

## Single Solid Phase Extraction Procedure Yields Base/Neutral and Acid Fractions, for HPLC Analysis of Pesticides in Groundwater

Investigators at the Nevada Division of Agriculture have developed a fast, solvent-conserving, automation-compatible solid phase extraction for concentrating pesticides from 400mL–2-liter samples of groundwater. Base/neutral and acidic pesticides are extracted simultaneously in an ENVI-Carb SPE tube, then are eluted as separate groups. Each fraction is concentrated and analyzed by reversed phase HPLC with diode array detection. Low background in the carbon-based adsorbent allows detection limits of 0.1–5.0µg/liter for 29 monitored compounds.

### Key Words:

- pesticides
- solid phase extraction
- sample preparation
- groundwater

Under the Federal Insecticide, Fungicide and Rodenticide Act (FIFRA), the United States Environmental Protection Agency requires states to develop State Management Plans (SMP) for pesticides considered to endanger groundwater. The Nevada Division of Agriculture laboratory (Reno, Nevada, USA) is to assist in monitoring groundwater for those pesticides used in Nevada.

In use at the Nevada laboratory, official methods of sample extraction proved time-consuming and required large amounts of expensive, hazardous solvent that then had to be disposed of. To better serve their needs, these investigators developed a solid phase extraction (SPE) method that is quick, uses minimal solvent, and can be automated. (The Nevada investigators use a Zymark® AutoTrace automated solid phase extraction workstation.) The investigators follow this method in monitoring water samples for 29 base/neutral and acidic pesticides.

The method calls for processing water samples through ENVI™-Carb SPE tubes. In investigations elsewhere, the carbon-based packing in ENVI-Carb tubes has provided superior recovery and less variability for polar analytes, such as acids and bases, while maintaining comparable results for less polar compounds (1-4). In the Nevada analysts' method, base/neutral pesticides and acidic pesticides adsorb to the packing in the ENVI-Carb tube, then are eluted as separate groups by using selective elution steps. Each extract is concentrated, then analyzed by reversed phase HPLC with diode array detection. Chromatograms for representative base/neutral and acidic pesticides, in Figures A and B, respectively, show the analytes are well resolved, and that low background helps to ensure reliable quantification. In developing the procedure, the analysts determined recovery rates for each

**Figure A. Base/Neutral Pesticides in Water, Using SPE/HPLC**

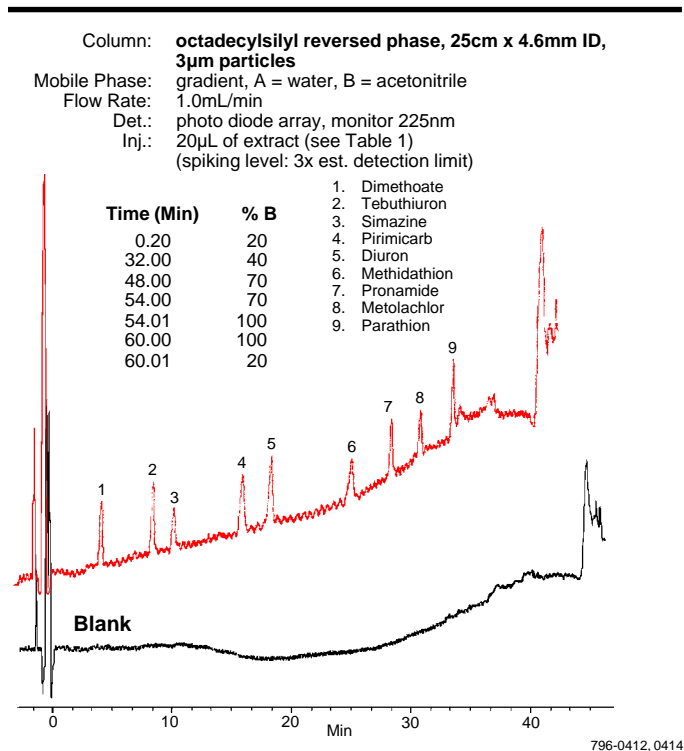


Figure provided by C.L. Ritland, S. Cohen, C.J. Mason, Nevada Div. Agriculture, Reno, Nevada, USA.

analyte at 3 times and at 10 times the estimated detection limit (EDL). Recovery rates at the 3x EDL level are summarized in Table 2. Detection limits for 29 monitored compounds ranged from 0.1 to 5.0µg/liter. Sensitivity can be increased by increasing the sample volume from 400mL to 2 liters. HPLC with diode array detection showed excellent linearity for all analytes, from 1x EDL to 50x EDL.

The results of this work show that automated solid phase extraction is a viable alternative to time-consuming liquid-liquid extractions for monitoring pesticides in water. The graphitized carbon-based adsorbent medium in ENVI-Carb SPE tubes and a selective elution process effectively separate base/neutral analytes from acidic analytes while ensuring high recovery rates.

**Figure B. Acidic Pesticides in Water, Using SPE/HPLC**

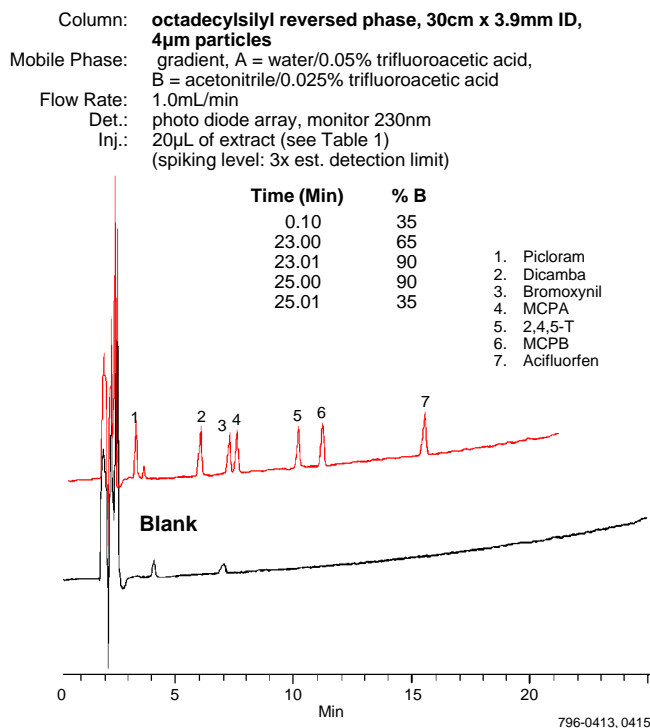


Figure provided by C.L. Ritland, S. Cohen, C.J. Mason, Nevada Div. Agriculture, Reno, Nevada, USA.

**Table 2. Recovery of Acidic and Base/Neutral Pesticides (3x Estimated Detection Limit)**

Analyte	Concentration (µg/L) EDL	Concentration (µg/L) in Sample	Recovery (%)	Mean % RSD
Dimethoate	5.0	15.0	89.2	8.0
Tebuthiuron	2.5	7.5	88.7	8.3
Simazine	0.1	0.3	86.2	7.4
Pirimicarb	1.3	3.7	83.4	9.4
Diuron	1.3	3.7	94.9	6.2
Methidathion	1.3	3.7	86.0	18.4
Pronamide	0.8	2.3	85.3	11.9
Metolachlor	0.8	2.3	88.4	18.8
Parathion	2.5	7.5	71.8	13.3
Bromacil	1.3	3.7	98.5	14.1
Cyanazine	0.2	0.6	93.9	18.9
Atrazine	0.2	0.6	95.1	17.4
Prometon	0.3	0.7	98.8	18.1
Chlorpropham	0.4	1.2	50.7	36.9
Alachlor	1.7	5.2	84.5	17.9
Dacthal	1.0	1.9	93.1	29.6
Picloram*	0.5	1.5	99.8	3.6
Dicamba*	1.3	3.7	128.8	12.0
Bromoxynil*	1.0	3.0	89.0	4.6
MCPA*	1.3	3.7	100.3	6.2
2,4,5-T*	1.5	4.5	102.0	6.2
MCPB*	1.2	3.6	93.8	5.4
Acifluorfen*	1.5	4.5	97.7	6.1
Bentazon	1.5	4.5	85.0	2.9
2,4-D	1.5	4.5	98.0	4.0
MCPP	1.3	3.7	92.3	3.2
2,4-DB	1.5	4.5	91.6	4.2
2,4,5-TP	1.5	4.5	95.7	4.0
Dinoseb	2.0	6.0	47.1	31.9

\*n = 8, all others n = 9.

**Table 1. Extracting Pesticides from Water**

- Condition ENVI-Carb SPE tube (3mL, 250mg packing) with 5mL methylene chloride:methanol, 4:1, followed by 2mL methanol, followed by 10mL 2% acetic acid.
- Pass 400mL water sample through tube.
- Rinse tube with 10mL water.
- Dry tube with clean nitrogen for 5 minutes.
- Recovery of Base/Neutral Pesticides**  
Pass 1mL methanol, then 2 x 2mL methylene chloride:methanol, 4:1 through tube.  
**Collect all 3 eluates (5mL total) in same recovery vessel.**
- Recovery of Acidic Pesticides**  
Pass 2 x 2mL methylene chloride:methanol, 3:2 (made basic with 0.016mol/liter KOH) through same tube.  
**Collect both eluates (4mL total) in same recovery vessel.**  
Add 350µL 2% trifluoroacetic acid in water (acids eluate only).
- Dry recovered eluates, then reconstitute each with 1mL methanol. Filter. Analyze by RP-HPLC.

**Acknowledgment**

Information and figures were provided by Christopher Ritland, Sharryn Cohen, and Christopher Mason, Nevada Division of Agriculture, 350 Capitol Hill Avenue, Reno, Nevada USA 89502.

**Note:** For information about extracting organochlorine, organo-phosphorus, nitrogen-containing, and carbamate pesticides from fruits and vegetables, request Bulletin 900.

**Ordering Information:**

Description	Cat. No.
<b>ENVI-Carb Solid Phase Extraction Tubes</b>	
3mL, 250mg packing, pk. of 54	<b>57088</b>
6mL, 250mg packing, pk. of 30	<b>57092</b>
6mL, 500mg packing, pk. of 30	<b>57094</b>

**AutoTrace SPE Tube Adaptors**

Prevent damage to packing bed in SPE tube	
3mL, pk. of 6	<b>57123</b>
6mL, pk. of 6	<b>57126</b>

**References**

- The Reporter*, Vol 15, No. 2, Supelco (1996).
- Application Note 27, Supelco (1994).
- DiCorcia, A. and M. Marchetti, *Anal. Chem.*, **63**: 580-585 (1991).
- DiCorcia, A., R. Samperi, A. Marcomini, and S. Stelluto, *Anal. Chem.*, **65**: 907-912 (1993).

References 3 and 4 not available from Supelco.

**Trademarks**

ENVI — Sigma-Aldrich Co.  
 Zymark — Zymark

For more information, or current prices, contact your nearest Supelco subsidiary listed below. To obtain further contact information, visit our website (www.sigma-aldrich.com), see the Supelco catalog, or contact Supelco, Bellefonte, PA 16823-0048 USA.

ARGENTINA · Sigma-Aldrich de Argentina, S.A. · Buenos Aires 1119 AUSTRALIA · Sigma-Aldrich Pty. Ltd. · Castle Hill NSW 2154 AUSTRIA · Sigma-Aldrich Handels GmbH · A-1110 Wien  
 BELGIUM · Sigma-Aldrich N.V./S.A. · B-2880 Bornem BRAZIL · Sigma-Aldrich Quimica Brasil Ltda. · 01239-010 São Paulo, SP CANADA · Sigma-Aldrich Canada, Ltd. · 2149 Winston Park Dr., Oakville, ON L6H 6J8  
 CZECH REPUBLIC · Sigma-Aldrich s.r.o. · 186 00 Praha 8 DENMARK · Sigma-Aldrich Denmark A/S · DK-2665 Vallensbaek Strand FINLAND · Sigma-Aldrich Finland/YA-Kemia Oy · FIN-00700 Helsinki  
 FRANCE · Sigma-Aldrich Chimie · 38297 Saint-Quentin-Fallavier Cedex GERMANY · Sigma-Aldrich Chemie GmbH · D-82041 Deisenhofen GREECE · Sigma-Aldrich (o.m.) Ltd. · Ilioupoli 16346, Athens  
 HUNGARY · Sigma-Aldrich Kft. · H-1067 Budapest INDIA · Sigma-Aldrich Co. · Bangalore 560 048 IRELAND · Sigma-Aldrich Ireland Ltd. · Dublin 24 ISRAEL · Sigma Israel Chemicals Ltd. · Rehovot 76100  
 ITALY · Sigma-Aldrich s.r.l. · 20151 Milano JAPAN · Sigma-Aldrich Japan K.K. · Chuo-ku, Tokyo 103 KOREA · Sigma-Aldrich Korea · Seoul MALAYSIA · Sigma-Aldrich (M) Sdn. Bhd. · Selangor  
 MEXICO · Sigma-Aldrich Química S.A. de C.V. · 50200 Toluca NETHERLANDS · Sigma-Aldrich Chemie BV · 3330 AA Zwijndrecht NORWAY · Sigma-Aldrich Norway · Torshov · N-0401 Oslo  
 POLAND · Sigma-Aldrich Sp. z o.o. · 61-663 Poznań PORTUGAL · Sigma-Aldrich Quimica, S.A. · Sintra 2710 RUSSIA · Sigma-Aldrich Russia · Moscow 103062 SINGAPORE · Sigma-Aldrich Pte. Ltd.  
 SOUTH AFRICA · Sigma-Aldrich (pty) Ltd. · Jet Park 1459 SPAIN · Sigma-Aldrich Quimica, S.A. · 28100 Alcobendas, Madrid SWEDEN · Sigma-Aldrich Sweden AB · 135 70 Stockholm  
 SWITZERLAND · Supelco · CH-9471 Buchs UNITED KINGDOM · Sigma-Aldrich Company Ltd. · Poole, Dorset BH12 4QH  
 UNITED STATES · Supelco · Supelco Park · Bellefonte, PA 16823-0048 · Phone 800-247-6628 or 814-359-3441 · Fax 800-447-3044 or 814-359-3044 · email:supelco@sial.com