

PRODUCT INFORMATION

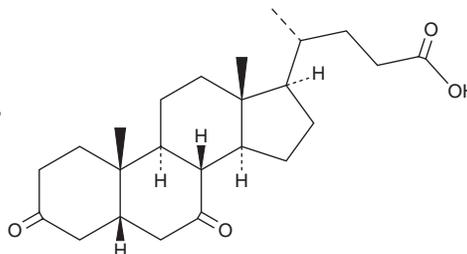


5 β -Cholanic Acid-3,7-dione

Item No. 43575

CAS Registry No.: 859-97-2
Formal Name: (5 β)-3,7-dioxo-cholan-24-oic acid
Synonyms: Dehydrochenodeoxycholic Acid,
3,7-Diketochenodeoxycholic Acid,
3,7-Dioxo-5 β -cholan-24-oic Acid

MF: C₂₄H₃₆O₄
FW: 388.5
Purity: \geq 98%
Supplied as: A 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

5 β -Cholanic acid-3,7-dione is supplied as a solid. A stock solution may be made by dissolving the 5 β -cholanic acid-3,7-dione in the solvent of choice, which should be purged with an inert gas. 5 β -Cholanic acid-3,7-dione is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

5 β -Cholanic acid-3,7-dione is a synthetic bile acid.¹ It decreases bile volume and bile levels of cholic acid and cholesterol in rabbits when administered at a dose of 5 mg/kg.

Reference

1. Hong, S.S., Kim, K.H., and Kim, W.J. Effect of bile acids on biliary excretion of cholesterol in rabbits. *Yonsei Med. J.* **14**, 109-115 (1973).

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, 07/15/2025

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