# Pesticide Residue Analyses of QuEChERS Extracts of Different Food Matrices Using an Online Robotic SPE **Clean-up Procedure Coupled to LC-MS/MS**

## F. Michael Hudson; Cristina C. Jacob; and Claudia P.B. Martins, Thermo Fisher Scientific, 355 River Oaks Pkwy, San Jose, CA 95134, USA

## **ABSTRACT**

**Purpose:** A simple and fast method for the clean-up of QuEChERS extracts, coupled with LC-MS/MS for multi-residue pesticide analysis in food matrices.

**Methods:** A robotic solid-phase extraction (SPE) using miniaturized cartridges on a CTC autosampler (TriPlus RSH with RTC) was coupled to Thermo Fisher Vanguish™ Flex Binary UHPLC Svstem interfaced with a Thermo Fisher TSQ Fortis<sup>™</sup> Triple Quadrupole Mass Spectrometer equipped with a Thermo Scientific<sup>™</sup> Heated Electrospray Ionization (HESI-II) probe operating in both positive and negative ionization modes.

**Results:** Improvements in analytical performance are here reported for an LC-MS/MS method for the multi-residue analysis of pesticides from food matrices after implementing a fully automated and online µSPE clean-up method for processing QuEChERS extracts.

## INTRODUCTION

Pesticides have been widely used to control pests and their residue analyses in crops and food products are routinely performed in regulatory and industrial laboratories. QuEChERS method is commonly used for the extraction of pesticide residues from food samples<sup>1</sup>. Despite its high pesticide extraction efficiency, fatty and pigmented matrix is often coextracted along with the compounds of interest when QuEChERS method is employed. This propensity for matrix coextraction can cause problems in the LC-MS analysis-such as ion suppression-and therefore clean-up is beneficial. We report herein a fully automated and online µSPE clean-up method of QuEChERS extracts with direct LC-MS/MS analysis.

## MATERIALS AND METHODS

#### **TriPlus RSH System Configuration**

• Triplus RSH configuration enables the use of dedicated syringes for μSPE extracts clean-up steps and LC/MS injection, and is also equipped with reservoirs for conditioning, washing and eluting solvents (Fig.1)

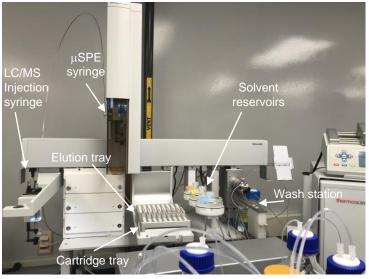


Figure 1. TriPlus RSH µSPE clean-up workflow setup

• μSPE Cartridges are sealed by a septum above the sorbent bed which allows the syringe to push sample extracts or solvents through the sorbent bed. Therefore, the syringe replaces the vacuum system of the classical SPE methodology with the added benefit of working at defined flow rates (Fig.

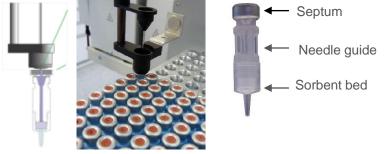


Figure 2. Cartridge needle transport and ITSP µSPE cartridges

• The small dimensions allows the scale-down of the μSPE clean-up workflow (low microliter level)

#### **uSPE workflow**

- (TriPlus RSH) with integrated injection to LC-MS/MS analysis (Table 1)
- are used as base sorbent
- Elution solvent 1:1 Acetonitrile/ Methanol + 100 mM ammonium formate (pH 5.8)
- pesticide standard mixture.
- allows continuous sample analysis within the cycle of the LC-MS system.

#### Table 1. QuEChERS extract automated clean-up workflow

1. Clean	Prep Syringe wit
2. Condition	$\mu SPE$ with 100 $\mu$
3. Move	$\mu SPE$ cartridge t
4. Load	100 μL QuEChE
5. Clean	The prep syringe
6. Elute	$\mu$ SPE cartridge v
7. Move	Used $\mu SPE$ cart
8. Change	To LC/MS injecti
9. Inject	To LC-MS/MS
10. Change	To prep syringe
11. Proceed	With prep-ahead Signal

#### LC-MS/MS system

equipped with a HESI ionization probe.

LC Conditions				
Analytical Column Thermo Accucore AQ, 2.1 x 100 mm, 2.6 $\mu m$	Gradient:	Time (min)	% Solvent B	
Flow rate 300 µL/min		0	0	
Column temperature 25 °C		0.5	0	
<b>Solvent A</b> Water containing 2mM ammonium formate, 2% methanol and 0.1% FA		7	70	
<b>Solvent B</b> Methanol containing 2mM ammonium formate, 2% water and 0.1%FA		9	100	
Injection volume 1 μL		12	100	
		12.1	0	
		15	0	

MS Source Parameters			
HESI Source: both positive and negative modes	Vaporizer temperature : 350 °C		
Spray Voltage: Positive 3.7 kV and Negative 2.5 kV	Cycle time :0.4 s		
Shealth Gas (Arb): 30	Q1 Resolution (FWHM): 0.7 Da		
Auxiliary Gas (Arb): 6	Q3 Resolution (FWHM): 1.2 Da		
Sweep Gas (Arb): 1	CID gas: 1.5 mTorr		
lon transfer tube temp.: 300 °C			

#### Data Analysis

The acquired data were processed using Thermo Scientific<sup>™</sup> TraceFinder<sup>™</sup> 4.1 software.

Robotic Solid-Phase Extraction (SPE) using miniaturized cartridges on a CTC autosampler

• ITSP SPE Cartridges QuEChERS Blend for LC; where C18/zirconia-coated silica and Carbon-X

• Blank matrix samples (Leek, avocado, and walnut) were purchased in local retail stores. The homogenized samples were then extracted with QuEChERS protocols and spiked with a 270

• The clean-up cycle is completed in less than 10 min. It is scheduled as a prep-ahead task which

- ith elution solvent
- μL elution solvent
- e to elution tray
- ERS extract onto µSPE cartridge
- e with 100 μL elution solvent
- tridge to cartridge tray
- tion syringe

e for next sample

ad for next extract sample upon Ready

#### Vanguis Flex Binary pump interfaced with a Fortis Triple Quadrupole Mass Spectrometer

## RESULTS

Leek QuEChERS extracts

- µSPE procedure resulted in clear extracts (efficient decolorization, Fig. 3)
- µSPE clean-up provided recoveries in the range 60-120% for most of the compounds (Fig.3)
- Precision <20% was achieve for most of the monitored compounds (Fig. 3)





Figure 3. Leek extract without clean-up and after µSPE clean-up

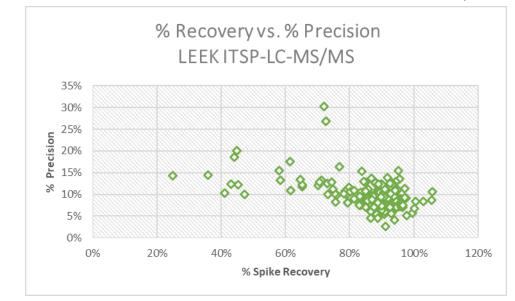


Figure 4. %Spike Recovery vs. % precision of Leek QuEChERS extracts

Matrix effects

78%

74%

73%

59%

59%

55%

53%

48%

47%

44%

44%

44%

40%

39%

39%

37%

36%

36%

33%

29%

33%

48% 32%

after mSPE

#### Avocado QuEChERS extracts

Table 2. Reduction od matrix effects for certain compounds

Selected Compounds

Brodifacoum

Etofenprox

Resmethrin

Fenpyroximat

Fenazaquin

Flufenoxuron

Bifenazate

chlorfluazuror

Clethodim

Profenophos

Tetramethrin

Tebufenpyrad

Penconazole

Tolfenpyrad

Hexythiazox

Furathiocarb

Propiconazole

Hexaconazole

Imibenconazole

Tebuconazole

Mexacarbate

Sulprofos

✓ High oil content and color which are removed by μSPE clean-up

Matrix effects

efore mSPE

90%

81%

79%

65%

65%

58%

94%

56%

54%

46%

40%

44%

46%

44%

44%

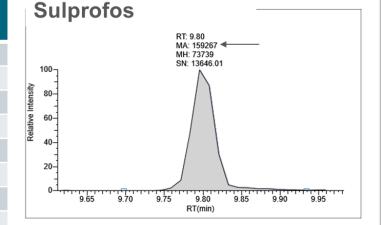
40%

48%

52%

49%

46%



### Sulprofos

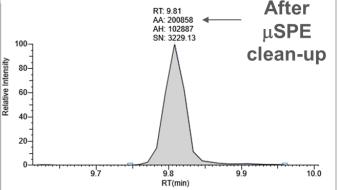
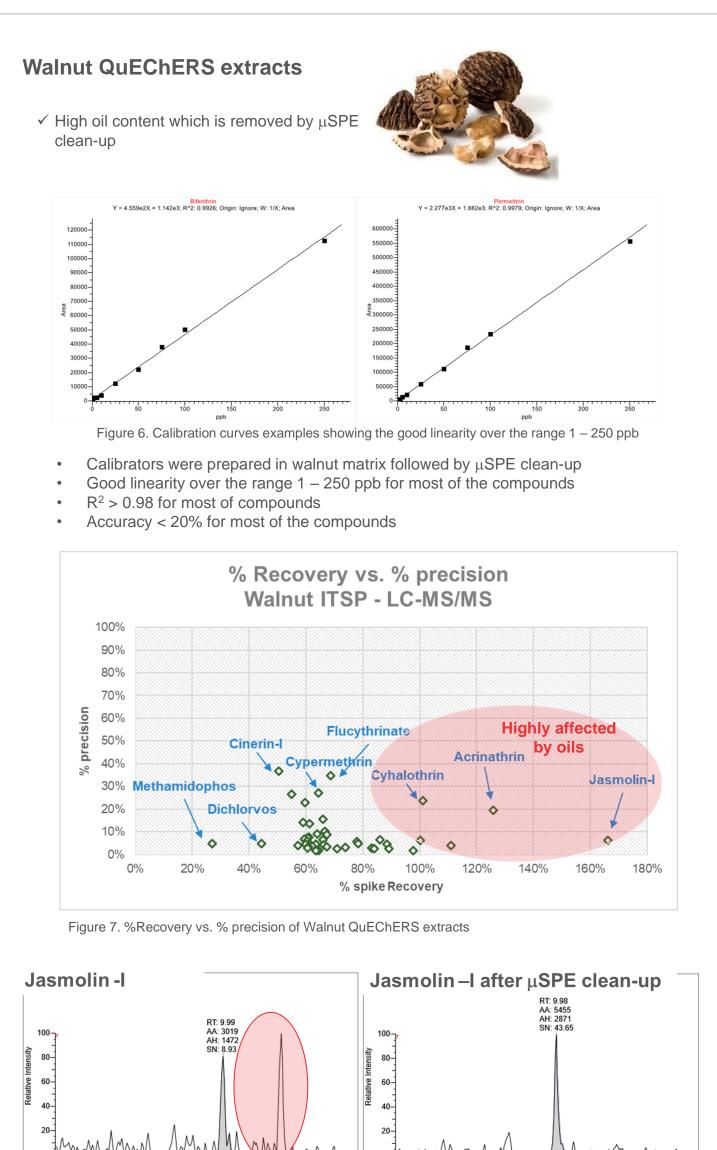


Figure 5. Sulprofos Extracted Ion Chromatograms showing better signal after  $\mu$ SPE clean-up.



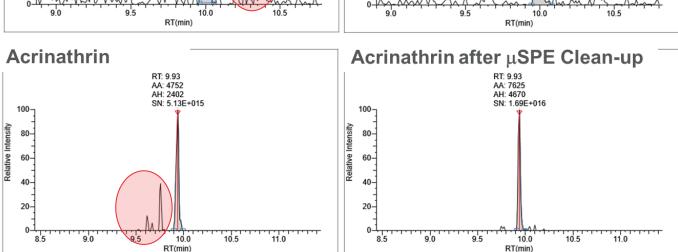


Figure 8. Extracted ion chromatograms showing the removal of non-polar matrix components

## CONCLUSIONS

Fully automated and online µSPE clean-up procedure with LC-MS/MS for the pesticide multi-residue analysis in food matrices was developed and evaluated.

- Average pesticides recoveries were in the range of 60 to 120% with RSDs <30% in different food</p> matrices (Leek, avocado and walnut)
- μSPE procedure improved removal of matrix components and analytical performance
- The developed μSPE workflow is reliable, robust and suitable for LC-MS/MS analyses
- The automated μSPE procedure uses the idle time between injections, therefore LC-MS/MS run times are not affected
- Lab efficiency is improved by removing the need for manual clean-up procedure

## REFERENCES

1. Anastassiades et al. Fast and easy multiresidue method employing acetonitrile extraction/partitioning and "dispersive solid-phase extraction" for the determination of pesticide residues in produce. J. Chromatogr. A. 2003. 1015, 163-184.

## ACKNOWLEDGEMENTS

We would like to thank Tom Flug and Brian Peat from CTC Analytics for the firmware, scripts, communications board, software, and training to operate the robotic liquid handler.

## **TRADEMARKS/LICENSING**

© 2019 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.

PO65484-EN0419S



