

Application

Note
140

Analysis of Fat Soluble Vitamins from Tablets, Using SPME/HPLC

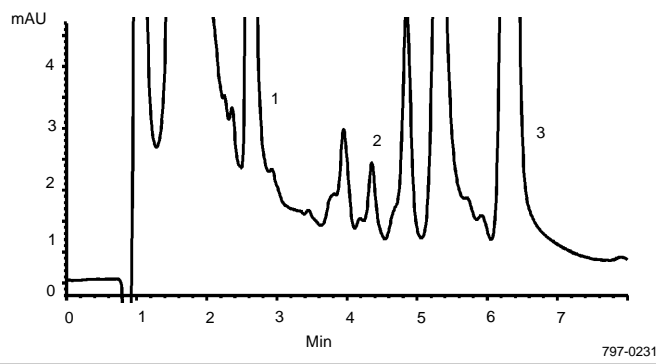
With the recent advent of the SPME/HPLC interface, solid phase microextraction (SPME) presents a novel and simple approach to concentrating and extracting vitamins A, D₃, and E from high potency tablets, for HPLC analysis.*

The best extraction for all three vitamins from a tablet was obtained with polydimethylsiloxane/divinylbenzene (PDMS/DVB) fiber (Figure A). We found that small adjustments to the solvent concentration shifted the equilibrium to such an extent that it considerably affected the fibers' extraction capabilities. Therefore, when determining your own sample extraction procedure, be sure to optimize it carefully.

Figure A. Fat Soluble Vitamins by SPME/HPLC

Sample: 2.5mL supernatant obtained from dissolving vitamin tablet in 1% acetic acid / 33% isopropanol
SPME Fiber: PDMS/DVB, 65µm
Cat. No.: 57311
Extraction: immersion (40 min), with stirring
Desorption: static, 2 min in mobile phase; dynamic, valve open during run
Column: SUPELCOSIL™ LC-8, 25cm x 4.6mm ID, 5µm particles
Cat. No.: 5-8297
Mobile Phase: acetonitrile:methanol:water (63:33:4)
Flow Rate: 2mL/min
Temp.: 35°C

1. Vitamin A
2. Vitamin D₃
3. Vitamin E



The best procedure for sampling vitamins A, D₃, and E from a generic brand high potency tablet is as follows: We sampled tablets which contained these vitamins at levels of 5000IU, 400IU, and 30IU, respectively. To prepare the vitamin tablet sample, we added 4mL of 1% acetic acid to one tablet weight of ground tablet powder. A low actinic container was flushed with nitrogen and shaken occasionally, then heated for 2 minutes in a 55°C water bath. After the container cooled to room temperature, 2mL of isopropanol was added, for a final volume of 6mL (33% isopropanol). The sample was gently shaken for 2 minutes, then

centrifuged. The vitamins were extracted from 2.5mL of the resulting supernatant using SPME, by immersing the exposed fiber for 40 minutes, with stirring.

The analytes are desorbed from the fiber in the SPME/HPLC interface. The fiber is exposed inside the desorption chamber for 2 minutes, for static desorption with mobile phase, before the valve is opened and dynamic desorption occurs during the run.

Ordering Information:

Description	Cat. No.
SPME Fibers for HPLC, pk. of 3	
50µm CWX/TPR [▲]	57315
65µm PDMS/DVB	57317
SPME Fibers Assortment Kit	
1 fiber each of 5-7315, 5-7317, 5-7301	57323-U
SPME Fiber Holder**	
HPLC holder	57331
SPME/HPLC Interface, includes 2 ferrules	
Valco® valve version	57350-U
Rheodyne® valve version	57353
SUPELCOSIL LC-8 Column	
25cm x 4.6mm ID, 5µm particles	58297
Fat Soluble Vitamins	
100mg, neat, unless otherwise noted. For additional vitamins refer to our catalog.	
Retinol acetate	46958
Retinol palmitate	46959-U
DL-α-Tocopherol	47783
rac-β-Tocopherol, 50mg/mL in 1mL hexane	46401
δ-Tocopherol	47784
γ-Tocopherol, 10mg	47785
DL-α-Tocopherol acetate	47786
Cholecalciferol (D ₃)	47763
Ergocalciferol (D ₂)	47768

*Solid phase microextraction technology licensed exclusively to Supelco. US patent pending; European patent #0523092.

**Initially you must order both holder and fiber assembly. Holder is reusable indefinitely.

*Carbowax®/templated resin.

Trademarks

Carbowax – Union Carbide Corp.
Rheodyne – Rheodyne, Inc.
SUPELCOSIL – Sigma-Aldrich Co.
Valco – Valco Instruments Co., Inc.

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 140

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.

BKK