

Quantification of tricyclic antidepressants in human plasma or serum by liquid chromatography-tandem mass spectrometry for clinical research

Authors

Claudio De Nardi, Thermo Fisher Scientific GmbH, Dreieich, Germany

Sergio Indelicato, Thermo Fisher Scientific, Les Ulis, France

Keywords

Tricyclic antidepressants, offline sample preparation, plasma, serum, mass spectrometry

Goal

Implementation of an analytical method for the quantification of 15 different tricyclic antidepressants in human plasma or serum on a Thermo Scientific™ TSQ Quantis™ triple quadrupole mass spectrometer.

Application benefits

- Simple offline sample preparation by protein precipitation
- 15 tricyclic antidepressants in a single quantitative method

Introduction

An analytical method for clinical research for the quantification of 15 tricyclic antidepressants in human plasma or serum is reported; the analysis includes amitriptyline, clomipramine, clozapine, desipramine, doxepin, imipramine, maprotiline, norclomipramine, norclozapine, nordoxepin, normaprotiline, nortrimipramine, nortriptyline, protriptyline, and trimipramine. Plasma or serum samples were extracted by offline internal standard addition and protein precipitation. Extracted samples were injected onto a Thermo Scientific™ Vanquish™ Flex Binary system connected to a Thermo Scientific™ TSQ Quantis™ triple quadrupole mass spectrometer with heated electrospray ionization. Detection was performed by selected-reaction monitoring (SRM) using 12 deuterated internal standards. Method performance was evaluated using the ClinMass® TDM Platform with the ClinMass Add-On Set for Tricyclic Antidepressants from RECIPE Chemicals + Instruments GmbH (Munich, Germany) in terms of linearity of response within the calibration ranges, accuracy, and intra- and inter-assay precision for each analyte.

Experimental

Target analytes

The analytes and corresponding concentration ranges covered by the calibrators used are reported in Table 1.

Table 1. Concentration ranges covered by calibrators.

Analyte	Concentration (ng/mL)
Amitriptyline	14.9–305
Clomipramine	18.8–393
Clozapine	54.9–1166
Desipramine	16.2–348
Doxepin	13.4–270
Imipramine	16.2–340
Maprotiline	21.2–434
Norclomipramine	20.7–431
Norclozapine	46.0–942
Nordoxepin	12.5–275
Normaprotiline	30.6–677
Nortrimipramine	14.1–315
Nortriptyline	16.4–344
Protriptyline	14.8–313
Trimipramine	16.4–345

Sample preparation

Reagents included four calibrators (including blank) and two controls from RECIPE, as well as 12 deuterated internal standards for quantification. Samples of 50 μ L of plasma or serum were protein precipitated using 100 μ L of precipitating solution containing the internal standards. Precipitated samples were vortex-mixed and centrifuged, and the supernatant was transferred to a clean plate or vial.

Liquid chromatography

Chromatographic separation was achieved using mobile phases and analytical column provided by RECIPE.

Details of the analytical method are reported in Table 2.

Total runtime was 4.0 minutes.

Table 2. Liquid chromatographic method description.

Gradient profile:

Time (min)	Flow Rate (mL/min)	A (%)	B (%)
0.00	0.8	85	15
0.05	0.8	85	15
0.06	0.8	70	30
2.10	0.8	70	30
2.11	0.8	62	38
2.70	0.8	62	38
3.00	0.8	25	75
3.35	0.8	25	75
3.40	0.8	85	15
4.00	0.8	85	15

Injection volume: 5 μ L

Column temp.: 40 $^{\circ}$ C

Mass spectrometry

Analytes and internal standards were detected by SRM on a TSQ Quantis triple quadrupole mass spectrometer with heated electrospray ionization operated in positive mode. A summary of the MS conditions is reported in Table 3. Two SRM transitions for each analyte were included in the acquisition method for quantification and confirmation, respectively.

Table 3. MS settings.

Source type:	Heated electrospray ionization (HESI)
Vaporizer temperature:	460 $^{\circ}$ C
Capillary temperature:	368 $^{\circ}$ C
Spray voltage (positive mode):	3500 V
Sheath gas:	50 AU
Sweep gas:	0 AU
Auxiliary gas:	15 AU
Data acquisition mode:	Selected-reaction monitoring (SRM)
Collision gas pressure:	2.0 mTorr
Cycle time:	0.300 s
Q1 mass resolution (FWMH):	0.7
Q3 mass resolution (FWMH):	0.7

Method evaluation

The method performance was evaluated in terms of linearity of response within the calibration ranges, accuracy, and intra- and inter-assay precision for each analyte. Analytical accuracy was evaluated in terms of percentage bias between nominal and average back-calculated concentrations using quality control samples at two different levels provided by RECIPE (MS9182 batch #1456), prepared and analyzed in replicates of five on three different days. Intra-assay precision was evaluated for each day on the same set of runs (control samples at two levels, replicates of five each day, three days) in terms of percentage coefficient of variation (%CV). Inter-assay precision was evaluated on the same controls including all the 15 replicates of the three days.

Data analysis

Data were acquired and processed using Thermo Scientific™ TraceFinder™ 4.1 software.

Results and discussion

The method proved to be linear in the calibration ranges covered by the calibrators. Representative chromatograms for the lowest calibrator for nordoxepin, trimipramine, and their internal standards are reported in Figure 1. Representative calibration curves for the same analytes are reported in Figure 2.

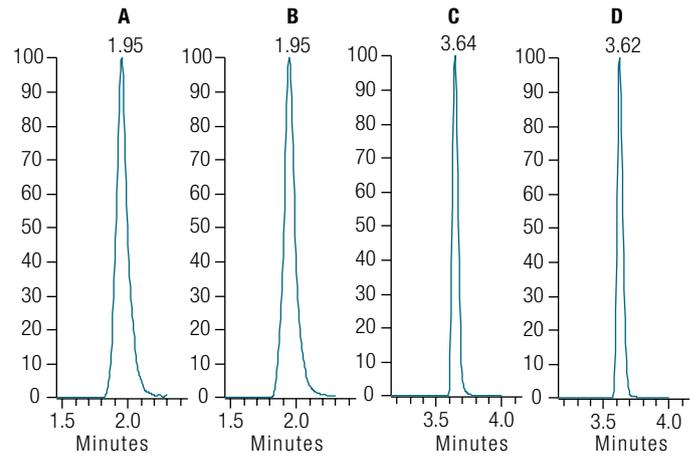


Figure 1. Representative chromatograms of the lowest calibrator for (A) nordoxepin, (B) d3-nordoxepin, (C) trimipramine, and (D) d3-trimipramine.

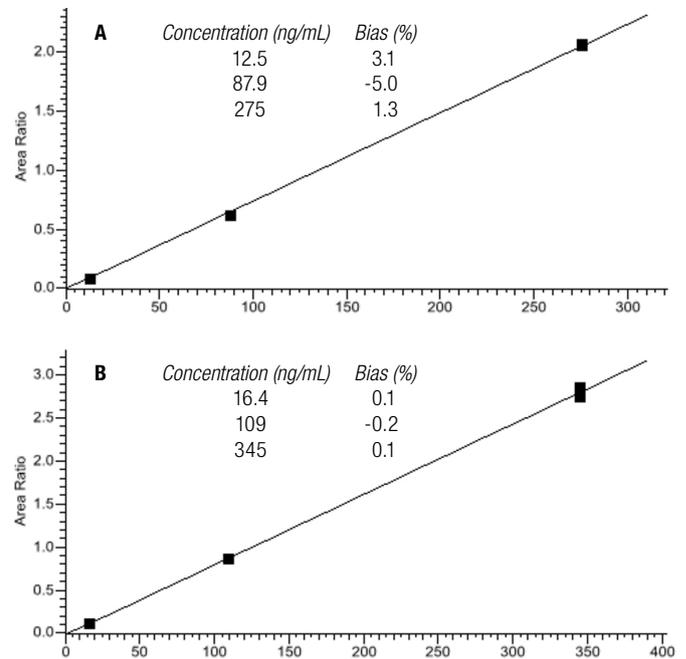


Figure 2. Representative calibration curves for (A) nordoxepin and (B) trimipramine – day 1.

The data demonstrated outstanding accuracy of the method with the percentage bias between nominal and average back-calculated concentration for the control samples ranging from -7.3% to 4.6%. Results are reported in Table 4.

The %CV for intra-assay precision was always below 7.7% for all the analytes (Table 5). The maximum %CV for inter-assay precision including all the analytes was 7.6% (Table 6).

Table 4. Analytical accuracy results for control MS9182 batch #1456.

Analyte	Control 1			Control 2		
	Nominal Concentration (ng/mL)	Average Calculated Concentration (ng/mL)	Bias (%)	Nominal Concentration (ng/mL)	Average Calculated Concentration (ng/mL)	Bias (%)
Amitriptyline	59.2	55.8	-5.8	135	125	-7.3
Clomipramine	73.9	71.1	-3.8	171	162	-5.5
Clozapine	217	209	-3.8	510	483	-5.3
Desipramine	64.3	62.3	-3.1	152	146	-3.8
Doxepin	50.8	49.5	-2.6	117	114	-3.0
Imipramine	63.9	65.4	2.4	148	145	-2.1
Maprotiline	82.7	79.1	-4.3	193	181	-6.1
Norclomipramine	79.9	78.1	-2.2	187	176	-5.7
Norclozapine	179	171.6	-4.1	418	395	-5.6
Nordoxepin	49.4	50.2	1.7	116	113	-2.7
Normaprotiline	121	125.6	3.8	280	293	4.6
Nortrimipramine	55.1	55.6	1.0	133	129	-2.9
Nortriptyline	64.5	64.0	-0.8	145	144	-0.9
Protriptyline	58.1	58.5	0.7	143	135	-5.5
Trimipramine	64.2	62.9	-2.0	155	146	-5.5

Table 5. Intra-assay precision results for control MS9182 batch #1456.

Analyte	Control 1						Control 2					
	Day 1		Day 2		Day 3		Day 1		Day 2		Day 3	
	Average Calculated Concentration (ng/mL)	CV (%)										
Amitriptyline	53.7	1.2	55.6	1.4	57.6	1.1	121	4.6	121	6.2	132	8.4
Clomipramine	69.5	3.1	70.5	3.0	73.1	2.2	159	4.9	158	7.9	168	7.4
Clozapine	202	3.5	206	3.6	216	2.6	465	3.5	488	2.3	491	6.5
Desipramine	61.7	2.6	63.3	4.6	61.8	2.4	143	2.7	149	4.1	146	6.7
Doxepin	47.9	1.3	49.5	2.8	50.6	2.9	108	4.0	114	3.8	118	6.2
Imipramine	67.1	2.0	62.5	4.4	67.0	6.2	140	5.1	147	7.8	147	7.8
Maprotiline	78.0	3.6	79.9	2.6	79.2	1.7	180	3.6	181	8.1	182	7.4
Norclomipramine	75.7	2.4	77.1	2.9	81.0	2.4	171	3.6	170	6.6	187	6.2
Norclozapine	170	1.6	172	3.0	173	1.6	400	6.2	385	5.8	400	6.3
Nordoxepin	49.4	2.9	49.2	2.8	51.9	2.5	111	3.6	111	5.2	117	6.0
Normaprotiline	130	2.8	120	5.0	128	2.1	310	2.3	279	3.5	293	3.3
Nortrimipramine	55.8	7.7	56.0	6.3	55.2	4.0	135	5.7	127	5.0	127	6.2
Nortriptyline	62.1	3.0	63.0	6.0	66.5	4.1	137	2.4	144	4.4	149	5.6
Protriptyline	57.1	6.7	58.1	4.1	60.0	1.6	135	2.2	130	6.0	140	6.7
Trimipramine	61.5	4.7	61.2	3.5	65.8	3.8	143	3.7	143	6.7	153	8.6

Table 6. Inter-assay precision results for control MS9182 batch #1456.

Analyte	Control 1		Control 2	
	Average Calculated Concentration (ng/mL)	CV (%)	Average Calculated Concentration (ng/mL)	CV (%)
Amitriptyline	55.8	3.1	125	7.6
Clomipramine	71.1	3.4	162	7.1
Clozapine	209	4.2	483	4.9
Desipramine	62.3	3.4	146	4.8
Doxepin	49.5	3.3	114	5.9
Imipramine	65.4	5.5	145	7.1
Maprotiline	79.1	2.6	181	6.4
Norclomipramine	78.1	3.8	176	7.1
Norclozapine	172	2.1	395	5.9
Nordoxepin	50.2	3.6	113	5.4
Normaprotiline	126	4.9	293	5.2
Nortrimipramine	55.6	5.6	129	6.1
Nortriptyline	64.0	5.3	144	5.6
Protriptyline	58.5	4.5	135	6.1
Trimipramine	62.9	5.1	146	7.2

Conclusions

A liquid chromatography-tandem mass spectrometry method for clinical research for the quantification of 15 different tricyclic antidepressants in human plasma or serum was implemented. The ClinMass TDM Platform with the ClinMass Add-On Set for Tricyclic Antidepressants from RECIPE was used. The method

was analytically evaluated on a Vanquish Flex Binary system connected to a TSQ Quantis triple quadrupole mass spectrometer. The method offers the quick and simple offline protein precipitation with concomitant internal standard addition. The described method meets research laboratory requirements in terms of sensitivity, linearity of response, accuracy, and precision.

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

Find out more at thermofisher.com/ClinicalResearchSolutions

ThermoFisher
SCIENTIFIC

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries unless otherwise specified. ClinMass is a registered trademarks of RECIPE Chemicals + Instruments GmbH. 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.

TN65128-EN 0118S