

Development of an LC-MS/MS Method for Measurement of a Steroid Panel in Serum for Clinical Research

Xiaolei Xie, Thermo Fisher Scientific, San Jose, CA, USA

ABSTRACT

Purpose: To develop a simple and fast LC-MS/MS method for quantification of 11 steroids in serum

Methods: A solid phase extraction (SPE) method was developed on a Thermo Scientific™ SOLAµ™ HRP 96-well plate to simultaneously extract 11 steroids from serum (**11-deoxycortisol, 17-OH progesterone, aldosterone, androstenedione, corticosterone, cortisol, cortisone, estradiol, estrone, progesterone and testosterone**). Extracted compounds were separated on a reverse phase column chromatographically followed by analysis on a Thermo Scientific™ TSQ Quantiva™ triple quadrupole mass spectrometer.

Results: The entire SPE process takes less than 20 minutes, and no pre-conditioning, evaporation or reconstitution is required. Comparing to conventional SPE method which involves those steps, our method is simpler and faster. The recovery rate ranged from 42% (aldosterone) to 95% (testosterone). In neat solution, lower limit of quantitation (LOQ) of androstenedione is 1 pg/mL. LOQ of testosterone is 2 pg/mL. LOQ of 11-deoxycortisol, 17-OH progesterone, cortisone, estradiol, estrone, and progesterone is 5 pg/mL. LOQ of aldosterone, corticosterone, and cortisol is 10 pg/mL.

INTRODUCTION

Over the past few years there has been a growing interest to use LC-MS/MS method to measure steroids in serum. LC-MS/MS offers the potential for more reliable measurement compared to other detection systems, such as immunoassays. However, the challenges are time-consuming sample pre-treatment, matrix interference and isobars separation. We developed a simple and fast SPE sample preparation method for an 11-steroid panel and evaluated the analytical performance on an LC-triple quadrupole MS/MS system.

MATERIALS AND METHODS

Sample Preparation

SPE was performed on a SOLAµ HRP 96-well plate. Calibrators/QCs were spiked into neat solution or charcoal stripped serum and mixed with water and methanol. The mixture was loaded directly onto the SPE plate; no preconditioning was required. After washing the plate with 30% methanol, the elution was performed with two volumes of 25 µL of methanol each. Eluates were diluted with 50 µL water. 50 µL of the diluted eluate was injected for LC-MS/MS analysis.

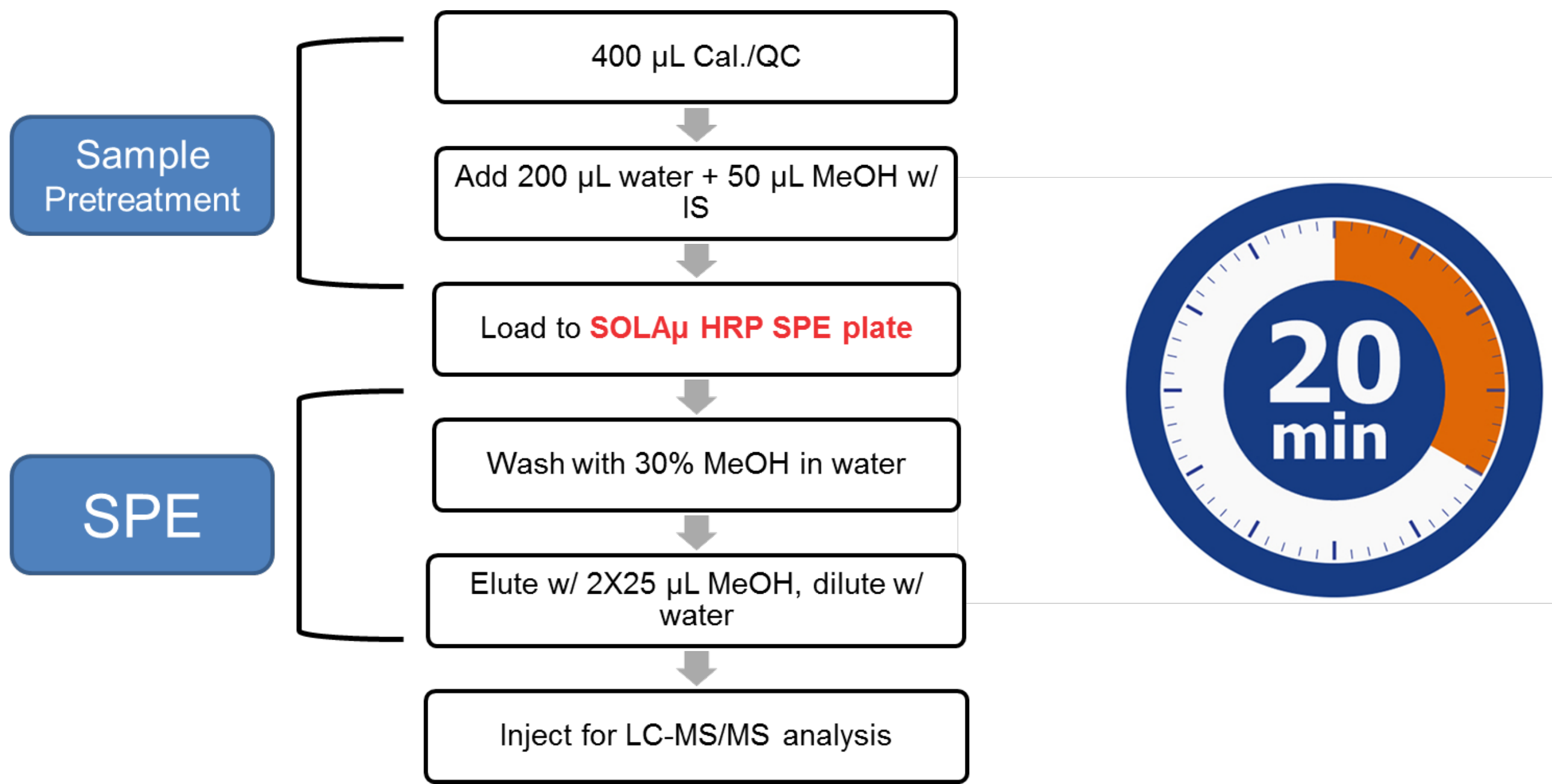


Figure 1. SPE process to extract 11 steroids from plasma.

Test Method(s)



Figure 2. Thermo Scientific™ UltiMate™ 3000 HPLC and TSQ Quantiva MS with heated ESI source

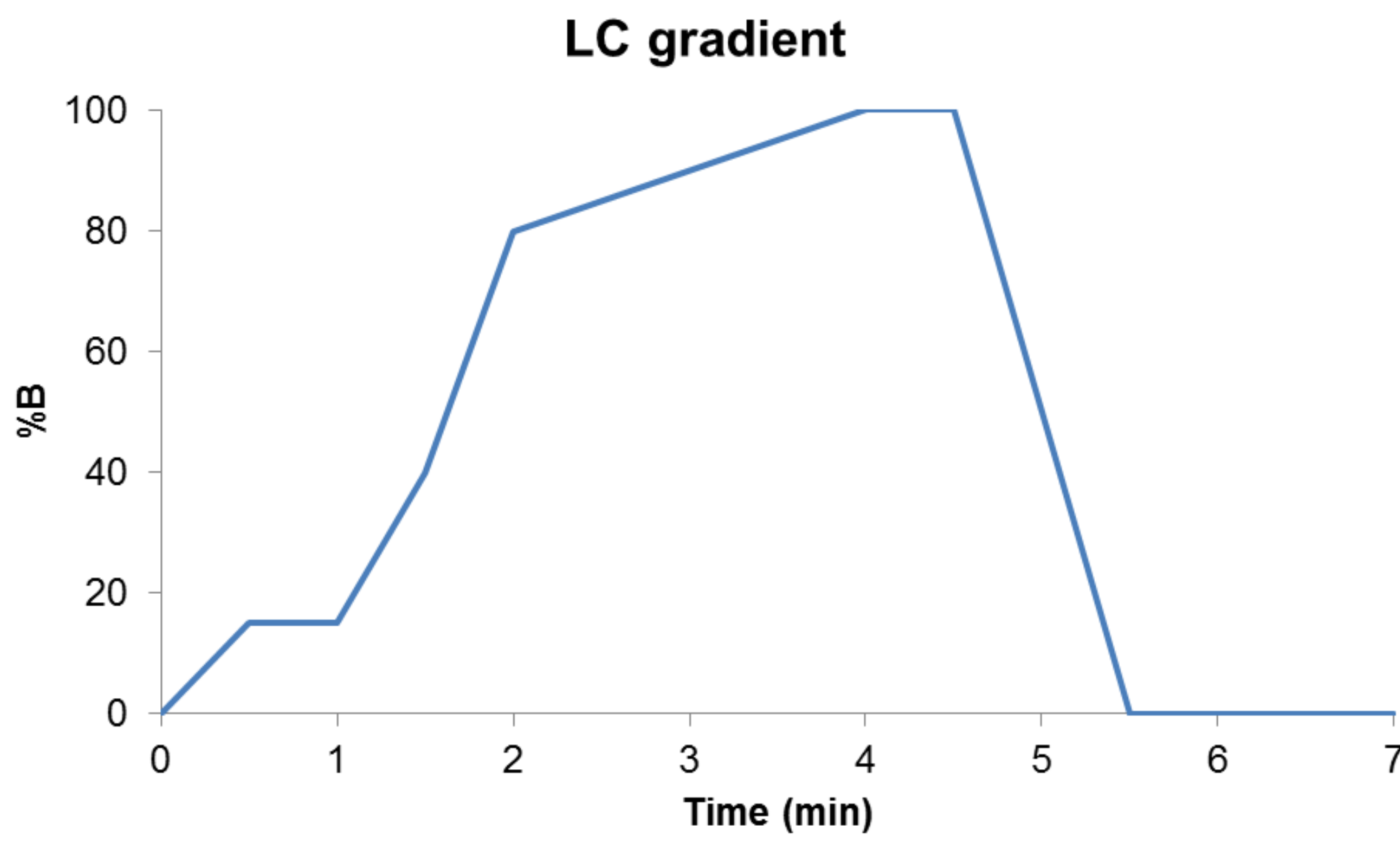


Figure 3. LC gradient (Column: Thermo Scientific™ Accucore™ RP-MS, 2.6 µm, 100 X 2.1 mm)

Compound	RT (min)	RT Window (min)	Polarity	Precursor (m/z)	Product (m/z)	Collision Energy (V)	RF Lens (V)
11-deoxycortisol	3.2	1	Positive	347.3	97	25	72
11-deoxycortisol	3.2	1	Positive	347.3	109	28	72
11-Deoxycortisol-d5	3.2	1	Positive	352.3	100	25	71
11-Deoxycortisol-d5	3.2	1	Positive	352.3	113	30	71
17-OH-P	3.4	1	Positive	331.2	97	24	69
17-OH-P	3.4	1	Positive	331.2	109.1	27	69
17OH-P-d8	3.4	1	Positive	339.3	100	26	70
17OH-P-d8	3.4	1	Positive	339.3	113	29	70
Aldosterone	3	1	Negative	359.3	189.1	17	69
Aldosterone	3	1	Negative	359.3	331.1	15	69
Androstenedione	3.3	1	Positive	287.2	97	23	64
Androstenedione	3.3	1	Positive	287.2	109.1	25	64
Androstenedione- ¹³ C3	3.3	1	Positive	290.2	100.1	22	65
Androstenedione- ¹³ C3	3.3	1	Positive	290.2	112	25	65
Corticosterone	3.2	1	Positive	347.3	121.1	25	68
Corticosterone	3.2	1	Positive	347.3	147.1	26	68
Cortisol	3	1	Positive	363.2	121.1	26	71
Cortisol	3	1	Positive	363.2	309.2	20	71
Cortisol-d4	3	1	Positive	367.3	121.1	26	72
Cortisol-d4	3	1	Positive	367.3	175.1	27	72
Cortisone	3	1	Positive	361.2	121.1	31	78
Cortisone	3	1	Positive	361.2	163	24	78
Estradiol	3.3	1	Negative	271.2	145	41	108
Estradiol	3.3	1	Negative	271.2	183	40	108
Estradiol-d5	3.3	1	Negative	276.1	147	40	115
Estradiol-d5	3.3	1	Negative	276.1	187	41	115
Estrone	3.4	1	Negative	269.2	145.1	40	90
Estrone	3.4	1	Negative	269.2	183.1	40	90
Estrone- ¹³ C3	3.3	1	Negative	272.2	146	54	106
Estrone- ¹³ C3	3.3	1	Negative	272.2	148.1	39	106
Progesterone	3.7	1	Positive	315.3	97.1	22	68
Progesterone	3.7	1	Positive	315.3	109.1	25	68
Progesterone-d9	3.7	1	Positive	324.5	100.1	22	67
Progesterone-d9	3.7	1	Positive	324.5	113	27	67
Testosterone	3.4	1	Positive	289.3	97	22	64
Testosterone	3.4	1	Positive	289.3	109.1	25	64
Testosterone- ¹³ C3	3.4	1	Positive	292.3	100	23	66
Testosterone- ¹³ C3	3.4	1	Positive	292.3	112	25	66

Figure 4. SRM transitions (two for each analyte/IS)

RESULTS

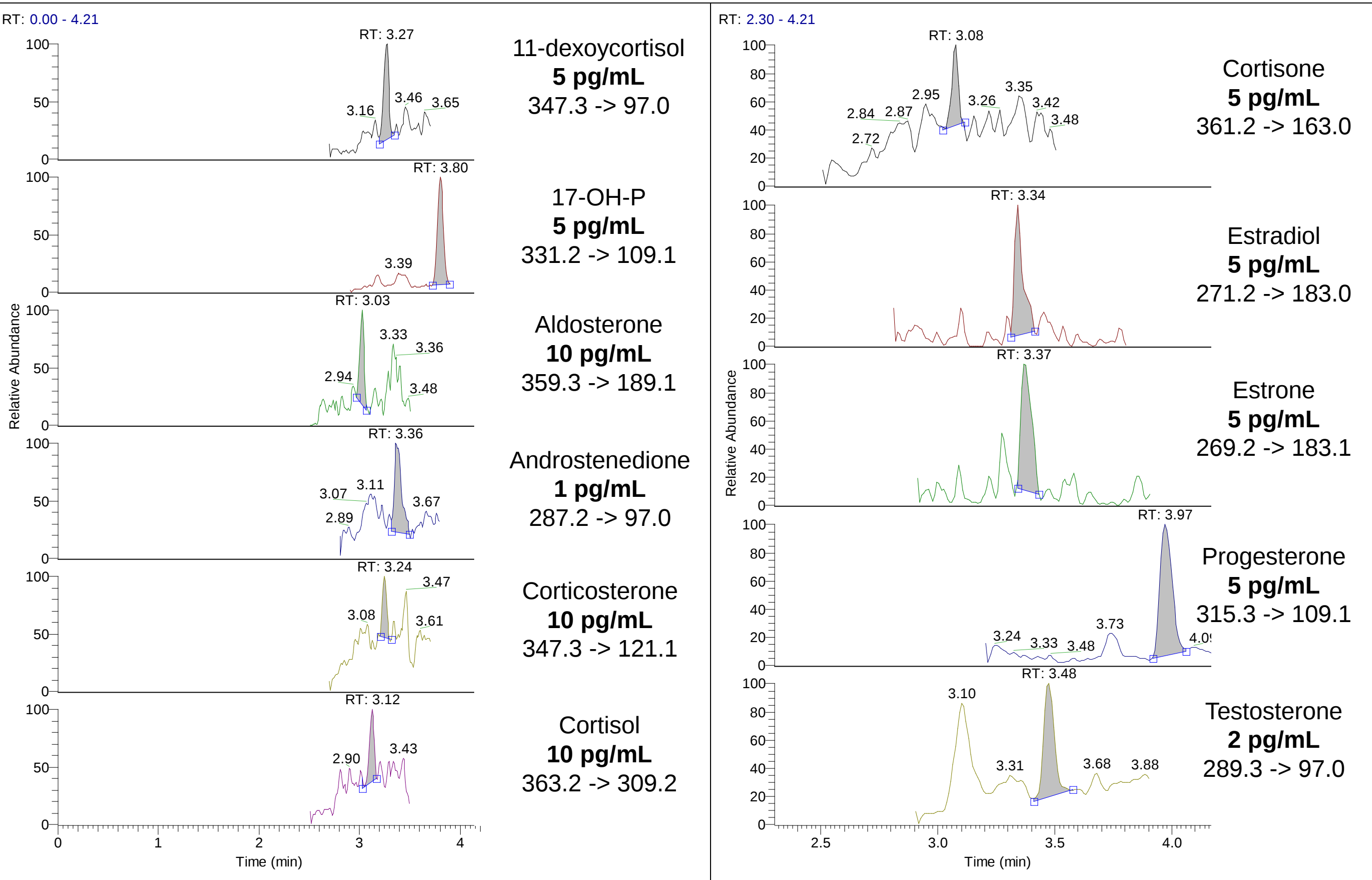


Figure 5. Chromatograms on the lower limit of quantitation level in neat solution

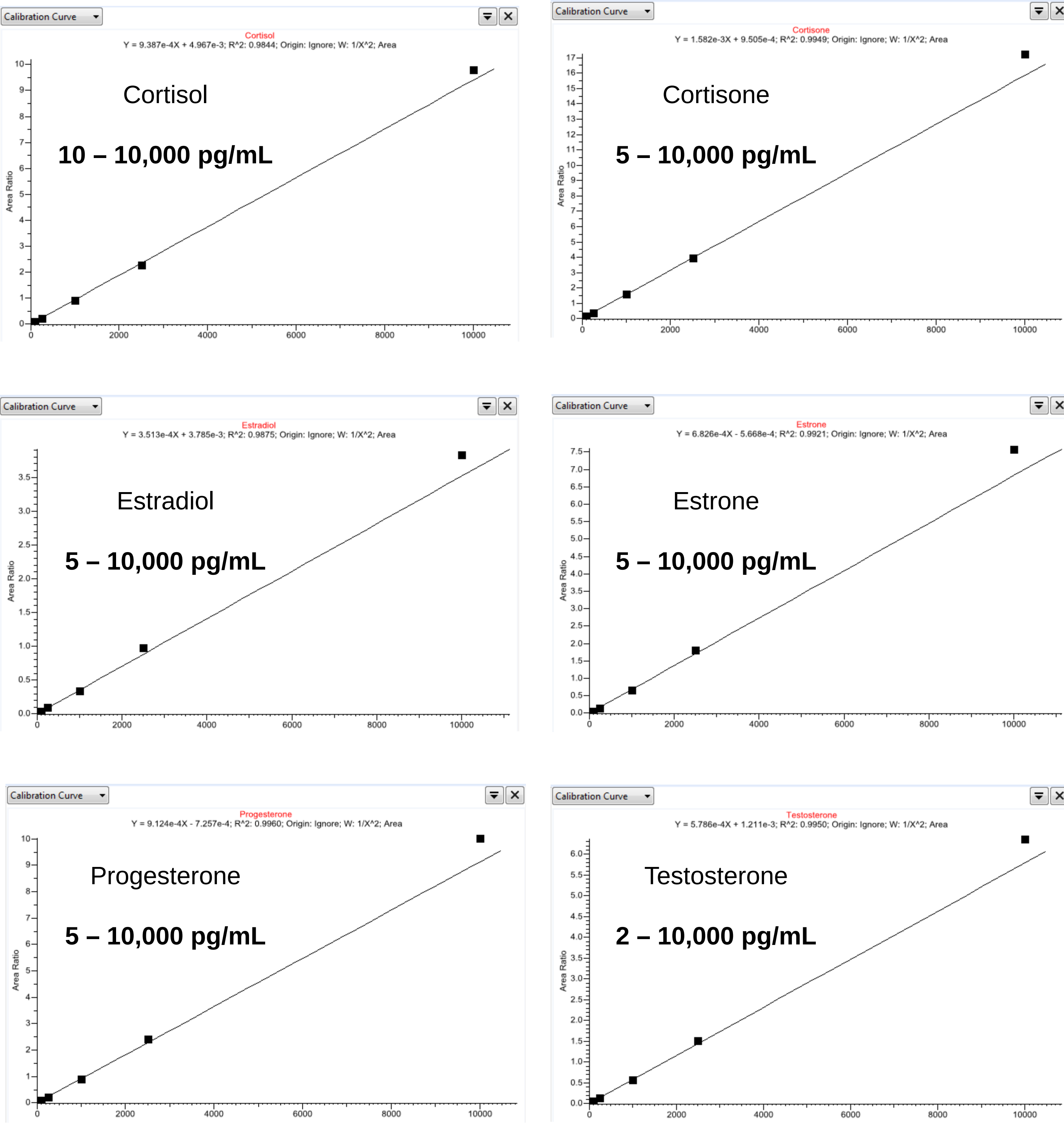
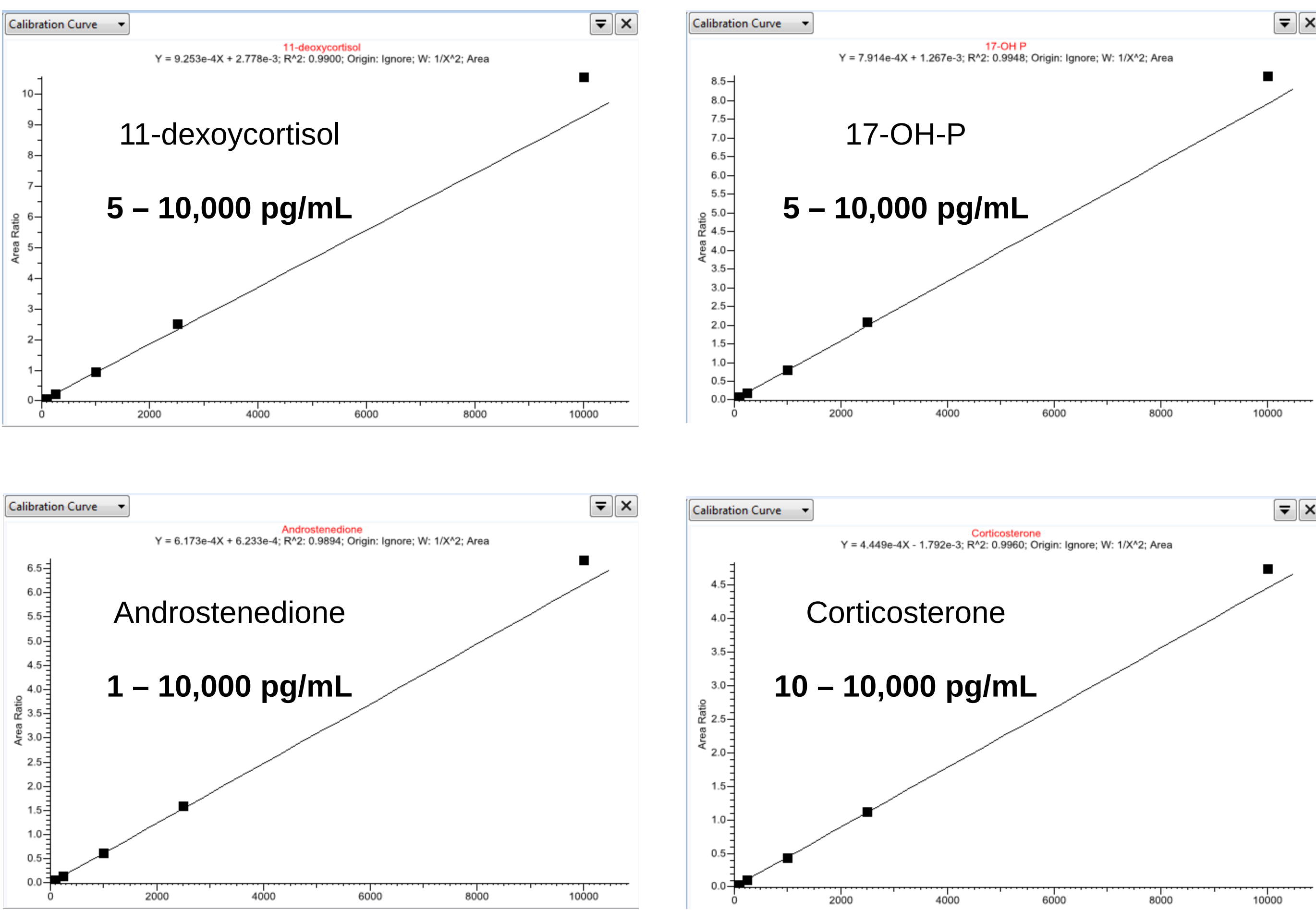


Figure 6. Linearity range in neat solution

Compound	SPE Recovery Rate (%)	Lower Limit of Quantitation (pg/mL)
11-deoxycortisol	63	5
11-Deoxycortisol-d5	64	
17-OH-P	84	5
17OH-P-d8	83	
Aldosterone	42	10
Androstenedione	93	1
Androstenedione- ¹³ C3	95	
Corticosterone	67	10
Cortisol	68	10
Cortisol-d4	70	
Cortisone	71	5
Estradiol	87	5
Estradiol-d5	88	
Estrone	89	5
Estrone- ¹³ C3	89	
Progesterone	84	5
Progesterone-d9	86	
Testosterone	95	2
Testosterone- ¹³ C3	97	

Figure 7. Recovery rate and lower limit of quantitation in neat solution

CONCLUSIONS

- A simple and fast SPE method (~20 min) was develop to simultaneously extract 11 steroids from serum with recovery rate ranged from 42% to 95%.
- A sensitive LC-MS/MS analytical method was demonstrated to analyze 11 steroids with lower limit of quantitation ranged from 1 pg/mL to 10 pg/mL in neat solution.

For Research Use Only. Not for use in diagnostic procedures.

TRADEMARKS/LICENSING

© 2017 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.