

Application

Note
78

Monitor Antioxidant Additives in Foods, Using HPLC (AOAC Method 983.15) or Capillary GC and a Supelco Reference Standards Kit

Supelco offers an antioxidant standards kit containing all 9 antioxidants listed in Association of Official Analytical Chemists (AOAC) Official Method 983.15: *Phenolic Antioxidants in Oils, Fats, and Butter Oil*, plus ethoxyquin, an additional commonly used antioxidant. These compounds are separated well, in less than 13 minutes, on a SUPELCOSIL LC-18 HPLC column, under the conditions described in the AOAC method. Several of the antioxidants also can be resolved on a SAC-5 capillary GC column.

Key Words:

- antioxidants • phenolic antioxidants • food preservatives

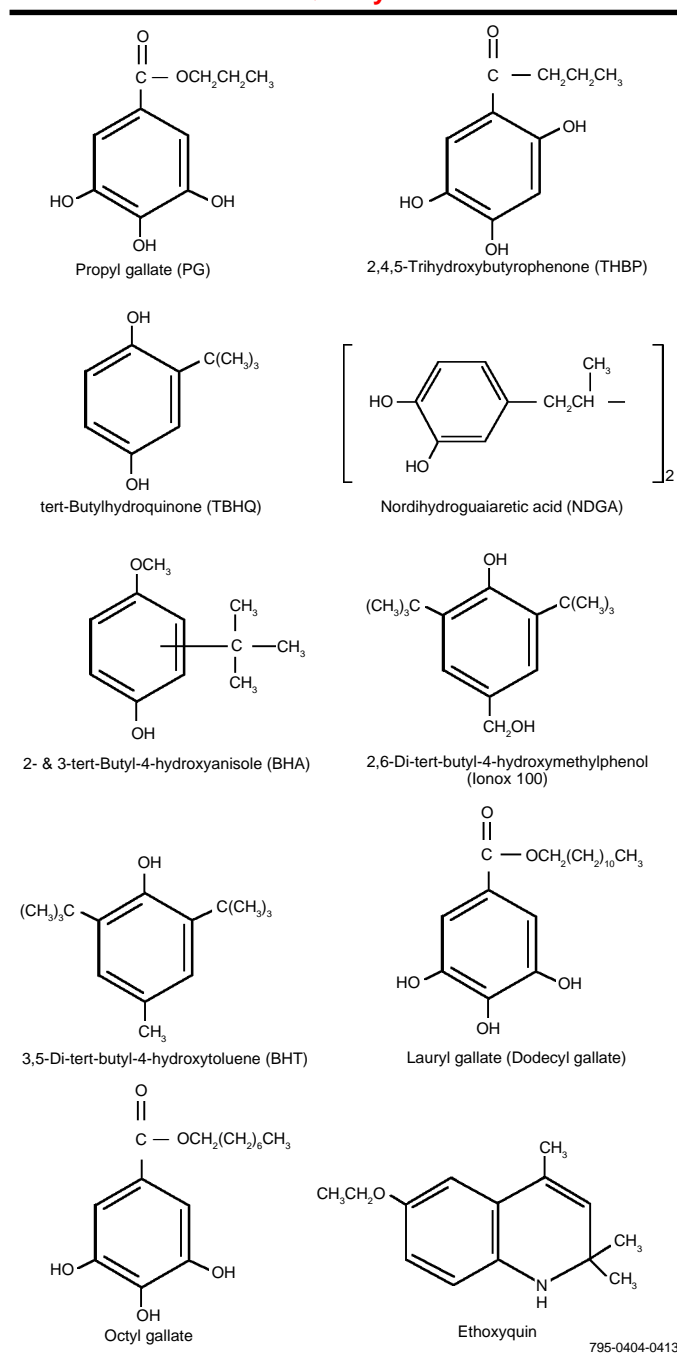
Rancidity caused by the oxidation of unsaturated fats in foods is a major problem for the food industry. With the trace metals present in foods acting as catalysts, the unsaturated fats react with oxygen, producing free radicals and peroxides. These compounds can destroy many vitamins, reducing the nutritional value of the food, destroy pigments, bleaching the food, and cause off-odors and off-flavors. While the portion of fat undergoing such changes can be small, the resulting off-odors are strong, and very much out of proportion to the amount of oil or fat involved.

Antioxidants are added to foods and other products to prevent rancidity. Although the exact mechanism(s) is (are) not clear, antioxidants react with free radicals and peroxides, slowing the development of rancidity. Certain other additives greatly enhance the effectiveness of antioxidants. Metal scavenger and chelating agents, such as citric acid and citrates, tie up the trace metals and greatly reduce their catalytic activity. Synergism between antioxidants also has been noted, and many commercial antioxidant mixes are formulated to contain mixtures of antioxidants. The most common of these mixtures contain both butyl-hydroxyanisole (BHA) and butyl-hydroxytoluene (BHT).

In order to assure consistent product quality, these additives must be monitored. In the past, food analysts had to go to several vendors to obtain good quality standards of all of the antioxidants they might be required to monitor. Now, Supelco offers an antioxidant standards kit, containing all 9 of the antioxidants listed in Association of Official Analytical Chemists (AOAC) Official Method 983.15: *Phenolic Antioxidants in Oils, Fats, and Butter Oil*, (Figure A) plus ethoxyquin, an additional antioxidant commonly used in spices and other food products, and in cosmetics.

AOAC Method 983.15 describes a linear gradient HPLC analysis for resolving the nine antioxidants shown in Figure A. Figure B shows a chromatogram generated on a 15cm x 4.6mm SUPELCOSIL™ LC-18 HPLC column, using the conditions described in the AOAC

Figure A. Phenolic Antioxidants Commonly Used to Preserve Food Quality



795-0404-0413

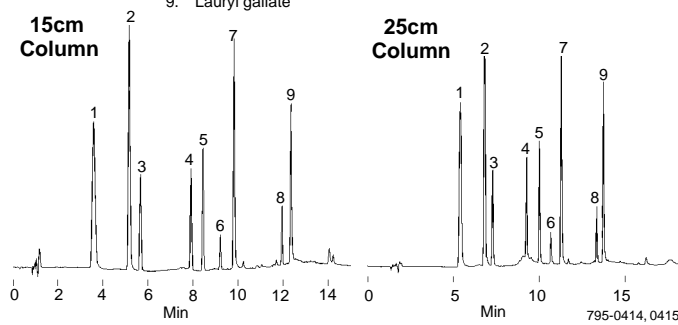
method. All 9 analytes were well resolved in less than 13 minutes. Ethoxyquin is not detected at the 280nm UV wavelength indicated by the AOAC method, and thus is not included here. Figure B also shows the analysis on a 25cm SUPELCOSIL LC-18 column. The analytes again are well resolved, and analysis time is only slightly prolonged.

In most food products, only 1 or 2 antioxidants are used. In some cases, gas chromatography can be used to monitor the antioxi-

Figure B. Antioxidants in AOAC Method 983.15 by HPLC

Column: **SUPELCOSIL LC-18, 15cm or 25cm x 4.6mm ID, 5µm particles**
 Cat. No.: **58230-U** (15cm column), **58298** (25cm column)
 Mobile Phase: A = 5% acetic acid in DI water
 B = acetonitrile:methanol, 50:50
 30% B to 100% B in 10 min, hold 10 min
 Flow Rate: 2mL/min
 Temp.: ambient
 Det.: UV, 280nm
 Inj.: 10µL, 20µg/mL each analyte

1. Propyl gallate
2. 2,4,5-Trihydroxybutyrophenone
3. tert-Butylhydroquinone
4. Nordihydroguaiaric acid
5. 2- & 3-tert-Butyl-4-hydroxyanisole
6. 2,6-Di-tert-butyl-4-hydroxymethylphenol (Ionox 100)
7. Octyl gallate
8. 3,5-Di-tert-butyl-4-hydroxytoluene
9. Lauryl gallate

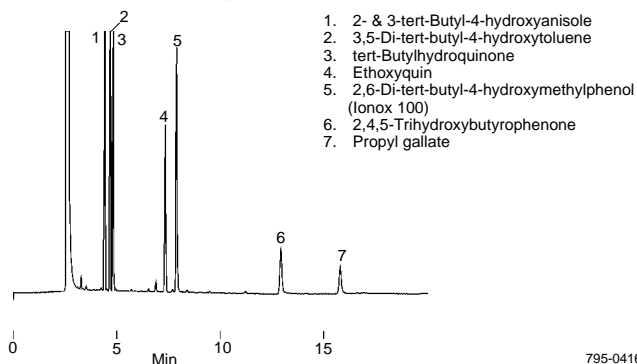


dants. (Many of the antioxidants in Figure A give a poor response to FID detection, however.) Figure C shows the chromatogram generated from a mix of 7 antioxidants, including ethoxyquin, separated on a 30m x 0.25mm x 0.25µm phase film SACTM-5 capillary GC column under isothermal (200°C) conditions. Resolution is good. Ethoxyquin was eluted at approximately 7 minutes. Note that at 200°C, an analysis time of more than 15 minutes was needed to elute propyl gallate. Lauryl gallate and octyl gallate cannot be monitored under these conditions – a very high oven temperature is required to elute these high boiling/low volatility compounds from the capillary column.

If your analyses include monitoring antioxidants, this standards kit should meet all your routine needs, and will prepare you for identifying unanticipated additional antioxidants in your samples. Note that the ethoxyquin standard is qualitative. Because of its highly reactive nature, we cannot assure a purity level of more than 70% for this compound. Each antioxidant is packaged neat, under nitrogen.

Figure C. Antioxidants with Lower Boiling Points Can Be Monitored by GC

Column: **SAC-5, 30m x 0.25mm, 0.25µm film**
 Cat. No.: **24156**
 Oven: 200°C
 Carrier: helium, 30cm/sec (set at 200°C)
 Det.: FID (250°C)
 Inj.: 2µL, 200µg/mL each analyte, split 100:1 (250°C)



1. 2- & 3-tert-Butyl-4-hydroxyanisole
2. 3,5-Di-tert-butyl-4-hydroxytoluene
3. tert-Butylhydroquinone
4. Ethoxyquin
5. 2,6-Di-tert-butyl-4-hydroxymethylphenol (Ionox 100)
6. 2,4,5-Trihydroxybutyrophenone
7. Propyl gallate

Ordering Information:

Phenolic Antioxidants Kit 47192

Individually packaged under nitrogen, neat, in quantities listed below.

Component	CAS No.	Qty. (mg)
Propyl gallate (PG)	121-79-9	500
2,4,5-Trihydroxybutyrophenone (THBP)	1421-63-2	500
tert-Butylhydroquinone (TBHQ)	1948-33-0	500
Nordihydroguaiaric acid (NDGA)	500-38-9	100
2- & 3-tert-Butyl-4-hydroxyanisole (BHA)	25013-16-5	500
2,6-Di-tert-butyl-4-hydroxymethylphenol (Ionox 100)	88-26-6	100
3,5-Di-tert-butyl-4-hydroxytoluene (BHT)	128-37-0	500
Lauryl gallate (Dodecyl gallate)	1166-52-5	500
Octyl gallate	1034-01-1	500
Ethoxyquin*	91-53-2	500

*Qualitative standard. Purity >70% cannot be assured.

SUPELCOSIL LC-18 HPLC Columns

Description	Cat. No.
15cm x 4.6mm, 5µm particles	58230-U
25cm x 4.6mm, 5µm particles	58298

SAC-5 Capillary GC Column

30m x 0.25mm, 0.25µm film	24156
---------------------------	--------------

Trademarks

SAC, SUPELCO, SUPELCOSIL – Supelco, Inc.

Reference

1. *Official Methods of Analysis* (16th ed.), Method 983.15 Association of Official Analytical Chemists, Arlington, VA USA (1995).

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

Nete 78

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-Aldrich 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

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.