

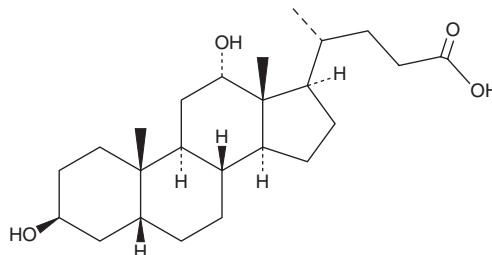
PRODUCT INFORMATION



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
Synonyms: 3 β -DCA, 3 β -Deoxycholic acid, EDCA, 3 β -Hydroxydeoxycholic acid
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

3-Epideoxycholic acid is a secondary bile acid formed *via* epimerization of deoxycholic acid (DCA; Item Nos. 20756 | 18231) by gut microbiota.¹ It has reduced detergent activity and is less active than DCA in inhibiting growth in a panel of seven gut bacteria species. 3-Epideoxycholic acid (10 and 20 μ M) increases the growth of Caco-2 colon cancer cells but to a lesser extent than DCA.² It increases the levels of Foxp3⁺CD4⁺ regulatory T cells (Tregs) in a co-culture of naïve T cells and dendritic cells.³

References

1. Devlin, A.S. and Fischbach, M.A. A biosynthetic pathway for a prominent class of microbiota-derived bile acids. *Nat. Chem. Biol.* **11**(9), 685-690 (2018).
2. Milovic, V., Teller, I.C., Faust, D., *et al.* Effects of deoxycholate on human colon cancer cells: Apoptosis or proliferation. *Eur. J. Clin. Invest.* **32**(1), 29-34 (2002).
3. Campbell, C., McKenney, P.T., Konstantinovskiy, D., *et al.* Bacterial metabolism of bile acids promotes generation of peripheral regulatory T cells. *Nature* **58**(7809), 475-479 (2020).

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, 10/21/2024

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