

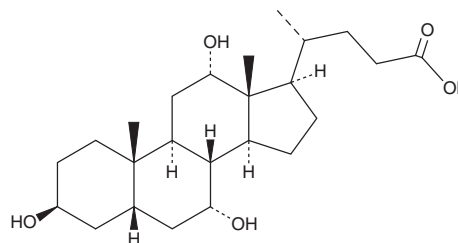
PRODUCT INFORMATION



3 β -Cholic Acid

Item No. 36447

CAS Registry No.: 3338-16-7
Formal Name: (5 β)-3 β ,7 α ,12 α -trihydroxy-cholan-24-oic acid
Synonym: 3 β -Epicholic Acid
MF: C₂₄H₄₀O₅
FW: 408.6
Purity: \geq 98%
Supplied as: A 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

3 β -Cholic acid is supplied as a solid. A stock solution may be made by dissolving the 3 β -cholic acid in the solvent of choice, which should be purged with an inert gas. 3 β -Cholic acid is slightly soluble in DMSO and ethanol (sonicated).

Description

3 β -Cholic acid is a bile acid and derivative of cholic acid (Item No. 20250) that has been found in human feces.¹

Reference

1. Eneroth, P., Gordon, B., and Sjövall, J. Characterization of trisubstituted cholanoic acids in human feces. *J. Lipid Res.* **7**(4), 524-530 (1966).

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, 03/03/2022

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