

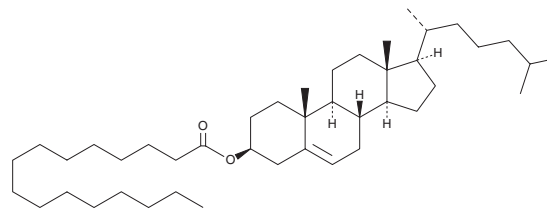
PRODUCT INFORMATION



Cholesteryl Palmitate

Item No. 26473

CAS Registry No.: 601-34-3
Formal Name: (3β)-cholest-5-en-3-ol, 3-hexadecanoate
Synonyms: CE(16:0), 16:0(CE), 16:0 Cholesterol ester, Cholesterol Palmitate, NSC 59692
MF: C₄₃H₇₆O₂
FW: 625.1
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

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

Description

Cholesteryl palmitate is a cholesterol ester. It is a major cholesterol ester found in human meibum and has been used to form a stable tear film to study amphiphilic block copolymers as surfactants for dry eyes.¹ Cholesteryl palmitate levels in amniotic fluid are decreased in pregnancies in which infants later developed respiratory distress syndrome and in patients with well-controlled diabetes.^{2,3} Levels of cholesteryl palmitate are increased in bronchoalveolar lavage fluid (BALF) from patients with chronic interstitial pneumonia.⁴

References

1. Shine, W.E. and McCulley, J.P. Polar lipids in human meibomian gland secretions. *Curr. Eye Res.* **26(2)**, 89-94 (2003).
2. Ludmir, J., Alvarez, J.G., Mennuti, M.T., *et al.* Cholesteryl palmitate as a predictor of fetal lung maturity. *Am. J. Obstet. Gynecol.* **157(1)**, 84-88 (1987).
3. Ludmir, J., Alvarez, J.G., Landon, M.B., *et al.* Amniotic fluid cholesteryl palmitate in pregnancies complicated by diabetes mellitus. *Obstet. Gynecol.* **73(3 Pt.1)**, 360-362 (1988).
4. Fukuhara, N., Tachihara, M., Tanino, Y., *et al.* Cholesteryl palmitate crystals in bronchoalveolar lavage fluid smears as a possible prognostic biomarker for chronic interstitial pneumonia: A preliminary study. *Respir. Investig.* **54(2)**, 109-115 (2016).

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