

Identify the Correct Heptachlor Epoxide Isomer by Capillary GC

Heptachlor epoxide exists in two isomer forms, each having specific behavior characteristics. Often analysts use the wrong isomer as a reference sample in their labwork. Supelco offers both isomers in high purity concentration form.

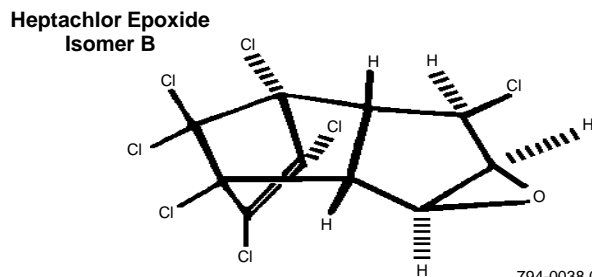
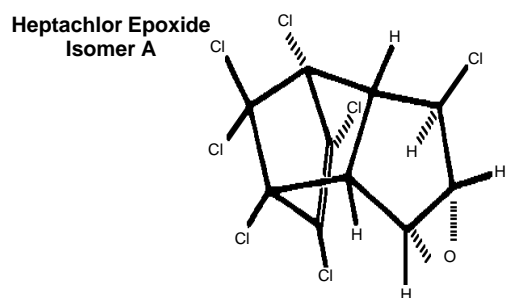
Key Words:

- heptachlor epoxide • isomer

For many decades, heptachlor was used routinely in agriculture as a pest control agent. Within the environment, heptachlor does not remain in its natural state, but metabolizes to form heptachlor epoxide. Heptachlor has been found to convert to heptachlor epoxide in animal species as well as in plants. Heptachlor epoxide is far more stable than heptachlor, and thus is more suitable for environmental testing (1).

Heptachlor epoxide exists in two isomeric forms (Figure A). Isomer A is readily synthesized and is commercially available from a variety of sources. It is water insoluble, and breaks down relatively quickly in nature. Isomer B is the more stable form of heptachlor epoxide. It is of greater concern to environmental scientists because of its ability to persist in the environment. This isomer is not generally available in the commercial market.

Figure A. Heptachlor Epoxide Isomers

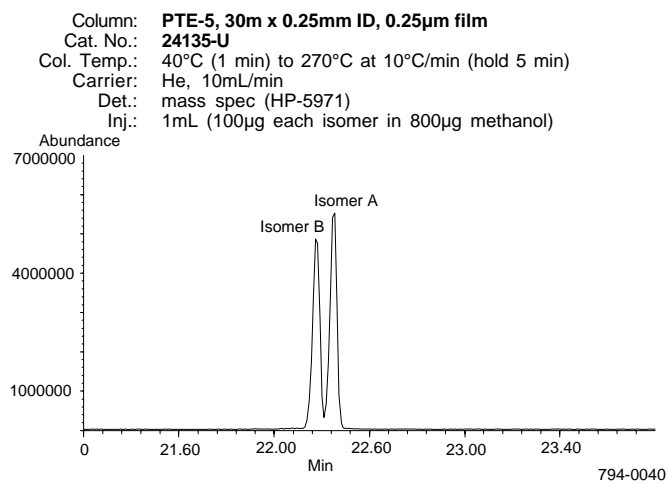


794-0038,0039

Analysts commonly and erroneously use isomer A as a reference sample in their environmental analyses, when isomer B actually is the compound of their concern. Realizing this common laboratory error, Supelco now offers both isomers in high purity concentration form.

Attempts to separate the two heptachlor epoxide isomers using packed column gas chromatography have not been successful. However, capillary GC completes this separation quite readily. We achieved the isomer separation using a 30m x 0.25mm ID x 0.25µm film PTE™ -5 fused silica capillary column, under conditions described in US Environmental Protection Agency Method 8270 for semivolatile organics. The total ion chromatogram (TIC) shows a 90% separation of the isomer peaks (Figure B).

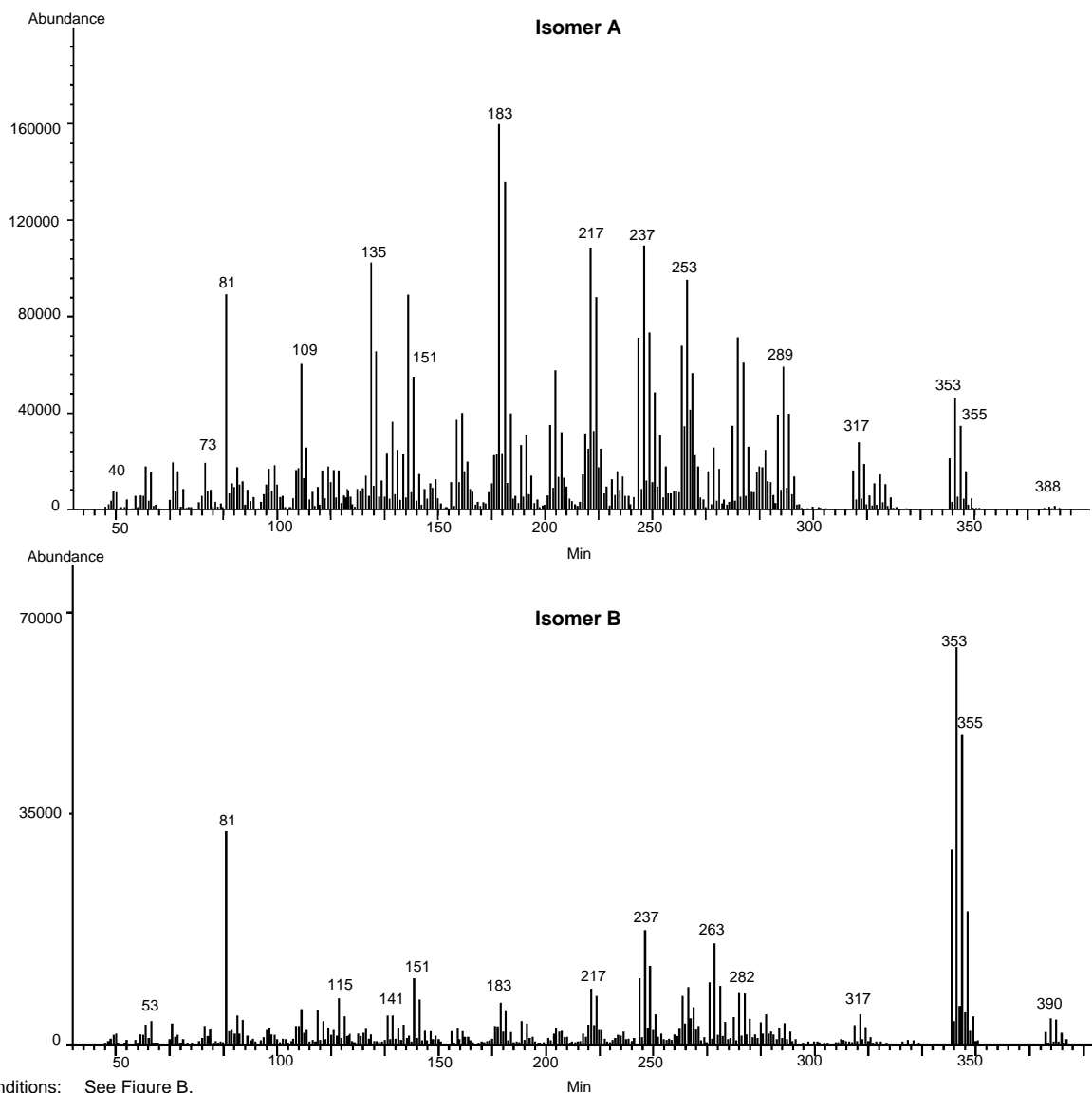
Figure B. Total Ion Chromatogram of Heptachlor Isomers A and B



Minor differences in the retention times of heptachlor epoxide between the analytical standard and the reference standard are sometimes observed. This is often the result of the reference standard isomer selected. If the system is calibrated using isomer A, a difference between an environmental sample retention, typically isomer B, and the calibration retention times is to be expected.

Further demonstration of the difference between the two heptachlor epoxide isomer forms is illustrated through consideration of the fragmentation patterns for these isomers. Fragmentation patterns were collected for the two peaks present in the TIC. The fragmentation patterns of the isomers are distinct and easily differentiated (Figure C). A search of the spectra at the National Institute of Standards and Testing (NIST) library identified the peak for isomer B, but not for isomer A. While it is possible that

Figure C. Fragmentation Patterns of Heptachlor Epoxide Isomers



794-0041,0042

isomer A could be found in the environment, it is more likely that this isomer is not readily observed.

Supelco has identified the important isomer B form of heptachlor epoxide, and made it available in both neat and single-analyte solution form.

If you are analyzing heptachlor epoxide, we strongly recommend that you examine the source of the reference material to determine its isomeric character. Then select the standard most appropriate for your analytical needs.

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

Ordering Information:

Description	Cat. No.
Heptachlor Epoxide Isomer A, 1mL 1000µg/mL in methanol	48198
Heptachlor Epoxide Isomer B, 1mL 1000µg/mL in methanol	40099
PTE-5 Fused Silica Capillary Column 30m x 0.25mm ID, 0.25µm film	24135-U

Reference

- Donnelly, Joseph R., G. Wayne Sovocool, and Richard L. Titus, Structures and Environmental Significance of Heptachlor Epoxide Isomers, Journal of AOAC International, Vol. 76, No. 5 (1993).
- Fused silica columns manufactured under HP US Pat. No. 4,293,415.
PTE is a trademark of Supelco, Inc.

Note 4

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 Vallengbaek 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. · Iloupoli 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

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.