

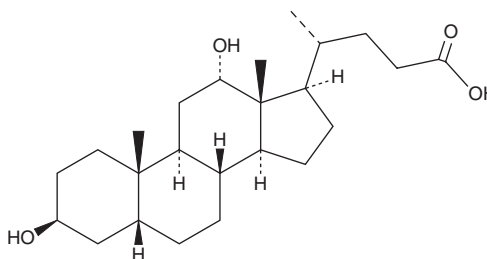
PRODUCT INFORMATION



3-Epideoxycholic Acid

Item No. 34984

CAS Registry No.: 570-63-8
Formal Name: (3 β ,5 β ,12 α)-3,12-dihydroxy-
cholan-24-oic acid
Synonym: EDCA
MF: C₂₄H₄₀O₄
FW: 392.6
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

3-Epideoxycholic acid (EDCA) is supplied as a solid. A stock solution may be made by dissolving the EDCA in the solvent of choice, which should be purged with an inert gas. EDCA is soluble in methanol.

Description

EDCA is a secondary bile acid and an epimer of deoxycholic acid (Item Nos. 20756 | 18231).¹ EDCA has been used as a steroidal host compound for the enantioselective purification of (2R,3S)-3-methyl-2-pentanol from its racemates *via* enclathration.² Fecal levels of EDCA are decreased in patients with pediatric non-alcoholic fatty liver disease (NAFLD).¹

References

1. Yu, J., Zhang, H., Chen, L., *et al.* Disease-associated gut microbiota reduces the profile of secondary bile acids in pediatric nonalcoholic fatty liver disease. *Front. Cell. Infect. Microbiol.* **11**, 698852 (2021).
2. Kato, K., Aburaya, K., Miyake, Y., *et al.* Excellent enantio-selective enclathration of (2R,3S)-3-methyl-2-pentanol in channel-like cavity of 3-epideoxycholic acid, interpreted by the four-location model for chiral recognition. *Chem. Commun. (Camb)* **23**, 2872-2873 (2003).

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, 12/13/2022

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