

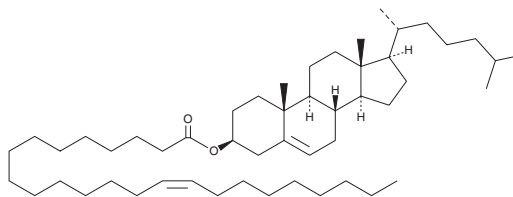
PRODUCT INFORMATION



Cholesteryl Nervonate

Item No. 26471

CAS Registry No.: 60758-73-8
Formal Name: cholest-5-en-3-ol, 3 β -[(15Z)-15-tetracosenoate]
Synonyms: C24:1-CE, 24:1 Cholesterol ester
MF: C₅₁H₉₀O₂
FW: 735.3
Purity: \geq 98%
Supplied as: A 50 mg/ml solution in chloroform
Storage: -20°C
Stability: \geq 4 years



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

Description

Cholesteryl nervonate is a cholesterol ester composed of cholesterol (Item No. 9003100) and nervonic acid (Item No. 13940).¹ It has been found in human meibum. Cholesteryl nervonate has been used in the generation of liposomes.²

References

1. Butovich, I.A. Cholesteryl esters as a depot for very long chain fatty acids in human meibum. *J. Lipid. Res.* **50(3)**, 501-513 (2009).
2. Miyazaki, Y., Hara-Hotta, H., Matsuama, T., *et al.* Hemolysis of phosphatidylcholine-containing erythrocytes by serratamic acid from *Serratia marcescens*. *Int. J. Biochem.* **25(7)**, 1033-1038 (1992).

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, 01/29/2026

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