



Food safety

Reducing running costs for GC-MS/MS analysis of pesticide residues using hydrogen carrier gas

Authors

Łukasz Rajski¹, Daniel Kutscher¹,
Adam Ladak²

¹Thermo Fisher Scientific, Bremen, Germany

²Thermo Fisher Scientific,
Hemel Hempstead, United Kingdom

Keywords

Helium saving, hydrogen carrier gas,
HeSaver-H₂Safer, TSQ 9610 Triple
Quadrupole GC-MS/MS, pesticides,
SRMs, timed-SRM, QuEChERS

Goal

This application note demonstrates the performance of the Thermo Scientific™ TSQ™ 9610 Triple Quadrupole GC-MS/MS system for the analysis of more than 180 pesticide residues using hydrogen carrier gas.

Introduction

Helium is the standard carrier gas for gas chromatography thanks to its high chromatographic efficiency and inertness. This is particularly important when considering applications involving mass spectrometric detection (GC-MS), where these characteristics ensure the best performance in terms of sensitivity. This has resulted in well-established, validated, and regulated GC and GC-MS methods using helium as the carrier gas. However, helium is a limited and non-renewable natural resource,¹ and recent price increases and supply issues caused by shortages have led laboratories and instrument manufacturers to investigate possible mitigation options, such as switching to alternative carrier gases or reducing the consumption of helium.

A possible alternative carrier gas, long established in regulated methods involving capillary GC combined with other detection systems such as flame ionization detection (FID), is hydrogen. Hydrogen has a diffusivity comparable to helium and is even preferable because its viscosity is about half, allowing lower inlet pressure for a given gas velocity, and maintaining the optimum separation efficiency at higher flow rates,

thus offering shorter analysis time. Additionally, hydrogen generators can create the required amount exactly when needed, offering laboratories a safe, steady, and renewable supply of hydrogen without the need to store a flammable gas. Therefore, using hydrogen instead of helium is an attractive solution because of its analytical and practical advantages; however, there are still important considerations before it can be used. Due to its flammability, a GC system using hydrogen needs to be equipped with a hydrogen sensor in the GC oven, capable of turning off the heating and the gas supply in case of detected leaks. Many existing validated methods are based on helium and must be re-optimized and re-validated in the case of carrier gas conversion. This is particularly important for GC-MS methods as the use of hydrogen affects the vacuum conditions in the ion source and therefore the ionization process, often with consequent decrease of the response. Despite these shortcomings, many GC and GC-MS applications are successfully run by using hydrogen, with clear advantages in terms of speed and cost of analysis.

A common application of gas chromatography coupled to triple quadrupole mass spectrometry is the analysis of pesticide residues in food samples. This is essential to monitor if pesticides are being applied on crops in compliance with regulations (label instructions) to protect the health of consumers. This application note demonstrates the performance of the Thermo Scientific™ TSQ™ 9610 Triple Quadrupole GC-MS/MS system equipped with a Thermo Scientific™ Advanced Electron Ionization (AEI) source for the analysis of more than 180 pesticide residues, many only accessible for analysis using GC-MS/MS.

Instrumentation

The Thermo Scientific™ TRACE™ 1610 GC connected to the TSQ 9610 mass spectrometer was equipped with a Thermo Scientific™ iConnect™ HeSaver-H₂Safer™ SSL Injector module. The HeSaver-H₂Safer technology decouples the gas used for pressurizing the inlet and splitting the sample from the actual carrier gas supplied only to the column for the gas chromatographic separation of the analytes. It can be used in conjunction with helium or hydrogen carrier gas and decreases the carrier gas consumption considerably. The pressurizing gas can be an inexpensive alternative (e.g., nitrogen or argon), replacing the valuable carrier gas for the split and purge flows, and it is efficiently prevented from entering the column during separation. The pressurizing gas has also the benefit to eliminate

the contact between the analytes and reactive carrier gas (hydrogen) in the hot injector, and consequently decreases the possibility of unwanted reactions. When the HeSaver-H₂Safer mode is used, a significant cost savings and prolonged cylinder lifetime can be realized when using helium, whereas the main advantage of the limited gas consumption for hydrogen is related to operational safety. The maximum amount of hydrogen flowing in the system is limited so that there is no risk to reach hazardous concentrations even in case of column breakage, and therefore a hydrogen sensor in the GC oven is not necessary.

The TSQ 9610 mass spectrometer was operated in timed-SRM mode. All the transitions were carefully optimized with Thermo Scientific™ AutoSRM software to obtain the highest possible sensitivity. The GC-MS method details are listed in Table 1, whereas the transitions, collision energies, and retention times can be found in Appendix 1. For the data acquisition and treatment, Thermo Scientific™ Chromeleon™ Chromatography Data System (CDS) software was used.

The compound identification criteria were taken from the DG SANTE guidance document.² According to this document, two ion transitions are required to identify a given pesticide. The chromatographic peaks of those transitions must fully overlap, and signal-to-noise ratio must be equal or greater than 3. The ion ratio should be within $\pm 30\%$ (relative) of average of calibration standards. The retention time shift should not be greater than ± 0.1 min. The aim was to check the performance at the default Maximum Residues Limit (MRL) of 0.01 mg/kg, applicable for baby food, without considering the MRLs mentioned in the Commission Directive 2006/125/EC of 5 December 2006.

A full list of transitions used with hydrogen as the carrier gas is included in Appendix 1. In a previously published method using helium, in general three transitions were monitored per analyte. However, the method presented here applied, in the great majority of cases, only two transitions. Moreover, the retention time window with hydrogen was set narrower than the window applied for helium (0.3 min vs. 0.7 min). The differences in the total transitions number and in the retention time window width influenced the dwell time applied per analyte, which could be extended. In the method with helium carrier gas, the shortest dwell time was 3.28 ms, whereas the current method with hydrogen carrier uses 8.48 ms, thus helping to increase the sensitivity and compensate for losses typically occurring with the use of hydrogen as the carrier gas.

Table 1. GC-MS method parameters

TRACE 1610 GC	
Injector	
Injector type	iConnect Split/Splitless (SSL) injector upgraded with the HeSaver-H ₂ Safer (P/N MI-190000-0022)
Temperature [°C]	260
Pressuring gas	Nitrogen
Split mode	Splitless
Split flow [mL/min]	50 (Nitrogen)
Splitless time [min]	1
Septum purge flow [mL/min]	5 (Nitrogen)
Vacuum compensation	On
Liner	Thermo Scientific™ LinerGOLD™ Single Taper Liner with Wool (P/N 453A1925-U)
Injection volume [μL]	1
Oven	
Analytical column	Thermo Scientific™ TraceGOLD™ TG-5SilMS (20 m × 0.18 mm × 0.18 μm) (P/N 26096-5780)
Carrier gas	Hydrogen
Carrier gas flow [mL/min]	1.2
Oven temperature program	
Ready delay [min]	1.3
Temperature 1 [°C]	40
Hold [min]	1.5
Rate [°C/min]	25
Temperature 2 [°C]	90
Hold [min]	1.5
Rate [°C/min]	25
Temperature 3 [°C]	180
Hold [min]	0
Rate [°C/min]	5
Temperature 4 [°C]	280
Hold [min]	0
Rate [°C/min]	10
Temperature 5 [°C]	300
Hold [min]	5
TSQ 9610 GC-MS/MS	
Ion source	Advanced Electron Ionization (AEI) Source
Transfer line temperature [°C]	300
Ion source temperature [°C]	320
Emission current [μA]	10
Electron energy [eV]	50

Sample preparation

During this study, two matrices (baby food and honey) were tested, and the results evaluated for a total of 181 pesticides.

Samples were prepared using the QuEChERS method.³ In short, a 10 g portion of the sample was weighed in a 50 mL PTFE centrifuge tube, subsequently 10 mL of acetonitrile was added and the samples were shaken in an automatic axial extractor for 7 min. Afterwards, 4 g of magnesium sulphate, 1 g of sodium chloride, 1 g of trisodium citrate dihydrate, and 0.5 g of disodium hydrogen citrate sesquihydrate were added (Thermo Scientific™ QuEChERS EN 15662 Method Extraction Kit, P/N S1-10-EN-KIT), and the samples were again shaken in the automatic axial extractor for 7 min. The extract was then centrifuged (3,700 rpm) for 5 min. A 5 mL volume of the supernatant was transferred to a 15 mL PTFE centrifuge tube containing 750 mg of MgSO₄ and 125 mg of PSA. The extract was shaken in a vortex for 30 s and centrifuged again (3,700 rpm) for 5 min. The extract was transferred to an amber vial and acidified with 10 μL formic acid 5% per mL of extract.

Consequences of switching the carrier gas type from helium to hydrogen

Substitution of helium with hydrogen has some important consequences for the chromatographic separation as well as for the MS detection. Maintaining the same column, carrier gas flow rate, and oven program, hydrogen provides shorter retention times than helium. That should be considered as an advantage because it improves the sample throughput. However, the retention time shift is not the same for all the analytes (in this study retention time shifts between 0.5 and 2 minutes were observed depending on the overall elution time of a given compound), so the new retention times must be confirmed first. This can be accomplished with additional injections of standard solutions with wider retention time windows, or in full scan MS mode.

Hydrogen provides a better chromatographic efficiency (often expressed as a higher number of the theoretical plates), which results in slightly narrower chromatographic peaks (compared to helium), however this had a negligible impact on the MS/MS method, as the number of the data points per chromatographic peak remained unchanged.

For some of the analytes, the presence of hydrogen affected the ionization process, and the spectrum obtained with hydrogen was different than the spectrum acquired with helium. Figure 1 shows a full mass spectrum for fipronil using both helium and hydrogen. Not only is the base peak different, but the relative intensities of the remaining major ions have also changed. This difference should not be considered a critical issue, as targeted methods used for pesticide residues analysis rely on compound identification based on retention time and ion ratios between compound specific transitions, and not on the fidelity to spectral libraries. Nevertheless, in some cases, a re-optimization of the SRM transitions must be performed to search for more sensitive alternatives. The sensitivity of the final method was sufficient to analyze the pesticides at 0.005 mg/kg. Figure 2 shows examples of the chromatographic peak obtained at this concentration level.

Linearity

Linearity was checked in the concentration range 0.005–0.500 mg/kg. According to the DG SANTE guidance document, the linearity evaluation is based on the back-calculated concentrations. To include a point into the calibration curve, its back-calculated concentration should not deviate from the true concentration by more than 20%. More than 97% of the 181 evaluated compounds showed a linear response within the investigated concentration range in both the tested matrices (more detailed information can be found in Appendix 2). Figure 3 highlights the linearity observed for chlorfenson in baby food and for triadimenol in honey, as a representative example.

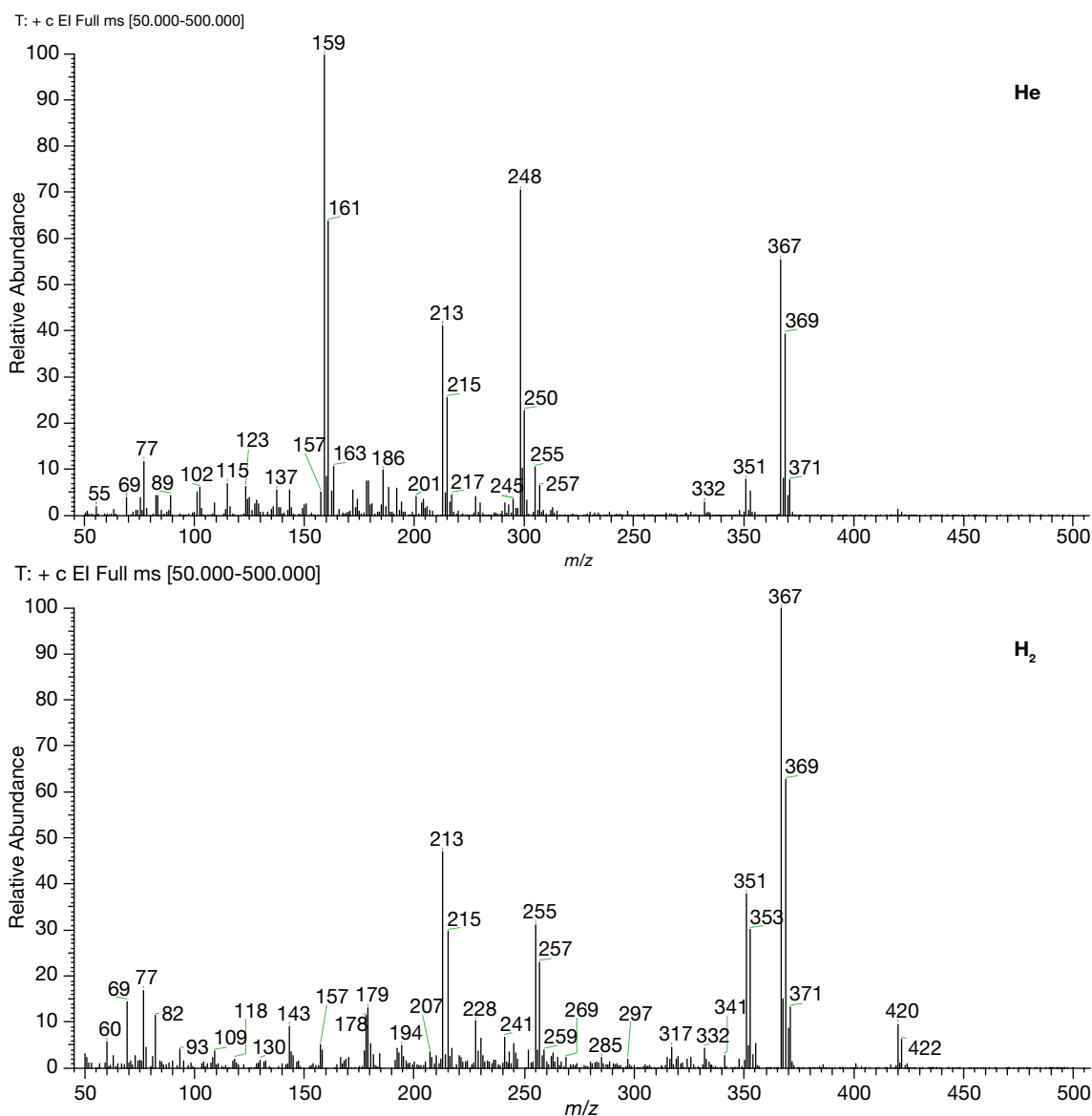


Figure 1. Fipronil, full scan spectra obtained with helium (top) and hydrogen (bottom) carrier gas

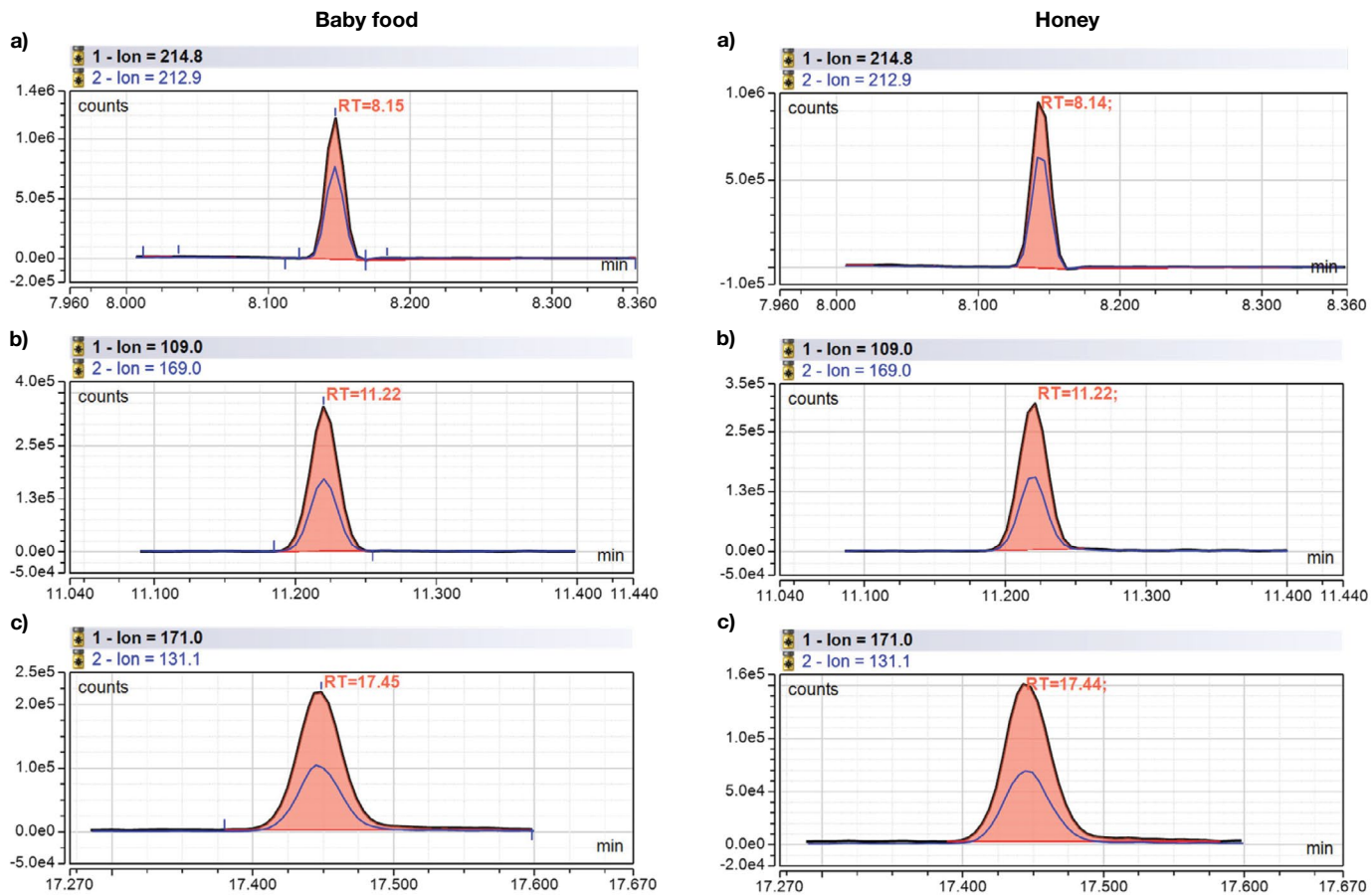


Figure 2. Examples of peaks obtained at 0.005 mg/kg in baby food and honey, using hydrogen as carrier gas: a) pentachlorobenzene, b) fenthion, c) tebufenpyrad

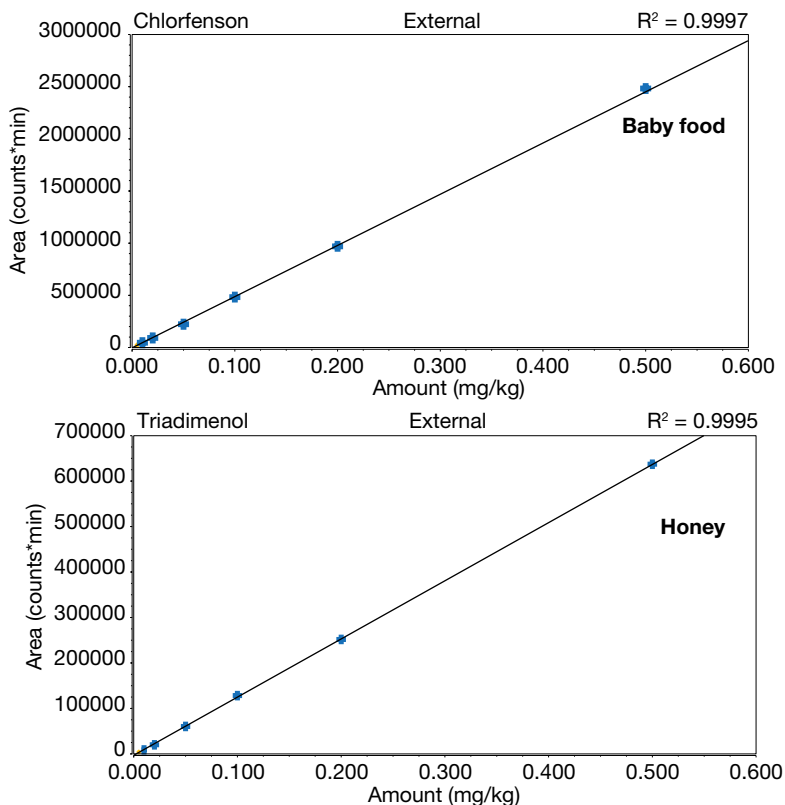


Figure 3. Calibration curve of chlorfensfen in baby food and triadimenol in honey, with hydrogen as carrier gas; concentration range 0.005–0.500 mg/kg

Repeatability

Repeatability was evaluated at two concentration levels: 0.01 mg/kg and 0.05 mg/kg. The samples were injected in quintuplicate and subsequently the relative standard deviation was calculated. In both matrices, 98% of the evaluated pesticides showed an RSD equal or lower than 10%. At 0.05 mg/kg the results were slightly better, with 99% of the analytes showing RSD \leq 10%. Although the SANTE guideline allows for RSD values of up to 20% for a given compound, it is important to consider that this number applies to the full workflow, including extraction and clean-up of the sample. Therefore, an analytical method must allow for some margin to always ensure compliance. Figure 4 shows the distribution of the RSD results (for more details see Appendix 2), including data obtained on baby food with helium carrier gas as reference (grey bar). As can be seen, the performance with hydrogen is highly comparable to the performance with helium.

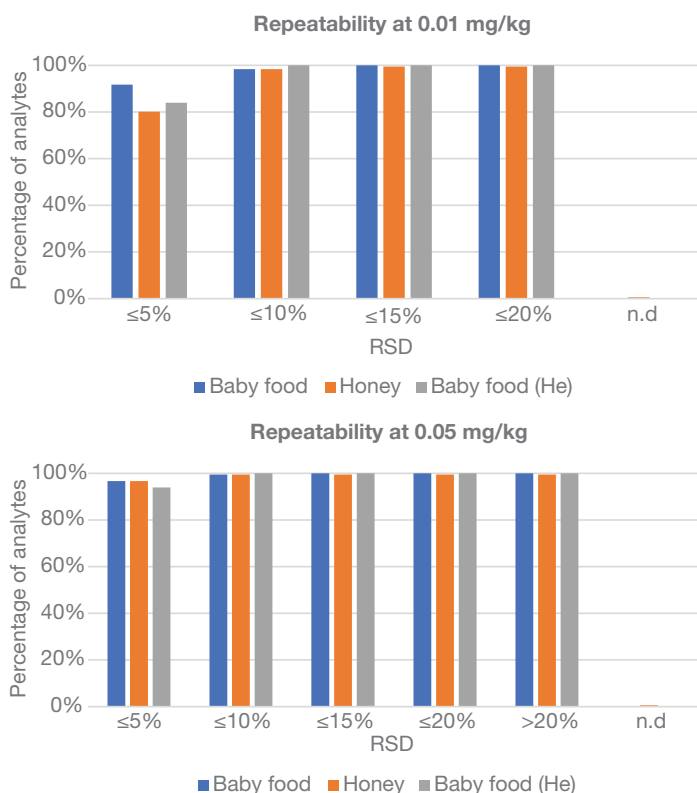


Figure 4. Repeatability (n = 5) of 181 pesticides in two matrices (baby food and honey) with hydrogen as carrier gas, compared with repeatability of pesticides in baby food using carrier helium

Implication on gas consumption

Although the data presented in this study clearly demonstrates compliance with SANTE guidelines, the use of hydrogen as carrier gas may not be the ideal solution in all cases. Optimum performance is often only achieved with helium. The use of the HeSaver-H₂Safer technology offers an alternative to the use of carrier hydrogen, especially for GC-MS methods. Thanks to the limited carrier gas consumption of a few mL/min, it offers the flexibility to maintain helium as the carrier gas with significant savings, mitigating helium shortage issues.

Helium gas consumption with a standard SSL injector and the GC-MS method described in this note is 1.3 liters per sample. However, this volume can be reduced to 0.32 liters per sample when the HeSaver-H₂Safer mode is used. Thus, the helium consumption with the HeSaver-H₂Safer mode would be four times lower than with the standard SSL injector. If the instrument was running 365 days per year and 24 hours per day, one helium cylinder would last almost 2 years. The [Thermo Scientific™ Helium Saver Calculator](#) tool⁴ offers an easy-to-use and intuitive interface to estimate helium consumption and cost impact on individual laboratories activities. GC parameters regarding column dimensions, carrier gas and split flow settings, as well as helium and nitrogen costs are adjustable to reflect a given laboratory's methodology and regional gas cost to provide estimates on helium cylinder lifetime and cost savings. It is worth highlighting that the use of the HeSaver-H₂Safer technology does not affect sensitivity or retention times and does not involve any method re-optimization when switching from a standard SSL injector.⁵

Conclusions

It has been demonstrated that the TSQ 9610 GC-MS/MS is capable of identifying and quantifying pesticides at 0.01 mg/kg concentration level when hydrogen is used as the carrier gas. The proposed method provides a viable alternative for laboratories performing the testing for pesticide residues in a wide range of food matrices and looking at ways to circumvent helium price increases or potential supply shortages. While the use of hydrogen may have a negative impact on detection sensitivity (and hence the ability to meet established detection limit requirements), the use of the AEI ion source together with a few changes in the choice of transitions can offset these and allow to achieve comparable performance at reduced running costs.

Key findings of the application are summarized in the following:

- The use of the HeSaver-H₂Safer technology allows for a safe and compliant use of hydrogen as an alternative carrier gas in GC/GC-MS applications without the need to install a hydrogen sensor and removes any risk of unwanted reactions with the sample in the hot SSL injector. Additionally, it offers the flexibility to maintain helium as the carrier gas with a significant savings, with the benefit to avoid method re-optimization while mitigating helium shortage issues.
- When used with hydrogen, the limited carrier gas consumption offered by the HeSaver-H₂Safer mode permits a very controlled system demand for hydrogen, making this solution ideal for laboratories working with hydrogen generators.
- Migrating the GC-MS/MS method from helium to hydrogen requires an adaptation of method parameters to address (i.e., retention time shift and different fragmentation patterns).
- All target pesticides met the regulatory RSD requirement of less than 20%. The relative standard deviation for N=5 repeats of each sample type at 0.01 mg/kg was equal or lower than 10% for 98% of the evaluated compounds.

- 97% of pesticides in baby food and 98% in honey showed a linear response in the concentration range 0.005–0.500 mg/kg.
- For laboratories requiring best detection sensitivity, or searching for non-target pesticides, the SSL upgraded to HeSaver-H₂Safer technology allows the use of carrier helium with a reduced consumption by a factor of four, maintaining the method.

References

1. The Science Of Helium And Why Global Supplies Are Running Low: [Short Wave : NPR](#)
2. Analytical Quality Control and Method Validation Procedures for Pesticide Residues Analysis In Food And Feed, SANTE 11312/2021.
3. Thermo Fisher Scientific, Application Note 000437: Confident analysis of ultra-trace pesticides residues in baby food using triple quadrupole GC-MS. <https://assets.thermofisher.com/TFS-Assets/CMD/Application-Notes/an-000437-gc-ms-pesticides-baby-food-an000437-en.pdf>
4. Thermo Fisher Scientific Helium Saver Calculator. <https://www.thermofisher.com/it/en/home/industrial/chromatography/chromatography-learning-center/chromatography-consumables-resources/chromatography-tools-calculators/helium-saver-calculator.html>
5. Thermo Fisher Scientific, Product Spotlight SL001922: Reduction of helium usage for the analysis of pesticide residues in botanical dietary supplements using GC-MS/MS with the HeSaver-H₂Safer technology. <https://assets.thermofisher.com/TFS-Assets/CMD/brochures/sl-001922-gc-ms-ms-hesaver-h2safer-botanical-pesticides-sl001922-en.pdf>

Appendix 1. List of the target compounds, retention times, and monitored transitions

Part 1

Pesticide	RT [min]	Window [min]	Precursor mass 1	Product mass 1	Collision energy 1	Precursor mass 2	Product mass 2	Collision energy 2	Precursor mass 3	Product mass 3	Collision energy 3
2,3,5,6-Tetrachloroaniline	8.72	0.3	230.8	157.9	18	230.8	159.8	18			
2,4'-Methoxychlor	15.65	0.3	121.1	78	20	197.1	165.1	15	227.1	121	10
3,4-Dichloroaniline	7.66	0.3	160.9	99	20	160.9	126	10			
4,4'-Methoxychlor olefin	15.19	0.3	238.1	152.1	34	238.1	223.1	10			
Acetochlor	10.35	0.3	146.1	130	24	223.1	132	20			
Acrinathrin	19.38	0.3	181	152	22	208.1	180.9	8	289	93.1	8
Alachlor	10.48	0.3	188.1	130	32	188.1	160.1	8			
Aldrin	11.16	0.3	254.9	219.9	20	262.7	192.9	32			
Allidochlor	6.67	0.3	132	56.1	8	134	56	8			
Anthraquinone	11.18	0.3	208.1	152	22	208.1	180.1	10			
Atrazine	9.44	0.3	200.1	122	8	215.1	58.1	10			
Azinphos-ethyl	18.95	0.3	132	77	12	160	77	16			
Azinphos-methyl	17.88	0.3	132	77	12	160	50.9	34	160	77	16
Benfluralin	8.98	0.3	292	160	20	292	264	8			
BHC, Alpha	9.13	0.3	180.9	144.9	14	216.9	181	8			
BHC, Beta	9.42	0.3	216.9	180.9	8	218.9	183	8			
BHC, delta	9.89	0.3	180.9	144.9	14	218.9	182.9	8			
BHC, gamma	9.53	0.3	180.9	144.9	12	218.9	183	8			
Bifenthrin	17.11	0.3	165.1	163.6	24	181	165.9	10	181	179	12
Biphenyl	7.38	0.3	154.1	115	26	155.1	154.1	12			
Bromfenvinphos-methyl	12.04	0.3	294.9	79.1	30	294.9	109	16			
Bromophos-ethyl	12.47	0.3	302.8	284.8	14	358.8	302.8	14			

Appendix 1. List of the target compounds, retention times, and monitored transitions

Part 2

Pesticide	RT [min]	Window [min]	Precursor mass 1	Product mass 1	Collision energy 1	Precursor mass 2	Product mass 2	Collision energy 2	Precursor mass 3	Product mass 3	Collision energy 3
Bromophos-methyl (Bromophos)	11.56	0.3	328.9	313.8	14	330.8	315.8	14			
Bromopropylate	16.91	0.3	184.9	156.9	12	340.8	185	14			
Bupirimate	13.6	0.3	273.1	108	14	273.1	193.2	8			
Carbophenothion	15.1	0.3	157	45	12	342	157	10			
Carfentrazon-ethyl	15.2	0.3	290	99.9	36	311.9	150.7	18	340.1	312.1	10
Chlorbenside	12.42	0.3	125	89	16	268	125	10			
Chlordane alpha-cis	12.71	0.3	271.9	236.9	14	374.8	265.9	20			
Chlordane gamma-trans	12.41	0.3	271.8	236.8	12	372.8	265.9	14			
Chlorfenapyr	13.83	0.3	248.9	112	24	327.9	246.9	14			
Chlorfenson	13.01	0.3	111	75	14	175	111	10			
Chlorobenzilate	14.19	0.3	139	111	12	251	139	14			
Chloroneb	8.11	0.3	190.9	113	14	193	53.1	32			
Chlorpropham	8.9	0.3	127	65	20	171	127	8	213	127	14
Chlorpyrifos-ethyl	11.17	0.3	198.9	171	14	313.9	257.9	12			
Chlorpyrifos-methyl	10.35	0.3	285.9	92.9	20	287.9	92.9	20			
Chlorthal-dimethyl (Dacthal)	11.25	0.3	300.9	222.9	22	300.9	272.9	12			
Chlorthiophos	14.53	0.3	268.9	205	14	324.9	268.9	12			
Chlozolate	12.01	0.3	186	145	14	188	147	14	259	187.9	12
Clomazone	9.47	0.3	125	89	16	204	107	18			
Coumaphos	20.11	0.3	209.9	119	22	209.9	182	10	226	163	18
Cycloate	8.79	0.3	154.1	55.1	18	154.1	83.1	8			
Cyhalothrin I (lambda)	18.85	0.3	180.9	152	22	197.1	141.1	10	207.9	180.9	8
Cyprodinil	11.83	0.3	224.1	197.1	20	224.1	208.1	18			
DDD p,p	14.35	0.3	235	165	20	235	199	14	237	165	20
DDD, o, p	13.45	0.3	235	165	20	237	165	20			
DDE o,p	12.54	0.3	246	176.1	28	248	176.1	30			
DDE p, p	13.32	0.3	246	176.1	28	315.9	246	14			
DDT o,p	14.4	0.3	212	151	30	212	176	25			
DDT p,p	15.38	0.3	235	165.1	22	236.8	165	22			
Deltamethrin	24.66	0.3	252.8	92.9	16	252.8	172	8			
Diazinon	9.71	0.3	137.1	84.1	12	199	92.9	14			
Dichlobenil	7.06	0.3	170.9	99.9	24	170.9	136	12			
Dichlorobenzophenone, 4, 4	11.42	0.3	139	74.9	26	141	113	10			
Dicloran (Bortran)	9.28	0.3	160	124.1	8	206	176	10			
Dieldrin	13.32	0.3	262.9	192.9	30	262.9	227.8	16			
Diphenamid	11.59	0.3	167.1	165.1	20	239.1	167.1	8			
Diphenylamine	8.73	0.3	168.1	167.1	14	169.1	168.1	12			
Disulfoton	9.82	0.3	88	59.8	6	142	81	10			
Edifenphos	15.12	0.3	172.9	109	8	310	109	26			
Endosulfan ether	10.14	0.3	238.9	204	12	240.9	206	14			
Endosulfan alpha	12.7	0.3	194.9	160	8	240.8	205.8	14			
Endosulfan beta	14.09	0.3	158.9	123	12	194.9	159	8			
Endosulfan sulfate	15.13	0.3	238.7	203.9	12	271.7	236.8	12			

Appendix 1. List of the target compounds, retention times, and monitored transitions

Part 3

Pesticide	RT [min]	Window [min]	Precursor mass 1	Product mass 1	Collision energy 1	Precursor mass 2	Product mass 2	Collision energy 2	Precursor mass 3	Product mass 3	Collision energy 3
Endrin	13.81	0.3	244.9	173	22	262.8	192.9	30	280.8	244.9	8
Endrin Aldehyde	14.48	0.3	249.8	214.9	24	278.9	242.9	10			
Endrin-Ketone	16.37	0.3	316.8	208.9	28	316.8	281	10			
EPN	16.85	0.3	157	77	22	169	77	22	169	141	8
Ethalfuralin	8.98	0.3	264	160	12	292	160	18	292	264	8
Ethion	14.46	0.3	230.9	128.9	22	230.9	174.9	12			
Etofenprox	22.24	0.3	163.1	107.1	16	163.1	135.1	10			
Fenarimol	18.68	0.3	139	74.9	26	139	111	14	219	107	10
Fenchlorfos	10.63	0.3	284.9	269.9	14	286.9	271.9	14			
Fenitrothion	10.87	0.3	277	109	14	277	260	6			
Fenpropathrin	17.33	0.3	181	126.8	28	181	151.9	22			
Fenson	11.51	0.3	141	50.9	30	141	77	8			
Fenthion	11.24	0.3	278	109	18	278	169	14			
Fenvalerate	23.25	0.7	125	89	18	167	125	10			
Fipronil	11.97	0.3	366.9	212.9	28	368.9	214.9	30			
Fluazifop-P-butyl	14.12	0.3	282.1	91.1	18	282.1	238.1	16			
Fluchloralin	9.74	0.3	306	264	8	326	63	12			
Fludioxonil	13.16	0.3	248	127	26	248	154	18			
Fluquinconazole	20.12	0.3	340	108.1	36	340	298	16			
Fluridone	22.51	0.7	328.1	189.1	38	328.1	258.8	24			
Flusilazole	13.53	0.3	233.1	151.9	14	233.1	164.9	16			
Flutolanil	13.11	0.3	173	95	28	281	173	10			
Flutriafol	12.86	0.3	123	75	24	219.1	95	34	219.1	123	12
Fluvalinate peak 1	23.65	1	250	55.1	16	250	199.9	18			
Fluvalinate peak 2	23.8	1	250	55.1	16	250	200	16			
Fonofos	9.64	0.3	137	109	6	246	109	14			
Heptachlor	10.56	0.3	271.8	236.8	12	273.8	238.8	14			
Hexachlorobenzene	9.17	0.3	283.8	213.8	30	283.8	248.8	16			
Hexazinone	15.55	0.3	171.1	71.1	14	171.1	85.1	12			
Iodofenfos	13	0.3	376.8	361.8	16	378.8	363.8	14			
Iprodione	16.72	0.3	314	245	10	315.7	247	10			
Isazophos	9.86	0.3	118.9	76	18	256.9	161.9	4			
Isodrin	11.7	0.3	146.9	111.1	10	192.9	123	28			
Isopropalin	11.73	0.3	280.1	180.2	10	280.1	238.2	8			
Lenacil	15.27	0.3	153	82.1	16	153	135.6	12			
Leptophos	17.86	0.3	171	77.1	18	171	124.3	10			
Linuron	10.96	0.3	187	124	20	248	61.1	8			
Malathion	11.07	0.3	127	99	6	158	125	6			
Metalaxyl	10.6	0.3	234.1	146.1	20	234.1	174.1	10			
Metazachlor	11.82	0.3	132.1	117.1	14	133.1	132.1	12	209	132.1	16
Methacrifos	8.07	0.3	124.9	47.1	12	207.9	180.1	6			
Methoxychlor	17.13	0.3	227.1	141.1	32	227.1	169.1	22			
Metolachlor	11.12	0.3	238.1	133.1	26	238.1	162.1	10			

Appendix 1. List of the target compounds, retention times, and monitored transitions

Part 4

Pesticide	RT [min]	Window [min]	Precursor mass 1	Product mass 1	Collision energy 1	Precursor mass 2	Product mass 2	Collision energy 2	Precursor mass 3	Product mass 3	Collision energy 3
Mevinphos	7.65	0.3	127	95	14	192	127	10			
MGK-264 A	11.63	0.3	164.1	80.1	24	164.1	93.1	10	164.1	98.1	10
MGK-264 B	11.88	0.3	164.1	67.1	6	164.1	98.1	12			
Mirex	18.19	0.3	272	236.8	14	273.8	238.8	14			
Myclobutanil	13.44	0.3	179	90	28	179	125	14			
N-(2,4-Dimethylphenyl) formamide	7.96	0.3	149.1	106.1	16	149.1	120.1	14	149.1	121.1	6
Nitralin	16.27	0.3	274	216	8	316.2	274	8			
Nitrofen	13.88	0.3	202	139	24	283	162	20			
Nonachlor-cis	14.26	0.3	406.8	299.9	14	410.8	301.8	14			
Nonachlor-trans	12.79	0.3	271.8	236.8	14	406.8	299.8	14			
Norflurazon	15.19	0.3	145	95	16	303	145	20			
Ortho-phenylphenol	8.17	0.3	170.1	115	34	170.1	141.1	22			
Oxadiazon	13.44	0.3	175	112	12	258	175	6			
Oxyfluorfen	13.62	0.3	252	146	30	300	223	14			
Paclobutrazol	12.6	0.3	138	103.1	14	236	125	12	236	167	10
Parathion (ethyl)	11.31	0.3	154.9	125	6	291	109	12			
Parathion-methyl	10.45	0.3	233	109	10	263	109	12			
Pebulate	7.86	0.3	128	72	6	128.1	57.1	8			
Penconazole	11.93	0.3	159	123	20	248.1	157	22			
Pendimethalin	11.82	0.3	252.1	161.1	14	252.1	162.1	8	252.1	191.3	8
Pentachloroaniline	10.12	0.3	262.9	191.9	20	266.9	193.9	20			
Pentachloroanisole	9.24	0.3	264.8	236.9	10	279.9	236.8	22			
Pentachlorobenzene	8.16	0.3	247.9	212.9	18	249.8	214.8	16			
Pentachlorobenzonitrile	9.51	0.3	272.9	237.9	16	274.8	239.9	18			
Pentachloroethoxyanisole	10.91	0.3	295.7	245.9	30	295.7	262.9	12			
Permethrin	20.2	0.6	183.1	165.1	10	184.1	91.1	20	184.1	141.1	15
Perthane (Ethylan)	14.01	0.3	223.1	167	12	223.1	179	20			
Phenothrin	18.01	0.3	123.1	41.1	24	123.1	79.1	14	123.1	81.1	8
Phorate	9.09	0.3	121	65	10	260	75	8			
Phosalone	17.85	0.3	182	74.8	30	182	111	14			
Phosmet	16.65	0.3	160	76.9	22	160	133	10			
Piperonyl butoxide	16.26	0.3	176.1	103.1	22	176.1	131.1	12			
Pirimiphos-ethyl	11.65	0.3	304.1	168.1	12	318.1	166.1	12			
Pirimiphos-methyl	10.86	0.3	233	151.1	6	290.1	125	20			
Pretilachlor	13.22	0.3	162	132.1	20	202.1	174.2	8	262.1	202.1	6
Prochloraz	20.27	0.3	180	69	14	180.1	138.1	12	308	70	12
Procymidone	12.22	0.3	283	68.1	24	283	96.1	8			
Prodiamine	10.93	0.3	275.1	255.1	8	321.1	203	10	321.1	279.1	6
Profenofos	13.23	0.3	336.9	266.9	12	338.9	268.9	14			
Profluralin	9.58	0.3	318.1	55	12	318.1	199.1	12	347.1	330.1	6
Propanil	10.29	0.3	161	90	24	161	99	24	217	161	8
Propargite	16.03	0.3	135.1	77.1	26	135.1	107.1	12	150.1	135.1	8
Propisochlor	10.56	0.3	162.1	120.1	12	162.1	144.1	8			

Appendix 1. List of the target compounds, retention times, and monitored transitions

Part 5

Pesticide	RT [min]	Window [min]	Precursor mass 1	Product mass 1	Collision energy 1	Precursor mass 2	Product mass 2	Collision energy 2	Precursor mass 3	Product mass 3	Collision energy 3
Propyzamide	9.66	0.3	172.9	109	24	172.9	145	14	174.9	147	14
Prothiofos	13.12	0.3	266.9	220.9	18	309	238.9	14			
Pyraclufos	19.41	0.3	194	138	18	360	194.1	12			
Pyrazophos	18.99	0.3	221	148.7	14	221	193.1	8			
Pyridaben	20.19	0.3	147.1	117.1	20	147.1	119.1	8	147.1	132.1	12
Pyridaphenthion	16.62	0.3	199	92.1	14	340	199.1	8			
Pyrimethanil	9.74	0.3	198.1	118	32	198.1	158.1	18			
Pyriproxyfen	18.31	0.3	136.1	96	10	226.1	186.1	12			
Quinalphos	12.15	0.3	146	118.1	10	157.1	102	22			
Quintozene	9.46	0.3	213.8	178.9	14	294.8	236.9	14			
Sulfotep	8.99	0.3	237.9	145.9	12	322	145.9	22			
Sulprofos	14.86	0.3	156	108	30	322	156.1	10			
Tebuconazole	15.78	0.3	125	89	16	250	125	20			
Tebufenpyrad	17.47	0.3	276.1	171	10	318.1	131.1	14			
Tecnazene	8.54	0.3	214.8	178.9	8	258.9	201	12			
Tefluthrin	9.91	0.3	177	127	14	177	137	16			
Terbacil	9.85	0.3	160	76	12	161	144	12			
Terbufos	9.6	0.3	231	128.9	20	231	175	10			
Terbutylazine	9.61	0.3	214.1	104.1	16	214.1	132.1	10			
Tetrachlorvinphos	12.62	0.3	328.9	109	18	330.9	109	18	332.9	109	14
Tetradifon	17.63	0.3	159	74.8	32	159	111	20			
Tetrahydrophthalimide (THPI)	8.03	0.3	151	79.9	6	151	122.1	8			
Tolclofos-methyl	10.46	0.3	265	219.9	20	265	250	12			
Triadimefon	11.37	0.3	208	111	20	208	126.7	12			
Triadimenol	12.24	0.3	128	65	18	128	100	10			
Triallate	9.94	0.3	86.1	43.3	6	268	183.9	18	268	226	12
Triazophos	14.86	0.3	161.1	134.1	8	257	162.1	6			
Tricyclazole	13.02	0.3	189	161	14	189	162	10			
Triflumizole	12.34	0.3	206	179	14	206	186	8			
Trifluralin	8.95	0.3	264	160	14	306.1	264.1	8			
Vinclozolin	10.43	0.3	186.8	124	18	212	172	14			

Appendix 2. Details of repeatability and linearity evaluation

Part 1

Pesticide	Repeatability (n = 5)						Linear range [mg/kg]	
	0.01 mg/kg			0.05 mg/kg			Baby food	Honey
	Baby food	Honey	Baby food (He)	Baby food	Honey	Baby food (He)		
2,3,5,6-Tetrachloroaniline	2%	4%	5%	4%	4%	5%	0.005–0.500	0.005–0.500
2,4'-Methoxychlor	6%	4%	5%	4%	4%	7%	0.005–0.500	0.005–0.500
3,4-Dichloroaniline	13%	-	9%	9%	-	8%	0.005–0.200	–
4,4'-Methoxychlor olefin	1%	1%	4%	1%	2%	4%	0.005–0.500	0.005–0.500
Acetochlor	4%	4%	3%	4%	2%	2%	0.005–0.500	0.005–0.500
Acrinathrin	5%	3%	6%	3%	1%	5%	0.005–0.500	0.005–0.500
Alachlor	3%	3%	5%	3%	1%	3%	0.005–0.500	0.005–0.500
Aldrin	3%	2%	3%	2%	4%	3%	0.005–0.500	0.005–0.500
Allidochlor	5%	4%	5%	4%	4%	4%	0.005–0.500	0.005–0.500
Anthraquinone	2%	3%	4%	3%	4%	3%	0.005–0.500	0.005–0.500
Atrazine	4%	5%	6%	5%	1%	2%	0.005–0.500	0.005–0.500
Azinphos-ethyl	1%	4%	7%	4%	5%	5%	0.005–0.500	0.005–0.500
Azinphos-methyl	5%	3%	6%	3%	3%	5%	0.005–0.200	0.005–0.500
Benfluralin	6%	3%	4%	3%	4%	3%	0.005–0.500	0.005–0.500
BHC, Alpha	2%	2%	4%	2%	3%	3%	0.005–0.500	0.005–0.500
BHC, Beta	2%	2%	4%	2%	2%	3%	0.005–0.500	0.005–0.500
BHC, delta	3%	1%	3%	1%	2%	2%	0.005–0.500	0.005–0.500
BHC, gamma	4%	2%	4%	2%	3%	3%	0.005–0.500	0.005–0.500
Bifenthrin	6%	2%	4%	2%	3%	4%	0.005–0.500	0.005–0.500
Biphenyl	4%	5%	4%	5%	2%	8%	0.005–0.500	0.005–0.500
Bromfenvinphos-methyl	3%	2%	5%	2%	2%	4%	0.005–0.500	0.005–0.500
Bromophos-ethyl	1%	1%	3%	1%	3%	3%	0.005–0.500	0.005–0.500
Bromophos-methyl (Bromophos)	2%	2%	4%	2%	3%	3%	0.005–0.500	0.005–0.500
Bromopropylate	3%	3%	5%	3%	3%	5%	0.005–0.500	0.005–0.500
Bupirimate	1%	2%	6%	2%	3%	4%	0.005–0.500	0.005–0.500
Carbophenothion	2%	2%	5%	2%	1%	4%	0.005–0.500	0.005–0.500
Carfentrazone-ethyl	4%	2%	5%	2%	2%	4%	0.005–0.500	0.005–0.500
Chlorbenside	2%	1%	4%	1%	2%	4%	0.005–0.500	0.005–0.500
Chlordane alpha-cis	2%	1%	5%	1%	3%	4%	0.005–0.500	0.005–0.500
Chlordane gamma-trans	1%	1%	5%	1%	3%	4%	0.005–0.500	0.005–0.500
Chlorfenapyr	5%	2%	3%	2%	5%	5%	0.005–0.500	0.005–0.500
Chlorfenson	2%	1%	5%	1%	1%	4%	0.005–0.500	0.005–0.500
Chlorobenzilate	3%	1%	4%	1%	3%	4%	0.005–0.500	0.005–0.500
Chloroneb	2%	5%	3%	5%	4%	5%	0.005–0.500	0.005–0.500
Chlorpropham	3%	2%	3%	2%	3%	2%	0.005–0.500	0.005–0.500
Chlorpyrifos-ethyl	2%	1%	3%	1%	3%	2%	0.005–0.500	0.005–0.500
Chlorpyrifos-methyl	4%	1%	4%	1%	2%	3%	0.005–0.500	0.005–0.500
Chlorthal-dimethyl (Dacthal)	1%	1%	6%	1%	3%	3%	0.005–0.500	0.005–0.500
Chlorthiophos	1%	1%	4%	1%	2%	4%	0.005–0.500	0.005–0.500
Chlozolate	3%	2%	4%	2%	2%	4%	0.005–0.500	0.005–0.500
Clomazone	3%	2%	4%	2%	2%	3%	0.005–0.500	0.005–0.500
Coumaphos	3%	5%	5%	5%	2%	4%	0.005–0.500	0.005–0.500
Cycloate	3%	4%	4%	4%	4%	6%	0.005–0.500	0.005–0.500

Appendix 2. Details of repeatability and linearity evaluation

Part 2

Pesticide	Repeatability (n = 5)						Linear range [mg/kg]	
	0.01 mg/kg			0.05 mg/kg			Baby food	Honey
	Baby food	Honey	Baby food (He)	Baby food	Honey	Baby food (He)		
Cyhalothrin I (lambda)	3%	4%	5%	4%	3%	5%	0.005–0.500	0.005–0.500
Cyprodinil	4%	2%	3%	2%	1%	4%	0.005–0.500	0.005–0.500
DDD p,p	5%	3%	3%	3%	4%	8%	0.005–0.500	0.005–0.500
DDD, o, p	4%	2%	3%	2%	3%	4%	0.005–0.500	0.005–0.500
DDE o,p	1%	3%	2%	3%	2%	3%	0.005–0.500	0.005–0.500
DDE p, p	1%	2%	4%	2%	2%	4%	0.005–0.500	0.005–0.500
DDT o,p	11%	12%	5%	12%	5%	3%	0.005–0.500	0.005–0.500
DDT p,p	14%	10%	6%	10%	6%	7%	0.005–0.500	0.005–0.500
Deltamethrin	4%	6%	4%	6%	2%	5%	0.005–0.500	0.005–0.500
Diazinon	3%	2%	3%	2%	2%	2%	0.005–0.500	0.005–0.500
Dichlobenil	4%	5%	5%	5%	5%	8%	0.005–0.500	0.005–0.500
Dichlorobenzophenone, 4, 4	1%	3%	2%	3%	4%	3%	0.005–0.500	0.005–0.500
Dicloran (Bortran)	3%	2%	2%	2%	2%	1%	0.005–0.500	0.005–0.500
Dieldrin	2%	1%	6%	1%	6%	4%	0.005–0.500	0.005–0.500
Diphenamid	3%	2%	4%	2%	2%	3%	0.005–0.500	0.005–0.500
Diphenylamine	3%	4%	5%	4%	2%	3%	0.005–0.500	0.005–0.500
Disulfoton	6%	2%	4%	2%	2%	1%	0.005–0.500	0.005–0.500
Edifenphos	2%	5%	3%	5%	2%	5%	0.005–0.500	0.005–0.500
Endosulfan ether	2%	3%	3%	3%	3%	3%	0.005–0.500	0.005–0.500
Endosulfan alpha	6%	1%	3%	1%	3%	3%	0.005–0.500	0.005–0.500
Endosulfan beta	3%	1%	3%	1%	2%	5%	0.005–0.500	0.005–0.500
Endosulfan sulfate	3%	2%	6%	2%	3%	4%	0.005–0.500	0.005–0.500
Endrin	4%	1%	6%	1%	3%	5%	0.005–0.500	0.005–0.500
Endrin Aldehyde	4%	1%	4%	1%	4%	4%	0.005–0.500	0.005–0.500
Endrin-Ketone	6%	4%	6%	4%	5%	5%	0.005–0.500	0.005–0.500
EPN	2%	3%	4%	3%	3%	5%	0.005–0.500	0.005–0.500
Ethalfuralin	3%	3%	5%	3%	4%	4%	0.005–0.500	0.005–0.500
Ethion	2%	2%	5%	2%	1%	4%	0.005–0.500	0.005–0.500
Etofenprox	3%	1%	5%	1%	1%	4%	0.005–0.500	0.005–0.500
Fenarimol	3%	3%	5%	3%	2%	4%	0.005–0.500	0.005–0.500
Fenchlorfos	2%	1%	3%	1%	2%	3%	0.005–0.500	0.005–0.500
Fenitrothion	5%	1%	4%	1%	2%	4%	0.005–0.500	0.005–0.500
Fenpropathrin	4%	1%	4%	1%	4%	4%	0.005–0.500	0.005–0.500
Fenson	2%	2%	4%	2%	2%	3%	0.005–0.500	0.005–0.500
Fenthion	4%	2%	3%	2%	1%	3%	0.005–0.500	0.005–0.500
Fenvalerate	2%	4%	6%	4%	2%	5%	0.005–0.500	0.005–0.500
Fipronil	5%	1%	5%	1%	2%	4%	0.005–0.500	0.005–0.500
Fluazifop-P-butyl	2%	2%	5%	2%	1%	4%	0.005–0.500	0.005–0.500
Fluchloralin	4%	2%	8%	2%	4%	3%	0.005–0.500	0.005–0.500
Fludioxonil	1%	1%	3%	1%	4%	4%	0.005–0.500	0.005–0.500
Fluquinconazole	4%	3%	5%	3%	2%	4%	0.005–0.500	0.005–0.500
Fluridone	4%	3%	4%	3%	5%	5%	0.005–0.500	0.005–0.500
Flusilazole	2%	4%	2%	4%	2%	4%	0.005–0.500	0.005–0.500

Appendix 2. Details of repeatability and linearity evaluation

Part 3

Pesticide	Repeatability (n = 5)						Linear range [mg/kg]	
	0.01 mg/kg			0.05 mg/kg			Baby food	Honey
	Baby food	Honey	Baby food (He)	Baby food	Honey	Baby food (He)		
Flutolanil	2%	1%	5%	1%	1%	3%	0.005–0.500	0.005–0.500
Flutriafol	4%	1%	5%	1%	3%	4%	0.005–0.500	0.005–0.500
Fluvalinate	3%	4%	4%	4%	1%	5%	0.005–0.500	0.005–0.500
Fonofos	3%	2%	3%	2%	2%	2%	0.005–0.500	0.005–0.500
Heptachlor	3%	3%	6%	3%	4%	3%	0.005–0.500	0.005–0.500
Hexachlorobenzene	3%	5%	6%	5%	5%	7%	0.005–0.500	0.005–0.500
Hexazinone	1%	1%	4%	1%	4%	5%	0.005–0.200	0.005–0.500
Iodofenfos	1%	1%	3%	1%	5%	4%	0.005–0.500	0.005–0.500
Iprodione	4%	3%	3%	3%	2%	4%	0.005–0.500	0.005–0.500
Isazophos	5%	3%	5%	3%	3%	2%	0.005–0.500	0.005–0.500
Isodrin	2%	2%	4%	2%	4%	3%	0.005–0.500	0.005–0.500
Isopropalin	3%	1%	3%	1%	3%	4%	0.005–0.500	0.005–0.500
Lenacil	5%	1%	4%	1%	3%	5%	0.005–0.500	0.005–0.500
Leptophos	1%	2%	4%	2%	2%	4%	0.005–0.500	0.005–0.500
Linuron	5%	2%	8%	2%	3%	5%	0.005–0.500	0.005–0.500
Malathion	3%	3%	4%	3%	1%	3%	0.005–0.500	0.005–0.500
Metaxalyl	3%	2%	5%	2%	3%	4%	0.005–0.500	0.005–0.500
Metazachlor	3%	2%	5%	2%	2%	4%	0.005–0.500	0.005–0.500
Methacrifos	3%	4%	3%	4%	3%	5%	0.005–0.500	0.005–0.500
Methoxychlor	9%	8%	5%	8%	6%	7%	0.005–0.500	0.005–0.500
Metolachlor	2%	1%	3%	1%	1%	3%	0.005–0.500	0.005–0.500
Mevinphos	3%	4%	4%	4%	3%	4%	0.005–0.500	0.005–0.500
MGK-264 A	4%	2%	2%	2%	1%	3%	0.005–0.500	0.005–0.500
MGK-264 B	3%	1%	3%	1%	1%	4%	0.005–0.500	0.005–0.500
Mirex	1%	1%	3%	1%	5%	5%	0.005–0.500	0.005–0.500
Myclobutanil	3%	1%	5%	1%	3%	4%	0.005–0.500	0.005–0.500
N-(2,4-Dimethylphenyl) formamide	5%	1%	3%	1%	5%	3%	0.005–0.500	0.005–0.500
Nitralin	2%	4%	7%	4%	2%	2%	0.005–0.500	0.005–0.500
Nitrofen	3%	1%	3%	1%	2%	5%	0.005–0.500	0.005–0.500
Nonachlor-cis	1%	1%	5%	1%	3%	5%	0.005–0.500	0.005–0.500
Nonachlor-trans	2%	2%	3%	2%	3%	3%	0.005–0.500	0.005–0.500
Norflurazon	2%	1%	3%	1%	3%	2%	0.005–0.500	0.005–0.500
Ortho-phenylphenol	3%	3%	5%	3%	3%	3%	0.005–0.500	0.005–0.500
Oxadiazon	1%	2%	5%	2%	2%	3%	0.005–0.500	0.005–0.500
Oxyfluorfen	1%	2%	5%	2%	2%	5%	0.005–0.500	0.005–0.500
Paclobutrazol	2%	2%	8%	2%	1%	4%	0.005–0.500	0.005–0.500
Parathion (ethyl)	5%	3%	2%	3%	3%	4%	0.005–0.500	0.005–0.500
Parathion-methyl	3%	2%	2%	2%	2%	4%	0.005–0.500	0.005–0.500
Pebulate	3%	5%	6%	5%	6%	9%	0.005–0.500	0.005–0.500
Penconazole	2%	1%	4%	1%	3%	3%	0.005–0.500	0.005–0.500
Pendimethalin	2%	1%	6%	1%	3%	4%	0.005–0.500	0.005–0.500
Pentachloroaniline	2%	4%	4%	4%	3%	2%	0.005–0.500	0.005–0.500
Pentachloroanisole	2%	4%	4%	4%	4%	5%	0.005–0.500	0.005–0.500

Appendix 2. Details of repeatability and linearity evaluation

Part 4

Pesticide	Repeatability (n = 5)						Linear range [mg/kg]	
	0.01 mg/kg			0.05 mg/kg			Baby food	Honey
	Baby food	Honey	Baby food (He)	Baby food	Honey	Baby food (He)		
Pentachlorobenzene	6%	8%	8%	8%	7%	7%	0.005–0.500	0.005–0.500
Pentachlorobenzonitrile	4%	4%	5%	4%	4%	2%	0.005–0.500	0.005–0.500
Pentachloroethoxybenzene	2%	4%	4%	4%	4%	3%	0.005–0.500	0.005–0.500
Permethrin	7%	2%	3%	2%	3%	4%	0.005–0.500	0.005–0.500
Perthane (Ethylan)	1%	1%	5%	1%	3%	4%	0.005–0.500	0.005–0.500
Phenothrin	5%	1%	5%	1%	2%	3%	0.005–0.500	0.005–0.500
Phorate	4%	3%	5%	3%	2%	2%	0.005–0.500	0.005–0.500
Phosalone	3%	4%	5%	4%	1%	5%	0.005–0.500	0.005–0.500
Phosmet	4%	5%	6%	5%	2%	5%	0.005–0.500	0.005–0.500
Piperonyl butoxide	1%	2%	5%	2%	3%	4%	0.005–0.500	0.005–0.500
Pirimiphos-ethyl	2%	1%	4%	1%	2%	3%	0.005–0.500	0.005–0.500
Pirimiphos-methyl	4%	2%	4%	2%	1%	2%	0.005–0.500	0.005–0.500
Pretilachlor	4%	1%	3%	1%	4%	4%	0.005–0.500	0.005–0.200
Prochloraz	5%	3%	8%	3%	5%	3%	0.005–0.500	0.005–0.500
Procymidone	6%	2%	5%	2%	1%	4%	0.005–0.500	0.005–0.500
Prodiamine	1%	3%	6%	3%	2%	5%	0.005–0.500	0.005–0.500
Profenofos	5%	2%	7%	2%	5%	5%	0.005–0.500	0.005–0.200
Profluralin	3%	1%	8%	1%	4%	4%	0.005–0.500	0.005–0.500
Propanil	2%	1%	4%	1%	3%	3%	0.005–0.500	0.005–0.500
Propargite	5%	2%	6%	2%	3%	4%	0.005–0.500	0.005–0.500
Propisochlor	3%	2%	4%	2%	2%	4%	0.005–0.500	0.005–0.500
Propyzamide	3%	2%	5%	2%	3%	3%	0.005–0.500	0.005–0.500
Prothiofos	3%	1%	4%	1%	1%	4%	0.005–0.500	0.005–0.500
Pyraclufos	2%	3%	5%	3%	4%	5%	0.005–0.500	0.005–0.500
Pyrazophos	7%	1%	5%	1%	1%	5%	0.005–0.200	0.005–0.500
Pyridaben	5%	4%	5%	4%	2%	4%	0.005–0.500	0.005–0.500
Pyridaphenthion	1%	3%	4%	3%	1%	5%	0.005–0.500	0.005–0.500
Pyrimethanil	3%	3%	2%	3%	2%	4%	0.005–0.500	0.005–0.500
Pyriproxyfen	3%	1%	5%	1%	2%	4%	0.005–0.500	0.005–0.500
Quinalphos	4%	1%	4%	1%	1%	4%	0.005–0.500	0.005–0.500
Quintozene	5%	3%	5%	3%	2%	5%	0.005–0.500	0.005–0.500
Sulfotep	5%	3%	2%	3%	2%	2%	0.005–0.500	0.005–0.500
Sulprofos	2%	1%	4%	1%	1%	4%	0.005–0.500	0.005–0.500
Tebuconazole	2%	2%	6%	2%	2%	4%	0.005–0.500	0.005–0.500
Tebufenpyrad	2%	2%	5%	2%	3%	4%	0.005–0.500	0.005–0.500
Tecnazene	3%	4%	4%	4%	4%	3%	0.005–0.500	0.005–0.500
Tefluthrin	2%	3%	3%	3%	3%	2%	0.005–0.500	0.005–0.500
Terbacil	4%	1%	5%	1%	1%	4%	0.005–0.500	0.005–0.500
Terbufos	4%	3%	3%	3%	1%	2%	0.005–0.500	0.005–0.500
Terbuthylazine	4%	2%	3%	2%	2%	2%	0.005–0.500	0.005–0.500
Tetrachlorvinphos	3%	3%	5%	3%	2%	4%	0.005–0.500	0.005–0.500
Tetradifon	3%	1%	4%	1%	2%	3%	0.005–0.500	0.005–0.500
Tetrahydrophthalimide (THPI)	6%	2%	3%	2%	2%	4%	0.005–0.500	0.005–0.500

Appendix 2. Details of repeatability and linearity evaluation

Part 5

Pesticide	Repeatability (n = 5)						Linear range [mg/kg]	
	0.01 mg/kg			0.05 mg/kg			Baby food	Honey
	Baby food	Honey	Baby food (He)	Baby food	Honey	Baby food (He)		
Tolclofos-methyl	3%	1%	2%	1%	2%	2%	0.005–0.500	0.005–0.500
Triadimefon	3%	0%	3%	0%	1%	3%	0.005–0.500	0.005–0.500
Triadimenol	3%	3%	5%	3%	3%	3%	0.005–0.500	0.005–0.500
Triallate	2%	2%	4%	2%	2%	3%	0.005–0.500	0.005–0.500
Triazophos	3%	3%	5%	3%	2%	4%	0.005–0.200	0.005–0.500
Tricyclazole	4%	1%	5%	1%	3%	5%	0.005–0.200	0.005–0.500
Triflumizole	2%	2%	3%	2%	2%	3%	0.005–0.500	0.005–0.500
Trifluralin	3%	3%	4%	3%	4%	5%	0.005–0.500	0.005–0.500
Vinclozolin	1%	4%	4%	4%	4%	3%	0.005–0.500	0.005–0.500

 Learn more at thermofisher.com/helium-saver