

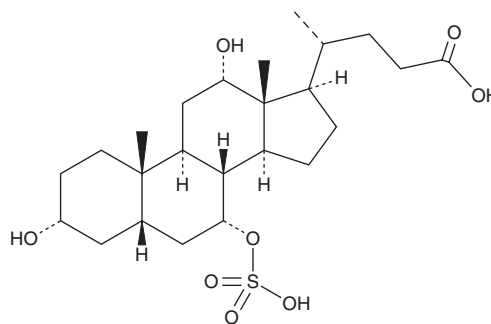
PRODUCT INFORMATION



Cholic Acid 7-sulfate

Item No. 9002532

CAS Registry No.: 60320-05-0
Formal Name: (3 α ,5 β ,7 α ,12 α)-3,12-dihydroxy-7-(sulfooxy)-cholan-24-oic acid
Synonyms: 7-Sulfocholate, 7-Sulfocholic Acid
MF: C₂₄H₄₀O₈S
FW: 488.6
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Cholic acid 7-sulfate is supplied as a crystalline solid. A stock solution may be made by dissolving the cholic acid 7-sulfate in the solvent of choice. Cholic acid 7-sulfate is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of cholic acid 7-sulfate in these solvents is approximately 0.3 and 1 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of cholic acid 7-sulfate can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of cholic acid 7-sulfate in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Cholic acid 7-sulfate is a metabolite of the primary bile acid cholic acid (Item No. 20250).^{1,2} It is produced by conjugation of a sulfate group with the hydroxy group at position 7 of cholic acid in the liver and gut. Fecal levels of cholic acid 7-sulfate are increased in male, but not female, mice fed a diet supplemented with cholic acid or chenodeoxycholic acid (Item No. 10011286).²

References

1. Alnouti, Y. Bile acid sulfation: A pathway of bile acid elimination and detoxification. *Toxicol. Sci.* **108**(2), 225-246 (2009).
2. Zhang, Y. and Klaassen, C.D. Effects of feeding bile acids and a bile acid sequestrant on hepatic bile acid composition in mice. *J. Lipid Res.* **51**(11), 3230-3242 (2010).

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, 04/11/2019

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