



Bond Elut Plexa Sample Preparation for LC/MS/MS Determination of Hormones in Pork

Application Note

Food Testing & Agriculture

Author

Chen-Hao (Andy) Zhai
Agilent Technologies Shanghai Ltd.

Introduction

Food safety is increasingly an important concern of people worldwide, because many chemicals added to food create potential hazards to human health. Hormones are a common food additive. Long term consumption of glucocorticoids can lead to hyperglycemia, hyperglycemia, osteoporosis, birth defects, and immune function decline. Other hormones, such as estrogen, androgen, and progesterone, are carcinogenic and can lead to breast cancer, ovarian cancer, and cell carcinoma. Many countries' regulations clearly define residual limits for these compounds in food. Maximum residue limits for these compounds can vary worldwide, but are generally in the low ppb concentration range. At these concentrations, the analysis of hormones in products such as meat is often very challenging due to the complexity of the sample.

In this application note, pork was prepared and analyzed for hormones at the ppb level, using a straightforward SPE methodology with Agilent Bond Elut Plexa polymeric SPE, efficient separation with an Agilent Poroshell 120 LC column, and sensitive detection with an Agilent 6460 Triple Quadrupole LC/MS system.



Agilent Technologies

Materials and Methods

LC conditions

| | |
|-------------------|--|
| Columns: | Agilent Poroshell 120 EC-C18, 2.1 × 50 mm, 2.7 μm (p/n 699775-902) |
| Eluent: | A, water; B, acetonitrile |
| Injection volume: | 10 μL |
| Flow rate: | 0.4 mL/min |
| Gradient: | 20% B, linear to 40% B in 5 min, linear to 90% B in 3 min, hold for 10 min |
| Temperature: | Ambient |
| Sample vials: | Agilent Certified Vials (p/n 5183-2072) |
| System: | Agilent 1260 Infinity LC |

MS conditions

| | |
|---------------------|--------------------------------------|
| Ionization mode: | ESI + Agilent Jet Stream |
| Gas temperature: | 325 °C |
| Gas flow: | 10 L/min |
| Nebulizer: | 50 psi |
| Sheath temperature: | 400 °C |
| Sheath gas flow: | 12 L/min |
| Capillary: | 4,500 V (ESI+), 3,500 V (ESI-) |
| Nozzle voltage: | 1,500 V (ESI+), 1,500 V (ESI-) |
| System: | Agilent 6460 Triple Quadrupole LC/MS |

The MRM transitions, fragmentors, and collision energies optimized for the hormones in this study are shown in Table 1.

Table 1. MRM transitions and other conditions for hormones.

| Compound name | Precursor ion | Product ion | Fragmentor | CE | Polarity |
|-----------------------------|---------------|-------------|------------|----|----------|
| Triamcinolone | 435.4 | 415.2 | 100 | 2 | Positive |
| | 435.4 | 397.2 | 100 | 5 | Positive |
| Nandrolone phenylpropionate | 407.4 | 257.2 | 150 | 10 | Positive |
| | 407.4 | 105.1 | 150 | 25 | Positive |
| Dexamethasone | 393.3 | 373.2 | 100 | 2 | Positive |
| | 393.3 | 355.1 | 100 | 2 | Positive |
| Methylprednisolone | 375.4 | 357.2 | 100 | 2 | Positive |
| | 375.4 | 161.2 | 100 | 15 | Positive |
| Hydrocortisone | 363.4 | 327.2 | 125 | 8 | Positive |
| | 363.4 | 121.1 | 125 | 20 | Positive |
| Prednisolone | 361.4 | 343.2 | 100 | 2 | Positive |
| | 361.4 | 147.2 | 100 | 20 | Positive |
| Prednisone | 359.3 | 341.2 | 125 | 2 | Positive |
| | 359.3 | 147.1 | 125 | 20 | Positive |
| Methyltestosterone | 303.4 | 109.1 | 125 | 25 | Positive |
| | 303.4 | 97.2 | 125 | 25 | Positive |
| Estriol | 287.3 | 171.2 | 125 | 30 | Negative |
| | 287.3 | 145.3 | 125 | 40 | Negative |
| Trenbolone | 271.4 | 253.2 | 150 | 15 | Positive |
| | 271.4 | 199.1 | 150 | 20 | Positive |
| Hexestrol | 269.3 | 134.1 | 125 | 4 | Negative |
| | 269.3 | 119 | 125 | 35 | Negative |
| Diethylstilbestrol | 267.3 | 251.2 | 150 | 15 | Negative |
| | 267.3 | 237.2 | 150 | 20 | Negative |

Sample preparation

To pretreat the sample, 5 g ground pork was placed into a 50 mL centrifuge tube, followed by 5 mL methanol and 20 mL water. The tube was then vortexed vigorously for 1 minute. Next, the tube was centrifuged for 5 minutes at 5,000 rpm at 4 °C, and all of the supernatant was transferred into another tube for SPE cleanup.

To extract the sample, the procedure shown in Figure 1 was used. Agilent Bond Elut Plexa cartridges (60 mg, 3 mL, p/n 12109603) were preconditioned with 3 mL methanol then 3 mL water. The extract (equivalent to 5 g sample) was passed through the cartridge at a rate of 1 mL/min. After the sample passed through completely, the cartridge was washed with 3 mL 35% methanol in water, and the entire effluent was discarded. The cartridge was dried under negative pressure (below 2.0 kPa) for 5 minutes. The sample was eluted with 5 mL of methanol. The eluate was collected and dried under nitrogen below 40 °C. The sample residue was then dissolved and brought to a constant volume of 1.0 mL using 20% methanol in water (v:v), filtered through a 0.2 μm filter membrane (Agilent Captiva Polyethersulfone, p/n 5190-5096), and analyzed by LC/MS/MS.

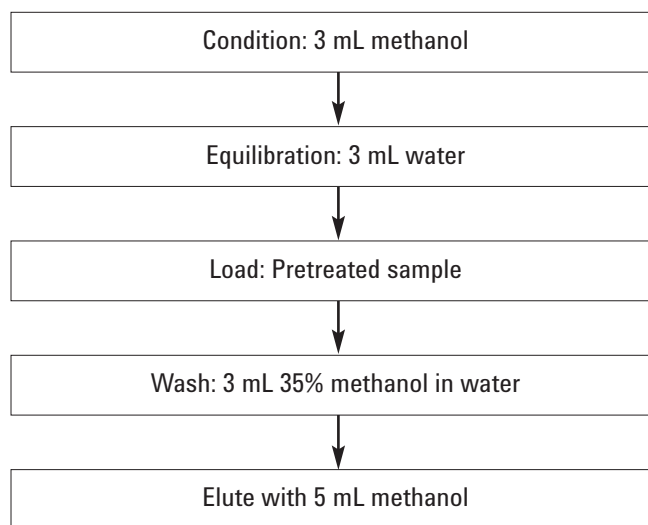


Figure 1. The SPE procedure for a pork sample.

Results and Discussion

Recovery was measured for each hormone at both low and high concentration levels (Table 2). Recovery was calculated by comparing the MRM peak area for samples spiked prior to SPE extraction, with the MRM peak area for samples spiked after SPE extraction (post spiked samples). Figure 2 shows a chromatogram obtained from the analysis of pork blank sample spiked with low levels of hormones. Figure 3 is the sample blank.

Table 2. Extraction recoveries of hormones from pork with SPE.

| Hormone | % Recovery (%RSD) n = 6 low level (1 ppb) | % Recovery (%RSD) n = 6 high level (10 ppb) |
|-----------------------------|---|---|
| Triamcinolone acetonide | 67.2 (5.4) | 78.3 (7.9) |
| Nandrolone phenylpropionate | 66.7 (9.8) | 70.3 (6.3) |
| Dexamethasone | 86.9 (4.5) | 91.4 (3.9) |
| Methylprednisolone | 95.8 (5.2) | 94.3 (8.3) |
| Hydrocortisone | 102.3 (7.1) | 98.7 (4.4) |
| Prednisolone | 82.8 (3.7) | 75.3 (3.9) |
| Prednisone | 77.2 (4.3) | 86.9 (1.4) |
| Methyltestosterone | 80.1 (1.7) | 87.6 (2.2) |
| Estriol | 53.7 (4.9) | 67.5 (8.7) |
| Trenbolone | 98.0 (3.5) | 103.5 (3.2) |
| Hexestrol | 34.7 (10.7) | 46.8 (8.4) |
| Diethylstilbestrol | 42.3 (5.1) | 36.2 (5.8) |

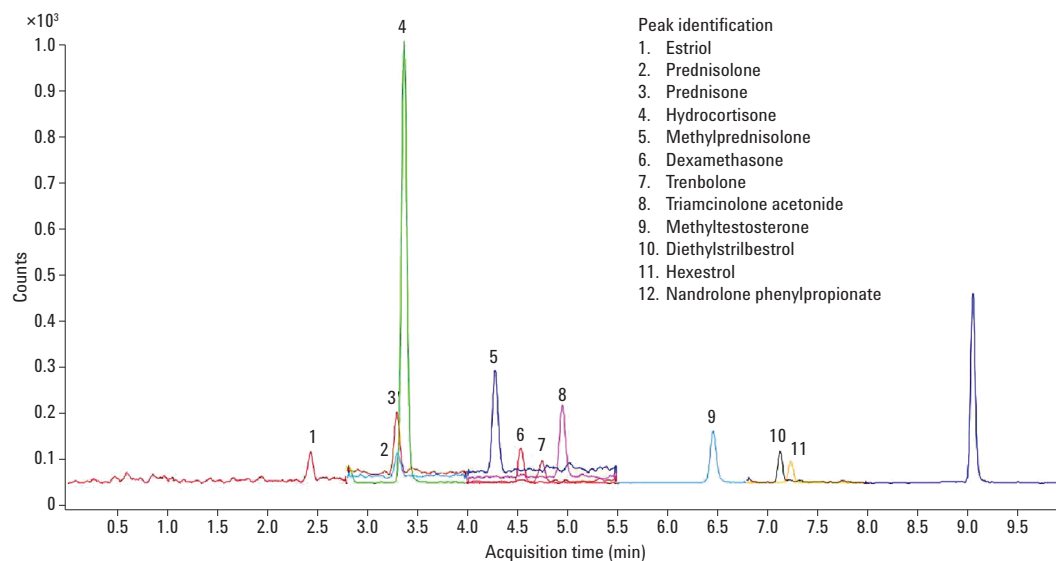


Figure 2. Chromatogram of hormones obtained from pork spiked with a low level sample,

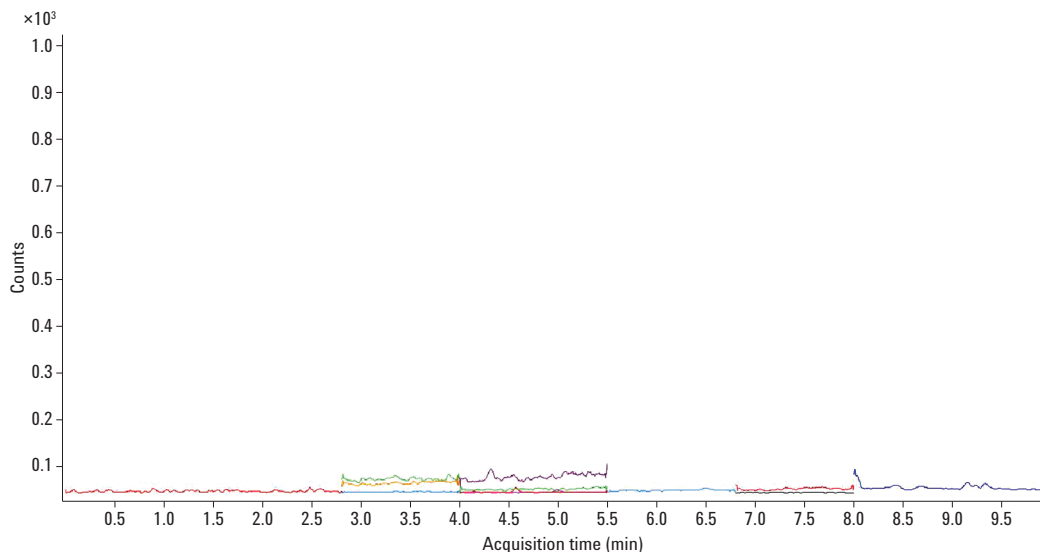


Figure 3. Chromatogram of hormones obtained from a pork blank sample.

Conclusions

Good recovery and reproducibility were obtained with Agilent Bond Elut Plexa SPE for most hormones in a pork matrix. However, for hexestrol and diethylstilbestrol, the results were influenced by matrix effects, and so isotope internal standards should be used for these two compounds to achieve better recoveries.

Bond Elut Plexa SPE combined with LC/MS/MS enables sensitive quantitation of hormones in meat samples at low ppb concentrations.

For More Information

These data represent typical results. For more information on our products and services, visit our Web site at www.agilent.com/chem.

www.agilent.com/chem

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc., 2013
 Printed in the USA
 November 21, 2013
 5991-3660EN



Agilent Technologies