



Ultra low level quantification of pesticides in baby foods using an advanced triple quadrupole GC-MS/MS system

Authors

Richard Law, Aaron Lamb, Paul Silcock, and Cristian Cojocariu
Thermo Fisher Scientific,
Runcorn, UK

Keywords

TSQ 9000, advanced electron ionization, AEI, baby food, sensitivity, robustness, pesticide residue, QuEChERS, triple quadrupole mass spectrometry, GC-MS/MS, Programmable Temperature Vaporization, food safety

Goal

The aim of the study was to assess the quantitative performance of the Thermo Scientific™ TSQ™ 9000 triple quadrupole GC-MS/MS system fitted with the Advanced Electron Ionization (AEI) source for the analysis of pesticide residues at ultra low levels in baby food.

Introduction

The detection and subsequent quantification of pesticides, contaminants, and other chemical residues are of paramount importance, especially when the food stuff is intended to be consumed by infants or young children. The maximum residue level (MRL) for the majority of pesticide-commodity combinations is set at the default level of 10 µg/kg.¹⁻³ However, the European Union (EU) has established LOD MRLs between 3–8 µg/kg for specific pesticides prohibited in baby foods.⁴ These pesticides and their metabolites may cause infants and young children (under worst-case intake conditions) to exceed the acceptable daily intake (ADI) values. The high sensitivity and selectivity of GC-MS/MS enables the detection and identification of residues of prohibited compounds, in compliance with the residue definitions, even when dealing with the diverse composition of multi-ingredient baby foods.

Also, the increased levels of selectivity and sensitivity provided by triple quadrupole instruments compared to single quadrupole instruments enabled analysts to adopt faster, less specific sample extraction procedures such as QuEChERS (quick, easy, cheap effective, rugged and safe).

The QuEChERS procedure has become the standard approach for sample preparation in many laboratories because of improvement in productivity.⁵ The method usually involves extraction with acetonitrile in the presence of various salts followed by dispersive solid phase extraction (dSPE) clean-up with a combination of PSA, C18, and carbon sorbents. The efficiency of the dSPE clean-up is limited so high concentrations of matrix-coextractives can remain in the final extract and cause system contamination. Also, use of acetonitrile (which has a high coefficient of expansion) limits the injection volume and hence the sensitivity of the method.

Taking all of these considerations together, it is evident that an ultra-sensitive, selective, reliable, and robust GC-MS/MS system is needed to address the challenge of routine high-throughput determination of pesticide residues at trace concentrations in baby foods. In this study, the quantitative performance of the Thermo Scientific™ TSQ™ 9000 triple quadrupole GC-MS/MS system was assessed for the analysis of more than 200 pesticides in baby food at ultra low concentrations (as low as 0.025 µg/kg). A complete evaluation of method performance included sample preparation, overall method suitability measured from pesticides recoveries, selectivity, sensitivity, linearity, and long-term robustness.

Experimental

Sample preparation

Samples of carrot/potato and apple/pear/banana baby food samples were extracted using the citrate-buffered QuEChERS protocol using Thermo Scientific™ HyperSep™ dispersive solid phase extraction (dSPE) products.

Homogenized sample (10 g) was extracted with acetonitrile (10 mL) followed by the addition of MgSO₄ (4 g), NaCl (1.0 g), disodium hydrogen citrate sesquihydrate (0.5 g), and trisodium citrate dihydrate (1.0 g). Dispersive solid phase extraction (dSPE) [MgSO₄ (150 mg), PSA (25 mg) and GCB (25 mg) per 1 mL of extract for carrot/potato and MgSO₄ (150 mg) and PSA (25 mg) for apple/pear/banana] was used for sample clean-up.

Preparation of matrix-matched calibrations

Immediately after dSPE clean-up, the final extracts (1 g sample/mL of acetonitrile) were acidified with 5% formic acid in acetonitrile and were spiked with a mixture of 211 pesticides at 14 concentrations spanning a range of 0.025–250 µg/kg. Robustness was tested using repeat injections of samples (carrot/potato) spiked at the 10 µg/kg level.

For method evaluation, samples of carrot/potato and apple/pear/banana baby food samples were each spiked at 1.0, 2.5, and 10.0 µg/kg (n = 6 for each concentration) before extraction, clean-up, and acidification were carried out as described above.

GC-MS/MS analysis

A TSQ 9000 triple quadrupole GC-MS/MS system equipped with a Thermo Scientific™ Advanced Electron Ionization (AEI) source and coupled with a Thermo Scientific™ TRACE™ 1310 GC system was used. The AEI source provides a highly efficient electron ionization of analytes and a more tightly focused ion beam that provides an unparalleled level of sensitivity.

Liquid injections of the sample extracts were performed using a Thermo Scientific™ TriPlus™ RSH™ autosampler, and chromatographic separation was achieved by a Thermo Scientific™ TraceGOLD™ TG-5SilMS 30 m × 0.25 mm I.D. × 0.25 µm film capillary column with 5 m integrated SafeGuard. Additional details of instrument parameters are displayed in Table 1.

Table 1. Gas chromatograph and mass spectrometer parameters.

| TRACE 1310 GC System Parameters | | | | |
|---------------------------------------|---|---------------|-------------------|-----------------|
| Injection Volume (µL) | 1 | | | |
| Liner | Siltek™ six baffle PTV liner (P/N 453T2120) | | | |
| Inlet (°C) | 70 | | | |
| Carrier Gas, (mL/min) | He, 1.2 | | | |
| Inlet Mode | Splitless (split flow 50 mL/min after 2 min) | | | |
| Column | TraceGOLD TG-5SiIMS with SafeGuard (30 m × 0.25 mm, 0.25 µm with 5 m integrated guard column (P/N 26096-1425) | | | |
| PTV Parameters | Rate (°C/s) | Temp. (°C) | Time (min) | Flow (mL/min) |
| Injection | - | 70 | 0.10 | - |
| Transfer | 5.0 | 300 | 2.00 | - |
| Cleaning | 14.5 | 320 | 5.00 | 75.0 |
| Oven Temperature Program | | | | |
| Ramp | RT (min) | Rate (°C/min) | Target Temp. (°C) | Hold Time (min) |
| Initial | 0 | - | 40 | 1.50 |
| 1 | 1.5 | 25.0 | 90 | 1.50 |
| 2 | 5.0 | 25.0 | 180 | 0.00 |
| 3 | 8.6 | 5.0 | 280 | 0.00 |
| Final | 28.6 | 10.0 | 300 | 5.00 |
| Run time | 35.6 | - | - | - |
| TSQ 9000 Mass Spectrometer Parameters | | | | |
| Transfer Line (°C) | 250 | | | |
| Ionization Type | EI | | | |
| Ion Source (°C) | 320 | | | |
| Acquisition Mode | timed-SRM | | | |
| Tuning parameters | AEI SmartTune | | | |
| Collision gas and pressure (psi) | Argon at 70 | | | |
| Peak Width (Da) | 0.7 (both Q1 and Q3) | | | |

Data processing

Data were acquired, processed, and reported using Thermo Scientific™ Chromeleon™ Chromatography Data System (CDS) software, which allows instrument control, method development, quantitative/qualitative analysis, and customizable reporting all within one package.

Data review is highly customizable, allowing the user to display the information required on screen in real time. Furthermore, the flexibility of Chromeleon CDS software ensures that SANTE³ compliance criteria can easily be flagged, tracked, and reported to the user's individual requirements.

Results and discussion

Compliance with EU SANTE criteria

The method performance was tested in accordance to the SANTE/11813/2017 guidance document, which requires that the following criteria are satisfied for identification of pesticide residues:

- I. A minimum of two product ions are detected for each pesticide with peak S/N >3 (or, in case noise is absent, a signal should be present in at least five subsequent scans) and with the mass resolution for precursor-ion isolation equal to or better than unit mass resolution.
- II. Retention time tolerance of ± 0.1 minutes compared with standards in the same sequence.
- III. Ion ratio within $\pm 30\%$ (relative) of the average of calibration standards from the same sequence.

Wherever SANTE compliance is referenced in this study, all three criteria have been met fully.

Recoveries

Pesticide recoveries were obtained from the QuEChERS extractions performed on the samples spiked before extraction. All detected compounds, at the three spiking levels in both matrices satisfied all SANTE requirements. More than 97% of the target pesticide residues at 1 $\mu\text{g}/\text{kg}$ had recoveries between 70% and 120%. An example of the recovery and precision data for the apple/pear/banana matrix spiked at the default MRL (10 $\mu\text{g}/\text{kg}$) is displayed in Figure 1. A full table of results can be found in Appendix B.

Chromatography and selectivity

Analysis of a large number of pesticides in a single injection requires careful optimization of parameters, especially when injecting acetonitrile. As acetonitrile is a low molecular weight low polarity solvent, it has a relatively high expansion volume and is insoluble in the low polarity phases normally used for routine pesticide analysis (this makes solvent focusing in a standard splitless type injection incredibly difficult). These issues can be addressed by using an optimized programmable temperature vaporisation (PTV) injection. Figure 2 shows an example of three pesticides eluting in the beginning (A–dichlobenil – 0.025 $\mu\text{g}/\text{kg}$), middle (B–dieldrin 0.5 $\mu\text{g}/\text{kg}$), and end (C–deltamethrin 0.05 $\mu\text{g}/\text{kg}$) of the chromatographic run in the lowest detectable standard in carrot and potato matrix, levels at which all compounds detected meet the SANTE requirements. Peak shapes were Gaussian and coefficient of determination (R^2) was >0.990 for all three compounds indicating good chromatography and excellent linear response.

Identification of all 210 component peaks was made using an in-house, commercially available Thermo Scientific SRM pesticide compound database (cdb). In addition to this, retention time alignment of target compounds can be easily performed using the Thermo Scientific RTA tool,⁸ eliminating the need for manual correction of compound retention times whenever column maintenance is performed. The cdb database contains >1000 compounds with >3700 unique SRM transitions. Due to the fast scanning speed of the EvoCell technology and the intelligent scheduling of the timed-SRM,⁹ it is possible to acquire data with several transitions per compound with minimal loss in sensitivity. This makes it simple to select the most optimal transitions in differing matrices to perform quantitation and qualification on, removing the need to develop matrix-matched SRM compound databases (Figure 3).

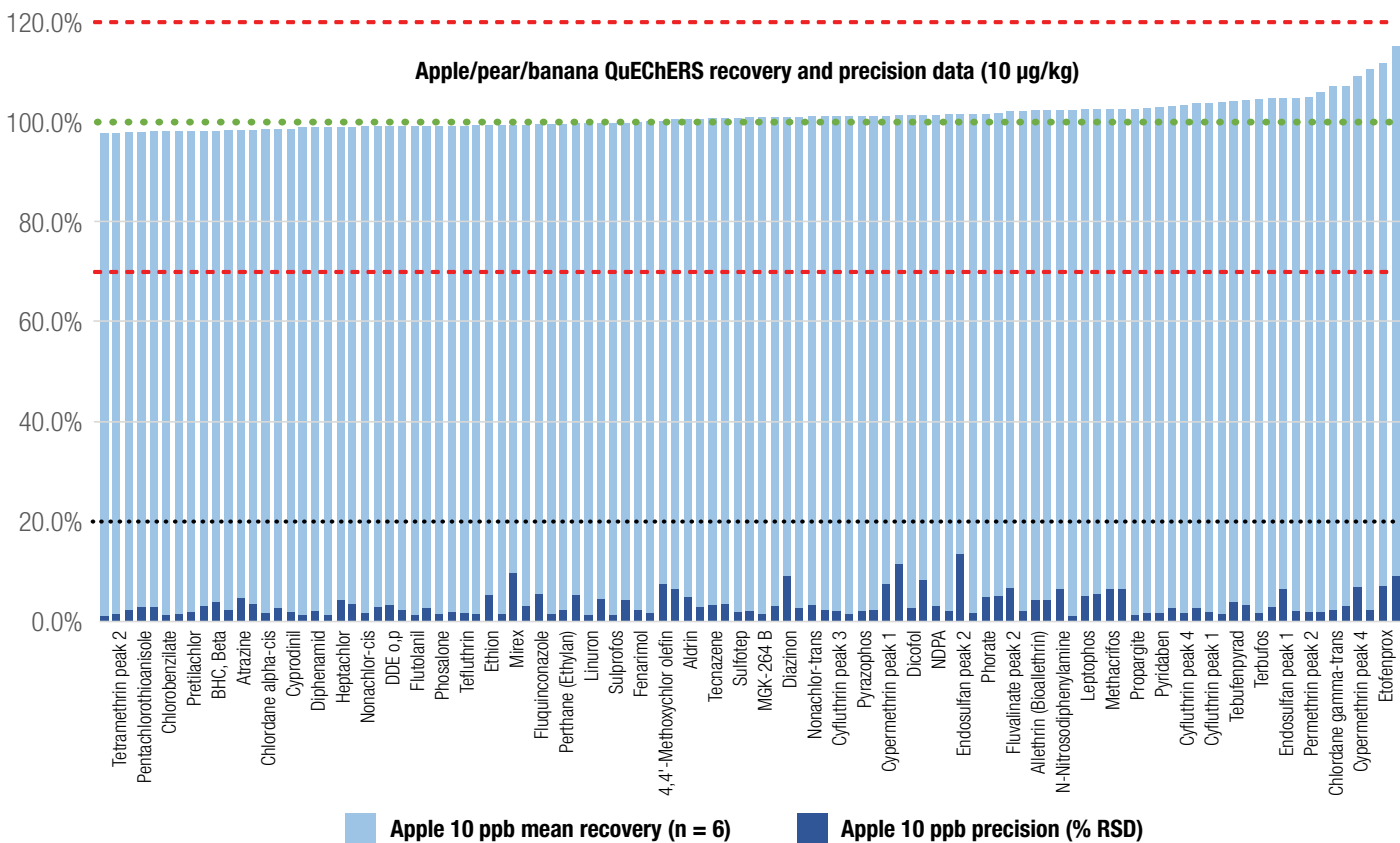
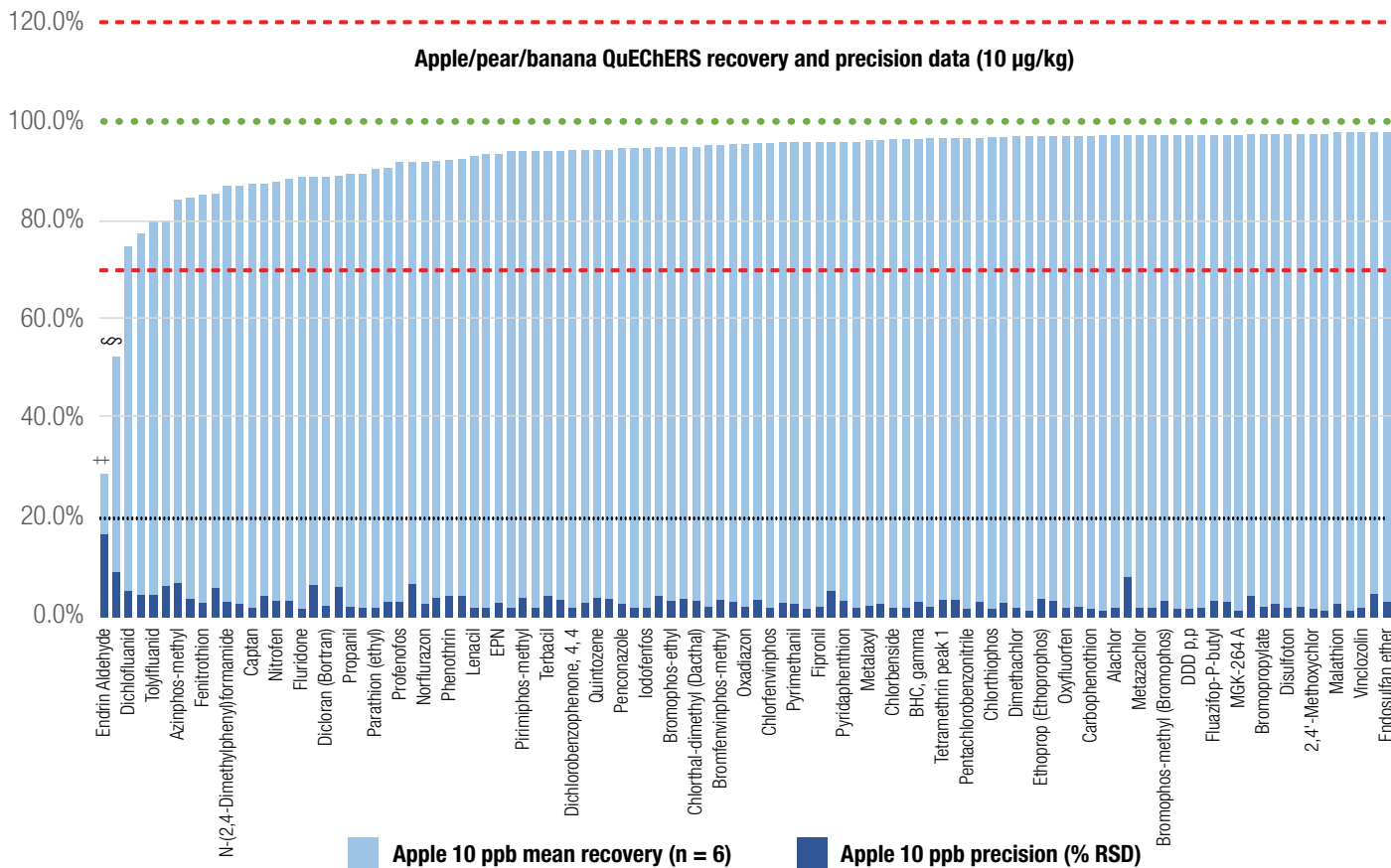


Figure 1. Recovery and precision data for apple/pear/banana extractions (n=6) at a concentration of 10 µg/kg. ‡ Endrin aldehyde recoveries were low, potentially due to reaction with PSA. § Recoveries of chlorothalonil, known to be problematic in QuEChERS extractions,⁷ were low.

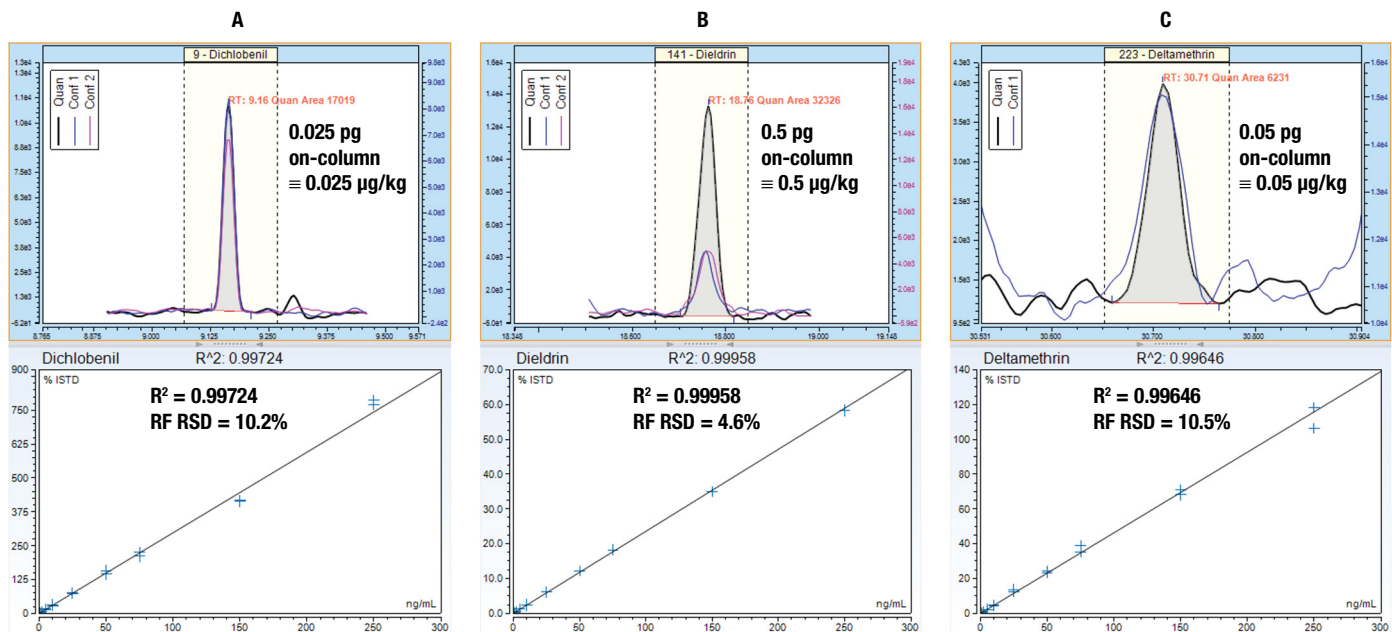


Figure 2. Example (A - Dichlobenil, B - dieldrin and C – deltamethrin) chromatographic peaks showing the lowest detectable matrix matched standard which meets SANTE requirements. The MRLs are 10 µg/kg, 3 µg/kg* and 10 µg/kg respectively. Calibration curves show duplicate injection at 14 discrete levels ranging from 0.025 pg to 250 pg on column. * Dieldrin is classed as a prohibited pesticide and 3 µg/kg considered to be the current limit of quantification, but is subject to regular review.⁴

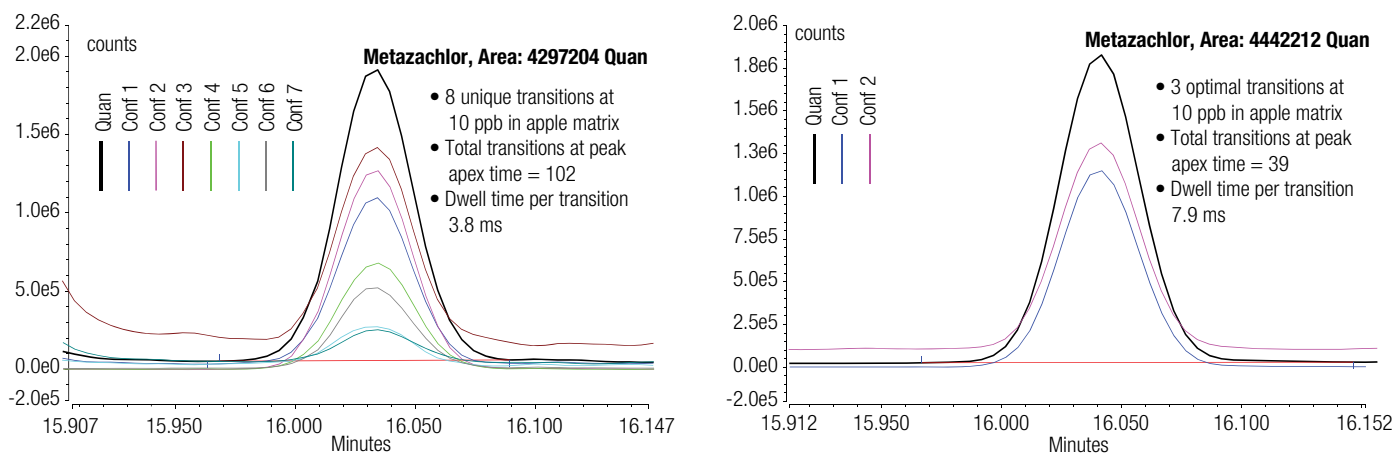


Figure 3. Comparison of Metazachlor SRM chromatographic peaks acquired using an injection containing 1317 unique transitions (left, 8 transitions) and an injection containing 663 (right, 3 transitions). No significant difference in the peak area for quantitation transition is observed indicating no loss in sensitivity.

Sensitivity and linearity

The TSQ 9000 AEI system easily met SANTE criteria (ion ratios $\pm 30\%$, etc.) at the default MRL of $10 \mu\text{g}/\text{kg}$ for all pesticides targeted. Moreover, over 90% of pesticides detected at $< 0.5 \mu\text{g}/\text{kg}$ meet the SANTE requirements and 10% of them meet SANTE criteria even at $0.025 \mu\text{g}/\text{kg}$ level (Figure 4). Resolution settings of 0.7 Daltons for Q1 and Q3 were used, ensuring the optimum combination of selectivity and sensitivity.

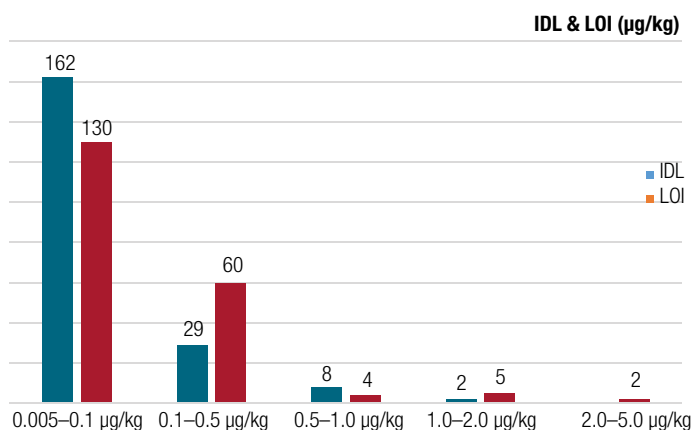


Figure 4. Number of target compounds satisfying the SANTE requirements, with over 90% below $0.5 \mu\text{g}/\text{kg}$, and over 60% below $0.1 \mu\text{g}/\text{kg}$ – 100 times lower than the default MRL [sample matrix – carrot/potato].

Over 90% of the target compounds had a Limit of Identification (LOI) (satisfying all SANTE requirements) below $0.5 \mu\text{g}/\text{kg}$, and over 60% below $0.1 \mu\text{g}/\text{kg}$.

System sensitivity, defined as instrumental detection limits (IDLs), was determined experimentally for each compound by performing $n=10$ replicate injections of the lowest matrix-matched standard of carrot and potato that met all SANTE criteria. Calculations of IDLs were then made using one-tailed student t -test at the 99% confidence interval for the corresponding degrees of freedom and taking into account the concentration and absolute peak area %RSD for each compound (Figures 5 and 6).

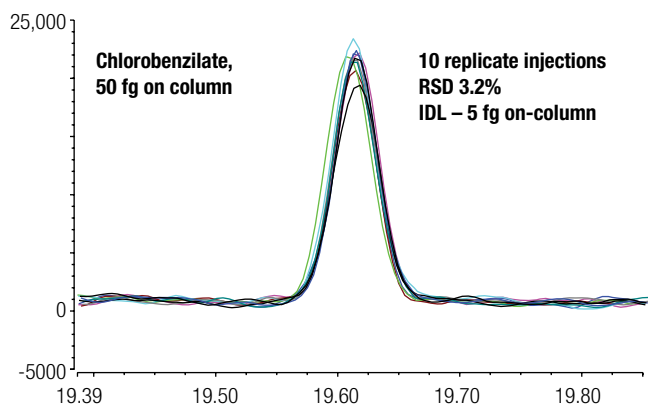
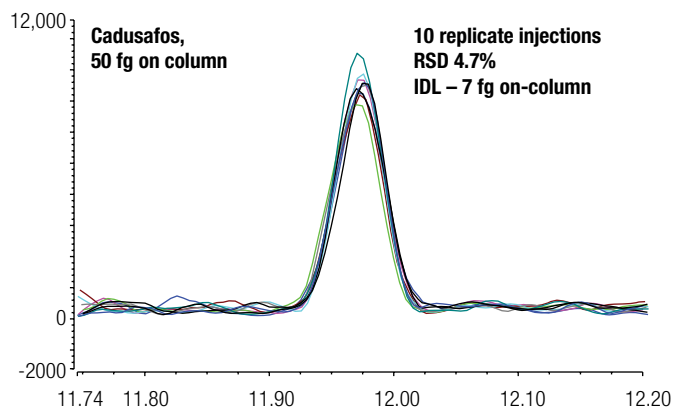


Figure 5. Example quantification SRM overlays of cadusafos and chlorobenzilate injected at the lowest level that met all SANTE criteria. Annotated are on column concentration, %RSD derived from absolute peak area response and calculated IDLs.

Fipronil and fipronil-desulfinyl, expressed as fipronil, have a multi-component MRL specified at $4 \mu\text{g}/\text{kg}$. Therefore, to satisfy the current regulations, each component must be identified at $2 \mu\text{g}/\text{kg}$. Figure 7 shows fipronil and fipronil-desulfinyl at concentrations of $0.2 \mu\text{g}/\text{kg}$, ten times lower than the requisite MRL, with back-calculated concentrations versus the linear calibration annotated.

Compound linearity was assessed by injecting matrix-matched standards in the range of 0.025 to $250 \mu\text{g}/\text{kg}$ in duplicate for both carrot/potato and apple/pear/banana. Both sets of linearity data showed $R^2 > 0.990$ and response factor (RF) % RSDs of $< 20\%$ for over 96% of component peaks indicating excellent linear response. Examples of linearity are shown in Figure 2 and in a comprehensive table provided in Appendix A.



Figure 6. Plot showing the calculated IDLs for all pesticides. IDLs ranged from ~5 fg (chlorobenzilate) to ~2.0 pg (bioallethrin) with >95% of compounds showing an IDL of less than 500 fg on column (equivalent to 0.5 µg/kg in sample extract). See Appendix A for tabulated data.

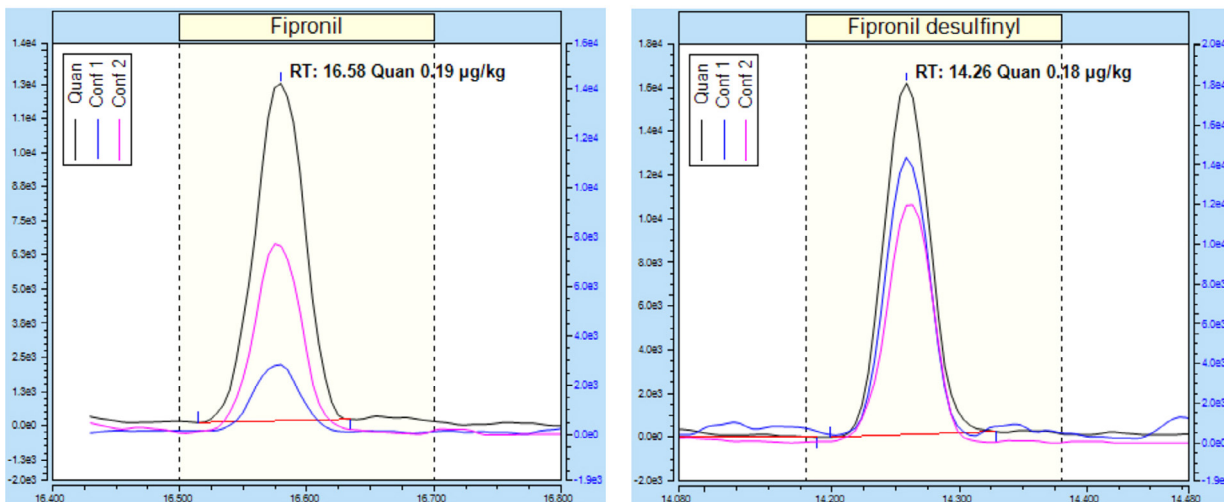


Figure 7. Fipronil and fipronil desulfinyl, at a concentration of 0.2 µg/kg equating to 0.4 µg/kg fipronil (sum), with SANTE compliance throughout.

AEI source robustness

The TSQ 9000 AEI system was set up as described in Table 1. After an initial source cleaning, repeat injections of a QuEChERS sample extract (1 g/mL carrot and potato) spiked at the default MRL (10 µg/kg) were made (Figure 8). Extracts resulting from the QuEChERS methodology contain many undesirable matrix co-extracted components which can easily contaminate the GC inlet, the chromatographic column and the MS ion source. To test the robustness of the AEI ion source

only (as far as reasonably practicable), after every 100 sample injections, the PTV liner was replaced along with the injector septum, approximately 10 cm was trimmed from the head of the guard column followed by automatic tuning of the system using the SmartTune feature. SmartTune uses the MS parameters established during the initial tuning on a clean source and intelligently assess the performance of the system, only re-tuning when MS performance has been compromised. No additional maintenance was performed.

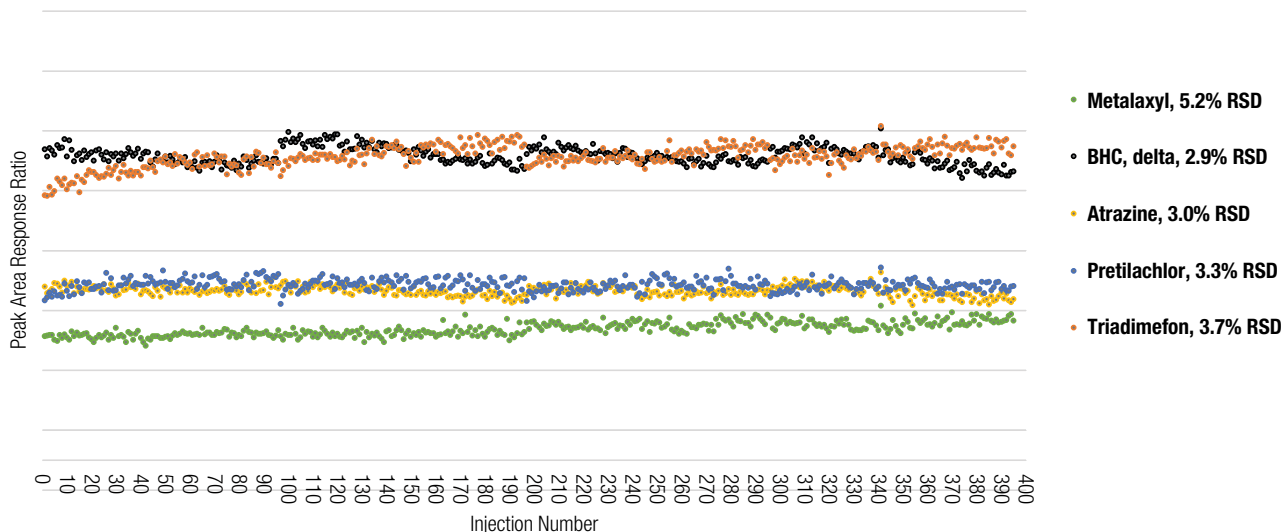


Figure 8A. Example of normalised peak area response for selected compounds across ~400 consecutive injections at the default MRL (10 µg/kg) in carrot/potato matrix.

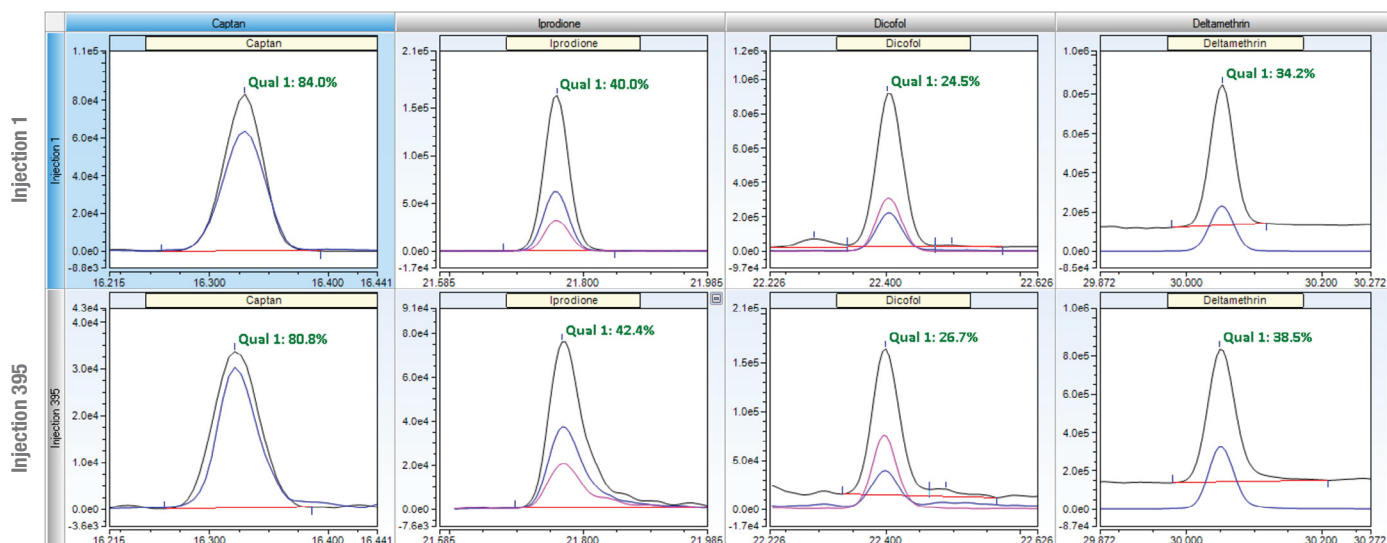


Figure 8B. Peak shapes, intensities and ion ratios of the primary qualifier ion for injection 1 (top row) and injection 395 (bottom row) for captan, iprodione, dicofol and deltamethrin.

Ion ratios at the default MRL were stable, Figure 9 shows pretilachlor ion qualifier ratios 1 and 2 in the first and last batches of injections. Ratios were well within the $\pm 30\%$ SANTE identification criteria.

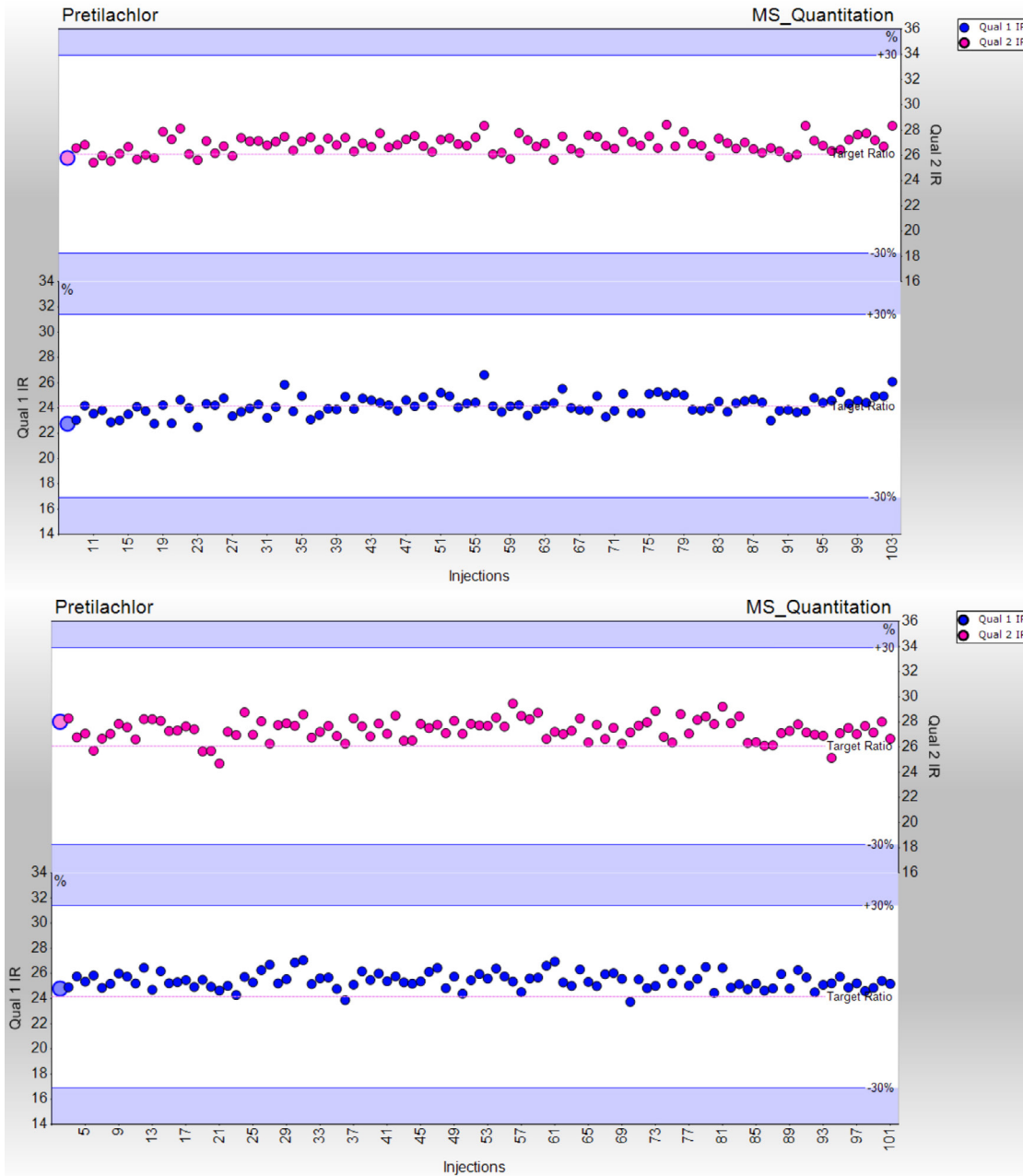


Figure 9. Pretilachlor Ion ratios of robustness injections 1-95 (top) and 295-395 (bottom).

Conclusions

In this work it has been demonstrated that by using QuEChERS with Thermo Scientific HyperSep dSPE products and a direct injection of acetonitrile extracts, the TSQ 9000 AEI system delivers outstanding quantitative performance for low-level pesticide residue analysis in baby food.

- QuEChERS extraction and subsequent clean-up of over 200 pesticides from replicate analysis (n=6 each at three concentrations) of each of two sample matrices, demonstrating excellent accuracy (recovery) and precision.
- Accurate, quantitative analysis of over 200 pesticides over up to five orders of magnitude (0.025–250 µg/kg), showing outstanding LODs and linear response.
- Robustness displayed over approximately 400 consecutive injections of sample matrix (1 g/mL), with SANTE compliance at the default MRL throughout.
- High sensitivity providing the real possibility to dilute the sample extract, thus limiting matrix contamination and system maintenance, leading to an increase in laboratory productivity.

The results of this study establish the TSQ 9000 triple quadrupole GC-MS/MS system, in combination with Chromeleon CDS software and HyperSep dSPE products, as the ideal solution for the routine analysis of pesticides in baby food, providing unprecedented sensitivity, robustness, ease of use, cost effectiveness, and reliability.

References

1. https://ec.europa.eu/food/plant/pesticides/max_residue_levels/guidelines_en
2. <http://eur-lex.europa.eu/legal-content/EN/TXT/?uri=URISERV:113002a>
3. SANTE/11813/2017. Guidance document on analytical quality control and method validation procedures for pesticides residues analysis in food and feed. Supersedes SANTE/11945/2015. Implemented by 01/01/2018. https://ec.europa.eu/food/sites/food/files/plant/docs/pesticides_mrl_guidelines_wrkdoc_2017-11813.pdf
4. <http://eur-lex.europa.eu/legal-content/EN/ALL/?uri=CELEX%3A32006L0125>
5. <http://quechers.cvua-stuttgart.de/>
6. <https://www.ncbi.nlm.nih.gov/pmc/articles/PMC4890451/>
7. Belmonte Valles, N.; Retamal, M.; Martinez-Uroz, M.A.; Mezcuca, M.; Fernandez-Alba, A.R.; de Kok, A. Determination of chlorothalonil in difficult-to-analyse vegetable matrices using various multiresidue methods. *Analyst*, 2012, 137(10), 2513–2520.
8. <https://assets.thermofisher.com/TFS-Assets/CMD/Application-Notes/an-10580-gc-ms-ms-ptv-pesticides-water-an10580-en.pdf>
9. <http://tools.thermofisher.com/content/sfs/posters/PN-10394-Fast-SRM-Transition-PN10394-EN.pdf>

Appendix A – Linearity data sets

Appendix A (Part 1). Linearity data sets.

| Chart Number | Compound Name | Apple/Pear/Banana Linearity | | | Carrot/Potato Linearity | | | IDL ₉₉ | |
|--------------|------------------------------|-----------------------------|-----------|-------------|-------------------------|-----------|-------------|-------------------|----------|
| | | R ² | RF RSD(%) | Range (ppb) | R ² | RF RSD(%) | Range (ppb) | pg on Column | IDL (fg) |
| 1 | 2,3,5,6-Tetrachloroaniline | 0.99931 | 6.80 | 0.1–250 | 0.99955 | 7.70 | 0.05–250 | 0.05 | 20 |
| 2 | 2,4'-Methoxychlor | 0.99987 | 3.50 | 0.1–250 | 0.99950 | 7.30 | 0.025–250 | 0.05 | 8 |
| 3 | 4,4'-Methoxychlor olefin | 0.99976 | 5.80 | 0.1–250 | 0.99932 | 4.20 | 0.05–250 | 0.05 | 31 |
| 4 | Acetochlor | 0.99972 | 3.60 | 0.2–250 | 0.99962 | 3.60 | 0.2–250 | 0.20 | 201 |
| 5 | Acrinathrin | 0.99963 | 2.70 | 0.2–250 | 0.99955 | 3.90 | 0.2–250 | 0.20 | 57 |
| 6 | Alachlor | 0.99965 | 5.50 | 0.1–250 | 0.99955 | 5.10 | 0.2–250 | 0.20 | 71 |
| 7 | Aldrin | 0.99959 | 8.00 | 0.1–250 | 0.99983 | 4.60 | 0.1–250 | 0.20 | 75 |
| 8 | Allethrin (Bioallethrin) | 0.99826 | 19.40 | 10–250 | 0.99888 | 7.50 | 5–250 | 5.00 | 2007 |
| 9 | Allidochlor | 0.99926 | 15.20 | 0.1–250 | 0.99631 | 6.30 | 0.2–250 | 0.20 | 145 |
| 10 | Anthraquinone | 0.99966 | 9.80 | 0.2–250 | 0.99988 | 17.60 | 0.025–250 | 0.05 | 27 |
| 11 | Atrazine | 0.99963 | 7.40 | 0.1–250 | 0.99990 | 6.70 | 0.05–250 | 0.05 | 19 |
| 12 | Azinphos-ethyl | 0.99465 | 9.40 | 0.2–250 | 0.99935 | 4.50 | 0.2–250 | 0.20 | 82 |
| 13 | Azinphos-methyl | 0.98165 | 16.90 | 1–250 | 0.99758 | 19.60 | 0.5–250 | 1.00 | 521 |
| 14 | BHC, Alpha | 0.99981 | 5.30 | 0.025–250 | 0.99949 | 6.60 | 0.025–250 | 0.05 | 15 |
| 15 | BHC, Beta | 0.99967 | 6.20 | 0.05–250 | 0.99985 | 8.80 | 0.025–250 | 0.05 | 15 |
| 16 | BHC, delta | 0.99971 | 4.10 | 0.05–250 | 0.99992 | 7.60 | 0.025–250 | 0.05 | 20 |
| 17 | BHC, gamma | 0.99970 | 7.30 | 0.05–250 | 0.99971 | 6.70 | 0.05–250 | 0.05 | 31 |
| 18 | Bifenthrin | 0.99989 | 4.20 | 0.5–250 | 0.99976 | 2.40 | 0.5–250 | 1.00 | 42 |
| 19 | Biphenyl | 0.99822 | 19.50 | 2–250 | 0.99573 | 14.50 | 5–250 | 5.00 | 865 |
| 20 | Bromfenvinphos | 0.99963 | 5.10 | 0.05–250 | 0.99960 | 7.60 | 0.025–250 | 0.05 | 31 |
| 21 | Bromfenvinphos-methyl | 0.99917 | 3.20 | 0.5–250 | 0.99971 | 3.60 | 0.1–250 | 0.20 | 33 |
| 22 | Bromophos-ethyl | 0.99946 | 3.30 | 0.1–250 | 0.99523 | 5.90 | 0.05–250 | 0.05 | 12 |
| 23 | Bromophos-methyl (Bromophos) | 0.99957 | 5.60 | 0.05–250 | 0.99848 | 5.90 | 0.05–250 | 0.05 | 24 |
| 24 | Bromopropylate | 0.99960 | 4.80 | 0.1–250 | 0.99806 | 5.40 | 0.1–250 | 0.20 | 61 |
| 25 | Bupirimate | 0.99947 | 10.10 | 0.05–250 | 0.99961 | 5.50 | 0.05–250 | 0.05 | 33 |
| 26 | Cadusafos | 0.99982 | 3.80 | 0.1–250 | 0.99952 | 6.40 | 0.025–250 | 0.05 | 7 |
| 27 | Captan | 0.98233 | 23.80 | 1–250 | 0.98303 | 16.60 | 0.5–250 | 1.00 | 733 |
| 28 | Carbophenothion | 0.99968 | 3.50 | 0.2–250 | 0.99970 | 4.40 | 0.1–250 | 0.20 | 30 |
| 29 | Carfentrazone-ethyl | 0.99929 | 6.10 | 0.2–250 | 0.99575 | 7.50 | 0.1–250 | 0.20 | 41 |
| 30 | Chlorbenside | 0.99981 | 5.50 | 0.025–250 | 0.99984 | 3.20 | 0.025–250 | 0.05 | 11 |
| 31 | Chlordane alpha-cis | 0.99875 | 8.10 | 0.05–250 | 0.98923 | 8.10 | 0.1–250 | 0.20 | 61 |
| 32 | Chlordane gamma-trans | 0.99949 | 6.50 | 0.05–250 | 0.99956 | 7.30 | 0.025–250 | 0.05 | 38 |
| 33 | Chlorfenapyr | 0.99979 | 6.10 | 0.2–250 | 0.99983 | 3.30 | 0.2–250 | 0.20 | 90 |
| 34 | Chlorfenson | 0.99986 | 3.60 | 0.025–250 | 0.99979 | 2.30 | 0.025–250 | 0.05 | 10 |
| 35 | Chlorfenvinphos | 0.99966 | 8.60 | 0.05–250 | 0.99987 | 5.70 | 0.025–250 | 0.05 | 16 |
| 36 | Chlorobenzilate | 0.99990 | 3.10 | 0.025–250 | 0.99976 | 3.70 | 0.025–250 | 0.05 | 5 |
| 37 | Chloroneb | 0.99962 | 4.70 | 0.5–250 | 0.99907 | 10.90 | 0.1–250 | 0.20 | 28 |
| 38 | Chlorothalonil | 0.99752 | 7.40 | 0.1–250 | 0.99635 | 18.40 | 0.05–250 | 0.05 | 24 |
| 39 | Chlorpropham | 0.99985 | 13.90 | 0.5–250 | 0.99981 | 12.60 | 2–250 | 5.00 | 166 |
| 40 | Chlorpyrifos-ethyl | 0.99948 | 15.00 | 0.1–250 | 0.99916 | 5.50 | 0.05–250 | 0.05 | 22 |

Appendix A (Part 2). Linearity data sets.

| Chart Number | Compound Name | Apple/Pear/Banana Linearity | | | Carrot/Potato Linearity | | | IDL ₉₉ | |
|--------------|-----------------------------|-----------------------------|-----------|-------------|-------------------------|-----------|-------------|-------------------|----------|
| | | R ² | RF RSD(%) | Range (ppb) | R ² | RF RSD(%) | Range (ppb) | pg on Column | IDL (fg) |
| 41 | Chlorpyrifos-methyl | 0.99969 | 12.30 | 0.1–250 | 0.99960 | 15.70 | 0.025–250 | 0.05 | 24 |
| 42 | Chlorthal-dimethyl | 0.99917 | 6.20 | 0.1–250 | 0.99912 | 5.80 | 0.025–250 | 0.05 | 20 |
| 43 | Chlorthiophos | 0.99976 | 7.20 | 0.1–250 | 0.99971 | 3.00 | 0.2–250 | 0.20 | 153 |
| 44 | Chlozolinate | 0.99981 | 7.90 | 0.2–250 | 0.99988 | 4.30 | 0.2–250 | 0.20 | 67 |
| 45 | Clomazone | 0.99983 | 4.30 | 0.025–250 | 0.99993 | 6.40 | 0.025–250 | 0.05 | 12 |
| 46 | Coumaphos | 0.99817 | 4.80 | 0.5–250 | 0.99966 | 8.50 | 0.2–250 | 0.20 | 78 |
| 47 | Cycloate | 0.99815 | 13.10 | 2–250 | 0.99890 | 4.80 | 2–250 | 5.00 | 789 |
| 48 | Cyhalothrin I (lambda) | 0.99966 | 7.50 | 0.2–250 | 0.99959 | 2.90 | 0.2–250 | 0.20 | 28 |
| 49 | Cyprodinil | 0.99983 | 5.00 | 0.1–250 | 0.99993 | 3.90 | 0.1–250 | 0.20 | 86 |
| 50 | DDD p,p | 0.99985 | 4.30 | 0.05–250 | 0.99992 | 4.00 | 0.025–250 | 0.05 | 22 |
| 51 | DDD, o, p | 0.99988 | 3.00 | 0.05–250 | 0.99985 | 7.50 | 0.05–250 | 0.05 | 20 |
| 52 | DDE o,p | 0.99974 | 3.80 | 0.025–250 | 0.99976 | 4.30 | 0.025–250 | 0.05 | 8 |
| 53 | DDE p, p | 0.99957 | 2.60 | 0.05–250 | 0.99989 | 4.20 | 0.05–250 | 0.05 | 9 |
| 54 | DDT o,p | 0.99988 | 5.70 | 0.05–250 | 0.99960 | 9.20 | 0.05–250 | 0.05 | 25 |
| 55 | DDT p,p | 0.99962 | 8.40 | 0.2–250 | 0.99937 | 14.40 | 0.1–250 | 0.20 | 41 |
| 56 | Deltamethrin | 0.99983 | 6.70 | 0.05–250 | 0.99646 | 11.00 | 0.05–250 | 0.05 | 22 |
| 57 | Diazinon | 0.99949 | 5.20 | 0.1–250 | 0.99906 | 5.00 | 0.1–250 | 0.20 | 33 |
| 58 | Dichlobenil | 0.99866 | 5.10 | 0.025–250 | 0.99724 | 10.20 | 0.025–250 | 0.05 | 5 |
| 59 | Dichlofuanid | 0.99949 | 6.40 | 0.2–250 | 0.99966 | 4.40 | 0.1–250 | 0.20 | 28 |
| 60 | Dichlorobenzophenone, 4, 4* | 0.99979 | 2.80 | 0.05–250 | 0.99957 | 7.50 | 0.025–250 | 0.05 | 17 |
| 61 | Dicloran (Bortran) | 0.99908 | 7.10 | 0.2–250 | 0.99801 | 7.00 | 0.2–250 | 0.20 | 58 |
| 62 | Dicofol* | 0.99272 | 13.50 | 0.5–250 | 0.98598 | 14.70 | 0.5–250 | 1.00 | 973 |
| 63 | Dieldrin | 0.99909 | 7.20 | 0.5–250 | 0.99958 | 4.60 | 0.5–250 | 1.00 | 162 |
| 64 | Dimethachlor | 0.99964 | 5.10 | 0.025–250 | 0.99968 | 4.60 | 0.025–250 | 0.05 | 10 |
| 65 | Dimethoate | 0.99903 | 6.30 | 0.2–250 | 0.99973 | 10.40 | 0.1–250 | 0.20 | 30 |
| 66 | Diphenamid | 0.99974 | 7.30 | 0.2–250 | 0.99974 | 5.20 | 0.2–250 | 0.20 | 62 |
| 67 | Diphenylamine | 0.99981 | 9.50 | 0.2–250 | 0.99931 | 17.00 | 0.1–250 | 0.20 | 28 |
| 68 | Disulfoton | 0.99982 | 6.80 | 0.2–250 | 0.99943 | 4.30 | 0.2–250 | 0.20 | 19 |
| 69 | Edifenphos | 0.99908 | 4.00 | 0.05–250 | 0.99967 | 11.50 | 0.025–250 | 0.05 | 10 |
| 70 | Endosulfan ether | 0.99983 | 5.70 | 0.05–250 | 0.99982 | 12.20 | 0.025–250 | 0.05 | 21 |
| 71 | Endosulfan peak 1 | 0.99963 | 4.50 | 0.2–250 | 0.99989 | 4.50 | 0.2–250 | 0.20 | 42 |
| 72 | Endosulfan peak 2 | 0.99982 | 5.20 | 0.5–250 | 0.99988 | 4.00 | 0.5–250 | 1.00 | 190 |
| 73 | Endosulfan sulfate | 0.99981 | 3.10 | 0.1–250 | 0.99980 | 3.60 | 0.05–250 | 0.05 | 20 |
| 74 | Endrin | 0.99981 | 3.70 | 0.5–250 | 0.99975 | 4.80 | 0.2–250 | 0.20 | 59 |
| 75 | Endrin Aldehyde | 0.99893 | 7.30 | 0.5–250 | 0.99786 | 9.30 | 0.5–250 | 1.00 | 209 |
| 76 | Endrin-Ketone | 0.99920 | 6.00 | 0.5–250 | 0.99872 | 5.80 | 0.5–250 | 1.00 | 353 |
| 77 | EPN | 0.99591 | 7.10 | 1–250 | 0.99334 | 14.00 | 1–250 | 1.00 | 302 |
| 78 | Ethion | 0.99981 | 3.60 | 0.1–250 | 0.99987 | 3.90 | 0.05–250 | 0.05 | 29 |
| 79 | Ethoprop (Ethoprophos) | 0.99975 | 6.10 | 0.05–250 | 0.99923 | 6.40 | 0.1–250 | 0.20 | 39 |
| 80 | Etofenprox | 0.99978 | 7.40 | 0.2–250 | 0.99992 | 3.70 | 0.2–250 | 0.20 | 42 |

* - 4,4-dichlorobenzophenone is a breakdown product of dicofol and therefore may be overestimated.

Appendix A (Part 3). Linearity data sets.

| Chart Number | Compound Name | Apple/Pear/Banana Linearity | | | Carrot/Potato Linearity | | | IDL ₉₉ | |
|--------------|-------------------------|-----------------------------|-----------|-------------|-------------------------|-----------|-------------|-------------------|----------|
| | | R ² | RF RSD(%) | Range (ppb) | R ² | RF RSD(%) | Range (ppb) | pg on Column | IDL (fg) |
| 81 | Etridiazole (Terrazole) | 0.99954 | 5.50 | 0.1–250 | 0.99599 | 16.30 | 0.025–250 | 0.05 | 18 |
| 82 | Fenamiphos | 0.99958 | 3.00 | 0.5–250 | 0.99848 | 3.70 | 0.5–250 | 1.00 | 147 |
| 83 | Fenarimol | 0.99984 | 3.20 | 0.2–250 | 0.99990 | 2.60 | 0.2–250 | 0.20 | 45 |
| 84 | Fenclorfos | 0.99964 | 3.90 | 0.05–250 | 0.99956 | 4.80 | 0.05–250 | 0.05 | 17 |
| 85 | Fenitrothion | 0.99241 | 13.50 | 0.1–250 | 0.99559 | 13.30 | 0.1–250 | 0.20 | 52 |
| 86 | Fenpropathrin | 0.99969 | 3.70 | 0.5–250 | 0.99976 | 5.00 | 1–250 | 1.00 | 147 |
| 87 | Fenson | 0.99989 | 4.40 | 0.05–250 | 0.99995 | 5.60 | 0.025–250 | 0.05 | 11 |
| 88 | Fenthion | 0.99959 | 10.20 | 0.05–250 | 0.99970 | 10.60 | 0.05–250 | 0.05 | 18 |
| 89 | Fenvalerate | 0.99992 | 3.10 | 0.1–250 | 0.99974 | 5.80 | 0.1–250 | 0.20 | 36 |
| 90 | Fipronil | 0.99923 | 5.10 | 0.1–250 | 0.99405 | 8.90 | 0.05–250 | 0.05 | 20 |
| 91 | Fipronil desulfinyl | 0.99826 | 5.60 | 0.05–250 | 0.98489 | 10.00 | 0.05–250 | 0.05 | 27 |
| 92 | Fluazifop-P-butyl | 0.99971 | 7.40 | 0.1–250 | 0.99976 | 10.50 | 0.1–250 | 0.20 | 45 |
| 93 | Fludioxonil | 0.99980 | 9.90 | 0.05–250 | 0.99951 | 8.00 | 0.025–250 | 0.05 | 29 |
| 94 | Fluquinconazole | 0.99953 | 7.70 | 0.05–250 | 0.99609 | 7.40 | 0.025–250 | 0.05 | 15 |
| 95 | Fluridone | 0.99890 | 13.60 | 0.2–250 | 0.99593 | 7.10 | 0.025–250 | 0.05 | 22 |
| 96 | Flusilazole | 0.99969 | 11.20 | 0.1–250 | 0.99982 | 9.00 | 0.1–250 | 0.20 | 66 |
| 97 | Flutolanil | 0.99966 | 5.80 | 0.05–250 | 0.99982 | 9.30 | 0.025–250 | 0.05 | 51 |
| 98 | Flutriafol | 0.99960 | 7.20 | 0.1–250 | 0.99989 | 9.50 | 0.025–250 | 0.05 | 24 |
| 99 | Folpet | 0.97866 | 23.90 | 0.5–250 | 0.98874 | 14.20 | 0.2–250 | 0.20 | 757 |
| 100 | Fonofos | 0.99970 | 4.80 | 0.05–250 | 0.99986 | 4.20 | 0.05–250 | 0.05 | 14 |
| 101 | Heptachlor | 0.99963 | 4.20 | 0.025–250 | 0.99976 | 6.40 | 0.025–250 | 0.05 | 7 |
| 102 | Hexachlorobenzene | 0.99862 | 7.80 | 0.025–250 | 0.99939 | 5.60 | 0.025–250 | 0.05 | 11 |
| 103 | Hexazinone | 0.99971 | 7.20 | 0.1–250 | 0.99983 | 5.50 | 0.05–250 | 0.05 | 18 |
| 104 | Iodofenfos | 0.99859 | 6.10 | 0.05–250 | 0.99012 | 11.60 | 0.05–250 | 0.05 | 19 |
| 105 | Iprodione | 0.99976 | 7.70 | 0.2–250 | 0.99536 | 20.80 | 0.1–250 | 0.20 | 80 |
| 106 | Isazophos | 0.99934 | 6.60 | 0.1–250 | 0.99945 | 12.10 | 0.1–250 | 0.20 | 46 |
| 107 | Isodrin | 0.99983 | 5.90 | 0.1–250 | 0.99992 | 6.20 | 0.1–250 | 0.20 | 26 |
| 108 | Lenacil | 0.99928 | 7.90 | 0.2–250 | 0.99971 | 5.30 | 0.1–250 | 0.20 | 83 |
| 109 | Leptophos | 0.99947 | 3.30 | 0.2–250 | 0.99909 | 3.80 | 0.2–250 | 0.20 | 36 |
| 110 | Linuron | 0.99913 | 8.00 | 0.5–250 | 0.99831 | 8.70 | 0.2–250 | 0.20 | 92 |
| 111 | Malathion | 0.99989 | 5.70 | 0.05–250 | 0.99972 | 6.10 | 0.025–250 | 0.05 | 12 |
| 112 | Metalaxyl | 0.99947 | 4.70 | 0.2–250 | 0.99985 | 21.80 | 0.1–250 | 0.20 | 54 |
| 113 | Metazachlor | 0.99958 | 3.80 | 0.1–250 | 0.99978 | 8.10 | 0.025–250 | 0.05 | 32 |
| 114 | Methacrifos | 0.99977 | 4.50 | 0.2–250 | 0.99819 | 4.80 | 0.2–250 | 0.20 | 91 |
| 115 | Methoxychlor | 0.99918 | 4.50 | 0.1–250 | 0.99921 | 5.40 | 0.1–250 | 0.20 | 38 |
| 116 | Metolachlor | 0.99978 | 4.20 | 0.05–250 | 0.99992 | 4.20 | 0.025–250 | 0.05 | 49 |
| 117 | Mevinphos | 0.99985 | 3.80 | 0.05–250 | 0.99937 | 4.60 | 0.1–250 | 0.20 | 31 |
| 118 | MGK-264 A | 0.99986 | 5.00 | 0.2–250 | 0.99966 | 4.00 | 0.2–250 | 0.20 | 51 |
| 119 | MGK-264 B | 0.99984 | 4.50 | 0.2–250 | 0.99974 | 4.40 | 0.2–250 | 0.20 | 65 |
| 120 | Mirex | 0.99980 | 4.60 | 0.025–250 | 0.99981 | 3.00 | 0.025–250 | 0.05 | 8 |

Appendix A (Part 4). Linearity data sets.

| Chart Number | Compound Name | Apple/Pear/Banana Linearity | | | Carrot/Potato Linearity | | | IDL ₉₉ | |
|--------------|----------------------------------|-----------------------------|-----------|-------------|-------------------------|-----------|-------------|-------------------|----------|
| | | R ² | RF RSD(%) | Range (ppb) | R ² | RF RSD(%) | Range (ppb) | pg on Column | IDL (fg) |
| 121 | Myclobutanil | 0.99978 | 1.90 | 0.1–250 | 0.99986 | 3.50 | 0.1–250 | 0.20 | 40 |
| 122 | N-(2,4-Dimethylphenyl) formamide | 0.99953 | 4.80 | 1–250 | 0.99982 | 8.30 | 1–250 | 1.00 | 106 |
| 123 | NDBA | 0.99866 | 18.00 | 0.5–250 | 0.99414 | 21.60 | 0.05–250 | 0.05 | 12 |
| 124 | NDEA | 0.99826 | 6.90 | 0.1–250 | 0.98989 | 9.80 | 0.2–250 | 0.20 | 74 |
| 125 | NDPA | 0.99865 | 8.70 | 0.5–250 | 0.99133 | 8.80 | 0.1–250 | 0.20 | 49 |
| 126 | NEMA | 0.99657 | 8.00 | 1–250 | 0.98500 | 12.80 | 0.2–250 | 0.20 | 87 |
| 127 | Nitrofen | 0.99590 | 13.50 | 0.05–250 | 0.99512 | 15.00 | 0.2–250 | 0.20 | 30 |
| 128 | N-Nitrosodiphenylamine | 0.99971 | 12.70 | 0.2–250 | 0.99931 | 17.00 | 0.1–250 | 0.20 | 28 |
| 129 | N-Nitrosomorpholine | 0.99759 | 18.50 | 0.5–250 | 0.99342 | 20.00 | 0.5–250 | 1.00 | 198 |
| 130 | N-Nitrosopiperidine | 0.99738 | 7.90 | 0.5–250 | 0.99368 | 14.20 | 0.5–250 | 1.00 | 217 |
| 131 | N-Nitrosopyrrolidine | 0.99830 | 14.70 | 1–250 | 0.99359 | 8.00 | 0.5–250 | 1.00 | 289 |
| 132 | Nonachlor-cis | 0.99591 | 7.20 | 0.2–250 | 0.97664 | 11.80 | 0.1–250 | 0.20 | 90 |
| 133 | Nonachlor-trans | 0.99924 | 4.70 | 0.1–250 | 0.99968 | 6.70 | 0.1–250 | 0.20 | 28 |
| 134 | Norflurazon | 0.99886 | 7.60 | 0.2–250 | 0.99937 | 5.10 | 0.05–250 | 0.05 | 23 |
| 135 | Ortho-phenylphenol | 0.99979 | 20.20 | 0.5–250 | 0.99957 | 14.70 | 0.5–250 | 1.00 | 102 |
| 136 | Oxadiazon | 0.99964 | 5.80 | 0.025–250 | 0.99970 | 14.10 | 0.1–250 | 0.20 | 22 |
| 137 | Oxyfluorfen | 0.99610 | 9.30 | 0.5–250 | 0.99445 | 18.80 | 0.5–250 | 1.00 | 73 |
| 138 | Paclobutrazol | 0.99977 | 5.40 | 0.05–250 | 0.99991 | 6.70 | 0.05–250 | 0.05 | 49 |
| 139 | Parathion (ethyl) | 0.99534 | 11.00 | 0.5–250 | 0.99395 | 19.40 | 0.5–250 | 1.00 | 158 |
| 140 | Parathion-methyl | 0.99478 | 12.00 | 0.05–250 | 0.99736 | 8.20 | 0.2–250 | 0.20 | 111 |
| 141 | Pebulate | 0.99885 | 14.30 | 0.5–250 | 0.99691 | 13.30 | 0.5–250 | 1.00 | 171 |
| 142 | Penconazole | 0.99985 | 6.30 | 0.05–250 | 0.99992 | 7.90 | 0.05–250 | 0.05 | 37 |
| 143 | Pentachloroaniline | 0.99973 | 6.10 | 0.1–250 | 0.99961 | 4.60 | 0.05–250 | 0.05 | 52 |
| 144 | Pentachloroanisole | 0.99939 | 5.30 | 0.05–250 | 0.99956 | 11.30 | 0.025–250 | 0.05 | 18 |
| 145 | Pentachlorobenzene | 0.99665 | 8.50 | 0.025–250 | 0.99765 | 11.40 | 0.025–250 | 0.05 | 12 |
| 146 | Pentachlorobenzonitrile | 0.99984 | 5.10 | 0.05–250 | 0.99973 | 8.40 | 0.025–250 | 0.05 | 23 |
| 147 | Pentachlorothioanisole | 0.99951 | 8.40 | 0.025–250 | 0.99973 | 6.00 | 0.05–250 | 0.05 | 22 |
| 148 | Perthane (Ethylan) | 0.99994 | 8.00 | 0.05–250 | 0.99982 | 4.50 | 0.1–250 | 0.20 | 30 |
| 149 | Permethrin peak 1 | 0.99971 | 9.90 | 1–250 | 0.99979 | 12.90 | 0.2–250 | 0.20 | 219 |
| 150 | Permethrin peak 2 | 0.99970 | 6.10 | 0.5–250 | 0.99979 | 5.80 | 0.5–250 | 1.00 | 48 |
| 151 | Phenothrin | 0.99950 | 17.90 | 1–250 | 0.99972 | 7.40 | 2–250 | 5.00 | 413 |
| 152 | Phorate | 0.99964 | 3.10 | 0.5–250 | 0.99910 | 13.50 | 0.025–250 | 0.05 | 18 |
| 153 | Phosalone | 0.99862 | 8.80 | 0.05–250 | 0.99982 | 8.90 | 0.05–250 | 0.05 | 18 |
| 154 | Phosmet | 0.99738 | 7.00 | 0.5–250 | 0.99916 | 24.50 | 0.2–250 | 0.20 | 54 |
| 155 | Piperonyl butoxide | 0.99977 | 6.50 | 0.1–250 | 0.99990 | 4.90 | 0.1–250 | 0.20 | 51 |
| 156 | Pirimiphos-ethyl | 0.99964 | 3.00 | 0.05–250 | 0.99967 | 5.30 | 0.025–250 | 0.05 | 21 |
| 157 | Pirimiphos-methyl | 0.99949 | 5.50 | 0.05–250 | 0.99949 | 4.80 | 0.025–250 | 0.05 | 18 |
| 158 | Pretilachlor | 0.99984 | 3.60 | 0.2–250 | 0.99989 | 2.40 | 0.2–250 | 0.20 | 44 |
| 159 | Prochloraz (parent) | 0.99749 | 14.30 | 1–250 | 0.99920 | 7.80 | 0.5–250 | 1.00 | 320 |
| 160 | Procymidone | 0.99991 | 4.00 | 0.1–250 | 0.99969 | 7.50 | 0.05–250 | 0.05 | 26 |

Appendix A (Part 5). Linearity data sets.

| Chart Number | Compound Name | Apple/Pear/Banana Linearity | | | Carrot/Potato Linearity | | | IDL ₉₉ | |
|--------------|------------------------------|-----------------------------|-----------|-------------|-------------------------|-----------|-------------|-------------------|----------|
| | | R ² | RF RSD(%) | Range (ppb) | R ² | RF RSD(%) | Range (ppb) | pg on Column | IDL (fg) |
| 161 | Profenofos | 0.99938 | 9.40 | 0.1–250 | 0.99654 | 10.40 | 0.1–250 | 0.20 | 53 |
| 162 | Propachlor | 0.99977 | 1.60 | 1–250 | 0.99926 | 15.30 | 0.025–250 | 0.05 | 13 |
| 163 | Propanil | 0.99925 | 12.30 | 0.025–250 | 0.99974 | 3.70 | 0.1–250 | 0.20 | 32 |
| 164 | Propargite | 0.99970 | 3.30 | 2–250 | 0.99881 | 19.40 | 2–250 | 5.00 | 1143 |
| 165 | Propisochlor | 0.99972 | 21.10 | 0.2–250 | 0.99953 | 5.80 | 0.5–250 | 1.00 | 284 |
| 166 | Propyzamide | 0.99966 | 4.50 | 0.1–250 | 0.99985 | 4.20 | 0.025–250 | 0.05 | 15 |
| 167 | Prothiofos | 0.99881 | 10.50 | 0.1–250 | 0.99842 | 5.90 | 0.1–250 | 0.20 | 28 |
| 168 | Pyraclufos | 0.99656 | 15.00 | 0.1–250 | 0.99920 | 20.60 | 0.05–250 | 0.05 | 28 |
| 169 | Pyrazophos | 0.99911 | 13.30 | 0.5–250 | 0.99979 | 16.60 | 0.2–250 | 0.20 | 163 |
| 170 | Pyridaben | 0.99987 | 3.10 | 0.2–250 | 0.99988 | 2.50 | 0.2–250 | 0.20 | 186 |
| 171 | Pyridaphenthion | 0.99941 | 3.80 | 0.2–250 | 0.99699 | 8.90 | 0.1–250 | 0.20 | 24 |
| 172 | Pyrimethanil | 0.99985 | 19.10 | 0.1–250 | 0.99971 | 8.40 | 0.05–250 | 0.05 | 23 |
| 173 | Pyriproxyfen | 0.99979 | 9.20 | 0.1–250 | 0.99990 | 2.30 | 0.2–250 | 0.20 | 24 |
| 174 | Quinalphos | 0.99926 | 10.80 | 0.5–250 | 0.99925 | 4.60 | 0.5–250 | 1.00 | 88 |
| 175 | Quintozene | 0.99912 | 10.20 | 0.2–250 | 0.99774 | 15.00 | 0.2–250 | 0.20 | 72 |
| 176 | Sulfotep | 0.99970 | 9.80 | 0.025–250 | 0.99962 | 9.40 | 0.025–250 | 0.05 | 26 |
| 177 | Sulprofos | 0.99986 | 3.30 | 0.05–250 | 0.99850 | 5.20 | 0.025–250 | 0.05 | 12 |
| 178 | Tebuconazole | 0.99983 | 9.10 | 0.5–250 | 0.99994 | 4.70 | 0.025–250 | 0.05 | 30 |
| 179 | Tebufenpyrad | 0.99980 | 4.20 | 0.05–250 | 0.99976 | 4.10 | 0.05–250 | 0.05 | 37 |
| 180 | Tecnazene | 0.99958 | 8.90 | 0.05–250 | 0.99815 | 13.90 | 0.025–250 | 0.05 | 15 |
| 181 | Tefluthrin | 0.99982 | 12.50 | 0.025–250 | 0.99944 | 6.60 | 0.025–250 | 0.05 | 34 |
| 182 | Terbacil | 0.99929 | 7.30 | 0.2–250 | 0.99974 | 5.40 | 0.1–250 | 0.20 | 95 |
| 183 | Terbufos | 0.99973 | 4.90 | 0.1–250 | 0.99978 | 5.00 | 0.05–250 | 0.05 | 13 |
| 184 | Terbuthylazine | 0.99967 | 8.60 | 0.2–250 | 0.99982 | 6.10 | 0.1–250 | 0.20 | 72 |
| 185 | Tetrachlorvinphos | 0.99941 | 8.20 | 0.05–250 | 0.99651 | 10.10 | 0.025–250 | 0.05 | 13 |
| 186 | Tetradifon | 0.99988 | 3.80 | 0.2–250 | 0.99990 | 17.50 | 0.025–250 | 0.05 | 17 |
| 187 | Tetrahydrophthalimide (THPI) | 0.99744 | 10.40 | 0.5–250 | 0.99985 | 6.00 | 0.5–250 | 1.00 | 67 |
| 188 | Tolclofos-methyl | 0.99985 | 4.70 | 0.05–250 | 0.99986 | 10.50 | 0.05–250 | 0.05 | 18 |
| 189 | Tolyfluanid | 0.99911 | 7.30 | 0.1–250 | 0.99952 | 7.90 | 0.1–250 | 0.20 | 27 |
| 190 | Triadimefon | 0.99965 | 8.80 | 0.05–250 | 0.99973 | 7.50 | 0.05–250 | 0.05 | 18 |
| 191 | Triadimenol | 0.99983 | 15.90 | 0.5–250 | 0.99973 | 11.20 | 0.5–250 | 1.00 | 116 |
| 192 | Triallate | 0.99983 | 2.20 | 0.1–250 | 0.99984 | 6.10 | 0.025–250 | 0.05 | 18 |
| 193 | Triazophos | 0.99937 | 6.30 | 0.1–250 | 0.99983 | 5.60 | 0.05–250 | 0.05 | 16 |
| 194 | Tricyclazole | 0.99883 | 9.50 | 2–250 | 0.99947 | 4.50 | 0.5–250 | 1.00 | 367 |
| 195 | Triflumizole | 0.99976 | 6.70 | 0.2–250 | 0.99978 | 6.00 | 0.5–250 | 1.00 | 147 |
| 196 | Vinclozolin | 0.99967 | 10.10 | 0.05–250 | 0.99969 | 5.50 | 0.05–250 | 0.05 | 22 |
| 197 | Tetramethrin peaks 1&2 | N/A | N/A | 0.5–250 | N/A | N/A | 0.5–250 | 5.00 | 929 |
| 198 | Resmethrin peaks 1&2 | N/A | N/A | 2–250 | N/A | N/A | 2–250 | 5.00 | 797 |
| 199 | Fluvalinate peaks 1&2 | N/A | N/A | 0.1–250 | N/A | N/A | 0.1–250 | 0.20 | 32 |
| 200 | Cypermethrin peaks I-IV | N/A | N/A | 1–250 | N/A | N/A | 1–250 | 1.00 | 214 |

Appendix A (Part 6). Linearity data sets.

| Chart Number | Compound Name | Apple/Pear/Banana Linearity | | | Carrot/Potato Linearity | | | IDL ₉₉ | |
|--------------|-----------------------|-----------------------------|-----------|-------------|-------------------------|-----------|-------------|-------------------|----------|
| | | R ² | RF RSD(%) | Range (ppb) | R ² | RF RSD(%) | Range (ppb) | pg on Column | IDL (fg) |
| 201 | Cyfluthrin peaks I-IV | N/A | N/A | 0.5–250 | N/A | N/A | 0.5–250 | 1.00 | 91 |
| N/A | Tetramethrin peak 1 | 0.99965 | 14.70 | N/A | 0.99968 | 16.50 | N/A | N/A | N/A |
| N/A | Tetramethrin peak 2 | 0.99974 | 18.80 | N/A | 0.99984 | 20.80 | N/A | N/A | N/A |
| N/A | Resmethrin peak 1 | 0.99950 | 13.60 | N/A | 0.99976 | 19.70 | N/A | N/A | N/A |
| N/A | Resmethrin peak 2 | 0.99983 | 6.60 | N/A | 0.99967 | 10.50 | N/A | N/A | N/A |
| N/A | Fluvalinate peak 1 | 0.99946 | 6.90 | N/A | 0.99936 | 5.50 | N/A | N/A | N/A |
| N/A | Fluvalinate peak 2 | 0.99933 | 10.20 | N/A | 0.99886 | 4.40 | N/A | N/A | N/A |
| N/A | Cypermethrin peak 1 | 0.99971 | 4.30 | N/A | 0.99989 | 13.40 | N/A | N/A | N/A |
| N/A | Cypermethrin peak 2 | 0.99987 | 2.40 | N/A | 0.99988 | 5.10 | N/A | N/A | N/A |
| N/A | Cypermethrin peak 3 | 0.99982 | 2.80 | N/A | 0.99975 | 3.10 | N/A | N/A | N/A |
| N/A | Cypermethrin peak 4 | 0.99981 | 2.70 | N/A | 0.99991 | 2.90 | N/A | N/A | N/A |
| N/A | Cyfluthrin peak 1 | 0.99967 | 2.50 | N/A | 0.99988 | 3.90 | N/A | N/A | N/A |
| N/A | Cyfluthrin peak 2 | 0.99905 | 4.40 | N/A | 0.99972 | 3.00 | N/A | N/A | N/A |
| N/A | Cyfluthrin peak 3 | 0.99837 | 10.70 | N/A | 0.99979 | 3.30 | N/A | N/A | N/A |
| N/A | Cyfluthrin peak 4 | 0.99723 | 8.10 | N/A | 0.99981 | 10.50 | N/A | N/A | N/A |

Appendix B – QuEChERS Recovery data

Appendix B (Part 1). QuEChERS Recovery data.

| Component Name | Carrot 1 µg/kg (n=6) | | Apple 1 µg/kg (n=6) | | Carrot 2.5 µg/kg (n=6) | | Apple 2.5 µg/kg (n=3) | | Carrot 10 µg/kg (n=6) | | Apple 10 µg/kg (n=6) | |
|------------------------------|----------------------|------------------|---------------------|------------------|------------------------|------------------|-----------------------|------------------|-----------------------|------------------|----------------------|------------------|
| | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) |
| 2,3,5,6-Tetrachloroaniline | 100.9% | 6.2% | 97.8% | 10.2% | 94.7% | 5.1% | 99.2% | 3.4% | 87.6% | 3.4% | 99.4% | 5.5% |
| 2,4'-Methoxychlor | 98.1% | 0.9% | 98.4% | 1.2% | 96.4% | 1.0% | 98.1% | 1.2% | 100.1% | 1.2% | 97.5% | 1.1% |
| 4,4'-Methoxychlor olefin | 99.9% | 2.8% | 106.7% | 3.5% | 98.8% | 2.6% | 99.4% | 1.6% | 101.0% | 1.1% | 100.2% | 1.6% |
| Acetochlor | 89.9% | 8.2% | 90.9% | 9.1% | 79.1% | 8.4% | 110.5% | 3.4% | 97.0% | 2.2% | 98.2% | 3.1% |
| Acrinathrin | 96.5% | 4.2% | 92.3% | 4.5% | 112.3% | 3.2% | 96.4% | 4.4% | 109.2% | 2.0% | 97.4% | 4.2% |
| Alachlor | 107.8% | 6.8% | 97.5% | 3.4% | 96.9% | 2.4% | 102.0% | 1.4% | 100.9% | 1.7% | 97.2% | 2.0% |
| Aldrin | 109.8% | 7.7% | 105.3% | 7.6% | 93.5% | 2.4% | 107.7% | 2.0% | 98.7% | 1.6% | 100.4% | 2.8% |
| Allethrin (Bioallethrin) | <LOQ | <LOQ | <LOQ | <LOQ | <LOQ | <LOQ | <LOQ | <LOQ | 105.7% | 1.2% | 102.1% | 2.2% |
| Allidochlor | 113.1% | 10.1% | 93.4% | 12.6% | 104.7% | 12.9% | 99.6% | 1.9% | 88.9% | 5.3% | 100.4% | 4.8% |
| Antraquinone | 36.3% | 27.8% | 102.4% | 14.5% | 39.6% | 6.5% | 94.6% | 3.5% | 34.1% | 47.7% | 95.8% | 2.7% |
| Atrazine | 111.1% | 4.1% | 94.4% | 6.0% | 90.9% | 4.0% | 98.7% | 5.0% | 98.7% | 0.9% | 98.4% | 2.4% |
| Azinphos-ethyl | 95.2% | 6.4% | 86.1% | 2.8% | 96.3% | 3.3% | 92.4% | 2.0% | 100.9% | 3.3% | 93.3% | 1.9% |
| Azinphos-methyl | 84.5% | 4.3% | 87.3% | 11.4% | 87.1% | 7.1% | 100.2% | 0.5% | 94.3% | 6.3% | 84.1% | 6.9% |
| BHC, Alpha | 102.0% | 3.6% | 95.0% | 5.6% | 94.3% | 3.6% | 104.2% | 3.2% | 94.1% | 1.9% | 99.8% | 4.2% |
| BHC, Beta | 98.3% | 2.3% | 96.3% | 7.7% | 94.8% | 0.9% | 96.7% | 6.3% | 97.7% | 1.0% | 98.2% | 1.9% |
| BHC, delta | 98.5% | 1.8% | 93.2% | 3.2% | 95.1% | 2.8% | 101.1% | 0.8% | 96.5% | 1.5% | 97.2% | 1.8% |
| BHC, gamma | 99.5% | 4.2% | 96.3% | 5.2% | 90.7% | 2.9% | 109.3% | 3.1% | 97.4% | 2.0% | 96.4% | 3.2% |
| Bifenthrin | 106.1% | 2.2% | 106.5% | 2.5% | 99.9% | 1.5% | 101.8% | 0.9% | 99.5% | 0.3% | 102.6% | 1.3% |
| Biphenyl | <LOQ | <LOQ | <LOQ | <LOQ | 86.4% | 10.5% | 118.1% | 3.0% | 79.5% | 7.7% | 111.6% | 7.1% |
| Bromfenvinphos | 111.2% | 3.2% | 97.0% | 2.7% | 91.8% | 1.9% | 103.6% | 2.8% | 97.3% | 1.7% | 94.6% | 2.0% |
| Bromfenvinphos-methyl | 110.5% | 3.4% | 87.8% | 5.1% | 92.3% | 2.5% | 106.4% | 2.3% | 98.4% | 0.9% | 95.2% | 3.6% |
| Bromophos-ethyl | 113.2% | 2.4% | 87.2% | 6.8% | 93.7% | 7.1% | 113.1% | 3.6% | 97.2% | 3.0% | 94.8% | 3.4% |
| Bromophos-methyl (Bromophos) | 104.0% | 4.3% | 90.0% | 7.6% | 93.2% | 2.9% | 114.8% | 2.5% | 95.9% | 2.6% | 97.3% | 3.2% |
| Bromopropylate | 101.2% | 14.2% | 97.2% | 5.3% | 96.1% | 2.0% | 111.1% | 2.6% | 97.1% | 2.3% | 97.4% | 2.3% |
| Bupirimate | 111.7% | 11.6% | 98.6% | 5.9% | 99.7% | 2.4% | 101.4% | 0.7% | 100.9% | 2.0% | 97.1% | 2.1% |
| Cadusafos | 101.9% | 3.4% | 101.4% | 3.8% | 96.4% | 3.8% | 101.9% | 1.8% | 97.6% | 1.7% | 101.4% | 3.1% |
| Captan | 85.0% | 16.5% | 80.4% | 9.1% | 96.7% | 11.9% | 78.8% | 2.4% | 97.6% | 5.6% | 87.5% | 4.2% |
| Carbophenothion | 103.2% | 4.3% | 99.3% | 3.6% | 95.6% | 1.5% | 96.6% | 1.9% | 99.2% | 1.0% | 97.1% | 1.5% |
| Carfentrazon-ethyl | 103.9% | 11.0% | 104.7% | 7.6% | 94.2% | 4.3% | 109.0% | 3.8% | 99.3% | 2.7% | 94.8% | 3.3% |
| Chlorbenside | 88.8% | 4.8% | 91.1% | 4.0% | 88.1% | 0.8% | 94.0% | 2.5% | 88.0% | 6.6% | 96.3% | 1.8% |
| Chlordane alpha-cis | 110.5% | 4.8% | 96.1% | 2.9% | 95.4% | 3.0% | 100.1% | 4.1% | 99.3% | 3.1% | 98.5% | 1.7% |
| Chlordane gamma-trans | 112.4% | 4.5% | 109.7% | 5.3% | 95.9% | 1.9% | 116.2% | 2.9% | 97.3% | 1.9% | 105.9% | 1.8% |
| Chlorfenapyr | 96.9% | 6.2% | 96.5% | 6.0% | 97.2% | 3.2% | 95.4% | 4.0% | 99.7% | 2.3% | 99.9% | 2.4% |
| Chlorfenoson | 97.9% | 2.5% | 95.3% | 3.5% | 92.9% | 1.8% | 94.8% | 0.4% | 93.1% | 2.2% | 94.4% | 3.7% |
| Chlorfenvinphos | 105.9% | 4.3% | 92.1% | 3.7% | 99.5% | 3.3% | 104.5% | 1.6% | 100.6% | 0.7% | 95.7% | 2.0% |
| Chlorobenzilate | 102.5% | 2.3% | 95.7% | 4.2% | 95.5% | 0.9% | 95.0% | 2.3% | 97.4% | 0.8% | 98.1% | 1.5% |
| Chloroneb | 102.6% | 4.3% | 97.6% | 5.5% | 99.9% | 6.3% | 105.6% | 4.8% | 89.0% | 4.6% | 102.0% | 6.6% |
| Chlorothalonil | 34.1% | 8.8% | 30.7% | 16.9% | 35.8% | 29.2% | 36.4% | 19.1% | 38.7% | 8.0% | 52.5% | 9.1% |
| Chlorpropham | 92.2% | 4.9% | 99.7% | 4.3% | 92.3% | 3.7% | 92.0% | 3.6% | 93.1% | 2.5% | 94.3% | 2.7% |
| Chlorpyrifos-ethyl | 100.1% | 5.6% | 99.5% | 6.2% | 94.1% | 2.1% | 101.7% | 1.2% | 96.7% | 1.2% | 97.1% | 1.2% |
| Chlorpyrifos-methyl | 113.9% | 2.8% | 88.1% | 5.5% | 93.3% | 5.8% | 104.9% | 2.5% | 95.3% | 1.9% | 98.0% | 2.7% |
| Chlorthal-dimethyl (Dacthal) | 117.1% | 4.7% | 96.9% | 4.0% | 93.1% | 1.5% | 109.9% | 5.4% | 98.0% | 1.4% | 94.8% | 3.8% |
| Chlorthiophos | 110.4% | 3.5% | 102.1% | 7.0% | 96.4% | 3.3% | 103.8% | 0.9% | 100.4% | 2.0% | 96.7% | 1.5% |
| Chlozolinate | 117.5% | 8.4% | 99.7% | 8.6% | 94.6% | 4.1% | 98.4% | 1.9% | 100.6% | 1.8% | 97.3% | 3.2% |
| Clomazone | 99.1% | 3.6% | 98.3% | 4.0% | 95.0% | 2.8% | 96.8% | 0.8% | 97.8% | 0.9% | 99.0% | 2.9% |
| Coumaphos | 98.6% | 6.7% | 93.0% | 3.5% | 88.5% | 5.6% | 91.4% | 2.3% | 94.9% | 3.7% | 92.5% | 4.2% |
| Cycloate | <LOQ | <LOQ | <LOQ | <LOQ | <LOQ | <LOQ | <LOQ | <LOQ | 114.7% | 6.3% | 115.2% | 9.0% |

Appendix B (Part 2). QuEChERS Recovery data.

| Component Name | Carrot 1 µg/kg (n=6) | | Apple 1 µg/kg (n=6) | | Carrot 2.5 µg/kg (n=6) | | Apple 2.5 µg/kg (n=3) | | Carrot 10 µg/kg (n=6) | | Apple 10 µg/kg (n=6) | |
|----------------------------|----------------------|------------------|---------------------|------------------|------------------------|------------------|-----------------------|------------------|-----------------------|------------------|----------------------|------------------|
| | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) |
| Cyfluthrin peak 1 | 109.7% | 3.2% | 103.2% | 7.3% | 106.0% | 3.1% | 96.1% | 16.7% | 106.3% | 2.0% | 103.6% | 2.6% |
| Cyfluthrin peak 2 | 95.5% | 3.5% | 98.5% | 2.5% | 104.1% | 1.7% | 103.0% | 1.4% | 105.7% | 2.6% | 101.2% | 2.3% |
| Cyfluthrin peak 3 | 107.8% | 4.8% | 105.7% | 7.2% | 103.3% | 3.5% | 97.5% | 2.5% | 104.7% | 2.2% | 101.1% | 2.1% |
| Cyfluthrin peak 4 | 110.3% | 4.9% | 98.3% | 6.4% | 100.5% | 2.4% | 99.9% | 1.8% | 105.3% | 2.6% | 103.3% | 2.7% |
| Cyhalothrin I (lambda) | 106.2% | 4.5% | 100.3% | 3.6% | 102.8% | 3.4% | 99.5% | 2.4% | 104.3% | 1.8% | 104.5% | 1.8% |
| Cypermethrin peak 1 | 80.2% | 3.6% | 94.8% | 2.4% | 94.8% | 4.0% | 101.6% | 1.2% | 101.1% | 2.8% | 101.2% | 2.2% |
| Cypermethrin peak 2 | 87.0% | 10.2% | 80.2% | 6.7% | 104.7% | 2.8% | 98.0% | 0.8% | 106.7% | 4.1% | 105.1% | 1.9% |
| Cypermethrin peak 3 | 102.0% | 5.6% | 99.0% | 4.9% | 109.1% | 1.3% | 101.9% | 2.7% | 105.6% | 3.6% | 107.0% | 2.4% |
| Cypermethrin peak 4 | 102.1% | 2.9% | 102.8% | 5.9% | 102.2% | 4.0% | 101.8% | 0.5% | 107.5% | 3.2% | 107.1% | 3.1% |
| Cyprodinil | 103.5% | 3.2% | 93.5% | 2.1% | 96.3% | 2.8% | 99.2% | 2.6% | 92.8% | 4.8% | 98.7% | 2.7% |
| DDD p,p | 101.1% | 1.9% | 102.0% | 2.4% | 96.6% | 1.2% | 99.2% | 0.8% | 99.1% | 1.3% | 97.3% | 1.5% |
| DDD, o, p | 104.1% | 2.2% | 104.0% | 4.1% | 95.5% | 1.8% | 101.4% | 1.3% | 98.5% | 1.3% | 96.3% | 1.6% |
| DDE o,p | 109.1% | 1.7% | 97.2% | 2.5% | 96.9% | 0.9% | 103.9% | 2.9% | 99.7% | 1.1% | 99.0% | 1.7% |
| DDE p, p | 106.9% | 1.7% | 99.3% | 4.0% | 96.2% | 1.5% | 97.8% | 2.4% | 98.8% | 1.1% | 97.5% | 1.5% |
| DDT o,p | 101.4% | 2.8% | 100.2% | 4.5% | 95.1% | 3.3% | 96.4% | 2.9% | 100.7% | 0.8% | 97.4% | 1.8% |
| DDT p,p | 97.9% | 2.6% | 97.0% | 0.8% | 95.4% | 2.0% | 97.3% | 2.1% | 100.7% | 0.8% | 96.6% | 1.5% |
| Deltamethrin | 77.2% | 5.2% | 95.2% | 7.0% | 114.9% | 6.0% | 100.0% | 8.5% | 110.6% | 2.3% | 101.2% | 7.4% |
| Diazinon | 102.0% | 3.5% | 103.9% | 4.4% | 102.3% | 2.9% | 97.4% | 4.0% | 99.0% | 2.1% | 100.9% | 3.1% |
| Dichlobenil | 97.2% | 6.4% | 94.6% | 10.5% | 98.9% | 7.3% | 103.4% | 5.6% | 86.1% | 5.6% | 102.5% | 6.5% |
| Dichlofuanid | 58.8% | 4.1% | 56.8% | 4.9% | 50.7% | 13.9% | 62.5% | 7.6% | 57.5% | 3.0% | 74.9% | 5.3% |
| Dichlorobenzophenone, 4, 4 | 115.3% | 2.7% | 99.1% | 4.1% | 97.3% | 1.8% | 95.2% | 1.4% | 98.0% | 2.7% | 94.2% | 1.7% |
| Dicloran (Bortran) | 86.6% | 6.0% | 84.9% | 6.8% | 86.6% | 3.8% | 91.7% | 4.1% | 86.4% | 3.7% | 88.8% | 2.4% |
| Dicofol | 34.9% | 15.0% | 69.0% | 12.1% | 113.2% | 5.2% | 92.3% | 9.6% | 105.9% | 15.6% | 101.3% | 11.5% |
| Dieldrin | 104.9% | 17.3% | 98.9% | 13.4% | 103.0% | 6.4% | 105.5% | 3.2% | 96.6% | 3.2% | 95.4% | 3.0% |
| Dimethachlor | 102.2% | 3.2% | 94.3% | 1.8% | 94.3% | 1.7% | 98.5% | 3.2% | 98.6% | 2.4% | 96.9% | 1.2% |
| Dimethoate | 97.7% | 10.4% | 71.9% | 8.5% | 87.0% | 10.1% | 77.3% | 1.3% | 94.9% | 1.8% | 85.3% | 5.8% |
| Diphenamid | 102.6% | 6.1% | 82.9% | 4.4% | 100.8% | 1.9% | 104.7% | 2.3% | 99.8% | 1.8% | 98.8% | 1.2% |
| Diphenylamine | 98.1% | 2.1% | 113.2% | 7.1% | 97.4% | 4.2% | 107.7% | 4.5% | 94.3% | 2.0% | 102.3% | 4.3% |
| Disulfoton | 96.5% | 4.0% | 94.3% | 8.8% | 95.8% | 1.5% | 100.1% | 5.1% | 98.3% | 1.6% | 97.4% | 2.5% |
| Edifenphos | 96.1% | 3.3% | 93.5% | 2.6% | 96.6% | 1.5% | 92.5% | 1.8% | 100.3% | 2.2% | 96.9% | 2.0% |
| Endosulfan ether | 105.3% | 5.6% | 98.8% | 3.0% | 95.7% | 2.6% | 101.5% | 3.0% | 94.5% | 1.5% | 97.8% | 3.0% |
| Endosulfan peak 1 | 101.7% | 6.6% | 107.1% | 6.2% | 94.8% | 2.8% | 116.7% | 4.6% | 99.0% | 2.1% | 104.7% | 2.8% |
| Endosulfan peak 2 | 97.9% | 6.1% | 95.6% | 5.4% | 100.7% | 5.3% | 98.2% | 2.9% | 98.0% | 1.8% | 101.6% | 2.1% |
| Endosulfan sulfate | 101.2% | 4.9% | 106.0% | 8.3% | 100.0% | 2.3% | 108.2% | 0.7% | 100.5% | 1.6% | 100.8% | 1.5% |
| Endrin | 105.6% | 3.5% | 93.4% | 10.7% | 99.3% | 2.1% | 99.7% | 1.2% | 98.0% | 1.5% | 96.8% | 2.8% |
| Endrin Aldehyde | 55.6% | 14.3% | 27.6% | 26.3% | 40.0% | 18.9% | 27.2% | 28.2% | 37.6% | 13.2% | 29.0% | 16.7% |
| Endrin-Ketone | 91.5% | 17.1% | 105.2% | 10.9% | 84.3% | 16.2% | 116.0% | 4.9% | 96.0% | 2.3% | 97.3% | 3.0% |
| EPN | 116.0% | 9.0% | 96.9% | 7.3% | 92.6% | 2.2% | 87.9% | 6.2% | 95.5% | 2.7% | 93.4% | 2.8% |
| Ethion | 107.0% | 3.7% | 98.3% | 2.3% | 96.8% | 1.4% | 97.8% | 1.4% | 100.0% | 1.2% | 99.3% | 1.5% |
| Ethoprop (Ethoprophos) | 97.2% | 4.3% | 91.4% | 4.9% | 96.5% | 2.6% | 96.0% | 3.8% | 95.0% | 3.0% | 97.0% | 3.8% |
| Etofenprox | 107.0% | 4.6% | 109.1% | 4.3% | 106.7% | 2.2% | 96.3% | 5.0% | 104.9% | 1.6% | 110.6% | 2.3% |
| Etridiazole (Terrazole) | 100.0% | 7.6% | 91.9% | 11.5% | 100.5% | 7.2% | 102.8% | 6.3% | 87.1% | 6.6% | 100.3% | 7.4% |
| Fenamiphos | 101.5% | 8.3% | 87.9% | 10.1% | 92.1% | 4.7% | 71.8% | 8.3% | 97.9% | 1.4% | 84.7% | 3.7% |
| Fenarimol | 107.9% | 3.7% | 101.3% | 1.6% | 104.5% | 1.4% | 102.5% | 1.7% | 100.3% | 0.8% | 99.8% | 1.1% |
| Fenchlorfos | 112.6% | 3.0% | 97.7% | 3.8% | 95.3% | 2.6% | 103.0% | 3.2% | 98.1% | 2.1% | 101.1% | 2.3% |
| Fenitrothion | 101.7% | 9.5% | 79.5% | 9.1% | 87.0% | 6.1% | 97.9% | 2.8% | 91.9% | 1.0% | 85.0% | 2.8% |
| Fenpropathrin | 97.2% | 6.0% | 80.4% | 5.8% | 102.5% | 6.0% | 97.5% | 2.7% | 98.1% | 1.3% | 103.4% | 1.6% |

Appendix B (Part 3). QuEChERS Recovery data.

| Component Name | Carrot 1 µg/kg (n=6) | | Apple 1 µg/kg (n=6) | | Carrot 2.5 µg/kg (n=6) | | Apple 2.5 µg/kg (n=3) | | Carrot 10 µg/kg (n=6) | | Apple 10 µg/kg (n=6) | |
|---------------------------------|----------------------|------------------|---------------------|------------------|------------------------|------------------|-----------------------|------------------|-----------------------|------------------|----------------------|------------------|
| | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) |
| Fenson | 98.9% | 2.5% | 95.5% | 3.8% | 96.5% | 2.2% | 96.0% | 2.0% | 98.8% | 1.1% | 97.0% | 2.0% |
| Fenthion | 109.0% | 4.5% | 92.3% | 2.3% | 100.4% | 2.7% | 108.7% | 1.9% | 99.6% | 2.2% | 95.1% | 2.2% |
| Fenvalerate | 101.3% | 2.9% | 102.6% | 4.2% | 110.1% | 2.6% | 104.6% | 2.6% | 110.3% | 1.8% | 104.2% | 4.0% |
| Fipronil | 103.5% | 7.8% | 83.8% | 8.6% | 94.9% | 3.7% | 110.6% | 4.4% | 100.5% | 2.8% | 95.9% | 3.3% |
| Fipronil desulfinyl | 101.5% | 8.4% | 92.2% | 8.3% | 95.2% | 5.4% | 108.4% | 2.1% | 99.5% | 1.8% | 98.5% | 3.6% |
| Fluazifop-P-butyl | 111.0% | 3.9% | 96.2% | 6.6% | 95.8% | 2.2% | 100.8% | 2.4% | 99.7% | 2.3% | 97.3% | 1.4% |
| Fludioxonil | 105.1% | 3.2% | 93.7% | 3.7% | 89.3% | 4.5% | 93.9% | 5.0% | 92.3% | 3.3% | 89.6% | 1.6% |
| Fluquinconazole | 103.1% | 16.6% | 102.9% | 7.0% | 94.3% | 3.0% | 91.1% | 3.9% | 98.2% | 2.3% | 99.4% | 3.0% |
| Fluridone | 100.7% | 17.2% | 85.6% | 11.3% | 92.1% | 4.2% | 107.2% | 0.4% | 100.9% | 2.3% | 88.7% | 1.5% |
| Flusilazole | 95.2% | 5.4% | 98.4% | 9.7% | 96.9% | 1.9% | 92.0% | 2.1% | 95.6% | 1.7% | 98.8% | 2.1% |
| Flutolanil | 99.0% | 2.7% | 95.1% | 1.9% | 95.0% | 0.9% | 96.7% | 0.6% | 97.0% | 0.8% | 99.1% | 1.6% |
| Flutriafol | 96.9% | 2.5% | 99.4% | 5.4% | 93.3% | 2.2% | 99.8% | 2.4% | 95.6% | 1.4% | 97.8% | 1.0% |
| Fluvalinate peak 1 | 100.3% | 5.9% | 115.1% | 18.5% | 118.9% | 3.6% | 98.5% | 3.5% | 120.0% | 2.4% | 100.6% | 3.2% |
| Fluvalinate peak 2 | 93.7% | 17.3% | 108.4% | 18.8% | 117.6% | 5.8% | 104.1% | 8.2% | 113.6% | 3.9% | 101.9% | 5.2% |
| Folpet | 90.2% | 11.4% | 72.2% | 4.1% | 86.6% | 10.8% | 70.7% | 7.1% | 97.0% | 4.2% | 77.3% | 4.4% |
| Fonofos | 105.9% | 3.9% | 99.9% | 4.6% | 99.5% | 2.1% | 105.4% | 4.8% | 97.1% | 2.7% | 101.3% | 2.5% |
| Heptachlor | 108.7% | 5.4% | 97.1% | 5.1% | 95.9% | 3.4% | 105.2% | 3.6% | 94.8% | 0.9% | 98.9% | 3.5% |
| Hexachlorobenzene | 95.1% | 7.1% | 94.6% | 8.3% | 86.7% | 2.9% | 109.1% | 4.6% | 75.4% | 9.7% | 99.3% | 5.4% |
| Hexazinone | 89.1% | 1.9% | 87.7% | 2.1% | 93.5% | 1.0% | 91.6% | 0.9% | 95.0% | 1.7% | 93.9% | 2.0% |
| Iodofenfos | 97.6% | 4.6% | 96.9% | 3.2% | 83.3% | 6.7% | 96.7% | 0.8% | 93.5% | 4.2% | 94.6% | 4.3% |
| Iprodione | 95.2% | 16.3% | 99.6% | 7.6% | 97.3% | 5.5% | 108.6% | 1.2% | 107.3% | 3.8% | 90.5% | 2.9% |
| Isazophos | 111.1% | 9.4% | 72.2% | 5.0% | 105.2% | 4.0% | 101.5% | 2.1% | 98.7% | 1.5% | 99.1% | 2.5% |
| Isodrin | 103.3% | 4.9% | 98.0% | 4.6% | 96.0% | 1.9% | 100.4% | 4.8% | 97.2% | 0.8% | 98.7% | 2.0% |
| Lenacil | 90.0% | 10.8% | 92.5% | 8.2% | 93.8% | 6.0% | 88.3% | 4.6% | 94.0% | 5.3% | 92.8% | 2.0% |
| Leptophos | 98.9% | 6.8% | 101.1% | 1.8% | 97.9% | 1.3% | 95.0% | 1.1% | 97.8% | 4.0% | 102.4% | 1.1% |
| Linuron | 104.4% | 9.9% | 82.4% | 3.2% | 92.9% | 4.7% | 97.2% | 7.8% | 90.9% | 3.6% | 99.7% | 4.5% |
| Malathion | 103.0% | 2.1% | 82.6% | 3.7% | 91.5% | 2.1% | 91.4% | 5.1% | 94.3% | 1.4% | 97.6% | 1.3% |
| Metalaxyl | 106.0% | 6.0% | 99.8% | 12.5% | 96.2% | 4.8% | 111.5% | 1.8% | 98.5% | 1.9% | 96.0% | 2.3% |
| Metazachlor | 98.5% | 5.0% | 99.8% | 4.9% | 97.4% | 1.3% | 100.4% | 0.2% | 98.7% | 0.9% | 97.2% | 1.6% |
| Methacrifos | 95.6% | 4.0% | 95.7% | 5.5% | 101.9% | 6.6% | 103.1% | 4.0% | 92.6% | 3.8% | 102.5% | 6.6% |
| Methoxychlor | 95.5% | 2.3% | 97.0% | 5.3% | 98.2% | 2.0% | 102.6% | 1.7% | 102.0% | 0.9% | 98.9% | 1.2% |
| Metolachlor | 98.9% | 2.9% | 93.3% | 2.4% | 94.2% | 1.1% | 96.7% | 3.1% | 97.8% | 1.5% | 97.3% | 1.2% |
| Mevinphos | 94.2% | 4.6% | 89.4% | 5.4% | 99.3% | 7.7% | 94.4% | 4.3% | 104.1% | 3.6% | 95.9% | 5.3% |
| MGK-264 A | 109.5% | 16.6% | 96.3% | 10.5% | 102.3% | 6.8% | 94.6% | 0.6% | 100.0% | 2.4% | 97.3% | 1.9% |
| MGK-264 B | 105.9% | 4.6% | 100.4% | 6.1% | 96.7% | 2.9% | 105.4% | 3.4% | 97.1% | 1.5% | 100.8% | 2.2% |
| Mirex | 97.6% | 4.0% | 103.6% | 2.2% | 93.9% | 1.8% | 102.0% | 1.0% | 102.5% | 1.7% | 99.3% | 1.4% |
| Myclobutanil | 100.6% | 4.5% | 97.6% | 5.4% | 96.7% | 2.7% | 101.5% | 2.8% | 98.1% | 2.0% | 96.5% | 2.1% |
| N-(2,4-Dimethylphenyl)formamide | 86.6% | 5.6% | 85.0% | 8.4% | 90.0% | 7.6% | 77.6% | 3.0% | 89.8% | 5.0% | 86.8% | 3.0% |
| NDBA | 110.7% | 7.6% | 110.9% | 10.8% | 105.2% | 9.0% | 116.5% | 7.8% | 90.0% | 7.3% | 104.8% | 6.4% |
| NDEA | 94.9% | 7.2% | 91.1% | 11.1% | 97.3% | 10.2% | 105.3% | 5.2% | 97.0% | 12.2% | 101.6% | 13.7% |
| NDPA | 97.2% | 8.1% | 74.8% | 13.5% | 97.3% | 7.7% | 90.5% | 11.7% | 85.4% | 9.4% | 101.3% | 8.4% |
| NEMA | 87.5% | 4.1% | 85.2% | 19.4% | 91.7% | 15.3% | 104.7% | 15.3% | 83.1% | 12.4% | 97.2% | 8.2% |
| Nitrofen | 109.8% | 9.3% | 97.4% | 2.2% | 90.9% | 3.3% | 93.7% | 1.7% | 92.8% | 1.5% | 87.9% | 3.2% |
| N-Nitrosodiphenylamine | 98.1% | 2.1% | 113.2% | 7.1% | 97.4% | 4.2% | 107.7% | 4.5% | 94.3% | 2.0% | 102.3% | 4.3% |
| N-Nitrosomorpholine | 84.9% | 8.2% | 80.7% | 10.3% | 85.1% | 8.4% | 80.5% | 3.5% | 81.8% | 6.9% | 91.8% | 6.8% |
| N-Nitrosopiperidine | 95.5% | 7.2% | 81.0% | 19.8% | 105.0% | 11.2% | 98.9% | 10.5% | 86.5% | 8.2% | 102.3% | 6.6% |
| N-Nitrosopyrrolidine | 95.4% | 11.0% | 114.0% | 13.8% | 101.6% | 6.4% | 91.4% | 14.2% | 82.5% | 6.1% | 99.4% | 9.7% |

Appendix B (Part 4). QuEChERS Recovery data.

| Component Name | Carrot 1 µg/kg (n=6) | | Apple 1 µg/kg (n=6) | | Carrot 2.5 µg/kg (n=6) | | Apple 2.5 µg/kg (n=3) | | Carrot 10 µg/kg (n=6) | | Apple 10 µg/kg (n=6) | |
|-------------------------|----------------------|------------------|---------------------|------------------|------------------------|------------------|-----------------------|------------------|-----------------------|------------------|----------------------|------------------|
| | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) | Mean Recovery | Precision RSD(%) |
| N-Nitrosopyrrolidine | 95.4% | 11.0% | 114.0% | 13.8% | 101.6% | 6.4% | 91.4% | 14.2% | 82.5% | 6.1% | 99.4% | 9.7% |
| Nonachlor-cis | 109.4% | 15.3% | 114.4% | 8.8% | 94.6% | 10.4% | 113.1% | 7.2% | 100.5% | 2.5% | 98.9% | 4.1% |
| Nonachlor-trans | 107.8% | 3.7% | 103.7% | 8.7% | 96.1% | 5.1% | 107.3% | 2.2% | 99.4% | 2.1% | 101.0% | 2.6% |
| Norflurazon | 100.4% | 7.8% | 83.9% | 17.3% | 93.7% | 3.5% | 96.7% | 1.8% | 95.5% | 1.8% | 91.8% | 2.7% |
| Ortho-phenylphenol | 90.6% | 5.1% | 115.5% | 3.5% | 82.3% | 6.2% | 102.6% | 2.1% | 90.9% | 3.2% | 99.7% | 5.4% |
| Oxadiazon | 100.3% | 4.3% | 97.7% | 4.6% | 95.2% | 1.9% | 92.9% | 1.1% | 96.9% | 2.1% | 95.5% | 3.6% |
| Oxyfluorfen | 102.4% | 15.7% | 113.2% | 8.9% | 97.7% | 7.1% | 101.6% | 4.5% | 95.6% | 2.0% | 97.0% | 3.2% |
| Paclobutrazol | 111.2% | 2.9% | 93.8% | 4.4% | 98.3% | 1.9% | 97.4% | 2.0% | 95.1% | 3.9% | 94.6% | 1.9% |
| Parathion (ethyl) | 118.4% | 7.4% | 111.1% | 5.7% | 96.0% | 4.0% | 111.5% | 4.5% | 92.0% | 3.4% | 90.4% | 1.8% |
| Parathion-methyl | 119.6% | 6.7% | 83.7% | 4.1% | 94.0% | 2.5% | 95.8% | 1.1% | 91.0% | 3.3% | 86.9% | 2.7% |
| Pebulate | 112.9% | 4.6% | 118.8% | 5.6% | 113.9% | 7.6% | 119.3% | 0.3% | 93.4% | 5.4% | 109.2% | 6.9% |
| Penconazole | 100.3% | 4.7% | 96.3% | 3.5% | 94.4% | 1.7% | 100.1% | 5.3% | 98.1% | 0.9% | 94.4% | 2.7% |
| Pentachloroaniline | 94.9% | 2.8% | 91.8% | 4.3% | 85.1% | 4.0% | 104.1% | 0.3% | 84.6% | 8.2% | 96.1% | 2.7% |
| Pentachloroanisole | 104.3% | 4.1% | 92.8% | 4.9% | 94.8% | 4.3% | 104.6% | 3.7% | 90.6% | 2.4% | 102.5% | 5.6% |
| Pentachlorobenzene | 100.1% | 4.0% | 94.5% | 6.3% | 95.8% | 5.6% | 117.9% | 6.9% | 83.2% | 4.4% | 100.9% | 8.9% |
| Pentachlorobenzonitrile | 102.5% | 5.7% | 91.4% | 3.3% | 95.2% | 6.4% | 103.3% | 1.9% | 87.7% | 2.7% | 96.6% | 3.5% |
| Pentachlorothioanisole | 106.7% | 6.3% | 97.4% | 3.2% | 86.8% | 3.4% | 106.5% | 0.2% | 82.8% | 9.7% | 98.0% | 2.3% |
| Permethrin peak 1 | 91.6% | 13.6% | 81.8% | 6.2% | 82.8% | 8.1% | 98.9% | 2.0% | 89.4% | 2.3% | 97.4% | 2.1% |
| Permethrin peak 2 | 97.6% | 7.0% | 98.7% | 3.2% | 96.2% | 2.6% | 98.9% | 0.3% | 100.4% | 2.5% | 104.8% | 2.0% |
| Perthane (Ethylan) | 109.8% | 3.1% | 101.1% | 2.4% | 96.6% | 0.7% | 103.1% | 0.8% | 101.0% | 0.8% | 99.4% | 1.4% |
| Phenothrin | <LOQ | <LOQ | <LOQ | <LOQ | 89.7% | 14.9% | 81.7% | 13.3% | 108.4% | 11.3% | 92.2% | 4.2% |
| Phorate | 116.6% | 15.9% | 103.1% | 7.0% | 102.5% | 4.8% | 103.3% | 3.0% | 98.9% | 3.0% | 101.7% | 4.8% |
| Phosalone | 100.2% | 3.7% | 97.2% | 4.6% | 98.4% | 1.8% | 70.4% | 13.3% | 100.4% | 2.1% | 99.1% | 1.2% |
| Phosmet | 85.2% | 3.8% | 89.6% | 6.9% | 89.3% | 3.6% | 89.7% | 2.2% | 95.3% | 3.3% | 88.7% | 6.4% |
| Piperonyl butoxide | 105.0% | 6.8% | 99.7% | 2.4% | 103.1% | 1.3% | 99.9% | 0.5% | 102.2% | 1.2% | 103.8% | 1.8% |
| Pirimiphos-ethyl | 104.9% | 4.3% | 107.5% | 4.0% | 100.2% | 2.0% | 102.9% | 4.2% | 100.0% | 2.5% | 98.1% | 2.8% |
| Pirimiphos-methyl | 117.2% | 5.1% | 97.0% | 4.8% | 93.6% | 1.9% | 110.7% | 4.2% | 99.1% | 1.6% | 94.0% | 4.0% |
| Pretilachlor | 103.3% | 4.1% | 95.3% | 5.0% | 94.6% | 2.0% | 98.4% | 5.2% | 99.4% | 1.0% | 98.1% | 1.2% |
| Prochloraz | 104.5% | 6.2% | 110.6% | 12.3% | 115.0% | 7.9% | 98.9% | 2.3% | 99.3% | 4.1% | 92.0% | 3.9% |
| Procymidone | 115.0% | 2.8% | 100.8% | 5.4% | 96.4% | 3.0% | 103.3% | 1.5% | 98.7% | 1.5% | 99.0% | 3.2% |
| Profenofos | 114.4% | 7.1% | 99.0% | 6.8% | 86.5% | 4.3% | 99.9% | 4.2% | 95.7% | 3.5% | 91.6% | 3.0% |
| Propachlor | 94.4% | 8.4% | 104.0% | 2.1% | 99.3% | 7.2% | 99.9% | 2.8% | 92.7% | 1.3% | 97.7% | 4.6% |
| Propanil | 89.6% | 7.5% | 78.8% | 10.2% | 91.3% | 5.6% | 91.6% | 2.3% | 95.1% | 4.2% | 89.5% | 2.1% |
| Propargite | <LOQ | <LOQ | <LOQ | <LOQ | 89.1% | 7.3% | 101.7% | 4.9% | 103.2% | 11.1% | 102.5% | 5.2% |
| Propisochlor | 97.5% | 7.0% | 99.2% | 3.6% | 101.7% | 3.2% | 99.2% | 1.5% | 101.3% | 3.0% | 99.1% | 1.6% |
| Propyzamide | 99.8% | 6.1% | 102.1% | 8.9% | 100.7% | 1.5% | 99.5% | 2.7% | 101.6% | 1.3% | 99.6% | 2.4% |
| Prothiofos | 115.8% | 4.8% | 92.3% | 2.1% | 92.1% | 3.5% | 105.4% | 4.7% | 96.0% | 1.2% | 98.2% | 4.0% |
| Pyraclufos | 96.9% | 8.3% | 89.6% | 7.2% | 92.3% | 3.6% | 87.6% | 1.8% | 94.1% | 6.6% | 94.1% | 4.1% |
| Pyrazophos | 110.6% | 2.6% | 96.0% | 3.0% | 93.7% | 3.8% | 97.5% | 1.9% | 101.9% | 3.1% | 101.1% | 1.4% |
| Pyridaben | 106.8% | 2.7% | 100.0% | 2.7% | 98.9% | 1.7% | 119.5% | 1.0% | 100.1% | 2.1% | 102.8% | 1.8% |
| Pyridaphenthion | 94.2% | 18.0% | 96.3% | 8.6% | 96.3% | 3.3% | 113.7% | 2.6% | 100.4% | 1.9% | 95.9% | 2.2% |
| Pyrimethanil | 118.4% | 7.6% | 80.4% | 13.3% | 117.0% | 9.1% | 104.7% | 9.1% | 101.3% | 2.7% | 95.8% | 2.8% |
| Pyriproxyfen | 103.9% | 4.5% | 105.7% | 1.4% | 99.5% | 2.9% | 80.0% | 7.3% | 101.2% | 1.3% | 102.9% | 1.6% |
| Quinalphos | 95.2% | 4.3% | 86.3% | 11.5% | 78.9% | 3.7% | 85.1% | 4.9% | 100.0% | 2.3% | 95.5% | 2.1% |
| Quintozene | 107.3% | 8.7% | 101.9% | 6.1% | 97.6% | 5.6% | 94.6% | 2.2% | 88.9% | 1.6% | 94.3% | 4.0% |
| Resmethrin peak 1 | <LOQ | <LOQ | <LOQ | <LOQ | 94.7% | 2.9% | 104.0% | 7.1% | 96.3% | 3.7% | 87.5% | 2.0% |
| Resmethrin peak 2 | <LOQ | <LOQ | <LOQ | <LOQ | 92.9% | 7.7% | 95.4% | 4.3% | 94.4% | 2.6% | 89.0% | 6.0% |

