

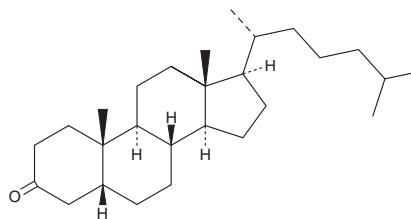
PRODUCT INFORMATION



5 β -Cholestan-3-one

Item No. 42284

CAS Registry No.: 601-53-6
Synonyms: Coprostanone, NSC 119031
MF: C₂₇H₄₆O
FW: 386.7
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 1 year



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

Laboratory Procedures

5 β -Cholestan-3-one is supplied as a solid. A stock solution may be made by dissolving the 5 β -cholestan-3-one in the solvent of choice, which should be purged with an inert gas. 5 β -Cholestan-3-one is sparingly soluble (1-10 mg/ml) in ethanol.

Description

5 β -Cholestan-3-one is an oxysterol and active metabolite of cholesterol (Item Nos. 9003100 | 39088) produced by gut microbiota.^{1,2} It induces apoptosis in primary dog gallbladder epithelial cells when used at a concentration of 1.4 mM.¹ Fecal levels of 5 β -cholestan-3-one are increased in patients with colon cancer or adenomatous polyps.²

References

1. Seo, D.W., Choi, H.S., Lee, S.P., *et al.* Oxysterols from human bile induce apoptosis of canine gallbladder epithelial cells in monolayer culture. *Am. J. Physiol. Gastrointest. Liver Physiol.* **287**(6), G1247-G1256 (2004).
2. Reddy, B.S. and Wynder, E.L. Metabolic epidemiology of colon cancer. Fecal bile acids and neutral sterols in colon cancer patients and patients with adenomatous polyps. *Cancer* **39**(6), 2533-2539 (1977).

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, 01/27/2025

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