

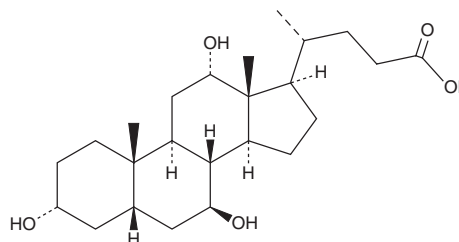
PRODUCT INFORMATION



Ursocholic Acid

Item No. 29883

CAS Registry No.: 2955-27-3
Formal Name: (3 α ,5 β ,7 β ,12 α)-3,7,12-trihydroxy-cholan-24-oic acid
Synonyms: 7 β -Hydroisocholic Acid, 7-Epichoic Acid, 3 α ,7 β ,12 α -Trihydroxy-5 β -cholanoic Acid, UCA
MF: C₂₄H₄₀O₅
FW: 408.6
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

Ursocholic acid is supplied as a solid. A stock solution may be made by dissolving the ursocholic acid in the solvent of choice, which should be purged with an inert gas. Ursocholic acid is slightly soluble in DMSO and ethanol.

Description

Ursocholic acid (UCA) is a bile acid and the 7 β -hydroxypimer of cholic acid (Item No. 20250).¹ Dietary administration of UCA (0.25%) decreases gallbladder cholesterol, phospholipid, and bile acid levels and the occurrence of gallstones in male mice fed a lithogenic diet.

Reference

1. Uchida, K., Akiyoshi, T., Igmi, H., *et al.* Differential effects of ursodeoxycholic acid and ursocholic acid on the formation of biliary cholesterol crystals in mice. *Lipids* **26**(7), 526-530 (1991).

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, 03/27/2025

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