

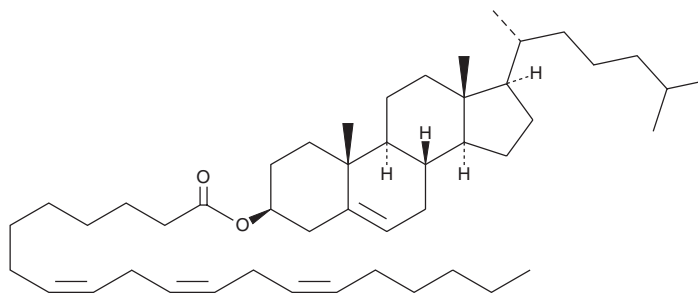
PRODUCT INFORMATION



Cholesteryl Homo- γ -Linolenate

Item No. 26469

CAS Registry No.: 7274-08-0
Formal Name: (3 β)-cholest-5-en-3-ol 3-(8Z,11Z,14Z)-8,11,14-eicosatrienoate
Synonyms: 20:3 (8Z,11Z,14Z) CE, 20:3 n-6 CE, 20:3 (8Z,11Z,14Z) Cholesterol Ester
MF: C₄₇H₇₈O₂
FW: 675.1
Purity: \geq 98%
Supplied as: An oil
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 homo- γ -linolenate is supplied as an oil. A stock solution may be made by dissolving the cholesteryl homo- γ -linolenate in the solvent of choice. Cholesteryl homo- γ -linolenate 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 homo- γ -linolenate is a cholesterol ester. It accumulates in the adrenal gland of rabbits fed a diet high in linolenic acid.¹ Levels of cholesteryl homo- γ -linolenate are decreased and positively correlate with cognitive decline in HIV-infected humans.²

References

1. Moore, J.H. and Williams, D.L. Dietary linoleic acid and the accumulation of homo-gamma-linolenic acid in the cholesterol esters of the adrenal gland of the rabbit. *Biochim Biophys. Acta.* **116(1)**, 181-183 (1966).
2. Bandaru, V.V., Mielke, M.M., Sacktor, N., *et al.* A lipid storage-like disorder contributes to cognitive decline in HIV-infected subjects. *Neurology* **81(17)**, 1492-1499 (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, 10/20/2022

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