TECHNICAL NOTE 66052

Quantification of 15 tricyclic antidepressants in human plasma by liquid chromatographytandem mass spectrometry for clinical research

Authors: Gaëtan Renoulin,¹ Mariana Barcenas,¹ and Claudio De Nardi²

¹Thermo Fisher Scientific, Les Ulis, France ²Thermo Fisher Scientific, Reinach, Switzerland Keywords: TraceFinder software, TSQ Quantis MS, Vanquish Flex UHPLC, LC-MS, mass spectrometry, offline sample preparation, plasma, tricyclic antidepressants

Application benefits

- Robust, sensitive hardware enables increased confidence in data
- Simple offline sample preparation by protein precipitation
- Fifteen tricyclic antidepressants in a single quantitative method

Goal

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

Introduction

Tricyclic antidepressants (TCAs) are a group of psychoactive drugs that are mainly used to alleviate symptoms of anxiety, endogenous depression, and pain.



In this report, an analytical method for clinical research for the quantification of 15 tricyclic antidepressants in human plasma in eight minutes is presented. The samples were prepared offline by protein precipitation. Extracted samples were chromatographically separated on a Thermo Scientific™ Vanquish™ Flex Binary UHPLC system. Detection was performed on a TSQ Quantis triple quadrupole mass spectrometer with heated electrospray ionization (HESI) operated in positive ionization mode. Method performance was evaluated using the ClinMass® Add-on Set for TCAs, including LC-MS/MS calibrators, controls, and internal standards from RECIPE Chemicals + Instruments GmbH (Munich, Germany) in terms of linearity of response, lower limit of quantification (LLOQ), accuracy, and intra- and interassay precision for all analytes.



Table 1. List of analytes, ranges of concentrations (MS9113 Batch #1389), and internal standards

Compound name	Formula	Concentration Range (μg/L)	Internal standard name	Formula
Amitriptyline	C ₂₀ H ₂₃ N	16.3–331	Amitriptyline-d ₃	C ₂₀ H ₂₀ D ₃ N
Clomipramine	$C_{19}H_{23}CIN_2$	18.3–361	Clomipramine-d ₃	$C_{19}H_{20}D_3CIN_2$
Clozapine	$C_{18}H_{19}CIN_4$	58.5-1202	Clozapine-d ₄	$C_{18}H15D_4CIN_4$
Desipramine	C ₁₈ H ₂₂ N ₂	18.1–377	Desipramine-d ₃	C ₁₈ H ₁₉ D3N ₂
Doxepin	$C_{19}H_{21}NO$	14.8–287	Doxepin-d ₃	$C_{19}H_{18}D_{3}NO$
Imipramine	C ₁₉ H ₂₄ N ₂	16.7–343	Imipramine-d₃	$C_{19}H_{21}D_3N_2$
Maprotiline	$C_{20}H_{23}N$	24.2-465	Maprotiline-d ₅	$C_{20}H_{20}D_3N$
Norclomipramine	$C_{18}H_{21}CIN_2$	21.1-416	Norclomipramine-d ₃	$C_{18}H_{18}D_3CIN_2$
Norclozapine	$C_{17}H_{17}CIN_4$	47.3-950	Norclozapine-d ₈	$C_{17}H_9D_8CIN_4$
Nordoxepin	C ₁₈ H ₁₉ NO	13.6–279	Nordoxepin-d ₃	$C_{18}H_{16}D_3NO$
Normaprotiline	$C_{19}H_{21}N$	33.4–725	Desipramine-d ₃	$C_{18}H_{19}D_{3}N_{2}$
Nortrimipramine	$C_{19}H_{24}N_2$	11.0–220	Imipramine-d ₃	$C_{19}H_{21}D_{3}N_{2}$
Nortriptyline	$C_{19}H_{21}N$	18.2–371	Nortriptyline-d ₃	$C_{19}H_{18}D_{3}N$
Protriptyline	C ₁₉ H ₂₁ N	16.9–344	Nortriptyline-d ₃	$C_{19}H_{18}D_{3}N$
Trimipramine	$C_{20}H_{26}N_2$	27.3-579	Trimipramine-d ₃	$C_{20}H_{23}D_3N_2$

Experimental

Target analytes

The complete list of analytes with their corresponding internal standards and the concentration ranges covered by the calibrators used are reported in Table 1.

Sample preparation

Reagents included four calibrators (including blank) and two controls from RECIPE, as well as an internal standard mix for quantitation. Samples of 50 μ L of plasma were protein precipitated using 100 μ L of acetonitrile containing the internal standards. Precipitated samples were vortex-mixed and centrifuged for 10 minutes. The supernatant was transferred to a clean vial.

Liquid chromatography

The supernatant was injected onto a Vanquish Flex Binary UHPLC system connected to a TSQ Quantis triple quadrupole mass spectrometer. Chromatographic separation was achieved by gradient elution on a Thermo Scientific™ Hypersil GOLD™ Phenyl (100 × 2.1 mm, 1,9 µm) column (P/N 25902-102130) kept at 40 °C.

Table 2. LC gradient profile

Time	Flow (mL/min)	%В
0.00	0.500	10.00
0.20	0.500	10.00
0.50	0.500	50.00
4.50	0.500	50.00
5.20	0.500	100.0
6.50	0.500	100.0
6.51	0.500	10.00
8.00	0.500	10.00

Mobile phases composition is the following:

- Mobile phase A: Water + 0.1% formic acid + 5 mM ammonium formate
- Mobile phase B: Methanol + 0.1% formic acid + 5 mM ammonium formate

Injection volume was 2 µL.

The LC gradient is described in Table 2. Total runtime was 8.0 minutes.

Mass spectrometry

Analytes and internal standards were detected by selected reaction monitoring (SRM) on a TSQ Quantis triple quadrupole mass spectrometer using a HESI source operated in positive ionization mode. A summary of the MS conditions is reported in Table 3. One confirming ion for each analyte was included in the acquisition method for confirmation (Table 4).

Method evaluation

The performance of the method was evaluated in terms of carryover, LLOQ, accuracy, and intra- and inter-assay precision for all analytes.

Carryover was calculated in terms of percentage ratio between peak area of the highest calibrator and a blank sample injected just after it.

Table 4. Description of the SRM parameters

Table 3. MS parameters

lon source parameters							
Heated Electrospray Source Ionization (HESI)							
3,500							
50							
10							
0							
300							
320							
Settings							
SRM							
SRM parameters							
0.3							
0.7							
1.2							
6							

Analyte/Internal standard	Retention	Quantification			Confirmation			RF lens
	(min)	Precursor ion (m/z)	Product ion (<i>m/z</i>)	Collision energy (V)	Precursor ion (m/z)	Product ion (<i>m/z</i>)	Collision energy (V)	(V)
Amitriptyline	5.6	278.2	105.1	23	278.2	233.2	17	122
Clomipramine	5.9	315.2	86.2	18	315.2	58.13	29	119
Clozapine	3.3	327.1	192.1	44	327.1	270.1	22	148
d ₃ -Amitriptyline	5.6	281.2	281.2	23	281.2	_	_	122
d ₃ -Clomipramine	5.9	318.2	318.2	18	318.2	_	_	119
d ₃ -Desipramine	5.1	270.2	270.2	16	270.2	_	_	101
d ₃ -Doxepin	3.4	283.2	283.2	23	283.2	_	_	123
d ₃ -Imipramine	5.1	284.2	284.2	28	284.2	_	_	111
d ₃ -Norclomipramine	5.9	304.2	304.2	17	304.2	_	_	112
d ₃ -Nordoxepin	3.4	269.2	269.2	22	269.2	_	_	117
d ₃ -Nortriptyline	5.6	267.2	267.2	14	267.2	_	_	111
d ₃ -Trimipramine	5.7	298.2	298.2	17	298.2	_	_	111
d ₄ -Clozapine	3.3	331.2	331.2	44	331.2	_	_	148
d ₅ -Maprotiline	5.4	283.2	283.2	19	283.2	_	_	128
d ₈ -Norclozapine	3.2	321.2	321.2	41	321.2	_	_	157
Desipramine	5.1	267.2	72.13	16	267.2	44.07	42	101
Doxepin	3.4	280.2	107	23	280.2	58.13	21	123
Imipramine	5.1	281.2	58.13	30	281.2	86.2	17	107
Maprotiline	5.4	278.2	250.2	25	278.2	178.1	39	128
Norclomipramine	5.9	301.2	72.13	17	301.2	44.07	49	112
Norclozapine	3.2	313.1	192.13	41	313.1	270.1	24	157
Nordoxepin	3.4	266.2	107.1	22	266.2	44.07	18	117
Normaprotiline	5.3	264.2	169.2	18	264.2	219.1	22	113
Nortrimipramine	5.7	281.2	44.07	28	281.2	55.05	30	108
Nortriptyline	5.6	264.2	233.2	14	264.2	105.1	20	111
Protriptyline	5.0	264.2	233.2	16	264.2	155.1	21	125
Trimipramine	5.7	295.2	100.1	17	295.2	58.13	28	111

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 #1389), prepared and analyzed in replicates of five on three different days. Intra-assay precision for each day was evaluated in terms of percentage coefficient of variation (%CV) using the controls at two different levels in replicates of five (n=5). Inter-assay precision was evaluated as the %CV on the full set of samples (control samples at two levels in replicates of five prepared and analyzed on three different days).

LLOQ was investigated by dilution of the lowest calibrator using blank matrix and was established as the lowest concentration with a mean accuracy and precision below than 20%.

Data analysis

Data were acquired and processed using Thermo Scientific™ TraceFinder™ 5.1 software.

Results and discussion

A linear interpolation with 1/x weighting was used for all analytes. The percentage bias between nominal and back-calculated concentration was always within ±10% for all the calibrators in all the runs. Representative chromatograms at the LLOQ for clozapine, norclomipramine, nortriptyline, trimipramine, and their corresponding internal standards are reported in Figure 1. Representative calibration curves for the same analytes are reported in Figure 2.

No carryover was observed for any of the analytes, with no signal detected in the blank injected immediately after the highest calibrator.

The data demonstrated good accuracy of the method with the percentage bias between nominal and average back-calculated concentration for the control samples ranging between -8.6% and 9.3% (Table 5). The %CV for intra-assay precision was always below 6.2% for all the analytes. The maximum %CV for inter-assay precision including all the analytes was 6.2%. Results for intra- and inter-assay precision are reported in Table 6.

LLOQs of all compounds were determined and reported in Table 7.

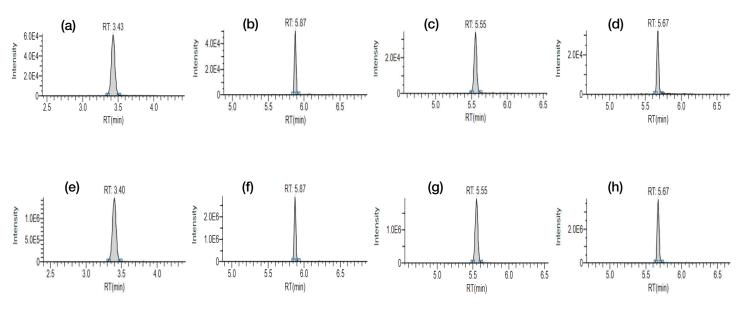


Figure 1. Representative chromatograms of the lower limits of quantification for (a) clozapine, (b) norclomipramine, (c) nortriptyline, (d) trimipramine, (e) d_3 -norclomipramine, (g) d_3 -nortriptyline, (h) d_3 -trimipramine

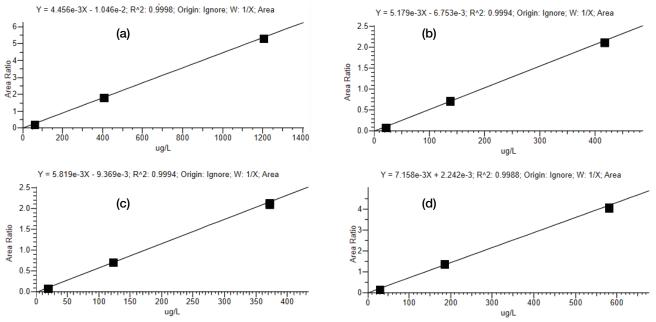


Figure 2. Representative calibration curves for (a) clozapine, (b) norclomipramine, (c) nortriptyline, (d) trimipramine

Table 5. Analytical accuracy results for control MS9182 Batch #1389

Analyte	Control	Nominal concentration (µg/L)	Average calculated concentration (µg/L)	Bias (%)
Amitriatulina	Level I	59.1	60.9	3.1
Amitriptyline	Level II	137	149	8.8
Clomipramine	Levell	67.3	69.3	3.0
Olompramme	Level II	161	165	2.7
Clozapine	Level I	229	228	-0.6
Olozapii le	Level II	529	540	2.1
Desipramine	Levell	68.1	66.9	-1.7
Desipramine	Level II	158	164	3.6
Dovonin	Levell	53.3	52.3	-1.9
Doxepin	Level II	123	128	4.3
landa ya ya in a	Levell	62.2	61.0	-1.9
Imipramine	Level II	143	152	6.1
Managatilia	Levell	86.6	83.0	-4.2
Maprotiline	Level II	199	200	0.4
Nevelemieromine	Levell	78.7	77.3	-1.7
Norclomipramine	Level II	180	183	1.9
Navalanasiaa	Levell	183	180	-1.5
Norclozapine	Level II	413	422	2.2
Nordovania	Levell	50.6	49.1	-3.0
Nordoxepin	Level II	118	121	2.5
N I a was a sawa tili a a	Level I	129	118	-8.6
Normaprotiline	Level II	297	295	-0.8
Nantainalana	Level I	40.4	39.9	-1.2
Nortrimipramine	Level II	90.8	98.2	8.1
N I a solution to allow a	Level I	69.1	69.0	-0.2
Nortriptyline	Level II	160	166	3.7
Duntalists die e	Level I	61.8	60.2	-2.5
Protriptyline	Level II	146	147	0.9
Tolorionomolor	Levell	102	108	6.2
Trimipramine	Level II	240	262	9.3
		5		

Table 6. Analytical intra- and inter-assay precision results for control MS9182 Batch #1389

	Intra-assay						Inter-assay		
	Control	Day 1		Day 2		Day 3			
Analyte		Average calculated concentration (µg/L)	CV (%)	Average calculated concentration (µg/L)	CV (%)	Average calculated concentration (µg/L)	CV (%)	Average calculated concentration (µg/L)	CV (%)
Amitriptyline	Level I	57.8	2.5	62.6	2.7	62.5	4.0	60.9	4.5
Aminiptyline	Level II	141	4.8	153	5.5	153	3.6	149	4.5
Clomipramine	Level I	66.2	2.2	70.7	3.5	71.0	2.9	69.3	3.8
Olompiamine	Level II	157	4.7	171	4.4	168	4.3	165	4.6
Clozapine	Level I	220	2.3	231	3.5	232	2.9	227	2.8
σιοΣαριπο	Level II	523	4.4	554	4.0	543	4.6	540	2.9
Desipramine	Level I	64.7	1.9	68.9	4.0	67.1	2.7	66.9	3.1
Desipramilie	Level II	158	5.3	171	4.8	163	3.6	164	4.0
Doxepin	Level I	50.5	3.1	52.7	3.1	53.7	2.6	52.3	3.1
Волорін	Level II	123	4.7	131	5.2	130	4.4	128	3.2
Imipramine	Level I	58.6	2.8	62.3	4.6	62.1	2.6	61.0	3.5
IIIIpramine	Level II	145	5.1	156	4.6	154	4.8	152	3.9
Maprotiline	Level I	80.4	2.5	85.7	3.5	82.8	2.6	83.0	3.2
Maprotiline	Level II	193	5.1	210	4.5	196	3.8	200	4.5
Norclomipramine	Level I	75.5	2.2	78.8	2.6	77.8	2.6	77.3	2.2
Nordiomplamile	Level II	179	4.3	189	4.7	182	3.2	183	2.7
Norclozapine	Level I	179	1.4	182	3.9	180	3.0	180	0.8
ΝοισιοΖαρίπο	Level II	419	3.9	431	3.4	417	3.6	422	1.7
Nordoxepin	Level I	47.4	3.5	50.1	2.4	49.7	2.0	49.1	3.0
Погаохерігі	Level II	118	5.2	125	4.5	120	4.3	121	3.3
Normaprotiline	Level I	113	5.4	122	4.7	119	3.4	118	3.9
Normaprodiline	Level II	284	4.6	316	5.2	285	5.0	295	6.2
Nortrimipramine	Level I	37.9	3.4	40.0	4.8	41.8	1.7	39.9	4.9
Northinipranine	Level II	93.7	6.2	100	5.7	101	4.1	98.2	4.0
Nortriptyling	Level I	66.7	3.1	69.9	4.0	70.3	3.2	69.0	2.9
Nortriptyline	Level II	159	5.1	172	5.4	166	4.1	166	4.0
Protriptyline	Level I	58.2	3.2	61.8	3.9	60.8	3.7	60.3	3.1
Promptyline	Level II	143	5.3	154	5.6	145	5.7	147	4.1
Triminramina	Level I	103	1.7	109	3.8	113	3.7	108	4.8
Trimipramine	Level II	249	5.3	267	4.8	270	4.3	262	4.4

thermoscientific

Table 7. LLOQs for all compounds

Analyte	LLOQ (μg/L)				
Amitriptyline	3.26				
Clomipramine	3.66				
Clozapine	11.7				
Desipramine	3.62				
Doxepin	2.96				
Imipramine	3.34				
Maprotiline	4.84				
Norclomipramine	4.22				
Norclozapine	2.37				
Nordoxepin	2.72				
Normaprotiline	18.2				
Nortrimipramine	16.9				
Nortriptyline	3.64				
Protriptyline	3.38				
Trimipramine	1.37				

Conclusion

A robust, reproducible, and sensitive liquid chromatography-tandem mass spectrometry method for clinical research for quantification of fifteen tricyclic antidepressants in human plasma was developed and implemented. The analytical method was validated on an Vanquish Flex Binary UHPLC system coupled to a TSQ Quantis triple quadrupole mass spectrometer. Method performance was evaluated using the ClinMass LC-MS/MS calibrators, controls, and internal standards. The method described here offers 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.

Find out more at thermofisher.com/ClinicalResearchApps



