

## Separating Homologs and Polymers by HPLC

Many products contain homologous series of components that must be accurately quantitated to ensure product uniformity. Molecular weight determines which method to use in separating homologs and polymers.

### Key Words:

- homologs • polymers • reversed phase HPLC
- normal phase HPLC • size exclusion chromatography

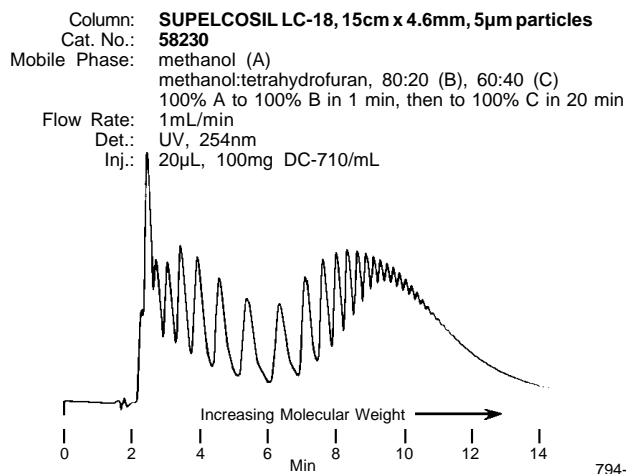
Homologs of many industrial polymers, with molecular weights of up to 10,000 daltons, can be separated by weight on reversed phase or normal phase HPLC columns. For example, Dow Corning® 710 fluid, a polymeric 50% phenyl silicone used as a lubricant or heat exchange fluid, contains a series of homologs of less than 10,000 daltons. The homologs are resolved according to increasing weight by using a 100Å pore reversed phase column, SUPELCOSIL™ LC-18 (octadecylsilyl bonded phase), and a mild gradient. The chromatogram shown in Figure A is typical of reversed phase separations of these homologs.

Normal phase HPLC also can be used to separate polymers with molecular weights of less than 10,000 daltons. This is shown for polyethoxylated octylphenol and nonylphenol surfactants in Figure B. The sample of Triton® X-100, containing an average of 9-10 ethoxy groups, was resolved to baseline on a 25cm SUPELCOSIL LC-Diol column. The surfactant containing 30 ethylene oxide units was analyzed within 40 minutes on the same column.

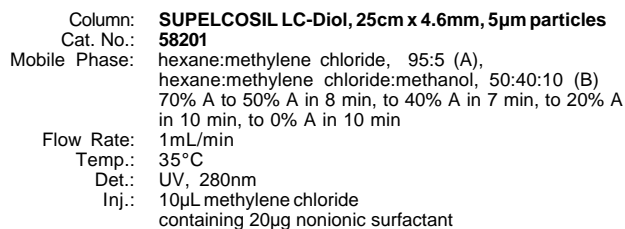
Homologs with molecular weights greater than 10,000 daltons are usually separated on high performance size exclusion chromatography (SEC) columns. The exclusion limits of SEC columns depend on the size of the pores in the packing particles, with larger pores admitting larger molecules into these particles. By combining, in series, columns having different pores sizes, analysts can expand the range of molecular weights that can be separated at one time. In Figure C, polystyrene standards having molecular weights from less than 1000 to over  $1 \times 10^6$  daltons were sharply separated on a 100Å pore and a 300Å pore column (trimethylsilyl bonded phase). Similarly, these SUPELCOSIL 100Å and 300Å pore columns in series (always install the smaller pore size column in front of the larger pore size column) will satisfactorily separate other polymers with molecular weights between  $10^3$  and  $10^6$  daltons. Figure D depicts a typical calibration curve for this pair of columns.

SUPELCOSIL size exclusion chromatography columns contain 5 micron spherical silica particles with 100Å or 300Å diameter pores. For SUPELCOSIL LC-1 and LC-301 columns, the silica is bonded with a trimethylsilyl bonded phase. The dimensions of these HPSEC columns, 25cm x 6.2mm, provide a good compromise between large sample capacity and the most rapid analysis time. Fittings for 1/16" OD tubing are included with each column.

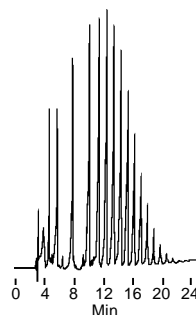
**Figure A. Phenyl Silicone Homologs (<10,000 daltons mol. wt.) by Reversed Phase HPLC**



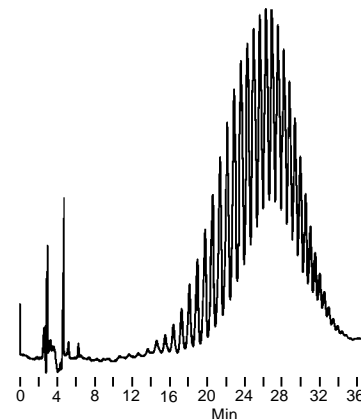
**Figure B. Nonionic Surfactants by Normal Phase HPLC**



**Triton X-100: 9-10 Ethoxy Groups (Octylphenol)**



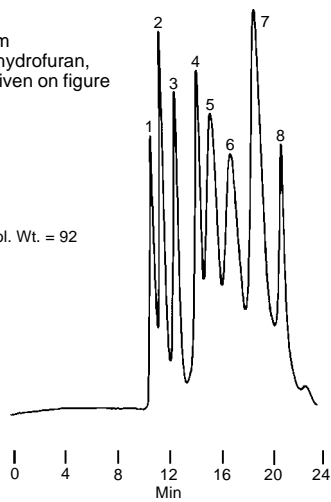
**Tergitol: 30 Ethoxy Groups (Nonylphenol)**



794-0875, 0876

**Figure C. Polystyrene Standards Using Size Exclusion Columns in Series**

Columns: **SUPELCO SIL LC-1, 25cm x 6.2mm, and SUPELCO SIL LC-301, 25cm x 6.2mm, in series**  
 Cat. Nos.: **58961**  
**58966**  
 Mobile Phase: tetrahydrofuran  
 Flow Rate: 1mL/min  
 Temp.: 45°C  
 Det.: UV, 254nm  
 Inj.: 10µL tetrahydrofuran, amounts given on figure

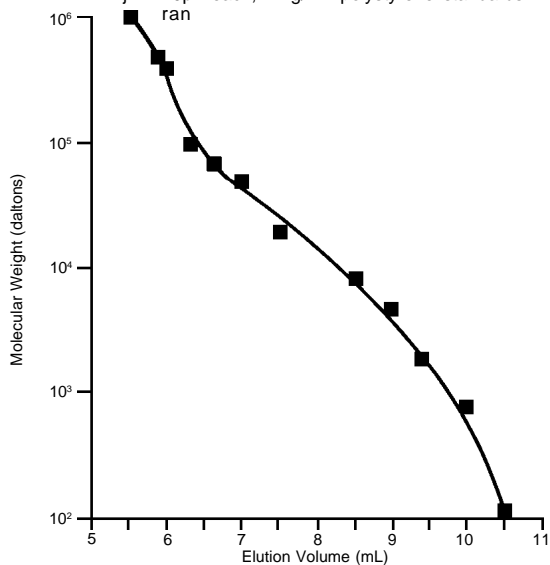


1. 1,800,000 (3.3µg)
2. 300,000 (3.3µg)
3. 100,000 (3.3µg)
4. 35,000 (5.0µg)
5. 17,500 (6.7µg)
6. 9,000 (6.7µg)
7. 2,000 (10.0µg)
8. Toluene (1.7µg) Mol. Wt. = 92

713-0781

**Figure D. 100Å Pore and 300Å Pore HPSEC Columns in Series Separate Homologs Having a Wide Molecular Weight Range**

Columns: **SUPELCO SIL LC-1, 25cm x 6.2mm, and SUPELCO SIL LC-301, 25cm x 6.2mm, in series**  
 Cat. Nos.: **58961**  
**58966**  
 Mobile Phase: tetrahydrofuran  
 Flow Rate: 1mL/min  
 Temp.: 45°C  
 Det.: UV, 254nm  
 Inj.: 20µL each, 2mg/mL polystyrene standards in tetrahydrofuran



794-0877

**Size Exclusion Columns (25cm x 6.2mm ID, 5µm packings)**

SUPELCO SIL Column	Pore Diameter	Cat. No.
<b>Silica</b>		
LC-Si	100Å	<b>58963</b>
LC-3Si	300Å	<b>58965</b>
<b>Trimethylsilyl Bonded Phase</b>		
LC-1	100Å	<b>58961</b>
LC-301	300Å	<b>58966</b>
<b>Diol Bonded Phase</b>		
LC-Diol	100Å	<b>58967</b>
LC-3Diol	300Å	<b>58968</b>

**SUPELCO SIL LC-18 and LC-Diol Columns (5µm packing, 100Å Pores)**

Dimensions	Cat. No.
<b>SUPELCO SIL LC-18 Modular Column</b>	
25cm x 4.6mm	<b>58298</b>
Supelguard LC-18 Columns, pk. of 2	<b>59564</b>
<b>SUPELCO SIL LC-Diol Modular Column</b>	
25cm x 4.6mm	<b>58201</b>
Supelguard LC-Diol Columns, pk. of 2	<b>59569</b>
<b>Cartridge Guard Column Holder</b>	<b>55205</b>
(connects guard column directly to modular column)	

**Mono-Disperse Polystyrene Standards (500mg bottles)**

Nominal Molecular Wt.	M <sub>w</sub> /M <sub>n</sub>	Cat. No.
800	1.30	<b>45700</b>
4,000	1.06	<b>45702</b>
35,000	1.06	<b>45705</b>
50,000	1.06	<b>45706</b>
100,000	1.06	<b>45707</b>
233,000	1.06	<b>45708</b>
300,000	1.06	<b>45709</b>
600,000	1.10	<b>45710</b>
900,000	1.10	<b>45711</b>

**Polystyrene Standards Kits, 250mg of each polymer**

<b>Low Molecular Weight Standard Kit</b>		<b>48937</b>
2500 MW	17,500 MW	
5000 MW	30,000 MW	
9000 MW	50,000 MW	
<b>High Molecular Weight Standard Kit</b>		<b>48938</b>
110,000 MW	600,000 MW	
220,000 MW	900,000 MW	
400,000 MW	1,800,000 MW	

**Trademarks**

Dow Corning — Dow Corning Corp.  
 SUPELCO SIL — Supelco, Inc.  
 Supelguard — Supelco, Inc.  
 Triton — Union Carbide

*Note 46*

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](http://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 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 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](mailto: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.