

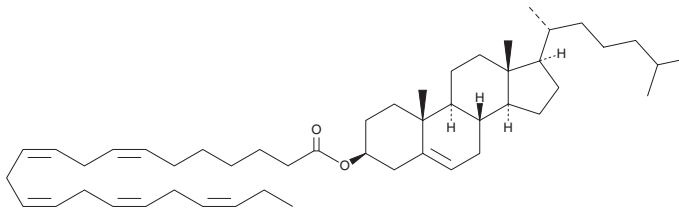
PRODUCT INFORMATION



Cholesteryl Docosapentaenoate

Item No. 22718

CAS Registry No.: 70110-49-5
Formal Name: cholest-5-en-3 β -ol,
3-[(7Z,10Z,13Z,16Z,19Z)-
7,10,13,16,19-docosapentaenoate]
MF: C₄₉H₇₈O₂
FW: 699.2
Purity: \geq 95%
Supplied as: A solution in chloroform
Storage: -20°C
Stability: \geq 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Cholesteryl docosapentaenoate is a cholesterol ester. It is elevated in liver, plasma, and skeletal muscle of acyl-CoA binding protein (ACBP) knockout mice compared with homozygous controls.¹

Reference

1. Gallego, S.F., Sporenger, R.R., Neess, D., *et al.* Quantitative lipidomics reveals age-dependent perturbations of whole-body lipid metabolism in ACBP deficient mice. *Biochim Biophys. Acta.* **1862(2)**, 145-155 (2017).

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, 11/28/2022

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