Determination of Pesticide Residues in Drinking Water Using Automated Solid-Phase Extraction and Gas Chromatography with Nitrogen Phosphorus Detection

Liu Qian, Zheng Hongguo and Yu Chongtian Thermo Fisher Scientific, People's Republic of China

Key Words

AutoTrace 280, SPE, TRACE 1310 GC, GC-NPD, Organophosphorous Pesticides, Chinese Method GB 5750.9-2006

Introduction

Organophosphorous pesticides have been widely used to prevent and treat plant diseases and pests in agricultural products and as a result, excessive residual pesticide content is frequently reported in drinking and surface water sources. A number of organophosphorous pesticides are acutely toxic to human and animal life, making the detection of residual organophosphorous pesticides a growing concern.

This application note addresses the detection of trace amounts of organophosphorus pesticides in water and optimizes the conditions for solid-phase extraction (SPE) and chromatographic analysis using the Thermo Scientific™ Dionex™ AutoTrace 280 Solid-Phase Extraction instrument and the Thermo Scientific™ TRACE™ 1300 Series gas chromatograph with a nitrogen-phosphorus detector (NPD). Eight different types of organophosphorus pesticide residues were detected in drinking water. The results show that this method achieves rapid, simple extraction with small quantities of solvent. In addition, the method has high sensitivity to the eight types of organophosphorus pesticides, a high recovery rate and good linear range.



Summary of the Method

The Chinese Ministry of the Environment Method GB 5750.9-2006 is used to extract and analyze organophosphorous pesticide residues in environmental waters. This method primarily uses liquid–liquid extraction (LLE) to remove the residues but automated SPE can be used in place of LLE to reduce the quantity of organic solvent used. A nitrogen–phosphorus detector (NPD) was used to selectively detect trace amounts of organophosphates in the water.

Sample Preparation

Methanol (2 mL) and NaCl (5 g) were added to 500 mL of tap water taken from the Beijing laboratory. These samples where then spiked with 50 μ L of 10 μ g/mL mixed standard solution. The sample was purified on a Thermo ScientificTM HyperSepTM Retain PEP SPE cartridge and eluted with ethyl acetate and dichloromethane. The collected liquid was concentrated to 1 mL by nitrogen blowing and analyzed using gas chromatography.



No.	Method	SPE Steps		
1	Add 2 mL of methanol and 5 g of NaCl to 500 mL of the sample water			
2	Spike with 50 μL of 10 μg/mL mixed standard solution			
3	Purify sample on a HyperSep Retain PEP SPE Cartridge (500 mg/6 mL, P/N 60107-206)			
4	Elute sample with ethyl acetate and dichloromethane			
5	Concentrate collected liquid to 1mL by nitrogen blowing and analyzed using gas chromatography	Collect		
Proces	Process six samples using the following procedure:			
1	Condition cartridge with 5.0 mL methanol (MeOH) into solvent waste			
2	Condition cartridge with 5.0 mL ethyl acetate (EtOAc) into solvent waste			
3	Condition cartridge with 5.0 mL dichloromethane (DCM) into solvent waste			
4	Condition cartridge with 5.0 mL ethyl acetate (EtOAc) into solvent waste			
5	Condition cartridge with 5.0 mL MeOH into solvent waste			
6	Load 550.0 mL of sample onto cartridge			
7	Load 550.0 mL of sample onto cartridge			
8	Dry cartridge with gas for 10 min			
9	Collect 4.0 mL fraction into sample tube using EtOAc			
10	Collect 4.0 mL fraction into sample tube using EtOAc			
11	Collect 4.0 mL fraction into sample tube using DCM			

Test Conditions				
Solvent 1	Water (reagent grade)			
Solvent 2	CH ₃ OH (methanol)			
Solvent 3	EtOAc (ethyl acetate)			
Solvent 4	DCM (dichloromethane)			
Condition flow	15.0 mL/min			
Load flow	10.0 mL/min			
Rinse flow	10.0 mL/min			
Elute flow	3.0 mL/min			
Condition air flow	15.0 mL/min			
Rinse air flow	10.0 mL/min			
Elute air flow	3.0 mL/min			

Table 1. Instrument parameters and settings.

landon mandal and	TRACE 1310 GC			
Instrument model and configuration	Instant Connect SSL injector			
Comigaration	AS 1310 autosampler			
chromatography column oven temperature program	TG-1701 MS, 30 m, 0.25 mm, 0.25 μm 80 °C (1 min), 20 °C/min to 250 °C (0 min), 5 °C/min			
Detector type and working parameters	Instant connect NPD: 300 °C Air: 60 mL/min Hydrogen: 2.3 mL/min Make-up gas: Nitrogen Add this info 15 mL/min			
Carrier gas type and flow rate	Nitrogen, constant current mode, flow rate: 1 mL/min			
Injection method and volume	Splitless, for 1 min, Injector temperature of 250 °C; sample volume: 1 mL			

No.	Extraction Procedure
1	Activate extraction cartridge with 5.0mL MeOH into the solvent waste channel
2	Activate extraction cartridge with 5.0 mL EtOAc into the solvent waste channel
3	Activate extraction cartridge with 5.0 mL DCM into solvent waste channel
4	Activate extraction cartridge with 5.0 mL EtOAc into solvent waste channel
5	Activate extraction cartridge with 5.0 mL MeOH into solvent waste channel
6	Activate extraction cartridge with 5.0 mL water into solvent waste channel
7	Load a 550.0 mL (10 mL/min) sample onto the extraction cartridge
8	Rinse the extraction cartridge with 5.0 mL of water into the waste channel
9	Dry the small extraction cartridge with nitrogen (15 mL/min for a period of 10 min)
10	Collect 4.0 mL of eluate from the extraction cartridge with EtOAc in the sample collection tube
11	Collect 4.0 mL of eluate from the extraction cartridge with EtOAc in the sample collection tube
12	Collect 4.0 mL of eluate from the extraction cartridge with EtOAc in the sample collection tube

Gas Chromatography Conditions				
Instrument model and configuration	TRACE 1310 GC Instant Connect SSL injector AS 1310 autosampler			
Chromatography column Oven temperature program	TG-1701 MS, 30 m, 0.25 mm, 0.25 μm 80 °C (1 min), 20 °C/min to 250 °C (0 min), 5 °C/min			
Detector type and working parameters	Instant Connect NPD: 300 °C Air: 60 mL/min Hydrogen: 2.3 mL/min Make-up gas: Nitrogen 15 mL/min			
Carrier gas type and flow rate	Nitrogen, constant current mode, flow rate: 1 mL/min			
Injection method and of volume	Splitless, for 1 min, injector temperature 250 °C; sample volume: 1 mL			

Results and Discussion

Optimization of Conditions for SPE Device

GB5749-2006 "Hygiene Standards for Domestic Drinking Water" requires that residual amounts of eight types of organophosphorus pesticide in drinking water be monitored. The HyperSep Retain PEP SPE cartridge (6 mL barrel with 500 mg sorbent material, part number 60107-206) was used for this method.

The elution efficiencies of solvents including methanol, ethyl acetate, dichloromethane and n-hexane prepared in different proportions on the organophosphorus pesticides retained on the HyperSep Retain PEP SPE cartridge were tested. The test results showed that carrying out elution with ethyl acetate/ethyl acetate/dichloromethane (4 mL/4 mL/2 mL) produced the highest elution efficiency with the smallest quantity of solvent. The boiling points of demeton and phorate were low and solubility in water was quite high and the recovery rate produced by the direct use of SPE was not ideal. However, adding NaCl to the water increased the recovery rate.

Extraction efficiencies were tested by adding 2, 5 and 10 g of NaCl to the water. The recovery rate was highest when adding 5 g NaCl to the water. This is due to the ability of NaCl to reduce the solubility of organic analytes in water, enabling greater retention during the extraction process and resulting in higher recovery rates.

Figure 1 shows the effect of adding different amounts of NaCl on phorate recovery rates. Another reason for low demeton recovery rates is the low boiling point, which makes it degrade easily during testing. Large quantities of demeton are lost easily during the concentration process. Therefore, lowering the the nitrogen blowing temperature to 30 °C during the concentration process and reducing the flow rate to prevent rapid degradation are recommended.

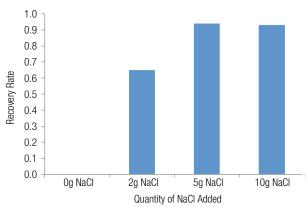


Figure 1. The effect of the quantity of NaCl added on recovery rates.

Standard Chromatogram and Spiked Chromatogram

See Figure 3 for the standard chromatograms of the eight organophosphate analytes. See Figure 2 for a blank sample that went through the entire preparation process and Figure 4 for chromatogram of the domestic drinking water sample spiked with the eight organophosphate species tested.

Linearity, Detection Limits and RSD

Calibration solutions were prepared in the following concentrations: 0.05, 0.1, 0.2, 0.5 and 1.0 μ g/L and analyzed the solutions using the instrument method specified in Table 1 to investigate linearity for each component within a concentration range of 0.05–1.0 μ g/mL. The test results showed a good linear relationship for the eight compounds between 0.05–1.0 μ g/mL. The linear correlation coefficients were all greater than 0.99 (see Table 2).

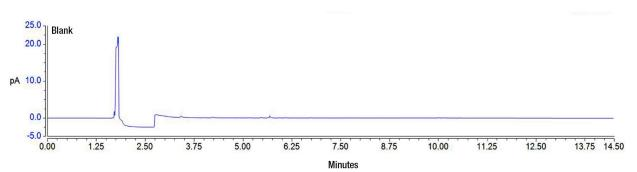


Figure 2. Blank sample chromatogram.

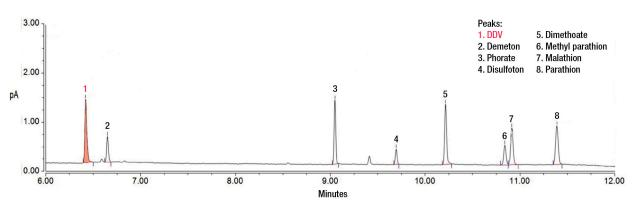


Figure 3. Chromatogram for 0.5 µg/mL mixed standard solution.

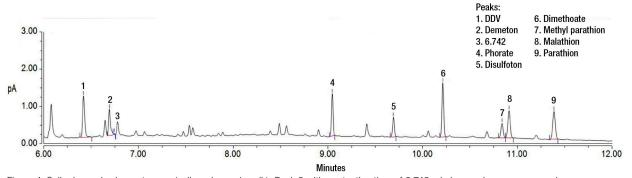


Figure 4. Spiked sample chromatogram (spike volume: $1 \mu g/L$). Peak 3 with a retention time of 6.742 min is an unknown compound.

Table 2. Linearity, detection limits and RSD Data (n=3).

Compound	Linear equation	R²	Spiked sample 0.2 µg/L Recovery rate/%	Spiked sample 0.4 µg/L Recovery rate/%	Spiked sample 1.0 µg/L Recovery rate/%	Detection limit/µg/L	RSD/% (n=6)
Dichlorvos	Y = (0.0415 + 6.0597)*X	0.9904	93.2	88.4	98.6	0.01	4.1
Demeton	Y = (0.0427 + 18.1974)*X	0.9948	83.3	88.5	90.7	0.05	8.7
Phorate	$Y = (0.0124 + 6.7178) \times X$	0.9936	98.7	91.5	95.4	0.05	6.3
Disulfoton	Y = (0.0081 + 14.7635)*X	0.9956	95.3	89.8	95.7	0.01	7.3
Dimethoate	Y = (0.0419 + 5.2029)*X	0.9909	88.5	89.3	99.8	0.01	2.4
Methyl parathion	Y = (0.0258 + 16.4662)*X	0.9933	94.0	90.9	89.3	0.05	3.8
Malathion	Y = (0.0266 + 8.1784) *X	0.9939	90.5	88.8	90.1	0.01	4.4
Parathion	Y = (0.0193 + 7.2690)*X	0.9943	93.2	95.7	96.4	0.01	3.5

A mixed standard solution was added to the water samples (spiked concentrations: 0.2, 0.4 and 1 μ g/L) to investigate the spike recovery rates of the eight pesticides. The test results showed that spike recovery rates for each component are between 83% and 100% and are in compliance with routine analysis detection requirements. At the same time, water samples were spiked with 0.2 μ g/L, 0.4 μ g/L and 2.0 μ g/L and analyzed six times in parallel. The average RSD value was between 2.4% and 8.7%, in compliance with stability requirements. In the experiment, the detection limit for each component was calculated using three times the signal-to-noise ratio. The detection limits for each component were 0.02–0.1 μ g/L (see Table 2).

Summary

The Dionex AutoTrace 280 SPE instrument can extract organic compounds from water samples ranging from 20 mL to 4 L. The Dionex AutoTrace 280 automatically carries out the activation, loading, rinsing and elution of the chromatography column, greatly reducing preprocessing times and reducing the time needed to manually load the samples (a 500 mL sample of water requires 50 min, so 5 hours are saved when processing six samples). In addition, the Dionex AutoTrace 280 offers excellent reproducibility, making it possible to carry out laboratory analysis and detection on largevolume water samples. Excellent sensitivity, repeatability and reproducibility in the analysis of organophosphate pesticides have been achieved using the innovative TRACE 1310 gas chromatograph. The NPD design uses ceramic beads that maintain excellent peak shape also for phosphorous specified. Its Instant Connect Split/Splitless injector and NPD detector are user-installable and, making the TRACE 1310 GC a versatile system that can be easily switched to other applications. By simply changing injector and detector modules, different methods can be run on the same GC platform in a matter of a few minutes.

Conclusion

This application note has established a method for determining eight different of organophosphate pesticides in water in accordance with the Chinese Ministry of the Environment Method GB 5750.9-2006 "Hygiene Standards for Domestic Drinking Water". The sensitivity of the method is low, 0.02–0.1 µg/L for each of the eight organophosphates. Recovery rates for the three spiked levels, 0.2, 0.4 and 1.0 µg/L, were between 83% and 100%. The extraction and analysis methods used here offer many advantages for the detection of organophosphates, including sensitivity, accuracy and a wide linear range, satisfying detection requirements for organophosphates in domestic drinking water.

References

- 1. Zhou Fang, Sun Cheng, Zhong Ming et al. Solid-phase extraction and capillary gas chromatography analysis of residual organophosphorus pesticides in water. *Environ. Pollut. Control.* **2005**, 27(2): 151-153.
- Maciej Tankiewicz, Jolanta Fenik, Marek Biziuk.
 Determination of organophosphorus and organonitrogen pesticides in water samples. *Trends Anal. Chem.* 2010, 29(9):1050-1064
- 3. Yao Chaoying, Yang Lili, Mu Yingfeng. Determination of demeton in water by gas chromatography[J]. *Sichuan Environment*. 2007, 26(4): 48-50.
- 4. Hygiene Standards for Domestic Drinking Water, Chinese Ministry of the Environment Method <GB5750-2006>. People's Republic of China.

www.thermoscientific.com/samplepreparation

©2014 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change Not all products are available in all countries. Please consult your local sales representative for details.

Africa +43 1 333 50 34 0 Australia +61 3 9757 4300 Austria +43 810 282 206 Belgium +32 53 73 42 41 Brazil +55 11 3731 5140

Canada +1 800 530 8447 **China** 800 810 5118 (free call domestic) 400 650 5118 Denmark +45 70 23 62 60 Europe-Other +43 1 333 50 34 0 Finland +358 9 3291 0200 France +33 1 60 92 48 00 Germany +49 6103 408 1014 India +91 22 6742 9494 Italy +39 02 950 591 Japan +81 6 6885 1213 Korea +82 2 3420 8600 Latin America +1 561 688 8700 Middle East +43 1 333 50 34 0 Netherlands +31 76 579 55 55 New Zealand +64 9 980 6700 Norway +46 8 556 468 00 Russia/CIS +43 1 333 50 34 0 Singapore +65 6289 1190 Sweden +46 8 556 468 00 Switzerland +41 61 716 77 00 Taiwan +886 2 8751 6655 UK/Ireland +44 1442 233555 USA +1 800 532 4752

ISO 9001:2008 Certified



A Thermo Fisher Scientific Brand