

Monitor Pesticides in Water Samples Using a Deactivated Reversed Phase HPLC Column

Because HPLC offers significant advantages compared to GC for many nonvolatile, thermally labile, and polar materials, this technique is becoming a preferred analytical method for monitoring many pesticides and their metabolites. A SUPELCOSM LC-ABZ column was used to separate a mixture of organochlorine and organophosphorous insecticides, triazines, phenylurea derivatives, carbamates, and other products in water samples. The SUPELCOSM LC-ABZ column, which has a unique deactivated stationary phase, provided excellent resolution and peak symmetry. Using this technique, samples of lake water, taken at varying depths, were found to contain traces of triazine herbicides — simazine, atrazine, and terbutylazine.

Key Words:

Herbicides • HPLC • Pesticides

United States Environmental Protection Agency (US EPA) methods for monitoring pesticides in water specify predominantly gas chromatography (GC) methods. Currently, there are only a few US EPA methods for high performance liquid chromatography (HPLC) analyses of pesticides, including Method 632 for urea and carbamate compounds, Method 629 for cyanazine, and the

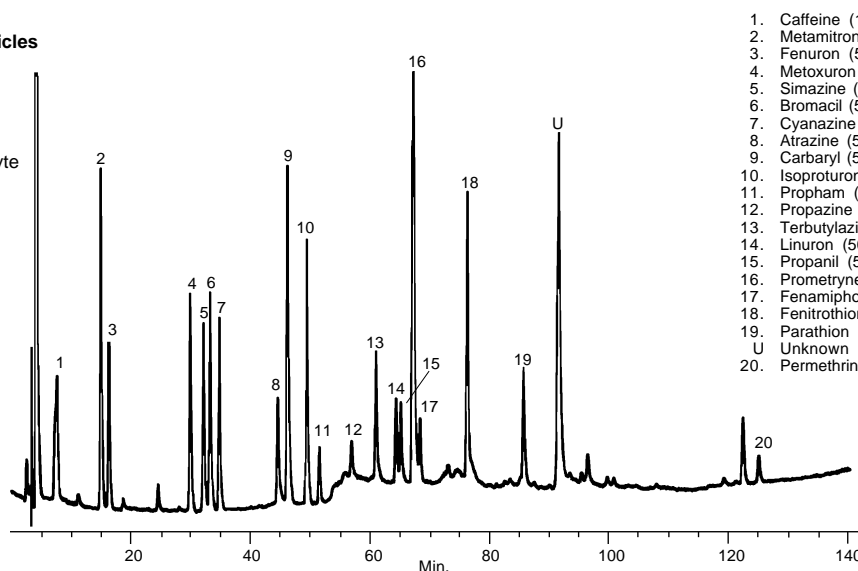
proposed Method 8318 for carbonyl. However, HPLC could be a useful method for a much broader range of pesticides, as demonstrated by investigators in Switzerland, using a SUPELCOSM LC-ABZ column.

With the trend toward biodegradable, more thermally labile, pesticides, HPLC is becoming a preferred analytical method for monitoring many insecticides and their metabolites, as well as herbicides, and plant growth regulators. HPLC offers significant advantages compared to GC for many nonvolatile, thermally labile, and polar materials. HPLC analysis minimizes the adsorptive loss encountered in GC analysis due to inlet and reactive column effects. Large volume injections of aqueous samples are possible using HPLC (e.g., sample screening, analysis of soluble but non-extractable compounds using on-column trace enrichment techniques). HPLC also allows the use of additional sensitive and selective detectors (UV/VIS variable wavelength, diode array, fluorescence, electrochemical, and others). The non-destructive nature of LC analysis allows the collection of the separated pesticides, and the use of post-column reactions to enhance sensitivity with LC detectors.

The SUPELCOSM LC-ABZ column is ideal for HPLC quantitation of pesticides and herbicides in water samples. This column features a unique surface deactivation process that provides well-resolved, symmetrical peaks, suitable for quantitative monitoring of these compounds.

Figure A. Pesticides and Herbicides Resolved on a SUPELCOSM LC-ABZ Column

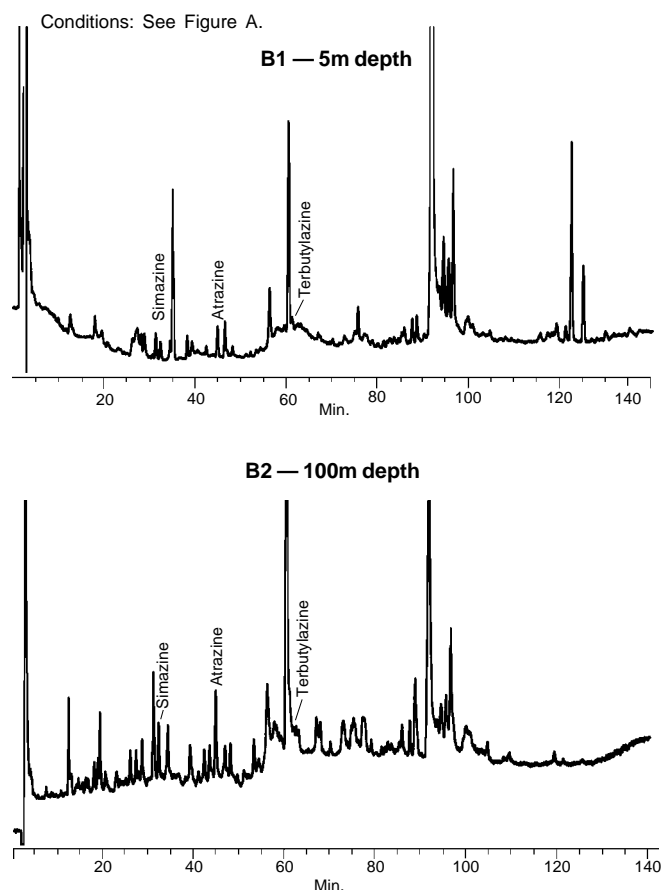
Column: SUPELCOSM LC-ABZ,
25cm x 4.6mm ID, 5µm particles
Mobile Phase: 10% to 90% acetonitrile in
water at 0.5%/min
Col. Temp.: 40°C
Flow Rate: 1mL/min
Det.: 225nm UV
Inj.: 50µL water, spiked with analyte
quantities shown on figure



794-0033

Investigators in Geneva, Switzerland have characterized the ability of the LC-ABZ column to separate mixtures of organochlorine and organophosphorous insecticides, triazines, phenylurea derivatives, carbamates, and related compounds in water samples. A mixture of twenty such compounds in water was separated with sharp, symmetrical peaks (Figure A). Of particular interest is the excellent recovery of carbaryl (peak 9), which is difficult to detect using GC because it thermally decomposes. The 70% recovery of metamitron (peak 2) using the HPLC method was far better than under GC conditions, where it thermally decomposes, producing two small, inconsistent peaks. Samples of water from Lake Geneva, taken at varying depths, were found to contain traces of triazine herbicides — simazine, atrazine, and terbutylazine — as well as unidentified impurities (Figure B).

Figure B. Herbicides Detected in Water Samples from Lake Geneva



Figures reproduced courtesy of Dr. Claude Corvi, Laboratoire Cantonal De Chimie, Geneva, Switzerland. 794-0034.0035

The choice of UV wavelength (225nm) in this separation was the best compromise for the twenty compounds. For example, carbaryl is not detected at 245nm, but is well detected at 218 or 225nm, and cyanazine and metoxuron maxima occur at approximately 220nm and 240nm, respectively.

Although EPA methods that specify HPLC do not exist for all pesticides and herbicides used throughout the US, this technique could be a useful, versatile method for various classes of compounds. We recommend the SUPELCOSIL LC-ABZ column for pesticides analysis because of its surface deactivation and superior performance over a wide range of analyte functionalities.

Ordering Information:

SUPELCOSIL LC-ABZ HPLC Columns

5µm particles, 100Å pores	
5cm x 4.6mm ID	59141
15cm x 4.6mm ID	59140-U
25cm x 4.6mm ID	59142

Guard Column

2cm x 4.6mm ID	59545-U
----------------	---------

Guard Column Kit

2cm x 4.6mm ID	59544-U
----------------	---------

Iso-Disc™ N-253 Syringe-Tip Filters, pk. of 50

25mm diameter, nylon, 0.2µm pores	59321
-----------------------------------	-------

Triazine Pesticides Standard

100µg/mL each component in methanol	
Ametryn	Propazine
Atrazine	Simazine
Prometryn	Terbutryn
Prometon	
	1mL
	48392

Neat Triazine Pesticides Kit

100mg each pesticide, individually packaged	
Ametryn	Propazine
Atrazine	Simazine
Prometryn	Terbutryn
Prometon	
	1mL
	49092

Iso-Disc and SUPELCOSIL are trademarks of Supelco, Inc.

Contact our Technical Service Department (phone 800-359-3041 or 814-359-3041, FAX 814-359-5468) for expert answers to your questions.

Acknowledgment

We are grateful to Dr. Claude Corvi and Dr. Sophal Khim-Heang of Laboratoire Cantonal De Chimie, Geneva, Switzerland for their work on the use of HPLC for environmental analyses.

Note 7

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 Quimica 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

H

Supelco is a member of the Sigma-Aldrich family. Supelco products are sold through Sigma-Aldrich, Inc. Sigma-Aldrich warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product for a particular use. Additional terms and conditions may apply. Please see the reverse side of the invoice or packing slip.

AQI