

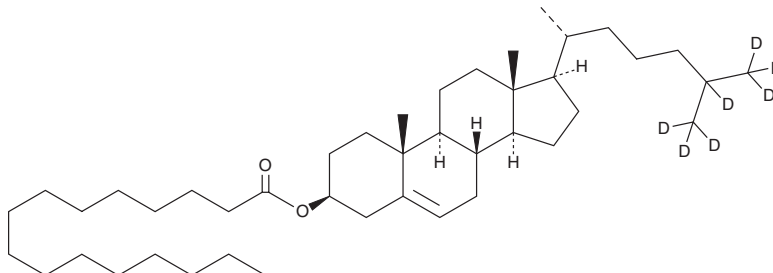
PRODUCT INFORMATION



Cholesteryl-d₇ Palmitate

Item No. 9000422

CAS Registry No.: 1416275-32-5
Formal Name: (3β)-cholest-5-en-25,26,26,26,27,27,27-d₇-3-ol, 3-hexadecanoate
Synonyms: 16:0 Cholesterol-d₇ ester, Cholesterol-d₇ Palmitate, 16:0(CE-d₇), CE-d₇(16:0)
MF: C₄₃H₆₉D₇O₂
FW: 632.1
Chemical Purity: ≥95% (Cholesteryl Palmitate)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₇); ≤1% d₀
Supplied as: A 10 mg/ml solution in chloroform
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-d₇ palmitate is intended for use as an internal standard for the quantification of cholesteryl palmitate (Item No. 26473) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

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, 11/24/2025

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