

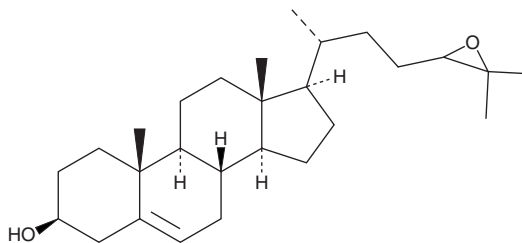
PRODUCT INFORMATION



24,25-epoxy Cholesterol

Item No. 25966

CAS Registry No.: 72542-49-5
Formal Name: (3 β)-24,25-epoxy-cholest-5-en-3-ol
Synonyms: (24R,S)-25-epoxy Cholesterol,
24,25-Oxidocholesterol
MF: C₂₇H₄₄O₂
FW: 400.6
Purity: $\geq 98\%$
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

24,25-epoxy Cholesterol is supplied as a crystalline solid. A stock solution may be made by dissolving the 24,25-epoxy cholesterol in the solvent of choice, which should be purged with an inert gas. 24,25-epoxy Cholesterol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 24,25-epoxy cholesterol in these solvents is approximately 3 mg/ml.

Description

24,25-epoxy Cholesterol is an oxysterol.¹ It inhibits the proliferation of HGC-27 gastric cancer cells when used at a concentration of 1 μ M, an effect that can be enhanced by knockdown of liver X receptor β (LXR β). Levels of 24,25-epoxy cholesterol are increased in the tumor tissues of patients with gastric cancer.

Reference

1. Guo, F., Hong, W., Yang, M., *et al.* Upregulation of 24(R/S),25-epoxycholesterol and 27-hydroxycholesterol suppresses the proliferation and migration of gastric cancer cells. *Biochem. Biophys. Res. Commun.* **504**(4), 892-898 (2018).

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, 04/07/2022

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