

# Parallel Reaction Monitoring and Selected Reaction Monitoring Exhibit Comparable Quantitative Performance in Clinical Research and Forensic Toxicology Applications

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

## Overview

### Purpose:

To evaluate the analytical performance of the parallel reaction monitoring (PRM) method on a Thermo Scientific™ benchtop high resolution quadrupole-Orbitrap™ mass spectrometer and draw a performance comparison to the selected reaction monitoring (SRM) method on a triple quadrupole mass spectrometer.

### Methods:

Linearity range, precision, and accuracy were compared on two applications, epi-vitamin D quantitation in plasma and barbiturate measurement in urine.

### Results:

We demonstrated comparable quantitative performance between PRM and SRM across multiple differing parameters including matrix, polarity, and ionization mode.

## Introduction

SRM, performed on triple quadrupole (QqQ) mass spectrometers (MS), has emerged as the MS “gold-standard” for targeted quantification in clinical research and forensic toxicology applications.

With the development of the hybrid quadrupole-Orbitrap MS instruments, an alternative quantitative method is PRM. In PRM, the third quadrupole of a triple quadrupole is essentially substituted with a high resolution accurate mass Thermo Scientific™ Orbitrap™ mass analyzer to permit the parallel detection of all target product ions in one, concerted high resolution mass analysis.

## Methods

### Sample Preparation

For the application of epi-vitamin D quantitation (25-OH vitamin D3, epi-25-OH vitamin D3, 25-OH vitamin D2), plasma samples were processed by protein precipitation followed by solid phase extraction.

For the application of barbiturate measurement, urine samples were diluted 20 fold. The five barbiturates are amobarbital (amo), butalbital, phenobarbital (pheno), pentobarbital (pento), secobarbital (seco).

After preparation, the same samples were divided equally to two aliquots.

### Liquid Chromatography

Identical front ends (autosampler, LC pump, and column heater) were used for both methods. They are Thermo Scientific™ Dionex™ UltiMate™ 3000RS LC system with an OAS autosampler.

For epi-vitamin D application, five minutes LC separation was run on Thermo Scientific™ Hypersil GOLD™ PFP column.

For barbiturates application, the LC column used was a Thermo Scientific™ Accucore™ C18 column. The total run time was six minutes.

### Mass Spectrometry

PRM was conducted on a Thermo Scientific™ Q Exactive™ Focus quadrupole-Orbitrap mass spectrometer (FIGURE 1).

SRM was conducted on Thermo Scientific™ TSQ Endura™ triple quadrupole mass spectrometer (FIGURE 1).

In PRM, a single precursor ion was selected in the quadrupole with an isolation width of 2.0 m/z and fragmented in the HCD cell using an optimized, compound-specific collision energy. The resulting MS/MS product ion spectrum was detected in the Orbitrap detector at a resolution of 35,000 (FWHM at m/z of 200). In SRM, the cycle time was 0.5 second. Q1 and Q3 resolution was set as 0.7 Da (FWHM).



Q Exactive Focus quadrupole-Orbitrap MS



TSQ Endura triple quadrupole MS

**FIGURE 1. Two mass spectrometers used in this study.**

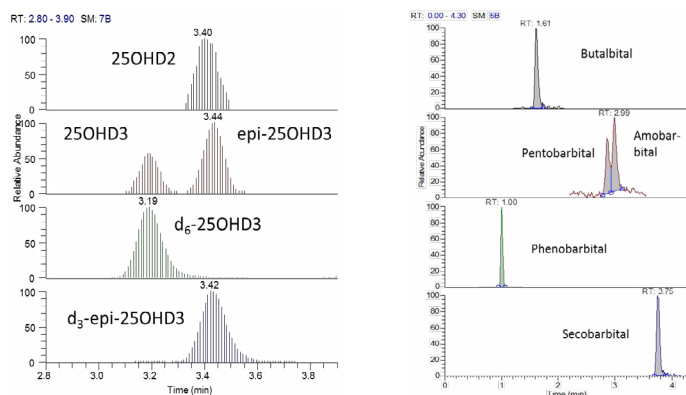
## Results

For epi-vitamin D analysis, both PRM and SRM were conducted in positive polarity with APCI probes.

For barbiturates analysis, both PRM and SRM were conducted in negative polarity with heated ESI probes.

For the PRM analysis of barbiturates, for each analyte and internal standard (IS), the precursor extract mass in the MS/MS spectrum was used for quantification and the most abundant fragment was used for confirmation. For epi-vitamin D analysis, a specific fragment from the MS/MS spectrum was selected as the quantifying ion. In both cases, the chromatograms were reconstructed with a mass accuracy of 5 ppm for quantification.

In SRM analysis of both sets of compound, two products ions were collected. One was for quantification and another one was for ion ratio confirmation. To ensure confident quantification, at least 15 scans were collected across each peak for both PRM and SRM analysis.



**FIGURE 2. Chromatographic separation of epi-vitamin D (left) and barbiturates (right).**

**TABLE 1. Linearity range (ng/mL)**

	25OH D2	25OH D3	epi-25OH D3
PRM	1-100	1-100	1-100
SRM	2-100	2-100	2-100

	Amo	Butalbital	Pento	Pheno	Seco
PRM	5-2000	5-2000	5-2000	25-2000	5-2000
SRM	10-2000	5-2000	10-2000	25-2000	5-2000

**TABLE 2. Inter-assay precision (% RSD, concentration ng/mL) using PRM (n=15). BQL, below quantitation limit.**

	25OH D2	25OH D3	epi-25OH D3
QC0	BQL	6.4 (8.4)	BQL
QC1	5.3 (6.0)	4.1 (14.4)	4.4 (6.6)
QC2	3.1 (15.0)	3.2 (23.4)	2.9 (15.6)
QC3	4.2 (50.0)	3.5 (58.4)	2.5 (50.6)

**TABLE 3. Inter-assay precision (% RSD, concentration is the same as above) using SRM (n=15).**

	25OH D2	25OH D3	epi-25OH D3
QC0	BQL	7.6	BQL
QC1	7.6	8.6	6.9
QC2	6.4	4.8	9.4
QC3	5.2	3.4	6.9

**TABLE 4. Inter-assay precision (% RSD) using PRM (n=15).**

	Amo	Butalbital	Pento	Pheno	Seco
Low QC (25 ng/mL)	4.6	4.5	6.5	5.5	5.1
Medium QC (100 ng/mL)	9.7	6.1	6.2	5.6	5.7
High QC (1000 ng/mL)	7.0	5.4	6.1	4.1	5.1

**TABLE 5. Inter-assay precision (% RSD) using SRM (n=15).**

	Amo	Butalbital	Pento	Pheno	Seco
Low QC	4.6	4.5	5.4	8.4	6.3
Medium QC	1.8	2.4	1.8	5.0	3.8
High QC	1.4	1.9	1.9	2.1	0.9

TABLE 6. Accuracy (% recovery, NIST standard concentration ng/mL) using PRM.

	25OH D2	25OH D3	epi-25OH D3
Level 1	BQL (0.54)	108 (28.8)	103 (1.84)
Level 2	BQL (0.81)	110 (18.1)	90.7 (1.29)
Level 3	97 (13.3)	104 (19.8)	114 (1.18)
Level 4	BQL (0.55)	106 (29.4)	90.9 (26.4)

TABLE 7. Accuracy (% recovery) using SRM.

	25OH D2	25OH D3	epi-25OH D3
Level 1	BQL	93.4	BQL
Level 2	BQL	102	BQL
Level 3	92.7	105	BQL
Level 4	BQL	98	100

Conclusions

- We demonstrated comparable quantitative performance between PRM (Q Exactive Focus quadrupole-Orbitrap MS) and SRM (TSQ Endura triple quadrupole MS).
- Q Exactive Focus MS is a versatile instrument. In addition to quantitation, it can also be used for screening, structure elucidation and bulk substance identification in clinical research and forensic toxicology applications.

www.thermofisher.com

©2016 Thermo Fisher Scientific Inc. All rights reserved. All 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  
**Brazil** +55 11 2730 3006  
**Canada** +1 800 530 8447  
**China** 800 810 5118 (free call domestic)  
400 650 5118

**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 6 6885 1213  
**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  
**Sweden** +46 8 556 468 00  
**Switzerland** +41 61 716 77 00  
**Taiwan** +886 2 8751 6655  
**UK/Ireland** +44 1442 233555  
**USA** +1 800 532 4752  
PN64701-EN 0616S

