

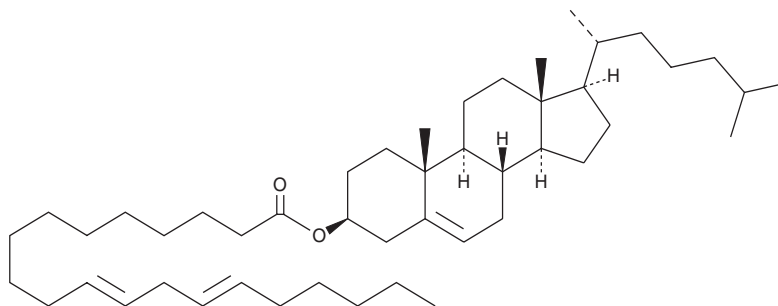
# PRODUCT INFORMATION



## Cholesteryl 11,14-Eicosadienoate

Item No. 26467

**CAS Registry No.:** 70110-48-4  
**Formal Name:** cholest-5-en-3 $\beta$ -ol,  
11,14-eicosadienoate  
**Synonyms:** 20:2 (11,14) CE,  
20:2 (11,14) Cholesterol ester  
**MF:** C<sub>47</sub>H<sub>80</sub>O<sub>2</sub>  
**FW:** 677.1  
**Purity:**  $\geq$ 95%  
**Supplied as:** A neat oil  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

Cholesteryl 11,14-eicosadienoate is supplied as a neat oil. A stock solution may be made by dissolving the cholesteryl 11,14-eicosadienoate in the solvent of choice. Cholesteryl 11,14-eicosadienoate is soluble in the organic solvent chloroform, which should be purged with an inert gas, at a concentration of approximately 10 mg/ml.

### Description

Cholesteryl 11,14-eicosadienoate is a cholesterol ester.<sup>1</sup> It has been used as a synthetic intermediate in the synthesis of cholesterol methyl esters.

### Reference

1. Tuckey, R.C. and Stevenson, P.M. Methanolysis of cholesteryl esters: Conditions for quantitative preparation of methyl esters. *Anal. Biochem.* **94**(2), 402-408 (1979).

**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, 12/06/2022

### 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