

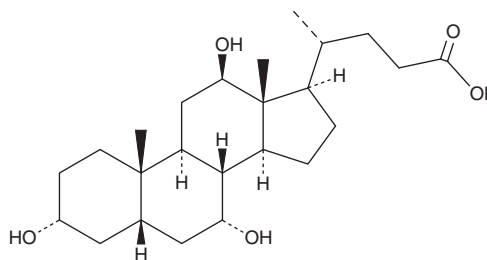
# PRODUCT INFORMATION



## 12 $\beta$ -Hydroxyisocholic Acid

Item No. 39638

**CAS Registry No.:** 71883-64-2  
**Formal Name:** (5 $\beta$ )-3 $\alpha$ ,7 $\alpha$ ,12 $\beta$ -trihydroxy-cholan-24-oic acid  
**Synonyms:** Lagocholic Acid, LCA  
**MF:** C<sub>24</sub>H<sub>40</sub>O<sub>5</sub>  
**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

12 $\beta$ -Hydroxyisocholic acid is supplied as a solid. A stock solution may be made by dissolving the 12 $\beta$ -hydroxyisocholic acid in the solvent of choice, which should be purged with an inert gas. 12 $\beta$ -Hydroxyisocholic acid is slightly soluble in DMF and methanol.

### Description

12 $\beta$ -Hydroxyisocholic acid is a bile acid and an isomer of cholic acid (Item No. 20250).<sup>1</sup> Hepatic levels of 12 $\beta$ -hydroxyisocholic acid are elevated in a mouse model of alcoholic liver damage.<sup>2</sup>

### References

1. Yamaga, N., Ikebuchi, J., Kohara, H., *et al.* Analysis of bile acids in urine specimens from healthy humans: Determination of several bile acids with  $\beta$ -hydroxyl and carbonyl groups. *J. Biochem.* **119(4)**, 725-730 (1996).
2. Yi, Z., Liu, X., Liang, L., *et al.* Antrodin A from *Antrodia camphorata* modulates the gut microbiome and liver metabolome in mice exposed to acute alcohol intake. *Food Funct.* **12(7)**, 2925-2937 (2021).

#### 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, 09/19/2023

#### 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