

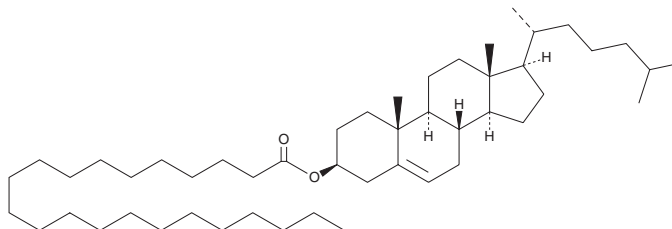
PRODUCT INFORMATION



Cholesteryl Behenate

Item No. 26466

CAS Registry No.: 61510-09-6
Formal Name: (3 β)-cholest-5-en-3-ol
3-docosanoate
Synonym: Cholesteryl Docosanoate
MF: C₄₉H₈₈O₂
FW: 709.2
Purity: \geq 98%
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 behenate is supplied as a crystalline solid. A stock solution may be made by dissolving the cholesteryl behenate in the solvent of choice. Cholesteryl behenate 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 behenate is a cholesterol ester.¹ It has been used as an internal standard for the quantification of cholesterol esters in human meibum samples.

Reference

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).

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, 09/30/2022

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