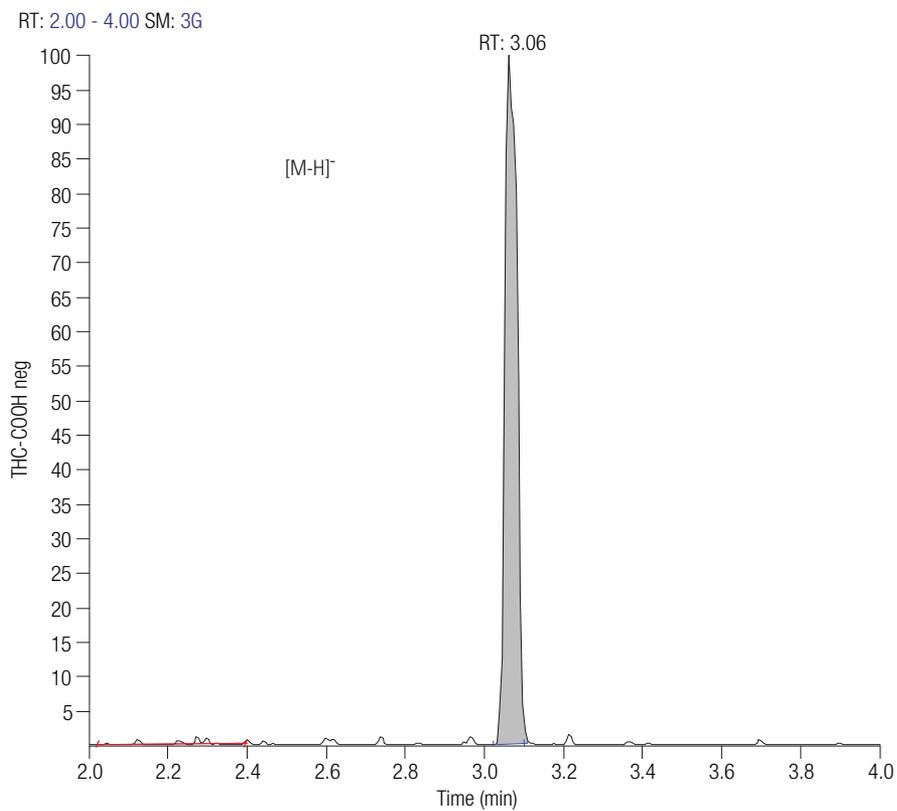


Figure 1. Ion chromatogram of 7.8 ng/mL urine calibration standard

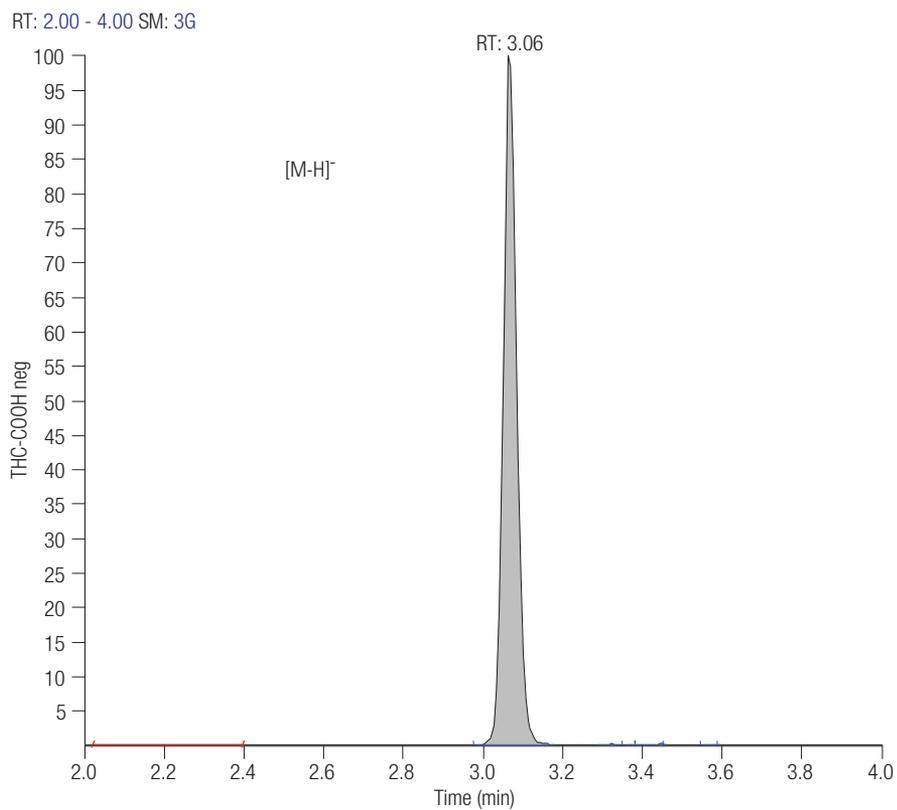


Figure 2. Ion chromatogram of 1000 ng/mL urine calibration standard

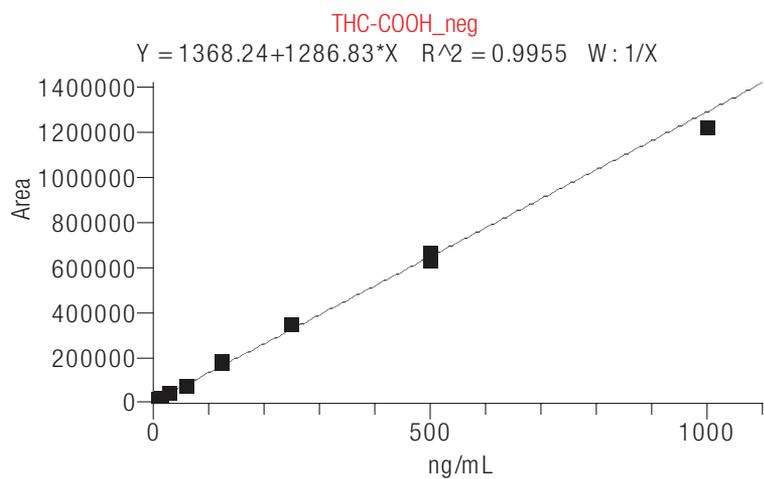


Figure 3. Calibration curve of THC-COOH in negative ionization mode

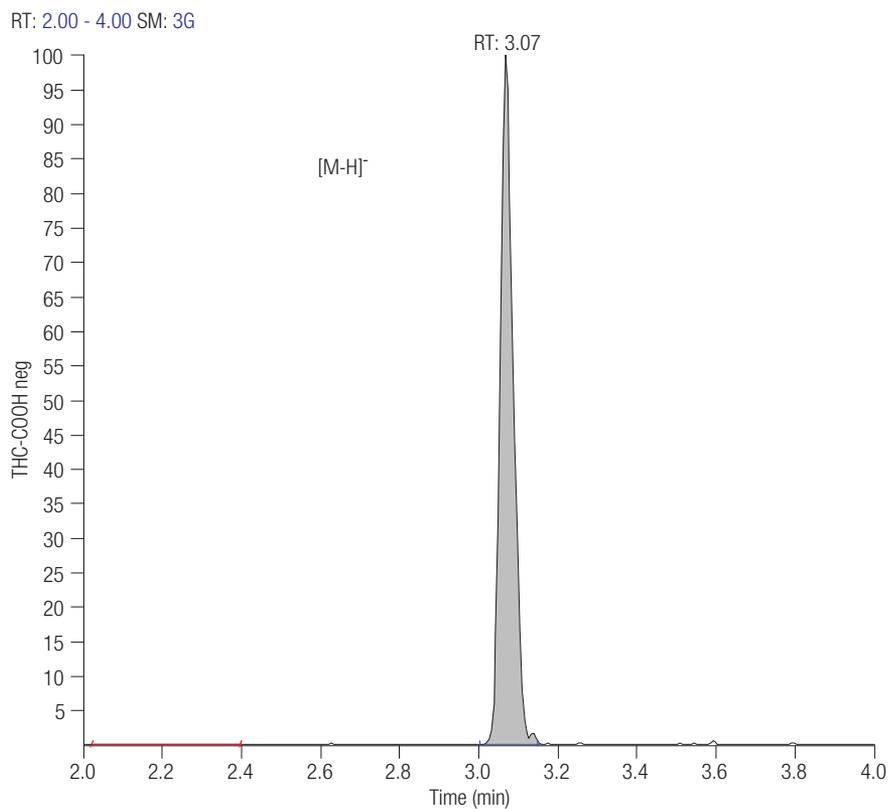


Figure 4. Ion chromatogram of urine sample containing 225 ng/mL. Sample was hydrolyzed and diluted 1:10 before the analysis.

RT: 2.00 - 4.00 SM: 3G

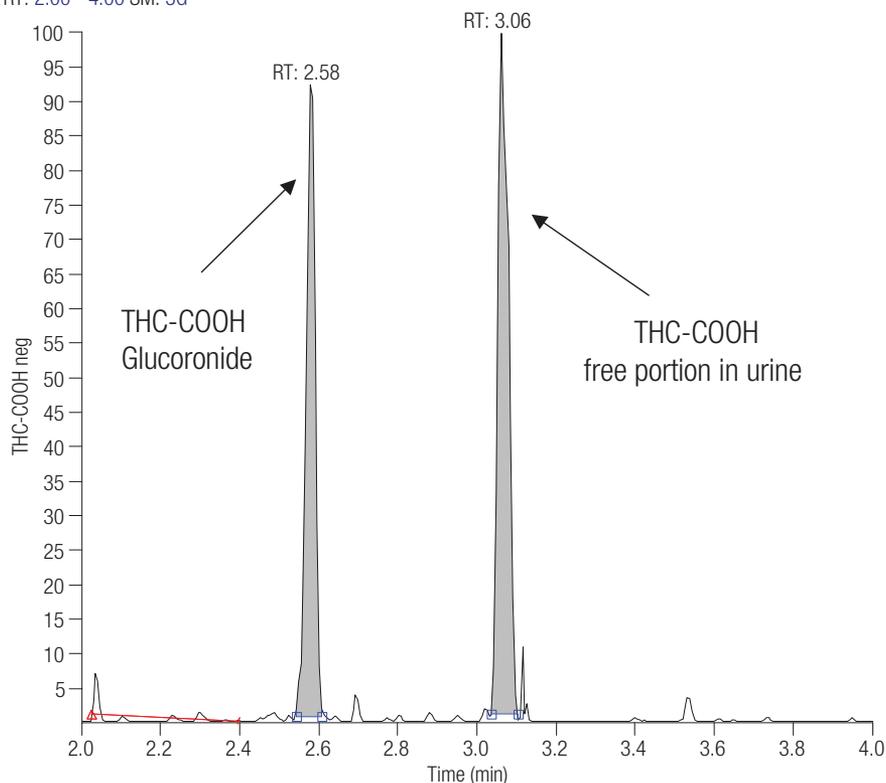


Figure 5. Ion chromatograms of urine sample containing 225 ng/mL. Sample was diluted 1:10 before the analysis; no hydrolysis was performed.

Conclusion

A robust 6-minute method for the quantification of THC-COOH with a dynamic range of 7.8-1000 ng/mL urine has been developed using the TSQ Quantum Access MAX mass spectrometer for forensic toxicology purposes.

www.thermoscientific.com

Legal Notices: ©2011 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



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

AN63443_E 10/11S