

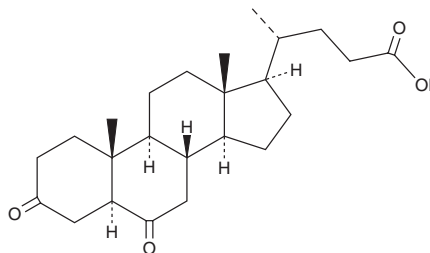
PRODUCT INFORMATION



5 α -Cholanic Acid-3,6-dione

Item No. 36184

CAS Registry No.: 6929-22-2
Formal Name: (5 α)-3,6-dioxo-cholan-24-oic acid
Synonym: NSC 18165
MF: C₂₄H₃₆O₄
FW: 388.5
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

5 α -Cholanic acid-3,6-dione is supplied as a solid. A stock solution may be made by dissolving the 5 α -cholanic acid-3,6-dione in the solvent of choice, which should be purged with an inert gas. 5 α -Cholanic acid-3,6-dione is soluble in chloroform.

Description

5 α -Cholanic acid-3,6-dione is a secondary bile acid.¹ Levels of 5 α -cholanic acid-3,6-dione are decreased in the large intestinal contents of germ-free mice.

Reference

1. Dempsey, J.L., Wang, D., Siginir, G., *et al.* Pharmacological activation of PXR and CAR downregulates distinct bile acid-metabolizing intestinal bacteria and alters bile acid homeostasis. *Toxicol. Sci.* **168**(1), 40-60 (2019).

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, 10/21/2022

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