

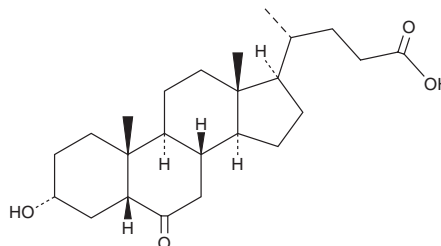
PRODUCT INFORMATION



6-keto Lithocholic Acid

Item No. 35373

CAS Registry No.: 2393-61-5
Formal Name: (3 α ,5 β)-3-hydroxy-6-oxo-cholan-24-oic acid
Synonyms: 5 β -Cholanic Acid-3 α -ol-6-one, 6-KLCA, 6-keto LCA, 6-oxo LCA, 6-oxo Lithocholic Acid, 6-keto Lithocholate, 6-oxo Lithocholate
MF: C₂₄H₃₈O₄
FW: 390.6
Purity: \geq 95%
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

6-keto Lithocholic acid is supplied as a solid. A stock solution may be made by dissolving the 6-keto lithocholic acid in the solvent of choice, which should be purged with an inert gas. 6-keto Lithocholic acid is sparingly soluble in dichloromethane and slightly soluble in methanol and DMSO.

Description

6-keto Lithocholic acid is a metabolite of the secondary bile acid lithocholic acid (Item Nos. 20253 | 31353).¹ It is formed from lithocholic acid by the cytochrome P450 (CYP) isoforms CYP3A1 and CYP3A2.

Reference

1. Deo, A.K. and Bandiera, S.M. Biotransformation of lithocholic acid by rat hepatic microsomes: Metabolite analysis by liquid chromatography/mass spectrometry. *Drug Metab. Dispos.* **36(2)**, 442-451 (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, 11/14/2022

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