

# Analysis of Multiclass Multiresidue Pesticides in Milk Using Agilent Captiva EMR—Lipid with LC/MS/MS and GC/MS/MS

## Authors

Xia Yang, Youjuan Zhang, and  
Zhiming Zhang  
Agilent Technologies, Inc.

## Abstract

This application note describes the analysis of multiclass, multiresidue pesticides in milk. The sample preparation method is based on Agilent QuEChERS extraction followed by Agilent Captiva EMR—Lipid cleanup and analysis. Two instrumental platforms were used: the Agilent 8890 gas chromatograph system coupled with the Agilent 7010B triple quadrupole mass spectrometry (GC/MS/MS) and the Agilent 1290 Infinity II LC system coupled with the Agilent 6470 triple quadrupole mass spectrometer (LC/MS/MS), respectively. The Captiva EMR—Lipid delivers highly efficient removal of matrix interferences in milk and allows detection of most pesticides down to 1 ng/mL. Results demonstrate the benefit of this workflow solution for multiresidue pesticide analysis in milk. For 171 pesticides studied, 98% of recoveries were within 60 to 120%, and over 95% achieve reproducibility of  $\leq 10\%$  RSD.

## Introduction

Milk is an important food in the diet, especially for infants and children. The presence of any contamination in milk is a common food safety concern, and great strides have been taken throughout the dairy industry to ensure the safety of milk. One of the main classes of contaminants in milk is pesticides, which can come from animals ingesting contaminated feed or water.<sup>1</sup> The maximum residue limits for pesticides in milk are often much lower than for general fruits and vegetables.<sup>2</sup> Therefore, the analysis of pesticides in milk requires a sample preparation method for better matrix removal and analytical instrument methods for increased sensitivity. The aim of this study is to develop a simple and efficient workflow for determining a wide range of pesticides that are broadly controlled in milk worldwide.

Captiva EMR–Lipid, which is a lipid removal product, combines size exclusion and hydrophobic interaction to selectively capture lipid hydrocarbon chains without the loss of target analytes. Compared with the traditional QuEChERS dispersive cleanup, Captiva EMR–Lipid cleanup provides simplified pass-through workflow and highly selective and efficient lipid removal. In this application note, 171 pesticides and related metabolites were tested using a simple workflow with QuEChERS extraction followed with Captiva EMR–Lipid cleanup by both LC/MS/MS and GC/MS/MS analysis.

## Experimental

### Chemicals and reagents

All reagents and solvents were HPLC or analytical grade. Acetonitrile (ACN) was from Honeywell (Muskegon, MI, USA). Formic acid (FA) was from J&K Scientific Ltd. (Beijing, China). The pesticides and metabolites standards were purchased from Alta (Tianjin, China).

### Solutions and standards

Individual pesticide stock solutions (100 µg/mL) in ACN were stored at –20 °C, and the mixed spiking solution (1 µg/mL) was prepared in ACN and stored at –20 °C. The 80/20 ACN/H<sub>2</sub>O elution solvent was prepared and stored at room temperature.

### Equipment and consumables

- Agilent Captiva EMR–Lipid, 6 mL cartridges, 600 mg (p/n 5190-1004)
- Agilent Vac Elut 20 manifold with collection rack for 13 × 100 mm test tubes (p/n 12234101)
- Agilent QuEChERS extraction kit EN 15662 method with 50 mL tubes with ceramic homogenizers (p/n 5982-5650CH)
- Agilent Bond Elut EMR–Lipid polish pouch, anhydrous MgSO<sub>4</sub> only (p/n 5982-0102)
- Agilent Bond Elut C18 cartridge, 500 mg, 6 mL, 40 µm (p/n 12102052)
- Agilent QuEChERS Dispersive Kit, fruits and vegetables with fats and waxes, AOAC method, 15 mL (p/n 5982-5158)

- SPEX SamplePrep 2010 Geno/Grinder (Metuchen, NJ, USA)
- Eppendorf Centrifuge 5810R (Hamburg, Germany)
- LC column and GC column, liner, etc.

### Instrument conditions

Instrument detection was carried out using both LC/MS/MS and GC/MS/MS platforms. The LC/MS/MS detection was performed on an Agilent 1290 Infinity II LC System. This consisted of the Agilent 1290 Infinity II high speed pump (G7120A), the Agilent 1290 Infinity II multisampler (G7167B), and the Agilent 1290 Infinity II multicolumn thermostat (G7116B). These were coupled with an 6470A triple quadrupole LC/MS (G6470A) with an Agilent Jet Stream electrospray ion source. The GC/MS/MS detection was on an Agilent 8890 GC with the Agilent 7010 triple quadrupole GC/MS. Agilent MassHunter workstation software was used for data acquisition and analysis.

Table 1 lists the GC/MS/MS method parameters for GC-amenable pesticides; Table 2 lists the LC/MS/MS method conditions for LC-amenable pesticides.

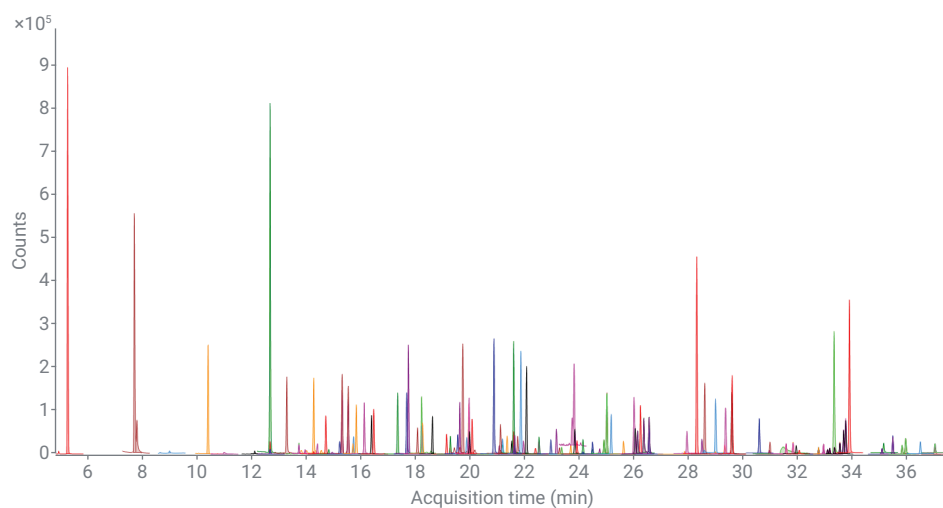
**Table 1.** GC/MS/MS method conditions.

| GC/MS/MS Parameter        | Setting                                                                                  |
|---------------------------|------------------------------------------------------------------------------------------|
| Column                    | Agilent HP-5ms UI, 30 m × 0.25 mm, 0.25 µm (p/n 19091S-433UI)                            |
| Carrier Gas               | Helium, 1.019 mL/min, constant flow                                                      |
| Injection Volume          | 1 µL                                                                                     |
| Inlet Liner               | Agilent Ultra Inert, splitless, single taper, glass wool (p/n 5190-2293)                 |
| Over Program              | 60 °C (1 min), 40 °C/min to 120 °C, then 5 °C/min to 310 °C                              |
| Inlet                     | Split/splitless; temperature: 280 °C<br>Splitless mode, purge flow 30 mL/min at 0.75 min |
| Run Time                  | 40.5 min                                                                                 |
| Transfer Line Temperature | 280 °C                                                                                   |
| Collision Cell EPC        | Quench Gas He, 2.25 mL/min ; Collision Gas N <sub>2</sub> , 1.5 mL/min                   |
| Source Temperature (HES)  | 280 °C                                                                                   |
| Quadrupole Temperature    | 150 °C                                                                                   |
| Data Monitoring           | dMRM                                                                                     |
| Gain Factor               | 10                                                                                       |
| Solvent Delay             | 3 min                                                                                    |

The MRM transitions and settings are listed in the appendix. Figures 1 and 2 show a typical GC/MS/MS MRM chromatogram and LC/MS/MS chromatogram for milk samples fortified with pesticides at the level of 50 ng/mL

**Table 2.** LC/MS/MS method conditions.

| LC/MS/MS Parameter                     | Setting                                                                                                                                                                               |           |                    |               |
|----------------------------------------|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|-----------|--------------------|---------------|
| Column                                 | Agilent ZORBAX Eclipse plus C18, 3.0 × 150 mm, 1.8 μm (p/n 959759-302)                                                                                                                |           |                    |               |
| Column Temperature                     | 45 °C                                                                                                                                                                                 |           |                    |               |
| Autosampler Temperature                | 10 °C                                                                                                                                                                                 |           |                    |               |
| Injection Volume                       | 2 μL                                                                                                                                                                                  |           |                    |               |
| Mobile Phase                           | A) Water, containing 4.5 mM ammonium formate, 0.5 mM ammonium floride, 0.1% formic acid<br>B) Methanol, containing 4.5 mM ammonium formate, 0.5 mM ammonium floride, 0.1% formic acid |           |                    |               |
| Gradient                               | Time (min)                                                                                                                                                                            | %B        | Flow rate (mL/min) |               |
|                                        | 0                                                                                                                                                                                     | 2         | 0.45               |               |
|                                        | 0.5                                                                                                                                                                                   | 2         | 0.45               |               |
|                                        | 1                                                                                                                                                                                     | 50        | 0.45               |               |
|                                        | 4                                                                                                                                                                                     | 65        | 0.45               |               |
|                                        | 16                                                                                                                                                                                    | 100       | 0.45               |               |
|                                        | 18                                                                                                                                                                                    | 100       | 0.45               |               |
|                                        | 18.1                                                                                                                                                                                  | 2         | 0.45               |               |
| 20                                     | 2                                                                                                                                                                                     | 0.45      |                    |               |
| Stop Time                              | 20 min                                                                                                                                                                                |           |                    |               |
| Source Parameters                      |                                                                                                                                                                                       |           |                    |               |
| Gas Temperature                        | 250 °C                                                                                                                                                                                |           |                    |               |
| Gas Flow                               | 11 L/min                                                                                                                                                                              |           |                    |               |
| Nebulizer                              | 40 psi                                                                                                                                                                                |           |                    |               |
| Sheath Gas Temperature                 | 350 °C                                                                                                                                                                                |           |                    |               |
| Sheath Gas Flow                        | 12 L/min                                                                                                                                                                              |           |                    |               |
| Capillary Voltage                      | +3,500                                                                                                                                                                                |           |                    |               |
| Nozzle Voltage                         | +300                                                                                                                                                                                  |           |                    |               |
| Time Segments                          |                                                                                                                                                                                       |           |                    |               |
| Agilent 1290 Infinity II binary system | Start time (min)                                                                                                                                                                      | Scan type | Div valve          | Delta EMV (+) |
|                                        | 0                                                                                                                                                                                     | DMRM      | To Waste           | 0             |
|                                        | 2.1                                                                                                                                                                                   | DMRM      | To MS              | 400           |
|                                        | 18                                                                                                                                                                                    | DMRM      | To Waste           | 0             |



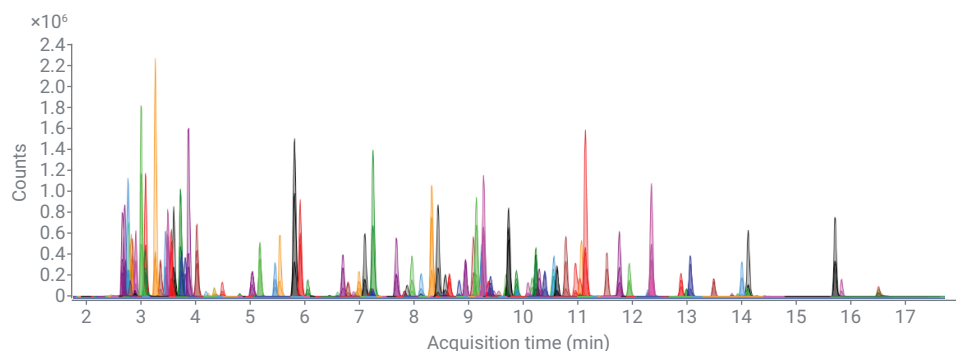
**Figure 1.** GC/MS/MS MRM chromatogram of pesticides in the fortified milk sample at the level of 50 ng/mL.

## Sample preparation

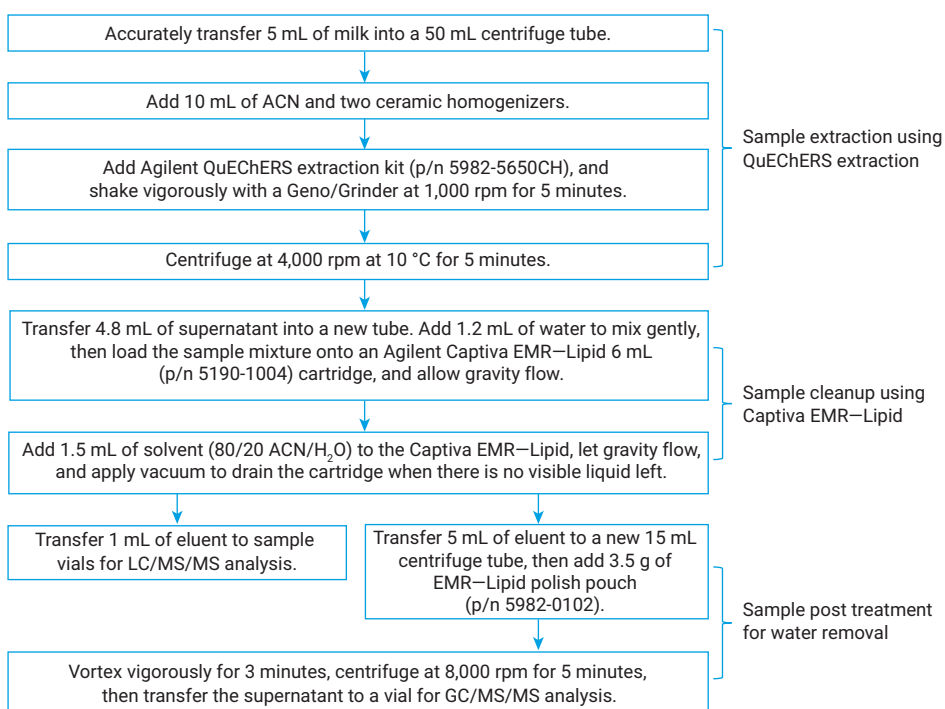
Milk was purchased from a local grocery store and prepared using the developed sample preparation method. Figure 3 shows the step-by-step procedure, featuring three major parts: sample extraction using the QuEChERS extraction kit, sample extract cleanup using Captiva EMR–Lipid, and post treatment for water removal using anhydrous  $\text{MgSO}_4$  salt partition specifically for GC/MS/MS analysis. The entire workflow introduced 3.125-fold and 2.5-fold dilution of the original sample concentration for LC/MS/MS and GC/MS/MS, respectively.

## Calibration standards and quality control (QC) samples

Prespiked QC samples were fortified by spiking the appropriate standard spiking solution into milk and vortexing, then equilibrating for five minutes. Matrix-matched calibration standards were prepared by spiking standard solution into the matrix blank, which was processed using the developed sample preparation workflow. The calibration standards correspond to 1, 2, 5, 10, 20, 50, 100, 200, and 500 ng/mL in milk. Four levels of QC samples were quantified against calibration curves at low level (5 and/or 10 ng/mL), midlevel (50 ng/mL), and high level (100 ng/mL) in milk with four replicates. Analyte identification, confirmation, and quantitation were determined from the retention times and MRM transitions.



**Figure 2.** LC/MS/MS MRM chromatogram of pesticides in the fortified milk sample at the level of 50 ng/mL.



**Figure 3.** The step-by-step procedure for milk sample preparation.

### Sample extraction and matrix removal investigation

The extraction method for a 5 mL milk sample was evaluated. First, a 15 mL ACN solvent extraction was compared to a QuEChERS extraction kit using 10 mL of ACN followed by the QuEChERS EN extraction. The cleanup step was also evaluated. The GC/MS full scan background profiles for sample extract without cleanup were compared to those for sample extract with the following cleanup techniques, including Captiva EMR–Lipid cleanup, Bond Elut C18 SPE cleanup, and Bond Elut QuEChERS dSPE cleanup.

## Results and discussion

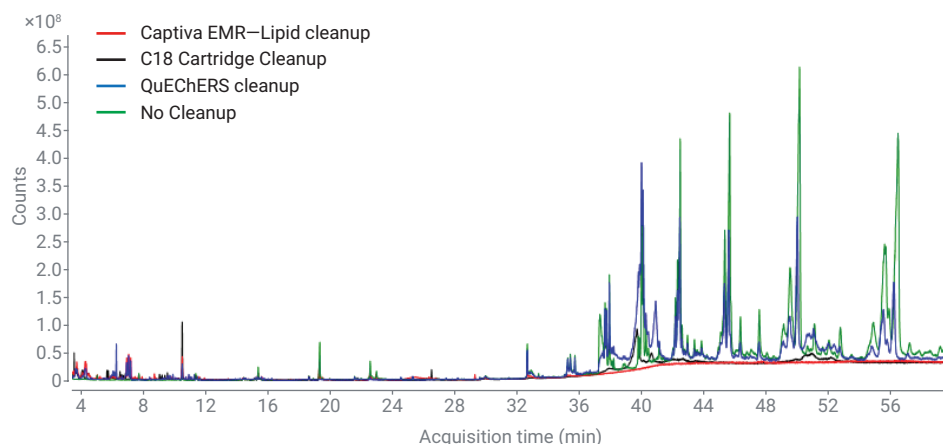
### Sample preparation optimization

Two extraction methods were evaluated: solvent extraction with ACN and QuEChERS extraction using ACN. Nonacidified ACN was used because it was reported that acidified ACN can cause degradation of some pesticides such as amitraz and semiamitraz.<sup>3</sup> The results from QuEChERS extraction showed better recoveries for many hydrophobic pesticides than results from ACN extraction only. The QuEChERS extraction method was selected for the milk sample extraction.

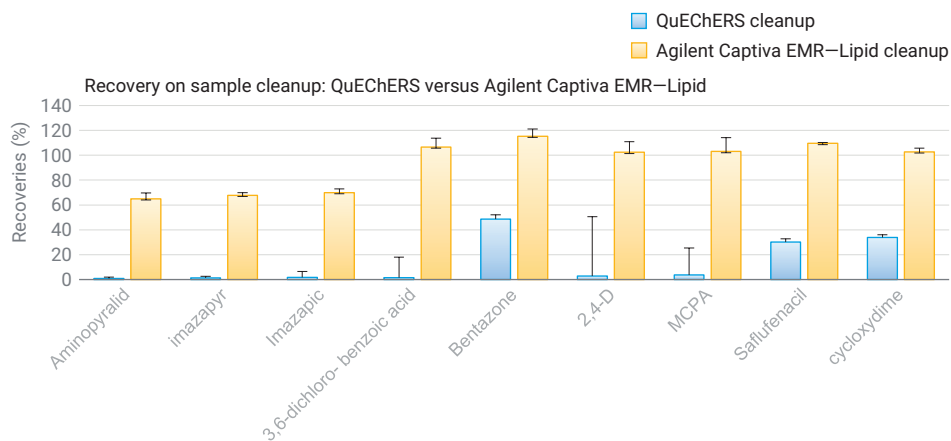
Matrix removal was investigated with a GC/MS full scan background profile comparison for sample extract without cleanup and with various cleanups by Captiva EMR–Lipid, Bond Elut C18 cartridge, and the QuEChERS dSPE kit. The overlapped full scan chromatograms are shown in Figure 4. By comparing the total integration peak area within

the entire retention window, Captiva EMR–Lipid cleanup was demonstrated for 99% of the matrix removal and significantly outperformed the other three cleanup methods. Cleanup with Captiva EMR–Lipid also removes late eluent interferences (as shown after 30 minutes), which are a significant contributor to instrument contamination.

The pesticide recoveries using various cleanup techniques were compared, and the results are shown in Figure 5. Results show that QuEChERS dispersive SPE cleanup caused significant loss of acidic pesticides, resulting from the use of basic sorbent PSA in the dispersive kit.



**Figure 4.** GC/MS full scan for the milk samples with different cleanup method.



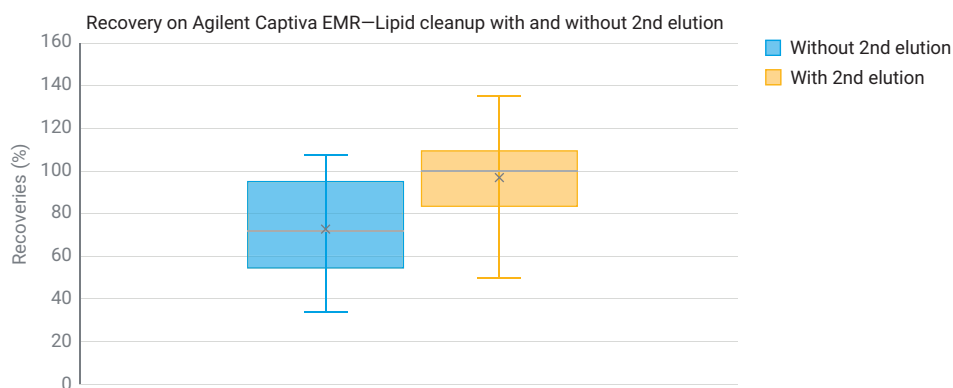
**Figure 5.** Comparison of the recoveries of acidic pesticides with Agilent QuEChERS dSPE cleanup and Agilent Captiva EMR–Lipid cleanup at 40 ng/mL spiking level by LC/MS/MS.

The recoveries for the Captiva EMR–Lipid cartridge cleanup step were subsequently studied for the total amount eluted with the two successive elution steps. From the box-and-whisker plot of the recovery data with and without 2nd elution in Figure 6, the middle line of the box represents the median recovery number, and the bottom and top lines of the box represent the median of the bottom half and top half recovery number, respectively. The whiskers (vertical lines) extend from the ends of the box to the minimum value and maximum value. Overall, the recoveries of the pesticides with 2nd elution were well improved. Therefore, a secondary elution with 1.5 mL of 80/20 ACN/water was applied after the initial 6 mL of sample mixture elution.

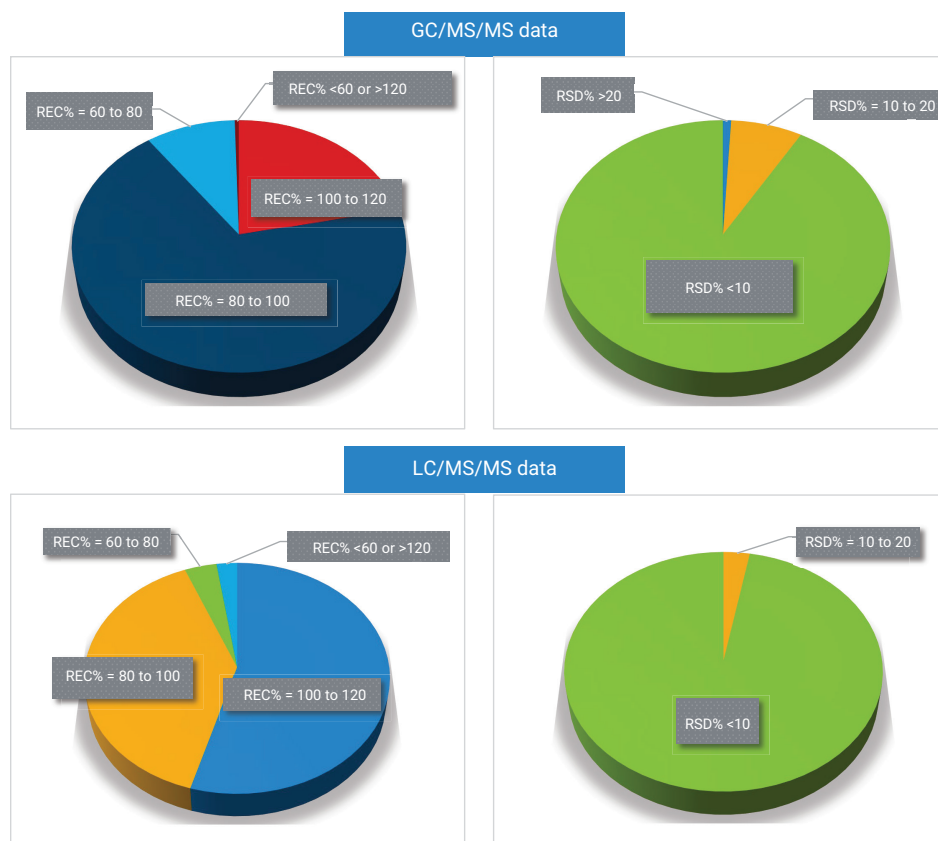
Introduction of water residue can be detrimental to a GC instrument, so the water in the eluent after Captiva EMR–Lipid cleanup must be removed. There are different ways to remove water before GC/MS/MS analysis that have been applied in various Agilent solutions.<sup>4,5</sup> For large panel pesticide analysis with a broad range of different chemical physical feature on analytes, MgSO<sub>4</sub> salting out was usually recommended and proved to be the appropriate method for water removal.

### Method validation

The quantitative method validation of GC/MS/MS and LC/MS/MS includes calibration curve linearity, analyte recovery, and precision at four spiking levels. The detailed quantitation results are shown in Tables 3 and 4, and the statistical summary results are shown in Figure 7. Acceptable recoveries (60% to 120%) were achieved for over 98% of analytes, and RSD values were less than 10% for over 95% of analytes. Of the 171 pesticides, 43 were verified with both GC/MS/MS and LC/MS/MS platform detection.



**Figure 6.** Comparison of with and without 2nd elution for the total pool of pesticides with LC/MS/MS and GC/MS/MS at 40 ppb spiking level in milk.



**Figure 7.** Statistical summary of the analysis of pesticides in milk method validation.

**Table 3.** Method quantitation results for 118 pesticides in milk with GC/MS/MS.

| Pesticide                       | Linearity Range (ng/mL) | R <sup>2</sup> | Low QC 5 ng/mL in Milk (n = 4) |      | Low QC 10 ng/mL in Milk (n = 4) |      | Mid QC 50 ng/mL in Milk (n = 4) |      | High QC 100 ng/mL in Milk (n = 4) |      |
|---------------------------------|-------------------------|----------------|--------------------------------|------|---------------------------------|------|---------------------------------|------|-----------------------------------|------|
|                                 |                         |                | Mean Recovery %                | RSD% | Mean Recovery %                 | RSD% | Mean Recovery %                 | RSD% | Mean Recovery %                   | RSD% |
| 2,4,6-Trichlorophenol           | 1 to 500                | 0.998          | 98.5                           | 5.2  | 101.2                           | 2.9  | 94.8                            | 7.2  | 92.0                              | 5.5  |
| Aldrin                          | 1 to 500                | 0.998          | 80.7                           | 6.2  | 78.4                            | 4.7  | 74.9                            | 5.6  | 71.2                              | 2.5  |
| Dichlorvos*                     | 5 to 500                | 0.998          | 116.6                          | 6.8  | 98.1                            | 9.1  | 96.3                            | 12.4 | 93.1                              | 4.8  |
| Phorate*                        | 2 to 500                | 0.993          | 96.7                           | 9.3  | 108.1                           | 4.2  | 99.8                            | 9.1  | 95.9                              | 2.6  |
| Methacrifos                     | 1 to 500                | 0.998          | 99.7                           | 5.3  | 110.1                           | 5.3  | 102.4                           | 8.1  | 97.8                              | 1.3  |
| Pentachloronitrobenzene         | 1 to 500                | 0.997          | 96.2                           | 7.9  | 94.2                            | 6.3  | 95.6                            | 3.8  | 85.4                              | 5.2  |
| Endosulfan I (alpha isomer)     | 2 to 500                | 0.997          | 85.0                           | 11.3 | 104.8                           | 4.0  | 95.3                            | 6.6  | 91.5                              | 1.8  |
| Acetamiprid*                    | 5 to 500                | 0.995          | 73.3                           | 20.7 | 69.8                            | 5.1  | 78.5                            | 4.1  | 83.6                              | 2.0  |
| Bentazone*                      | 10 to 500               | 0.994          | 98.8                           | 8.1  | 120.4                           | 8.4  | 103.3                           | 5.3  | 104.8                             | 5.6  |
| Chlorantraniliprole*            | 10 to 500               | 0.992          | 80.0                           | 11.9 | 110.6                           | 7.0  | 84.0                            | 6.2  | 94.7                              | 2.3  |
| Cyromazine*                     | 2 to 500                | 0.990          | 64.1                           | 10.3 | 61.4                            | 9.3  | 60.0                            | 5.0  | 55.0                              | 3.9  |
| Famoxadone*                     | 5 to 500                | 0.996          | 104.2                          | 20.2 | 105.8                           | 9.8  | 92.2                            | 2.0  | 98.2                              | 3.4  |
| Methamidophos*                  | 1 to 500                | 0.994          | 89.5                           | 7.8  | 91.2                            | 14.4 | 80.6                            | 13.2 | 81.9                              | 1.1  |
| Spirodiclofen*                  | 10 to 500               | 0.997          | —                              | —    | 83.3                            | 22.1 | 92.8                            | 7.0  | 97.0                              | 1.9  |
| Oxamyl*                         | 5 to 500                | 0.993          | 92.0                           | 12.0 | 106.7                           | 3.3  | 98.5                            | 4.0  | 99.0                              | 3.1  |
| Prochloraz*                     | 2 to 500                | 0.998          | 100.5                          | 14.8 | 80.5                            | 7.2  | 89.7                            | 6.7  | 99.3                              | 5.5  |
| Azinphos-ethyl                  | 1 to 500                | 0.993          | 98.4                           | 8.8  | 96.6                            | 2.1  | 85.4                            | 5.1  | 94.6                              | 0.7  |
| Captan                          | 50 to 500               | 0.999          | —                              | —    | —                               | —    | 70.4                            | 14.4 | 72.2                              | 5.6  |
| Hexachlorobenzene               | 2 to 500                | 1.000          | 93.4                           | 9.5  | 69.9                            | 12.6 | 79.6                            | 6.1  | 64.6                              | 1.2  |
| Fenamidone                      | 1 to 500                | 0.991          | 83.4                           | 10.8 | 86.2                            | 4.2  | 79.9                            | 1.3  | 87.8                              | 3.3  |
| Ethoprophos*                    | 1 to 500                | 0.993          | 101.6                          | 6.1  | 105.4                           | 2.3  | 100.9                           | 11.1 | 96.3                              | 0.5  |
| Azinphos-methyl                 | 5 to 500                | 0.995          | —                              | —    | 95.6                            | 6.5  | 73.4                            | 9.7  | 89.9                              | 2.7  |
| Phosmet                         | 1 to 500                | 0.995          | 96.3                           | 15.1 | 90.0                            | 3.0  | 72.2                            | 7.5  | 87.1                              | 1.9  |
| Azoxystrobin*                   | 5 to 500                | 0.995          | 75.3                           | 16.9 | 92.1                            | 11.8 | 92.0                            | 4.4  | 101.9                             | 1.7  |
| Bifenthrin*                     | 1 to 500                | 0.990          | 73.2                           | 12.6 | 76.8                            | 5.1  | 70.3                            | 2.8  | 71.4                              | 3.5  |
| Bitertanol*                     | 5 to 500                | 0.998          | 89.4                           | 4.6  | 90.6                            | 9.5  | 90.5                            | 8.2  | 94.5                              | 1.2  |
| Boscalid*                       | 1 to 500                | 0.994          | 99.0                           | 7.7  | 96.8                            | 1.4  | 90.0                            | 5.6  | 98.6                              | 1.2  |
| Buprofezin*                     | 10 to 500               | 0.995          | —                              | —    | 90.8                            | 7.2  | 87.1                            | 7.9  | 86.2                              | 3.8  |
| Carbaryl*                       | 2 to 500                | 0.991          | 96.4                           | 12.2 | 105.3                           | 4.9  | 99.0                            | 5.8  | 100.4                             | 1.4  |
| Chinomethionate (Oxythioquinox) | 1 to 500                | 0.995          | 88.8                           | 7.9  | 89.8                            | 1.6  | 88.3                            | 5.6  | 88.8                              | 1.6  |
| Chlordane-cis                   | 1 to 500                | 0.998          | 87.4                           | 7.7  | 90.4                            | 1.8  | 88.6                            | 5.1  | 86.6                              | 1.5  |
| Chlordane-oxy                   | 1 to 500                | 0.998          | 88.9                           | 4.8  | 93.4                            | 3.3  | 93.0                            | 5.2  | 88.1                              | 2.1  |
| Chlordane-trans                 | 1 to 500                | 0.998          | 91.3                           | 2.7  | 89.4                            | 0.4  | 90.8                            | 4.2  | 83.6                              | 0.9  |
| Chlorfenvinphos                 | 1 to 500                | 0.994          | 88.5                           | 14.3 | 91.5                            | 4.4  | 80.7                            | 10.9 | 94.3                              | 2.5  |
| Chlorpropham                    | 1 to 500                | 1.000          | 104.8                          | 3.2  | 105.7                           | 3.5  | 105.9                           | 3.7  | 99.0                              | 3.5  |
| Chlorpyrifos*                   | 1 to 500                | 0.993          | 90.7                           | 9.6  | 98.5                            | 3.8  | 98.1                            | 8.1  | 94.1                              | 1.4  |
| Chlorpyrifos-methyl*            | 1 to 500                | 0.990          | 94.3                           | 5.9  | 97.6                            | 4.2  | 98.4                            | 6.1  | 96.9                              | 2.1  |
| Clofentezine*                   | 1 to 500                | 1.000          | 111.5                          | 1.8  | 106.4                           | 5.1  | 107.7                           | 1.4  | 107.8                             | 1.4  |
| Coumaphos                       | 5 to 500                | 0.992          | 83.6                           | 10.1 | 87.4                            | 4.9  | 85.7                            | 7.4  | 96.6                              | 2.0  |
| Cyfluthrin-1                    | 2 to 500                | 0.998          | 101.9                          | 5.0  | 89.1                            | 6.8  | 83.0                            | 5.2  | 88.8                              | 3.4  |
| Cyfluthrin-2                    | 2 to 500                | 0.996          | 94.4                           | 6.9  | 92.0                            | 4.5  | 88.8                            | 5.7  | 91.9                              | 2.5  |
| Cyfluthrin-3                    | 5 to 500                | 0.998          | 94.5                           | 7.8  | 88.3                            | 6.4  | 85.4                            | 5.0  | 93.8                              | 1.4  |
| Cyfluthrin-4                    | 5 to 500                | 0.996          | 106.4                          | 5.6  | 86.9                            | 9.7  | 89.1                            | 2.9  | 94.4                              | 0.9  |
| Cypermethrin-1                  | 5 to 500                | 0.997          | —                              | —    | 92.3                            | 7.7  | 85.7                            | 6.9  | 91.9                              | 0.5  |

| Pesticide                   | Linearity Range (ng/mL) | R <sup>2</sup> | Low QC 5 ng/mL in Milk (n = 4) |      | Low QC 10 ng/mL in Milk (n = 4) |      | Mid QC 50 ng/mL in Milk (n = 4) |      | High QC 100 ng/mL in Milk (n = 4) |      |
|-----------------------------|-------------------------|----------------|--------------------------------|------|---------------------------------|------|---------------------------------|------|-----------------------------------|------|
|                             |                         |                | Mean Recovery %                | RSD% | Mean Recovery %                 | RSD% | Mean Recovery %                 | RSD% | Mean Recovery %                   | RSD% |
| Cypermethrin-2              | 5 to 500                | 0.997          | —                              | —    | 91.0                            | 5.6  | 91.4                            | 5.9  | 97.8                              | 1.9  |
| Cypermethrin-3              | 2 to 500                | 0.996          | —                              | —    | 97.5                            | 3.5  | 84.0                            | 5.6  | 89.3                              | 2.0  |
| Cypermethrin-4              | 2 to 500                | 0.996          | —                              | —    | 93.2                            | 2.9  | 86.6                            | 5.5  | 91.5                              | 1.4  |
| Cyprodinil*                 | 1 to 500                | 0.993          | 87.8                           | 6.4  | 93.2                            | 2.3  | 89.5                            | 7.2  | 93.7                              | 1.1  |
| DDD-o,p'                    | 1 to 500                | 1.000          | 85.1                           | 7.6  | 83.5                            | 4.2  | 81.9                            | 5.8  | 79.8                              | 1.2  |
| DDD-p,p'                    | 1 to 500                | 1.000          | 81.9                           | 7.9  | 78.9                            | 3.3  | 80.3                            | 7.9  | 77.8                              | 2.2  |
| DDE-p,p'                    | 1 to 500                | 0.999          | 105.9                          | 5.3  | 84.6                            | 2.6  | 80.9                            | 6.6  | 80.6                              | 2.7  |
| DDT-o,p'                    | 1 to 500                | 0.997          | 80.0                           | 7.4  | 76.5                            | 2.9  | 75.3                            | 5.8  | 73.0                              | 1.6  |
| DDT-p,p'                    | 1 to 500                | 0.996          | 81.2                           | 11.3 | 72.0                            | 3.3  | 68.7                            | 5.3  | 67.6                              | 1.7  |
| Deltamethrin                | 2 to 500                | 0.993          | 92.3                           | 8.8  | 95.2                            | 1.1  | 90.5                            | 2.9  | 92.8                              | 2.5  |
| Demeton-S-methyl            | 5 to 500                | 0.991          | 95.0                           | 11.6 | 111.0                           | 7.6  | 90.5                            | 11.0 | 94.4                              | 3.3  |
| Diazinon*                   | 1 to 500                | 0.995          | 101.1                          | 5.8  | 107.0                           | 4.1  | 101.1                           | 7.2  | 98.6                              | 1.4  |
| Dichlofenthion              | 1 to 500                | 0.993          | 99.7                           | 5.8  | 105.3                           | 2.8  | 98.2                            | 7.0  | 95.8                              | 1.5  |
| Dichloran                   | 1 to 500                | 0.995          | 105.9                          | 5.2  | 106.9                           | 4.4  | 101.0                           | 5.9  | 99.3                              | 0.9  |
| Dieldrin                    | 2 to 500                | 0.999          | 92.3                           | 8.5  | 90.2                            | 5.1  | 90.3                            | 5.3  | 87.6                              | 2.1  |
| Difenoconazole I            | 1 to 500                | 0.994          | 100.5                          | 12.7 | 99.4                            | 6.4  | 92.8                            | 4.5  | 97.2                              | 1.2  |
| Difenoconazole II           | 1 to 500                | 0.992          | 98.0                           | 8.2  | 106.6                           | 2.4  | 93.4                            | 3.7  | 98.4                              | 1.2  |
| Dimethipin                  | 1 to 500                | 0.997          | 95.9                           | 4.7  | 105.1                           | 4.1  | 108.3                           | 3.8  | 103.6                             | 0.7  |
| Dimethoate*                 | 2 to 500                | 0.996          | 101.9                          | 7.4  | 104.0                           | 2.5  | 90.1                            | 9.7  | 99.0                              | 2.1  |
| Diphenylamine*              | 1 to 500                | 0.998          | 101.5                          | 3.5  | 103.6                           | 4.4  | 99.9                            | 6.9  | 96.4                              | 1.0  |
| Endosulfan II (beta isomer) | 1 to 500                | 0.997          | 93.0                           | 4.1  | 91.0                            | 3.7  | 92.3                            | 4.5  | 95.2                              | 0.7  |
| Endosulfan sulfate          | 1 to 500                | 0.993          | 88.5                           | 7.1  | 86.5                            | 1.1  | 83.6                            | 5.8  | 90.9                              | 1.6  |
| Endrin                      | 2 to 500                | 0.993          | 89.3                           | 9.2  | 91.3                            | 2.6  | 86.1                            | 8.1  | 88.2                              | 2.0  |
| Ethion                      | 1 to 500                | 0.996          | 90.7                           | 9.1  | 92.5                            | 1.3  | 89.3                            | 6.3  | 94.5                              | 0.9  |
| Ethofenprox                 | 1 to 500                | 0.992          | 92.3                           | 6.7  | 84.6                            | 3.2  | 79.5                            | 7.3  | 81.7                              | 1.1  |
| Fenamiphos sulfone*         | 10 to 500               | 0.992          | 77.9                           | 16.2 | 83.9                            | 2.1  | 74.6                            | 10.1 | 82.0                              | 2.6  |
| Fenitrothion                | 1 to 500                | 0.996          | 101.7                          | 7.6  | 101.9                           | 3.1  | 97.7                            | 7.4  | 102.3                             | 0.5  |
| Fenpropathrin               | 2 to 500                | 0.993          | 82.2                           | 13.9 | 85.4                            | 6.8  | 79.2                            | 2.3  | 87.3                              | 3.9  |
| Fenpropimorph*              | 1 to 500                | 0.994          | 78.6                           | 4.2  | 78.9                            | 3.8  | 79.7                            | 7.3  | 82.3                              | 2.0  |
| Fensulfothion               | 2 to 500                | 0.995          | 97.3                           | 4.9  | 91.8                            | 1.9  | 83.8                            | 5.9  | 97.0                              | 0.7  |
| Fenthion                    | 1 to 500                | 0.992          | 93.9                           | 9.0  | 102.2                           | 2.4  | 97.2                            | 7.3  | 99.8                              | 1.3  |
| Fenvalerate I               | 1 to 500                | 0.994          | 102.5                          | 6.4  | 92.8                            | 3.2  | 87.3                            | 7.0  | 90.0                              | 1.5  |
| Fenvalerate II              | 1 to 500                | 0.994          | 98.8                           | 8.4  | 92.1                            | 2.5  | 89.0                            | 5.2  | 92.5                              | 1.3  |
| Fipronil*                   | 1 to 500                | 0.994          | 103.5                          | 5.6  | 101.8                           | 2.4  | 97.5                            | 4.2  | 101.8                             | 1.5  |
| Fipronil sulfide*           | 1 to 500                | 0.995          | 90.4                           | 5.2  | 96.1                            | 1.7  | 97.1                            | 4.7  | 102.9                             | 1.2  |
| Fipronil sulfone            | 1 to 500                | 0.995          | 99.0                           | 12.4 | 102.7                           | 2.3  | 92.6                            | 6.1  | 100.7                             | 1.3  |
| Flusilazole*                | 1 to 500                | 0.991          | 90.9                           | 9.8  | 97.3                            | 2.1  | 90.4                            | 6.5  | 100.9                             | 1.2  |
| HCH- <i>alpha</i>           | 1 to 500                | 0.999          | 100.2                          | 5.6  | 104.6                           | 2.7  | 104.0                           | 7.8  | 92.4                              | 2.8  |
| HCH- <i>beta</i>            | 1 to 500                | 0.999          | 102.2                          | 6.4  | 104.9                           | 4.2  | 106.1                           | 4.8  | 97.3                              | 1.1  |
| HCH- <i>delta</i>           | 1 to 500                | 0.998          | 101.8                          | 3.4  | 104.5                           | 3.0  | 104.5                           | 3.9  | 98.6                              | 0.5  |
| HCH- <i>gamma</i>           | 1 to 500                | 0.999          | 102.0                          | 3.3  | 101.3                           | 5.3  | 101.9                           | 3.9  | 95.7                              | 1.6  |
| Heptachlor                  | 1 to 500                | 0.996          | 93.7                           | 2.2  | 95.4                            | 5.7  | 94.3                            | 4.6  | 93.6                              | 0.9  |
| Heptachlor exo-epoxide      | 2 to 500                | 0.994          | 92.7                           | 8.3  | 91.8                            | 4.0  | 93.5                            | 7.0  | 88.1                              | 1.9  |
| Isopyrzazam*                | 2 to 500                | 0.991          | 101.3                          | 7.6  | 91.5                            | 4.7  | 84.4                            | 6.5  | 93.3                              | 0.4  |
| Malathion                   | 1 to 500                | 0.995          | 92.1                           | 1.6  | 98.0                            | 2.9  | 94.5                            | 7.2  | 99.2                              | 1.8  |



| Pesticide               | Linearity Range (ng/mL) | R <sup>2</sup> | Low QC 5 ng/mL in Milk (n = 4) |      | Low QC 10 ng/mL in Milk (n = 4) |      | Mid QC 50 ng/mL in Milk (n = 4) |      | High QC 100 ng/mL in Milk (n = 4) |      |
|-------------------------|-------------------------|----------------|--------------------------------|------|---------------------------------|------|---------------------------------|------|-----------------------------------|------|
|                         |                         |                | Mean Recovery %                | RSD% | Mean Recovery %                 | RSD% | Mean Recovery %                 | RSD% | Mean Recovery %                   | RSD% |
| Mecarbam                | 5 to 500                | 0.996          | 103.6                          | 8.3  | 96.8                            | 7.5  | 94.2                            | 7.4  | 98.7                              | 2.5  |
| Methidathion*           | 1 to 500                | 0.991          | 99.7                           | 11.2 | 98.9                            | 2.2  | 90.8                            | 6.9  | 98.8                              | 2.0  |
| Metrafenone*            | 5 to 500                | 0.994          | 95.7                           | 9.3  | 92.8                            | 2.9  | 85.4                            | 6.4  | 94.1                              | 0.6  |
| Dicrotofos              | 5 to 500                | 0.997          | 86.2                           | 11.0 | 96.3                            | 9.5  | 69.3                            | 22.0 | 85.2                              | 5.1  |
| Parathion               | 1 to 500                | 0.990          | 103.4                          | 8.6  | 102.3                           | 1.0  | 101.5                           | 7.3  | 100.3                             | 1.0  |
| Permethrin, (1R)-cis-   | 5 to 500                | 0.996          | 83.5                           | 7.1  | 82.0                            | 5.2  | 77.9                            | 6.3  | 77.5                              | 1.7  |
| Permethrin, (1R)-trans- | 5 to 500                | 0.995          | 92.3                           | 9.0  | 82.0                            | 4.2  | 80.2                            | 8.8  | 80.2                              | 2.6  |
| Phenthoate              | 1 to 500                | 0.996          | 90.9                           | 2.4  | 98.1                            | 3.6  | 95.6                            | 6.1  | 98.3                              | 1.9  |
| Phorate sulfone*        | 2 to 500                | 0.992          | 106.5                          | 7.1  | 94.1                            | 2.8  | 96.3                            | 8.0  | 103.0                             | 5.0  |
| Phosalone               | 1 to 500                | 0.991          | 91.0                           | 9.9  | 91.6                            | 2.2  | 83.9                            | 3.6  | 93.4                              | 1.6  |
| Pirimicarb              | 1 to 500                | 0.993          | 102.8                          | 6.8  | 103.9                           | 2.1  | 100.1                           | 7.3  | 99.8                              | 0.8  |
| Pirimiphos-methyl*      | 1 to 500                | 0.994          | 84.7                           | 12.1 | 98.8                            | 4.1  | 93.7                            | 7.7  | 94.6                              | 2.8  |
| Profenofos*             | 1 to 500                | 0.993          | 82.9                           | 13.1 | 83.7                            | 4.6  | 73.6                            | 9.9  | 82.2                              | 2.9  |
| Propanil                | 1 to 500                | 0.994          | 94.7                           | 5.5  | 105.9                           | 2.5  | 98.0                            | 6.3  | 102.6                             | 1.0  |
| Propiconazole I         | 5 to 500                | 0.995          | 113.5                          | 11.1 | 101.5                           | 2.3  | 88.6                            | 6.7  | 88.7                              | 4.0  |
| Prothiofos              | 1 to 500                | 0.991          | 82.8                           | 7.2  | 83.2                            | 2.6  | 80.5                            | 6.5  | 79.4                              | 1.9  |
| Pyraclostrobin*         | 10 to 500               | 0.991          | —                              | —    | 92.8                            | 12.4 | 87.1                            | 5.0  | 101.5                             | 1.3  |
| Pyrimethanil*           | 1 to 500                | 0.996          | 101.4                          | 4.9  | 104.4                           | 2.9  | 100.9                           | 7.1  | 98.6                              | 1.3  |
| Pyriproxyfen            | 1 to 500                | 0.997          | 80.9                           | 12.2 | 81.1                            | 4.3  | 74.2                            | 5.0  | 83.5                              | 2.1  |
| Quinalphos              | 1 to 500                | 0.995          | 93.9                           | 7.6  | 94.1                            | 2.3  | 94.9                            | 7.0  | 97.3                              | 1.3  |
| Quinoxifen*             | 1 to 500                | 0.992          | 87.0                           | 7.1  | 88.9                            | 1.0  | 83.8                            | 5.1  | 89.6                              | 1.1  |
| Ronnel                  | 1 to 500                | 0.991          | 94.5                           | 8.0  | 101.5                           | 3.1  | 95.1                            | 7.6  | 94.3                              | 0.5  |
| Sulfoxaflor*            | 2 to 500                | 0.998          | 102.6                          | 4.0  | 105.8                           | 6.5  | 105.3                           | 5.2  | 100.7                             | 2.4  |
| Terbufos                | 1 to 500                | 0.995          | 103.5                          | 9.0  | 104.8                           | 1.0  | 98.0                            | 8.4  | 100.0                             | 1.2  |
| Terbufos sulfone*       | 1 to 500                | 0.996          | 98.7                           | 5.8  | 104.2                           | 4.9  | 96.9                            | 7.3  | 95.6                              | 3.0  |
| Tetradifon              | 1 to 500                | 0.997          | 93.1                           | 8.2  | 93.3                            | 4.0  | 85.1                            | 2.1  | 89.3                              | 2.5  |
| Thiabendazole*          | 1 to 500                | 0.994          | 85.5                           | 6.5  | 96.8                            | 3.1  | 88.3                            | 6.8  | 95.8                              | 1.1  |
| Triadimefon*            | 1 to 500                | 0.993          | 94.2                           | 5.7  | 104.8                           | 2.7  | 100.0                           | 7.2  | 100.9                             | 1.7  |
| Triazophos              | 2 to 500                | 0.990          | 93.6                           | 8.3  | 97.3                            | 0.5  | 87.4                            | 6.3  | 97.6                              | 0.5  |

\* Pesticides validated with two instrumental platforms of GC/MS/MS and LC/MS/MS.

**Table 4.** Method quantitation results for 95 pesticides in milk with LC/MS/MS.

| Name                          | Linearity Range (ng/mL) | R <sup>2</sup> | Low QC 5 ng/mL in Milk (n = 4) |      | Low QC 10 ng/mL in Milk (n = 4) |      | Mid QC 50 ng/mL in Milk (n = 4) |      | High QC 100 ng/mL in Milk (n = 4) |      |
|-------------------------------|-------------------------|----------------|--------------------------------|------|---------------------------------|------|---------------------------------|------|-----------------------------------|------|
|                               |                         |                | Mean Recovery %                | RSD% | Mean Recovery %                 | RSD% | Mean Recovery %                 | RSD% | Mean Recovery %                   | RSD% |
| 2,4-D                         | 1 to 500                | 0.999          | 83.1                           | 10.8 | 87.1                            | 16.0 | 94.7                            | 2.5  | 100.3                             | 3.1  |
| 2-Chlorobenzoic acid          | 20 to 500               | 0.995          | —                              | —    | —                               | —    | 98.0                            | 14.5 | 109.7                             | 14.1 |
| 3,6-dichloro- benzoic acid    | 1 to 500                | 0.996          | 92.2                           | 9.6  | 96.6                            | 4.7  | 103.7                           | 2.4  | 105.3                             | 1.9  |
| 4-Hydroxy chlorothalonil      | 1 to 500                | 0.998          | 97.2                           | 16.2 | 102.0                           | 7.8  | 95.0                            | 5.6  | 99.6                              | 1.9  |
| Acetamiprid*                  | 1 to 200                | 0.998          | 103.8                          | 7.3  | 108.5                           | 3.0  | 113.2                           | 0.7  | 110.3                             | 0.8  |
| Aldicarb                      | 1 to 200                | 0.999          | 96.3                           | 5.1  | 102.3                           | 4.9  | 114.4                           | 2.2  | 115.0                             | 1.8  |
| Aldicarb-sulfone (Aldoxycarb) | 1 to 200                | 0.999          | 105.1                          | 2.8  | 104.6                           | 4.7  | 110.7                           | 3.6  | 110.5                             | 2.5  |
| Aldicarb-sulfoxide            | 1 to 500                | 1.000          | 92.6                           | 5.2  | 98.3                            | 3.2  | 102.9                           | 1.1  | 105.7                             | 0.7  |
| Aminopyralid                  | 1 to 100                | 0.994          | 69.7                           | 10.5 | 85.1                            | 6.2  | 87.2                            | 1.7  | 82.9                              | 1.9  |
| Amitraz                       | 1 to 500                | 0.997          | 69.7                           | 9.8  | 66.0                            | 3.0  | 64.2                            | 3.5  | 63.7                              | 1.9  |
| Azoxystrobin*                 | 1 to 500                | 0.999          | 95.4                           | 4.6  | 97.7                            | 1.2  | 99.7                            | 2.1  | 102.4                             | 1.9  |
| Bentazone*                    | 1 to 500                | 1.000          | 91.0                           | 8.6  | 93.1                            | 6.7  | 103.2                           | 2.7  | 105.6                             | 3.0  |
| benzovindiflupyr              | 1 to 500                | 0.995          | 101.6                          | 8.9  | 107.2                           | 4.4  | 119.2                           | 3.1  | 121.7                             | 1.5  |
| Bifenazate                    | 1 to 200                | 0.999          | 99.2                           | 5.9  | 104.9                           | 2.5  | 110.7                           | 0.4  | 110.9                             | 1.0  |
| Bifenazate metabolite B       | 1 to 200                | 0.998          | 91.2                           | 6.9  | 91.5                            | 2.8  | 96.0                            | 1.6  | 101.3                             | 0.5  |
| Bifenthrin*                   | 1 to 500                | 1.000          | 71.9                           | 8.5  | 81.1                            | 2.3  | 84.1                            | 4.7  | 80.2                              | 7.1  |
| Bitertanol*                   | 1 to 500                | 0.997          | 101.9                          | 5.3  | 106.8                           | 2.4  | 115.6                           | 0.7  | 116.2                             | 1.6  |
| Boscalid*                     | 1 to 500                | 0.994          | 102.5                          | 7.2  | 109.7                           | 1.6  | 117.8                           | 1.6  | 117.1                             | 0.4  |
| Buprofezin*                   | 1 to 500                | 0.998          | 89.7                           | 7.3  | 94.3                            | 3.3  | 98.6                            | 1.3  | 102.3                             | 0.2  |
| Carbaryl*                     | 1 to 500                | 0.999          | 96.5                           | 5.1  | 102.3                           | 4.1  | 107.8                           | 1.2  | 108.3                             | 1.3  |
| Carbendazim                   | 1 to 200                | 0.999          | 94.3                           | 6.5  | 94.9                            | 1.7  | 100.1                           | 1.4  | 106.9                             | 0.3  |
| Chlorantraniliprole*          | 1 to 500                | 0.997          | 97.3                           | 8.2  | 103.1                           | 4.4  | 113.7                           | 0.8  | 116.0                             | 1.5  |
| Chlorpyrifos*                 | 1 to 500                | 0.999          | 85.3                           | 2.7  | 91.1                            | 5.6  | 95.9                            | 1.6  | 99.1                              | 0.5  |
| Chlorpyrifos-methyl*          | 2 to 500                | 0.998          | 89.7                           | 8.0  | 98.7                            | 7.2  | 102.9                           | 2.5  | 105.8                             | 2.8  |
| Clofentezine*                 | 1 to 500                | 0.999          | 94.8                           | 9.9  | 98.2                            | 2.6  | 104.5                           | 1.7  | 106.6                             | 1.2  |
| Clothianidin                  | 1 to 500                | 0.994          | 103.8                          | 5.2  | 116.7                           | 3.5  | 120.1                           | 0.7  | 111.5                             | 1.4  |
| Cycloxydime                   | 1 to 500                | 1.000          | 88.9                           | 7.1  | 89.5                            | 3.0  | 94.5                            | 0.4  | 96.7                              | 2.6  |
| Cyhalothrin                   | 1 to 200                | 0.998          | 91.1                           | 7.2  | 103.1                           | 5.2  | 125.9                           | 1.4  | 126.3                             | 1.6  |
| Cypermethrin                  | 1 to 500                | 0.995          | 89.8                           | 4.8  | 100.6                           | 2.6  | 110.1                           | 3.3  | 110.3                             | 1.4  |
| Cyprodinil*                   | 1 to 500                | 0.998          | 89.5                           | 6.7  | 95.7                            | 2.0  | 107.9                           | 2.6  | 107.6                             | 1.6  |
| Cyromazine*                   | 1 to 500                | 0.999          | 55.9                           | 9.7  | 58.1                            | 5.4  | 64.1                            | 1.7  | 64.9                              | 1.2  |
| Diazinon*                     | 1 to 500                | 1.000          | 92.5                           | 3.9  | 97.7                            | 3.6  | 103.5                           | 2.7  | 105.1                             | 0.4  |
| Dicamba                       | 5 to 500                | 0.997          | —                              | —    | 85.3                            | 9.0  | 82.9                            | 10.7 | 90.0                              | 11.1 |
| Dichlorvos*                   | 5 to 500                | 0.999          | —                              | —    | 101.4                           | 8.5  | 112.0                           | 2.6  | 115.2                             | 3.6  |
| Difenoconazole                | 1 to 500                | 0.999          | 93.0                           | 5.1  | 98.4                            | 4.0  | 106.5                           | 1.1  | 107.9                             | 1.6  |
| Diffubenzuron                 | 1 to 200                | 0.996          | 100.2                          | 8.4  | 107.6                           | 2.7  | 114.8                           | 1.7  | 111.3                             | 0.8  |
| Dimethoate*                   | 1 to 500                | 0.998          | 98.4                           | 5.9  | 100.9                           | 3.1  | 107.3                           | 1.3  | 110.4                             | 0.9  |
| Dinotefuran                   | 1 to 500                | 0.998          | 96.5                           | 4.1  | 101.9                           | 4.1  | 111.0                           | 2.2  | 110.3                             | 1.3  |
| Dinotefuran metabolite UF     | 1 to 200                | 1.000          | 88.2                           | 4.6  | 91.4                            | 2.9  | 96.4                            | 1.2  | 98.8                              | 0.3  |
| Diphenylamine*                | 20 to 500               | 0.999          | —                              | —    | —                               | —    | 108.2                           | 7.5  | 123.6                             | 2.5  |
| Ethoprophos*                  | 1 to 500                | 0.998          | 98.3                           | 5.9  | 104.5                           | 3.7  | 111.8                           | 2.0  | 112.6                             | 1.2  |
| Etofenprox                    | 1 to 500                | 0.998          | 75.1                           | 3.0  | 85.9                            | 3.1  | 85.8                            | 3.6  | 78.8                              | 13.1 |
| Famoxadone*                   | 1 to 200                | 0.995          | 102.1                          | 7.3  | 114.0                           | 5.5  | 125.6                           | 3.0  | 118.1                             | 2.5  |

| Name                                                                             | Linearity Range (ng/mL) | R <sup>2</sup> | Low QC 5 ng/mL in Milk (n = 4) |      | Low QC 10 ng/mL in Milk (n = 4) |      | Mid QC 50 ng/mL in Milk (n = 4) |      | High QC 100 ng/mL in Milk (n = 4) |      |
|----------------------------------------------------------------------------------|-------------------------|----------------|--------------------------------|------|---------------------------------|------|---------------------------------|------|-----------------------------------|------|
|                                                                                  |                         |                | Mean Recovery %                | RSD% | Mean Recovery %                 | RSD% | Mean Recovery %                 | RSD% | Mean Recovery %                   | RSD% |
| Fenamiphos                                                                       | 1 to 500                | 0.994          | 96.1                           | 4.2  | 103.8                           | 1.9  | 110.2                           | 2.0  | 114.8                             | 1.1  |
| Fenamiphos sulfone*                                                              | 1 to 200                | 0.998          | 105.9                          | 3.6  | 110.9                           | 3.8  | 115.3                           | 1.1  | 111.4                             | 1.2  |
| Fenamiphos sulfoxide                                                             | 1 to 500                | 0.998          | 101.4                          | 4.1  | 108.4                           | 2.4  | 117.9                           | 1.2  | 114.7                             | 1.1  |
| Fenpropimorph*                                                                   | 1 to 500                | 0.999          | 77.6                           | 0.7  | 79.8                            | 2.8  | 86.1                            | 3.8  | 85.7                              | 0.7  |
| Fenvalerate                                                                      | 1 to 500                | 0.997          | 89.9                           | 10.1 | 107.0                           | 5.5  | 108.4                           | 3.3  | 108.5                             | 2.4  |
| Fipronil Desulfinyl                                                              | 1 to 500                | 1.000          | 90.9                           | 7.8  | 94.8                            | 3.2  | 100.9                           | 1.6  | 105.6                             | 0.9  |
| Fipronil*                                                                        | 1 to 500                | 0.999          | 93.2                           | 2.4  | 98.2                            | 5.9  | 104.0                           | 1.9  | 107.6                             | 1.8  |
| Fipronil sulfide*                                                                | 1 to 500                | 0.999          | 89.9                           | 8.5  | 89.1                            | 5.6  | 101.4                           | 2.4  | 106.9                             | 1.0  |
| Fipronil-sulfone                                                                 | 1 to 500                | 1.000          | 89.8                           | 2.6  | 96.2                            | 9.4  | 101.3                           | 3.7  | 107.3                             | 2.0  |
| Flubendiamide                                                                    | 1 to 500                | 0.998          | 91.3                           | 2.3  | 90.0                            | 6.0  | 97.4                            | 2.6  | 100.2                             | 2.9  |
| Fluopicolide                                                                     | 1 to 500                | 0.995          | 104.6                          | 5.1  | 109.8                           | 3.7  | 117.9                           | 2.2  | 118.4                             | 2.1  |
| Flusilazole*                                                                     | 1 to 500                | 0.995          | 98.0                           | 9.4  | 107.1                           | 4.2  | 115.1                           | 2.4  | 114.6                             | 2.4  |
| Hexythiazox                                                                      | 1 to 500                | 1.000          | 91.5                           | 5.1  | 93.5                            | 2.7  | 99.5                            | 1.1  | 99.2                              | 1.3  |
| Imazapyr                                                                         | 1 to 500                | 0.999          | 76.0                           | 5.5  | 81.9                            | 3.0  | 89.2                            | 2.1  | 88.7                              | 1.1  |
| Imidacloprid                                                                     | 1 to 200                | 0.998          | 98.4                           | 6.4  | 103.5                           | 3.6  | 107.1                           | 0.9  | 108.1                             | 0.9  |
| Isopyrazam*                                                                      | 1 to 500                | 1.000          | 94.2                           | 7.8  | 96.9                            | 1.5  | 102.3                           | 0.7  | 104.9                             | 1.0  |
| Kresoxim-methyl acid                                                             | 1 to 200                | 0.994          | 94.2                           | 6.2  | 103.7                           | 6.4  | 112.2                           | 3.1  | 110.7                             | 0.7  |
| MCPA                                                                             | 1 to 500                | 0.999          | 82.9                           | 9.1  | 90.5                            | 6.8  | 98.0                            | 3.8  | 101.6                             | 2.6  |
| Metaflumizone                                                                    | 1 to 200                | 0.996          | 100.8                          | 6.7  | 112.7                           | 6.2  | 122.5                           | 2.4  | 113.8                             | 1.7  |
| Methamidophos*                                                                   | 1 to 500                | 0.998          | 92.9                           | 5.4  | 98.3                            | 2.7  | 104.4                           | 1.1  | 106.3                             | 0.5  |
| Methidathion*                                                                    | 1 to 500                | 0.999          | 98.0                           | 3.8  | 103.2                           | 2.1  | 108.5                           | 1.1  | 109.0                             | 1.8  |
| Metrafenone*                                                                     | 1 to 500                | 1.000          | 92.6                           | 6.0  | 93.4                            | 4.0  | 99.6                            | 2.1  | 101.5                             | 0.6  |
| Novaluron                                                                        | 1 to 500                | 0.998          | 101.1                          | 6.7  | 104.7                           | 1.3  | 111.9                           | 1.4  | 113.5                             | 1.0  |
| Oxamyl*                                                                          | 1 to 500                | 1.000          | 99.2                           | 5.2  | 103.7                           | 3.1  | 108.3                           | 0.5  | 108.7                             | 1.3  |
| Penthiopyrad                                                                     | 1 to 500                | 0.998          | 94.3                           | 4.8  | 98.5                            | 3.4  | 105.1                           | 2.0  | 107.8                             | 0.9  |
| Penthiopyrad metabolite (3-(Trifluoromethyl)-1-methyl-1H-pyrazole-4-carboxamide) | 1 to 200                | 0.999          | 94.9                           | 5.2  | 97.9                            | 2.9  | 102.6                           | 1.3  | 108.5                             | 0.6  |
| Phorate*                                                                         | 1 to 500                | 0.999          | 89.5                           | 11.8 | 97.3                            | 1.8  | 107.1                           | 1.1  | 109.0                             | 1.6  |
| Phorate Sulfone*                                                                 | 1 to 500                | 0.996          | 95.3                           | 2.6  | 102.8                           | 2.2  | 111.7                           | 1.4  | 113.2                             | 0.5  |
| Phorate Sulfoxide                                                                | 1 to 500                | 0.999          | 96.8                           | 5.0  | 99.8                            | 3.0  | 106.0                           | 2.1  | 109.1                             | 1.3  |
| Pirimiphos-methyl*                                                               | 1 to 500                | 0.999          | 92.9                           | 4.4  | 95.8                            | 2.3  | 100.7                           | 2.6  | 101.8                             | 1.1  |
| Prochloraz*                                                                      | 1 to 500                | 1.000          | 92.2                           | 8.5  | 94.0                            | 2.7  | 100.8                           | 1.6  | 102.5                             | 1.0  |
| Profenofos*                                                                      | 1 to 500                | 0.998          | 94.6                           | 6.8  | 101.2                           | 1.1  | 108.0                           | 1.5  | 108.7                             | 0.4  |
| Propargite                                                                       | 1 to 500                | 1.000          | 93.5                           | 9.6  | 98.5                            | 2.7  | 107.7                           | 2.9  | 107.4                             | 1.5  |
| Propiconazole                                                                    | 1 to 500                | 0.997          | 99.5                           | 6.1  | 102.8                           | 3.5  | 113.4                           | 3.9  | 113.2                             | 0.8  |
| Prothioconazole-desthio                                                          | 1 to 500                | 0.999          | 95.1                           | 6.8  | 96.6                            | 2.7  | 99.0                            | 1.6  | 100.5                             | 1.1  |
| Pyraclostrobin*                                                                  | 1 to 500                | 1.000          | 94.8                           | 6.6  | 98.4                            | 2.5  | 104.0                           | 1.4  | 105.3                             | 1.1  |
| Pyrimethanil*                                                                    | 1 to 500                | 0.999          | 95.2                           | 6.1  | 99.9                            | 4.1  | 108.8                           | 1.1  | 109.8                             | 0.9  |
| Quinoxifen*                                                                      | 1 to 500                | 0.999          | 83.1                           | 9.5  | 85.9                            | 1.6  | 90.2                            | 1.6  | 92.1                              | 0.8  |
| Saflufenacil                                                                     | 1 to 200                | 0.998          | 104.2                          | 3.5  | 110.8                           | 2.1  | 118.1                           | 1.3  | 114.7                             | 1.6  |
| Semiamitraz                                                                      | 1 to 500                | 0.996          | 73.7                           | 9.8  | 87.6                            | 5.3  | 95.9                            | 2.3  | 97.4                              | 2.1  |
| Spirodiclofen*                                                                   | 1 to 500                | 1.000          | 89.2                           | 5.4  | 92.9                            | 3.1  | 96.9                            | 1.8  | 98.8                              | 0.7  |
| Spirotetramat                                                                    | 1 to 500                | 0.994          | 91.7                           | 7.0  | 107.9                           | 4.1  | 112.3                           | 3.7  | 112.2                             | 3.2  |

| Name                                       | Linearity Range (ng/mL) | R <sup>2</sup> | Low QC 5 ng/mL in Milk (n = 4) |      | Low QC 10 ng/mL in Milk (n = 4) |      | Mid QC 50 ng/mL in Milk (n = 4) |      | High QC 100 ng/mL in Milk (n = 4) |      |
|--------------------------------------------|-------------------------|----------------|--------------------------------|------|---------------------------------|------|---------------------------------|------|-----------------------------------|------|
|                                            |                         |                | Mean Recovery %                | RSD% | Mean Recovery %                 | RSD% | Mean Recovery %                 | RSD% | Mean Recovery %                   | RSD% |
| Spirotetramet metabolite BYI08330-cis-enol | 1 to 200                | 1.000          | 95.9                           | 5.6  | 100.6                           | 3.2  | 109.2                           | 0.9  | 109.5                             | 0.6  |
| Sulfoxaflor*                               | 1 to 500                | 0.998          | 100.2                          | 5.2  | 105.5                           | 3.9  | 112.0                           | 0.7  | 113.4                             | 1.0  |
| Tebufenozide                               | 1 to 500                | 0.998          | 102.4                          | 8.4  | 109.3                           | 5.9  | 116.2                           | 6.2  | 115.3                             | 1.9  |
| Terbufos Sulfone*                          | 1 to 500                | 1.000          | 90.5                           | 3.0  | 95.9                            | 0.9  | 102.4                           | 2.3  | 106.6                             | 1.1  |
| Terbufos-sulfoxide                         | 1 to 500                | 0.993          | 95.4                           | 6.9  | 100.8                           | 2.5  | 107.8                           | 1.6  | 113.6                             | 0.3  |
| Thiabendazole*                             | 1 to 500                | 0.996          | 96.0                           | 6.7  | 104.6                           | 4.9  | 112.8                           | 1.4  | 113.1                             | 2.3  |
| Thiacloprid                                | 1 to 200                | 1.000          | 97.9                           | 5.5  | 101.1                           | 4.0  | 105.6                           | 1.2  | 109.0                             | 1.5  |
| Thiamethoxam                               | 1 to 200                | 0.999          | 99.8                           | 6.4  | 108.7                           | 3.6  | 116.1                           | 0.9  | 117.9                             | 1.7  |
| Triadimefon*                               | 1 to 200                | 0.999          | 96.6                           | 2.3  | 102.9                           | 3.4  | 109.8                           | 1.5  | 108.3                             | 0.7  |
| Triadimenol                                | 1 to 500                | 0.999          | 99.9                           | 7.0  | 104.4                           | 3.0  | 108.2                           | 0.6  | 108.0                             | 1.2  |
| Triforine                                  | 1 to 200                | 0.999          | 92.7                           | 5.2  | 102.9                           | 3.8  | 114.6                           | 2.4  | 112.6                             | 1.3  |

\* Pesticides validated with two instrumental platforms of GC/MS/MS and LC/MS/MS.

## Conclusion

A rapid, reliable, and robust workflow using QuEChERS extraction followed by Agilent Captiva EMR—Lipid cartridge cleanup was developed and verified for the analysis of 171 pesticide multiresidues in milk using LC/MS/MS and GC/MS/MS. The sample preparation procedure was investigated and optimized for extraction method, matrix removal, elution efficiency, and water removal. Captiva EMR—Lipid cleanup provides the best matrix removal efficiency and the least instrumental contamination. The total workflow delivers satisfactory quantitation results, and demonstrates an efficient solution for the analysis of multiclass multiresidue pesticides in milk.

## References

1. Abou Donia, M. A. *et al.* Chemical Composition of Raw Milk and the Accumulation of Pesticide Residues in Milk Products. *Global Veterinaria* **2010**, 4(1), 6.
2. GB 2763-2019 National Food Safety Standard—Maximum Residue Limits for Pesticides in Food: <https://www.codeofchina.com/standard/GB2763-2019.html>.
3. Yang, X. Analysis of Formamidine Pesticides and Metabolites in Pork and Porcine Liver Using Agilent Captiva EMR—Lipid and LC/MS/MS. *Agilent Technologies application note*, publication number 5994-0357EN, **2019**.
4. Zhao, L. Determination of Multiclass, Multiresidue Pesticides in Olive Oils by Captiva EMR—Lipid Cleanup and GC/MS/MS. *Agilent Technologies application note*, publication number 5994-0405EN, **2018**.

## Appendix

### MRM parameters setting of GC/MS/MS and LC/MS/MS

| Compound Name                   | GC/MS/MS MRM Setting |                  |                      |
|---------------------------------|----------------------|------------------|----------------------|
|                                 | RT (min)             | Quant Transition | Qualifier Transition |
| 2,4,6-Trichlorophenol           | 7.73                 | 131.8 -> 97.0    | 96.9 -> 62.0         |
| Acetamidiprid                   | 27.87                | 126.0 -> 73.0    | 152.0 -> 116.1       |
| Aldrin                          | 19.57                | 262.9 -> 192.9   | 254.9 -> 220.0       |
| Azinphos-ethyl                  | 30.62                | 132.0 -> 77.1    | 160.0 -> 77.1        |
| Azinphos-methyl                 | 29.35                | 160.0 -> 77.0    | 160.0 -> 132.1       |
| Azoxystrobin                    | 37.06                | 344.1 -> 329.0   | 344.1 -> 171.9       |
| Bentazone                       | 20.36                | 119.0 -> 92.0    | 198.0 -> 119.0       |
| Bifenthrin                      | 28.33                | 181.2 -> 165.2   | 181.2 -> 166.2       |
| Bitertanol                      | 31.51                | 170.1 -> 141.1   | 170.1 -> 115.0       |
| Boscalid                        | 33.36                | 140.0 -> 112.0   | 140.0 -> 76.0        |
| Buprofezin                      | 23.76                | 104.0 -> 51.0    | 104.0 -> 77.0        |
| Captan                          | 21.42                | 151.0 -> 80.0    | 149.0 -> 79.1        |
| Carbaryl                        | 18.25                | 144.1 -> 116.1   | 144.1 -> 89.0        |
| Chinomethionate (Oxythioquinox) | 21.89                | 233.9 -> 206.1   | 206.0 -> 148.1       |
| Chlorantraniliprole             | 28.34                | 277.8 -> 215.0   | 277.8 -> 248.8       |
| Chlordane-cis                   | 22.55                | 271.8 -> 236.9   | 372.8 -> 265.9       |
| Chlordane-oxy                   | 21.14                | 114.9 -> 51.1    | 114.9 -> 87.0        |
| Chlordane-trans                 | 21.99                | 271.7 -> 236.9   | 372.8 -> 265.8       |
| Chlorfenvinphos                 | 21.55                | 266.9 -> 159.1   | 322.8 -> 266.8       |
| Chlorpropham                    | 13.31                | 153.0 -> 90.0    | 153.0 -> 125.1       |
| Chlorpyrifos                    | 19.99                | 198.9 -> 171.0   | 196.9 -> 169.0       |
| Chlorpyrifos-methyl             | 18.10                | 285.9 -> 93.0    | 287.9 -> 92.9        |
| Clofentezine                    | 5.28                 | 136.7 -> 102.0   | 138.7 -> 102.0       |
| Coumaphos                       | 31.97                | 210.0 -> 182.0   | 361.9 -> 109.0       |
| Cyfluthrin-1                    | 32.79                | 226.0 -> 206.0   | 198.9 -> 170.1       |
| Cyfluthrin-2                    | 32.97                | 226.0 -> 206.0   | 198.9 -> 170.1       |
| Cyfluthrin-3                    | 33.12                | 226.0 -> 206.0   | 198.9 -> 170.1       |
| Cyfluthrin-4                    | 33.20                | 226.0 -> 206.0   | 198.9 -> 170.1       |
| Cypermethrin-1                  | 33.11                | 163.0 -> 91.0    | 163.0 -> 127.0       |
| Cypermethrin-2                  | 33.20                | 163.0 -> 91.0    | 163.0 -> 127.0       |
| Cypermethrin-3                  | 33.37                | 163.0 -> 127.0   | 163.0 -> 91.0        |
| Cypermethrin-4                  | 33.56                | 163.0 -> 91.0    | 163.0 -> 127.0       |
| Cyprodinil                      | 20.90                | 225.2 -> 224.3   | 224.2 -> 208.2       |
| Cyromazine                      | 15.47                | 151.0 -> 109.0   | 165.9 -> 151.0       |
| DDD- <i>o,p'</i>                | 23.72                | 235.0 -> 165.2   | 237.0 -> 165.2       |
| DDD- <i>p,p'</i>                | 24.93                | 234.9 -> 165.1   | 236.9 -> 165.2       |
| DDE- <i>p,p'</i>                | 23.42                | 246.1 -> 176.2   | 315.8 -> 246.0       |
| DDT- <i>o,p'</i>                | 25.04                | 235.0 -> 165.2   | 237.0 -> 165.2       |
| DDT- <i>p,p'</i>                | 26.27                | 235.0 -> 165.2   | 237.0 -> 165.2       |
| Deltamethrin                    | 36.52                | 252.9 -> 93.0    | 250.7 -> 172.0       |
| Demeton-S-methyl                | 12.70                | 88.0 -> 60.0     | 142.0 -> 78.9        |
| Diazinon                        | 16.42                | 137.1 -> 84.0    | 137.1 -> 54.0        |

| Compound Name                       | GC/MS/MS MRM Setting |                  |                      |
|-------------------------------------|----------------------|------------------|----------------------|
|                                     | RT (min)             | Quant Transition | Qualifier Transition |
| Dichlofenthion                      | 17.76                | 278.9 -> 222.9   | 222.9 -> 204.9       |
| Dichloran                           | 14.74                | 206.1 -> 176.0   | 160.1 -> 124.1       |
| Dichlorvos                          | 6.13                 | 109.0 -> 79.0    | 184.9 -> 93.0        |
| Dicrotofos                          | 13.75                | 127.0 -> 109.0   | 127.0 -> 95.0        |
| Dieldrin                            | 23.38                | 262.9 -> 193.0   | 277.0 -> 241.0       |
| Difenoconazole I                    | 35.85                | 322.8 -> 264.8   | 264.9 -> 202.0       |
| Difenoconazole II                   | 35.98                | 322.8 -> 264.8   | 264.9 -> 202.0       |
| Dimethipin                          | 15.25                | 118.0 -> 58.0    | 124.0 -> 76.0        |
| Dimethoate                          | 14.85                | 87.0 -> 46.0     | 142.9 -> 111.0       |
| Diphenylamine                       | 12.70                | 169.0 -> 168.2   | 168.0 -> 167.2       |
| Endosulfan I ( <i>alpha</i> isomer) | 22.42                | 194.9 -> 159.0   | 194.9 -> 125.0       |
| Endosulfan II ( <i>beta</i> isomer) | 24.51                | 206.9 -> 172.0   | 194.9 -> 124.9       |
| Endosulfan sulfate                  | 26.03                | 271.9 -> 237.0   | 273.8 -> 238.9       |
| Endrin                              | 24.16                | 262.8 -> 193.0   | 244.8 -> 173.0       |
| Ethion                              | 25.19                | 230.9 -> 129.0   | 230.9 -> 175.0       |
| Ethofenprox                         | 33.92                | 163.0 -> 107.1   | 163.0 -> 135.1       |
| Ethoprophos                         | 12.99                | 157.9 -> 97.0    | 157.9 -> 114.0       |
| Famoxadone                          | 37.06                | 197.0 -> 115.0   | 223.9 -> 196.2       |
| Fenamidone                          | 28.62                | 238.0 -> 237.2   | 268.0 -> 180.2       |
| Fenamiphos sulfone                  | 27.89                | 319.8 -> 292.0   | 171.0 -> 107.0       |
| Fenitrothion                        | 19.17                | 277.0 -> 260.1   | 277.0 -> 109.0       |
| Fenpropathrin                       | 28.52                | 181.1 -> 152.1   | 207.9 -> 181.0       |
| Fenpropimorph                       | 19.98                | 128.1 -> 70.1    | 128.1 -> 110.1       |
| Fensulfothion                       | 24.77                | 291.8 -> 156.0   | 291.8 -> 108.8       |
| Fenthion                            | 19.90                | 278.0 -> 109.0   | 278.0 -> 169.0       |
| Fenvalerate I                       | 35.11                | 167.0 -> 125.1   | 224.9 -> 119.0       |
| Fenvalerate II                      | 35.51                | 167.0 -> 125.1   | 224.9 -> 119.0       |
| Fipronil                            | 21.64                | 366.8 -> 212.8   | 368.8 -> 214.8       |
| Fipronil sulfide                    | 21.38                | 351.0 -> 254.9   | 420.0 -> 350.9       |
| Fipronil sulfone                    | 23.96                | 382.8 -> 254.9   | 384.8 -> 256.8       |
| Flusilazole                         | 23.86                | 233.0 -> 165.1   | 233.0 -> 91.0        |
| HCH- <i>alpha</i>                   | 14.30                | 216.9 -> 181.0   | 218.9 -> 183.0       |
| HCH- <i>beta</i>                    | 15.34                | 181.0 -> 145.0   | 216.9 -> 181.1       |
| HCH- <i>delta</i>                   | 16.50                | 181.1 -> 145.1   | 217.0 -> 181.1       |
| HCH- <i>gamma</i>                   | 15.56                | 181.0 -> 145.0   | 216.9 -> 181.0       |
| Heptachlor                          | 18.28                | 271.7 -> 236.9   | 273.7 -> 238.9       |
| Heptachlor exo-epoxide              | 21.10                | 352.8 -> 262.9   | 354.8 -> 264.9       |
| Hexachlorobenzene                   | 14.56                | 283.8 -> 213.9   | 283.8 -> 248.8       |
| Isoprazam                           | 31.01                | 159.0 -> 42.1    | 159.0 -> 139.0       |
| Malathion                           | 19.65                | 126.9 -> 99.0    | 172.9 -> 99.0        |
| Mecarbam                            | 21.63                | 158.9 -> 131.0   | 130.9 -> 74.0        |

| Compound Name           | GC/MS/MS MRM Setting |                  |                      |
|-------------------------|----------------------|------------------|----------------------|
|                         | RT (min)             | Quant Transition | Qualifier Transition |
| Methacrifos             | 10.43                | 207.9 -> 180.1   | 207.9 -> 93.0        |
| Methamidophos           | 5.84                 | 141.0 -> 95.0    | 141.0 -> 79.0        |
| Methidathion            | 22.09                | 144.9 -> 85.0    | 144.9 -> 58.1        |
| Metrafenone             | 30.98                | 208.9 -> 166.0   | 394.8 -> 364.8       |
| Oxamyl                  | 11.02                | 162.0 -> 114.9   | 98.0 -> 58.0         |
| Parathion               | 20.01                | 139.0 -> 109.0   | 290.9 -> 109.0       |
| Pentachloronitrobenzene | 15.76                | 295.0 -> 237.0   | 236.9 -> 142.9       |
| Permethrin, (1R)-cis-   | 31.61                | 183.1 -> 168.1   | 183.1 -> 153.0       |
| Permethrin, (1R)-trans- | 31.85                | 183.1 -> 168.1   | 183.1 -> 153.0       |
| Phenthoate              | 21.66                | 273.7 -> 121.0   | 273.7 -> 124.9       |
| Phorate                 | 14.20                | 260.0 -> 75.0    | 230.9 -> 128.9       |
| Phorate sulfone         | 19.76                | 124.9 -> 96.9    | 153.0 -> 97.0        |
| Phosalone               | 29.38                | 182.0 -> 111.0   | 182.0 -> 102.1       |
| Phosmet                 | 27.97                | 160.0 -> 77.1    | 160.0 -> 133.1       |
| Pirimicarb              | 17.37                | 166.0 -> 55.1    | 238.0 -> 166.2       |
| Pirimiphos-methyl       | 19.30                | 290.0 -> 125.0   | 232.9 -> 151.0       |
| Prochloraz              | 32.09                | 195.9 -> 96.9    | 180.0 -> 138.0       |

| Compound Name    | GC/MS/MS MRM Setting |                  |                      |
|------------------|----------------------|------------------|----------------------|
|                  | RT (min)             | Quant Transition | Qualifier Transition |
| Profenofos       | 23.30                | 207.9 -> 63.0    | 338.8 -> 268.7       |
| Propanil         | 17.70                | 161.0 -> 99.0    | 161.0 -> 90.0        |
| Propiconazole I  | 26.16                | 172.9 -> 145.0   | 172.9 -> 74.0        |
| Prothiofos       | 23.19                | 266.9 -> 239.0   | 308.9 -> 238.9       |
| Pyraclostrobin   | 35.18                | 132.0 -> 104.0   | 132.0 -> 77.1        |
| Pyrimethanil     | 16.15                | 198.0 -> 118.1   | 198.0 -> 183.1       |
| Pyriproxyfen     | 29.61                | 136.1 -> 78.1    | 136.1 -> 96.0        |
| Quinalphos       | 21.63                | 146.0 -> 118.0   | 146.0 -> 91.0        |
| Quinoxifen       | 26.03                | 271.9 -> 237.1   | 237.0 -> 208.1       |
| Ronnel           | 18.64                | 285.0 -> 269.9   | 286.9 -> 272.0       |
| Spirodiclofen    | 31.55                | 109.1 -> 81.1    | 109.1 -> 79.1        |
| Sulfoxaflor      | 12.70                | 173.7 -> 104.1   | 173.7 -> 154.0       |
| Terbufos         | 15.86                | 230.9 -> 175.0   | 230.9 -> 129.0       |
| Terbufos sulfone | 21.22                | 153.0 -> 97.0    | 198.9 -> 96.9        |
| Tetradifon       | 29.02                | 158.9 -> 131.0   | 226.9 -> 199.0       |
| Thiabendazole    | 21.22                | 201.0 -> 174.0   | 201.9 -> 175.0       |
| Triadimefon      | 20.10                | 208.0 -> 181.1   | 208.0 -> 111.0       |
| Triazophos       | 25.64                | 161.2 -> 134.2   | 161.2 -> 106.1       |

| Compound Name                 | LC/MS/MS MRM Setting |               |             |            |                  |          |
|-------------------------------|----------------------|---------------|-------------|------------|------------------|----------|
|                               | RT (min)             | Precursor Ion | Product Ion | Fragmentor | Collision Energy | Polarity |
| 2,4-D                         | 5.79                 | 219.0         | 161.0       | 90         | 15               | -        |
|                               |                      | 221.0         | 163.0       |            | 15               |          |
| 2-Chlorobenzoic acid          | 4.19                 | 155.0         | 111.0       | 65         | 4                | -        |
|                               |                      | 155.0         | 35.1        |            | 10               |          |
| 3,6-Dichloro- benzoic acid    | 4.03                 | 204.9         | 160.9       | 80         | 10               | -        |
|                               |                      | 204.9         | 125.0       |            | 22               |          |
| 4-Hydroxy chlorothalonil      | 4.89                 | 244.9         | 181.9       | 146        | 34               | -        |
|                               |                      | 244.9         | 174.9       |            | 30               |          |
| Acetamiprid                   | 3.60                 | 223.0         | 126.1       | 80         | 18               | -        |
|                               |                      | 223.0         | 99.0        |            | 44               |          |
| Aldicarb                      | 4.52                 | 208.0         | 116.0       | 65         | 6                | +        |
|                               |                      | 208.0         | 89.1        |            | 10               |          |
| Aldicarb-sulfone (Aldoxycarb) | 2.92                 | 223.1         | 86.1        | 80         | 8                | +        |
|                               |                      | 223.1         | 76.0        |            | 0                |          |
| Aldicarb-sulfoxide            | 2.85                 | 207.1         | 131.9       | 65         | 2                | +        |
|                               |                      | 207.1         | 105.2       |            | 4                |          |
| Aminopyralid                  | 2.77                 | 207.0         | 161.0       | 90         | 20               | +        |
|                               |                      | 207.0         | 134.0       |            | 32               |          |
| Amitraz                       | 14.30                | 294.2         | 163.0       | 90         | 12               | +        |
|                               |                      | 294.2         | 122.0       |            | 32               |          |
| Azoxystrobin                  | 7.11                 | 404.0         | 372.2       | 105        | 10               | +        |
|                               |                      | 404.0         | 344.0       |            | 24               |          |
| Bentazone                     | 4.66                 | 239.0         | 197.0       | 105        | 17               | -        |
|                               |                      | 239.0         | 132.0       |            | 25               |          |
| Benzovindiflupyr              | 9.48                 | 398.1         | 378.0       | 90         | 14               | +        |
|                               |                      | 398.1         | 342.1       |            | 18               |          |

| Compound Name           | LC/MS/MS MRM Setting |               |             |            |                  |          |
|-------------------------|----------------------|---------------|-------------|------------|------------------|----------|
|                         | RT (min)             | Precursor Ion | Product Ion | Fragmentor | Collision Energy | Polarity |
| Bifenazate              | 8.32                 | 301.1         | 198.2       | 90         | 5                | +        |
|                         |                      | 301.1         | 170.1       |            | 18               |          |
| Bifenazate metabolite B | 10.78                | 299.1         | 213.0       | 85         | 10               | +        |
|                         |                      | 299.1         | 197.0       |            | 18               |          |
| Bifenthrin              | 15.90                | 440.2         | 198.2       | 100        | 4                | +        |
|                         |                      | 440.2         | 181.0       |            | 14               |          |
| Bitertanol              | 10.55                | 338.1         | 269.2       | 70         | 2                | +        |
|                         |                      | 338.1         | 251.2       |            | 6                |          |
| Boscalid                | 7.65                 | 343.2         | 307.2       | 140        | 20               | +        |
|                         |                      | 343.2         | 271.0       |            | 35               |          |
| Buprofezin              | 12.42                | 306.1         | 201.2       | 105        | 9                | +        |
|                         |                      | 306.1         | 116.0       |            | 15               |          |
| Carbaryl                | 5.42                 | 202.0         | 145.0       | 65         | 2                | +        |
|                         |                      | 202.0         | 127.1       |            | 28               |          |
| Carbendazim             | 3.40                 | 192.1         | 160.1       | 105        | 16               | +        |
|                         |                      | 192.1         | 132.1       |            | 32               |          |
| Chlorantraniliprole     | 6.76                 | 483.9         | 452.9       | 105        | 16               | +        |
|                         |                      | 483.9         | 286.0       |            | 12               |          |
| Chlorpyrifos            | 13.05                | 351.9         | 199.9       | 100        | 15               | +        |
|                         |                      | 349.9         | 197.9       |            | 22               |          |
| Chlorpyrifos-methyl     | 11.08                | 322.0         | 290.0       | 110        | 10               | +        |
|                         |                      | 322.0         | 125.0       |            | 25               |          |
| Clofentezine            | 10.55                | 303.0         | 138.0       | 80         | 15               | +        |
|                         |                      | 303.0         | 102.1       | 110        | 40               |          |
| Clothianidin            | 3.42                 | 250.0         | 169.0       | 90         | 8                | +        |
|                         |                      | 250.0         | 131.9       |            | 8                |          |
| Cycloxydime             | 11.33                | 326.2         | 280.1       | 90         | 14               | +        |
|                         |                      | 326.2         | 180.0       |            | 22               |          |
| Cyhalothrin             | 13.99                | 467.1         | 450.1       | 90         | 4                | +        |
|                         |                      | 467.1         | 225.0       |            | 12               |          |
| Cypermethrin            | 14.18                | 435.1         | 193.0       | 100        | 10               | +        |
|                         |                      | 433.1         | 191.0       |            | 12               |          |
| Cyprodinil              | 9.51                 | 226.1         | 108.1       | 140        | 24               | +        |
|                         |                      | 226.1         | 93.1        |            | 40               |          |
| Cyromazine              | 2.64                 | 167.0         | 125.0       | 120        | 20               | +        |
|                         |                      | 167.0         | 108.0       |            | 15               |          |
| Diazinon                | 10.37                | 305.1         | 169.0       | 105        | 26               | +        |
|                         |                      | 305.1         | 153.1       |            | 24               |          |
| Dicamba                 | 4.20                 | 221.0         | 177.0       | 60         | 2                | -        |
|                         |                      | 219.0         | 175.0       |            | 2                |          |
| Dichlorvos              | 5.04                 | 221.0         | 109.0       | 100        | 12               | +        |
|                         |                      | 221.0         | 95.1        |            | 32               |          |
| Difenoconazole          | 11.08                | 406.0         | 251.1       | 120        | 25               | +        |
|                         |                      | 406.0         | 188.1       |            | 46               |          |
| Diflubenzuron           | 9.09                 | 311.0         | 158.0       | 80         | 18               | +        |
|                         |                      | 311.0         | 141.0       |            | 46               |          |
| Dimethoate              | 3.70                 | 230.0         | 199.0       | 70         | 4                | +        |
|                         |                      | 230.0         | 171.0       |            | 12               |          |

| Compound Name             | LC/MS/MS MRM Setting |               |             |            |                  |          |
|---------------------------|----------------------|---------------|-------------|------------|------------------|----------|
|                           | RT (min)             | Precursor Ion | Product Ion | Fragmentor | Collision Energy | Polarity |
| Dinotefuran               | 2.82                 | 203.1         | 157.0       | 85         | 2                | +        |
|                           |                      | 203.1         | 129.0       |            | 6                |          |
| Dinotefuran metabolite UF | 2.83                 | 159.1         | 102.1       | 85         | 10               | +        |
|                           |                      | 159.1         | 67.1        |            | 22               |          |
| Diphenylamine             | 8.42                 | 170.1         | 93.1        | 115        | 32               | +        |
|                           |                      | 170.1         | 92.6        |            | 24               |          |
| Ethoprophos               | 9.02                 | 243.0         | 173.1       | 90         | 12               | +        |
|                           |                      | 243.0         | 131.0       |            | 16               |          |
| Etofenprox                | 15.79                | 394.2         | 359.2       | 100        | 12               | +        |
|                           |                      | 394.2         | 177.2       |            | 13               |          |
| Famoxadone                | 10.14                | 392.0         | 331.1       | 75         | 6                | +        |
|                           |                      | 392.0         | 238.1       |            | 18               |          |
| Fenamiphos                | 9.30                 | 304.1         | 234.0       | 120        | 16               | +        |
|                           |                      | 304.1         | 217.1       |            | 20               |          |
| Fenamiphos sulfone        | 5.15                 | 336.1         | 308.0       | 110        | 12               | +        |
|                           |                      | 336.1         | 266.0       |            | 16               |          |
| Fenamiphos sulfoxide      | 5.02                 | 320.1         | 233.0       | 115        | 20               | +        |
|                           |                      | 320.1         | 171.1       |            | 16               |          |
| Fenpropimorph             | 6.99                 | 304.0         | 147.0       | 120        | 30               | +        |
|                           |                      | 304.0         | 130.0       |            | 30               |          |
| Fenvalerate               | 14.80                | 437.2         | 181.0       | 90         | 38               | +        |
|                           |                      | 437.2         | 167.0       |            | 14               |          |
| Fipronil Desulfinyl       | 8.80                 | 387.0         | 351.0       | 90         | 14               | -        |
|                           |                      | 387.0         | 331.0       |            | 34               |          |
| Fipronil                  | 9.24                 | 435.0         | 330.0       | 70         | 12               | -        |
|                           |                      | 435.0         | 250.0       |            | 28               |          |
| Fipronil sulfide          | 9.50                 | 419.2         | 383.2       | 80         | 18               | -        |
|                           |                      | 419.2         | 262.0       |            | 13               |          |
| Fipronil-sulfone          | 9.92                 | 450.9         | 415.0       | 100        | 9                | -        |
|                           |                      | 450.9         | 282.0       |            | 25               |          |
| Flubendiamide             | 9.61                 | 681.0         | 274.0       | 101        | 14               | -        |
|                           |                      | 681.0         | 271.9       |            | 14               |          |
| Fluopicolide              | 8.02                 | 385.0         | 147.0       | 112        | 60               | +        |
|                           |                      | 382.9         | 172.9       | 110        | 32               |          |
| Flusilazole               | 9.26                 | 316.0         | 247.2       | 120        | 15               | +        |
|                           |                      | 316.0         | 219.1       |            | 26               |          |
| Hexythiazox               | 13.11                | 353.0         | 228.1       | 90         | 12               | +        |
|                           |                      | 353.0         | 194.1       |            | 24               |          |
| Imazapyr                  | 3.47                 | 262.1         | 217.0       | 121        | 22               | +        |
|                           |                      | 262.1         | 69.2        |            | 30               |          |
| Imidacloprid              | 3.36                 | 256.0         | 209.1       | 80         | 12               | +        |
|                           |                      | 256.0         | 175.1       |            | 22               |          |
| Isopyrazam                | 11.17                | 360.0         | 320.0       | 75         | 10               | +        |
|                           |                      | 360.0         | 244.1       |            | 20               |          |
| Kresoxim-methyl acid      | 7.44                 | 300.1         | 253.0       | 85         | 2                | +        |
|                           |                      | 300.1         | 222.0       |            | 10               |          |
| MCPA                      | 6.10                 | 201.0         | 143.0       | 100        | 15               | -        |
|                           |                      | 199.0         | 141.0       |            | 15               |          |



| Compound Name           | LC/MS/MS MRM Setting |               |             |            |                  |          |
|-------------------------|----------------------|---------------|-------------|------------|------------------|----------|
|                         | RT (min)             | Precursor Ion | Product Ion | Fragmentor | Collision Energy | Polarity |
| Metaflumizone           | 12.29                | 507.0         | 287.1       | 150        | 20               | +        |
|                         |                      | 507.0         | 267.0       |            | 40               |          |
| Methamidophos           | 2.75                 | 142.0         | 125.0       | 85         | 10               | +        |
|                         |                      | 142.0         | 94.1        |            | 12               |          |
| Methidathion            | 6.71                 | 302.9         | 145.0       | 55         | 6                | +        |
|                         |                      | 302.9         | 85.1        |            | 20               |          |
| Metrafenone             | 10.38                | 409.1         | 226.9       | 106        | 22               | +        |
|                         |                      | 409.1         | 209.0       |            | 14               |          |
| Novaluron               | 11.55                | 493.1         | 310.0       | 90         | 16               | +        |
|                         |                      | 493.1         | 158.1       |            | 18               |          |
| Oxamyl                  | 2.94                 | 237.1         | 220.0       | 50         | 2                | +        |
|                         |                      | 237.1         | 90.0        |            | 2                |          |
| Penthiopyrad            | 9.36                 | 360.1         | 276.0       | 90         | 14               | +        |
|                         |                      | 360.1         | 256.0       |            | 22               |          |
| Penthiopyrad metabolite | 3.05                 | 194.1         | 174.0       | 80         | 6                | +        |
|                         |                      | 194.1         | 154.0       |            | 14               |          |
| Phorate                 | 10.69                | 261.0         | 199.0       | 60         | 2                | +        |
|                         |                      | 261.0         | 75.1        |            | 6                |          |
| Phorate Sulfone         | 6.07                 | 293.0         | 171.0       | 80         | 6                | +        |
|                         |                      | 293.0         | 143.0       |            | 15               |          |
| Phorate Sulfoxide       | 5.92                 | 277.0         | 199.0       | 80         | 5                | +        |
|                         |                      | 277.0         | 171.0       |            | 12               |          |
| Pirimiphos-methyl       | 10.72                | 306.0         | 164.2       | 130        | 24               | +        |
|                         |                      | 306.0         | 136.1       |            | 32               |          |
| Prochloraz              | 10.42                | 376.0         | 308.0       | 70         | 4                | +        |
|                         |                      | 376.0         | 265.9       |            | 12               |          |
| Profenofos              | 11.99                | 374.9         | 347.0       | 120        | 9                | +        |
|                         |                      | 373.0         | 344.9       |            | 10               |          |
| Propargite              | 13.55                | 368.1         | 81.0        | 82         | 44               | +        |
|                         |                      | 368.0         | 231.1       |            | 1                |          |
| Propiconazole           | 10.29                | 342.0         | 158.9       | 115        | 36               | +        |
|                         |                      | 342.0         | 123.1       |            | 56               |          |
| Prothioconazole-desthio | 9.10                 | 312.1         | 125.0       | 111        | 46               | +        |
|                         |                      | 312.1         | 70.0        |            | 26               |          |
| Pyraclostrobin          | 10.44                | 388.0         | 194.1       | 95         | 6                | +        |
|                         |                      | 388.0         | 163.1       |            | 20               |          |
| Pyrimethanil            | 7.57                 | 200.1         | 167.9       | 120        | 28               | +        |
|                         |                      | 200.1         | 107.1       |            | 20               |          |
| Quinoxifen              | 12.89                | 308.0         | 214.1       | 120        | 40               | +        |
|                         |                      | 308.0         | 197.0       |            | 36               |          |
| Saflufenacil            | 7.00                 | 501.1         | 459.0       | 160        | 8                | +        |
|                         |                      | 501.1         | 349.0       |            | 24               |          |
| Semiamitraz             | 3.14                 | 163.1         | 122.0       | 116        | 10               | +        |
|                         |                      | 163.1         | 107.0       |            | 18               |          |
| Spirodiclofen           | 14.10                | 411.1         | 313.0       | 110        | 6                | +        |
|                         |                      | 411.1         | 213.1       |            | 36               |          |
| Spirotetramat           | 8.69                 | 374.1         | 330.3       | 117        | 10               | +        |
|                         |                      | 374.1         | 302.2       |            | 14               |          |

| Compound Name            | LC/MS/MS MRM Setting |               |             |            |                  |          |
|--------------------------|----------------------|---------------|-------------|------------|------------------|----------|
|                          | RT (min)             | Precursor Ion | Product Ion | Fragmentor | Collision Energy | Polarity |
| Spirotetramat metabolite | 5.65                 | 302.2         | 270.1       | 131        | 22               | +        |
|                          |                      | 302.2         | 216.0       |            | 34               |          |
| Sulfoxaflor              | 3.81                 | 278.2         | 174.1       | 95         | 8                | +        |
|                          |                      | 278.2         | 154.1       |            | 32               |          |
| Tebufenozide             | 9.40                 | 353.0         | 297.2       | 95         | 5                | +        |
|                          |                      | 353.0         | 133.1       |            | 15               |          |
| Terbufos Sulfone         | 7.20                 | 321.0         | 171.0       | 80         | 5                | +        |
|                          |                      | 321.0         | 142.9       |            | 20               |          |
| Terbufos-sulfoxide       | 7.30                 | 305.1         | 186.9       | 50         | 10               | +        |
|                          |                      | 305.1         | 96.9        |            | 50               |          |
| Thiabendazole            | 3.68                 | 202.0         | 175.1       | 151        | 22               | +        |
|                          |                      | 202.0         | 131.1       |            | 30               |          |
| Thiacloprid              | 3.86                 | 253.0         | 126.0       | 100        | 20               | +        |
|                          |                      | 253.0         | 99.1        |            | 44               |          |
| Thiamethoxam             | 3.11                 | 292.0         | 211.1       | 85         | 10               | +        |
|                          |                      | 292.0         | 181.1       |            | 28               |          |
| Triadimefon              | 8.13                 | 294.1         | 225.0       | 111        | 2                | +        |
|                          |                      | 294.1         | 197.0       |            | 6                |          |
| Triadimenol              | 8.50                 | 298.1         | 70.1        | 70         | 10               | +        |
|                          |                      | 296.1         | 227.2       |            | 6                |          |
| Triforine                | 6.76                 | 434.9         | 389.9       | 110        | 12               | +        |
|                          |                      | 434.9         | 214.9       |            | 28               |          |

[www.agilent.com/chem](http://www.agilent.com/chem)

DE.4167708333

This information is subject to change without notice.

© Agilent Technologies, Inc. 2020  
Printed in the USA, June 24, 2020  
5994-2038EN