

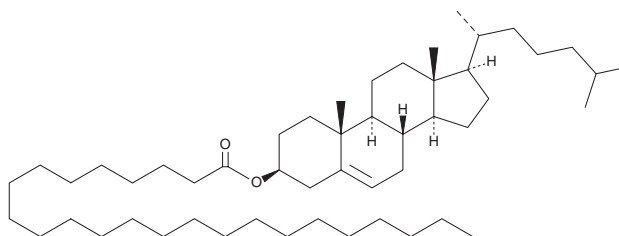
PRODUCT INFORMATION



Cholesteryl Lignocerate

Item No. 26470

CAS Registry No.: 73024-96-1
Formal Name: cholest-5-en-(3 β)-ol, 3-tetracosanoate
Synonyms: 24:0 CE, 24:0 Cholesterol ester
MF: C₅₁H₉₂O₂
FW: 737.3
Purity: \geq 98%
Supplied as: A 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 lignocerate is supplied as a solid. A stock solution may be made by dissolving the cholesteryl lignocerate in the solvent of choice. Cholesteryl lignocerate 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 lignocerate is a cholesterol ester that has been found in human meibum.¹ Cholesteryl lignocerate can be hydrolyzed by cellular extracts from cultured human skin fibroblasts isolated from healthy individuals but not patients with cholesteryl ester storage disease (CESD) or Wolman disease.²

References

1. Chen, J., Green, K.B., and Nichols, K.K. Quantitative profiling of major neutral lipid classes in human meibum by direct infusion electrospray ionization mass spectrometry. *Invest. Ophthalmol. Vis. Sci.* **54**(8), 5730-5753 (2013).
2. Michels, V.V. and Beaudet, A.L. Cholesteryl lignocerate hydrolysis in adrenoleukodystrophy. *Pediatr. Res.* **14**(1), 21-23 (1980).

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, 12/19/2022

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