

PRODUCT INFORMATION



Desmosteryl Palmitate

Item No. 34900

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 palmitate

Synonym: 16:0 Desmosteryl ester

MF: C₄₃H₇₄O₂

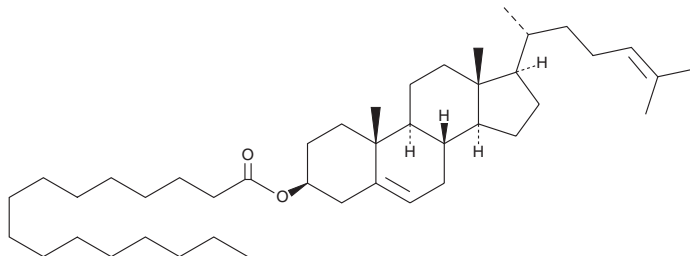
FW: 623.0

Purity: ≥95%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

Desmosteryl palmitate is supplied as a crystalline solid. A stock solution may be made by dissolving the desmosteryl palmitate in the solvent of choice, which should be purged with an inert gas. Desmosteryl palmitate is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

Description

Desmosteryl palmitate 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, 09/30/2021

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