

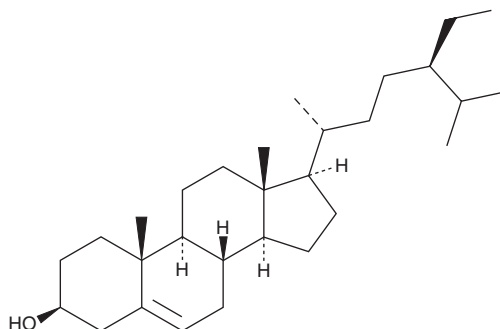
PRODUCT INFORMATION



24 α -ethyl Cholesterol

Item No. 37386

CAS Registry No.: 83-46-5
Formal Name: stigmaster-5-en-3 β -ol
Synonyms: Cupreol, α -Dihydrofucosterol, NSC 8096, NSC 18173, NSC 49083, Rhamnol, β -Sitosterol, SKF 14463, 22,23-dihydro Stigmasterol
MF: C₂₉H₅₀O
FW: 414.7
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/Unknown sp.



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

Laboratory Procedures

24 α -ethyl Cholesterol is supplied as a solid. A stock solution may be made by dissolving the 24 α -ethyl cholesterol in the solvent of choice, which should be purged with an inert gas. 24 α -ethyl Cholesterol is soluble in the organic solvents chloroform and ethanol. 24 α -ethyl Cholesterol is soluble in chloroform at a concentration of approximately 20 mg/ml and is slightly soluble in ethanol.

Description

24 α -ethyl Cholesterol is a phytosterol that has been found in *S. plebian* and has diverse biological activities.¹⁻⁵ It protects lard from thermal oxidation when used at concentrations of 0.02 and 0.04%.¹ 24 α -ethyl Cholesterol (100 μ M) induces glucose uptake and lipolysis in primary rat differentiated adipocytes and adipogenesis in primary rat preadipocytes.² It is cytotoxic to MCF-7 breast cancer cells (IC₅₀ = 264.83 μ M).³ 24 α -ethyl Cholesterol increases protein levels of the glucocorticoid receptor in human bronchial smooth muscle cells cultured with TGF- β 1 and prevents pulmonary fibrosis in a mouse model of allergic asthma induced by ovalbumin.⁴ Dietary administration of 24 α -ethyl cholesterol (0.5%) reduces cholesterol-induced increases in liver weight and hepatic levels of cholesterol (Item No. 9003100) in rabbits.⁵ 24 α -ethyl Cholesterol has also been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vitro* and *in vivo*.⁶ Formulations containing 24 α -ethyl cholesterol have been used as dietary supplements.

References

1. Weng, X.C. and Weng, W. *Food Chem.* **71**(4), 489-493 (2000).
2. Chai, J.W., Lim, S.L., Kanthimathi, M.S., et al. *Genes Nutr.* **6**(2), 181-188 (2011).
3. Hadrian, E., Sari, A.P., Mayanti, T., et al. *Indones. J. Chem.* **23**(1), 200-209 (2023).
4. Xu, J., Yang, L., and Lin, T. *Pulm. Pharmacol. Ther.* **78**, 102183 (2023).
5. Ikeda, I., Kawasaki, A., Samezima, K., et al. *J. Nutr. Sci. Vitaminol. (Tokyo)* **27**(3), 243-251 (1981).
6. Medjmedj, A., Ngalle-Loth, A., Clemençon, R., et al. *Nanomaterials (Basel)* **12**(14), 2446 (2022).

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, 07/25/2023

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