

Quantification of 36 benzodiazepines in human plasma or serum by LC-HRAM(MS) for clinical research

Authors: Claudio De Nardi¹, Katharina Kern², Steffen Peters²

¹Thermo Fisher Scientific GmbH, Dreieich, Germany

²RECIPE Chemicals + Instruments GmbH, Munich, Germany

Keywords: Benzodiazepines, offline sample preparation, plasma, serum, Orbitrap, high-resolution mass spectrometry

Application benefits

- Increased accuracy of method by implementation of a comprehensive ClinMass[®] kit for sample preparation
- High-resolution mass spectrometry for improved selectivity
- Robust, sensitive hardware enables increased confidence in data
- Simple offline sample preparation by protein precipitation
- 36 benzodiazepines in a single quantitative method

Goal

Implementation of an analytical method for the quantification of 36 benzodiazepines in human plasma or serum on a Thermo Scientific™ Q Exactive™ Plus hybrid quadrupole-Orbitrap™ mass spectrometer.



Introduction

Benzodiazepines (BZP) are a group of psychoactive drugs with a broad range of therapeutic effects. These controlled substances are some of the most frequently prescribed medications worldwide and are frequently prone to abuse by those without valid medical prescriptions.

An analytical method for clinical research for the quantification of 36 benzodiazepines in human plasma or serum is reported in this study. A comprehensive list of analytes, corresponding internal standards and concentration ranges covered by the method are reported in Table 1. Most reported LC-MS analyses of the above-mentioned benzodiazepines involve triple quadrupole mass spectrometers, traditionally used for targeted, sensitive

quantitation assays. However, in this report we present LC-MS data acquired using high-resolution accurate-mass (HRAM) mass spectrometry leveraging Orbitrap technology. This report demonstrates the capability of HRAM for routine quantitation analyses in addition to its use for performing in-depth qualitative investigations.

Table 1. Analytes, internal standards, and concentration ranges

| Analyte | Internal standard | Conc. range (ng/mL) |
|------------------------|-------------------------------------|---------------------|
| 3-Hydroxybromazepam | 7-Aminoflunitrazepam-D ₇ | 16.6–258 |
| 7-Aminoclonazepam | 7-Aminoclonazepam-D ₄ | 5.02–70.3 |
| 7-Aminoflunitrazepam | 7-Aminoflunitrazepam-D ₇ | 5.36–76.5 |
| 7-Aminonitrazepam | 7-Aminoclonazepam-D ₄ | 21.6–304 |
| alpha-OH-Alprazolam | alpha-OH-Alprazolam-D ₅ | 5.74–81.1 |
| alpha-OH-Midazolam | alpha-OH-Midazolam-D ₄ | 10.6–135 |
| alpha-OH-Triazolam | alpha-OH-Triazolam-D ₄ | 3.92–63.5 |
| Alprazolam | Alprazolam-D ₅ | 5.14–76.4 |
| Bromazepam | alpha-OH-Alprazolam-D ₅ | 21.3–319 |
| Brotizolam | Temazepam-D ₅ | 2.00–30.0 |
| Chlordiazepoxide | Chlordiazepoxide-D ₅ | 260–3512 |
| Clobazam | Nordiazepam-D ₅ | 27.7–460 |
| Clonazepam | Clonazepam-D ₄ | 4.94–72.3 |
| Demoxepam | 7-Aminoflunitrazepam-D ₇ | 231–3524 |
| Desalkylflurazepam | Zolpidem-D ₆ | 17.1–274 |
| Desmethylflunitrazepam | Oxazepam-D ₅ | 4.78–84.3 |
| Diazepam | Diazepam-D ₅ | 128–1733 |
| Estazolam | Estazolam-D ₅ | 44.6–679 |
| Flunitrazepam | Flunitrazepam-D ₇ | 5.25–76.6 |
| Flurazepam | Midazolam-D ₄ | 10.4–155 |
| Lorazepam | Lorazepam-D ₄ | 19.2–276 |
| Lormetazepam | Midazolam-D ₄ | 1.97–27.2 |
| Medazepam | Prazepam-D ₅ | 75.3–1047 |
| Midazolam | Midazolam-D ₄ | 20.5–300 |
| Nitrazepam | Nitrazepam-D ₅ | 18.9–293 |
| Norclobazam | Estazolam-D ₅ | 240–4281 |
| Nordiazepam | Nordiazepam-D ₅ | 116–1653 |
| Oxazepam | Oxazepam-D ₅ | 117–1808 |
| Prazepam | Prazepam-D ₅ | 89.4–1259 |
| Temazepam | Temazepam-D ₅ | 115–1587 |
| Tetrazepam | Prazepam-D ₅ | 45.0–640 |
| Trazodone | Midazolam-D ₄ | 180–2396 |
| Triazolam | Triazolam-D ₄ | 3.46–43.8 |
| Zaleplone | 7-Aminoflunitrazepam-D ₇ | 17.3–253 |
| Zolpidem | Zolpidem-D ₆ | 36.6–603 |
| Zopiclone | Chlordiazepoxide-D ₅ | 16.5–279 |

Plasma or serum samples were processed by offline internal standard addition and protein precipitation. Extracted samples were injected onto a Thermo Scientific™ Vanquish™ Flex Duo UHPLC system connected to a Q Exactive Plus hybrid quadrupole-Orbitrap mass spectrometer with heated electrospray ionization (H-ESI II). Detection was performed by full scan coupled to data-dependent fragmentation (fullMS-ddMS²) using 21 deuterated internal standards. The full scan experiment was used for quantification, and fragmentation for confirmation. Method performance was evaluated using the ClinMass® TDM Platform with the ClinMass® Add-On Set for Benzodiazepines from RECIPE® Chemicals + Instruments GmbH (Munich, Germany) in terms of linearity of response within the calibration ranges, lower limit of quantification (LLOQ), carryover, accuracy, trueness of measurements, and intra- and inter-assay precision for each analyte.

Experimental

Target analytes

The concentration ranges covered by the calibrators (MS6013 batch #1069) used are reported in Table 1.

Sample preparation

Reagents included four calibrators (including blank) and two controls from RECIPE (MS6082 batch #1267), as well as 20 deuterated internal standards for the quantification. Samples of 50 µL of plasma or serum were protein-precipitated using 100 µL of precipitating solution containing the internal standards. Precipitated samples were vortex-mixed and centrifuged. 100 µL of supernatant were transferred to a clean plate or vial and diluted with 100 µL of a dilution solution provided with the kit (MS9022).

Liquid chromatography

A Vanquish Duo UHPLC system, a dual-channel instrument configured for both LC-only and online SPE applications (Figure 1), was used for chromatographic separation. The LC-only channel was used in this case, utilizing mobile phases and an analytical column provided by RECIPE. Details of the analytical method are reported in Table 2. Total runtime was 6.8 minutes.

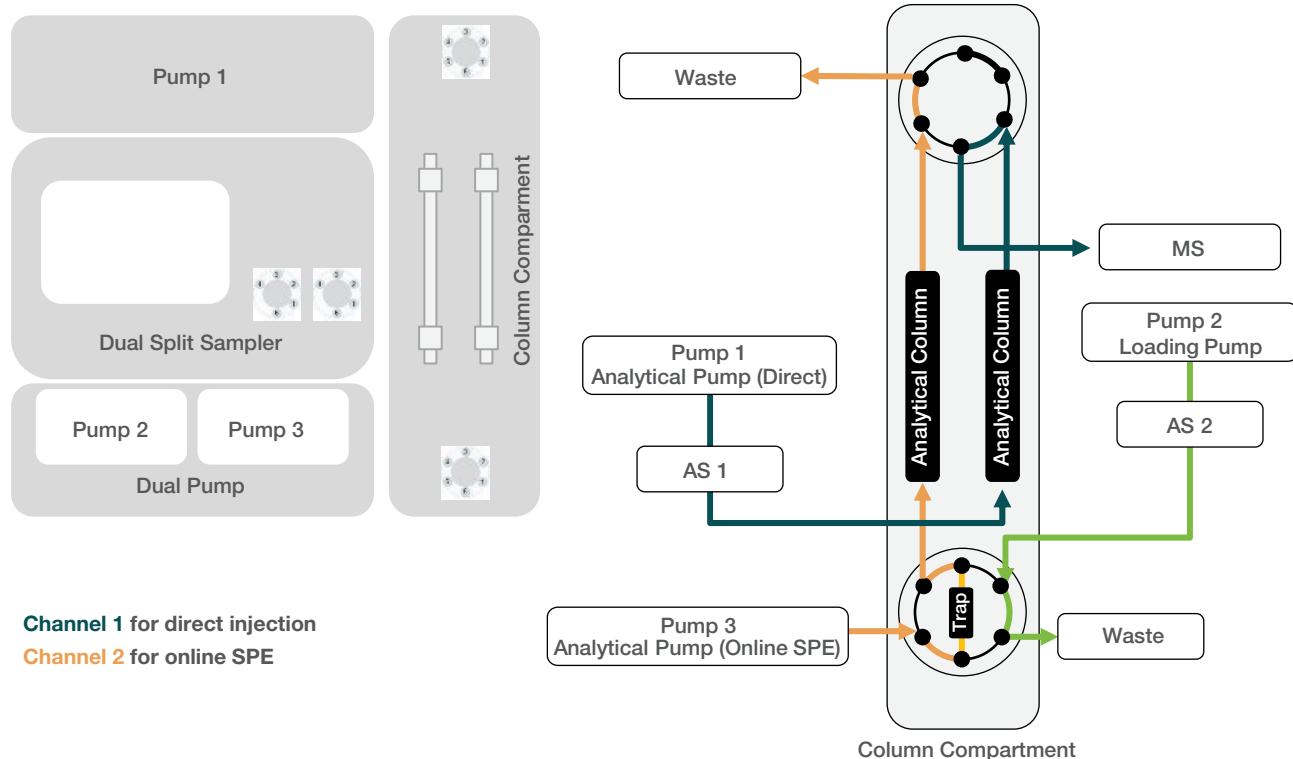


Figure 1. Schematic representation of the Vanquish Duo UHPLC system setup

Table 2. Liquid chromatography method description

| Gradient profile | | | | |
|--------------------|--------------------|-----------------|-------|--|
| Time (min) | Flow rate (mL/min) | A (%) | B (%) | |
| 0.00 | 0.6 | 90 | 10 | |
| 0.21 | 0.6 | 90 | 10 | |
| 0.30 | 0.6 | 76 | 24 | |
| 3.00 | 0.6 | 74 | 26 | |
| 6.20 | 0.6 | 40 | 60 | |
| 6.21 | 0.6 | 15 | 85 | |
| 6.40 | 0.6 | 15 | 85 | |
| 6.41 | 0.6 | 90 | 10 | |
| 6.80 | 0.6 | 90 | 10 | |
| Other parameters | | | | |
| Injection volume | | 15 μ L | | |
| Column temperature | | 40 $^{\circ}$ C | | |

Mass spectrometry

Analytes and internal standards were detected by FullIMS-ddMS² mode on a Q Exactive Plus hybrid quadrupole-Orbitrap mass spectrometer with heated electrospray ionization (H-ESI II) operated in positive ion mode. A summary of the MS conditions is reported in Table 3.

Table 3. MS settings

| Parameter | Value |
|---|---|
| Source type | Heated electrospray ionization (H-ESI II) |
| Vaporizer temperature | 450 $^{\circ}$ C |
| Capillary temperature | 275 $^{\circ}$ C |
| Spray voltage (positive mode) | 3000 V |
| Sheath gas | 58 AU |
| Sweep gas | 0 AU |
| Auxiliary gas | 16 AU |
| S-Lens RF level | 60 |
| Data acquisition mode | FullIMS-ddMS ² |
| FullIMS resolution @ m/z 200 | 70,000 |
| FullIMS scan range | 110 – 400 m/z |
| ddMS ² resolution @ m/z 200 | 17,500 |
| ddMS ² isolation window | 2.0 m/z |
| Stepped normalized collision energy (NCE) | 15, 25, 35 |

Method evaluation

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

To determine the LLOQ, a four-step serial dilution of the lowest calibrator using blank matrix was performed; a full set of calibrators (three levels), diluted calibrators (four levels), and controls (two levels) were extracted in replicates of five (n=5), injected in a single batch, and all used for the linear interpolation. The LLOQ was set as the lowest level that could be determined with a CV < 20% across the entire batch of samples.

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 back-calculated concentrations at two different levels using the quality control samples provided by RECIPE, prepared and analyzed in replicates of five on three different days.

Trueness of measurement was also evaluated as percentage bias using certified external quality controls (GTFCh BZF1-2019 A and B, Instand 876, 877 and 878 3-2019 Probes 31 and 32) prepared and analyzed in replicates of five on a single day.

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).

Data analysis

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

Results and discussion

A linear response (quadratic for trazodone) with 1/x weighting was obtained for all the analytes not only in the calibration range covered by the calibrators, but also down to a LLOQ reported in Table 4. The percentage bias between nominal and back-calculated concentration was always within $\pm 10\%$ for all the calibrators ($\pm 15\%$

for the lowest calibrator) in all the runs. Representative chromatograms for the LLOQ for midazolam, prazepam, and the corresponding internal standards are depicted in Figure 2. Representative calibration curves for the same analytes in the concentration range covered by the kit (three calibrators) are shown in Figure 3.

No significant carryover was registered, with no peak detected in the blank sample injected immediately after the highest calibrator.

Table 4. Analytes and corresponding LLOQ

| Analyte | LLOQ (ng/mL) |
|------------------------|--------------|
| 3-Hydroxybromazepam | 8.30 |
| 7-Aminoclonazepam | 0.251 |
| 7-Aminoflunitrazepam | 0.268 |
| 7-Aminonitrazepam | 1.08 |
| alpha-OH-Alprazolam | 2.87 |
| alpha-OH-Midazolam | 0.53 |
| alpha-OH-Triazolam | 1.96 |
| Alprazolam | 0.514 |
| Bromazepam | 10.7 |
| Brotizolam | 2.00 |
| Chlordiazepoxide | 13.0 |
| Clobazam | 13.9 |
| Clonazepam | 2.47 |
| Demoxepam | 11.6 |
| Desalkylflurazepam | 0.855 |
| Desmethylflunitrazepam | 2.39 |
| Diazepam | 6.40 |
| Estazolam | 2.23 |
| Flunitrazepam | 0.525 |
| Flurazepam | 0.52 |
| Lorazepam | 1.92 |
| Lormetazepam | 0.985 |
| Medazepam | 3.77 |
| Midazolam | 1.03 |
| Nitrazepam | 0.945 |
| Norclobazam | 120 |
| Nordiazepam | 5.80 |
| Oxazepam | 5.85 |
| Prazepam | 4.47 |
| Temazepam | 5.75 |
| Tetrazepam | 2.25 |
| Trazodone | 9.00 |
| Triazolam | 1.73 |
| Zaleplone | 0.865 |
| Zolpidem | 18.3 |
| Zopiclone | 1.65 |

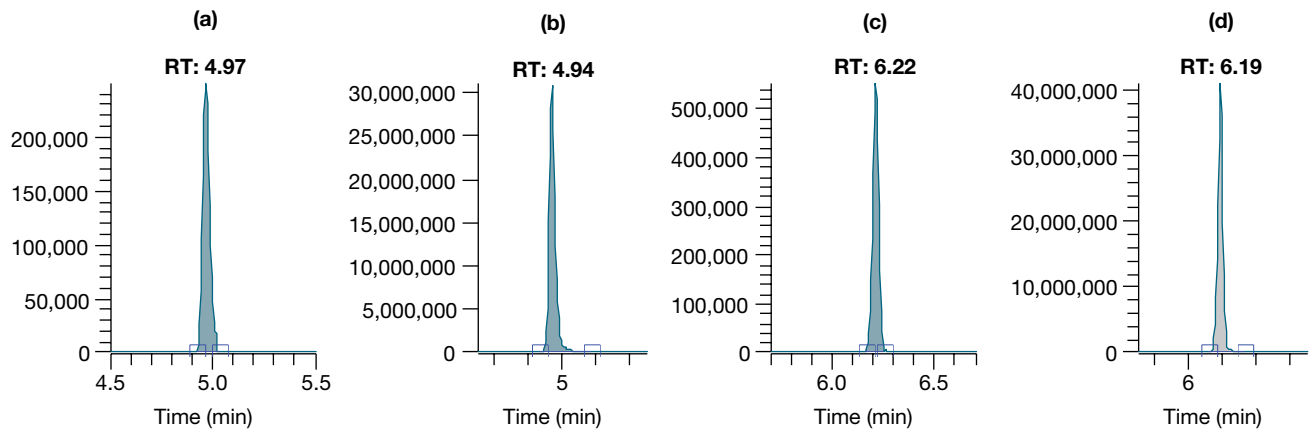


Figure 2. Representative chromatograms of the LLOQ for (a) midazolam, (b) midazolam-D₄, (c) prazepam, and (d) prazepam-D₅

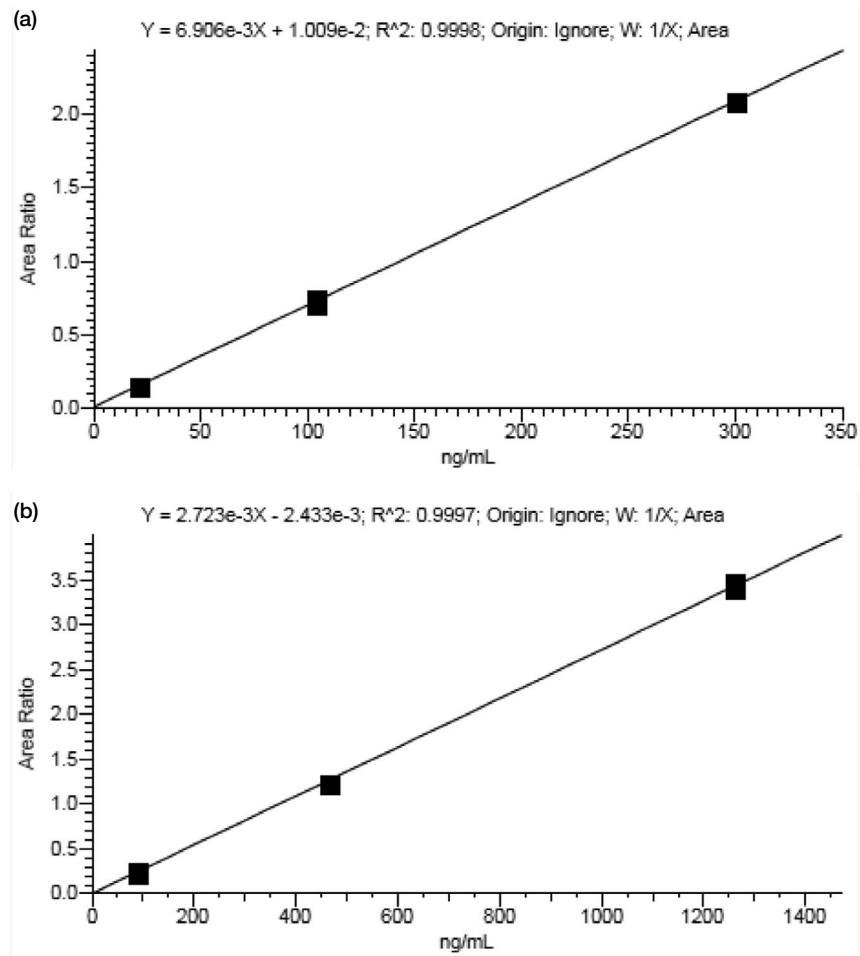


Figure 3. Representative calibration curves for (a) midazolam and (b) prazepam

The data presented in this report demonstrate the outstanding accuracy of the method with the percentage bias between nominal and average back-calculated concentration for the used control samples ranging between -13.2% and 8.6% (Table 5).

Table 5. Analytical accuracy results for controls MS6082 batch #1267

| 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 (%) |
| 3-Hydroxybromazepam | 42.4 | 45.1 | 6.4 | 144 | 156 | 8.6 |
| 7-Aminoclonazepam | 14.3 | 14.5 | 1.6 | 49 | 47.8 | -1.4 |
| 7-Aminoflunitrazepam | 15.0 | 15.7 | 4.7 | 50 | 50.8 | 2.7 |
| 7-Aminonitrazepam | 63.5 | 65.2 | 2.7 | 211 | 207 | -1.8 |
| alpha-OH-Alprazolam | 16.2 | 16.3 | 0.7 | 54.1 | 54.7 | 1.1 |
| alpha-OH-Midazolam | 55.0 | 53.7 | -2.3 | 177 | 174 | -1.5 |
| alpha-OH-Triazolam | 15.7 | 15.3 | -2.2 | 51.9 | 52.6 | 1.3 |
| Alprazolam | 16.9 | 15.5 | -8.2 | 57.2 | 52.3 | -8.5 |
| Bromazepam | 93.3 | 88.1 | -5.6 | 305 | 306 | 0.4 |
| Brotizolam | 4.56 | 4.12 | -9.6 | 15.4 | 14.0 | -8.9 |
| Chlordiazepoxide | 631 | 635 | 0.6 | 2053 | 1998 | -2.7 |
| Clobazam | 89.5 | 87.8 | -1.9 | 292 | 289 | -0.9 |
| Clonazepam | 8.13 | 7.90 | -2.8 | 61.3 | 58.9 | -4.0 |
| Demoxepam | 646 | 657 | 1.7 | 2189 | 2255 | 3.0 |
| Desalkylflurazepam | 29.9 | 29.5 | -1.4 | 101 | 103 | 1.7 |
| Desmethylflunitrazepam | 14.6 | 15.0 | 2.7 | 50.9 | 52.1 | 2.3 |
| Diazepam | 290 | 297 | 2.6 | 939 | 955 | 1.7 |
| Estazolam | 127 | 134 | 5.5 | 425 | 444 | 4.4 |
| Flunitrazepam | 16.2 | 15.2 | -6.2 | 54.2 | 51.1 | -5.6 |
| Flurazepam | 62.3 | 59.4 | -4.7 | 199 | 193 | -3.1 |
| Lorazepam | 65.6 | 56.9 | -13.2 | 203 | 186 | -8.3 |
| Lormetazepam | 5.65 | 5.38 | -4.8 | 18.3 | 18.2 | -0.5 |
| Medazepam | 274 | 288 | 5.2 | 838 | 879 | 4.9 |
| Midazolam | 30.2 | 30.0 | -0.6 | 78.9 | 81.1 | 2.8 |
| Nitrazepam | 45.3 | 41.3 | -8.9 | 148 | 140 | -5.6 |
| Norclobazam | 771 | 799 | 3.6 | 2733 | 2709 | -0.9 |
| Nordiazepam | 244 | 222 | -9.2 | 782 | 719 | -8.0 |
| Oxazepam | 360 | 355 | -1.3 | 1205 | 1197 | -0.6 |
| Prazepam | 271 | 259 | -4.6 | 866 | 842 | -2.7 |
| Temazepam | 202 | 185 | -8.4 | 562 | 501 | -10.8 |
| Tetrazepam | 126 | 131 | 4.2 | 418 | 427 | 2.2 |
| Trazodone | 509 | 496 | -2.5 | 1664 | 1509 | -9.3 |
| Triazolam | 7.55 | 7.96 | 5.5 | 24.1 | 25.0 | 3.8 |
| Zaleplone | 24.7 | 25.6 | 3.8 | 83.2 | 88.0 | 5.8 |
| Zolpidem | 128 | 124 | -3.1 | 426 | 409 | -4.0 |
| Zopiclone | 18.9 | 18.3 | -3.0 | 66.2 | 66.1 | -0.2 |

Excellent results were obtained also from the evaluation of trueness of measurement, with a percentage bias between -8.7% and 15.2% (Table 6 and Table 7).

Table 6. Analytical accuracy results for controls GTFCh BZF1-2019 and Instand 876 3-2019

| Analyte | GTFCh BZF1-2019 | | | | | | Instand 876 3-2019 | | | | | |
|------------------------|-----------------------|----------------------------------|----------|-----------------------|----------------------------------|----------|-----------------------|----------------------------------|----------|-----------------------|----------------------------------|----------|
| | Probe A | | | Probe B | | | Probe 31 | | | Probe 32 | | |
| | Nominal conc. (ng/mL) | Average calculated conc. (ng/mL) | Bias (%) | Nominal conc. (ng/mL) | Average calculated conc. (ng/mL) | Bias (%) | Nominal conc. (ng/mL) | Average calculated conc. (ng/mL) | Bias (%) | Nominal conc. (ng/mL) | Average calculated conc. (ng/mL) | Bias (%) |
| 7-Aminoflunitrazepam | N/A | N/A | N/A | 17.6 | 19.3 | 8.9 | N/A | N/A | N/A | N/A | N/A | N/A |
| Alprazolam | 46.0 | 47.3 | 2.8 | N/A | N/A | N/A | 32.1 | 32.8 | 2.0 | 10.7 | 10.9 | 1.4 |
| Bromazepam | 146 | 165 | 11.3 | N/A | N/A | N/A | N/A | N/A | N/A | N/A | N/A | N/A |
| Chlordiazepoxide | N/A | N/A | N/A | N/A | N/A | N/A | 2248 | 2417 | 7.0 | 634 | 704 | 9.9 |
| Clonazepam | 59.2 | 63.5 | 6.8 | N/A | N/A | N/A | N/A | N/A | N/A | N/A | N/A | N/A |
| Desalkylflurazepam | N/A | N/A | N/A | N/A | N/A | N/A | 68.4 | 77.3 | 11.5 | 38.6 | 36.8 | -5.0 |
| Desmethylflunitrazepam | N/A | N/A | N/A | 20.5 | 21.1 | 2.8 | N/A | N/A | N/A | N/A | N/A | N/A |
| Diazepam | 368 | 396 | 7.1 | N/A | N/A | N/A | N/A | N/A | N/A | N/A | N/A | N/A |
| Flunitrazepam | N/A | N/A | N/A | 14.8 | 15.1 | 2.2 | N/A | N/A | N/A | N/A | N/A | N/A |
| Lorazepam | N/A | N/A | N/A | 119 | 119 | 0.1 | N/A | N/A | N/A | N/A | N/A | N/A |
| Lormetazepam | N/A | N/A | N/A | 15.5 | 15.7 | 1.3 | N/A | N/A | N/A | N/A | N/A | N/A |
| Midazolam | N/A | N/A | N/A | 91.5 | 94.7 | 3.3 | N/A | N/A | N/A | N/A | N/A | N/A |
| Nitrazepam | 109 | 105 | -3.5 | N/A | N/A | N/A | N/A | N/A | N/A | N/A | N/A | N/A |
| Nordiazepam | 416 | 413 | -0.7 | N/A | N/A | N/A | 236 | 225 | -5.1 | 321 | 318 | -0.9 |
| Oxazepam | N/A | N/A | N/A | 292 | 289 | -0.9 | 639 | 667 | 4.2 | 1259 | 1343 | 6.2 |
| Trazodone | N/A | N/A | N/A | N/A | N/A | N/A | 661 | 657 | -0.6 | 1379 | 1275 | -8.2 |
| Zolpidem | 108 | 109 | 0.6 | N/A | N/A | N/A | N/A | N/A | N/A | N/A | N/A | N/A |
| Zopiclone | N/A | N/A | N/A | 59.1 | 69.7 | 15.2 | N/A | N/A | N/A | N/A | N/A | N/A |

Table 7. Analytical accuracy results for controls Instand 877 and 878 3-2019

| Analyte | Instand 877 3-2019 | | | | | | Instand 878 3-2019 | | | | | |
|---------------|-----------------------|----------------------------------|----------|-----------------------|----------------------------------|----------|-----------------------|----------------------------------|----------|-----------------------|----------------------------------|----------|
| | Probe 31 | | | Probe 32 | | | Probe 31 | | | Probe 32 | | |
| | Nominal conc. (ng/mL) | Average calculated conc. (ng/mL) | Bias (%) | Nominal conc. (ng/mL) | Average calculated conc. (ng/mL) | Bias (%) | Nominal conc. (ng/mL) | Average calculated conc. (ng/mL) | Bias (%) | Nominal conc. (ng/mL) | Average calculated conc. (ng/mL) | Bias (%) |
| Bromazepam | N/A | N/A | N/A | N/A | N/A | N/A | 167 | 186 | 10.3 | 106 | 112 | 5.0 |
| Clobazam | N/A | N/A | N/A | N/A | N/A | N/A | N/A | N/A | N/A | 85.4 | 83.2 | -2.6 |
| Clonazepam | 91.3 | 98.3 | 7.1 | 13.2 | 12.7 | -3.6 | N/A | N/A | N/A | N/A | N/A | N/A |
| Flunitrazepam | 43.7 | 44.8 | 2.5 | 30.4 | 31.1 | 2.1 | N/A | N/A | N/A | N/A | N/A | N/A |
| Flurazepam | 317 | 320 | 0.8 | 286 | 274 | -4.4 | N/A | N/A | N/A | N/A | N/A | N/A |
| Lorazepam | N/A | N/A | N/A | N/A | N/A | N/A | 21.4 | 22.1 | 3.2 | 12.5 | 13.4 | 6.5 |
| Nitrazepam | 94.6 | 88.3 | -7.2 | 151 | 139 | -8.7 | N/A | N/A | N/A | N/A | N/A | N/A |
| Temazepam | 613 | 591 | -3.7 | 104 | 103 | -0.9 | N/A | N/A | N/A | N/A | N/A | N/A |

The %CV for intra-assay precision was always below 9.1%.
The maximum %CV for inter-assay precision was 8.5%.
Results for intra- and inter-assay precision are reported in
Table 8 and Table 9, respectively.

Table 8. Intra-assay precision results for control MS6082 batch #1267

| Analyte | Control 1 | | | | | | Control 2 | | | | | |
|-----------------------|--|--------|--|--------|--|--------|--|--------|--|--------|--|--------|
| | Day 1 | | Day 2 | | Day 3 | | Day 1 | | Day 2 | | Day 3 | |
| | Average calculated concentration (ng/mL) | CV (%) | Average calculated concentration (ng/mL) | CV (%) | Average calculated concentration (ng/mL) | CV (%) | Average calculated concentration (ng/mL) | CV (%) | Average calculated concentration (ng/mL) | CV (%) | Average calculated concentration (ng/mL) | CV (%) |
| 3-Hydroxybromazepam | 45.1 | 2.6 | 45.8 | 2.2 | 44.5 | 2.9 | 154 | 2.1 | 155 | 2.6 | 159 | 2.0 |
| 7-Aminoclonazepam | 14.5 | 5.1 | 14.1 | 7.3 | 15.0 | 5.2 | 47.8 | 3.1 | 45.8 | 3.2 | 49.8 | 5.3 |
| 7-Aminoflunitrazepam | 15.2 | 2.8 | 16.0 | 2.3 | 15.8 | 2.5 | 48.5 | 2.4 | 52.1 | 4.2 | 51.9 | 3.5 |
| 7-Aminonitrazepam | 65.7 | 3.1 | 63.4 | 8.5 | 66.6 | 4.3 | 215 | 3.1 | 192 | 3.4 | 215 | 3.5 |
| alpha-OH-Alprazolam | 16.1 | 2.6 | 16.2 | 4.3 | 16.6 | 3.5 | 52.6 | 3.0 | 54.5 | 2.6 | 57.1 | 3.1 |
| alpha-OH-Midazolam | 55.1 | 4.2 | 52.2 | 2.6 | 54.2 | 3.0 | 179 | 2.1 | 173 | 0.6 | 171 | 1.4 |
| alpha-OH-Triazolam | 15.0 | 4.4 | 15.1 | 4.4 | 15.9 | 4.3 | 49.8 | 1.9 | 53.7 | 3.4 | 54.3 | 3.6 |
| Alprazolam | 15.9 | 2.7 | 15.4 | 2.0 | 15.4 | 3.5 | 53.8 | 2.7 | 51.7 | 2.6 | 51.4 | 1.8 |
| Bromazepam | 87.2 | 2.7 | 89.3 | 1.9 | 87.6 | 3.6 | 299 | 4.3 | 309 | 4.4 | 311 | 6.5 |
| Brotizolam | 4.11 | 4.6 | 4.01 | 4.4 | 4.25 | 4.3 | 13.6 | 1.6 | 14.6 | 6.0 | 13.8 | 2.7 |
| Chlordiazepoxide | 641 | 4.3 | 623 | 4.9 | 641 | 3.8 | 1979 | 1.7 | 1984 | 1.0 | 2029 | 2.2 |
| Clobazam | 90.5 | 4.7 | 84.2 | 7.5 | 89.2 | 4.4 | 295 | 4.1 | 277 | 3.6 | 297 | 2.2 |
| Clonazepam | 7.93 | 7.0 | 8.25 | 8.1 | 7.53 | 9.1 | 59.0 | 2.2 | 58.6 | 3.9 | 58.9 | 4.3 |
| Demoxepam | 660 | 2.0 | 669 | 6.7 | 643 | 3.6 | 2137 | 3.5 | 2366 | 3.2 | 2261 | 2.1 |
| Desalkylflurazepam | 29.9 | 3.6 | 29.6 | 4.3 | 29.1 | 2.9 | 100 | 5.0 | 105 | 3.8 | 103 | 2.9 |
| Desmethyflunitrazepam | 15.8 | 6.4 | 15.0 | 4.8 | 14.3 | 3.9 | 53.2 | 2.6 | 53.7 | 2.1 | 49.4 | 2.4 |
| Diazepam | 300 | 2.4 | 297 | 2.2 | 297 | 1.9 | 963 | 1.8 | 950 | 2.3 | 951 | 2.2 |
| Estazolam | 136 | 1.6 | 131 | 2.4 | 136 | 2.7 | 436 | 1.5 | 441 | 2.3 | 454 | 3.4 |
| Flunitrazepam | 15.4 | 3.2 | 15.3 | 2.7 | 14.9 | 3.7 | 50.5 | 3.8 | 51.7 | 3.0 | 51.2 | 3.7 |
| Flurazepam | 59.7 | 4.0 | 56.2 | 4.9 | 62.3 | 6.7 | 196 | 3.9 | 187 | 2.5 | 196 | 1.9 |
| Lorazepam | 56.5 | 0.9 | 57.2 | 2.1 | 57.0 | 1.5 | 181 | 1.7 | 190 | 2.5 | 187 | 2.3 |
| Lormetazepam | 5.65 | 1.6 | 5.30 | 4.7 | 5.23 | 5.1 | 17.8 | 3.1 | 18.0 | 4.4 | 18.9 | 5.4 |
| Medazepam | 286 | 4.7 | 274 | 7.7 | 304 | 1.6 | 883 | 1.3 | 857 | 3.2 | 897 | 3.7 |
| Midazolam | 29.8 | 5.2 | 29.5 | 4.7 | 30.7 | 5.9 | 81.9 | 2.4 | 80.6 | 3.7 | 80.8 | 2.2 |
| Nitrazepam | 41.2 | 4.2 | 41.6 | 4.4 | 40.9 | 4.8 | 142 | 5.2 | 145 | 1.6 | 133 | 1.8 |
| Norclobazam | 790 | 6.1 | 812 | 2.8 | 793 | 2.8 | 2811 | 5.4 | 2770 | 4.2 | 2546 | 4.4 |
| Nordiazepam | 220 | 6.2 | 227 | 4.7 | 218 | 3.0 | 715 | 2.4 | 719 | 1.8 | 724 | 2.2 |
| Oxazepam | 355 | 2.6 | 355 | 3.2 | 356 | 3.8 | 1212 | 5.3 | 1194 | 2.6 | 1186 | 2.3 |
| Prazepam | 264 | 1.3 | 258 | 2.1 | 255 | 2.0 | 837 | 1.8 | 858 | 1.5 | 832 | 1.5 |
| Temazepam | 181 | 0.4 | 183 | 2.5 | 190 | 3.0 | 497 | 1.4 | 500 | 2.9 | 507 | 3.3 |
| Tetrazepam | 131 | 2.2 | 130 | 2.4 | 134 | 3.4 | 420 | 2.2 | 430 | 3.4 | 432 | 1.4 |
| Trazodone | 496 | 3.2 | 483 | 1.6 | 509 | 5.3 | 1490 | 2.1 | 1490 | 3.5 | 1548 | 4.5 |
| Triazolam | 7.90 | 2.0 | 8.15 | 3.9 | 7.83 | 5.4 | 25.4 | 3.0 | 25.0 | 2.3 | 24.6 | 2.6 |
| Zaleplone | 24.5 | 5.0 | 26.9 | 3.8 | 25.3 | 3.0 | 84.9 | 5.8 | 86.7 | 3.6 | 92.4 | 2.3 |
| Zolpidem | 125 | 2.8 | 124 | 5.1 | 123 | 2.2 | 408 | 2.3 | 411 | 3.3 | 409 | 2.2 |
| Zopiclone | 19.2 | 1.3 | 17.4 | 7.2 | 18.6 | 2.4 | 67.4 | 3.4 | 64.6 | 3.7 | 66.3 | 2.9 |

Table 9. Inter-assay precision results for control MS6082 batch #1267

| Analyte | Control 1 | | Control 2 | |
|------------------------|--|--------|--|--------|
| | Average calculated concentration (ng/mL) | CV (%) | Average calculated concentration (ng/mL) | CV (%) |
| 3-Hydroxybromazepam | 45.1 | 2.7 | 156 | 2.6 |
| 7-Aminoclonazepam | 14.5 | 6.3 | 47.8 | 5.2 |
| 7-Aminoflunitrazepam | 15.7 | 3.1 | 50.8 | 4.7 |
| 7-Aminonitrazepam | 65.2 | 5.8 | 207 | 6.3 |
| alpha-OH-Alprazolam | 16.3 | 3.5 | 54.7 | 4.4 |
| alpha-OH-Midazolam | 53.7 | 3.8 | 174 | 2.4 |
| alpha-OH-Triazolam | 15.3 | 4.8 | 52.6 | 4.9 |
| Alprazolam | 15.5 | 3.0 | 52.3 | 3.1 |
| Bromazepam | 88.1 | 2.8 | 306 | 5.1 |
| Brotizolam | 4.12 | 4.8 | 14.0 | 4.9 |
| Chlordiazepoxide | 635 | 4.2 | 1998 | 2.0 |
| Clobazam | 87.8 | 6.2 | 289 | 4.4 |
| Clonazepam | 7.90 | 8.5 | 58.9 | 3.4 |
| Demoxepam | 657 | 4.7 | 2255 | 5.1 |
| Desalkylflurazepam | 29.5 | 3.5 | 103 | 4.2 |
| Desmethylflunitrazepam | 15.0 | 6.2 | 52.1 | 4.4 |
| Diazepam | 297 | 2.0 | 955 | 2.1 |
| Estazolam | 134 | 2.7 | 444 | 3.0 |
| Flunitrazepam | 15.2 | 3.2 | 51.1 | 3.4 |
| Flurazepam | 59.4 | 6.8 | 193 | 3.6 |
| Lorazepam | 56.9 | 1.6 | 186 | 2.9 |
| Lormetazepam | 5.38 | 5.2 | 18.2 | 4.9 |
| Medazepam | 288 | 6.6 | 879 | 3.3 |
| Midazolam | 30.0 | 5.2 | 81.1 | 2.8 |
| Nitrazepam | 41.3 | 4.2 | 140 | 4.9 |
| Norclobazam | 799 | 3.9 | 2709 | 6.2 |
| Nordiazepam | 222 | 4.7 | 719 | 2.1 |
| Oxazepam | 355 | 3.0 | 1197 | 3.5 |
| Prazepam | 259 | 2.2 | 842 | 2.0 |
| Temazepam | 185 | 3.1 | 501 | 2.6 |
| Tetrazepam | 131 | 2.9 | 427 | 2.6 |
| Trazodone | 496 | 4.2 | 1509 | 3.8 |
| Triazolam | 7.96 | 4.2 | 25.0 | 2.8 |
| Zaleplone | 25.6 | 5.4 | 88.0 | 5.3 |
| Zolpidem | 124 | 3.4 | 409 | 2.5 |
| Zopiclone | 18.3 | 5.9 | 66.1 | 3.6 |

Conclusions

An LC-HRAM-based method (a Vanquish Duo UHPLC system connected to a Q Exactive Plus hybrid quadrupole-Orbitrap MS) is reported here, demonstrating the power of Orbitrap technology in performing accurate qualitative analyses and routine quantitation with high efficiency. This method for clinical research was developed and implemented for the quantification of 36 benzodiazepines in human plasma or serum. The ClinMass TDM Platform with the ClinMass Add-On Set for Benzodiazepines from RECIPE was used. The method incorporates a quick and simple offline protein precipitation step 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/clinicalresearch