

Poster Reprint

ASMS 2023
Poster number MP 216

Automated sample preparation using CTC PAL3 to analyze >570 pesticides in an orange by the combination of GC/MS/MS and LC/MS/MS techniques

Zhiming, Zhang; Ge, Meng; Zhicong, Wang; Jianguo, Ji;
Chunxiao, Wang, Lorna De-Leoz

Agilent Technologies, Inc., Shanghai, P.R. China

Introduction

Automated sample preparation can not only increase testing efficiency, but also reduces dependency on in-person activity. For example many lab chemists had to stay at home due to COVID-19 pandemic, resulting in a lot of sample testing been delayed due to the shortage of chemist availability. Compared with manual work, an automated workflow, especially coupled online to mass spectrometry, is less manpower reliant and a best choice to tackle this kind of situation.

The manual sample preparation process for QuEChERS includes solvent extraction, salting and cleaning, with multi-time of shaking/vortex and centrifugation. This process avoids dependency on chemist time and increases the precision of the results due to reducing human error.

An automated workflow was developed to analyze multi-class pesticides in orange with the combination of GC/MS/MS and LC/MS/MS techniques. Sample preparation was automatically done with CTC PAL3 on the basis of QuEChERS, which ensured consistent and high quality results. Worth mention, using scripts, the sample preparation was done in overlapped mode, avoiding additional sample prep overhead time.

Validations were done at three different spiking levels (10µg/Kg, 20µg/Kg, 50µg/Kg) with 5 replicates for each level. The validation results for automatic sample preparation and manual sample preparation were summarized and compared.

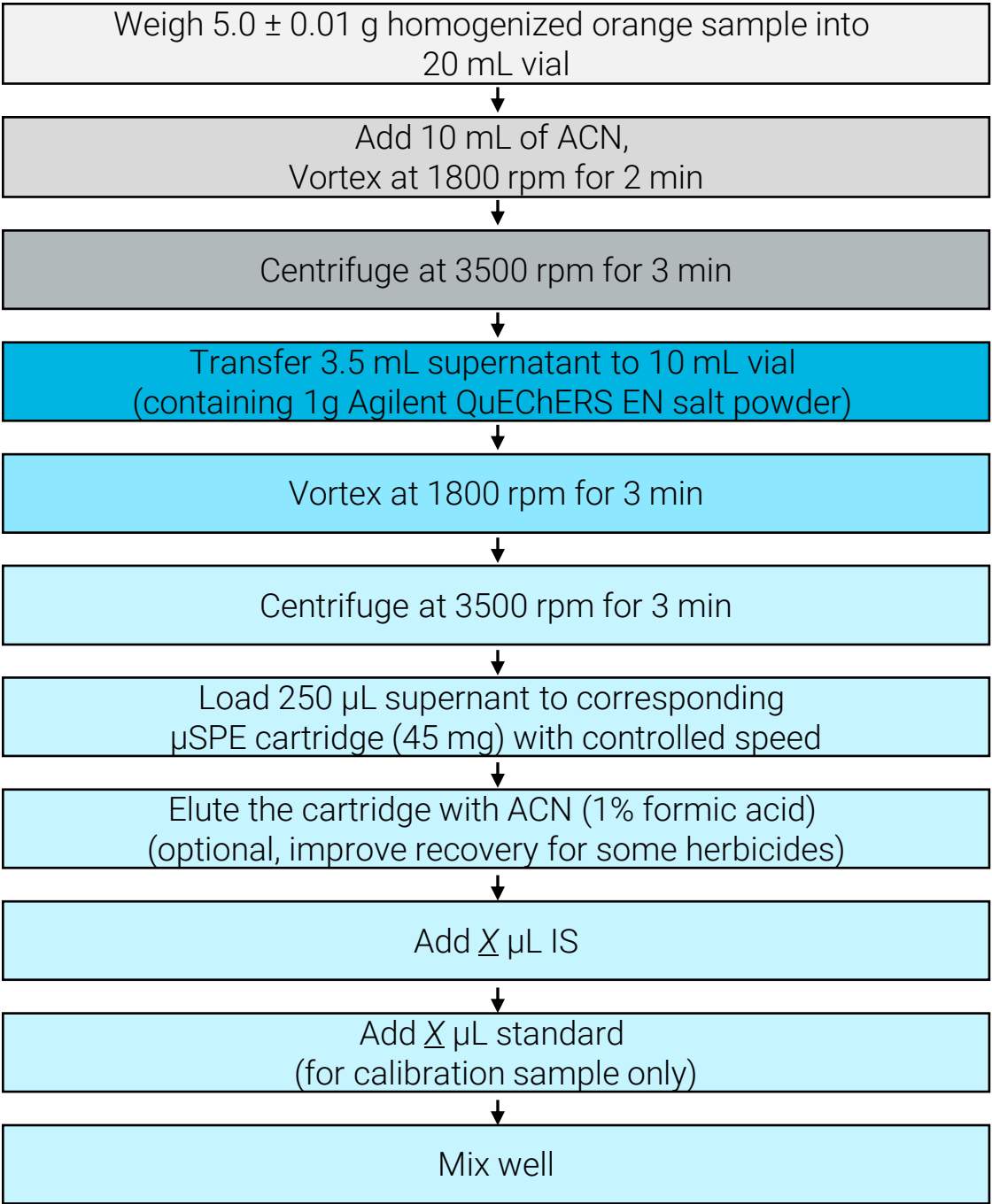


1290Infinity II LC + 6470

Figure 1. Instruments used in the experiment

Experimental

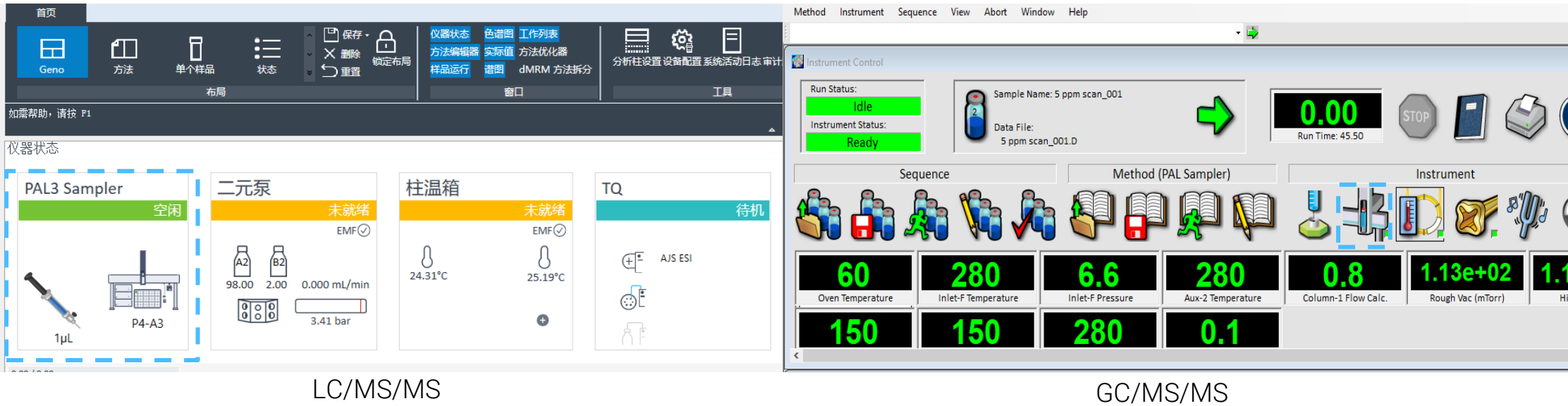
Automatic sample Prep.



Partial parameters of PAL3 auto sample prep.

⌵ Conditioning Conditioning Solvent Source none Conditioning Solvent Index 2 Conditioning Solvent Volume 100 µL Conditioning Solvent Fill Speed 10 µL/s	⌵ Internal Standard Internal Standard Source none Internal Standard Index 54 Internal Standard Volume 7.7 µL Internal Standard Fill Speed 1 µL/s
⌵ Sample µSPE µSPE Sample Load Volume 257 µL µSPE Sample Fill Speed 5 µL/s	⌵ Solvent Addition Solvent Source none Solvent Index 1 Solvent Volume 16.45 µL Solvent Fill Speed 5 µL/s
⌵ Elution Elution Solvent Source none Elution Solvent Index 2 Elution Volume 100 µL Elution Solvent Fill Speed 5 µL/s	⌵ Target Standard Target Standard Source none Target Standard Index 52 Target Standard Volume 3.85 µL Target Standard Fill Speed 1 µL/s
⌵ Protectants Protectant Source none Protectant Index 3 Protectant Volume 0 µL Protectant Fill Speed 5 µL/s	⌵ Mixing Mix Cycles 5 Mix Volume 500 µL
	⌵ Injection

Instrument control panel



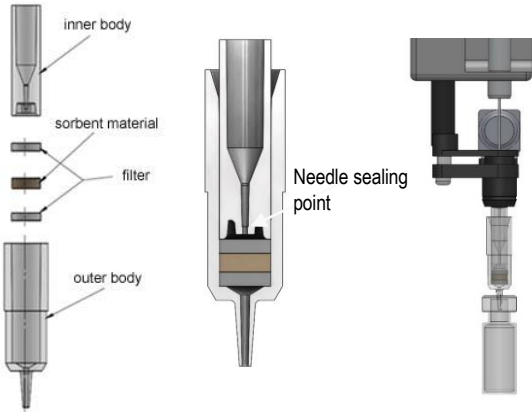
Picture: liquid tools



Picture: centrifuge

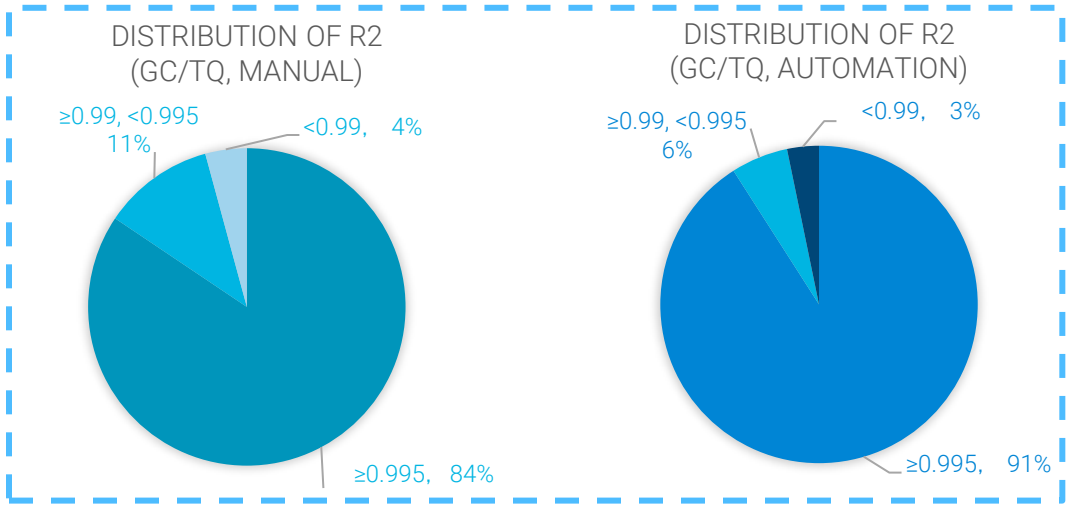
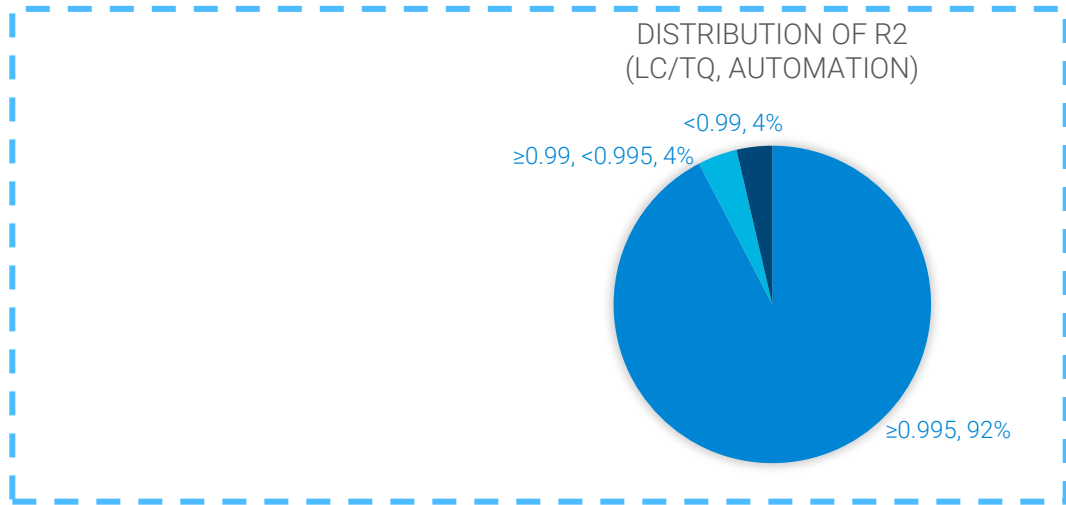


Picture: vortex

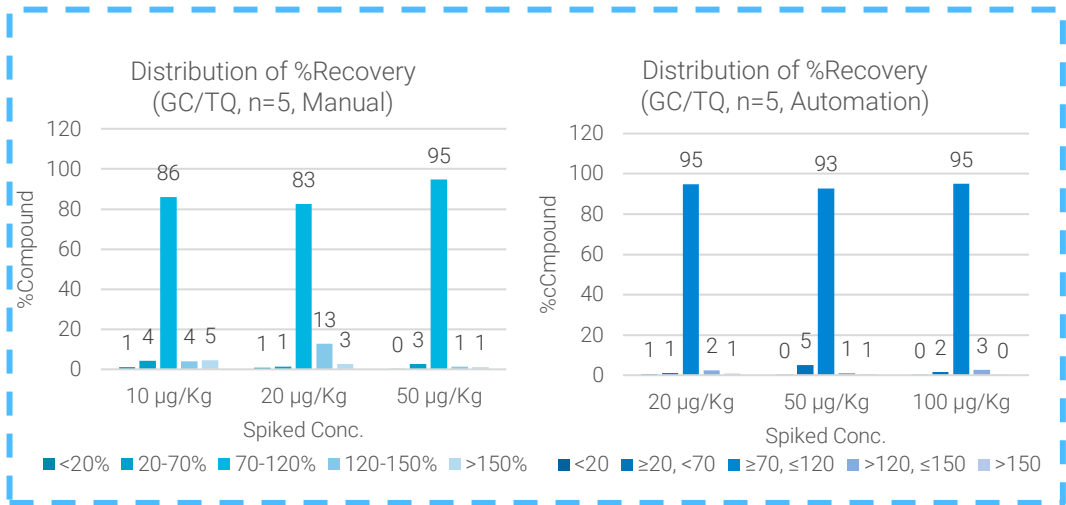
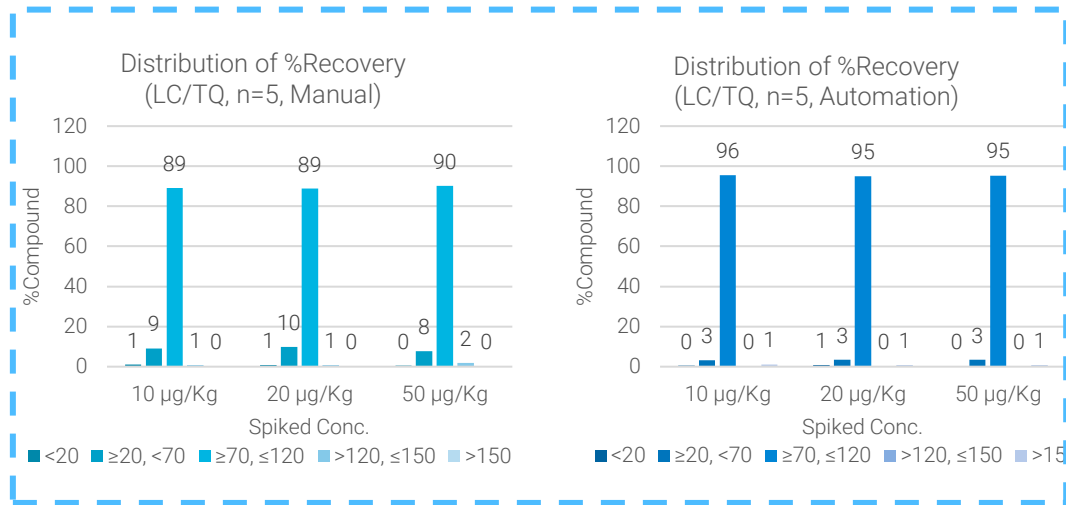


Picture: Design of μSPE

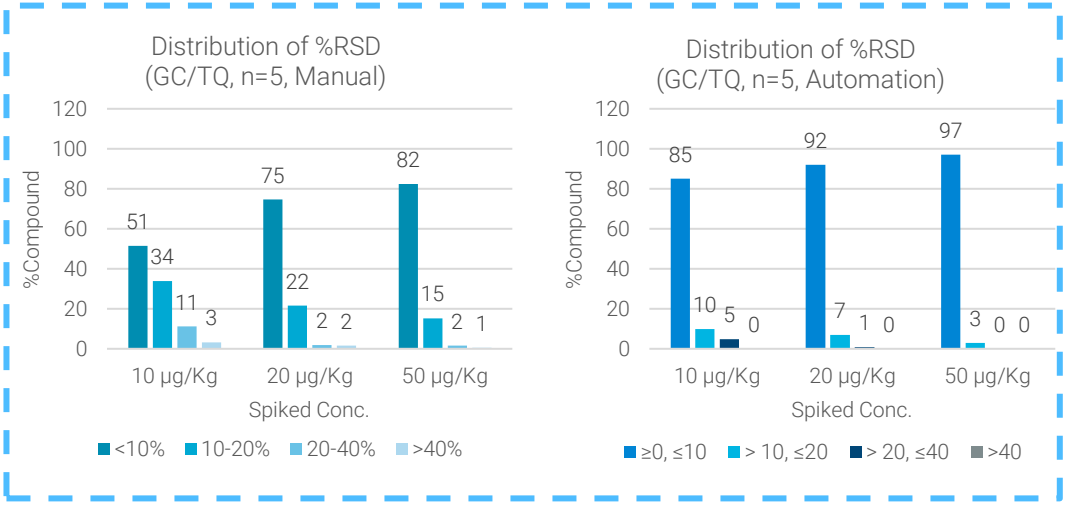
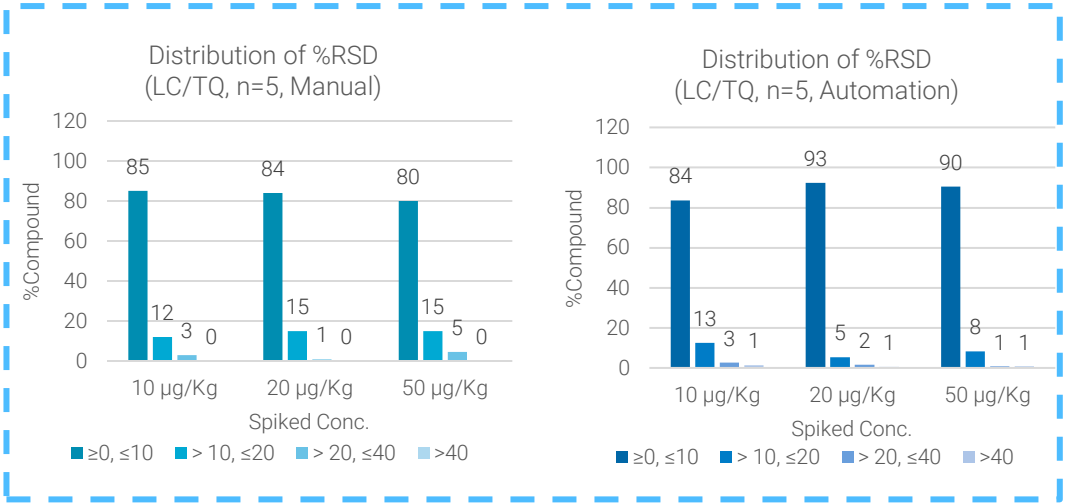
Linearity (manual vs. automation)



Accuracy (manual vs. automation)



Accuracy (manual vs. automation)



Conclusions

Full automation, high efficiency, low cost

This is the first time to develop a whole automatic sample preparation workflow for analyzing > 570 pesticides in orange by the combination of GC/MS/MS and LC/MS/MS techniques.

In general, automatic workflow has better performance than manual workflow by comparison of linearity, accuracy and repeatability of validation results done by manual and automated sample preparation.

References

- GB2763-2021: National food safety standard- Maximum residue limits for pesticides in food.
- GB23200.121-2021: National food safety standard- Determination of 331 pesticides and metabolites residues in foods of plant origin-LC/MS/MS
- GB23200.121-2018: National food safety standard- Determination of 208 pesticides and metabolites residues in foods of plant origin-GC/MS/MS