

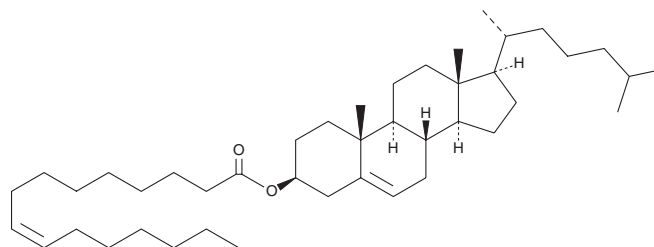
PRODUCT INFORMATION



Cholesteryl Palmitoleate

Item No. 26474

CAS Registry No.: 16711-66-3
Formal Name: (3 β)-cholest-5-en-3-ol 3-[(9Z)-9-hexadecenoate]
Synonyms: 16:1(9Z) CE, CE(16:1), 16:1(9Z) Cholesterol ester, Cholesterol Palmitoleate
MF: C₄₃H₇₄O₂
FW: 623.1
Purity: \geq 95%
Supplied as: A low melting solid
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 palmitoleate is supplied as a neat oil. A stock solution may be made by dissolving the cholesteryl palmitoleate in the solvent of choice, which should be purged with an inert gas. Cholesteryl palmitoleate is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

Description

Cholesteryl palmitoleate is a cholesterol ester.¹⁻³ Plasma levels of cholesteryl palmitoleate are increased in *ApoE*^{-/-} mice exposed to cigarette smoke and in pediatric patients with biliary atresia.^{1,2} Cholesteryl palmitoleate has been used as a standard for the identification of cholesterol esters in human meibomian gland secretions.³

References

1. Boué, S., Tarasov, K., Jänis, M., *et al.* Modulation of atherogenic lipidome by cigarette smoke in apolipoprotein E-deficient mice. *Atherosclerosis* **225**(2), 328-334 (2012).
2. Robberecht, E., Koletzko, B., and Christophe, A. Several mechanisms contribute to the abnormal fatty acid composition of serum phospholipids and cholesterol esters in cholestatic children with extrahepatic biliary atresia. *Prostaglandins Leukot. Essent. Fatty Acids* **56**(3), 199-204 (1997).
3. Butovich, I.A. Fatty acid composition of cholesteryl esters of human meibomian gland secretions. *Steroids* **75**(10), 726-733 (2010).

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, 12/07/2022

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