thermo scientific



Quantification of Catecholamines and Metanephrines in Urine Using the Thermo Scientific[™] TSQ Endura[™] Mass Spectrometer for Research Use

ABSTRACT

Purpose: A rapid analytical method for the quantification of adrenaline, noradrenaline, metanephrine, normetanephrine and dopamine in urine, involving a simple sample preparation procedure, has been developed and analytically validated.

Methods: The method involves a liquid liquid extraction (LLE) of compounds of interest from urine followed by injection onto a Thermo Scientific[™] UltiMate[™] 3000 system; mass spectrometric detection is performed by Single Reaction Monitoring (SRM) on a Thermo Scientific[™] TSQ Endura[™] triple quadrupole mass spectrometer using heated electrospray ionization in positive mode.

Results: The method was analytically validated using charcoal stripped urine spiked with the compounds of interest for lower limit of quantification, linearity range, accuracy, intra- and inter-assay precision and matrix effect evaluation.

INTRODUCTION

Adrenaline (also called epinephrine), noradrenaline (also called norepinephrine) and dopamine are generally called catecholamines. Catecholamines are produced by enzymatic decarboxylation of the amino acid tyrosine in the organism, which is derived from dietary sources as well as synthesis from phenylalanine. They act as neuromodulators in the central nervous system being messengers and play a key role in the blood circulation. Catecholamines are produced mainly by the chromaffin cells of the adrenal medulla and the postganglionic fibers of the sympathetic nervous system. Metanephrine and normetanephrine are produced by the degradation of catecholamines. They are eliminated in urine as free and conjugated metabolites.

Metabolism of catecholamines is a two steps process: they are first methylated by catecholamine-O-methyltransferase (COMT) producing the metanephrines and afterwards the deamination by monoamine oxidase (MAO) produces the homovanilic acid (HVA) and the vanillylmandelic acid (VMA). While COMT is not present in sympathetic cells but is present in adrenal medulla cells, metanephrines are specific markers of catecholamines in chromaffin cells and their tumors.

Liquid chromatography coupled with tandem mass spectrometry is the gold standard technique for the determination and quantification of these neurotransmitters in urine, due to its analytical sensitivity. The use of labeled internal standard increases the accuracy of the quantification. To reduce the matrix effect, methods presented in literature usually involve an extraction step such as SPE (solid phase extraction) which is complex to set-up. A simple liquid-liquid extraction procedure, using a complexing reagent, has been developed in the reported analytical method for research use. Target analytes and corresponding internal standards are reported in Figure 1.

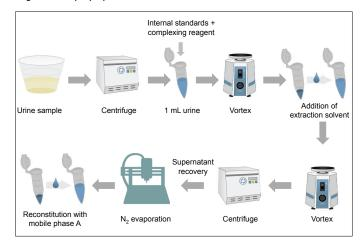
MATERIALS AND METHODS

Sample Preparation

Sample clean-up is performed using a simple LLE procedure with ethyl acetate as the extraction solvent and 2-aminoethyl diphenylboronate as complexing reagent. The diphenyl boronate forms a stable negatively charged complex with cis-hydroxyl groups of catecholamines, that shows a strong affinity for a polar solvent, when operating in alkali media¹.

50 µL of internal standard solution and 1600 µL of complexing reagent solution at the concentration of 2 g/L were added to 1.0 mL of centrifuged urine in a centrifugation tube. pH was adjusted to 9.5 using acetic acid. The tube was vortex-mixed for 60 seconds and 1.5 mL of ethyl acetate was added. After vortex-mixing and centrifugation, 800 µL of supernatant was recovered, evaporated to dryness using a flow of nitrogen and finally reconstituted with 200 µL of mobile phase A. Sample preparation procedure is summarized in Figure 2.

Figure 2. Sample preparation scheme



Thermo Fisher

Liquid chromatography

A 10.0-minute gradient elution was performed using a Thermo Scientific[™] UltiMate[™] 3000RS LC system. Mobile phases consisted of 2 mM ammonium formate in water with 0.1% formic acid and acetonitrile with 0.1% formic acid. The chromatographic separation was achieved using a 2.1 × 150 mm (3µ) Thermo Scientific[™] Acclaim[™] Mixed Mode WCX-1 column at 40°C. Injection volume was10 µL. Further details of the chromatographic method are reported in Table 1.

Table 1. HPLC settings

| Mobile phases | A: 2 mM ammo water + 0.019 | | B: acetonitrile + 0.1% formic acid | | |
|------------------|-------------------------------|-----------------------|------------------------------------|-------|--|
| | Time (min) | Flow Rate (mL/min) | A (%) | В (%) | |
| | 0.0 | 0.40 | 98 | 2 | |
| | 3.5 | 0.40 | 98 | 2 | |
| Gradient Profile | 5.0 | 0.40 | 10 | 90 | |
| | 5.5 | 0.50 | 10 | 90 | |
| | 7.0 | 0.50 | 10 | 90 | |
| | 7.5 | 0.40 | 98 | 2 | |
| | 10.0 | 0.40 | 98 | 2 | |

Mass spectrometry

Target analytes and internal standards were detected by scheduled SRM on a TSQ Endura triple quadrupole mass spectrometer with heated electrospray ionization operating in positive mode. The MS settings are reported in Table 2. A 0.5-minute acquisition time window was used for each analyte and two SRM transitions were included in the acquisition method for quantification and confirmation. Details of the SRM transitions for compounds and internal standards, together with the corresponding source fragmentation, RF lens and collision energy values, are reported in Table 3.

Table 3. SRM settings for target compounds.

| Compound | Precursor Ion (m/z) | Source Fragm. (V) | RF Lens (V) | Product Ion (m/z) | Collision Energy (V) | lon Type |
|--------------------|---------------------------|-------------------------|-------------------|-------------------------|----------------------------|-------------|
| Dopamine | 154.06 | 0 | 142 | 137.05 | 10 | Quan |
| Dopamine | 154.00 | U | 142 | 91.03 | 24 | Confirming |
| Dopamine-D4 | 158.06 | 0 | 80 | 140.99 | 10 | Quan |
| Dopannie-D4 | 156.00 | | 80 | 94.04 | 22 | Confirming |
| Eninophrino | 166.06 | 28.6 | 153 | 107.05 | 20 | Quan |
| Epinephrine | 100.00 | 20.0 | 155 | 135.00 | 15 | Confirming |
| Epinephrine-D6 | 172.06 | 28.6 | 172 | 157.07 | 20 | Quan |
| Ebiliebili lile-Do | | | 172 | 111.05 | 21 | Confirming |
| | 180.06 | 0 | 116 | 165.04 | 17 | Quan |
| Metanephrine | | | | 148.04 | 18 | Confirming |
| Metanephrine-D3 | 183.06 | 0 | 123 | 151.07 | 19 | Quan |
| Metallepilline-D5 | | | | 168.13 | 18 | Confirming |
| Norepinephrine | 152.06 | 4.1 | 93 | 135.00 | 13 | Quan |
| Norepinepinnie | | | | 107.07 | 18 | Confirming |
| Norepinephrine-D6 | 158.06 | 4 | 136 | 111.05 | 18 | Quan |
| Norepinepinine-D6 | | | | 139.07 | 14 | Confirming |
| Normetanephrine | 166.06 | 10.2 | 96 | 121.07 | 17 | Quan |
| | | | 86 | 149.07 | 13 | Confirming |
| Normatananhrina D2 | rine-D3 169.06 | 20.4 | 133 | 137.04 | 17 | Quan |
| Normetanephrine-D3 | | | 133 | 109.11 | 20 | Confirming |

Method evaluation

The analytical method performance was evaluated by defining the following parameters: limit of quantification, linearity range, accuracy, intra- and inter-assay precision and matrix effect for each analyte. Eight calibration levels containing all analytes were prepared by spiking charcoal stripped urine with the proper volumes of standard solutions of the compounds of interest. Each calibrator was extracted and analyzed in 5 replicates to evaluate sensitivity and linearity. A maximum bias percentage between nominal and back-calculated concentration of 15% was set as acceptance criterion for all the calibrators (20% for the lowest). Controls at 4 levels containing all the analytes were also extracted and analyzed in triplicate for three consecutive days, in order to evaluate the precision of the method. A maximum percentage bias between nominal and back-calculated concentration of 20% was set as acceptance criterion for all the control samples. Accuracy for the assay was evaluated in terms of trueness of measurement, measuring the bias percentage between nominal and avarage back-calculated concentration for each calibrator level.

Intra-assay precision was evaluated as the percentage RSD (%RSD) using the controls in replicates of three (n=3) analyzed in one batch.

Inter-assay precision was evaluated on the same controls in replicates of nine (n=9) prepared and analyzed on three different days.

Matrix effect was measured for each analyte as the percentage ratio between analyte / internal standard area ratio in matrix samples and in water at the same concentration for three replicates.

Data Analysis

Data were acquired and processed using Thermo Scientific $^{\rm TM}$ TraceFinder $^{\rm TM}$ 4.1 software.

RESULTS

The assay proved to be linear in the tested calibration range (0.4 - 900 ng/mL) for all the analytes of interest, with a lower limit of quantification (LLOQ) of 0.4 ng/mL and a correlation factor (R²) always above 0.998. The concentrations of prepared calibrators and controls and a calibration performance summary for all the target analytes are reported in Table 4 and 5, respectively. Representative chromatograms for the LLOQ (0.4 ng/mL) for all compounds together with the corresponding calibration curves are reported in Figure 3. The %RSD values for all the controls for each compound for intra-(n=3) and inter-assay (n=9) precision are reported in Table 6 and 7 respectively. Matrix effect values are reported in Table 8.

Table 4. Concentration of calibrators and controls in charcoal stripped urine

| Calibrator ID | Concentration (ng/mL) |
|---------------|-----------------------|
| Cal-0 | 0 |
| Cal-1 | 0.4 |
| Cal-2 | 1.2 |
| Cal-3 | 3.7 |
| Cal-4 | 11.1 |
| Cal-5 | 33.3 |
| Cal-6 | 100 |
| Cal-7 | 300 |
| Cal-8 | 900 |
| Ctrl-1 | 1 |
| Ctrl-50 | 50 |
| Ctrl-200 | 200 |
| Ctrl-800 | 800 |

Table 5. Calibration Performances Summary

| Target Compound | Internal Standard | Curve type | Origin | Weighting | R ² | LOQ (ng/mL) |
|-----------------|---|---------------|--------|-----------|----------------|----------------|
| Metanephrine | Metanephrine- ² H ₃ | Linear | Ignore | 1/X | 0.9991 | 0.4 |
| Normetanephrine | Normetanephrine- ${}^{2}H_{3}$ | Linear | Ignore | 1/X | 0.9992 | 0.4 |
| Dopamine | Dopamine- ² H ₄ | Linear | Ignore | 1/X | 0.9991 | 0.4 |
| Epinephrine | Epinephrine- ² H ₆ | Linear | Ignore | 1/X | 0.9983 | 0.4 |
| Norepinephrine | Norepinephrine- ² H ₆ | Linear | Ignore | 1/X | 0.9979 | 0.4 |

Figure 3. Chromatograms at LLOQ level (0.4 ng/mL) and calibration curves for metanephrine (a), normetanephrine (b), epinephrine (C), norepinephrine (d) and dopamine (e)

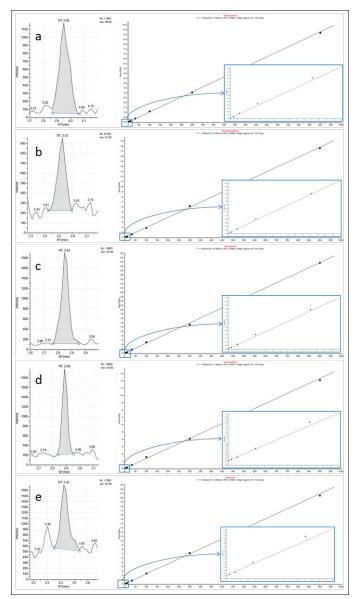


Table 6. %RSD for intra-assay (n=3) precison

| Target Compound | Ctrl-1 | Ctrl-50 | Ctrl-200 | Ctrl-800 |
|-----------------|--------|---------|----------|----------|
| Metanephrine | 8.34 | 3.12 | 4.44 | 2.09 |
| Normetanephrine | 3.67 | 9.89 | 2.33 | 0.45 |
| Epinephrine | 14.90 | 7.51 | 3.69 | 0.66 |
| Norepinephrine | 2.91 | 10.59 | 2.09 | 1.68 |
| Dopamine | 12.37 | 5.60 | 1.03 | 3.90 |

Table 7. %RSD for inter-assay (n=9) precison

| Target Compound | Ctrl-1 | Ctrl-50 | Ctrl-200 | Ctrl-800 |
|-----------------|--------|---------|----------|----------|
| Metanephrine | 7.82 | 6.02 | 3.32 | 2.89 |
| Normetanephrine | 4.82 | 9.71 | 2.80 | 2.45 |
| Epinephrine | 12.91 | 8.75 | 2.91 | 2.17 |
| Norepinephrine | 11.38 | 8.10 | 2.71 | 2.10 |
| Dopamine | 14.55 | 3.96 | 3.11 | 2.98 |

Table 8. Matrix effect

| Target Compound | Cal 3 (%) | Cal 4 (%) | Cal 5 (%) | Cal 6 (%) | Cal 7 (%) |
|-----------------|--------------|--------------|--------------|--------------|--------------|
| Epinephrine | 127.0 | 90.9 | 86.5 | 95.5 | 105.5 |
| Norepinephrine | 115.0 | 109.5 | 95.4 | 93.4 | 112.3 |
| Metanephrine | 130.0 | 85.4 | 86.5 | 88.7 | 102.8 |
| Normetanephrine | 123.6 | 96.1 | 85.7 | 86.3 | 105.4 |
| Dopamine | 187.5 | 203.5 | 86.0 | 83.7 | 87.0 |

CONCLUSIONS

A liquid chromatography tandem mass spectrometry method used for clinical research was developed for the quantification of metanephrine, normetanephrine, epinephrine, norepinephrine and dopamine in urine and analytically validated on a Thermo Scientific UltiMate 3000 system connected to a TSQ Endura triple quadrupole mass spectrometer. This analytical method, based on a simple LLE for sample clean-up, meets research laboratory requirements in terms of sensitivity, linearity of response, accuracy and intra- and inter-assay precision.

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

REFERENCES

1. D. Talwar et al. / J. Chromatogr. B 769 (2002) 341-349

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

Find out more at www.thermofisher.com

©2017 Thermo Fisher Scientific Inc. All rights reserved. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific 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 representatives for details. **PN65046-EN 0817S**

