

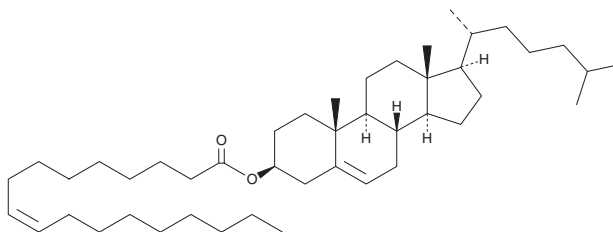
PRODUCT INFORMATION



Cholesteryl Oleate

Item No. 26472

CAS Registry No.: 303-43-5
Formal Name: cholest-5-en-3 β -ol, 3-[(9Z)-9-octadecenoate]
Synonyms: CE(18:1), 18:1(9Z) CE,
18:1(9Z) Cholesterol ester, NSC 18186
MF: C₄₅H₇₈O₂
FW: 651.1
Purity: \geq 95%
Supplied as: A crystalline 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 oleate is supplied as a crystalline solid. A stock solution may be made by dissolving the cholesteryl oleate in the solvent of choice, which should be purged with an inert gas. Cholesteryl oleate is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

Description

Cholesteryl oleate is a cholesterol ester.^{1,2} It has been used in the generation of nanoparticles and liposomes for drug delivery.

References

1. Abd-Rabou, A.A., Bharali, D.J., and Mousa, S.A. Taribavirin and 5-fluorouracil-loaded pegylated-lipid nanoparticle synthesis, p38 docking, and antiproliferative effects on MCF-7 breast cancer. *Pharm. Res.* **35(4)**, 76 (2018).
2. Wang, J., Kang, Y.-X., Pan, W., *et al.* Enhancement of anti-inflammatory activity of curcumin using phosphatidylserine-containing nanoparticles in cultured macrophages. *Int. J. Mol. Sci.* **17(5)**, E969 (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, 12/05/2022

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