

Determination of Pesticides and Persistent Organic Pollutants in Honey by Accelerated Solvent Extraction and GC-MS/MS

Aaron Kettle and Fabrizio Galbiati, Thermo Fisher Scientific, 1214 Oakmead Parkway, Sunnyvale, CA, USA, 94085

INTRODUCTION

Honey is a natural product that is widely used for both nutritional and medicinal purposes. It is generally considered a natural and healthy product of animal origin, free of impurities. However, honeybees are subject to a number of viral, bacterial, fungal, and parasitic diseases and infestations. Insecticides, fungicides, and acaricides are used to protect colonies against infestations from hive beetle and parasites. Many pollutants in the environment can also contaminate the bees themselves in addition to their pollen, honey, and other bee products. Pollutants such as organochlorine pesticides (OCs), polychlorobiphenyls (PCBs), organophosphates (OPs), and polybromodiphenylethers (PBDEs) are a particular threat due to their environmental persistence and ability to bioaccumulate in the food chain. Due to the potential toxicity, a comprehensive workflow method for the extraction and analysis of these environmental pollutants is of growing importance to ensure the health and safety of bees and their honey.

Among the available extraction techniques, accelerated solvent extraction (ASE) offers shorter extraction times and reduced solvent consumption. ASE uses high temperatures combined with high pressure. A high temperature allows a higher rate of extraction due to a reduction in viscosity and surface tension, and increases the solubility and diffusion rate into the sample. The method reported here is applicable for the extraction and analysis of four different classes of compounds (6 PCBs, 7 PBDEs, 16 OCs, and 19 OPs) in honey using ASE and GC-MS/MS.

MATERIALS AND METHODS

Sample Collection and Preparation

Beekeepers from three different Italian regions: Calabria, South Italy (14 samples); Trentino Alto Adige, North Italy (18 samples) and Lombardia, North Italy (27 samples) provided 59 organic honey samples, as summarized in Table 1. All samples were stored at -20 °C prior to analysis to prevent matrix decomposition.

Table 1. Areas of Origin for Organic Honey Samples.

| Number of Samples | Area of Origin | Botanical Source | Potential Environmental Contamination Sources |
|-------------------|----------------------------|---------------------|---|
| 27 | Lombardia (North of Italy) | Multifloral | Industrialized Area (OCs, PCBs, PBDEs) |
| 14 | Trentino (North of Italy) | Multifloral | Intensive Apple Orchard (Pesticides) |
| 18 | Calabria (South of Italy) | Citrus (Monofloral) | Intensive Citrus Orchard (Pesticides) |

Working solutions were prepared by diluting the stock solution in hexane for pesticides and then stored at -40 °C. Mixed compound calibration solution, in hexane, was prepared daily from the stock solutions (10 µg/mL) and the appropriate volume was used as a spiking solution.

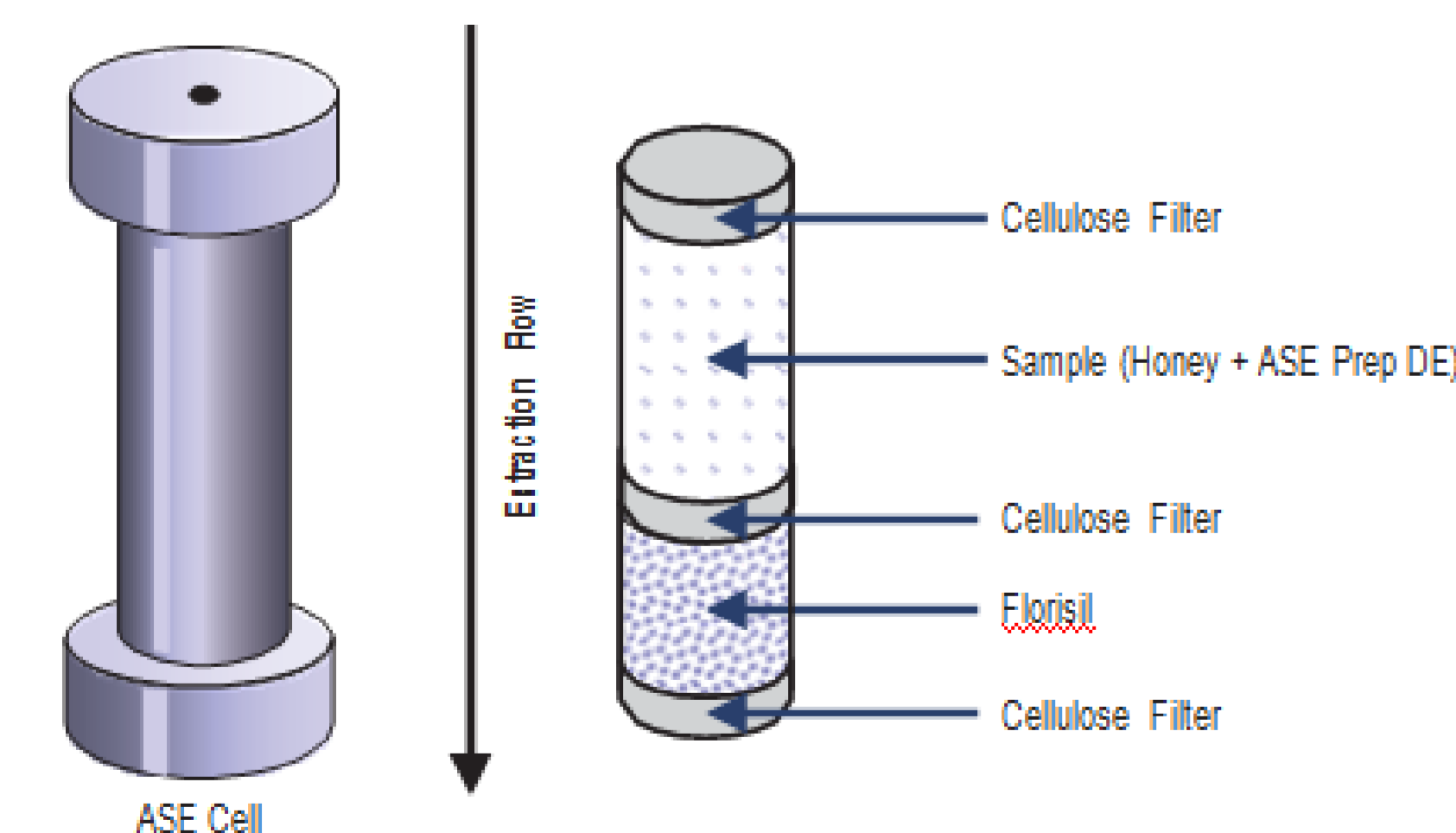
The extractions were carried out using a Thermo Scientific™ Dionex™ ASE™ 350 Accelerated Solvent Extractor shown in Figure 1, equipped with 34 mL stainless steel extraction cells. The extracts were collected in 60 mL vials, treated with sodium sulfate and directly concentrated in a 2 mL autosampler glass vial using a Thermo Scientific™ Rocket™ Evaporator system (Figure 1).

Figure 1. Dionex ASE 350 Accelerated Solvent Extractor and Rocket Evaporator.



A cellulose filter was placed in the bottom of a 34 mL extraction cell (Figure 2), followed by 5 g of activated Florisil and another cellulose filter. A 2 g sample of honey was homogenized with an equal weight of Thermo Scientific™ Dionex™ ASE™ Prep DE dispersant, sodium sulfate and transferred into the cell. One mL of isooctane solution containing the three internal standards was added. The remaining empty volume was filled with Dionex ASE Prep DE dispersant. The extractor was programmed according to the conditions reported in Table 2. The extracts were collected in 60 mL vials and treated with sodium sulfate to remove any possible water. After filtration, the organic phase was concentrated to dryness using the Rocket Evaporator system, dissolved in 200 µL of isooctane, and submitted to analysis by GC-MS/MS.

Figure 2. Extraction Cell Schematic.



| Parameter | Setting |
|--|-----------------------------------|
| Table 2. Dionex ASE 350 Accelerated Solvent Extractor Extraction Method. | |
| Solvent | n-Hexane/Ethyl Acetate (4:1, v/v) |
| Temperature | 80 °C |
| Pressure | 1500 psi |
| Static Cycles | 3 |
| Static Cycle Time | 3 min |
| Rinse Volume | 90% |
| Purge Time | 90 s |
| Extraction Time per Sample | ~15 min |
| Solvent Used per Sample | ~50 mL |

Analytical Methods

The samples were analyzed using a Thermo Scientific™ TRACE™ 1310 gas chromatograph equipped split/splitless injector, a fused-silica capillary column (Rt-5MS Crossbond-5% diphenyl 95% dimethylpolysiloxane, 35 m × 0.25 mm × 0.25 µm) and a Thermo Scientific™ TSQ™ 8000 Triple Quadrupole GC-MS/MS (Figure 3). The method conditions for the gas chromatograph and mass spectrometer are listed in Tables 3 and 4.

Figure 3. TSQ 8000 Triple Quadrupole GC-MS/MS.



Table 3. GC and Injector Conditions.

| Parameter | Setting |
|----------------------|---|
| Injector Type | PTV, Splitless |
| Injector Temperature | 250 °C |
| Liner | 2 x 2.75 x 120 mm |
| Injected Volume | 1 µL |
| Splitless Time | 0.5 min |
| Splitflow | 20 mL/min |
| GC Column | Rt-5MS (35m x 0.25 mm x 0.25 µm) |
| Carrier Gas | Helium, 99.999% purity |
| Flow Rate | 1.0 mL/min, constant |
| Initial Temperature | 80 °C (3 min) 10 °C/min to 170 °C 3 °C/min to 190 °C 2 °C/min to 240 °C 3 °C/min to 280 °C 10 °C/min to 310 °C |
| Final Temperature | 310 °C (5 min) |

Table 4. Mass Spectrometer Parameters.

| Parameter | Setting |
|--------------------|-------------|
| Source Temperature | 270 °C |
| Ionization | EI |
| Electron Energy | 70 eV |
| Emission Current | 50 µA |
| Q2 Gas Pressure | 1.5 mTorr |
| Collision Energy | 10 to 30 eV |
| Q1 Peak Width FWHM | 0.7 Da |
| Q3 Peak Width FWHM | 0.7 Da |

RESULTS

A multiresidue method for the analysis of organic contaminants and pesticides was developed. The ASE extraction with inline cleanup was necessary for the removal of interfering substances (e.g., carbohydrates) from honey samples. Florisil proved to be very efficient for the cleanup of different foods as well as honey samples. The proposed method was optimized for the multiresidual analysis of 59 pesticides and persistent organic pollutants (POPs). Total ion current chromatograms (GC-MS/MS) of blank honey samples spiked with investigated compounds and a naturally contaminated sample are shown in Figures 4 and 5.

Figure 4. Total Ion Current (GC-MS/MS) Chromatogram of a Spiked Honey Sample (10 ng/g).

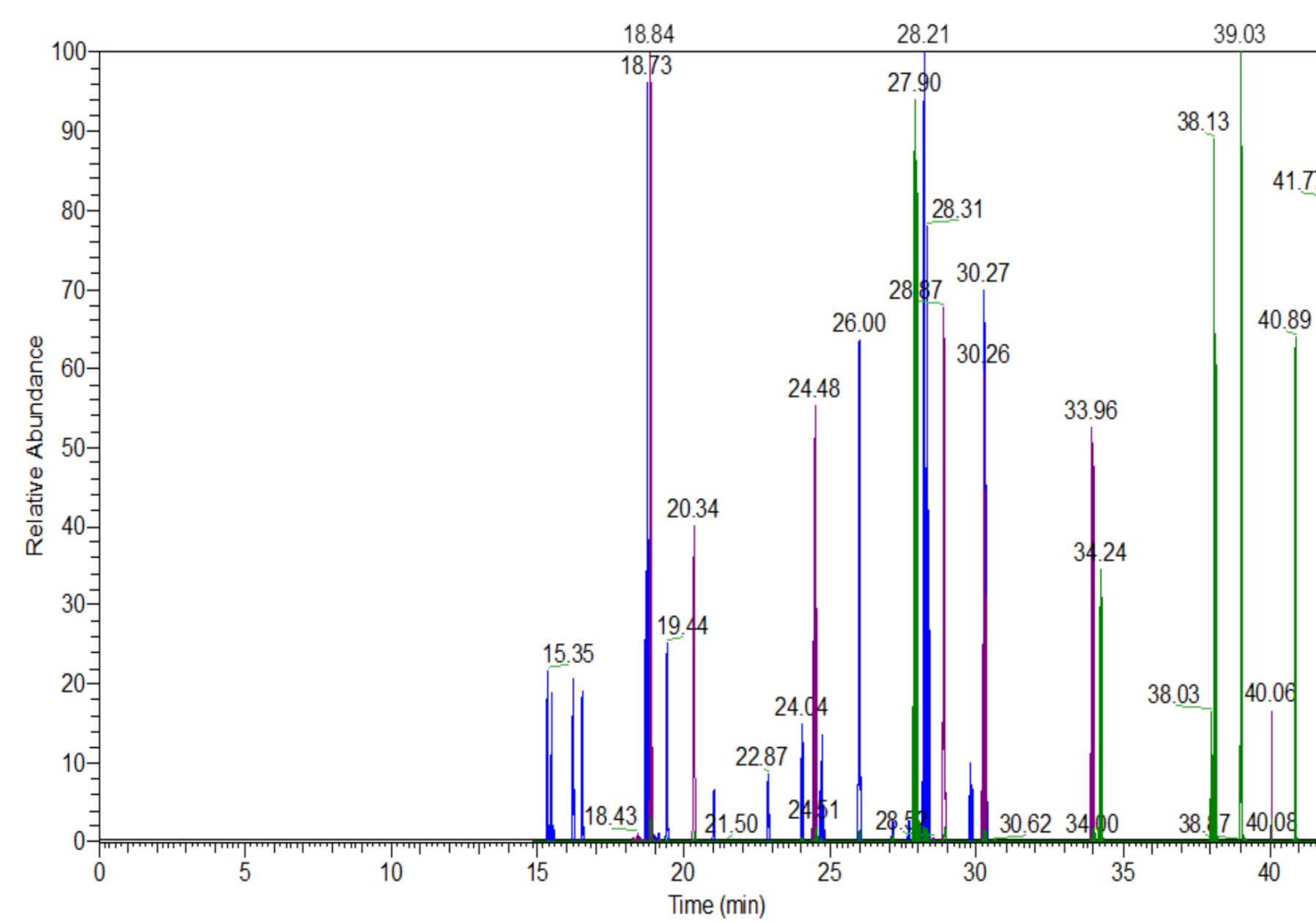
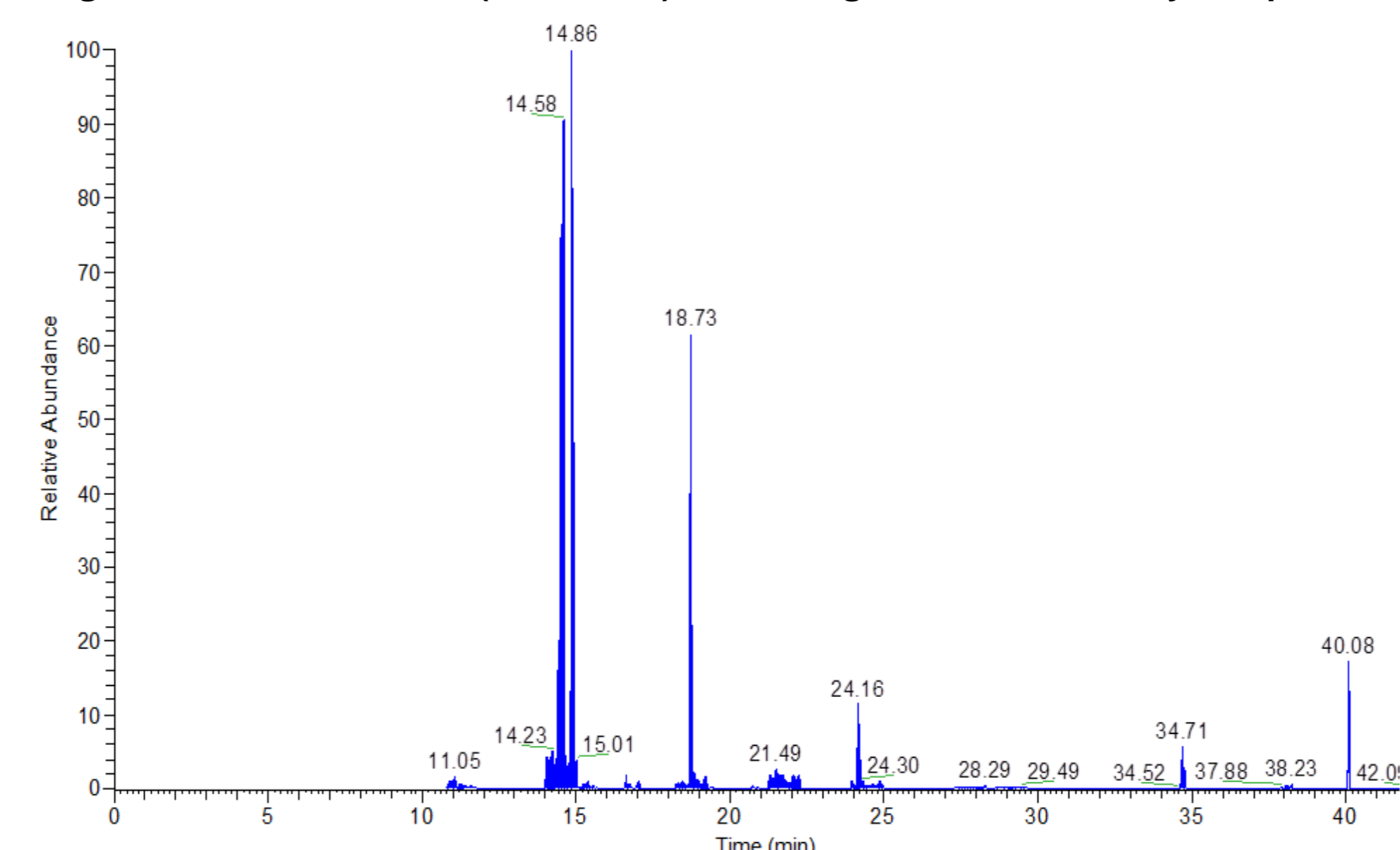


Figure 5. Total Ion Current (GC-MS/MS) Chromatogram of a Raw Honey Sample.



The method showed good linearity with determination coefficients equal to or higher than 0.99 for all of the compounds investigated. The method also showed good repeatability, demonstrating that it is effective for monitoring compounds belonging to different chemical classes (Table 5). The recoveries ranged from 97 to 102% for PCBs and PBDEs, from 75 to 95% for OCs, from 75 to 97% for OPs and from 75% to 102% for the additional agrochemicals. The CVs ranged from 4 to 14%. The one-step ASE method, using Florisil as an interference retainer, is rapid, cost-effective, and minimizes waste generation compared to the classic methods. Combining the extraction and the two clean-up steps (i.e., GPC and SPE) in a single ASE step reduced laboratory time by half. At present, this is the first ASE application using an inline clean-up step to screen for the presence of different pesticides and organic contaminants in honey.

| Contaminants | LOD (ng/g) | LOQ (ng/g) | Recovery % (RSD) | Coefficient of Determination (r ²) |
|---|------------|------------|------------------|--|
| Polychlorobiphenyls (PCBs) | | | | |
| PCB 28 | 0.08 | 0.24 | 102 (7) | 0.9994 |
| PCB 52 | 0.07 | 0.21 | 103 (7) | 0.9999 |
| PCB 101 | 0.04 | 0.12 | 97 (4) | 0.9999 |
| PCB 138 | 0.05 | 0.15 | 105 (4) | 0.9999 |
| PCB 153 | 0.02 | 0.06 | 102 (4) | 0.9999 |
| PCB 180 | 0.06 | 0.18 | 98 (9) | 0.9999 |
| Polybrominated Diphenyl Ethers (PBDEs) | | | | |
| PBDE 28 | 0.01 | 0.03 | 100 (9) | 0.9991 |
| PBDE 33 | 0.02 | 0.06 | 98 (9) | 0.9999 |
| PBDE 47 | 0.02 | 0.06 | 97 (8) | 0.9996 |
| PBDE 99 | 0.03 | 0.09 | 102 (7) | 0.9998 |
| PBDE 100 | 0.01 | 0.03 | 103 (7) | 0.9998 |
| PBDE 153 | 0.03 | 0.09 | 97 (10) | 0.9992 |
| PBDE 154 | 0.02 | 0.06 | 100 (12) | 0.9999 |
| Organochlorine Pesticides (OCs) | | | | |
| α-HCH | 0.99 | 2.97 | 78 (10) | 0.9959 |
| Hexachlorobenzene | 1.26 | 3.78 | 80 (12) | 0.9945 |
| β-HCH | 1.17 | 3.51 | 85 (12) | 0.9995 |
| Lindane (γ-HCH) | 0.79 | 2.39 | 96 (10) | 0.9985 |
| Heptachlor | 0.95 | 2.84 | 93 (12) | 0.9996 |
| Aldrin | 0.85 | 2.55 | 75 (14) | 0.9991 |
| Heptachlor Epoxide | 0.91 | 2.73 | 77 (14) | 0.9994 |
| trans-Chlordane | 1.48 | 4.44 | 92 (10) | 0.9993 |
| Endosulfan I | 1.13 | 3.38 | 80 (13) | 0.9992 |
| pp'-DDE | 0.85 | 2.55 | 97 (12) | 0.9994 |
| Endrin | 0.99 | 2.98 | 88 (11) | 0.9998 |
| Endosulfan II | 1.14 | 3.42 | 90 (10) | 0.9993 |
| pp'-DDD | 0.91 | 2.74 | 87 (14) | 0.9986 |
| op-DDT | 0.94 | 2.83 | 82 (14) | 0.9963 |
| Endosulfan Sulfate | 1.07 | 3.22 | 85 (12) | 0.9921 |
| pp'-DDT | 0.91 | 2.74 | 95 (12) | 0.9992 |
| Organophosphorus (OPs) | | | | |
| Mevinphos | 0.75 | 2.25 | 75 (12) | 0.9996 |
| Ethoprophos | 0.44 | 1.32 | 86 (10) | 0.9991 |
| Dichlorvos | 0.33 | 0.99 | 93 (10) | 0.9997 |
| Phorate | 0.52 | 1.56 | 75 (13) | 0.9993 |
| Demephron (-O and -S) | 1.12 | 3.36 | 77 (14) | 0.9992 |
| Diazinon | 1.10 | 3.30 | 90 (10) | 0.9994 |
| Disulfoton | 0.90 | 2.70 | 80 (14) | 0.9998 |
| Parathion-methyl | 0.83 | 2.49 | 97 (8) | 0.9993 |
| Fenchlorphos | 1.12 | 3.36 | 88 (11) | 0.9986 |
| Chlorpyrifos | 0.95 | 2.85 | 90 (9) | 0.9963 |
| Fenthion | 0.78 | 2.34 | 87 (12) | 0.9996 |
| Trichlorfon | 0.98 | 2.94 | 85 (14) | 0.9998 |
| Tetrachlorvinphos | 1.12 | 3.36 | 85 (12) | 0.9998 |
| Prothiofos | 0.75 | 2.25 | 92 (12) | 0.9992 |
| Terbuphos | 0.68 | 2.04 | 90 (12) | 0.9963 |
| Fensulfothion | 1.09 | 3.27 | 88 (10) | 0.9921 |
| Sulprofos | 0.98 | 2.94 | 87 (10) | 0.9992 |
| Azinphos-methyl | 1.07 | 3.21 | 87 (12) | 0.9978 |
| Coumaphos | 0.78 | 2.34 | 84 (14) | 0.9962 |

CONCLUSIONS

An analytical method was developed and successfully applied to evaluate pesticides and POP residues in organic honey samples produced in three different Italian regions that are characterized by different contamination sources. The method proved to be simple and rapid, requiring small sample sizes, and minimizing solvent consumption, due to the ASE with an inline cleanup step. MS/MS detection provided both quantitative information and the confirmation of POP residues in honey, confirming the one-step ASE method as a valid alternative to classical extraction methods.

REFERENCES

- Thermo Fisher Scientific Customer Application Note (CAN) 125: Determination of Pesticides and Persistent Organic Pollutants in Honey by Accelerated Solvent Extraction and GC-MS/MS.
- Chiesa, L.M.; Labella, G.F.; Panseri, S.; Britti, D.; Galbiati, F.; Villa, R.; Arioli, F. Accelerated solvent extraction by using an 'in-line' clean-up approach for multiresidue analysis of pesticides in organic honey *Food Additives & Contaminants: Part A*. 2017, 34, 809-818.

ACKNOWLEDGEMENTS

We would like to thank Professor Luca Chiesa from the University of Milan for the fruitful collaboration and Dr. Sara Panseri and her team who performed the work on the honey samples.

TRADEMARKS/LICENSING

© 2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.