

PRODUCT INFORMATION



Desmosteryl Linoleate

Item No. 35423

Formal Name: (3S,8S,9S,10R,13R,14S,17R)-10,13-dimethyl-17-((R)-6-methylhept-5-en-2-yl)-2,3,4,7,8,9,10,11,12,13,14,15,16,17-tetradecahydro-1H-cyclopenta[a]phenanthren-3-yl (9Z,12E)-octadeca-9,12-dienoate

Synonyms: DE(18:2), 18:2(9Z,12Z) DE, 18:2(9Z,12Z) Desmosterol ester, Desmosteryl Linoleic Acid

MF: C₄₅H₇₄O₂

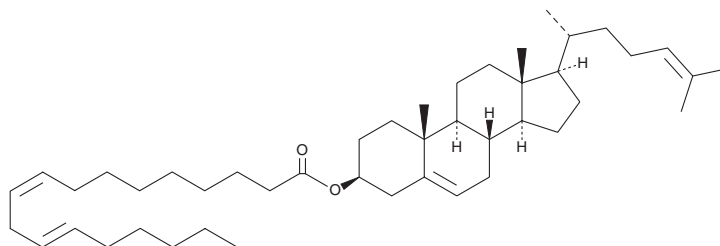
FW: 647.1

Purity: ≥98%

Supplied as: A solution in heptane

Storage: -20°C

Stability: ≥1 year



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

Description

Desmosteryl linoleate is an ester form of desmosterol (24-dehydro cholesterol; Item No. 14943), a precursor in the biosynthesis of cholesterol (Item No. 9003100), that is converted to cholesterol by Δ^{24} -dehydro cholesterol reductase (DHCR24).¹

Reference

1. Clayton, P.T. Disorders of cholesterol biosynthesis. *Arch. Dis. Child.* **78(2)**, 185-189 (1998).

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, 11/02/2021

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