

Application

Note
99

SPME/HPLC: A Rapid and Sensitive Analysis of Polynuclear Aromatic Hydrocarbons in Water

An SPME/HPLC interface enables HPLC analysts to take advantage of the time and solvent savings offered by SPME. Peaks for 16 polynuclear aromatic hydrocarbons were sharp and symmetric, confirming efficient transfer from the SPME fiber to the HPLC column. Relative to peaks for directly injected analytes, peaks for most of the extracted analytes were significantly larger, indicating a concentrating effect that can increase the sensitivity of the analysis for trace-level analytes.

Key Words:

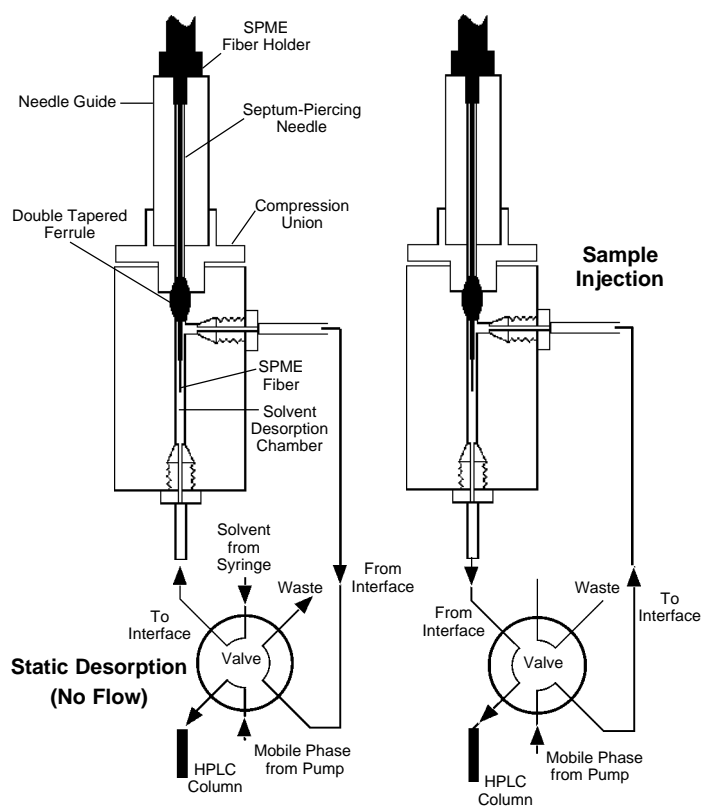
- polynuclear aromatic hydrocarbons • sample extraction
- solid phase microextraction • SPME

Solid phase microextraction – SPME* – has rapidly been established among the practical alternatives for sample preparation for gas chromatography. Polynuclear aromatic hydrocarbons and other weakly volatile or thermally labile compounds are more effectively monitored by HPLC but, initially, there was no simple way to introduce analytes extracted by SPME onto an HPLC column.

An SPME/HPLC interface developed at Supelco enables HPLC analysts to take advantage of the time and solvent savings offered by SPME. Easily installed and removed, the interface consists of a six-port injection valve and a desorption chamber (Figure A) that replaces the injection valve in the HPLC system. The SPME fiber is introduced into the desorption chamber under ambient pressure when the injection valve is in the “load” position. After the fiber is inserted through the ferrule, the unit is made leak-tight by closing a sealing clamp and compressing the ferrule against the SPME needle. All surfaces which contact the SPME fiber or the mobile phase are stainless steel or VESPEL®. Within the interface, mobile phase contacts the SPME fiber, removes the adsorbed analytes, and delivers them to the column for separation. Analytes can be removed in a moving stream of mobile phase (dynamic desorption) or, when analytes are more strongly adsorbed to the fiber, the fiber can be soaked in mobile phase before the injection is made (static desorption).

Analysts at the University of Waterloo (Ontario, Canada) analyzed polynuclear aromatic hydrocarbons, using SPME/HPLC (1). Chemists in Supelco laboratories also have evaluated the practicality of the technique for monitoring PAHs in water, using SPME fibers with 7 μ m, 30 μ m, or 100 μ m polydimethylsiloxane coating to extract the 16 PAHs listed in US Environmental Protection Agency methods 610 and 8310. All but three of the analytes were present at a concentration of 100ppb or 200ppb. Equilibration was fastest

Figure A. SPME/HPLC Interface: Desorption Chamber



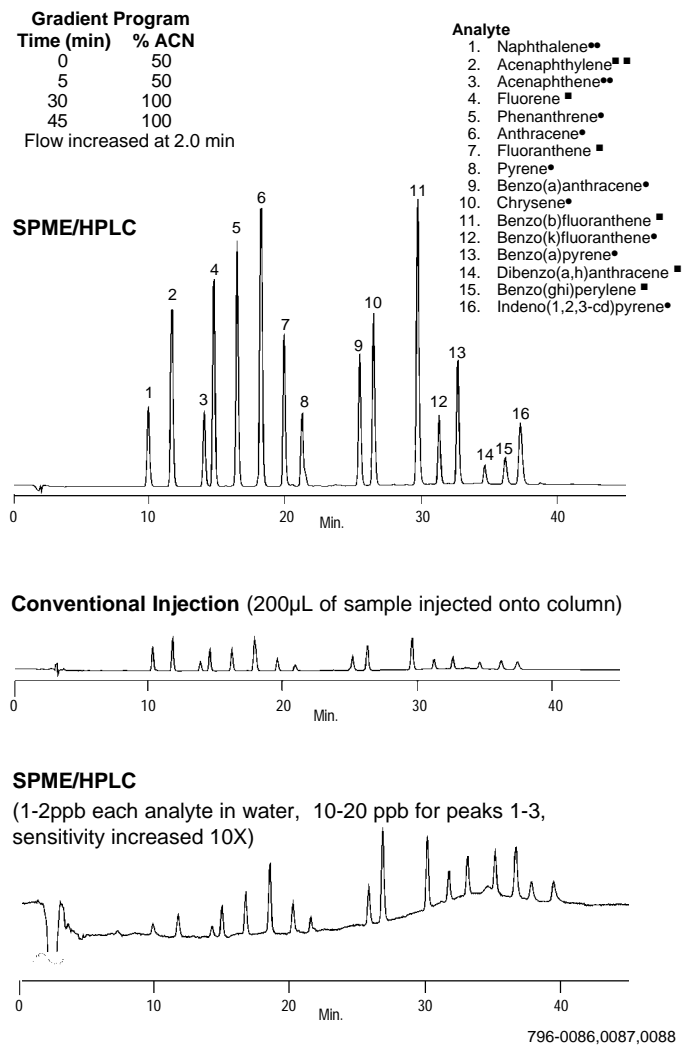
96-0068, 796-0032, 0033

with the 7 μ m coating, but the 100 μ m coating provided the highest extraction levels (i.e., greatest sensitivity). After a 30-minute immersion in the sample, the fiber was introduced into the SPME/HPLC interface, and with the injection valve in the “load” position, the desorption reservoir was filled with 200 μ L of 40:60 acetonitrile:water. The fiber was soaked in the solvent for 2 minutes, then the analysts switched the valve to the “inject” position and began integration. During the first 2 minutes, the analytes were delivered to the column at a low flow rate (0.2mL/min), to minimize band broadening on the column. After 2 minutes, the valve was returned to the “load” position and the desorption chamber was flushed with 500 μ L of acetonitrile to eliminate the possibility of analyte carry-over to the next extraction. The desorption chamber was flushed and refilled with 500 μ L of 40:60 acetonitrile:water, the fiber was removed, and the flow

Figure B. PAHs Effectively Extracted from Water

Sample: 5µL PAH mix (Cat. No. 48743) in 5mL water (SPME) or acetonitrile:water, 40:60 (conventional injection)
 •100ng/mL ■200ng/mL ••1000ng/mL ■■2000ng/mL

SPME Fiber: 100µm polydimethylsiloxane
 Cat. No.: 57301
 Extraction: immersion, 30 min (rapid stirring)
 Desorption: static, 200µL acetonitrile:water, 40:60, 2 min
 Column: **SUPELCOSIL LC-PAH, 15cm x 4.6mm ID, 5µm particles**
 Cat. No.: 58318
 Mobile Phase: acetonitrile:water gradient (see program)
 Flow Rate: 0-2 min: 0.2mL/min
 2-45 min: 1.0mL/min
 Det.: UV, 254nm



rate was increased to 1mL/min. A 1-minute air drying period was sufficient to prepare the fiber for the next extraction.

The analytes were best separated on a 15cm SUPELCOSIL™ LC-PAH column. Sharp, symmetric peaks indicate efficient transfer of the analytes from the SPME fiber to the HPLC column (Figure B). Relative to peaks for directly injected analytes, peaks for the extracted analytes were significantly larger, indicating a concentrating effect that can increase the sensitivity of the analysis for trace-level analytes. Subsequent extraction of the same analytes at 1-2ppb each clearly demonstrates the effectiveness and sensitivity of the extraction technique (Figure B).

If your HPLC analyses call for time-consuming and expensive solvent-based extractions, SPME may very well be a better approach. In addition to analyses of PAHs, investigators have shown that SPME can be effective for monitoring drugs and drug metabolites in biological fluids (for an example – amphetamines in urine – request Application Note 83). The SPME/HPLC interface will accelerate the development of many new environmental, pharmacological, and food and beverage applications.

Ordering Information:

Description	Cat. No.
SPME/HPLC Interface with Valco® Valve	57350-U
SPME/HPLC Interface with Rheodyne® Valve	57353
SPME Holder	
Initially you must order both holder and fiber assembly. Holder is reusable indefinitely.	
For HPLC (also for Varian 8100/8200 AutoSampler)	57331
SPME Fiber Assembly	
100µm polydimethylsiloxane, pk. of 3	
For HPLC (also for Varian 8100/8200 AutoSampler)	57301
SUPELCOSIL LC-PAH Column	
15cm x 4.6mm ID, 5µm particles	58318
EPA 610 Polynuclear Aromatic Hydrocarbons Mix	
16 analytes at concentrations indicated on Figure B in methanol:methylene chloride, 50:50.	
1mL	48743
To see our complete selection of SPME fibers, refer to the Supelco catalog, or call our Technical Service chemists.	

Reference

1. Chen, J. and J. Pawliszyn, *Anal. Chem.* **67**: 2530-2533 (1995).
 Reference not available from Supelco. For information, contact the authors at The Guelph-Waterloo Centre for Graduate Work in Chemistry, University of Waterloo, Waterloo, Ontario, Canada N2L 3G1.

Trademarks

Rheodyne — Rheodyne, Inc.
 SUPELCOSIL — Sigma-Aldrich Co.
 Valco — Valco Instruments Co., Inc.
 VESPEL — E.I. du Pont de Nemours & Co., Inc.
 *Technology licensed exclusively to Supelco. US patent no. 5,691,206; European patent #0523092.

Contact our Technical Service Department
 (phone 800-359-3041 or 814-359-3041, FAX 800-359-3044 or 814-359-5468) for expert answers to your questions.

Note 99

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 Química 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 Química, 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 Química, 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

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.

ASF