

## 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

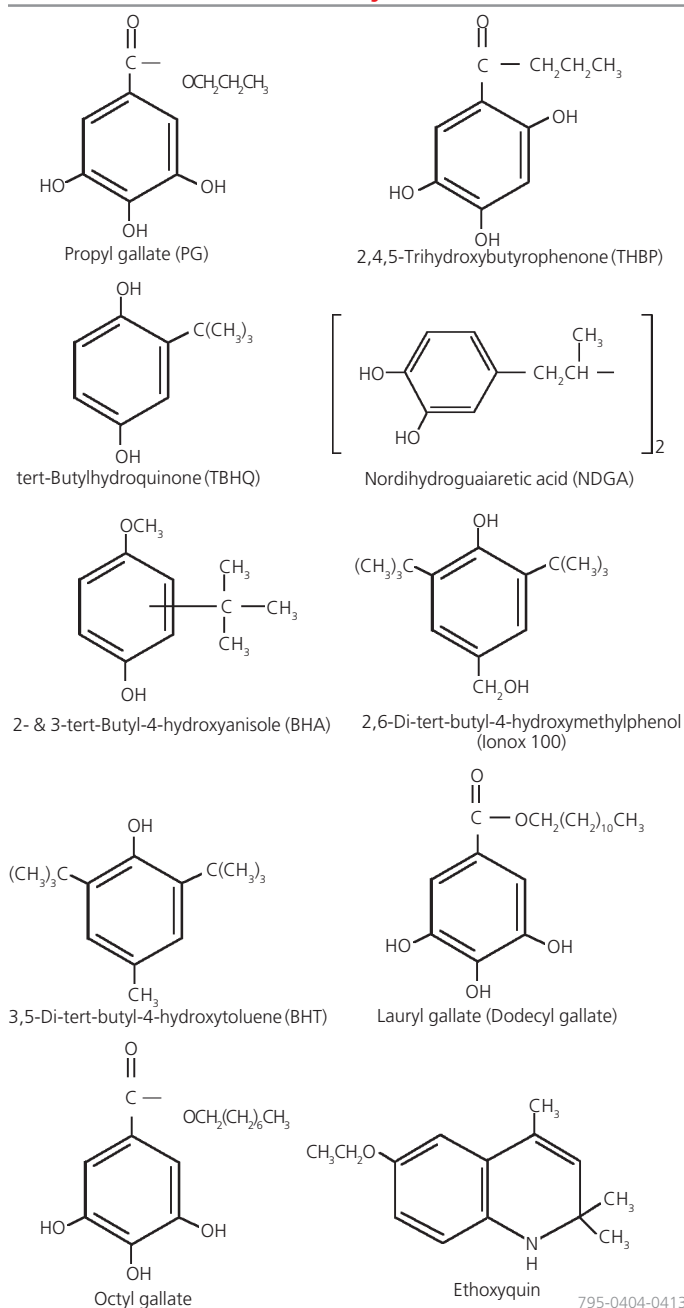
Rancidity caused by the oxidation of unsaturated fats in foods is a major problem for the food industry. With the trace metals present in foods acting as catalysts, the unsaturated fats react with oxygen, producing free radicals and peroxides. These compounds can destroy many vitamins, reducing the nutritional 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

**Figure A. Phenolic Antioxidants Commonly Used to Preserve Food Quality**



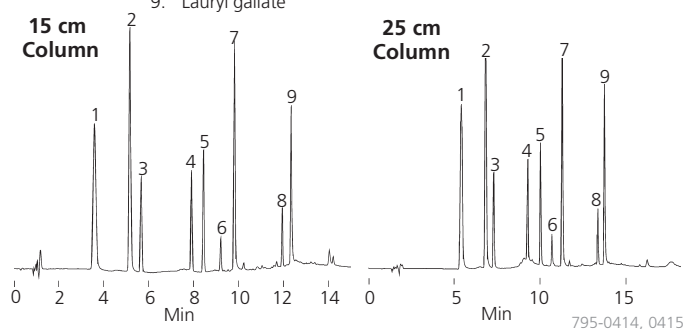
795-0404-0413

a chromatogram generated on a 15 cm x 4.6 mm SUPELCOSIL™ LC-18 HPLC column, using the conditions described in the AOAC method. All 9 analytes were well resolved in less than 13 minutes. Ethoxyquin is not detected at the 280 nm UV wavelength indicated by the AOAC method, and thus is not included here. Figure B also shows the analysis on a 25 cm SUPELCOSIL LC-18 column. The analytes again are well resolved, and analysis time is only slightly prolonged.

**Figure B. Antioxidants in AOAC Method 983.15 by HPLC**

column: SUPELCOSIL LC-18, 15 cm or 25 cm x 4.6 mm I.D., 5 µm particles (**58230-U** (15 cm column), **58298** (25 cm column))  
 mobile phase A: 5% acetic acid in DI water  
 mobile phase B: acetonitrile:methanol, 50:50  
 30% B to 100% B in 10 min., hold 10 min  
 flow rate: 2 mL/min.  
 temp.: ambient  
 det.: UV, 280 nm  
 injection: 10 µL, 20 µg/mL each analyte

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

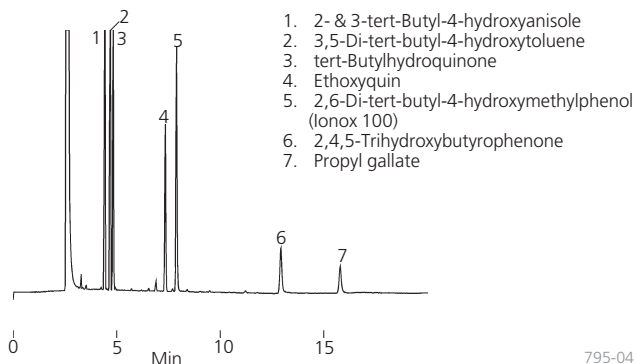


In most food products, only 1 or 2 antioxidants are used. In some cases, gas chromatography can be used to monitor the antioxidants. (Many of the antioxidants in Figure A give a poor response to FID detection, however.) Figure C shows the chromatogram generated from a mix of 7 antioxidants, including ethoxyquin, separated on a 30 m x 0.25 mm x 0.25 µm phase film SAC™-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, 30 m x 0.25 mm, 0.25 µm film (24156)  
 oven: 200 °C  
 det.: FID (250 °C)  
 carrier gas: helium, 30 cm/sec. (set at 200 °C)  
 injection: 2 µL, 200 µg/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 (lonox 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       |
| Nordihydroguaiaretic acid (NDGA)                    | 500-38-9   | 100       |
| 2- & 3-tert-Butyl-4-hydroxyanisole (BHA)            | 25013-16-5 | 500       |
| 2,6-Di-tert-butyl-4-hydroxymethylphenol (lonox 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.       |
|--------------------------------|----------------|
| 15 cm x 4.6 mm, 5 µm particles | <b>58230-U</b> |
| 25 cm x 4.6 mm, 5 µm particles | <b>58298</b>   |

**SAC-5 Capillary GC Column**

|                              |              |
|------------------------------|--------------|
| 30 m x 0.25 mm, 0.25 µm film | <b>24156</b> |
|------------------------------|--------------|

**Trademarks**

SAC, SUPELCO, SUPELCOSIL – Sigma-Aldrich Co.

**Reference**

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

Order/Customer Service 800-247-6628, 800-325-3010  
 Fax 800-325-5052 • E-mail supelco@sial.com

Technical Service 800-359-3041, 814-359-3041  
 Fax 800-359-3044, 814-359-5468 • E-mail techservice@sial.com

SUPELCO • 595 North Harrison Road, Bellefonte, PA 16823-0048 • 814-359-3441

We are committed to the success of our Customers, Employees and Shareholders through leadership in Life Science, High Technology and Service.  
 The SIGMA-ALDRICH Family SIGMA ALDRICH Fluka Riedel-deHaën SUPELCO ISOTEK

©2004 Sigma-Aldrich Co. Printed in USA. Supelco brand products are sold through Sigma-Aldrich, Inc. Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip. Supelco is a registered trademark of Sigma-Aldrich Co.  
 ARM T395078

**SUPELCO**  
 sigma-aldrich.com/supelco