

# PRODUCT INFORMATION



## Oxidized Lipid HPLC Mixture

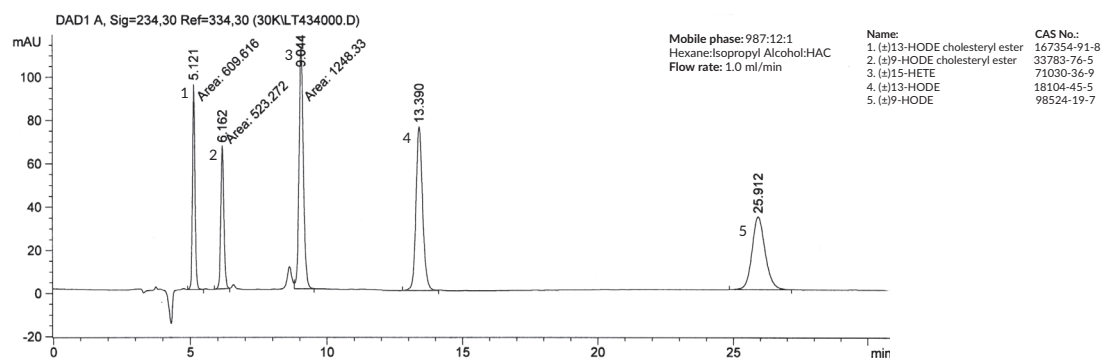
Item No. 34004

**Purity:** ≥98% for each compound  
**Supplied as:** A solution in ethanol (5 µg of each compound)  
**Storage:** -20°C  
**Stability:** ≥1 year

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

The oxidized lipid HPLC mixture contains a combination of oxidized fatty acids found in atherosclerotic plaques supplied in ethanol. This HPLC mixture contains 5 µg of the following oxidized fatty acids: (±)9-HODE, (±)13-HODE, (±)9-HODE cholesteryl ester, (±)13-HODE cholesteryl ester, and (±)15-HETE.



#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/08/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM