

Application

Note
65

ENVI-18 SPE Tube Ensures Low Background for Monitoring Organic Compounds in Drinking Water by EPA Method 525

ENVI-18 solid phase extraction tubes help ensure rapid, economical, reliable sample preparation and provide excellent analyte recovery and clean chromatograms in the screening or monitoring of drinking water samples.

Key Words:

- ENVI-18 SPE tubes • organic contaminants
- drinking water • EPA Method 525

Based on an ongoing review of sample preparation techniques, the US Environmental Protection Agency (EPA) has designated solid phase extraction cartridges for the cleanup and concentration of organic contaminants from drinking water samples (1,2).

ENVI™-18 SPE tubes are compatible with the large sampled volume (1 liter) and low pH (pH 2) specified in EPA Method 525. They can consistently accept large samples and low pH without loss of phase from the silica support. EPA precautions warn that, if phase is stripped from the extraction tube packing at pH 2, the chromatographic analysis can be complicated by high background (Figure A). High background can obscure compounds of

interest. In contrast, the specially prepared packing in ENVI-18 tubes meets the background specifications set by the EPA for extraction analysis by GC-MS (Figure B). ENVI-18 tubes also ensure consistent, high compound recovery .

Supelco has also developed procedures for extracting several classes of environmental contaminants from water samples — phthalate esters, organochlorine pesticides and triazine herbicides, and polynuclear aromatic hydrocarbons — using drinking water samples spiked with contaminants (Tables 1 and 2).

To significantly reduce preparation time, sample volumes were kept at 250mL. These smaller samples can also reduce the costs and inconveniences of sampling, transporting, and storing larger

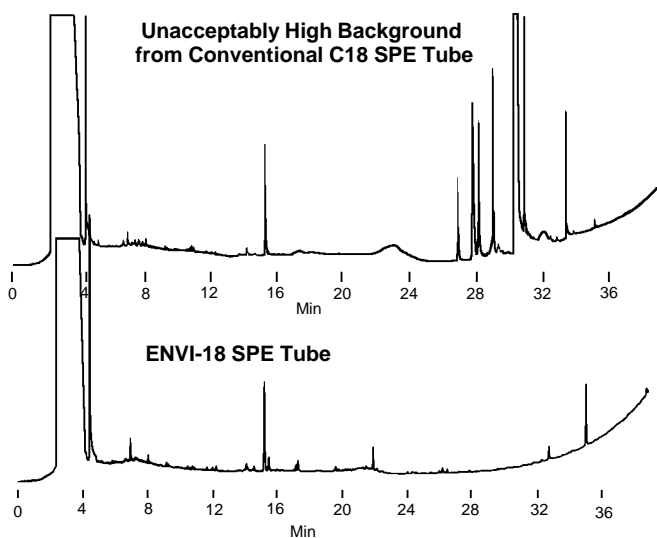
Figure B. Total Ion Chromatograms from Extracted Blank SPE Tubes

Conditions for Figure B2

Sample: 1 liter water
Extraction Tube: Supelclean ENVI-18, 6mL, 0.5g packing
Cat. No.: 57064
Column: PTE-5 fused silica, 30m x 0.25mm ID, 0.25µm film
Cat. No.: 24135-U
Col. Temp.: 40°C (1 min) rapidly to 160°C (3 min) to 320°C at 6°C/min
Carrier: helium, 33cm/sec
Inj.: see Figure A

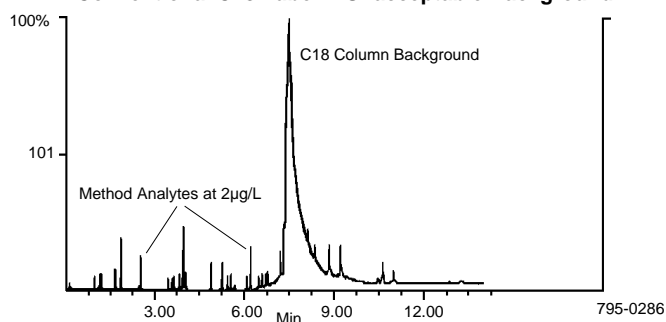
Figure A. GC-FID Chromatograms: ENVI-18 Phase Is Not Stripped from the Support at pH 2

Sample: 1 liter water
Extraction Tube: Supelclean ENVI-18, 6mL, 0.5g packing
Cat. No.: 57064
Column: PTE-5 QTM fused silica, 15m x 0.53mm ID, 0.50µm film
Cat. No.: 25355
Inj.: 1µL

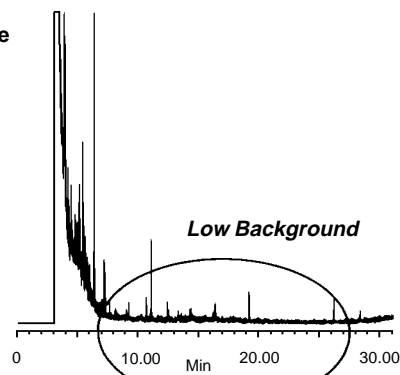


795-0284, 0285

B1 — Conventional C18 Tube — Unacceptable Background*



B2 — ENVI-18 Tube



*Figure B1 adapted from Reference 1.

795-0287

samples. Optimized methods for these smaller sample volumes ensure that elution volumes are reduced – yet compounds of interest can be easily analyzed and quantified without excessive solvent consumption or prolonged eluate concentration times.

Table 1. Procedures for Extracting Organic Compounds from Drinking Water, Using ENVI-18 SPE Tubes (500mg packing bed)

Phthalates	
Conditioning:	2 x 6mL methylene chloride:methanol (1:1) 6mL methanol 6mL deionized water
Sample Addition:	250mL water, unadjusted pH passed slowly through tube (10mL/min)
Drying:	5-10 min
Elution:	2 x 1.5mL methylene chloride:methanol (1:1)* concentrate extract to 1mL
Total Time:	30-40 min
Pesticides/Herbicides	
Conditioning:	2 x 6mL hexane:ethyl ether (1:1) 6mL methanol 6mL deionized water
Sample Addition:	250mL water, unadjusted pH passed slowly through tube (10mL/min)
Drying:	5-10 min (or use sulfate tube)
Elution:	2 x 1.5mL hexane:ethyl ether (1:1)* concentrate extract to 2mL
Total Time:	30-40 min
Polynuclear Aromatic Hydrocarbons	
Conditioning:	2 x 6mL toluene:methanol (10:1) 6mL methanol 6mL deionized water
Sample Addition:	250mL water, unadjusted pH passed slowly through tube (10mL/min)
Drying:	5-10 min
Elution:	2 x 1mL toluene:methanol (10:1)
Total Time:	30-40 min

* Extract can be passed through a bed of sodium sulfate, to remove traces of water, before concentration.

With our extraction procedures, recovery exceeded 90% for almost every compound spiked into the water samples (Table 2). We noted significant recovery variation in only one case. In analyses of PAHs, recovery of acenaphthylene appeared to depend on the drying step before elution. If the packing bed was slightly moist, recovery was 90% or greater. If the bed was allowed to dry completely, recovery was reduced.

Excellent analyte recovery and very clean chromatograms can be obtained by using ENVI-18 SPE tubes. If you are screening drinking water samples according to EPA Method 525 – or monitoring phthalates, organochlorine pesticides, triazine herbicides, or PAHs in drinking water – Supelco's ENVI-18 SPE tubes will help ensure rapid, economical, reliable sample preparation.

Reference 1 available from: NTIS, 5285 Port Royal, Springfield, VA USA 22161 (order number PB 89-220-461).

Fused silica columns manufactured under HP US Pat. No. 4,293,415.

ENVI and PTE are trademarks of Sigma-Aldrich Co.

Table 2. Recovery of Organic Compounds

Compound	Recovery (mean % ± CV)**
Phthalates (20µg/L)	
Dimethyl phthalate	93 ± 5.9
Diethyl phthalate	97 ± 4.6
Di-n-butyl phthalate	97 ± 6.8
Benzyl butyl phthalate	100 ± 5.4
Bis(2-ethylhexyl)adipate	99 ± 11
Bis(2-ethylhexyl)phthalate	111 ± 8.3
Pesticides (8µg/L)	
Hexachlorobenzene	87 ± 11
Lindane	99 ± 13
Heptachlor	96 ± 12
Aldrin	94 ± 13
Heptachlor epoxide	98 ± 13
Endrin	93 ± 11
Methoxychlor	110 ± 13
Simazine	89 ± 5.1
Atrazine	100 ± 5.1
Polynuclear Aromatic Hydrocarbons (20µg/L)	
Acenaphthylene	66 ± 28
Fluorene	95 ± 7.5
Phenanthrene	95 ± 8.0
Anthracene	95 ± 11
Pyrene	97 ± 7.1
Benzo(a)anthracene	94 ± 8.3
Chrysene	101 ± 8.8
Benzo(a)fluoranthene	103 ± 10
Benzo(k)fluoranthene	103 ± 9.4
Benzo(a)pyrene	94 ± 11
Indeno(1,2,3-cd)pyrene	102 ± 10
Dibenzo(a,h)anthracene	104 ± 9.7
Benzo(ghi)perylene	101 ± 11

**Mean for 5 extractions.

Ordering Information:

Description	Cat. No.
ENVI-18 Solid Phase Extraction Tubes	
100mg (1mL), box of 100	57062
500mg (3mL), box of 50	57063
500mg (6mL), box of 30	57064
1000mg (6mL), box of 30	57065-U
Capillary GC Columns	
PTE™-5 column	
30m x 0.25mm ID x 0.25µm film	24135-U
PTE-5 QTM column	
15m x 0.53mm ID x 0.50µm film	25355
Chemical Standards	
Phthalate Esters Mix 525***	48223
Polynuclear Aromatic Hydrocarbons Mix 525***	48953-U

***Components are listed in Table 2.

References

- Method 525. Determination of Organic Compounds in Drinking Water by Liquid-Solid Extraction and Capillary Column Gas Chromatography/Mass Spectrometry (Revision 2.1). Environmental Monitoring Systems Laboratory, Office of Research and Development, U.S. Environmental Protection Agency, Cincinnati, OH USA 45268.
- National Primary Drinking Water Regulations; Analytical Techniques 40 CFR Parts 141 and 143 (Final Rule), Federal Register 53 (No. 33), 5142-5147 (Feb. 19, 1988).

Note 65

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.

ARA