



# FAST, WIDE-RANGE SCREENING OF BANNED OR NOT REGULATED VETERINARY DRUGS IN URINE BY LIQUID CHROMATOGRAPHY COUPLED TO HIGH RESOLUTION MASS SPECTROMETRY

León N.<sup>1</sup>, Yusá V.<sup>1</sup>, Roca M.<sup>1</sup>, Igualada C.<sup>1</sup>, Martins C.<sup>2</sup>

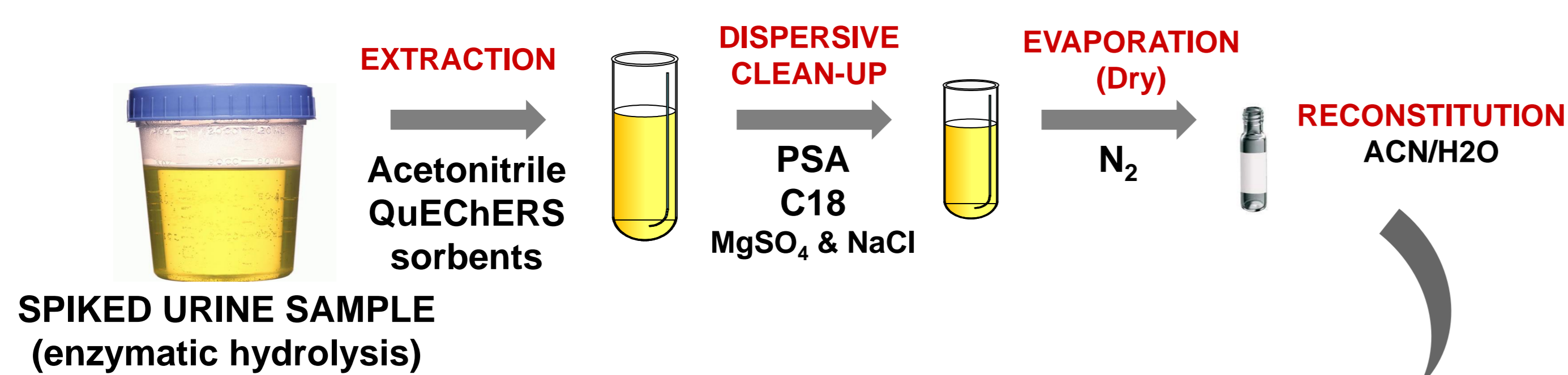
<sup>1</sup>Center for Public Health Research (CSISP), Valencia (Spain), <sup>2</sup>Thermo Fisher Scientific, Barcelona (Spain)

## INTRODUCTION

- ✓ The use of pharmacologically active substances such as hormones, thyrostats, many  $\beta$ -agonists and other drugs, is prohibited in stock farming
- ✓ Low reporting levels in urine must be analytically achieved to successfully monitor their administration
- ✓ LC-MS/MS is the methodology of choice for confirmatory purposes but for multi-class and multi-component analysis this approach requires extensive compound-dependent parameter optimization
- ✓ Full Scan approaches using high resolution mass spectrometry (HRMS) can provide a rapid wide-range screening of veterinary drugs overcoming the limitations of LC-MS/MS analysis

➔ The aim of this study was to develop a UHPLC-HRMS method for the multi-residue screening of around 100 veterinary drugs banned or not regulated in bovine urine

## EXPERIMENTAL



### Chromatographic conditions

System: Autosample: Accela Open AS

UHPLC: Accela 1250 pump

Mobile phase: Eluent A: HAc 0.1%

Eluent B: Acetonitrile 0.1% HAc

Column: Hypersil Gold aQ (100 x 2.1 mm, 1.9  $\mu$ )

Elution mode: Gradient

Flow rate: 400  $\mu$ L/min

Injection volume: 10  $\mu$ L

Time (min)	0	5	8	8.5	10	10.5*
% A	100	50	50	5	5	100
% B	0	50	50	95	95	0

\* Equilibration time: 3min

### Mass conditions

System: Source: HESI II

Analyzer: Orbitrap Exactive™



Spray voltage: 3 kV

Sheath gas flow rate: 8

Auxiliary gas flow rate: 0

Skimmer voltage: 38 V

Heater temperature: 30 °C

Capillary temperature: 250 °C

Tube lens voltage: 120 V

Automatic gain control : 10<sup>6</sup>(Balanced)

Resolving power (FWHM): 50.000 (High)

Mass range (m/z): 150-2000

### ➔ Acquisition functions

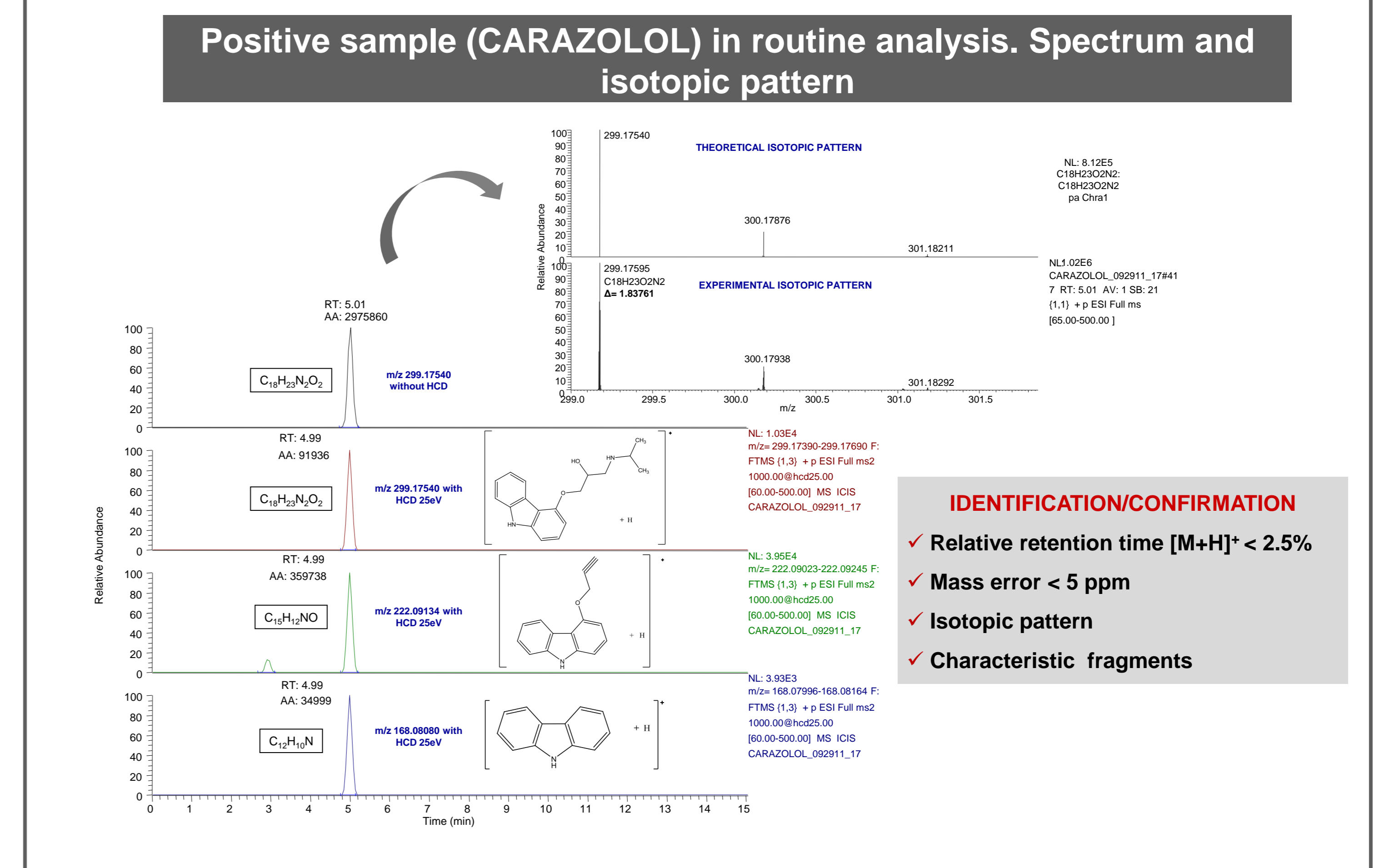
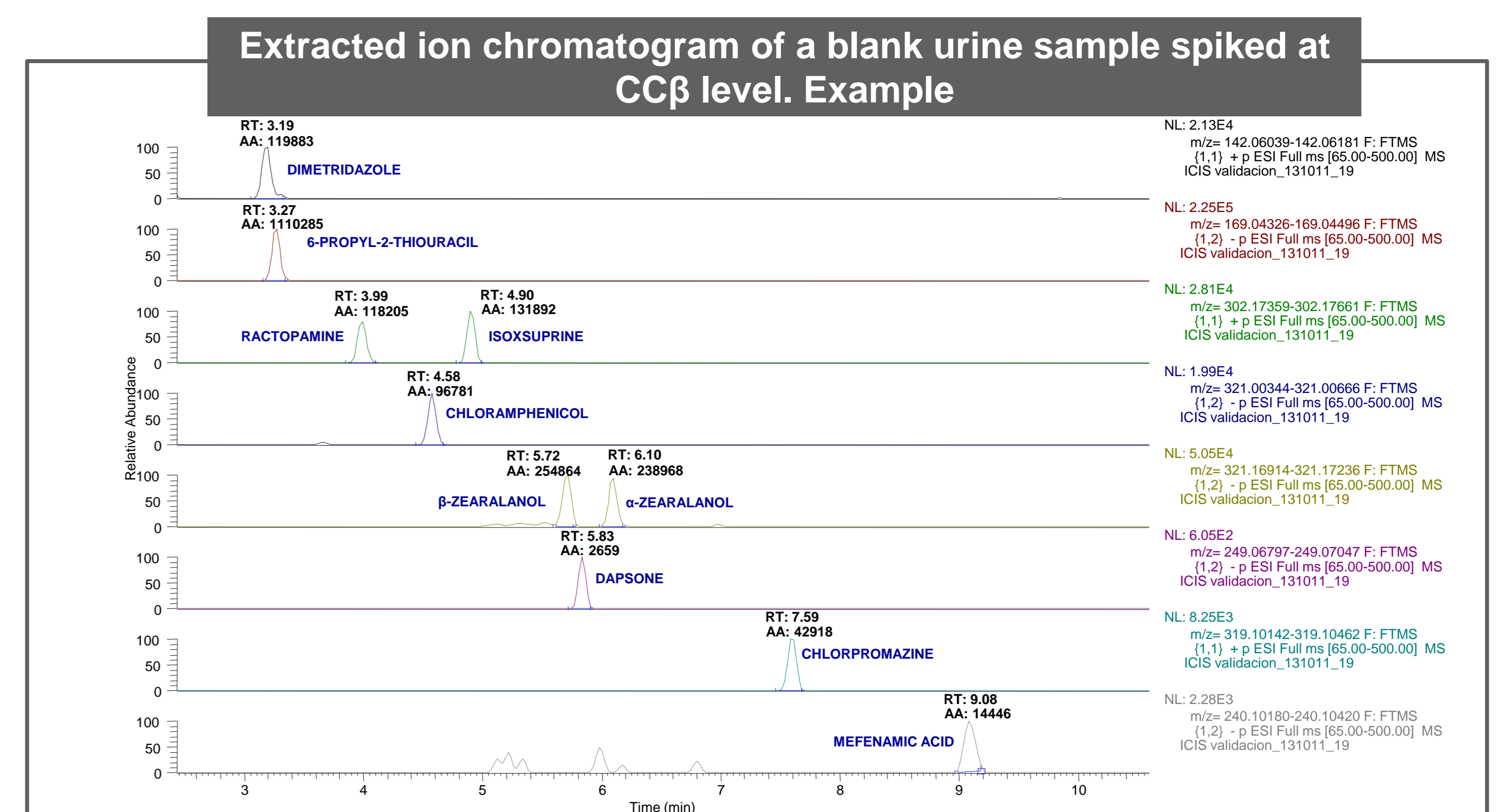
- ✓ Polarity switching: Positive and negative mode
- ✓ Pseudo MS/MS to obtain fragments for confirmation (HCD)

## RESULTS ➔ About 100 compounds

- ✦  **$\beta$ -agonists:** clenbuterol, brombuterol, ractopamine, zilpaterol, isoxuprine, tulobuterol, chorprenaline, cimaterol, ritodrine, terbutaline, metaproterenol, salbutamol, salmeterol, fenoterol, cimbuterol, chrolbrombuterol, mabuterol, mapenterol, hydroxymethylclenbuterol, clenpenterol, clenispenterol, clenicyclohexerol, clenhexerol, formoterol, clenbuterol-d9 and ractopamine-d5
- ✦ **Steroids:**  $\alpha$ -nortestosterone,  $\beta$ -nortestosterone,  $\alpha$ -boldenone,  $\beta$ -boldenone,  $\alpha$ -testosterone, hydroxystanozolol,  $\alpha$ -trenbolone,  $\beta$ -trenbolone, stanozolol, 16 $\beta$ -hydroxystanozolol, 3'-hydroxystanozolol, 4 $\alpha$ -hydroxystanozolol, 4 $\beta$ -hydroxystanozolol, androstenedione, boldione, methyltestosterone,  $\beta$ -nortestosterone-d3,  $\beta$ -boldenone-d3, methyltestosterone-d3, 16 $\beta$ -hydroxystanozolol-d3 and  $\beta$ -trenbolone-d3
- ✦ **RALs:** zearalanone,  $\alpha$ -zearalenol,  $\beta$ -zearalenol,  $\alpha$ -zearalanol,  $\beta$ -zearalanol, zearalenone,  $\alpha$ -zearalanol-d4 and  $\beta$ -zearalanol-d4
- ✦ **Stilbenes:** diethylstilbestrol, dienestrol, hexestrol, diethylstilbestrol-d6, dienestrol-d2 and hexestrol-d4
- ✦ **Nitroimidazoles:** dimetridazole, metronidazole, ronidazole, ipronidazole, HMMNI, ipronidazole-OH, metronidazole-OH, dimetridazole-d3, ronidazole-d3 and ipronidazole-d3
- ✦ **Corticosteroids:** dexamethasone, betamethasone, prednisolone, methylprednisolone, flumethasone and dexamethasone-d4
- ✦ **AINES:** phenylbutazone, oxyphenylbutazone, naproxen, mefenamic acid and phenylbutazone-d10.
- ✦ **Thyrostats:** phenylthiouracil, mercaptobenzimidazole, propylthiouracil, thiouracil, tapazol, methylthiouracil
- ✦ **Sedatives:** chlorpromazine, propionylpromazine, acepromazine, carazolol, azaperone, azaperol, atenolol, haloperidol, xylazine, haloperidol metabolite II and chlorpromazine-d6 and atenolol-d7
- ✦ **Anphenicols:** chloramphenicol, thianphenicol, florphenicol, florphenicol amine, chloramphenicol-d4
- ✦ **Others:** dapsone, monoacetyl dapsone...

## ✓ Database ➔ Characterization of compounds

Elemental composition	Theoretical accurate mass (5 decimals)
Polarity (+/-)	Expected Retention time (min)
Adducts (H <sup>+</sup> , CH <sub>3</sub> COO <sup>-</sup> )	Fragments



## CONCLUSIONS

- ➔ The developed method is simple, fast and wide-range for the screening of veterinary drugs banned or not regulated in bovine urine.
- ➔ The performance characteristics have been determined according to the EU criteria for screening methods (2002/657/EC) and the guideline of the CRLs (2010). The validation was achieved analyzing 20 bovine urine blanks from different animals spiked at CC $\beta$  levels.
- ➔ In all the cases CC $\beta$  levels were equal or lower than action limits or recommended by EU laboratories.
- ➔ From our point of view, the use of high-resolution liquid chromatography combined with high resolution mass spectrometry is a powerful and reliable tool for identification and confirmation in multi-residue analysis. The information obtained from the combination of exact mass at high resolution (R  $\geq$  20000), isotopic pattern and fragments is highly specific, and should be taking into account in future for the establishment of new confirmation criteria in revision of the Decision 2002/657/EC, according to these new technologies.