Application Note: 482

Detection of Glucocorticoid Residues in Animalderived Food by HPLC-MS/MS

Chen Dongmei¹, Fei Liu², and Charles Yang³

¹National Reference Laboratory of Veterinary Drug Residues, Huazhong Agricultural University, Wuhan, China;

²Thermo Fisher Scientific, Shanghai, China; ³Thermo Fisher Scientific, San Jose, CA, USA

Key Words

- Food residue analysis
- Veterinary drugs
- TSQ Quantum
- SRM (Selective Reaction Monitoring)

Introduction

Glucocorticoids are a class of steroid hormones, produced by the adrenal cortex, that affect metabolism and have immunosuppressive and anti-inflammatory properties. The common chemical structure of these compounds is shown in Figure 1. The chemical structures of eight glucocorticoid drugs are listed in Table 1.

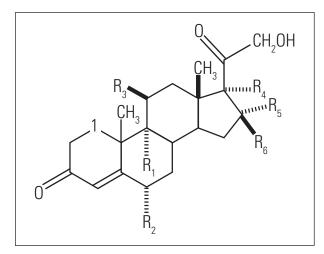


Figure 1. Common chemical structure of glucocorticoids

Table 1. Glucocorticoid drugs and their functional groups

	R ₁	R_2	R_3	R_4	R_5	R_6
Dexamethasone	-F	-H	-0H	-OH	-CH ₃	-H
Betamethasone	-F	-H	-0H	-0H	-H	-CH ₃
Prednisone	-H	-H	=0	-0H	-H	-H
Prednisolone	-H	-H	-0H	-0H	-H	-H
Methylprednisolone	-H	-CH ₃	-0H	-OH	-H	-H
Beclomethasone	-CI	-H	-0H	-OH	-H	-CH ₃
Hydrocortisone	-H	-H	-0H	-OH	-H	-H
Fludrocortisone	-F	-H	-OH	-OH	-H	-H

Synthetic glucocorticoids are often administered to livestock to treat inflammatory diseases, disorders of the musculoskeletal system, and other diseases common to farm animals. They can also be used to improve feed intake and stimulate growth in livestock. However, glucocorticoid residues in animal-derived food may be

harmful to humans. Therefore, the use of certain substances with a hormonal or thyrostatic action as growth-promoting additives has been banned by the European Commission. Dexamethasone, betamethasone, prednisolone, and methylprednisolone are the only glucocorticoids permitted for therapeutic purposes in livestock. Maximum residue levels (MRLs) have been established for these compounds in tissues and in milk intended for human consumption by the EC2-5 and other countries. Table 2 lists the MRLs for glucocorticoids that were set by the European Union. In China, the Institute for the Control of Agrochemicals, Ministry of Agriculture, has established MRLs of 0.75 $\mu g/kg$ for dexamethasone in the muscles, livers, and kidneys of cattle, pigs, and horses and 10 $\mu g/kg$ for hydrocortisone in milk.

Table 2. European Union MRLs for glucocorticoids

Substance	Animal Species	MRL	Target Tissues
Dexamethasone	Cattle	0.3 μg/kg	Milk
	Cattle, Pig, Horse	0.75 μg/kg 2.0 μg/kg 0.75 μg/kg	Muscle Liver Kidney
Betamethasone	Cattle	0.75 µg/kg 2.0 µg/kg 0.75 µg/kg 0.3 µg/kg	Muscle Liver Kidney Milk
	Pig	0.75 μg/kg 2.0 μg/kg 0.75 μg/kg	Muscle Liver Kidney
Prednisolone	Cattle	4 μg/kg 4 μg/kg 10 μg/kg 10 μg/kg 6 μg/kg	Muscle Fat Liver Kidney Milk
Methylprednisolone	Cattle	10 µg/kg 10 µg/kg 10 µg/kg 10 µg/kg	Muscle Fat Liver Kidney

Goal

To develop an effective LC-MS/MS method to detect eight gluococorticoid residues in the liver and muscles of pig, cattle, and lamb, as well as in chicken, eggs and milk.



Experimental Conditions

Sample Preparation

A 2 g tissue sample was weighed into a 50 mL centrifuge tube and 15 mL of ethyl acetate was added. The sample was vortexed for 3 minutes and then centrifuged at 4000 rpm for 10 minutes. The supernatant was transferred into another 50 mL centrifuge tube. The sample was extracted again with 10 mL sodium hydroxide (0.1 M) and 15 mL ethyl acetate. The two supernatants were combined and evaporated to dryness at 40 °C. The residue was dissolved in 1.0 mL of ethyl acetate and 5.0 mL of n-hexane before being loaded to the silica extraction cartridges (SPE column). The SPE column was previously conditioned with 5 mL of n-hexane. The column was washed with 5 mL of n-hexane. After drying, the analytes were eluted with 5 mL n-hexane:acetone (40:60, v/v) and evaporated to dryness under nitrogen at 40 °C. The residues were reconstituted in 1.0 mL of 20% acetonitrile aqueous solution. The resulting solutions were vortexed for 1 minute and then centrifuged at 10,000 rpm for 10 minutes. The upper clear solutions were transferred to another sample vial for LC-MS/MS analysis.

The glucocorticoid standards were obtained from Sigma (Sigma Chemical Company, St. Louis, MO). All other chemicals were HPLC grade.

LC

HPLC analysis was performed using the Thermo Scientific Surveyor HPLC system. Each 10 μL of sample was injected into a Thermo Scientific Hypersil GOLD column (150 x 2.1 mm, 5 μm). A gradient LC method was used and mobile phases A (water with 0.1% formic acid) and B (acetonitrile) were at a flow rate of 250 $\mu L/min$ together. Table 3 illustrates the gradient LC method.

Table 3. Gradient details

Retention	Time (min)	A (%)	B (%)
	0	70	30
	18	40	60
2	23	40	60
2	3.1	70	30
2	28	70	30

MS

MS analysis was performed on a Thermo Scientific TSQ Quantum triple stage quadrupole mass spectrometer with an electrospray ionization (ESI) probe. The MS conditions were as follows:

Ion source polarity: Negative ion mode

Spray voltage: 3500 V
Sheath gas pressure (N₂): 40 units
Auxiliary gas pressure (N₂): 5 units
Capillary temperature: 350 °C
Collision gas pressure (Ar): 1.5 mTorr
Q1/Q3 Peak Resolution: 0.7 FWHM

The SRM transitions that were monitored are summarized in Table 4.

Table 4. SRM transitions

	SRM Transitions (m/z)		
Drug	Quantitative	Qualitative	
Prednisone	403.0 → 326.9	$403.0 \rightarrow 357.2$	
Prednisolone	$405.0 \rightarrow 329.0$	$405.0 \rightarrow 358.8$	
Hydrocortisone	407.1 → 331.1	407.1 → 361.1	
Methylprednisolone	419.0 → 342.9	$419.0 \rightarrow 373.1$	
Dexamethasone	437.0 → 361.0	$437.0 \rightarrow 391.5$	
Betamethasone	$437.0 \rightarrow 361.0$	$437.0 \rightarrow 391.5$	
Beclomethasone	$453.0 \rightarrow 406.9$	$453.0 \rightarrow 376.8$	
Fludrocortisone acetate	467.1 → 420.8	467.1 → 349.2	

Results and Discussion

Figure 2 displays the SRM chromatograms for the eight glucocorticoids. The limit of quantitation (LOQ) for prednisone, prednisolone, dexamethasone, betamethasone, and methylprednisolone in milk is 0.1 μ g/L. The LOQ of these residues in muscle, eggs and livers is 0.2 μ g/kg. For fludrocortisone acetate and beclomethasone, the LOQ is 0.2 μ g/L in milk, and 0.2 μ g/kg in eggs, and livers. For hydrocortisone, the LOQ in milk is 0.2 μ g/L and in eggs and livers is 0.5 μ g/kg. These LOQs easily meet the specified MRLs of the European Union and China.

The extraction recovery of glucocorticoids from the muscle and livers of cattle, pigs, and lamb; chicken; eggs; and milk is between 60% and 110%. This achieves the minimum detection requirements. Thus, the qualification method is accurate and reproducible.

The method validation data is summarized in Table 5. Linearity of the method was assumed because the R^2 values were greater than 0.99 for the linear regression equations (1/x weighted).

Table 5. Linearity and dynamic range

Drug	Dynamic Range (μg/L, μg/kg)	Equation	R²
Prednisone	0.2 ~ 50	y = 106952 x + 5288	0.9992
Prednisolone	0.2 ~ 50	y = 100276 x - 2420.7	0.9954
Hydrocortisone	0.8 ~ 200	y = 85500 x + 6127.5	0.9947
Methylprednisolone	0.2 ~ 50	y = 98154 x - 5291.5	0.9995
Dexamethasone	0.2 ~ 50	y = 106501 x - 8596.1	0.9991
Betamethasone	0.2 ~ 50	y = 116052 x - 17056	0.9986
Beclomethasone	0.4 ~ 100	y = 52195 x - 6621.2	0.9911
Fludrocortisone acetate	e 0.4 ~ 100	y = 33779 x - 2193.9	0.9992

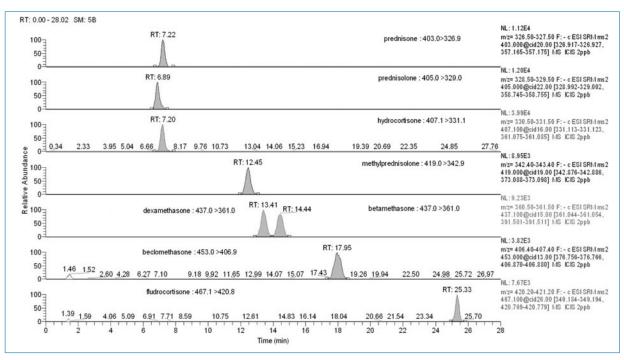


Figure 2. SRM chromatograms for the eight glucocorticoids.

Conclusion

The LC-MS/MS method described above is able to detect glucocorticoid residues with high recoveries and enable their accurate quantification. The sensitivity, extraction recovery, and reproducibility of this method meet international regulation and detection requirements.

References

- European Community, Council Directive 96/22/EC concerning the prohibition on the use in stockfarming of certain substances having a hormonal or thyrostatic action and of β-agonists, (Official Journal of the European Communities Official Journal L 125, 23 May 1996, pp. 3-9).
- European Community, Commission Regulation (EC) No 508/1999 of 4
 March 1999 amending Annexes I to IV of Council Regulation (EEC) No
 2377/90 laying down a Community procedure for the establishment of
 maximum residue limits of veterinary medicinal products in foodstuffs of
 animal origin, (Official Journal of the European Communities Official
 Journal L 060, of 9 March 1999, pp. 16-52).
- 3. European Community, Commission Regulation (EC) No 2593/1999 of 8 December 1999 amending Annexes I, II and III of Council Regulation (EEC) No 2377/90 laying down a Community procedure for the establishment of maximum residue limits of veterinary medicinal products in foodstuffs of animal origin, (Official Journal of the European Communities Official Journal L 315, of 8 December 1999, pp. 26-31).
- 4. European Community, Commission Regulation (EC) No 2535/2000 of 17 November 2000 amending Annex I of Council Regulation (EEC) No 2377/90 laying down a Community procedure for the establishment of maximum residue limits of veterinary medicinal products in foodstuffs of animal origin, (Official Journal of the European Communities Official Journal L 291, of 17 November 2000, pp. 9-10).
- 5. European Community, Commission Regulation (EC) No 77/2002 of 17 January 2002 amending Annexes I and III to Council Regulation (EEC) No 2377/90 laying down a Community procedure for the establishment of maximum residue limits of veterinary medicinal products in foodstuffs of animal origin, (Official Journal of the European Communities Official Journal L 16, of 17 January 2002, pp. 9-11).

In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.

Africa-Other

Australia +61 2 8844 9500

Austria

+43 1 333 50 34 0

Belgium +32 2 482 30 30

Canada +1 800 530 8447

China +86 10 841<u>9 3588</u>

Denmark +45 70 23 62 60

Europe-Other

Finland/Norway/

Sweden +46 8 556 468 00

France +33 1 60 92 48 00

Germany +49 6103 408 1014

India +91 22 6742 9434

Italy +39 02 950 591

Japan +81 45 453 9100

Latin America +1 608 276 5659

Middle East +43 1 333 50 34 0

Netherlands +31 76 579 55 55

South Africa

Spain +34 914 845 965

Switzerland

UK

+44 1442 233555 USA

+1 800 532 4752

www.thermo.com

Legal Notices

©2009 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific Inc. 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.

View additional Thermo Scientific LC/MS application notes at: www.thermo.com/appnotes



San Jose, CA USA is ISO Certified.

AN62970 E 10/09S

