



Screening and quantitation of pesticide residues in rice using LC-(HESI)-MS/MS

Authors

Ramiz M.R. Azad and
Dasharath Oulkar
Customer Solution Center,
Ghaziabad,
Thermo Fisher Scientific, India

Keywords

TraceFinder, pesticide residues,
rice, QuEChERS, LC-MS/MS,
TSQ Quantis

Goal

The objective of this work was to develop a screening solution followed by quantitation of 160 pesticide residues in rice matrix using liquid chromatography-triple quadrupole mass spectrometry. The optimized method was validated in accordance with the SANTE guidelines and in compliance with the requirements of FSSAI and the European Commission (EC) MRLs.

Introduction

The Central Insecticide Board and Registration Committee (CIBRC) has few chemicals registered for rice.¹ The European Commission (EC) and FSSAI have set the maximum residue levels (MRLs) for many pesticides in rice at 0.01 mg/kg, but the MRLs for fipronil and fipronil sulfone are set at 0.005 mg/kg.^{2,3} The QuEChERS (Quick, Easy, Cheap, Effective, Rugged and Safe) method has been adopted for pesticide residue extraction in most food samples.⁴ Additionally, the instrument method plays an important role in delivering accurate and precise results to meet the regulatory requirements.

The aim of this work was the optimization and method validation of a multi-residue method for pesticides in rice by using LC-MS/MS with a Thermo Scientific™ TSQ Quantis™ triple quadrupole mass spectrometer. Sample

extraction was carried out using the acetic acid buffered version of the QuEChERS.⁴ The data acquisition and processing was carried out by using Thermo Scientific™ TraceFinder™ software. The optimized method was validated according to the SANTE/11813/ 2017 guidelines.⁵ This method was applied to real samples to demonstrate the application of streamlined workflow in compliance with the EU and FSSAI MRL requirements.

Experimental

Chemicals and apparatus

- Acetonitrile, Optima™ LC/MS Grade, Fisher Chemical
- Methanol, Optima™ LC/MS Grade, Fisher Chemical
- Water, Optima™ LC/MS Grade, Fisher Chemical
- Formic acid, Fisher Chemical
- Acetic acid, Fisher Chemical
- Ammonium formate, LC-MS Grade, Fisher Chemical
- Anhydrous magnesium sulfate, Thermo Fisher Scientific
- Sodium acetate, LR Grade, Fisher Chemical
- Certified reference materials
- Analytical balance (Aczet, CY2202, San Diego, CA) and precision balance (Aczet, CY205C, San Diego, CA)
- Vortex mixer (Thermo Scientific, P/N 88880017TS, also known as 88880017)
- Refrigerated centrifuge (Thermo Scientific™ Sorvall™ ST8 ventilated benchtop centrifuge)
- Variable volume micropipettes (Thermo Scientific)
- QuEChERS Salts (2007.01) Mylar Pouch 6 g magnesium sulfate (anhydrous), 1.5 g sodium acetate 50 pk Thermo Scientific™ (P/N 60105-341)

LC-MS/MS analysis

The Thermo Scientific™ Vanquish™ Flex Binary UHPLC system was coupled with the TSQ Quantis triple quadrupole mass spectrometer, which included the Heated Electrospray Ionization (HESI) source. The instrument was operated using an electrospray source in both positive/negative mode. The detailed conditions are given in Table 1. Multiple reaction monitoring (MRM) conditions were optimized for each pesticide during infusion (Appendix, Table 2).

Table 1. LC-MS/MS instrument conditions

| Liquid chromatography method | | | | |
|-------------------------------------|---|------------------|-----------|--------------|
| Instrumentation: | Vanquish Flex Binary UHPLC | | | |
| Column: | Thermo Scientific™ Hypersil GOLD™ (100 mm × 2.1 mm × 1.9 μm) (P/N 25002-102130) | | | |
| Sample compartment temp.: | 10 °C | | | |
| Column oven temp.: | 25 °C | | | |
| Mobile phase A: | 2 mM ammonium formate + 0.1% formic acid in water: acetonitrile (90:10, v/v) | | | |
| Mobile phase B: | 2 mM ammonium formate + 0.1% formic acid in water: acetonitrile (10:90, v/v) | | | |
| Total run time: | 18.0 min | | | |
| Gradient program: | <i>Time (min)</i> | <i>Flow Rate</i> | <i>%B</i> | <i>Curve</i> |
| | 0.0 | 0.400 | 1 | 5 |
| | 1.5 | 0.400 | 1 | 5 |
| | 5.0 | 0.400 | 5 | 5 |
| | 8.5 | 0.400 | 95 | 5 |
| | 13.5 | 0.400 | 95 | 5 |
| | 14.0 | 0.400 | 1 | 5 |
| | 18.0 | 0.400 | 1 | 5 |
| Mass spectrometry method | | | | |
| Instrumentation: | TSQ Quantis triple quadrupole tandem mass spectrometer | | | |
| Method type: | Time-based Selective-Reaction Monitoring (T-SRM) | | | |
| Ion source type: | H-ESI | | | |
| Polarity: | Positive/Negative switching | | | |
| Spray voltage: | Static Positive: 3500 V Negative: 2500 V | | | |
| Sheath gas: | 50 Arb | | | |
| Aux gas: | 10 Arb | | | |
| Sweep gas: | 1 Arb | | | |
| Ion transfer tube temp.: | 325 °C | | | |
| Vaporizer temp.: | 350 °C | | | |

Sample preparation

The rice grain matrix purchased from a local market was homogenized using a heavy-duty homogenizer to reduce the particle size to approximately 200 to 500 μm . The QuEChERS method was used for extraction as described below without cleanup.

Sample extraction

- Weigh 5 g homogenized sample into a 50 mL extraction tube (Note sample spiking at this step).
- Add 15 mL of HPLC grade water (containing 1% acetic acid) and leave the sample to soak for 10 min.
- Add 15 mL acetonitrile.
- Shake vigorously for 1 minute on a vortex mixer at 2500 rpm.
- Add 6 g anhydrous MgSO_4 and 1.5 g sodium acetate to the tube and again mix vigorously for 1 minute on a vortex mixer at 2500 rpm.
- Centrifuge with 5000 rpm for 5 min at ambient conditions.
- Filter the acetonitrile supernatant through a 0.2 μm PTFE membrane filter and dilute with HPLC grade water (1:1) before analyzing by LC-MS/MS.

Data acquisition and processing

The data acquisition and processing were carried out using TraceFinder software version 4.1. The data was acquired in Timed-SRM mode, which includes two or more transitions per analyte from the compound database (CDB). For data processing, the ion ratio ($\pm 30\%$), retention time (± 0.1 min), linearity (>0.99 with

residuals ± 20), recovery (70–120%) and precision ($\pm 20\%$) were set as user-defined filters as per SANTE guideline criteria.⁵

Results and discussion

Sample preparation

Rice is a dry powder and complex matrix that contains carbohydrates (9%) and proteins (5%). Rice has close to neutral pH (6.2–6.7), therefore acidification was maintained by using 1% acetic acid in water. It has very low moisture content, so water (1% acetic acid) was used, which is required for liquid-liquid partitioning with acetonitrile. Acidification improved the stability of base-sensitive compounds during extraction. No cleanup was applied, and the extract was diluted 6-fold as per the defined protocol to minimize the matrix effect without losses of target analytes. The final diluted extract, e.g. 0.01 mg/kg, corresponds to 0.0016 mg/kg.

LC-MS/MS analysis

Total LC-MS/MS method conditions were optimized for pesticide residues analysis, showing excellent sensitivity for 160 compounds. The total ion chromatogram (TIC) is shown in Figure 1. The optimized liquid chromatographic method offered excellent separation for the target analytes (spinetoram and spinosad D presented in Figure 2) and the absence of an isobaric interference from the matrix. In this method, the dwell time was automatically optimized in the range of 2–10 ms per transitions, which offered ≥ 12 points per peak. For early eluting compounds like 3-OH-carbofuran, this offered more than 12 points per peak as shown in Figure 3. This optimized instrument conditions provided excellent selectivity, repeatability, and reproducibility.

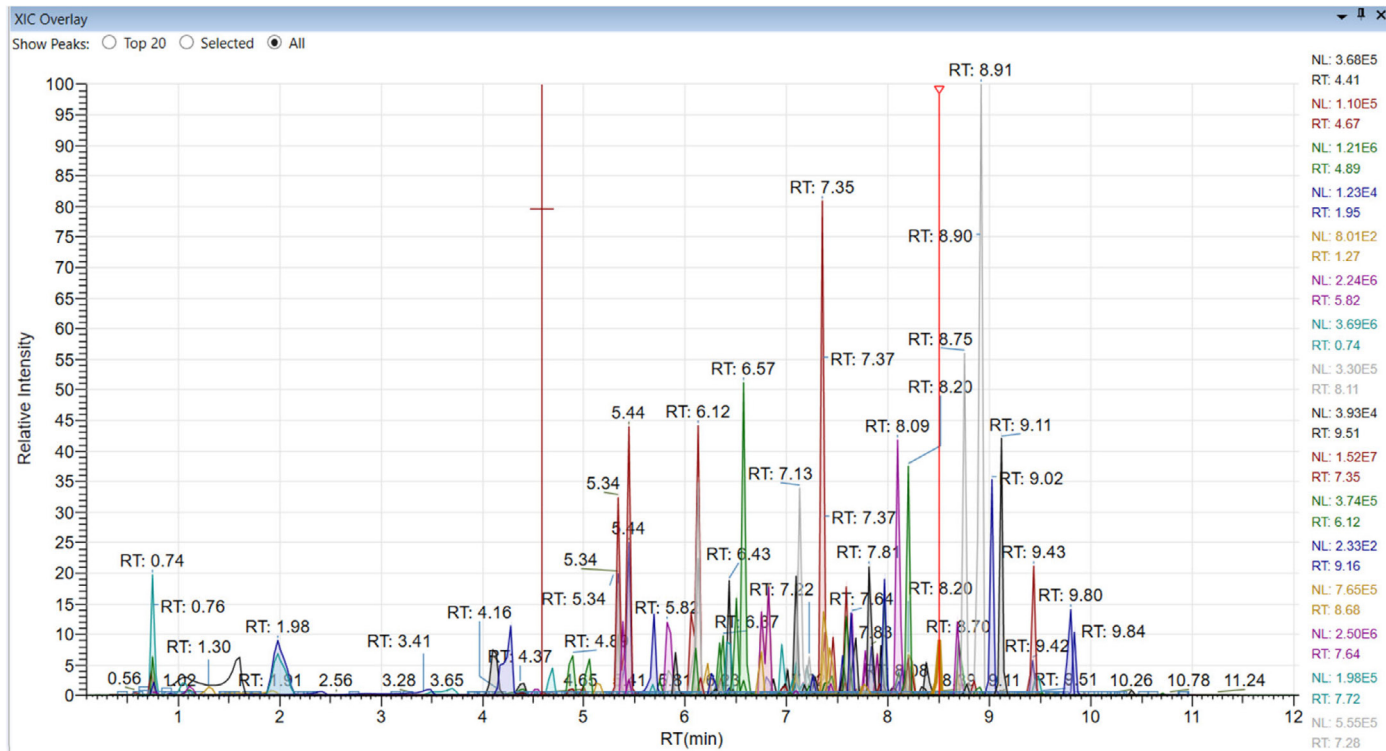


Figure 1. Extracted ion chromatogram for 160 compounds overlaid in a single window

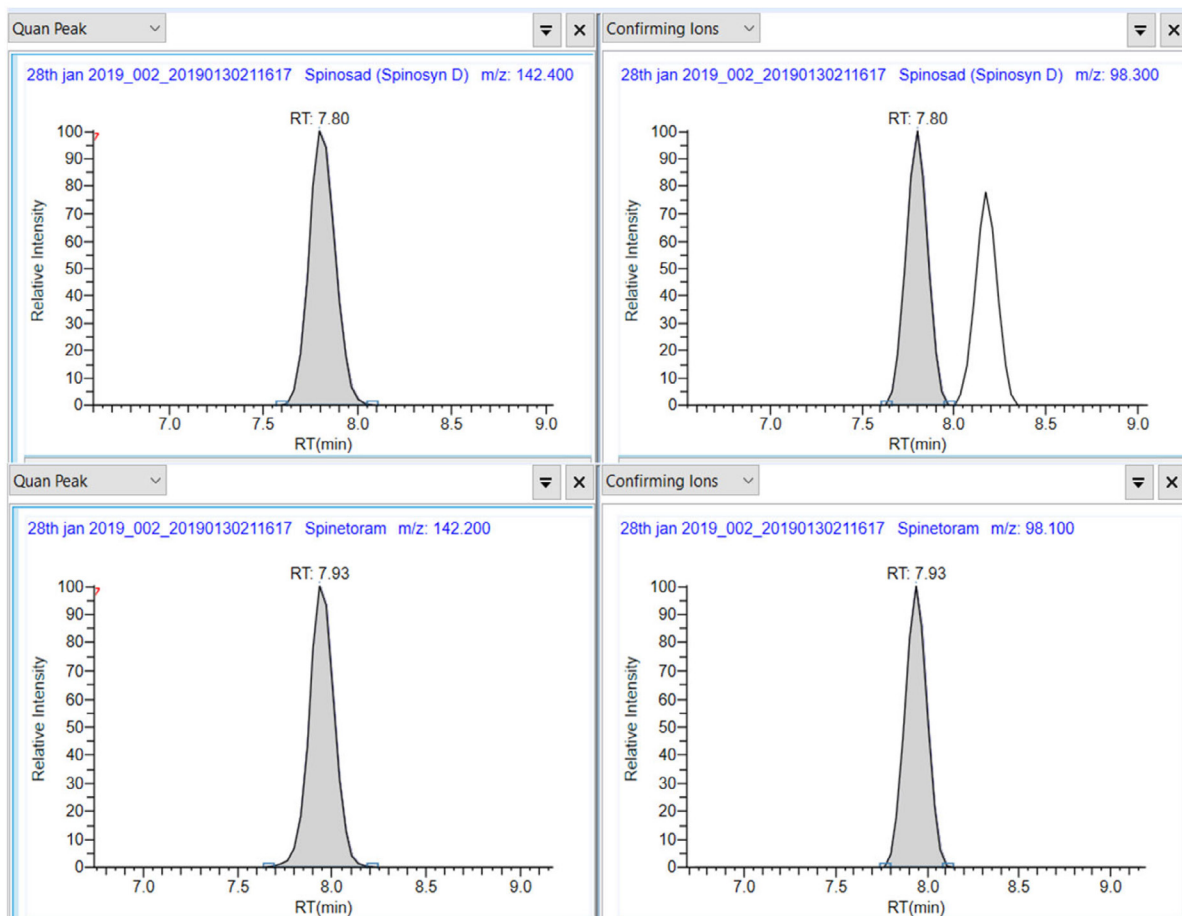
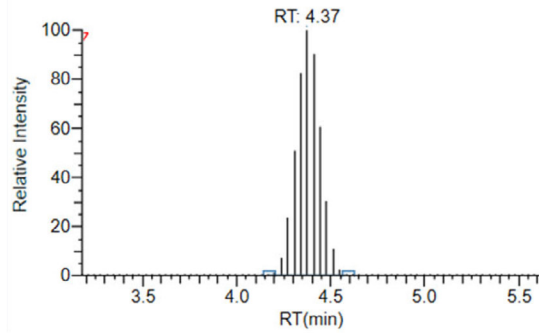


Figure 2. Effect of optimized gradient program on separation of isobaric compounds

28th jan 2019_008 3-Hydroxycarbofuran m/z: 181.000



28th jan 2019_008 3-Hydroxycarbofuran m/z: 163.000

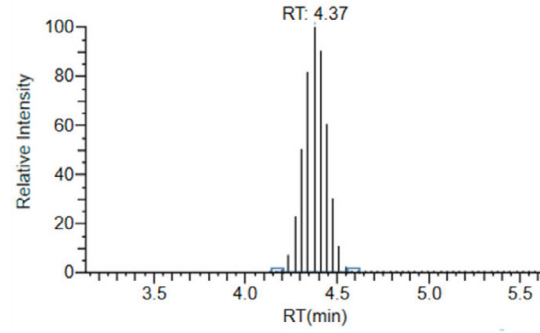


Figure 3. Impact of optimized dwell time on the data points per peak

Identification and quantitation

Parameters set in the TraceFinder Master Method for data processing included two transitions per analyte, retention time, correlation coefficient, and calibration graph residuals. Based on these parameters, the data was processed automatically with flagging. These color-coded flags indicated whether results passed or failed the acceptance criteria set in the processing method. The results passed SANTE guideline criteria as shown by the green colored flags (Figure 4).

In Figure 3, an identification of 3-hydroxycarbofuran in rice was demonstrated with two transitions 238.1→181.0 (quantitative) and 238.1→163.0 (confirmatory) at the same retention time (4.37 min, ±0.1), with observed ion ratio of 66.0% (45.21%–83.96%) in rice in comparison with the neat standard.

The linearity for 3-hydroxycarbofuran provided a correlation coefficient >0.999 with <15% residuals in compliance with the SANTE guidelines.

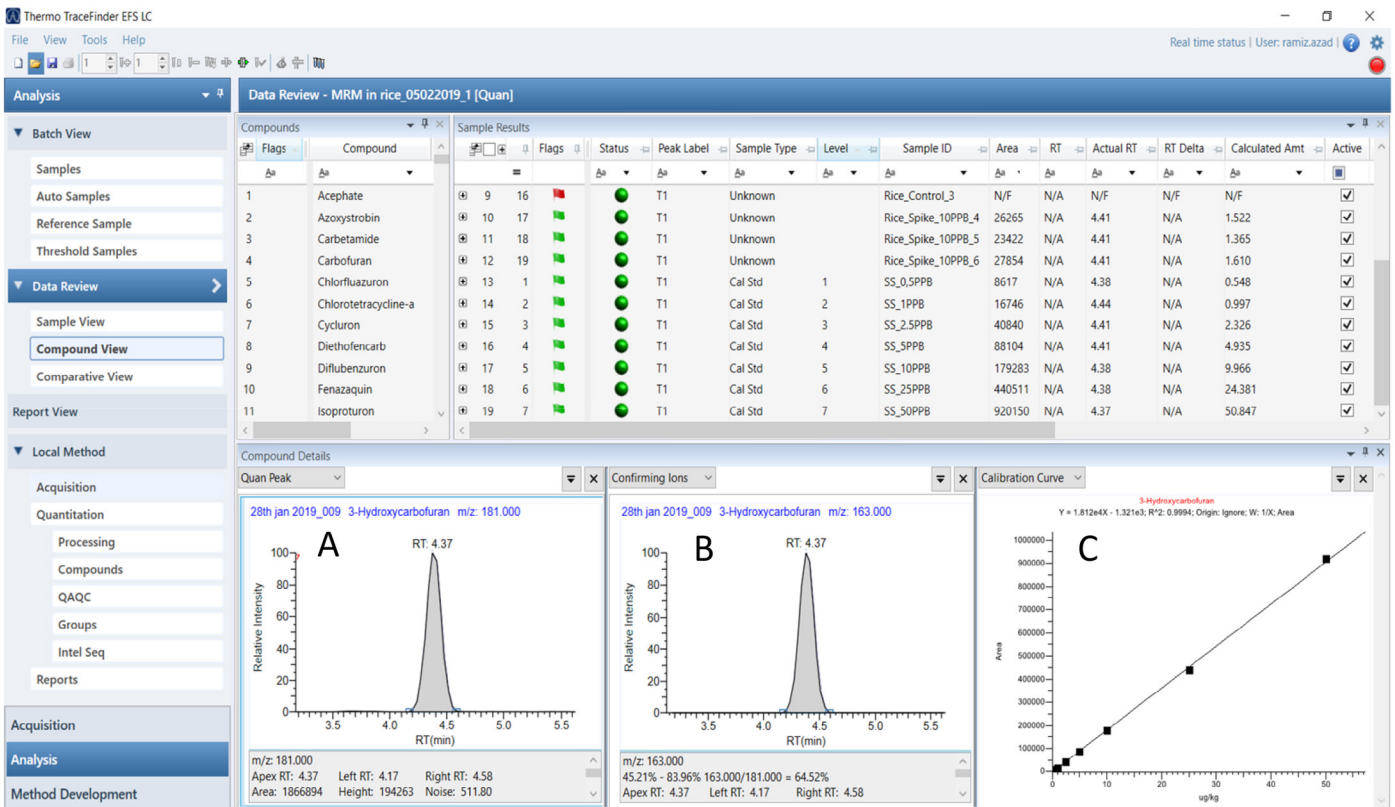


Figure 4. (A) Extracted ion chromatogram for quantifier ion of 3-hydroxycarbofuran, (B) identification based on the selectivity of confirmatory ions and confirmed with ion ratio, and (C) calibration curve

Method performance

The instrument detection limit was 0.05 ng/mL for 95% of the 160 target pesticides (Figure 5). Calibration standards were linear over the range of 0.0005 to 0.1 mg/kg with the correlation coefficient (>0.99) and residuals $<20\%$ for all the target analytes in both solvent and rice matrix. Because of the dilution of rice extracts, the LOQs in rice matrix were 0.01 mg/kg with acceptable recoveries (70%–120%) and precision ($<20\%$). The validation experiment was carried out by analysis of six replicates spiked at 0.01 (LOQ) and 0.05 mg/kg. Average recoveries were in the range of 76% to 116% with $<15\%$ RSD (Appendix, Table 3), which are in compliance with the SANTE criteria (recovery 70%–120% and precision $<20\%$).⁵ The repeatability of results obtained by analysis of an extended sequence (n=50 injection) was $<10\%$ for the area and $<\pm 0.05$ min retention time for acetamiprid and thiabendazole in Figures 6 and 7 (Figure 7 next page). This reveals that the optimized method offered excellent repeatability in results.

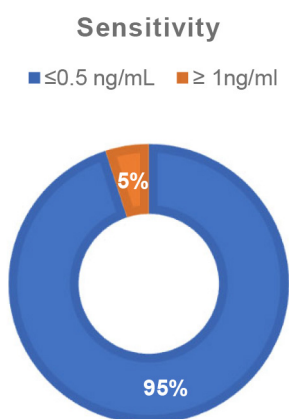


Figure 5. Sensitivity for the target list of analytes in the solvent standard

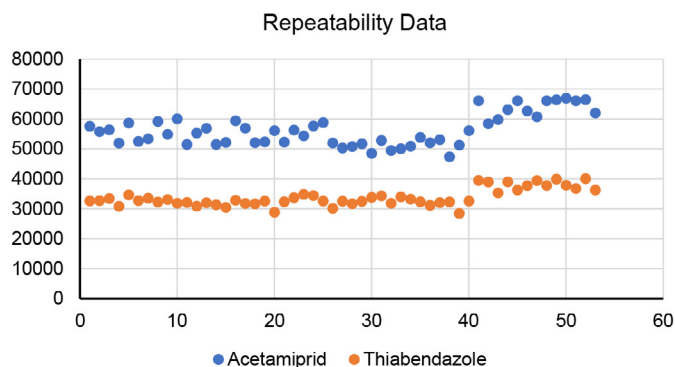


Figure 6. Area repeatability for acetamiprid and thiabendazole (n=54)

Conclusion

The developed method provides a solution for the simultaneous screening and quantitation of pesticide residues (160 pesticides) in rice by using a Thermo Scientific LC-HESI-MS/MS system. Use of the QuEChERS method for extraction followed by LC-MS/MS analysis could increase the overall high throughput of the commercial food testing laboratory. By following this approach at least 50 injections (standards, samples, blank) could be completed in a 24 h cycle. The optimized method results showed that LC separations in combination with SRM windows allowed maintaining the number of transitions monitored in single injection by auto-optimized dwell time without compromising data quality. This validated method data meets the SANTE guidelines.⁵ Also, this method complies with the EU and FSSAI MRL requirements by achieving an excellent lower limit of quantitation (LOQ).

References

1. Insecticides / Pesticides Registered under section 9(3) of the Insecticides Act, 1968 for use in the Country (as on 31/12/2018) <http://ppqs.gov.in/insecticides-pesticides-registered-under-section-93-insecticides-act-1968-use-country-31122018>
2. FSSAI Manual for food safety, 17th Edition, 2017 (THE FOOD SAFETY AND STANDARDS ACT, 2006)
3. EU Pesticides Database. <http://ec.europa.eu/food/plant/pesticides/eu-pesticides-database/public/?event=product.resultat&language=EN&selectedID=237>
4. Lehotay, S.J. Determination of Pesticide Residues in Foods by Acetonitrile Extraction and Partitioning with Magnesium Sulfate: Collaborative Study. *J. AOAC Int.* **2007**, *90*(2), 485–520.
5. SANTE guidelines. https://ec.europa.eu/food/sites/food/files/plant/docs/pesticides_mrl_guidelines_wrkdoc_2017-11813.pdf

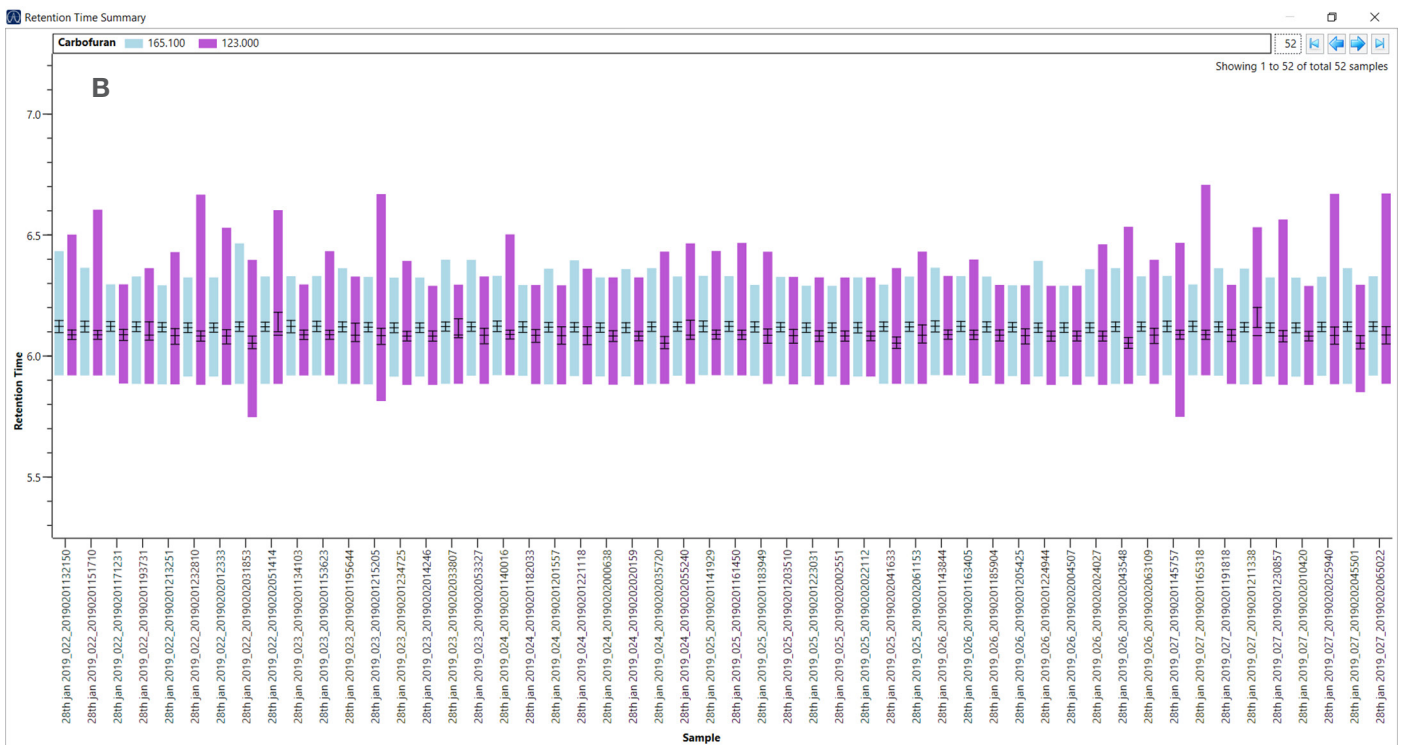
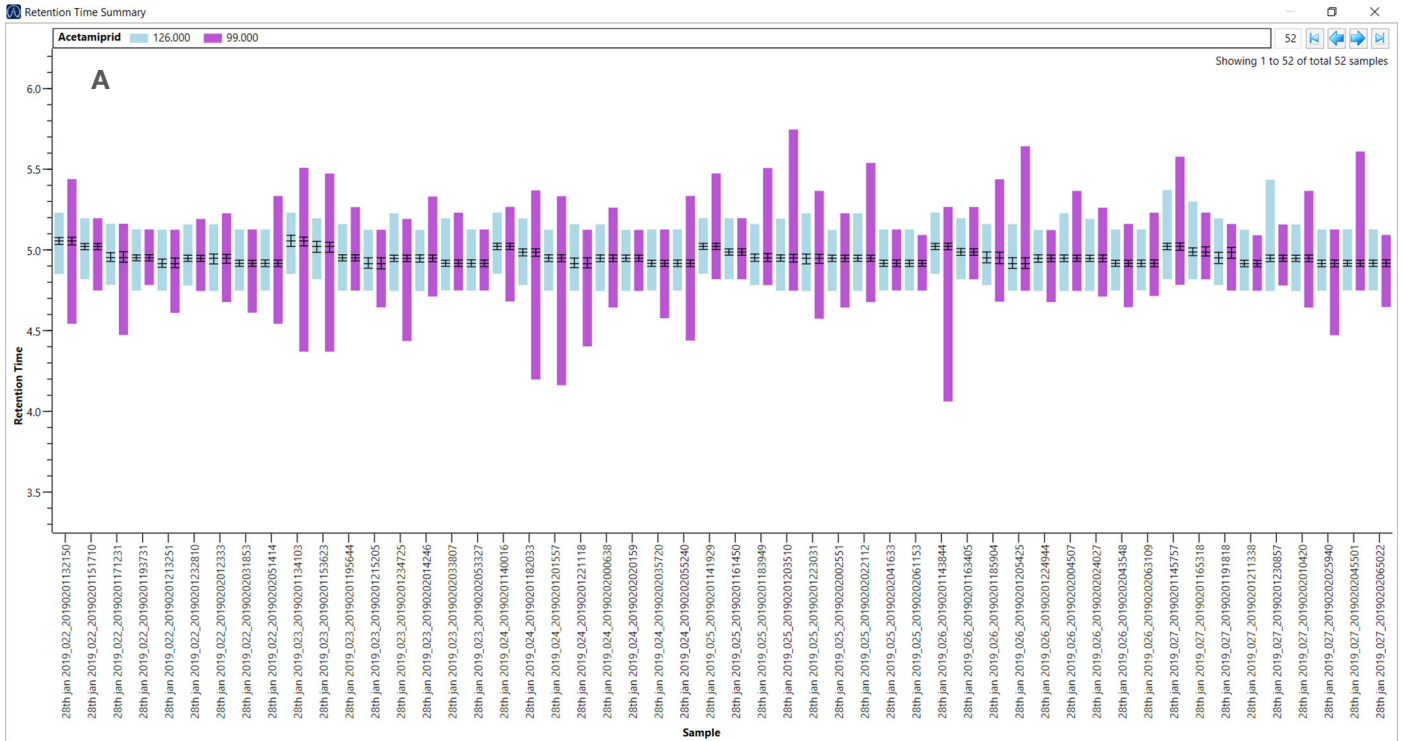


Figure 7. Retention time repeatability for acetamidrid (A) and carbofuran (B) (n=52)

Appendix

Table 2A. List of pesticides with MRM transitions used

| Compound | RT (min) | Polarity | Precursor (m/z) | Product (m/z) | CE (V) |
|---------------------|----------|----------|-----------------|---------------|--------|
| 3-Hydroxycarbofuran | 3.89 | Positive | 238.1 | 163 | 19 |
| 3-Hydroxycarbofuran | 3.89 | Positive | 238.1 | 181 | 15 |
| Acephate | 4.17 | Positive | 184.1 | 49 | 35 |
| Acephate | 4.17 | Positive | 184.1 | 143 | 11 |
| Acetamiprid | 4.51 | Positive | 223 | 99 | 53 |
| Acetamiprid | 4.51 | Positive | 223 | 126 | 29 |
| Aldicarb sulfone | 1.46 | Positive | 240.1 | 86.2 | 28 |
| Aldicarb sulfone | 1.46 | Positive | 240.1 | 148.2 | 19 |
| Aldicarb sulfoxide | 0.98 | Positive | 207.1 | 89.1 | 19 |
| Aldicarb sulfoxide | 0.98 | Positive | 207.1 | 132.1 | 9 |
| Ametryn | 5.75 | Positive | 228.1 | 96 | 35 |
| Ametryn | 5.75 | Positive | 228.1 | 186.1 | 25 |
| Aminocarb | 0.94 | Positive | 209.1 | 137.1 | 33 |
| Aminocarb | 0.94 | Positive | 209.1 | 152 | 19 |
| Amitraz | 7.71 | Positive | 294.2 | 91.2 | 57 |
| Amitraz | 7.71 | Positive | 294.2 | 148.3 | 22 |
| Avermectin Ba | 8.95 | Positive | 890.5 | 305 | 35 |
| Avermectin Ba | 8.95 | Positive | 890.5 | 567.5 | 17 |
| Azoxystrobin | 6.96 | Positive | 404.1 | 344.1 | 33 |
| Azoxystrobin | 6.96 | Positive | 404.1 | 372.1 | 19 |
| Benalaxyl | 7.72 | Positive | 326.2 | 148.1 | 29 |
| Benalaxyl | 7.72 | Positive | 326.2 | 294.1 | 15 |
| Bendiocarb | 5.98 | Positive | 224.1 | 109 | 33 |
| Bendiocarb | 5.98 | Positive | 224.1 | 167.1 | 15 |
| Benzoximate | 8.1 | Positive | 364 | 105 | 31 |
| Benzoximate | 8.1 | Positive | 364 | 199 | 11 |
| Bifenazate | 7.64 | Positive | 301.1 | 170.1 | 27 |
| Bifenazate | 7.64 | Positive | 301.1 | 198.1 | 13 |
| Bitertanol | 7.35 | Positive | 338.2 | 70 | 29 |
| Bitertanol | 7.35 | Positive | 338.2 | 269.2 | 13 |
| Boscalid | 7.07 | Positive | 343 | 140 | 25 |
| Boscalid | 7.07 | Positive | 343 | 307 | 27 |
| Bromucanozole | 7.04 | Positive | 378 | 70 | 47 |
| Bromucanozole | 7.04 | Positive | 378 | 159 | 37 |
| Bupirimate | 7.13 | Positive | 317 | 108 | 35 |
| Bupirimate | 7.13 | Positive | 317 | 159.1 | 33 |
| Buprofezin | 8.28 | Positive | 306.2 | 116.2 | 23 |
| Buprofezin | 8.28 | Positive | 306.2 | 201.1 | 17 |
| Butafenacil | 7.6 | Positive | 492.1 | 331 | 27 |
| Butafenacil | 7.6 | Positive | 492.1 | 349 | 19 |

Continued on next page

Table 2B. List of pesticides with MRM transitions used (continued from previous page)

| Compound | RT (min) | Polarity | Precursor (m/z) | Product (m/z) | CE (V) |
|----------------------|-----------------|-----------------|------------------------|----------------------|---------------|
| Butoxycarboxim | 1.87 | Positive | 223.1 | 106 | 13 |
| Butoxycarboxim | 1.87 | Positive | 223.1 | 159 | 11 |
| Carbendazim | 1.64 | Positive | 192.2 | 132.1 | 41 |
| Carbendazim | 1.64 | Positive | 192.2 | 160.2 | 25 |
| Carbetamide | 5.02 | Positive | 237.1 | 118.1 | 17 |
| Carbetamide | 5.02 | Positive | 237.1 | 192 | 13 |
| Carbofuran | 5.98 | Positive | 222.1 | 123 | 29 |
| Carbofuran | 5.98 | Positive | 222.1 | 165.1 | 17 |
| Carboxin | 5.98 | Positive | 236.1 | 87 | 33 |
| Carboxin | 5.98 | Positive | 236.1 | 143 | 21 |
| Chlorantraniliprole | 6.47 | Positive | 484 | 285.9 | 17 |
| Chlorantraniliprole | 6.47 | Positive | 484 | 452.9 | 21 |
| Chlorotoluron | 5.97 | Positive | 213.1 | 46.2 | 35 |
| Chlorotoluron | 5.97 | Positive | 213.1 | 72.2 | 31 |
| Chloroxuron | 6.99 | Positive | 291.1 | 72.4 | 47 |
| Chloroxuron | 6.99 | Positive | 291.1 | 218.1 | 33 |
| Clethodim | 8.27 | Positive | 360.101 | 164 | 29 |
| Clethodim | 8.27 | Positive | 360.101 | 268.1 | 17 |
| Clothianidin | 3.86 | Positive | 250 | 132 | 21 |
| Clothianidin | 3.86 | Positive | 250 | 169 | 19 |
| Cyazofamid | 7.72 | Positive | 325.2 | 108 | 18 |
| Cyazofamid | 7.72 | Positive | 325.2 | 261.2 | 14 |
| Cycluron | 6.24 | Positive | 199.1 | 89 | 21 |
| Cycluron | 6.24 | Positive | 199.1 | 89.1 | 21 |
| Cyproconazole Isomer | 6.79 | Positive | 292 | 70 | 31 |
| Cyproconazole Isomer | 6.79 | Positive | 292 | 125 | 29 |
| Cyprodinil | 7.29 | Positive | 226 | 77 | 61 |
| Cyprodinil | 7.29 | Positive | 226 | 93 | 47 |
| Desmedipham | 6.73 | Positive | 318.1 | 154 | 35 |
| Desmedipham | 6.73 | Positive | 318.1 | 182 | 19 |
| Diclobutrazol | 7.54 | Positive | 328.2 | 59.1 | 48 |
| Diclobutrazol | 7.54 | Positive | 328.2 | 70.2 | 48 |
| Diclotophos | 2.78 | Positive | 238.1 | 112.1 | 17 |
| Diclotophos | 2.78 | Positive | 238.1 | 193 | 13 |
| Diethofencarb | 6.75 | Positive | 268.1 | 124 | 43 |
| Diethofencarb | 6.75 | Positive | 268.1 | 226.1 | 13 |
| Difenoconazole | 7.76 | Positive | 406.1 | 251.1 | 41 |
| Difenoconazole | 7.76 | Positive | 408.2 | 253.1 | 31 |
| Dimethoate | 4.15 | Positive | 230 | 125 | 29 |
| Dimethoate | 4.15 | Positive | 230 | 199 | 13 |
| Dimethomorph Isomer | 6.38 | Positive | 388.1 | 165.1 | 45 |
| Dimethomorph Isomer | 6.38 | Positive | 388.1 | 301 | 29 |

Continued on next page

Table 2C. List of pesticides with MRM transitions used (continued from previous page)

| Compound | RT (min) | Polarity | Precursor (m/z) | Product (m/z) | CE (V) |
|------------------------|----------|----------|-----------------|---------------|--------|
| Dimethomorph Isomer _1 | 6.6 | Positive | 388.101 | 165.1 | 45 |
| Dimethomorph Isomer _1 | 6.6 | Positive | 388.101 | 301 | 29 |
| Dimoxystrobin | 7.41 | Positive | 327.1 | 116 | 29 |
| Dimoxystrobin | 7.41 | Positive | 327.1 | 205 | 23 |
| Diniconazole | 7.61 | Positive | 326.2 | 70.2 | 50 |
| Diniconazole | 7.61 | Positive | 326.2 | 159 | 43 |
| Dinotefuran | 1.07 | Positive | 203.1 | 129.2 | 17 |
| Dinotefuran | 1.07 | Positive | 203.1 | 157.2 | 11 |
| Dioxacarb | 4.11 | Positive | 224.1 | 123 | 21 |
| Dioxacarb | 4.11 | Positive | 224.1 | 167 | 11 |
| Diuron | 6.21 | Positive | 233.1 | 72 | 37 |
| Diuron | 6.21 | Positive | 235.1 | 72.1 | 33 |
| Doramectin | 9.31 | Positive | 916.6 | 331.4 | 33 |
| Doramectin | 9.31 | Positive | 916.6 | 593.5 | 19 |
| Emamectin-benzoate b1a | 8.54 | Positive | 886.5 | 82.1 | 65 |
| Emamectin-benzoate b1a | 8.54 | Positive | 886.5 | 158.1 | 49 |
| Epoxiconazole | 7.1 | Positive | 330 | 101.1 | 65 |
| Epoxiconazole | 7.1 | Positive | 330 | 121.1 | 49 |
| Eprinomectin | 8.65 | Positive | 914.6 | 154.2 | 49 |
| Eprinomectin | 8.65 | Positive | 914.6 | 186.2 | 25 |
| Etaconazole | 7.17 | Positive | 328.1 | 159 | 31 |
| Etaconazole | 7.17 | Positive | 328.1 | 205 | 23 |
| Ethirimol | 4.2 | Positive | 210.2 | 98.1 | 39 |
| Ethirimol | 4.2 | Positive | 210.2 | 140.1 | 31 |
| Etoxazole | 8.9 | Positive | 360.1 | 57.2 | 51 |
| Etoxazole | 8.9 | Positive | 360.1 | 141 | 59 |
| Fenamidone | 6.96 | Positive | 312.1 | 92 | 35 |
| Fenamidone | 6.96 | Positive | 312.1 | 236.1 | 21 |
| Fenarimol | 6.9 | Positive | 331 | 81 | 55 |
| Fenarimol | 6.9 | Positive | 331 | 268 | 35 |
| Fenazaquin | 9.69 | Positive | 307.1 | 147 | 25 |
| Fenazaquin | 9.69 | Positive | 307.1 | 161.1 | 27 |
| Fenbuconazole | 7.31 | Positive | 337 | 70 | 39 |
| Fenbuconazole | 7.31 | Positive | 337 | 124.9 | 55 |
| Fenpropimorph | 7.23 | Positive | 304 | 117 | 65 |
| Fenpropimorph | 7.23 | Positive | 304 | 147 | 39 |
| Fenpyroximate | 8.92 | Positive | 422 | 135.1 | 53 |
| Fenpyroximate | 8.92 | Positive | 422 | 366.1 | 23 |
| Fenuron | 3.8 | Positive | 165.1 | 46 | 29 |
| Fenuron | 3.8 | Positive | 165.1 | 72.1 | 45 |
| Fipronil | 7.7 | Negative | 435 | 250 | 35 |
| Fipronil | 7.7 | Negative | 435 | 330 | 20 |

Continued on next page

Table 2D. List of pesticides with MRM transitions used (continued from previous page)

| Compound | RT (min) | Polarity | Precursor (m/z) | Product (m/z) | CE (V) |
|-----------------|----------|----------|-----------------|---------------|--------|
| Fluazinam | 8.63 | Positive | 465 | 91 | 43 |
| Fluazinam | 8.63 | Positive | 465 | 148.9 | 43 |
| Flubendiamide | 7.63 | Positive | 683.1 | 274.1 | 41 |
| Flubendiamide | 7.63 | Positive | 683.1 | 408 | 9 |
| Flufenacet | 7.42 | Positive | 364.1 | 152.1 | 27 |
| Flufenacet | 7.42 | Positive | 364.1 | 194.2 | 17 |
| Fluometuron | 5.95 | Positive | 233.1 | 46 | 37 |
| Fluometuron | 5.95 | Positive | 233.1 | 72.1 | 35 |
| Fluoxastrobin | 7.46 | Positive | 459.2 | 188 | 47 |
| Fluoxastrobin | 7.46 | Positive | 459.2 | 427.2 | 23 |
| Flusilazole | 7.34 | Positive | 316.1 | 165.1 | 37 |
| Flusilazole | 7.34 | Positive | 316.1 | 247.1 | 21 |
| Flutolanil | 7.38 | Positive | 324.1 | 242.1 | 35 |
| Flutolanil | 7.38 | Positive | 324.1 | 262.1 | 31 |
| Forchlorfenuron | 6.05 | Positive | 248 | 93.1 | 49 |
| Forchlorfenuron | 6.05 | Positive | 248 | 129.1 | 25 |
| Formetanate HCl | 1.05 | Positive | 222.1 | 120 | 37 |
| Formetanate HCl | 1.05 | Positive | 222.1 | 165 | 23 |
| Fuberidazole | 2.5 | Positive | 185 | 65 | 57 |
| Fuberidazole | 2.5 | Positive | 185 | 157 | 37 |
| Furalaxyl | 6.76 | Positive | 302.1 | 95 | 39 |
| Furalaxyl | 6.76 | Positive | 302.1 | 242.1 | 21 |
| Furathiocarb | 8.43 | Positive | 383.1 | 195.1 | 25 |
| Furathiocarb | 8.43 | Positive | 383.1 | 252.1 | 17 |
| Hexaconazole | 7.44 | Positive | 314.1 | 70 | 55 |
| Hexaconazole | 7.44 | Positive | 314.1 | 159 | 33 |
| Hexafluomuron | 7.49 | Positive | 461.1 | 141.1 | 57 |
| Hexafluomuron | 7.49 | Positive | 461.1 | 158.2 | 23 |
| Hexythiazox | 8.67 | Positive | 353.1 | 168 | 37 |
| Hexythiazox | 8.67 | Positive | 353.1 | 228 | 19 |
| Imidacloprid | 4.09 | Positive | 256 | 175.1 | 25 |
| Imidacloprid | 4.09 | Positive | 256 | 209.1 | 21 |
| Indoxacarb | 8.15 | Positive | 528 | 203 | 47 |
| Indoxacarb | 8.15 | Positive | 528 | 218 | 35 |
| Ipconazole | 7.68 | Positive | 334.2 | 70 | 37 |
| Ipconazole | 7.68 | Positive | 334.2 | 125 | 47 |
| Iprovalicarb | 6.83 | Positive | 321.2 | 119 | 47 |
| Iprovalicarb | 6.83 | Positive | 321.2 | 203.1 | 13 |
| Isoprocarb | 6.2 | Positive | 194.1 | 95 | 21 |
| Isoprocarb | 6.2 | Positive | 194.1 | 137 | 13 |
| Isoproturon | 6.14 | Positive | 207.2 | 46.1 | 35 |
| Isoproturon | 6.14 | Positive | 207.2 | 72.1 | 29 |

Continued on next page

Table 2E. List of pesticides with MRM transitions used (continued from previous page)

| Compound | RT (min) | Polarity | Precursor (m/z) | Product (m/z) | CE (V) |
|-----------------------|-----------------|-----------------|------------------------|----------------------|---------------|
| Ivermectin | 9.78 | Positive | 892.6 | 307.3 | 31 |
| Ivermectin | 9.78 | Positive | 892.6 | 569.5 | 19 |
| Linuron | 6.85 | Positive | 249.1 | 160 | 25 |
| Linuron | 6.85 | Positive | 249.1 | 182.1 | 21 |
| (Monceren) Pencycuron | 7.95 | Positive | 329.1 | 125 | 31 |
| (Monceren) Pencycuron | 7.95 | Positive | 329.1 | 218.1 | 23 |
| Mandipropamid | 7.08 | Positive | 412.1 | 328.1 | 19 |
| Mandipropamid | 7.08 | Positive | 412.1 | 356.1 | 15 |
| Mefenacet | 7.06 | Positive | 299 | 120.1 | 35 |
| Mefenacet | 7.06 | Positive | 299 | 148.1 | 21 |
| Mepanipyrim | 7.36 | Positive | 224 | 77 | 55 |
| Mepanipyrim | 7.36 | Positive | 224 | 106 | 35 |
| Mepronil | 7.26 | Positive | 270.1 | 119.1 | 31 |
| Mepronil | 7.26 | Positive | 270.1 | 228 | 21 |
| Metalaxyl | 6.02 | Positive | 280.1 | 192.2 | 25 |
| Metalaxyl | 6.02 | Positive | 280.1 | 220.2 | 19 |
| Metconazole | 7.51 | Positive | 320.1 | 70 | 43 |
| Metconazole | 7.51 | Positive | 320.1 | 125 | 53 |
| Methabenzthiazuron | 5.98 | Positive | 222.1 | 150.3 | 45 |
| Methabenzthiazuron | 5.98 | Positive | 222.1 | 165.2 | 21 |
| Methamidophos | 0.63 | Positive | 142 | 94 | 19 |
| Methamidophos | 0.63 | Positive | 142 | 125 | 19 |
| Methiocarb | 6.71 | Positive | 226.1 | 121.1 | 27 |
| Methiocarb | 6.71 | Positive | 226.1 | 169.1 | 13 |
| Methoprotryne | 5.72 | Positive | 272.2 | 198 | 31 |
| Methoprotryne | 5.72 | Positive | 272.2 | 240.2 | 27 |
| Methoxyfenozide | 7.25 | Positive | 369.1 | 149.1 | 21 |
| Methoxyfenozide | 7.25 | Positive | 369.1 | 313.2 | 11 |
| Metobromuron | 6.28 | Positive | 259 | 148.2 | 21 |
| Metobromuron | 6.28 | Positive | 259 | 170.2 | 25 |
| Metribuzin | 5.28 | Positive | 215.1 | 84.1 | 31 |
| Metribuzin | 5.28 | Positive | 215.1 | 187.1 | 25 |
| Mevinphos Isomer | 3.91 | Positive | 225.101 | 127.1 | 21 |
| Mevinphos Isomer | 3.91 | Positive | 225.101 | 193.2 | 11 |
| Mevinphos Isomer_1 | 4.55 | Positive | 225.102 | 127.1 | 21 |
| Mevinphos Isomer_1 | 4.55 | Positive | 225.102 | 193.2 | 11 |
| Mexacarbate | 1.88 | Positive | 223.2 | 151 | 31 |
| Mexacarbate | 1.88 | Positive | 223.2 | 159.1 | 21 |
| Monocrotophos | 1.94 | Positive | 224.1 | 98 | 17 |
| Monocrotophos | 1.94 | Positive | 224.1 | 127.1 | 21 |
| Monolinuron | 6.07 | Positive | 215.1 | 99 | 47 |

Continued on next page

Table 2F. List of pesticides with MRM transitions used (continued from previous page)

| Compound | RT (min) | Polarity | Precursor (m/z) | Product (m/z) | CE (V) |
|--------------------|----------|----------|-----------------|---------------|--------|
| Monolinuron | 6.07 | Positive | 215.1 | 126.1 | 23 |
| Moxidectin | 9.74 | Positive | 640.4 | 498.5 | 17 |
| Moxidectin | 9.74 | Positive | 640.4 | 528.5 | 13 |
| Myclobutanil | 6.99 | Positive | 289 | 70 | 41 |
| Myclobutanil | 6.99 | Positive | 289 | 125 | 39 |
| Nitenpyram | 2.22 | Positive | 271 | 126 | 35 |
| Nitenpyram | 2.22 | Positive | 271 | 225.2 | 17 |
| Nuarimol | 6.72 | Positive | 315 | 81 | 49 |
| Nuarimol | 6.72 | Positive | 315 | 251.1 | 37 |
| Omethoate | 0.85 | Positive | 214 | 124.9 | 31 |
| Omethoate | 0.85 | Positive | 214 | 182.8 | 17 |
| Oxadixyl | 5.26 | Positive | 279.1 | 132.1 | 43 |
| Oxadixyl | 5.26 | Positive | 279.1 | 219.1 | 15 |
| Paclobutrazol | 6.65 | Positive | 294 | 70 | 49 |
| Paclobutrazol | 6.65 | Positive | 294 | 125 | 41 |
| Penconazole | 7.57 | Positive | 284.1 | 70 | 37 |
| Penconazole | 7.57 | Positive | 284.1 | 159 | 35 |
| Phenmedipham | 6.72 | Positive | 301.2 | 107.9 | 44 |
| Phenmedipham | 6.72 | Positive | 301.2 | 168 | 12 |
| Picoxystrobin | 7.62 | Positive | 368 | 145 | 29 |
| Picoxystrobin | 7.62 | Positive | 368 | 205 | 13 |
| Piperonyl butoxide | 8.37 | Positive | 356.2 | 119.1 | 47 |
| Piperonyl butoxide | 8.37 | Positive | 356.2 | 177.2 | 13 |
| Pirimicarb | 3.89 | Positive | 239.2 | 72.1 | 33 |
| Pirimicarb | 3.89 | Positive | 239.2 | 182.1 | 21 |
| Prochloraz | 7.56 | Positive | 376 | 70 | 43 |
| Prochloraz | 7.56 | Positive | 376 | 308 | 15 |
| Promecarb | 6.85 | Positive | 208.1 | 109 | 21 |
| Promecarb | 6.85 | Positive | 208.1 | 151 | 13 |
| Prometon | 5.32 | Positive | 226.1 | 86 | 39 |
| Prometon | 5.32 | Positive | 226.1 | 142 | 33 |
| Prometryne | 6.41 | Positive | 242.2 | 158.1 | 33 |
| Prometryne | 6.41 | Positive | 242.2 | 200.1 | 25 |
| Propamocarb | 1.13 | Positive | 189.2 | 102 | 25 |
| Propamocarb | 1.13 | Positive | 189.2 | 144 | 19 |
| Propargite | 8.91 | Positive | 368.2 | 175.1 | 23 |
| Propargite | 8.91 | Positive | 368.2 | 231.1 | 15 |
| Propham | 6.08 | Positive | 180.1 | 120 | 23 |
| Propham | 6.08 | Positive | 180.1 | 138 | 11 |
| Propiconazole | 7.65 | Positive | 342.1 | 69 | 39 |
| Propiconazole | 7.65 | Positive | 342.1 | 159 | 31 |
| Propoxur | 5.54 | Positive | 210.1 | 111 | 19 |

Continued on next page

Table 2G. List of pesticides with MRM transitions used (continued from previous page)

| Compound | RT (min) | Polarity | Precursor (m/z) | Product (m/z) | CE (V) |
|-----------------------|----------|----------|-----------------|---------------|--------|
| Propoxur | 5.54 | Positive | 210.1 | 168.1 | 11 |
| Pyracarbolid | 5.78 | Positive | 218.1 | 97 | 37 |
| Pyracarbolid | 5.78 | Positive | 218.1 | 125 | 25 |
| Pyraclostrobin | 7.83 | Positive | 388 | 163 | 31 |
| Pyraclostrobin | 7.83 | Positive | 388 | 194 | 17 |
| Pyridaben | 9.23 | Positive | 365 | 147 | 33 |
| Pyridaben | 9.23 | Positive | 365 | 309 | 19 |
| Pyrimethanil | 6.24 | Positive | 200 | 82 | 37 |
| Pyrimethanil | 6.24 | Positive | 200 | 107 | 33 |
| Pyriproxyfen | 8.57 | Positive | 322 | 96 | 21 |
| Pyriproxyfen | 8.57 | Positive | 322 | 185 | 31 |
| Quinoxifen | 8.91 | Positive | 308.1 | 162.1 | 63 |
| Quinoxifen | 8.91 | Positive | 308.1 | 197.1 | 45 |
| Rotenone | 7.42 | Positive | 395.1 | 192.1 | 33 |
| Rotenone | 7.42 | Positive | 395.1 | 213.1 | 31 |
| Secbumeton | 5.38 | Positive | 226.2 | 100 | 37 |
| Secbumeton | 5.38 | Positive | 226.2 | 170.1 | 25 |
| Siduron | 6.61 | Positive | 233.3 | 94 | 31 |
| Siduron | 6.61 | Positive | 233.3 | 137.2 | 23 |
| Simetryn | 5.04 | Positive | 214 | 124 | 29 |
| Simetryn | 5.04 | Positive | 214 | 144 | 29 |
| Spinetoram | 8.44 | Positive | 748.5 | 98.1 | 65 |
| Spinetoram | 8.44 | Positive | 748.5 | 142.2 | 43 |
| Spinosad (Spinosyn A) | 7.85 | Positive | 732.5 | 98.1 | 65 |
| Spinosad (Spinosyn A) | 7.85 | Positive | 732.5 | 142.2 | 39 |
| Spinosad (Spinosyn D) | 8.24 | Positive | 746.8 | 98.3 | 65 |
| Spinosad (Spinosyn D) | 8.24 | Positive | 746.8 | 142.4 | 41 |
| Spirodiclofen | 9.91 | Positive | 411.3 | 71.3 | 31 |
| Spirodiclofen | 9.91 | Positive | 411.3 | 313.3 | 17 |
| Spiromesifen | 9.21 | Positive | 371.2 | 255.2 | 31 |
| Spiromesifen | 9.21 | Positive | 371.2 | 273.2 | 11 |
| Spirotetramat | 6.8 | Positive | 374.2 | 302.2 | 27 |
| Spirotetramat | 6.8 | Positive | 374.2 | 330.2 | 23 |
| Spiroxamine | 7.3 | Positive | 298.2 | 100.1 | 43 |
| Spiroxamine | 7.3 | Positive | 298.2 | 144.2 | 29 |
| Sulfentrazone | 5.93 | Positive | 387 | 146 | 59 |
| Sulfentrazone | 5.93 | Positive | 387 | 307.1 | 29 |
| Tebufenozide | 7.59 | Positive | 353.2 | 133 | 23 |
| Tebufenozide | 7.59 | Positive | 353.2 | 297.2 | 11 |
| Tebufenpyrad | 8.28 | Positive | 334 | 117 | 47 |
| Tebufenpyrad | 8.28 | Positive | 334 | 145 | 37 |
| Tebuthiuron | 5.05 | Positive | 229.1 | 116.1 | 37 |

Continued on next page

Table 2H. List of pesticides with MRM transitions used (continued from previous page)

| Compound | RT (min) | Polarity | Precursor (m/z) | Product (m/z) | CE (V) |
|--------------------|----------|----------|-----------------|---------------|--------|
| Tebuthiuron | 5.05 | Positive | 229.1 | 172.4 | 25 |
| Terbumeton | 5.38 | Positive | 226.1 | 100 | 41 |
| Terbumeton | 5.38 | Positive | 226.1 | 170.1 | 23 |
| Terbutryn | 6.51 | Positive | 242.1 | 68.1 | 61 |
| Terbutryn | 6.51 | Positive | 242.1 | 186.1 | 25 |
| Tetraconazole | 7.18 | Positive | 372.1 | 70 | 47 |
| Tetraconazole | 7.18 | Positive | 372.1 | 159 | 35 |
| Thiabendazole | 2.41 | Positive | 202.1 | 131.2 | 45 |
| Thiabendazole | 2.41 | Positive | 202.1 | 175.1 | 35 |
| Thiacloprid | 5.03 | Positive | 253 | 99 | 59 |
| Thiacloprid | 5.03 | Positive | 253 | 126 | 29 |
| Thiamethoxam | 2.86 | Positive | 292 | 181 | 31 |
| Thiamethoxam | 2.86 | Positive | 292 | 211 | 17 |
| Thidiazuron | 5.28 | Positive | 221.1 | 102.1 | 23 |
| Thidiazuron | 5.28 | Positive | 221.1 | 127.9 | 23 |
| Thiophanate-methyl | 5.5 | Positive | 343 | 151.1 | 31 |
| Thiophanate-methyl | 5.5 | Positive | 343 | 311 | 17 |
| Triadimefon | 7.06 | Positive | 294 | 197.1 | 21 |
| Triadimefon | 7.06 | Positive | 294 | 225 | 19 |
| Triadimenol | 6.65 | Positive | 296.1 | 70 | 33 |
| Triadimenol | 6.65 | Positive | 296.1 | 227.1 | 17 |
| Trichlorfon | 5.03 | Positive | 256.9 | 109.1 | 25 |
| Trichlorfon | 5.03 | Positive | 256.9 | 127 | 23 |
| Tricyclazole | 5 | Positive | 190 | 136 | 39 |
| Tricyclazole | 5 | Positive | 190 | 163 | 33 |
| Trifloxystrobin | 8.2 | Positive | 409 | 186 | 21 |
| Trifloxystrobin | 8.2 | Positive | 409 | 206 | 19 |
| Triflumizole | 7.85 | Positive | 346.1 | 73 | 27 |
| Triflumizole | 7.85 | Positive | 346.1 | 278.1 | 17 |
| Vamidothion | 3.86 | Positive | 288 | 118 | 37 |
| Vamidothion | 3.86 | Positive | 288 | 146 | 17 |

Table 3A. Method validation data (Linearity, LOQ, % recovery, and precision)

| Sr. No. | Compound | R ² | LOQ (mg/kg) | 0.010 mg/kg | | 0.050 mg/kg | |
|---------|------------------------|----------------|-------------|-------------|-------|-------------|-------|
| | | | | % Rec | %RSD | % Rec | %RSD |
| 1 | 3-Hydroxycarbofuran | 0.9992 | 0.01 | 88 | 4.79 | 89 | 3.32 |
| 2 | Acephate | 0.9994 | 0.01 | 101 | 11.66 | 93 | 7.89 |
| 3 | Acetamiprid | 0.9993 | 0.01 | 95 | 1.55 | 90 | 2.15 |
| 4 | Aldicarb sulfone | 0.9991 | 0.01 | 100 | 8.59 | 93 | 6.41 |
| 5 | Aldicarb sulfoxide | 0.9933 | 0.05 | ND | ND | 101 | 3.76 |
| 6 | Ametryn | 0.9984 | 0.01 | 103 | 5.59 | 101 | 7.63 |
| 7 | Aminocarb | 0.9993 | 0.01 | 119 | 3.28 | 118 | 3.4 |
| 8 | Amitraz | 0.9994 | 0.01 | 108 | 10.71 | 102 | 15.01 |
| 9 | Avermectin B1a | 0.9947 | 0.01 | 89 | 17.09 | 112 | 11.89 |
| 10 | Azoxystrobin | 0.9991 | 0.01 | 71 | 19.1 | 76 | 7.71 |
| 11 | Benalaxyl | 0.9992 | 0.01 | 92 | 19.04 | 99 | 7.54 |
| 12 | Bendiocarb | 0.9991 | 0.01 | 96 | 14.82 | 101 | 5.34 |
| 13 | Benzoximate | 0.9978 | 0.01 | 77 | 13.86 | 110 | 19.14 |
| 14 | Bifenazate | 0.9986 | 0.01 | 91 | 17.7 | 107 | 5.12 |
| 15 | Bitertanol | 0.9991 | 0.01 | 74 | 16.95 | 84 | 3.71 |
| 16 | Boscalid | 0.9925 | 0.05 | ND | ND | 100 | 8.34 |
| 17 | Bromucanazole Isomer 1 | 0.9973 | 0.01 | 74 | 11.04 | 90 | 11.32 |
| 18 | Bromucanazole Isomer 2 | 0.9946 | 0.01 | 87 | 9.06 | 86 | 18.16 |
| 19 | Bupirimate | 0.9996 | 0.01 | 82 | 8.52 | 95 | 3.47 |
| 20 | Buprofezin | 0.9933 | 0.01 | 73 | 17.16 | 119 | 11.56 |
| 21 | Butafenacil | 0.9941 | 0.01 | 89 | 15.07 | 105 | 15.96 |
| 22 | Butoxycarboxim | 0.9997 | 0.01 | 92 | 1.16 | 94 | 1.59 |
| 23 | Carbendazim | 0.9968 | 0.01 | 164 | 2.77 | 244 | 1.4 |
| 24 | Carbetamide | 0.9992 | 0.01 | 87 | 8.51 | 109 | 6.07 |
| 25 | Carbofuran | 0.9989 | 0.01 | 83 | 4.49 | 83 | 6.53 |
| 26 | Carboxin | 0.9962 | 0.01 | 94 | 8.94 | 85 | 3.51 |
| 27 | Chlorantraniliprole | 0.9967 | 0.01 | 118 | 14.16 | 90 | 9.13 |
| 28 | Chlorotoluron | 0.9918 | 0.01 | 93 | 8.24 | 85 | 7.27 |
| 29 | Chloroxuron | 0.9939 | 0.01 | 91 | 15.79 | 97 | 9.22 |
| 30 | Clethodim | 0.9942 | 0.01 | 84 | 7.72 | 81 | 12.66 |
| 31 | Clothianidin | 0.9989 | 0.01 | 78 | 17.93 | 92 | 5.82 |
| 32 | Cyazofamid | 0.9953 | 0.01 | 82 | 16.75 | 92 | 14.09 |
| 33 | Cycluron | 0.9986 | 0.01 | 80 | 5.29 | 85 | 3.49 |
| 34 | Cyproconazole Isomer 1 | 0.999 | 0.01 | 90 | 13.56 | 95 | 5.41 |
| 35 | Cyproconazole Isomer 2 | 0.999 | 0.01 | 88 | 6.94 | 95 | 5.29 |
| 36 | Cyprodinil | 0.9977 | 0.01 | 84 | 12.94 | 81 | 9.31 |
| 37 | Desmedipham | 0.9943 | 0.01 | 78 | 6.67 | 79 | 17.01 |
| 38 | Diclobutrazol | 0.9942 | 0.01 | 85 | 11.09 | 79 | 13.62 |
| 39 | Diclotophos | 0.9971 | 0.01 | 100 | 18.15 | 95 | 16.84 |
| 40 | Diethofencarb | 0.9997 | 0.05 | ND | ND | 95 | 19.61 |

Continued on next page. Those analytes without MRL values were considered as 0.01 mg/kg (default MRL set at LOQ)

Table 3B. Method validation data (Linearity, LOQ, % recovery, and precision; continued from previous page)

| Sr. No. | Compound | R ² | LOQ (mg/kg) | 0.010 mg/kg | | 0.050 mg/kg | |
|---------|------------------------|----------------|-------------|-------------|-------|-------------|-------|
| | | | | % Rec | %RSD | % Rec | %RSD |
| 41 | Difenoconazole | 0.9981 | 0.01 | 117 | 10.97 | 101 | 7.3 |
| 42 | Dimethoate | 0.9991 | 0.01 | 87 | 3.12 | 84 | 1.34 |
| 43 | Dimethomorph Isomer 1 | 0.9975 | 0.01 | 77 | 11.99 | 82 | 15.07 |
| 44 | Dimethomorph Isomer 2 | 0.9986 | 0.01 | 80 | 12.26 | 84 | 12.55 |
| 45 | Dimoxystrobin | 0.9958 | 0.01 | 103 | 15.13 | 113 | 10.33 |
| 46 | Diniconazole | 0.9917 | 0.01 | 110 | 14.11 | 85 | 11.01 |
| 47 | Dinotefuran | 0.9981 | 0.01 | 99 | 19.28 | 91 | 4.24 |
| 48 | Dioxacarb | 0.9982 | 0.01 | 91 | 5.73 | 83 | 5.43 |
| 49 | Diuron | 0.9979 | 0.01 | 99 | 14.7 | 85 | 8.01 |
| 50 | Doramectin | 0.9975 | 0.01 | 117 | 8.69 | 117 | 7.9 |
| 51 | Emamectin-benzoate b1a | 0.9942 | 0.01 | 86 | 9.43 | 79 | 11.45 |
| 52 | Epoxiconazole | 0.9956 | 0.01 | 109 | 17.8 | 102 | 8.28 |
| 53 | Eprinomectin | 0.9923 | 0.01 | 120 | 14.45 | 117 | 7.75 |
| 54 | Etaconazole | 0.9973 | 0.01 | 95 | 9.16 | 108 | 4.98 |
| 55 | Ethirimol | 0.9987 | 0.01 | 77 | 10.19 | 87 | 4.37 |
| 56 | Etoxazole | 0.9977 | 0.01 | 94 | 10.05 | 97 | 10.01 |
| 57 | Fenamidone | 0.9936 | 0.01 | 97 | 13.4 | 94 | 19.98 |
| 58 | Fenarimol | 0.9948 | 0.01 | 70 | 11.45 | 106 | 17.26 |
| 59 | Fenzaquin | 0.9967 | 0.01 | 75 | 7.99 | 85 | 12 |
| 60 | Fenbuconazole | 0.9941 | 0.01 | 101 | 15.97 | 100 | 14.81 |
| 61 | Fenpropimorph | 0.9913 | 0.01 | 95 | 16.68 | 71 | 4.21 |
| 62 | Fenpyroximate | 0.9982 | 0.01 | 75 | 6.53 | 89 | 9.19 |
| 63 | Fenuron | 0.9987 | 0.01 | 86 | 6.92 | 88 | 2.66 |
| 64 | Fipronil | 0.9949 | 0.0025 | 89 | 17.37 | 105 | 17.8 |
| 65 | Fluazinam† | 0.9973 | 0.01 | 115 | 14.08 | 116 | 12.36 |
| 66 | Flubendiamide | 0.9959 | 0.01 | 112 | 18.06 | 120 | 3.34 |
| 67 | Flufenacet | 0.9999 | 0.01 | 115 | 4.67 | 119 | 2.93 |
| 68 | Fluometuron | 0.9942 | 0.01 | 94 | 8.56 | 79 | 6.77 |
| 69 | Fluoxastrobin | 0.9999 | 0.01 | 96 | 13.56 | 99 | 14.47 |
| 70 | Flusilazole | 0.9975 | 0.01 | 114 | 14.2 | 102 | 19.46 |
| 71 | Flutolanil | 0.9944 | 0.01 | 77 | 19.85 | 114 | 14.81 |
| 72 | Forchlorfenuron | 0.9975 | 0.01 | 90 | 11.16 | 71 | 10.03 |
| 73 | Formetanate HCl | 0.9985 | 0.01 | 106 | 4.94 | 110 | 2.49 |
| 74 | Fuberidazole | 0.999 | 0.01 | 91 | 1.62 | 91 | 1.64 |
| 75 | Furalaxyl | 0.9992 | 0.01 | 97 | 17.52 | 119 | 4.19 |
| 76 | Furathiocarb | 0.9954 | 0.01 | 79 | 6.2 | 88 | 10.4 |
| 77 | Hexafluomuron | 0.9938 | 0.05 | ND | ND | 108 | 12.19 |
| 78 | Hexaconazole | 0.9947 | 0.01 | 91 | 12.82 | 74 | 12.29 |
| 79 | Hexythiazox | 0.9947 | 0.01 | 83 | 11.82 | 83.71 | 19.71 |
| 80 | Imidacloprid | 0.9993 | 0.01 | 114 | 6.07 | 110 | 4.92 |

Continued on next page. Those analytes without MRL values were considered as 0.01 mg/kg (default MRL set at LOQ)

Table 3C. Method validation data (Linearity, LOQ, % recovery, and precision; continued from previous page)

| Sr. No. | Compound | R ² | LOQ (mg/kg) | 0.010 mg/kg | | 0.050 mg/kg | |
|---------|-----------------------|----------------|-------------|-------------|-------|-------------|-------|
| | | | | % Rec | %RSD | % Rec | %RSD |
| 81 | Indoxacarb | 0.9969 | 0.01 | 116 | 14.52 | 105 | 17.95 |
| 82 | Ipconazole | 0.9955 | 0.01 | 98 | 11.5 | 114 | 16.95 |
| 83 | lprovalicarb | 0.995 | 0.01 | 111 | 12.8 | 119 | 6.99 |
| 84 | Isoprocarb | 0.9928 | 0.01 | 89 | 18.67 | 120 | 12.11 |
| 85 | Isoproturon | 0.9974 | 0.01 | 101 | 5.71 | 93 | 9.63 |
| 86 | Ivermectin | 0.9985 | 0.01 | 119 | 8.66 | 110 | 5.24 |
| 87 | Linuron | 0.9925 | 0.01 | 83 | 12.62 | 79 | 7.65 |
| 88 | (Monceren) Pencycuron | 0.9999 | 0.01 | 113 | 10.71 | 110 | 1.37 |
| 89 | Mandipropamid | 0.9975 | 0.01 | 103 | 18.35 | 120 | 19.38 |
| 90 | Mefenacet | 0.9927 | 0.01 | 117 | 14.89 | 118 | 1.63 |
| 91 | Mepanipyrim | 0.9936 | 0.01 | 86 | 16.23 | 82 | 16.03 |
| 92 | Mepronil | 0.9966 | 0.01 | 91 | 13.52 | 81 | 11.19 |
| 93 | Metalaxyl | 0.9968 | 0.01 | 106 | 5.86 | 120 | 11.61 |
| 94 | Metconazole | 0.996 | 0.01 | 93 | 10.18 | 99 | 11.55 |
| 95 | Methabenzthiazuron | 0.9971 | 0.01 | 97 | 4.09 | 104 | 4.88 |
| 96 | Methamidophos | 0.9944 | 0.01 | 97 | 5.29 | 90 | 3.93 |
| 97 | Methiocarb | 0.9934 | 0.01 | 117 | 7.97 | 79 | 5.65 |
| 98 | Methoprotryne | 0.9992 | 0.01 | 93 | 8.69 | 118 | 2.67 |
| 99 | Methoxyfenozide | 0.9905 | 0.01 | 110 | 14.66 | 117 | 16.96 |
| 100 | Metobromuron | 0.9936 | 0.01 | 101 | 10.46 | 116 | 8.2 |
| 101 | Metribuzin | 0.9951 | 0.01 | 100 | 5.99 | 83 | 7.54 |
| 102 | Mevinphos Isomer 1 | 0.9982 | 0.01 | 90 | 7.47 | 87 | 4.37 |
| 103 | Mevinphos Isomer 2 | 0.9988 | 0.01 | 85 | 8.44 | 88 | 6.16 |
| 104 | Mexacarbate | 0.9997 | 0.01 | 91 | 1.61 | 93 | 1.65 |
| 105 | Monocrotophos | 0.9985 | 0.01 | 118 | 9.42 | 118 | 2.11 |
| 106 | Monolinuron | 0.9979 | 0.01 | 98 | 9.65 | 107 | 4.73 |
| 107 | Moxidectin | 0.9979 | 0.01 | 104 | 6.15 | 113 | 5.8 |
| 108 | Myclobutanil | 0.994 | 0.01 | 84 | 19.98 | 86 | 11.16 |
| 109 | Nitenpyram | 0.9924 | 0.01 | 119 | 7.91 | 120 | 3.44 |
| 110 | Nuarimol | 0.9978 | 0.05 | ND | ND | 19.45 | 94.54 |
| 111 | Omethoate | 0.9951 | 0.01 | 114 | 5.1 | 114 | 3.87 |
| 112 | Oxadixyl | 0.9985 | 0.01 | 86 | 3.68 | 98 | 5.56 |
| 113 | Paclobutrazol | 0.9954 | 0.01 | 96 | 17.77 | 76 | 3.61 |
| 114 | Penconazole | 0.9975 | 0.01 | 83 | 11.76 | 76 | 15.87 |
| 115 | Phenmedipham | 0.9918 | 0.01 | 98 | 16.74 | 106 | 5.52 |
| 116 | Picoxystrobin | 0.9962 | 0.01 | 91 | 15.02 | 88 | 8.86 |
| 117 | Piperonyl butoxide | 0.9951 | 0.01 | 99 | 14.39 | 107 | 5.99 |
| 118 | Pirimicarb | 0.9993 | 0.01 | 89 | 1.79 | 94 | 1.69 |
| 119 | Prochloraz | 0.9969 | 0.01 | 95 | 7.55 | 92 | 5.94 |
| 120 | Promecarb | 0.9938 | 0.01 | 87 | 4.8 | 88 | 3.13 |

Continued on next page. Those analytes without MRL values were considered as 0.01 mg/kg (default MRL set at LOQ)

Table 3D. Method validation data (Linearity, LOQ, % recovery, and precision; continued from previous page)

| Sr. No. | Compound | R ² | LOQ (mg/kg) | 0.010 mg/kg | | 0.050 mg/kg | |
|---------|-----------------------|----------------|-------------|-------------|-------|-------------|-------|
| | | | | % Rec | %RSD | % Rec | %RSD |
| 121 | Prometon | 0.9993 | 0.01 | 79 | 7.66 | 100 | 4.61 |
| 122 | Prometryne | 0.9991 | 0.01 | 94 | 3.97 | 98 | 1.9 |
| 123 | Propamocarb | 0.9986 | 0.01 | 86 | 4.91 | 82 | 3.17 |
| 124 | Propargite | 0.9958 | 0.05 | ND | ND | 93 | 15.98 |
| 125 | Propiconazole | 0.9947 | 0.01 | 110 | 16.5 | 94 | 15.46 |
| 126 | Propoxur | 0.992 | 0.05 | ND | ND | 96 | 7.65 |
| 127 | Pyracarbolid | 0.9957 | 0.01 | 76 | 7 | 81 | 4.02 |
| 128 | Pyraclostrobin | 0.9972 | 0.01 | 107 | 13.18 | 113 | 18.22 |
| 129 | Pyridaben | 0.9978 | 0.01 | 79 | 12.33 | 93 | 7.27 |
| 130 | Pyrimethanil | 0.9985 | 0.01 | 80 | 6.44 | 86 | 5.28 |
| 131 | Pyriproxyfen | 0.9931 | 0.01 | 76 | 11.98 | 84 | 11.86 |
| 132 | Quinoxifen | 0.9973 | 0.01 | 84 | 15.61 | 94 | 6.53 |
| 133 | Rotenone | 0.9978 | 0.01 | 104 | 18.1 | 88 | 18.8 |
| 134 | Seccubumeton | 0.999 | 0.01 | 87 | 5.41 | 90 | 2.16 |
| 135 | Siduron | 0.9938 | 0.01 | 101 | 8.39 | 94 | 10.02 |
| 136 | Simetryn | 0.9996 | 0.01 | 93 | 10.76 | 98 | 5 |
| 137 | Spinetoram | 0.996 | 0.01 | 85 | 9.39 | 90 | 8.9 |
| 138 | Spinosad (Spinosyn A) | 0.9963 | 0.01 | 92 | 11.51 | 102 | 15.45 |
| 139 | Spinosad (Spinosyn D) | 0.996 | 0.01 | 104 | 10.32 | 104 | 5.26 |
| 140 | Spirodiclofen | 0.9926 | 0.01 | 104 | 10.33 | 107 | 7.48 |
| 141 | Spiromesifen | 0.9968 | 0.01 | 104 | 10.33 | 99 | 6.15 |
| 142 | Spirotetramat | 0.9961 | 0.01 | 91 | 10.04 | 102 | 14.87 |
| 143 | Spiroxamine | 0.9983 | 0.01 | 88 | 4.19 | 94 | 1.1 |
| 144 | Tebufenozide | 0.9924 | 0.01 | 89 | 16.12 | 100 | 17.95 |
| 145 | Tebufenpyrad | 0.9929 | 0.01 | 80 | 13.74 | 98 | 10.26 |
| 146 | Tebuthiuron | 0.9968 | 0.01 | 83 | 8.23 | 120 | 3.29 |
| 147 | Terbumeton | 0.9991 | 0.01 | 86 | 4.69 | 93 | 1.65 |
| 148 | Terbutryn | 0.9994 | 0.01 | 87 | 2.87 | 99 | 2.52 |
| 149 | Tetraconazole | 0.9954 | 0.01 | 100 | 9.28 | 93 | 8.94 |
| 150 | Thiabendazole | 0.9996 | 0.01 | 91 | 3.52 | 88 | 0.56 |
| 151 | Thiacloprid | 0.9984 | 0.01 | 76 | 5.05 | 106 | 9.01 |
| 152 | Thiamethoxam | 0.9994 | 0.01 | 94 | 4.3 | 95 | 5.22 |
| 153 | Thidiazuron | 0.9973 | 0.01 | 101 | 10.26 | 74 | 12.76 |
| 154 | Thiophanate-methyl | 0.9972 | 0.01 | 91 | 7.15 | 91 | 6.27 |
| 155 | Triadimefon | 0.9958 | 0.01 | 73 | 14.39 | 90 | 16.51 |
| 156 | Triadimenol | 0.9952 | 0.01 | 101 | 6.91 | 78 | 6.77 |
| 157 | Tricyclazole | 0.9993 | 0.01 | 101 | 15.27 | 113 | 15.56 |
| 158 | Trifloxystrobin | 0.9947 | 0.01 | 77 | 15.52 | 87 | 16.37 |
| 159 | Triflumizole | 0.9942 | 0.01 | 86 | 7.8 | 82 | 7.1 |
| 160 | Vamidotion | 0.9994 | 0.01 | 87 | 3.51 | 86 | 2.68 |

Those analytes without MRL values were considered as 0.01 mg/kg (default MRL set at LOQ)

thermo**scientific**

Find out more at [thermofisher.com](https://www.thermofisher.com)

©2019 Thermo Fisher Scientific Inc. All rights reserved. All 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. **AN72961-EN 0519S**

ThermoFisher
S C I E N T I F I C