

Evaluation of Benchtop Quadrupole Orbitrap Ultra-High-Resolution Mass Spectrometer in Rapid Quantitative Analysis of Immunosuppressant Drugs in Blood Samples

Mindy Gao, Marta Kozak
Thermo Fisher Scientific, San Jose, CA, USA

Overview

Purpose: To evaluate a benchtop quadrupole Orbitrap mass spectrometer for rapid quantitative analysis of immunosuppressant drugs in human blood samples for clinical research.

Methods: Blood samples were processed by protein precipitation, and analytes were chromatographically separated with a 3-minute LC gradient method. The mass spectrometer collected high resolution MS/MS spectra for each analyte, and the most abundant fragment was used for quantitation. The method was evaluated by obtaining limits of quantitation, intra- and inter-assay precisions, and investigating matrix effects. Cross-correlation between this method and a method from a collaborator's research laboratory was determined.

Results: We developed a fast and cost efficient method for analysis of immunosuppressant drugs in blood. The method performance met clinical research laboratory criteria. Implementation of high resolution mass spectrometers in clinical research laboratories has advantages over conventionally used triple quadrupole instruments because of their versatile application range: screening, quantitation, and structure elucidation.

Introduction

Clinical research labs commonly use selective and cost-efficient LC-MS techniques for the analysis of immunosuppressant drugs. Usually, the quantitative method is developed on a triple quadrupole mass spectrometer. Here, we evaluated a high-throughput method implemented on a Thermo Scientific™ Q Exactive Focus™ hybrid quadrupole-Orbitrap mass spectrometer for improved selectivity when compared to the conventional LC-MS platform.

Methods

Calibrators

Calibrators containing cyclosporin A, everolimus, sirolimus, and tacrolimus in whole blood were purchased from Chromsystems Instruments & Chemicals GmbH (TABLE 1).

TABLE 1. Concentrations of calibrators.

Analyte	Cal 1	Cal 2	Cal 3	Cal 4	Cal 5	Cal 6
	Concentration (ng/mL)					
Tacrolimus	2.35	5.96	11.9	17.6	24.6	42.2
Sirolimus	2.27	6.12	11.9	18.4	27.9	46.1
Everolimus	2.27	5.94	11.7	18.1	25.1	45.4
Cyclosporin A	26.1	123	294	485	764	942

Sample Preparation

•100 μ L of blood + 400 μ L of precipitation solvent containing 150 μ L of 0.1 M ZnSO₄ in water and 250 μ L of internal standard (10 ng/mL ascomycin and 250 ng/mL cyclosporin D) in methanol

- Vortex, store at approx. 4 °C for 10 min
- Vortex, centrifuge for 10 min at 12,000 rpm
- Transfer supernatant into HPLC injection vial
- Inject 20 μ L into LC-MS system

Liquid Chromatography

LC system

Pump: Thermo Scientific™ Dionex™ UltiMate™ 3000

Autosampler: Thermo Scientific™ OAS-3X00TXRS

Mobile phase A: 0.1% formic acid, 10 mM ammonium formate in water

Mobile phase B: 0.1% formic acid, 10 mM ammonium formate in methanol

Mobile phase C: acetonitrile/isopropyl alcohol/acetone 45:45:10 v/v/v

Mass Spectrometry

Thermo Scientific™ Q Exactive™ Focus Orbitrap mass spectrometer with HESI ionization source

Data acquisition method: parallel reaction monitoring (PRM) experiment collecting MS/MS spectra for each analyte (TABLE 3)

Precursor isolation width: 1 amu

Resolution: 17.5K

Data Analysis

The most abundant fragment from MS/MS spectrum was selected for quantification (TABLE 3, FIGURE 1). Thermo Scientific™ TraceFinder™ software version 3.2 was used for data acquisition and data processing.

TABLE 2. HPLC gradient method.

Time (min)	Flow rate (mL/min)	%A	%B	%C
0.00	0.5	70	30	0
0.25	0.5	70	30	0
0.50	0.5	0	100	0
1.80	0.5	0	100	0
1.81	0.5	0	0	100
2.00	0.5	0	0	100
2.01	1.5	0	0	100
2.29	1.5	0	0	100
2.30	1.5	0	100	0
2.31	1.5	70	30	0
3.00	1.5	70	30	0

TABLE 3. Mass spectrometer scan parameters and quantification fragments

Analyte	Precursor Mass [m/z]	Quantitation Fragment [m/z]	Collision Energy (eV)	Polarity
Cyclosporin D (IS)	1233.9	1216.8633	24	Positive
Cyclosporin A	1219.9	1202.8478	20	Positive
Everolimus	975.6	908.5494	20	Positive
Sirolimus	931.6	864.5242	22	Positive
Tacrolimus	821.5	768.4666	22	Positive
Ascomycin (IS)	809.5	756.4670	24	Positive

Method Performance Evaluation

Quality Control (QC) samples containing cyclosporin A, everolimus, sirolimus, and tacrolimus in whole blood were purchased from Chromsystems Instruments & Chemical GmbH (TABLE 4).

TABLE 4. QC sample concentrations.

Analyte	QC1	QC2	QC3	QC4
Tacrolimus	2.6	7.3	16.7	34.2
Sirolimus	2.9	10.1	20.4	38.5
Everolimus	2.3	4.4	8.5	28.8
Cyclosporin A	53.0	276	514	1111

Limit of quantitation (LOQ) were defined as the lowest concentrations that had back-calculated values within 20% of nominal and RSD for 5 replicates within 20%. FIGURE 1 shows chromatograms of the lowest calibration standards.

Method accuracy was evaluated by obtaining % difference from nominal for quality control samples.

Method precision was evaluated by analyzing five replicates of each QC sample in three different days (TABLE 5 and TABLE 6).

System robustness and reproducibility were evaluated by analyzing five replicate injections of each calibration standard containing cyclosporin A, everolimus, sirolimus, and tacrolimus (TABLE 7).

Matrix effects were evaluated by spiking 10 ng/mL of everolimus, sirolimus, tacrolimus, and 60 ng/mL of cyclosporin A into whole blood samples from five different donors in two replicates, and to three solvent blanks. Each sample was injected twice. Relative and absolute % recoveries were calculated against data collected for spiked solvent (TABLE 8). A total of 97 donor blood samples containing tacrolimus and cyclosporin A provided by collaborator lab were analyzed. Correlation between data generated with collaborator clinical research method using triple quadrupole mass spectrometer and method on Q Exactive Focus was calculated (FIGURE 4)

Results

FIGURE 1. MS/MS spectra for analytes and internal standards.

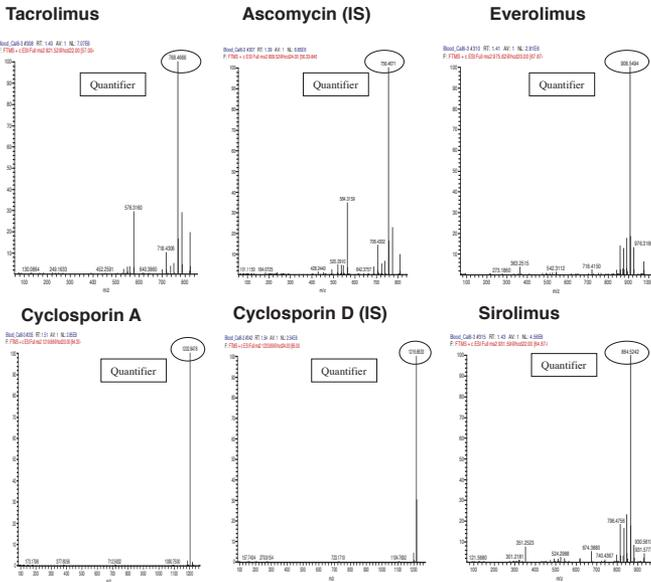


FIGURE 2. Chromatogram of lowest calibration standard (LOQ) reconstructed with a mass accuracy of 5 ppm

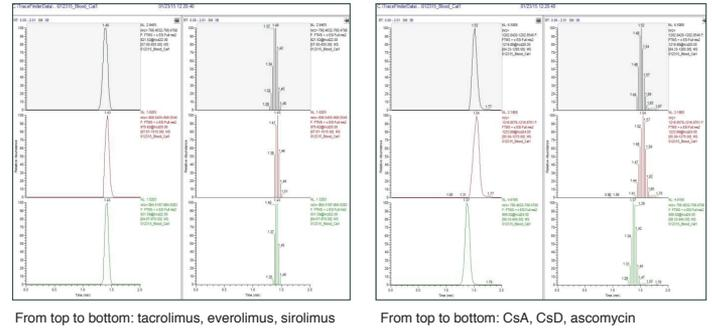


TABLE 5. Intra-assay precision and accuracy. Five replicates of QC samples containing cyclosporin A, tacrolimus, sirolimus, and everolimus were processed and analyzed in three batches.

Analyte	QC1	QC2	QC3	QC4
Tacrolimus	4.12 – 5.01	3.68 – 5.64	2.65 – 4.19	3.88 – 5.12
Sirolimus	2.27 – 5.49	3.40 – 4.48	3.21 – 6.06	4.28 – 4.96
Everolimus	3.50 – 4.59	2.04 – 4.29	3.24 – 5.67	2.12 – 5.25
Cyclosporin A	2.21 – 4.21	2.76 – 4.88	3.86 – 6.85	4.58 – 5.08

Analyte	Day-1	Day-2	Day-3
Tacrolimus	7.73% to -14.7%	4.23% to -11.2%	8.77% to -9.38%
Sirolimus	5.91% to -8.76%	9.41% to -9.48%	9.58% to -13.3%
Everolimus	8.34% to -6.49%	7.24% to -9.17%	10.2% to -6.62%
Cyclosporin A	8.82% to -16.8%	10.5% to -12.0%	11.7% to -13.5%

FIGURE 3. Calibration curves and lowest calibration standard peaks.

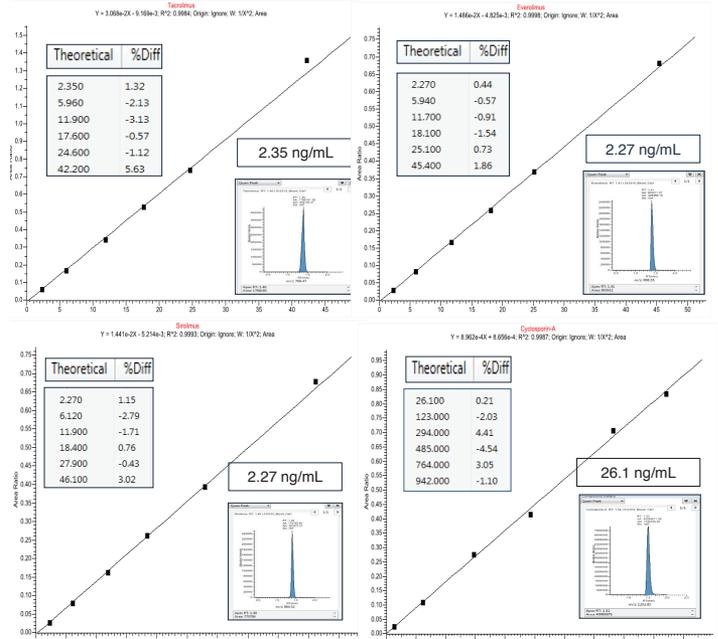


TABLE 6. Inter-assay precision and accuracy. Five replicates of QC samples were processed and analyzed in three batches.

Analyte	QC1	QC2	QC3	QC4
Tacrolimus	4.65	4.55	3.77	4.21
Sirolimus	4.16	3.71	4.22	4.24
Everolimus	5.92	4.47	4.46	3.88
Cyclosporin A	3.46	3.5	5.11	4.59

Analyte	QC1	QC2	QC3	QC4
Tacrolimus	101	94.2	95.2	99.8
Sirolimus	99.9	98.2	99.1	102
Everolimus	96.6	96.3	101	102
Cyclosporin A	101	94.4	94.7	104

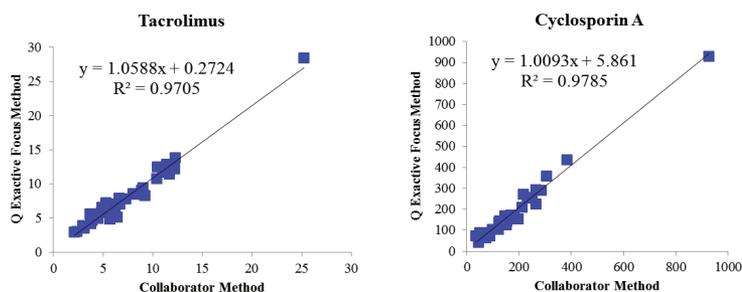
TABLE 7. System robustness and reproducibility. Five replicate injections of each calibration standard were performed to demonstrate the system robustness and reproducibility.

Analyte	Std-1	Std-2	Std-3	Std-4	Std-5	Std-6
%RSD						
Tacrolimus	3.65	3.03	3.05	3.50	3.06	4.39
Sirolimus	10.3	4.28	3.23	2.46	1.64	3.42
Everolimus	5.57	1.98	2.34	4.78	3.58	4.20
Cyclosporin A	3.93	2.81	0.97	1.17	3.82	1.83
Mean %Recovery						
Tacrolimus	102	96.2	101	96.2	101	104
Sirolimus	101	96.3	106	95.3	103	98.2
Everolimus	100	96.9	107	95.5	103	97.6
Cyclosporin A	99.8	101	102	96.6	101	99.9

TABLE 8. Matrix effects.

Analyte	Blood-1	Blood-2	Blood-3	Blood-4	Blood-5
Absolute %Recovery					
Tacrolimus	96.9	107	109	107	108
Sirolimus	94.3	93.2	103	103	104
Everolimus	104	98.6	112	109	108
Cyclosporin A	90.8	91.5	101	101	82.0
Relative (IS ratio) %Recovery					
Tacrolimus	85.1	92.9	97.0	94.8	93.7
Sirolimus	82.8	80.9	92.1	90.8	90.3
Everolimus	91.6	85.5	99.9	96.2	94.0
Cyclosporin A	98.5	95.9	96.3	101	85.5

FIGURE 4. Methods cross-correlation. Samples from collaborator lab that contain cyclosporin A and tacrolimus.



Conclusion

- We demonstrated a fast, cost-efficient, highly specific method for the analysis of immunosuppressant drugs on a high-resolution Orbitrap mass spectrometer.
- Implementation of a high-resolution instrument resulted in chromatograms containing just analyte peak, which significantly improves accuracy of data processing and reduces data review time.
- Method meets clinical research requirements for analysis of immunosuppressant drugs in whole blood samples.

For research use only. Not for use in diagnostic procedures

www.thermofisher.com

©2016 Thermo Fisher Scientific Inc. All rights reserved. Cerilliant is a trademark of Sigma-Aldrich Inc. 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 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.

Africa +43 1 333 50 34 0
Australia +61 3 9757 4300
Austria +43 810 282 206
Belgium +32 53 73 42 41
Canada +1 800 530 8447
China 800 810 5118 (free call domestic)
 400 650 5118
 PNG4481-EN 0816S

Denmark +45 70 23 62 60
Europe-Other +43 1 333 50 34 0
Finland +358 10 3292 200
France +33 1 60 92 48 00
Germany +49 6103 408 1014
India +91 22 6742 9494
Italy +39 02 950 591

Japan +81 45 453 9100
Korea +82 2 3420 8600
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
Norway +46 8 556 468 00

Russia/CIS +43 1 333 50 34 0
Singapore +65 6289 1190
Spain +34 914 845 965
Sweden +46 8 556 468 00
Switzerland +41 61 716 77 00
UK +44 1442 233555
USA +1 800 532 4752

Thermo
 SCIENTIFIC

A Thermo Fisher Scientific Brand