Quantification of 28 neuroleptics in human plasma by LC-HRAM-MS for clinical research

Author: Mariana Barcenas Thermo Fisher Scientific, Les Ulis, France

Keywords: Orbitrap Exploris 120 MS, Vanquish Flex UHPLC, TraceFinder software, neuroleptics, antipsychotic drugs, offline sample preparation, plasma, high-resolution mass spectrometry

Application benefits

- Accurate results, simple sample preparation, and rapid quantitation
- High-resolution mass spectrometry for improved selectivity
- Robust, sensitive LC and MS platforms enable increased confidence in data

Goal

Implementation of an analytical method for the quantification of 28 neuroleptics in human plasma on a Thermo Scientific[™] Orbitrap Exploris[™] 120 mass spectrometer connected to a Thermo Scientific[™] Vanquish[™] Binary Flex UHPLC system.

Introduction

Neuroleptics are a class of psychoactive drugs that are used to treat psychoses, a symptom of schizophrenia. Since antipsychotic drugs are very powerful and can cause



adverse side effects, they are usually administered at low daily dosages. While most antipsychotic drugs are found in plasma in the low ng/mL range, some are found in very low concentrations in brain tissue. High performance liquid chromatography (HPLC) with UV detection is widely used for measurement of these drugs. However, identification and quantitation of antipsychotics at low concentrations require technology that can offer high resolution (to separately identify every analyte with confidence) and high sensitivity. The use of high-resolution accurate-mass (HRAM) mass spectrometers coupled to UHPLC systems can offer the necessary selectivity and specificity while providing the required sensitivity.



In this study, a clinical research analytical method for the quantification of 28 neuroleptics in human plasma is reported. Plasma samples were extracted by internal standard addition and protein precipitation. Method performance was evaluated using calibrators and controls from RECIPE Chemicals + Instruments GmbH (Munich, Germany) in terms of linearity of response within the calibration ranges, lower limit of quantification (LLOQ), carryover, accuracy, and intra- and inter-assay precision for the entire panel of analytes.

Experimental

Target analytes

The list of analytes, their corresponding internal standards, and concentration ranges covered by the calibrators (MS9313 batch #1368) are reported in Table 1.

Table 1. Analytes, internal standards, and concentration ranges covered by calibrators

Analyte			Concentration (ng/mL)			
	Internal standard	Retention time (min)	ы	L2	L3	
Amisulpride	Amisulpride-d ₅	2.46	37.1	255	751	
Aripiprazole	Aripiprazole-d ₈	3.66	54.8	362	1129	
Chlorpromazine	Chlorpromazine-d ₆	4.19	18.5	126	375	
Chlorprothixene	Chlorprothixene-d ₆	4.22	17.7	125	378	
Clozapine	Clozapine-d ₄	3.26	57.1	406	1223	
Dehydro-Aripiprazole	Dehydro-Aripiprazole-d ₈	3.59	10.5	70.9	217	
Desmethylolanzapine	Olanzapine-d ₃	2.29	6.75	49.3	146	
Flupentixol	Flupenthixol-d ₄	4.33	0.648	4.56	13.2	
Fluphenazine	Fluphenazine-d ₈	4.31	0.590	4.12	12.7	
Haloperidol	Haloperidol-d ₄	3.40	0.630	4.28	12.5	
Levomepromazine	Levomepromazine-d ₃	3.84	8.54	60.2	182	
Melperone	Melperone-d ₄	2.95	11.0	72.4	220	
Norclozapine	Norclozapine-d ₈	3.19	45.3	323	961	
Norquetiapine	Quetiapine-d ₈	3.38	8.71	61.3	183	
Olanzapine	Olanzapine-d ₃	2.31	7.14	50.2	145	
Paliperidone	Paliperidone-d ₄	2.96	6.70	45.3	135	
Perazine	Perazine-d ₈	3.87	23.9	160	487	
Pipamperone	Pipamperone-d ₁₀	2.59	30.9	218	636	
Promethazine	Promethazine-d ₆	3.58	20.8	139	416	
Prothipendyl	Prothipendyl-d ₆	3.43	1.08	6.93	22.0	
Quetiapine	Quetiapine-d ₈	3.39	30.0	200	571	
Risperidone	Risperidone-d ₄	3.01	6.51	44.0	139	
Sertindole	Sertindole-d ₄	4.27	10.3	71.0	222	
Sulpiride	Sulpiride-d ₃	1.90	54.7	382	1154	
Thioridazine	Thioridazine-d ₃	4.28	19.1	142	428	
Ziprasidone	Ziprasidone-d ₈	3.21	16.8	116	335	
Zotepine	Sertindole-d ₄	4.25	9.21	59.1	189	
Zuclopenthixol	Zuclopenthixol-d ₄	4.30	3.85	26.4	81.8	

Sample preparation

Reagents included four calibrators (including blank) and two controls (MS9382 batch #1279) from RECIPE, as well as internal standards for quantification. Samples of 50 μ L of plasma were protein precipitated using 100 μ L of acetonitrile containing the internal standard. Precipitated samples were vortex-mixed and centrifuged. The supernatant was diluted 10-fold using mobile phase A and transferred to a clean vial.

Liquid chromatography

LC separation was performed on a Thermo Scientific[™] Vanquish[™] Flex Binary UHPLC system using the following mobile phases:

- Mobile phase A: 0.1% formic acid in water
- Mobile phase B: 0.1% formic acid in methanol

Chromatographic separation was achieved by gradient elution on a Thermo Scientific[™] Hypersil GOLD[™] 2.1 × 50 mm (1.9 µm) analytical column (P/N 25002-052130) maintained at 40 °C at a flow rate of 0.4 mL/min. Total run time was 5.5 minutes. The chromatographic conditions are given in Table 2.

Table 2. LC gradient profile

Time (min)	Flow (mL/min)	В (%)			
0.00	0.4	2			
0.80	0.4	2			
2.50	0.4	55			
3.40	0.4	55			
3.50	0.4	98			
4.50	0.4	98			
4.51	0.4	2			
5.50	0.4	2			
(Other parameters				
Column temperatu	40 °C				
Injection volume	Injection volume				

Mass spectrometry

Detection was performed on an Orbitrap Exploris 120 mass spectrometer equipped with a heated electrospray ionization (HESI) ion source operated in positive ionization mode. Data was acquired in Full MS-ddMS² mode using a resolution of 60,000 (FWHM) at m/z 200 on a scan range of m/z 100 to 500. The ion source conditions and the mass spectrometer settings are presented in Tables 3 and 4, respectively.

Table 3. Ion source settings

Parameter	Value
Source type	Heated Electrospray Ionization (HESI)
Vaporizer temperature	350 °C
lon transfer tube	300 °C
Capillary voltage	3,500 V
Sheath gas	40 AU
Sweep gas	0 AU
Auxiliary gas	10 AU

Table 4. MS settings

Parameter	Value
Resolution at <i>m/z</i> 200	60,000
Scan range	100–500
AGC target	Standard
RF lens	70%
Maximum injection time mode	Auto
Data type	Profile
Polarity	Positive
Source fragmentation	Off
Mild trapping	Off

Method evaluation

The parameters used to evaluate the performance of the method included linearity of response, LLOQ, intraand inter-assay accuracy and precision, and carryover for all the analytes. Carryover was calculated in terms of percentage ratio between peak area of the highest calibrator and a blank sample injected immediately after it. Analytical accuracy was evaluated in terms of percentage bias between nominal and average backcalculated concentrations using quality control samples at two different levels provided by RECIPE, 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). The LLOQ was investigated by dilution of the lowest calibrator with blank matrix and was established as the lowest concentration with a mean accuracy and precision better than 20%.

Data analysis

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

Results and discussion

A summary of the LLOQs obtained for all the analytes is presented in Table 5. A linear regression with 1/× weighting down to the LLOQ value was obtained for all the analytes, except for norclozapine where a quadratic regression was used. The percentage bias between nominal and backcalculated concentration was always within ±10% for all the calibrators in all the runs.

Table 5. Lower limit of quantitation

Assel	Concentration (ng/mL)				
Analyte	Lowest calibrator	LLOQ			
Amisulpride	37.1	3.71			
Aripiprazole	54.8	5.48			
Chlorpromazine	18.5	3.70			
Chlorprothixene	17.7	3.54			
Clozapine	57.1	5.71			
Dehydro-Aripiprazole	10.5	5.25			
Desmethylolanzapine	6.75	2.25			
Flupentixol	0.648	0.648			
Fluphenazine	0.590	0.590			
Haloperidol	0.630	0.630			
Levomeprazine	8.54	2.85			
Melperone	11.0	2.20			
Norclozapine	45.3	9.06			
Norquetiapine	8.71	2.90			
Olanzapine	7.14	2.38			
Paliperidone	6.70	2.23			
Perazine	23.9	7.97			
Pipamperone	30.9	6.18			
Promethazine	20.8	6.93			
Prothipendyl	1.08	1.08			
Quetiapine	30.0	3.00			
Risperidone	6.51	1.30			
Sertindole	10.3	5.15			
Sulpiride	54.7	5.47			
Thioridazine	19.1	3.82			
Ziprasidone	16.8	3.36			
Zotepine	9.21	4.61			
Zuclopenthixol	3.85	1.93			

Representative chromatograms for the LLOQ for clozapine, quetiapine, and perazine, and their corresponding internal standards are reported in Figure 1. Representative calibration curves for the same analytes are reported in Figure 2.

No significant carryover was observed for any of the analytes.

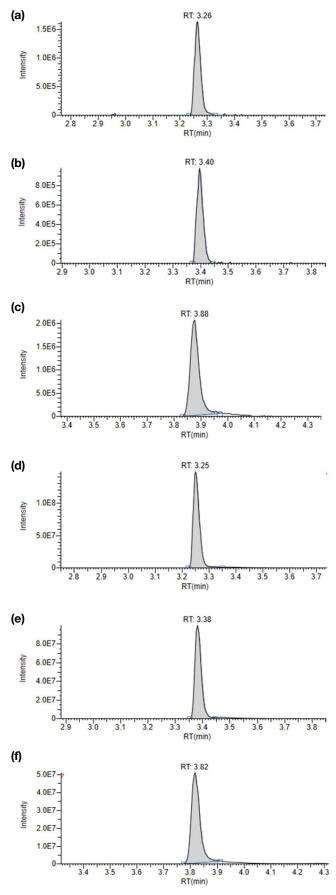


Figure 1. Representative chromatograms of the LLOQ for (a) clozapine, (b) quetiapine, (c) perazine, (d) clozapine- d_4 , (e) quetiapine- d_8 , and (f) perazine- d_8

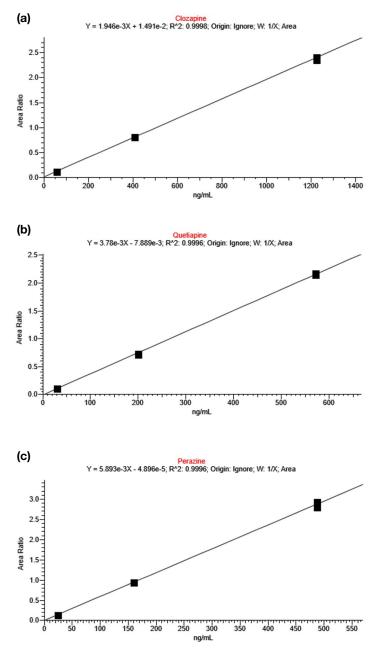


Figure 2. Representative calibration curves for (a) clozapine, (b) quetiapine, and (c) perazine

The data demonstrated good accuracy of the method with the percentage bias between nominal and average backcalculated concentration for the control samples ranging between -8.0% and 5.1% (Table 6). The %CV for intraassay precision was always below 8% for all the analytes, reported in Table 7. The maximum %CV for inter-assay precision including all the analytes was 7.9% (Table 8).

Table 6. Analytical accuracy results for control MS9382 batch #1279

		CTRL1	CTRL2			
Analyte	Nominal concentration (ng/L)	Average calculated concentration (ng/L)	Bias %	Nominal concentration (ng/L)	Average calculated concentration (ng/L)	Bias %
Amisulpride	135	136	0.9	317	299	-5.7
Aripiprazole	212	222	4.8	488	449	-8.0
Chlorpromazine	74.3	75.3	1.3	172	161	-6.1
Chlorprothixene	72.9	73.5	0.9	170	157	-7.7
Clozapine	235	247	5.1	551	513	-7.0
Dehydro-Aripiprazole	39.3	40.7	3.6	91.7	86.8	-5.3
Desmethylolanzapine	29.7	28.8	-2.9	68.5	63.1	-7.9
Flupentixol	2.69	2.60	-3.2	6.12	5.72	-6.6
Fluphenazine	2.44	2.48	1.8	5.55	5.28	-4.8
Haloperidol	2.56	2.61	2.0	6.04	5.65	-6.4
Levomepromazine	52.1	53.2	2.1	122	112	-8.1
Melperone	41.1	42.6	3.7	93.4	87.0	-6.8
Norclozapine	184	193	4.8	424	404	-4.7
Norquetiapine	70.2	73.1	4.1	168	156	-7.3
Olanzapine	29.9	29.2	-2.3	70.2	66.2	-5.7
Paliperidone	26.3	26.7	1.5	61.7	57.9	-6.2
Perazine	94.5	93.9	-0.7	217	203	-6.5
Pipamperone	125	126	1.1	288	270	-6.2
Promethazine	22.9	23.0	0.5	51.7	48.3	-6.7
Prothipendyl	11.9	12.1	1.7	28.1	26.2	-6.9
Quetiapine	142	143	0.8	322	300	-6.8
Riperidone	25.8	27.1	4.9	59.9	55.3	-7.6
Sertindole	43.4	44.3	2.0	102	95.5	-6.4
Sulpiride	225	231	2.6	529	495	-6.5
Thioridazine	80.6	84.4	4.7	194	180	-7.1
Ziprasidone	64.7	64.4	-0.4	154	144	-6.2
Zotepine	38.4	38.2	-0.6	87.9	81.3	-7.5
Zuclopenthixol	16.6	16.4	-1.0	39.0	36.6	-6.0

Table 7 (part 1). Analytical intra-assay precision results for control MS9382 (CTRL1) batch #1279

	CTRL1							
	Day 1		Day 2		Day 3			
Analyte	Average calculated concentration (ng/L)	CV %	Average calculated concentration (ng/L)	CV %	Average calculated concentration (ng/L)	CV %		
Amisulpride	135	5.6	144	1.4	129	3.1		
Aripiprazole	218	5.2	225	1.9	224	1.1		
Chlorpromazine	75.6	4.0	76.8	0.8	73.4	2.0		
Chlorprothixene	73.9	3.3	75.8	1.5	71.0	2.1		
Clozapine	245	3.2	255	0.6	241	1.9		
Dehydro-Aripiprazole	40.6	3.7	41.9	3.5	39.7	4.3		
Desmethylolanzapine	27.4	3.3	31.6	2.4	27.4	1.7		
Flupentixol	2.51	4.7	2.82	4.7	2.53	2.4		
Fluphenazine	2.41	5.6	2.62	1.5	2.43	3.1		
Haloperidol	2.56	5.2	2.70	2.4	2.58	3.2		
Levomepromazine	51.5	7.9	55.5	2.6	52.6	2.8		
Melperone	42.3	4.2	43.9	0.7	41.6	2.7		
Norclozapine	196	1.9	197	0.7	185	2.1		
Norquetiapine	70.4	7.4	73.6	1.3	75.2	2.8		
Olanzapine	27.9	3.6	32.3	1.0	27.5	1.2		
Paliperidone	25.3	5.2	28.6	1.3	25.9	1.8		
Perazine	92.4	4.0	98.3	1.8	90.9	3.3		
Pipamperone	128	6.1	129	4.5	122	5.4		
Promethazine	22.7	4.4	24.5	1.5	21.8	1.7		
Prothipendyl	11.8	6.3	12.5	2.2	12.1	4.2		
Quetiapine	141	5.4	147	0.8	141	2.8		
Riperidone	26.4	3.8	27.5	1.2	27.4	4.5		
Sertindole	43.2	3.7	46.5	1.3	43.1	1.5		
Sulpiride	223	7.2	240	0.9	229	4.0		
Thioridazine	83.6	6.0	85.7	1.9	83.8	2.1		
Ziprasidone	62.9	4.8	67.0	0.8	63.4	2.0		
Zotepine	37.1	3.7	40.8	1.4	36.6	1.8		
Zuclopenthixol	16.1	3.9	17.5	6.9	15.7	2.3		

Table 7 (part 2). Analytical intra-assay precision results for control MS9382 (CTRL2) batch #1279

	CTRL2							
	Day 1		Day 2		Day 3			
Analyte	Average calculated concentration (ng/L)	CV %	Average calculated concentration (ng/L)	CV %	Average calculated concentration (ng/L)	CV %		
Amisulpride	299	2.2	296	1.5	302	1.8		
Aripiprazole	441	0.5	454	1.1	451	1.1		
Chlorpromazine	163	1.6	163	1.0	158	1.2		
Chlorprothixene	160	1.7	156	1.5	155	0.7		
Clozapine	520	2.0	505	1.7	513	1.4		
Dehydro-Aripiprazole	85.0	3.1	90.8	4.4	84.5	1.2		
Desmethylolanzapine	63.9	1.1	62.8	1.2	62.7	1.3		
Flupentixol	5.77	2.0	5.84	1.8	5.54	1.3		
Fluphenazine	5.15	2.5	5.56	3.1	5.14	1.3		
Haloperidol	5.67	3.4	5.69	2.0	5.60	1.3		
Levomepromazine	113	2.6	112	1.0	112	1.0		
Melperone	86.8	1.4	86.3	1.3	87.9	2.1		
Norclozapine	423	3.9	393	1.4	397	1.5		
Norquetiapine	153	0.7	155	1.5	159	2.5		
Olanzapine	66.5	2.5	65.3	1.2	66.7	2.5		
Paliperidone	57.9	2.4	57.7	1.6	58.0	2.5		
Perazine	201	1.3	205	2.2	203	1.8		
Pipamperone	270	2.6	271	1.9	270	1.0		
Promethazine	48.0	1.2	48.3	1.3	48.5	2.4		
Prothipendyl	26.0	2.7	26.2	0.7	26.4	1.7		
Quetiapine	302	1.1	301	0.7	297	0.9		
Riperidone	54.6	0.7	56.0	1.1	55.4	1.7		
Sertindole	94.6	2.1	96.0	2.4	95.7	1.9		
Sulpiride	498	2.3	494	1.1	492	1.3		
Thioridazine	183	2.9	179	1.2	179	1.8		
Ziprasidone	143	2.1	144	2.1	147	2.0		
Zotepine	80.6	1.4	81.6	0.7	81.6	1.1		
Zuclopenthixol	36.3	1.7	37.2	2.3	36.5	1.1		

thermo scientific

	CTRL1		CTRL2	CTRL2		
Analyte	Average calculated concentration (ng/L)	CV %	Average calculated concentration (ng/L)	CV %		
Amisulpride	136	6.0	299	1.9		
Aripiprazole	222	3.3	449	1.5		
Chlorpromazine	75.3	3.1	161	1.8		
Chlorprothixene	73.5	3.6	157	1.9		
Clozapine	247	3.1	513	2.0		
Dehydro-Aripiprazole	40.7	4.2	86.8	4.5		
Desmethylolanzapine	28.8	7.5	63.1	1.4		
Flupentixol	2.60	6.7	5.72	2.8		
Fluphenazine	2.48	5.2	5.28	4.4		
Haloperidol	2.61	4.3	5.65	2.3		
Levomepromazine	53.2	5.7	112	1.6		
Melperone	42.6	3.6	87.0	1.7		
Norclozapine	193	3.5	404	4.1		
Norquetiapine	73.1	5.0	156	2.2		
Olanzapine	29.2	7.9	66.2	2.2		
Paliperidone	26.7	6.3	57.9	2.1		
Perazine	93.9	4.5	203	1.9		
Pipamperone	126	5.7	270	1.8		
Promethazine	23.0	5.7	48.3	1.7		
Prothipendyl	12.1	4.8	26.2	1.9		
Quetiapine	143	3.9	300	1.2		
Riperidone	27.1	3.7	55.3	1.5		
Sertindole	44.3	4.3	95.5	2.1		
Sulpiride	231	5.3	495	1.6		
Thioridazine	84.4	3.7	180	2.2		
Ziprasidone	64.4	4.0	144	2.3		
Zotepine	38.2	5.6	81.3	1.2		
Zuclopenthixol	16.4	6.8	36.6	2.0		

Table 8. Analytical inter-assay precision results for control MS9382 batch #1279

Conclusions

A robust, reproducible, and sensitive liquid chromatography-high resolution mass spectrometry method for clinical research for quantification of 28 neuroleptics in human plasma was developed, implemented, and analytically validated. Sample preparation is based on a rapid 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

For General Lab Use only. Not for use in diagnostic procedures. © 2021 Thermo Fisher Scientific Inc. All rights reserved. ClinMass is a registered trademark of RECIPE Chemicals + Instruments GmbH. 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 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. TN66065-EN 0821S

