

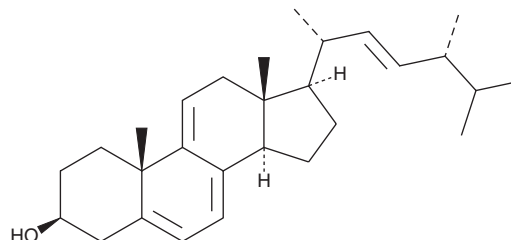
PRODUCT INFORMATION



Dehydroergosterol

Item No. 19709

CAS Registry No.: 516-85-8
Formal Name: ergosta-5,7,9(11),22E-tetraen-3 β -ol
Synonyms: DHE, 9,11-dehydro Ergosterol
MF: C₂₈H₄₂O
FW: 394.6
Purity: \geq 98%
UV/Vis.: λ_{max} : 311, 324 nm
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

Dehydroergosterol is supplied as a crystalline solid. A stock solution may be made by dissolving the dehydroergosterol in the solvent of choice, which should be purged with an inert gas. Dehydroergosterol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of dehydroergosterol in these solvents is approximately 20, 0.1, and 2 mg/ml, respectively.

Dehydroergosterol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, dehydroergosterol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Dehydroergosterol has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Dehydroergosterol is a naturally occurring, fluorescent analog of cholesterol (ex/em = 324/375 nm) that mimics the properties of cholesterol in cell membranes.¹ It is readily bound by cholesterol-binding proteins and has been used for real-time probing of the sterol environment and to elucidate intracellular sterol trafficking in living organisms.¹

Reference

1. McIntosh, A.L., Atshaves, B.P., Huang, H., *et al.* Fluorescence techniques using dehydroergosterol to study cholesterol trafficking. *Lipids* **43**(12), 1185-1208 (2008).

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, 12/20/2022

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