

High-Throughput LC-MS/MS Measurement of Pregnenolone in Human Blood Serum for Research Purposes

Joe Di Bussolo¹, Marta Kozak¹, Ian White², Emily Herman², Raidiri Castillo³, Amit Shah³ and Hashim Othman³; ¹Thermo Fisher Scientific, West Chester, PA; ²West Chester University of PA, West Chester, PA; ³Bio-Reference Laboratories, Elmwood Park, NJ

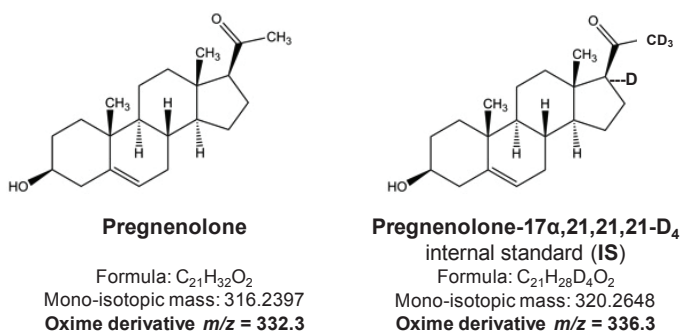
ABSTRACT

We developed an LC-MS/MS method for researchers to accurately measure pregnenolone in blood serum from 10 to 500 ng/dL with a throughput of at least 12 injections per hour. This was done by liquid-liquid extraction (LLE) of blood serum followed by derivatization with hydroxylamine and quantitative analysis using a 4-channel ultra high-performance liquid chromatography (UHPLC) system coupled to a tandem mass spectrometer (MS/MS). Oxime derivatives of pregnenolone and its internal standard were separated from sample matrix components by gradient elution through a UHPLC column packed with solid-core silica particles having an alkyl-aromatic bonded phase on its surfaces. Analytes were eluted to a heated electro-spray ionization (HESI) probe of a triple-quadrupole mass spectrometer where selected-reaction monitoring (SRM) detected the analytes. The desired analytical range of 10 to 500 ng/dL was achieved with inter- and intra-batch reproducibility of less than 8%, carryover less than 0.2% and acceptable correlation of specimen results with those from a reference laboratory. Throughputs of 13 to 52 injections/hour were achieved.

INTRODUCTION

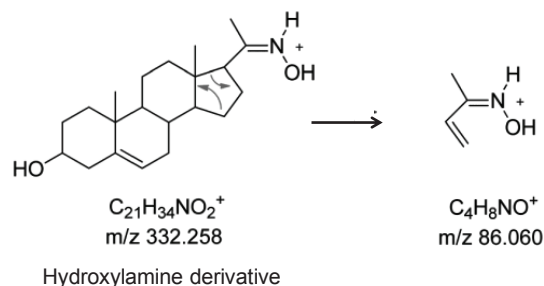
Pregnenolone is a biosynthetic precursor to other steroids such as corticosteroids, androgens and estrogens. It is converted to progesterone by 3-beta-hydroxysteroid dehydrogenase or to 17-OH-pregnenolone by 17-alpha-hydroxylase. Researchers investigating anomalies of these enzymes, which result in steroid hormone imbalances, need to measure pregnenolone within an analytical range of 10 to 500 ng/dL of blood serum. The chemical structures of pregnenolone and its deuterated internal standard (IS) are shown in Figure 1. Since pregnenolone does not ionize well in either atmospheric-pressure chemical-ionization (APCI) or electro-spray ionization (ES) sources, derivatization with hydroxylamine was necessary to achieve the desired lower-limit of quantitation (LOQ). The resulting positive-ion oximes (Figure 2) were quantitated via UHPLC-MS/MS using pregnenolone-D4 as the IS.

Figure 1. Chemical structures of analyte and IS



Reactions with hydroxylamine form positive-ion oxime

Figure 2. Pregnenolone oxime fragmentation



MATERIALS AND METHODS

Sample Preparation

200 μ L aliquots of fresh blood serum specimens, as well as calibrators and quality control specimens (QCs), were spiked with pregnenolone-D4 intern standard (IS) before being subjected to liquid-liquid extraction with 1 mL *m*-butyl ether (MTBE). After drying the ether extracts by heated nitrogen flow residue of each was reacted with hydroxylamine to form positive-ion oxime derivatives. After drying the derivative preparations, the residue of each was reconstituted with water and methanol (1:1) to a total volume of 200 μ L and μ L injections were made into the LC-MS/MS system.

Test Methods

Using one or more channels of a Thermo Scientific™ Transcend™ LX-4 sy chromatographic segregation of the steroid oximes from unwanted sample components was achieved by gradient elution through a Thermo Scientific Accucore™ Phenyl-X column (2.6 μ m, 50 x 2.1 mm), which was heated to Chromatographic conditions are described in Figure 3. The Thermo Scientific TSQ Endura™ triple-quadrupole mass spectrometer was used with a heated electro-spray ionization (HESI) probe. Ion source and MS/MS conditions are described in Figure 4.

Instrument Control & Data Analysis

Thermo Scientific™ TraceFinder™ with Aria™ MX software was used to control the Transcend LX-4 and Endura MS/MS systems, submit batches to desired channels as well as for analyzing data and reporting results.

Figure 3. Liquid chromatography parameters

Column: Accucore Phenyl-X, 2.6 μ, 50 x 2.1 mm, @40°C						
Solvent A: Water + 0.1% Formic Acid						
Solvent B: Methanol						
Step	Start	Sec	Flow	Gradient	%A	%B
1	0.00	10	0.5	Step	50	50
2	0.17	20	0.5	Ramp	20	80
3	0.50	120	0.5	Ramp	-	100
4	2.50	60	0.5	Step	-	100
5	3.50	60	0.5	Step	50	50
Start data: 1.0 min Duration: 1.0 min						
Total run time: 4.5 min						

Figure 4. Mass spectrometer HESI & SRM parameters

Ion Source: HESI, +3500 V, vaporizer temp: 400°C				
SRM Transitions: Q1 & Q3 resolutions: 0.7				
Analyte	Q1	Q2	CE	RF
Pregnenolone oxime (Quan)	332.25	86.25	27	95
Pregnenolone oxime (Confirm)	332.25	300.25	22	95
Pregnenolone-D4 oxime (Quan)	336.30	90.30	27	95
Pregnenolone-D4 oxime (Confirm)	336.30	304.25	22	95

RESULTS

Quantitation Reliability

Typical results from calibrators, QCs and specimens are show in Figure 5. Ion ratios (Confirming/Quan transitions) averaged 66% for pregnenolone and 50 % for its IS.

IS peak areas among calibrators & QC averaged 17,400 with RSD of 12%. IS peak areas in specimens ranged from 10,600 to 14,700 and showed an average recovery of 65%. Carryover was less than 0.2%. Inter- and intra-batch precisions were less than 5 and 8% coefficient of variation (CV), respectively (Tables 1a & 1b).

Figure 5. Typical pregnenolone quantitative results

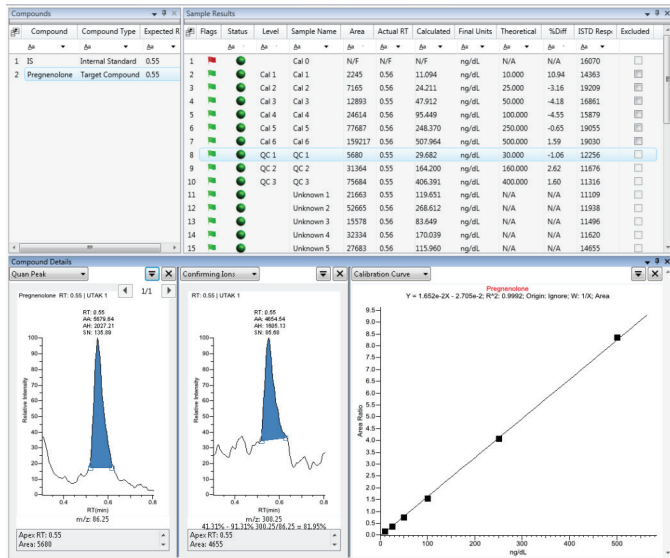


Table 1. Inter- and intra-batch precisions

Injection	Pregnenolone (ng/dL)	Date	Injection	Pregnenolone (ng/dL)
1	110	Day 1, 7/13/15	1	42
2	104		2	41
3	102		3	40
4	108		4	39
5	100		5	38
6	105	Day 2, 7/14/15	1	41
7	109		2	40
8	103		3	41
9	104		4	38
10	109		5	40
11	95	Day 3, 7/15/15	1	37
12	94		2	36
13	107		3	36
14	105		4	37
15	101		5	36
16	105	Day 4, 7/16/15	1	35
17	100		2	42
18	99		3	43
19	96		4	33
20	100		5	32
Mean	102.80		Mean	38.35
STD	4.69		SD	3.07
CV %	4.6		CV %	7.99

a. Inter-batch

b. Intra-batch

Throughput

Single-channel throughput was 13 injections per hour. When multi-channelled across 2, 3 or 4 channels, the throughput increased to 26, 39 and 52 injections per hour, respectively. Pregnenolone batches were also multi-channelled with estrogen batches which utilized the same MS source conditions.

Accuracy Assessment

Comparison of 40 specimen results ranging from 13 to 130 ng/dL with reference-lab results showed differences that averaged 3.8% and ranged from -20.0 to 24.7% (Table 2). Only 5 of 40 test samples (7.5%), deviated 20% or more. Deming regression analysis showed R² = 0.9706, slope = 1.080, intercept = -1.8 and standard error of estimate = 7.3%.

Table 2. Reference lab results comparison

Test Sample	Current Method	Reference Lab	Difference %	Test Sample	Current Method	Reference Lab	Difference %
1	116	93	24.73	21	27	30	-10.00
2	85	78	8.97	22	49	52	-5.77
3	86	73	17.81	23	66	73	-9.59
4	13	14	-7.14	24	22	24	-8.33
5	74	68	8.82	25	55	60	-8.33
6	32	32	0.00	26	110	121	-9.09
7	43	42	2.38	27	20	25	-20.00
8	17	16	6.25	28	35	30	16.67
9	85	69	23.19	29	55	64	-14.06
10	16	16	0.00	30	77	62	24.19
11	51	46	10.87	31	36	31	16.13
12	20	19	5.26	32	90	90	0.00
13	131	118	11.02	33	38	33	15.15
14	52	48	8.33	34	33	30	10.00
15	52	65	-20.00	35	53	45	17.78
16	28	30	-6.67	36	64	56	14.29
17	52	48	8.33	37	18	17	5.88
18	83	86	-3.49	38	37	36	2.78
19	13	13	0.00	39	50	50	0.00
20	28	27	3.70	40	42	37	13.51

CONCLUSIONS

Pregnenolone can be accurately measured in blood serum by this method which achieved:

- Analytical range from 10 to 500 ng/dL
- Throughputs of 13, 26 or 52 injections per hour from a 1-, 2- or 4-channel system
- Inter- & intra-batch precisions less than 8% and carryover less than 0.2%
- Multi-channeling with other methods utilizing the same HESI source

REFERENCES

1. Kushnir MM, et al. Development and Performance Evaluation of a Tandem Mass Spectrometry Assay for 4 Adrenal Steroids. Clin Chem 2006; 52:1559–1567.

ACKNOWLEDGEMENTS

The authors thank Dr. Lori Vermeulen, Dean of the College of Arts & Sciences of West Chester University of Pennsylvania and Dr. Fred Monson, Director of the Center for Microanalysis & Imaging Research & Training, for the use of space and resources needed for some of this work. We also thank Dr. Stephen Zimmiski, Director, Pharmaceutical Product Development Program, for providing student internship opportunities for this project.

TRADEMARKS/LICENSING

© 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.

For research use only. Not for use in diagnostic procedures.

- Multi-channeling with other methods utilizing the same HESI source

REFERENCES

1. Nelson, RE, et al. Liquid chromatography–tandem mass spectrometry assay for simultaneous measurement of estradiol and estrone in human plasma. Clin Chem, 2004, 50:373-384.
2. Kushnir, MM, et al. High-sensitivity tandem mass spectrometry assay for serum estrone and estradiol. Am J Clin Pathol, 2008, 129:530-539.

ACKNOWLEDGEMENTS

The authors thank Dr. Lori Vermeulen, Dean of the College of Arts & Sciences of West Chester University of Pennsylvania and Dr. Fred Monson, Director of the Center for Microanalysis & Imaging Research & Training, for the use of space and resources needed for some of this work. We also thank Dr. Stephen Zimmiski, Director, Pharmaceutical Product Development Program, for providing student internship opportunities for this project.

For research use only. Not for use in diagnostic procedures.

Find out more at thermofisher.com

ThermoFisher
SCIENTIFIC

batch precisions less than 0.7% and carryover less than 0.2%
For researching with other methods utilizing the same HESI source

and error of estimate = 7.3%.

lab results comparison

Reference Lab	Difference %	Test Sample	Current Method	Reference Lab	Difference %
93	24.73	21	27	30	-10.00
78	8.97	22	49	52	-5.77
73	17.81	23	66	73	-9.59
14	-7.14	24	22	24	-8.33
68	8.82	25	55	60	-8.33
32	0.00	26	110	121	-9.09
42	2.38	27	20	25	-20.00
16	6.25	28	35	30	16.67
69	23.19	29	55	64	-14.06
16	0.00	30	77	62	24.19
46	10.87	31	36	31	16.13
19	5.26	32	90	90	0.00
118	11.02	33	38	33	15.15
48	8.33	34	33	30	10.00
65	-20.00	35	53	45	17.78
30	-6.67	36	64	56	14.29
48	8.33	37	18	17	5.88
86	-3.49	38	37	36	2.78
27	3.70	40	42	37	13.51

ICES

et al. Development and Performance Evaluation of a Tandem
 netry Assay for 4 Adrenal Steroids. Clin Chem 2006; 52:1559–

LEDGEMENTS

ank Dr. Lori Vermeulen, Dean of the College of Arts & Sciences of
 iversity of Pennsylvania and Dr. Fred Monson, Director of the
 analysis & Imaging Research & Training, for the use of space
 eeded for some of this work. We also thank Dr. Stephen
 r, Pharmaceutical Product Development Program, for providing
 p opportunities for this project.

ARKS/LICENSING

Fisher Scientific Inc. All rights reserved. All trademarks are the
 no Fisher Scientific and its subsidiaries. This information is not
 urage use of these products in any manner that might infringe
 roperty rights of others.

ONS

be accurately measured in blood serum by this method

from 10 to 500 ng/dL

13, 26 or 52 injections per hour from a 1-, 2- or 4-channel

batch precisions less than 8% and carryover less than 0.2%
 with other methods utilizing the same HESI source

ICES

al. Development and Performance Evaluation of a Tandem
 etry Assay for 4 Adrenal Steroids. Clin Chem 2006; 52:1559–

EDGEMENTS

For research use only. Not for use in diagnostic procedures.

Find out more at thermofisher.com

