

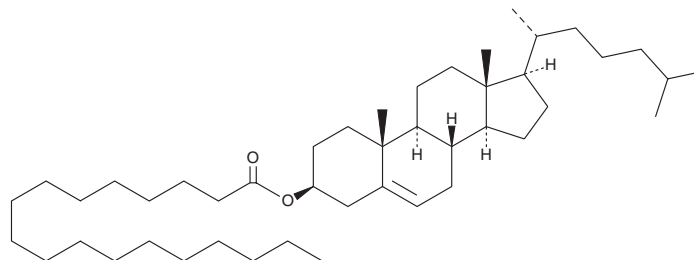
PRODUCT INFORMATION



Cholesteryl Stearate

Item No. 26475

CAS Registry No.: 35602-69-8
Formal Name: (3 β)-cholest-5-en-3-ol,
3-octadecanoate
Synonyms: 18:0 CE, 18:0 Cholesterol ester,
CE(18:0), Cholesterol Stearate,
NSC 59693
MF: C₄₅H₈₀O₂
FW: 653.1
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 stearate is supplied as a crystalline solid. A stock solution may be made by dissolving the cholesteryl stearate in the solvent of choice. Cholesteryl stearate 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 stearate is a cholesterol ester.¹ It has been used as an internal standard for the quantification of sterol esters in various plant tissues by quadropole time-of-flight (Q-TOF) MS. Cholesteryl stearate levels are decreased in coronary arteries with atherosclerotic plaques, isolated fatty streaks, and isolated atherosclerotic plaques compared with non-atherosclerotic coronary arteries.²

References

1. Wewer, V. and Dörmann, P. Determination of sterol lipids in plant tissues by gas chromatography and Q-TOF mass spectrometry. *Methods Mol. Biol.* **1153**, 115-133 (2014).
2. Ylä-Herttuala, S. Enrichment of long-chain ω 9 and ω 6 fatty acids in arterial cholesteryl esters in the early phase of atherogenesis. *Prog. Lipid Res.* **25**, 475-478 (1986).

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