TECHNICAL NOTE 66064

Quantification of eight antimycotics in human plasma using an Orbitrap Exploris 120 high-resolution accurate mass mass spectrometer for clinical research

Authors: Gaëtan Renoulin¹, Claudio De Nardi²

¹Thermo Fisher Scientific, Les Ulis, France

²Thermo Fisher Scientific, Reinach,
Switzerland

Keywords: Orbitrap Exploris 120 MS, TraceFinder software, Vanquish Flex UHPLC, antifungals, antimycotics, offline sample preparation, plasma, high-resolution mass spectrometry

Application benefits

- High-resolution accurate mass mass spectrometry for improved selectivity
- Simple offline sample preparation by protein precipitation
- Eight antimycotics drugs in a single 3.5-minute quantitative method

Goal

Implementation of an analytical method for the quantification of eight antimycotics in human plasma on a Thermo Scientific™ Orbitrap Exploris™ 120 mass spectrometer.

Introduction

Antifungals, also known as antimycotics, typically refer to a class of pharmaceutical fungicides used in the relief and prevention of mycosis, ranging from athlete's foot to ringworm to serious infections, such as cryptococcal meningitis. Voriconazole, posaconazole, fluconazole,



ketoconazole, and other similar antimycotics are used to address life-threatening fungal infections along with prevention of infections in immunocompromised individuals. The narrow therapeutic drug monitoring (TDM) research ranges of these antifungal agents, in addition to other complications, could lead to very different drug exposure from the same dosage regimen, and therefore, very different individual outcomes. Analytical methods to quantify such antimycotics were traditionally performed using high-performance liquid chromatography (HPLC) coupled with UV detectors. However, these methods require complicated extraction procedures and timeconsuming chromatography. LC-MS based methods are known for their superior selectivity and often result in significant reduction of the time spent on complicated sample preparation procedures and chromatography.



In this report, a clinical research analytical method for the quantification of eight antimycotics in human plasma in 3.5 minutes is presented. Samples were prepared by protein precipitation followed by chromatographic separation on a Thermo Scientific™ Vanquish™ Flex Binary UHPLC system. Detection was performed on an Orbitrap Exploris 120 mass spectrometer with heated electrospray ionization (HESI) operated in positive ionization mode. Method performance was evaluated using the ClinMass® TDM Platform with the ClinMass Add-On Set for Antimycotics from RECIPE Chemicals + Instruments GmbH (Munich, Germany) in terms of linearity of response, carryover, accuracy, and intra- and inter-assay precision for all analytes.

Experimental

Target analytes

The complete list of analytes and corresponding internal standards is reported in Table 1. The concentration ranges covered by the calibrators used are reported in Table 2.

Sample preparation

Reagents included four calibrators (MS9613 batch #1369) including blank and two controls from RECIPE (MS9682 batch #1369), as well as eight stable-isotope-labeled internal standards for the quantification.

Samples of 50 μ L of plasma or serum were protein precipitated using 100 μ L of precipitating solution containing the internal standards. Precipitated samples were vortex-mixed (30 seconds) and centrifuged (5 minutes at 10,000 \times g). The supernatant was diluted by ten by using a solution provide by RECIPE (MS9022). Finally, the sample was transfered to a clean vial.

Liquid chromatography

A Vanquish Flex Binary UHPLC system was used for chromatographic separation, utilizing mobile phases and an analytical column provided by RECIPE. Details of the analytical gradient are reported in Table 3. Total runtime was 3.5 minutes.

Table 1. List of analytes and internal standards

Compound name	Formula	Expected mass (m/z)	Internal standard name	Formula	Expected mass (m/z)
5-Fluorocytosine	C ₄ H ₄ FN ₃ O	130.0411	¹³ C- ¹⁵ N ₂ -5-Fluorocytosine	¹³ C ¹⁵ N ₂ C ₃ H ₄ FNO	133.0385
Fluconazole	C ₁₃ H ₁₂ F ₂ N ₆ O	307.1113	d ₄ -Fluconazole	$C_{13}H_8D_4F_2N_6O$	311.1365
Isavuconazole	C ₂₂ H ₁₇ F ₂ N ₅ OS	438.5078	¹³ C-d ₄ -Isavuconazole	¹³ CC ₂₁ H ₁₃ D ₄ F ₂ N ₅ OS	443.1480
Itraconazole	C ₃₅ H ₃₈ Cl ₂ N ₈ O ₄	705.2466	d ₅ -Itraconazole	C ₃₅ H ₃₃ D ₅ Cl ₂ N ₈ O ₄	710.2780
Ketoconazole	$C_{26}H_{28}CI_2N_4O_4$	531.1560	d ₈ -Ketoconazole	$C_{26}H_{20}D_8CI_2N_4O_4$	539.2063
OH-Itraconazole	C ₃₅ H ₃₈ Cl ₂ N ₈ O ₅	721.2415	d ₅ -OH-Itraconazole	C ₃₅ H ₃₃ D ₅ Cl ₂ N ₈ O ₅	726.2729
Posaconazole	$C_{37}H_{42}F_2N_8O_4$	701.3370	d ₄ -Posaconazole	$C_{37}H_{38}D_4F_2N_8O_4$	705.3621
Voriconazole	C ₁₆ H ₁₄ F ₃ N ₅ O	350.1223	d ₃ -Voriconazole	C ₁₆ H ₁₁ D ₃ F ₃ N ₅ O	353.1412

Table 2. Concentration ranges covered by the calibrators (MS9613 batch #1369)

Compound name	Concentration range (mg/L)
5-Fluorocytosine	4.90–108
Fluconazole	0.622-13.5
Isavuconazole	0.481–10.8
Itraconazole	0.146-3.11
Ketoconazole	0.430-8.88
OH-Itraconazole	0.171-3.55
Posaconazole	0.233-4.90
Voriconazole	0.275-5.96

Table 3. LC gradient profile

0.6	0	
0.6		
0.0	0	
0.6	30	
0.6	60	
0.6	98	
0.6	98	
0.6	0	
0.6	0	
ther parameters		
Injection volume		
	40 °C	
	0.6 0.6 0.6 0.6 0.6	

Mass spectrometry

Analytes and internal standards were detected by Full Scan – data-dependent MS² (ddMS²) on an Orbitrap Exploris 120 mass spectrometer using a HESI source operated in positive ionization mode. The Full Scan experiment was used for quantification and ddMS² for confirmation based on ion ratio. A summary of the MS conditions is reported in Table 4.

Table 4. MS parameters

lon source parameters						
Source type	Heated Electrospray Source Ionization (H-ESI)					
Spray voltage - Positive (V)	3,750					
Sheath gas (Arb)	55					
Aux gas (Arb)	10					
Sweep gas (Arb)	2					
Ion transfer tube temp. (°C)	320					
Vaporizer temp (°C)	450					
Settings						
Mild trapping	No					
Internal mass calibration	RunStart EASY-IC™					
Data acquisition mode	Full Scan - ddMS					
Full Scan parameters						
Resolution (at m/z 200)	60,000					
Scan range (m/z)	70-1400					
Expected LC peak width(s)	6					
RF lens (%)	70					
AGC target	Standard (1e6)					
Maximum injection time mode	Auto					
Polarity	Positive					
Data-Dependent MS ² scan properties						
Isolation window (m/z)	2					
Collision energy type	Normalized					
HCD collision energy (%)	30					
Resolution (at m/z 200)	15,000					
Scan range mode	Auto					

Method evaluation

The method performance was evaluated in terms of linearity of response within the calibration ranges, LLOQ, carryover, accuracy, intra- and inter-assay precision for all analytes. Carryover was calculated in terms of percentage ratio between the peak area of the highest calibrator and a subsequent blank sample injection. Analytical accuracy was evaluated in terms of percentage bias between nominal and average calculated concentrations using quality control samples at two different levels provided by RECIPE (MS9682 batch #1369). The LLOQ was evaluated by diluting the lowest calibrator down to 20-fold using blank matrix. LLOQs were determined as the lowest concentration with inaccuracy and precision below 20%.

Quality control samples were prepared and analyzed in replicates of five over three different days. Intra-assay precision for each day was evaluated in terms of percentage coefficient of variation (%CV) using the controls at two different levels in replicates of five (n=5). Inter-assay precision was evaluated as the %CV on the full set of samples (control samples at two levels in replicates of five prepared and analyzed on three different days, n=15).

Data analysis

Data were acquired and processed using Thermo Scientific™ TraceFinder™ 5.1 software.

Results and discussion

A linear response with 1/x weighting was obtained for all the analytes. The percentage bias between nominal and back-calculated concentration was always within $\pm 10\%$ for all the calibrators in all the runs. Chromatograms for the lowest calibrator for representative analytes and their internal standards are reported in Figure 1. Representative calibration curves are reported in Figure 2.

No significant carryover was observed for any of the analytes, with no signal detected in the blank injection just after the highest calibrator.

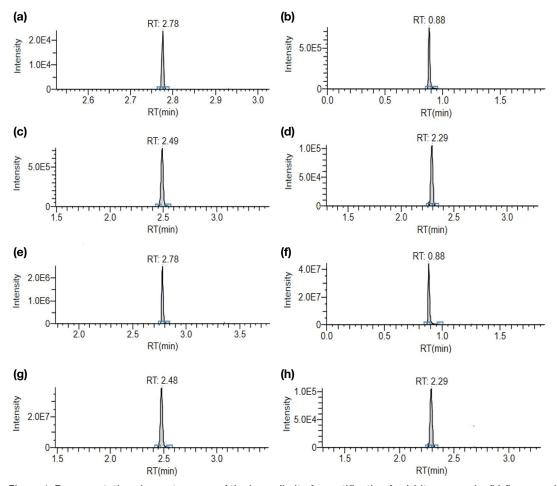


Figure 1. Representative chromatograms of the lower limit of quantification for (a) itraconazole, (b) fluconazole, (c) isavuconazole, (d) OH-itraconazole, (e) d_s -itraconazole, (f) d_4 -fluconazole, (g) 13 C- d_4 -isavuconazole, (h) d_s -OH-itraconazole

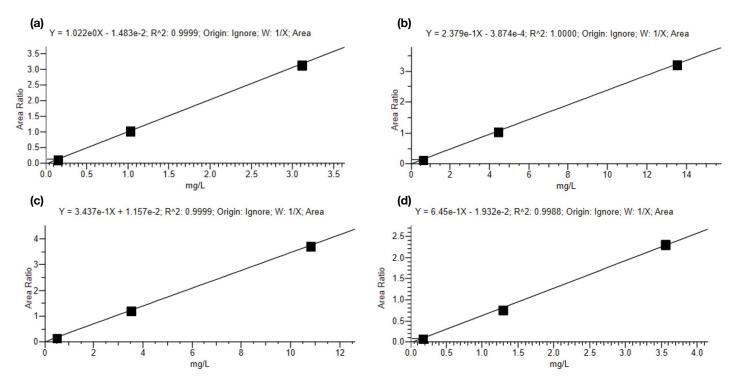


Figure 2. Representative calibration curves for (a) itraconazole, (b) fluconazole, (c) isavuconazole, (d) OH-itraconazole

The data demonstrated good accuracy of the method with the percentage bias between nominal and average backcalculated concentration for the used control samples ranging between -4.3% and 6.0% (Table 5). The %CV for intra-assay precision was always below 9.3% for all the analytes. The maximum %CV for inter-assay precision including all the analytes was 7.0%. Results for intra- and inter-assay precision are reported in Table 6.

Table 5. Analytical accuracy results for control MS9682 batch #1369

Compound name	Control	Nominal concentration (mg/L)	Average calculated concentration (mg/L)	Bias (%)
5-Fluorocytosine	Level I	19.9	20.4	2.4
5-Fluorocytosine	Level II	46.7	48.8	4.5
Fluconazole	Level I	2.43	2.58	6.0
Flucoriazole	Level II	5.79	6.03	4.2
Isavuconazole	Level I	1.95	2.05	5.0
Isavucoriazoie	Level II	4.59	4.82	4.9
Itraconazole	Level I	0.590	0.565	-4.3
Itracoriazoie	Level II	1.31	1.36	4.2
Ketoconazole	Level I	1.71	1.80	5.2
Keloconazoie	Level II	3.90	4.12	5.7
OH-Itraconazole	Level I	0.678	0.679	0.2
On-itracoriazole	Level II	1.60	1.58	-1.2
Posaconazole	Level I	0.909	0.934	2.8
r usacunazule	Level II	2.19	2.18	-0.6
Voriconazole	Level I	1.10	1.14	3.4
VONGONAZOIE	Level II	2.59	2.69	4.0

Table 6. Analytical intra- and inter-assay precision results for control MS9682 batch #1369

				Intra-assay				Inter-assay	
		Day 1		Day 2		Day 3			
Compound name	Control	Average calculated concentration (mg/L)	CV (%)	Average calculated concentration (mg/L)	CV (%)	Average calculated concentration (mg/L)	CV (%)	Average calculated concentration (mg/L)	CV (%)
5-Fluorocytosine	Level I	21.5	2.5	19.7	4.7	20.0	4.2	20.4	4.7
5-Fluorocytosine	Level II	51.9	5.0	47.7	7.5	46.7	6.5	48.8	5.6
Fluconazole	Level I	2.68	1.9	2.51	4.9	2.54	4.0	2.58	3.5
Fluconazole	Level II	6.23	1.8	5.97	8.2	5.90	5.9	6.03	2.8
1	Level I	2.12	1.7	2.01	4.6	2.01	4.3	2.05	2.9
Isavuconazole	Level II	4.96	1.8	4.78	8.4	4.71	5.9	4.82	2.6
11	Level I	0.585	3.5	0.554	2.3	0.554	5.9	0.565	3.2
Itraconazole	Level II	1.43	3.3	1.32	7.2	1.34	4.1	1.36	4.4
IZ-L	Level I	1.87	2.6	1.76	5.2	1.77	4.1	1.80	3.2
Ketoconazole	Level II	4.25	1.9	4.10	9.3	4.02	6.7	4.12	2.8
	Level I	0.733	6.1	0.658	4.6	0.646	7.9	0.679	7.0
OH-Itraconazole	Level II	1.66	1.1	1.55	9.3	1.53	5.5	1.58	4.5
December	Level I	0.96	2.3	0.93	4.4	0.91	5.6	0.93	2.9
Posaconazole	Level II	2.23	2.7	2.17	7.8	2.13	4.8	2.18	2.5
Variance	Level I	1.19	2.6	1.12	4.4	1.12	3.7	1.14	3.7
Voriconazole	Level II	2.75	1.7	2.64	8.0	2.61	5.5	2.67	2.8

thermoscientific

LLOQs of all compounds were determined and reported in Table 7.

Table 7. LLOQs for all compounds

Compound name	LLOQ (mg/L)
5-Fluorocytosine	0.490
Fluconazole	0.0622
Isavuconazole	0.0481
Itraconazole	0.146
Ketoconazole	0.0430
OH-Itraconazole	0.171
Posaconazole	0.0233
Voriconazole	0.0275

Conclusions

A robust, reproducible, and sensitive liquid chromatography-HRAM mass spectrometry method for clinical research for quantification of eight antimycotics in human plasma was developed and implemented.

Sample preparation consisted of a simple offline protein precipitation with concomitant internal standard addition. The method was evaluated on an Vanquish Flex UHPLC system coupled to an Orbitrap Exploris 120 mass spectrometer. Method performance was evaluated using the ClinMass TDM Platform with the ClinMass Add-On Set for Antimycotics from RECIPE. The described method meets research laboratory requirements in terms of sensitivity, linearity of response, accuracy, and precision.

Find out more at thermofisher.com/ClinicalResearchApps

For General Lab Use only. Not for use in diagnostic procedures. © 2021 Thermo Fisher Scientific Inc. All rights reserved. ClinMass is a registered trademark of RECIPE Chemicals + Instruments GmbH. All other trademarks are the property of Thermo Fisher Scientific 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. TN66064-EN 0721S

