

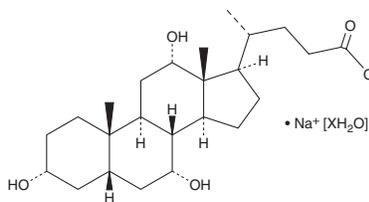
PRODUCT INFORMATION



Cholic Acid (sodium salt hydrate)

Item No. 44008

CAS Registry No.: 206986-87-0
Formal Name: (5 β)-3 α ,7,12 α -trihydroxy-cholan-24-oic acid, sodium salt hydrate
Synonyms: CA, Cholalic Acid, Cholalin, NSC 6135
MF: C₂₄H₃₉O₅ • Na [XH₂O]
FW: 488.6
Purity: \geq 98%
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

Cholic acid (sodium salt hydrate) is supplied as a solid. A stock solution may be made by dissolving the cholic acid (sodium salt hydrate) in the solvent of choice, which should be purged with an inert gas. Cholic acid (sodium salt hydrate) is soluble (\geq 10 mg/ml) in ethanol and DMSO.

Description

Cholic acid is a primary bile acid.¹ It is formed from cholesterol *via* a multistep process catalyzed by the cytochrome P450 (CYP) isoforms CYP7A1, CYP8B1, and CYP27A1. Cholic acid is conjugated to glycine or taurine by bile acid-CoA:amino acid N-acyltransferase (BAAT) to produce glycocholic acid (GCA; Item No. 20276) and taurocholic acid (TCA; Item No. 16215), respectively, in the liver, and is transformed into the secondary bile acid deoxycholic acid (DCA; Item Nos. 20756 | 18231) by intestinal microbiota.¹⁻³ It induces *C. difficile* colony formation in an agar dilution assay when used at a concentration of 0.1% w/v.⁴ Dietary administration of cholic acid (0.4% w/w) increases serum cholesterol levels, biliary phospholipid secretion, and fecal DCA levels in rats.⁵

References

- Šarenac, T.M. and Mikov, M. Bile acid synthesis: From nature to the chemical modification and synthesis and their applications as drugs and nutrients. *Front. Pharmacol.* **9**, 939 (2018).
- Hunt, M.C., Siponen, M.I., and Alexson, S.E.H. The emerging role of acyl-CoA thioesterases and acyltransferases in regulating peroxisomal lipid metabolism. *Biochim. Biophys. Acta* **1822(9)**, 1397-1410 (2012).
- Staley, C., Weingarden, A.R., Khoruts, A., *et al.* Interaction of gut microbiota with bile acid metabolism and its influence on disease states. *Appl. Microbiol. Biotechnol.* **101(1)**, 47-64 (2017).
- Sorg, J.A. and Sonenshein, A.L. Bile salts and glycine as cogerminants for *Clostridium difficile* spores. *J. Bacteriol.* **190(7)**, 2505-2512 (2008).
- Uchida, K., Nomura, Y., and Takeuchi, N. Effects of cholic acid, chenodeoxycholic acid, and their related bile acids on cholesterol, phospholipid, and bile acid levels in serum, liver, bile, and feces of rats. *J. Biochem.* **87(1)**, 187-194 (1980).

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, 08/04/2025

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