

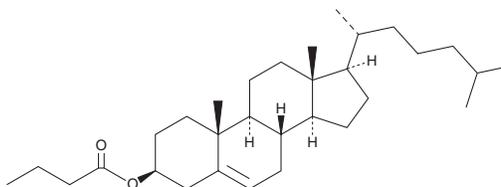
PRODUCT INFORMATION



Cholesteryl Butyrate

Item No. 40950

CAS Registry No.: 521-13-1
Formal Name: cholest-5-en-3 β -ol, 3-butanoate
Synonyms: Chol-Butyrate, Cholesterol *n*-butyrate
MF: C₃₁H₅₂O₂
FW: 456.8
Purity: \geq 95%
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 butyrate is supplied as a solid. A stock solution may be made by dissolving the cholesteryl butyrate in the solvent of choice, which should be purged with an inert gas. Cholesteryl butyrate is sparingly soluble (1-10 mg/ml) in chloroform.

Description

Cholesteryl butyrate is a cholesterol ester that has been used with other lipids in the generation of lipid nanoparticles (LNPs).¹ LNPs containing cholesteryl butyrate decrease proliferation, induce cell death, as well as decrease c-Myc levels in cancer cells.

Reference

1. Serpe, L., Laurora, S., Pizzimenti, S., *et al.* Cholesteryl butyrate solid lipid nanoparticles as a butyric acid pro-drug: Effects on cell proliferation, cell-cycle distribution and c-myc expression in human leukemic cells. *Anticancer Drugs* **15(5)**, 525-536 (2004).

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, 05/28/2024

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