thermoscientific



A quantitative determination of pesticide residues in chili powder using GC-MS/MS

Authors

Sarvendra P. Singh, Subodh K. Budakoti, and Dasharath P. Oulkar Customer Solution Center, Thermo Fisher Scientific, Ghaziabad, India

Keywords

Quantitation, pesticides, chili powder, TraceFinder, GC-MS/MS, QuEChERS, TSQ 9000, ExtractaBrite

Goal

The objective of the experiments described here was to set up a complete analytical solution to enable commercial food testing laboratories to analyze pesticide residues in chili powder, in compliance with the requirements of the Food Safety and Standards Authority of India (FSSAI).

Introduction

Chili (*Capsicum annum* Linn.), an essential spice in Indian cuisine, is grown by farmers on a large scale, i.e., about 800,000 hectares in India. The states of Andhra Pradesh and Karnataka account for a substantial share of chili production in comparison with the other states of India. The chili crop is vulnerable to a multitude of pests such as chili thrips, scirtothriops, dorsalis hood, and yellow mite. Consequently pesticides are widely used to maintain crop quality and yield. To date, relatively few pesticides have been registered under the Central Insecticide Board and Registration Committee (CIBRC) for use on chili.¹ Current agriculture practices rely on various pesticides applications to control the insect and pest attack, but these are harmful to human health. The frequent use of these chemicals can result in the residues in the harvested crop,² which is a food safety concern. Therefore, monitoring of pesticide residues in chili is necessary. Effective residue monitoring requires appropriate sample preparation for optimum extraction efficiency, accuracy, and precision.



Goon et al.³ have reported the QuEChERS method for the extraction of pesticides residue analysis in spices including chili followed by LC-MS analysis. This technique can also be applied for GC-MS/MS amenable pesticides.

Endosulfan-
$$\alpha$$
Endosulfan- β

Endosulfan-sulfate

Dimethoate

Figure 1. Structure of, endosulfan- α , endosulfan- β , endosulfan-sulfate, dicofol, and dimethoate

The aim of this work was to develop and demonstrate the sample preparation and determination of targeted pesticide residues in chili powder. Data were acquired using the Thermo Scientific™ TSQ™ 9000 GC-MS/MS (gas chromatograph-triple quadrupole mass spectrometer). The optimized method was validated at LOQ and LOQ × 2 in terms of the accuracy and precision

as per the SANTE/11813/2017 guidelines and assessed for FSSAI MRL compliance. This method was also applied to the analysis of real samples for comprehensive method evaluation.

Experimental

Chemicals, reagents, and apparatus

- Acetonitrile, Optima[™] LC/MS Grade, Fisher Chemical (P/N 514 L-16923 U)
- Anhydrous Magnesium Sulfate, Thermo Scientific (P/N 80020-415-500)
- Sodium Acetate, Thermo Scientific (P/N 80020-424)
- C18 Octadecyl Endcapped, Thermo Scientific (P/N 80020-413-100)
- PSA (Primary, secondary amine), Thermo Scientific (P/N 80020-416-100)
- GCB (Graphitized carbon black), Thermo Scientific (P/N 80020-417-100)
- The reference standards (dimethoate, dicofol, endosulfan-alfa, endosulfan beta, and endosulfan sulfate)
- Sample handling equipment: weighing balance (analytical and precision), benchtop centrifuge for 50 mL and 15 mL tubes, vortex mixer, micro-pipettes

Standard solutions preparations

Individual stock standard solutions were purchased from Sigma-Aldrich, Bangalore, India, and prepared by dissolving 10 mg of each in 10 mL acetonitrile corresponding to 1000 μ g/mL. These were used to prepare a mixed working solution of the pesticides at 10 μ g/mL in acetonitrile. The working standard solutions given in Table 1 were prepared by serial dilution of the 10 μ g/mL standard.

Table 1. Matrix-matched calibration standards preparation

Working standard (µg/mL)	Volume from working standard (μL)	Acetonitrile (μL)	Final concentration (mg/kg)	Total volume (μL)
4.000	50	950	0.200	1000
3.000	50	950	0.150	1000
2.000	50	950	0.100	1000
1.500	50	950	0.075	1000
1.000	50	950	0.050	1000
0.500	50	950	0.025	1000
0.200	50	950	0.010	1000
0.100	50	950	0.005	1000

Sample preparation

Chili powder (finely homogenized powder) was purchased from a local retail outlet.

Extraction

- Weigh 2 g homogenized chili powder sample into a 50 mL extraction tube. Note: Spike sample at this step and wait for 10 min.
- Add 15 mL of 1% acetic acid in the water, shake, and soak for 10 min.
- Add 15 mL acetonitrile and mix vigorously for 1 min on a vortex mixer at 2500 rpm.
- Add 6 g anhydrous MgSO₄ and 1.5 gm of sodium acetate, again mix vigorously for 1 min on a vortex mixer.
- Centrifuge at 4000 rpm for 5 min at ambient temperature.

Dispersive solid phase extraction (dSPE) cleanup

- Decant 5 mL supernatant into 15 mL centrifuge tube and add 750 mg MgSO4, 250 mg PSA, 250 mg C18, and 50 mg of GCB.
- Vortex for 30 s and centrifuge at 10,000 rpm for 5 min.
- Collect 2.5 mL cleaned supernatant in the separate glass tube.
- Evaporate, to near dryness, under a gentle stream of nitrogen gas and then reconstitute with 1 mL acetonitrile.
- Inject 2 µL to the GC-MS/MS through the Thermo Scientific™ TriPlus™ RSH™ Autosampler.

Preparation of matrix-matched calibration standards

The blank matrix was analyzed for assuring the incurred residues of target analytes. After assuring the absence of target analytes, the same matrix was used for the further study. Following the above protocol, 2.5 mL cleaned blank matrix extract was evaporated and used to prepare matrix match calibration standards at the concentrations given in Table 1.

Recovery and precision

The mixed working standards were spiked in chili powder before extraction at 0.025 and 0.050 mg/kg (n=6 for each).

GC-MS/MS analysis

Chromatographic separation and data acquisition were carried out using the Thermo Scientific™ TRACE™ 1310 gas chromatograph coupled with a Thermo Scientific™ TSQ™ 9000 triple quadrupole mass spectrometer. The optimized instrument parameters for each compound are given in Table 2, while the compound-dependent parameters, i.e., selected reaction monitoring (SRM) parameters, are presented in Table 3.

Table 2A. Instrument parameters (GC)

Gas chromatogra	aphy method				
Instrumentation:	TRACE 1310 GC with TSQ 9000 triple quadrupole and				
	TriPlus RSH autosampler				
Column:	Thermo Scientific™ TraceGOLD™				
	TG-5MS GC column				
	$(30 \text{ m} \times 0.25 \text{ mm i.d.} \times 0.25 \mu\text{m})$				
	P/N 26096-1420				
Liner:	Siltek [™] six baffle PTV liner				
India at a m	(P/N 453T2120)				
Injector:	Programmed Temperature Vaporizing Injector (PTV)				
Injector mode:	Splitless (PTV)				
Splitless time:	2 min				
Injection volume:	2 μL				
PTV program:	90 °C for 0.1 min;				
	14.5 °C/min, 90-300 °C;				
	2.0 min (hold);				
	14.5 °C/min, 300–320 °C;				
	5 min (hold). Flow 75.0 mL/min				
Cleaning phase:	On				
Flow control mode:	Constant Flow				
Column flow:	1.20 mL/min				
Carrier gas					
and purity:	Helium (99.999%)				
Purge flow:	5.00 mL/min				
Split flow:	50.00 mL/min				
Post-cycle temp.:	Cool Down				
Total run time:	24.83 min				
GC oven program:	70 °C, 1 min (hold);				
	30 °C/min, 70–150 °C;				
	5 °C/min, 150–200 °C;				
	10 °C/min, 200–280 °C;				
	30 °C/min, 280-300 °C, 5 min (hold)				

Table 2B. Instrument parameters (MS)

Mass spectrometry method					
Instrumentation:	TSQ 9000 triple quadrupole				
	mass spectrometer with				
	ExtractaBrite ion source				
Method type:	Acquisition-General (SRM mode)				
MS transfer line					
temperature:	310 °C				
lon source					
temperature:	250 °C				
lonization:	El (Electron Ionization)				

Data acquisition and processing

Thermo Scientific™ TraceFinder™ 4.1 software was used for instrument control, data acquisition, and processing, data review, and reporting. The data was acquired in SRM mode with a minimum of two SRM transitions per analyte. The data processing included user-defined criteria as; two transitions per analytes, retention time (±0.1 min) and ion ratio (±30%) for identification, and confirmations as per the SANTE guidelines.⁴

Results and discussion

Sample preparation

Pesticide residue analysis in chili powder is challenging due to the matrix complexity. The chili powder contains red colored pigments (alkaloids) that can quickly contaminate the GC liner and column. The use of dSPE cleanup with GCB material significantly removed the amount of pigment, reducing the intensity of color more than 40% in comparison with the raw extract. After the cleanup, the red coloration of the extract became a pale vellow. This cleanup step reduced the maintenance of the liner and column therefore reducing the cost per sample. Still, matrix enhancement (>20%) was observed compared to solvent standards (Figure 2), hence the need for matrixmatched calibration standards in the range of 0.005 to 0.20 mg/kg for accurate quantitation. Also, the final acetonitrile extract carried 0.33 g/mL (3× dilution). So, the lower calibration level of 0.005 mg/kg (actual level) was used to cover the range of final diluted (3x) acetonitrile extract (0.33 g matrix/mL).

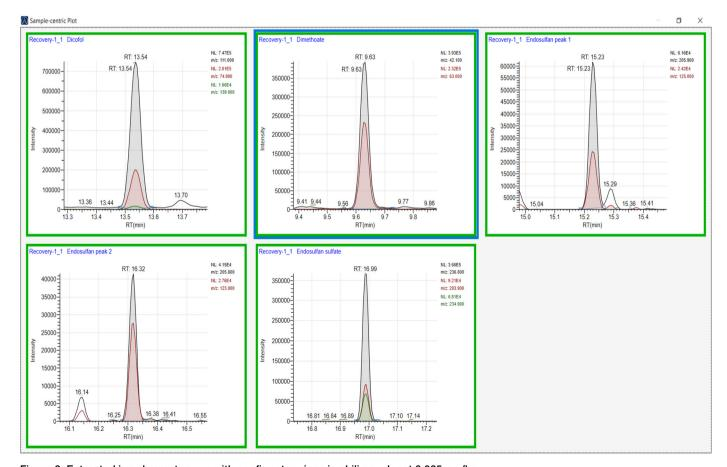


Figure 2. Extracted ion chromatogram with confirmatory ions in chili powder at 0.025 mg/kg $\,$

GC-MS/MS analysis

The optimized instrumental conditions offered Gaussian peak shapes for all target analytes for the spiked concentration, i.e., 0.025 mg/kg (Figure 2). The QuChERS method for extraction was used and the solvent exchange was avoided, i.e. from acetonitrile to any non-polar solvent, hence directly acetonitrile. Because of the high expansion volume, a 2 μ L injection volume was preferred in the PTV injector, which offered symmetrical peak shapes without losing the target analyte's peak quality.

Method performance

Due to the matrix enhancement, the matrix-matched calibration curve was plotted in the range of 0.005 to

0.20 mg/kg (Figure 3). This curve offered linearity with $R^2 > 0.99$ and within 15% residuals by following the linear curve fitting and the 1/x weighting factor. Average recoveries were within 72% to 104% with <6.2% RSD at 0.025 (LOQ) and 0.05 mg/kg (n=6). For the identification, both the ions should overlap (±0.1 min retention time window) on the same retention time as per the SANTE guidelines, which has been demonstrated in Figure 2 and the retention time repeatability in Figure 4. The overall optimized method provided good ion ratios (±30%) presented in Figure 5. Recoveries and precision presented in Table 3 were within the acceptance criteria as per SANTE/11813/2017.⁴ Hence, the described method can be confidently used for the routine analysis of pesticides in chili powder.

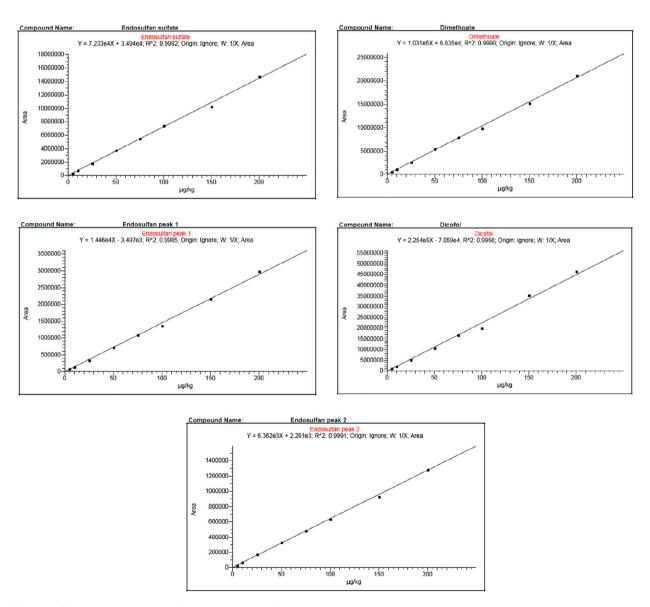


Figure 3. The matrix-matched calibration standards linearity for target analytes

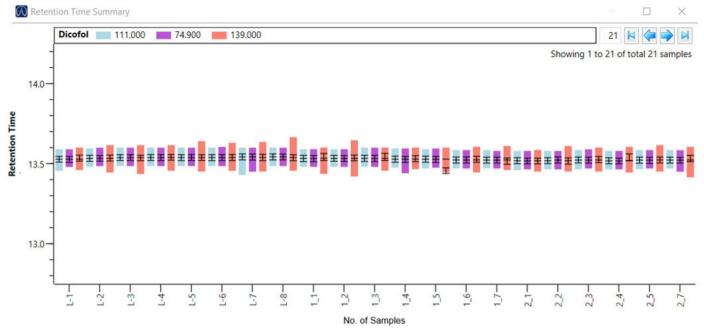


Figure 4. Retention time stability of dicofol in chili powder as assessed across n=25 consecutive matrix injections in a batch

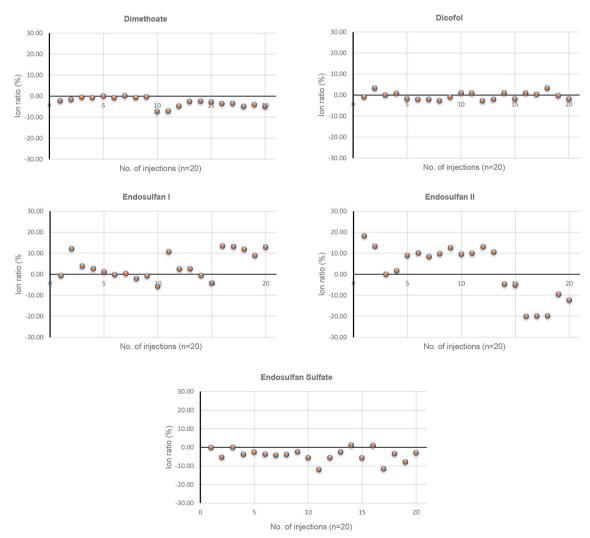


Figure 5. Ion ratio % difference observed range (±30%) for the above target analytes in chili powder spiked at 0.025 mg/kg in replicate injections (n=20)

Table 3. SRM transitions, retention time, R2, LOQ, recovery and precision, and FSSAI MRLs for target analytes

							% Recovery (%RSD)		FSSAI
Name of compound	RT (min)	Q1 (<i>m/z</i>)	Q3 (m/z)	CE (V)	R²	LOQ (mg/kg)	0.025 (mg/kg)	0.05 (mg/kg)	MRLs (mg/kg)
		87	42	10			00	00	
Dimethoate	9.63	93	63	8	0.999	0.025	90 (2.3)	96 (6.2)	5.00
		125	79	8			(2.0)	(0.2)	
	13.54	139	111	12	0.995	0.025	104 (2.4)	07	10.00
Dicofol		111	75	12				97 (3.8)	
		251	139	12				(0.0)	
		241	206	14			00	72	
Endosulfan-alpha	15.24	195	125	22	0.998	0.025	90 (5.5)	(2.7)	
		195	159	8			(0.0)	(=)	
Endosulfan-beta	16.31	159	123	12	0.999	0.025	85 (5.3)	06	0.025*
		195	159	8				86 (6.0)	
		241	206	12				(0.0)	
Endosulfan-sulfate	16.99	272	237	12	0.998	0.025	85 (3.5)	75	
		239	204	12				(3.5)	
		272	235	12			(0.0)	(0.0)	

Q1=precursor ion, Q3=product ion, RT=retention time, CE=collision energy

Conclusion

The developed method successful complies with the MRL requirements of the FSSAI and the method performance criteria of the EU SANTE guidelines for the determination of pesticides in chili powder by GC-MS/MS. The optimized method demonstrated that the LOQ (0.025 mg/kg) is much lower than the MRLs, except for endosulfan, without compromising data quality. Even at the LOQ level, identification and confirmation with retention time, ion ratio, recoveries (70-120%) and precision (<20%) offered by the method are within acceptance criteria of SANTE. This method was fast, allowing a batch of 10 samples to be analyzed within eight hours. This method can be implemented for routine pesticide residue analysis and target analytes sensitivity meeting the FSSAI MRLs requirement (The Food Safety Standard Act, 2006), in the chili powder.

References

- Government of India, Ministry of Agriculture and Farmers Welfare, Department of Agriculture, Cooperation & Farmers Welfare, Directorate of Plant Protection, Quarantine and Storage, List of Pesticides which are Banned, Refused Registration And Restricted in Use as on 19.03.2019. http://ppqs.gov.in/divisions/cib-rc/registered-products
- 2. C. Sivanandha reddy, S. Kulavardhana reddy and Gopireddy Venkata Subba Reddy, Int. J. Eng. Sci Invention (IJESI) ISSN (Online): 2319–6734, (2018), 2319–6726.
- 3. Arnab Goon, Zareen Khan, Dasharath Oulkar, Raviraj Shinde, Suresh Gaikwad, Kaushik Banerjee, *J. Chromatogr.*, *A*, **2018**, *1532*,105–111.
- Guidance document on analytical quality control and method validation procedures for pesticide residues and analysis in food and feed. https://ec.europa.eu/food/sites/food/ files/plant/docs/pesticides_mrl_guidelines_wrkdoc_2017-11813.pdf. (Accessed April 24, 2019)
- 10.1.04. AOAC Official Method 2007.01 Pesticide Residues in Foods by Acetonitrile Extraction and Partitioning with Magnesium Sulfate. 2007 AOAC international.
- Food Safety and Standards Authority of India, The Food Safety and Standards Regulations. https://www.fssai.gov.in/home/fss-legislation/fss-regulations.html. (Accessed April 24, 2019)

Find out more at thermofisher.com



^{*} MRL set at LOQ