

Quantitation of 12 Benzodiazepines and Metabolites in Urine Using Ultrahigh Resolution LC-MS for Forensic Toxicology Use

Kent Johnson, Fortes Laboratories, Portland, OR; Marta Kozak, Thermo Fisher Scientific, San Jose, CA

Goal

To demonstrate the quantitation of 12 benzodiazepines in urine using a liquid chromatography-mass spectrometry (LC-MS) method and ultrahigh resolution with the Thermo Scientific Exactive benchtop mass spectrometer for forensic analysis.

Experimental

Standards and Samples Preparation

Calibration standards were prepared by spiking blank urine with 12 benzodiazepines (lorazepam, nordiazepam, oxazepam, temazepam, hydroxytriazolam, 7-aminoclonazepam, 7-aminonitrazepam, hydroxyalprazolam, 7-aminoflunitrazepam, desalkylflurazepam, diazepam, and 2-hydroxyethylflurazepam) to final concentrations ranging from 10 ng/mL to 2,000 ng/mL.

Calibration standards and urine samples were spiked with internal standards (10 deuterated benzodiazepines), hydrolyzed and processed using a solid phase extraction (SPE) procedure.

Third party QC samples containing 6 benzodiazepines were processed and analyzed to obtain method accuracy and precision.

HPLC

HPLC analysis was performed using a Thermo Scientific Accela liquid chromatography system with a Thermo Scientific Hypersil GOLD PFP column (50 x 2.1 mm; 5 µm). A processed sample of 5 µL was analyzed with a 6-minute gradient method.

Mass Spectrometry

MS analysis was carried out on an Exactive™ benchtop LC-MS instrument with an electrospray ionization (ESI) source. Full scan data with resolution of 100,000 (FWHM) was acquired.

Results and Discussion

Figure 1 displays 6 of the 12 selected benzodiazepines at 10 ng/mL and internal standards. Chromatograms for compound detection and quantitation are reconstructed with a mass tolerance of 5 ppm.

Figure 2 shows the calibration curve for this set. Data results for the other six benzodiazepines are available upon request.

Conclusion

The Exactive benchtop LC-MS instrument provides excellent quantitative analysis of 12 benzodiazepines, from 10 ng/mL to 5000 ng/mL in urine, using ultrahigh resolution full scan data acquisition in a 6-minute method. The accuracy, precision, LOQ, and linearity range of the method meet the demands of today's forensic toxicology laboratories.

Method Performance Summary

Target Analytes	Benzodiazepines
Matrix	Urine
Limit of Quantitation (LOQ)	10 ng/mL
Recovery	> 85%
Assay Linearity	10 ng/mL – 5000 ng/mL
Precision (%CV)	< 4%
Carryover at Lower Limit of Quantitation (LLOQ)	< 1%
Sample Volume	2 mL
Analysis Time	6 minutes

Analyte	Mean Conc.(ng/mL)	% Recovery	%RSD
Oxazepam	248	99.3	1
Nordiazepam	234	93.5	1.4
Temazepam	218	87.1	4
Desalkylflurazepam	214	85.7	4
Lorazepam	227	90.8	0.4
Hydroxyalprazolam	255	102	0.4

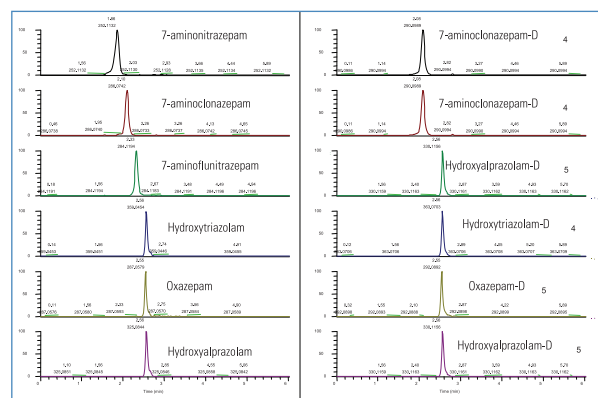


Figure 1: Chromatograms of 6 of the 12 selected benzodiazepines at 10 ng/mL and internal standards.

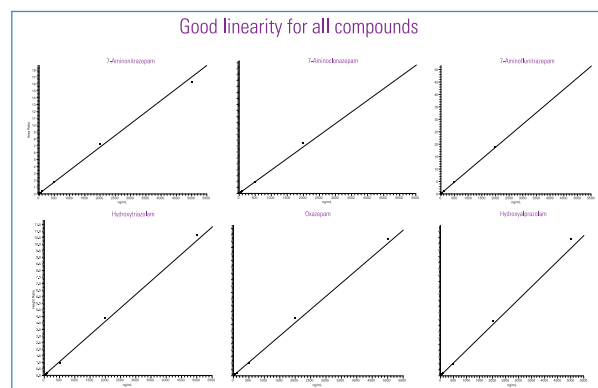


Figure 2: Calibration curves (10-5000 ng/mL) for all analytes

Key Words

- Exactive
- Accela HPLC
- Pain Management
- Forensic Toxicology

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.thermoscientific.com

AN63221b_E 06/16S

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.