

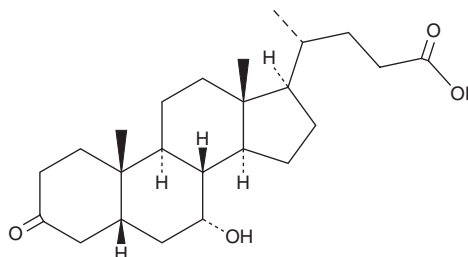
PRODUCT INFORMATION



7 α -hydroxy-3-oxo-5 β -Cholan-24-oic Acid

Item No. 38159

CAS Registry No.: 4185-00-6
Formal Name: (5 β)-7 α -hydroxy-3-oxo-cholan-24-oic acid
Synonyms: 3-keto-7 α -hydroxy-5 β -Cholanic Acid,
3-Oxochenodeoxycholic Acid
MF: C₂₄H₃₈O₄
FW: 390.6
Purity: \geq 95%
UV/Vis.: λ_{max} : 279 nm
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

7 α -hydroxy-3-oxo-5 β -Cholan-24-oic acid is supplied as a solid. A stock solution may be made by dissolving the 7 α -hydroxy-3-oxo-5 β -cholan-24-oic acid in the solvent of choice, which should be purged with an inert gas. 7 α -hydroxy-3-oxo-5 β -Cholan-24-oic acid is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide. The solubility of 7 α -hydroxy-3-oxo-5 β -cholan-24-oic acid in these solvents is approximately 10 mg/ml.

Description

7 α -hydroxy-3-oxo-5 β -Cholan-24-oic acid is a metabolite of chenodeoxycholic acid (Item Nos. 35346 | 10011286).¹ It is formed from chenodeoxycholic acid primarily by the cytochrome P450 (CYP) isoform CYP3A4 but also by CYP3A5 and other P450 enzymes at lower rates.

Reference

1. Deo, A.K. and Bandiera, S.M. Identification of human hepatic cytochrome p450 enzymes involved in the biotransformation of cholic and chenodeoxycholic acid. *Drug Metab. Dispos.* **36(10)**, 1983-1991 (2008).

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, 08/14/2023

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