

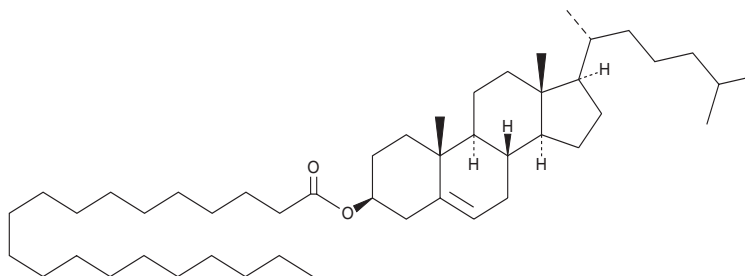
PRODUCT INFORMATION



Cholesteryl Arachidate

Item No. 26465

CAS Registry No.: 2573-03-7
Formal Name: (3 β)-cholest-5-en-3-ol
3-eicosanoate
Synonyms: CE(20:0), Cholesterol Arachidate,
Cholesteryl Eicosanoate,
20:0 Cholesteryl ester
MF: C₄₇H₈₄O₂
FW: 681.2
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 arachidate is supplied as a crystalline solid. A stock solution may be made by dissolving the cholesteryl arachidate in the solvent of choice. Cholesteryl arachidate 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 arachidate is a cholesterol ester that has been found in human meibum.^{1,2} It has been used as an internal standard for the quantification of fatty acids in *B. napus*.³

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. 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).
3. Madey, E., Nowack, L.M., and Thompson, J.E. Isolation and characterization of lipid in phloem sap of canola. *Planta* **4(4)**, 625-634 (2002).

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