

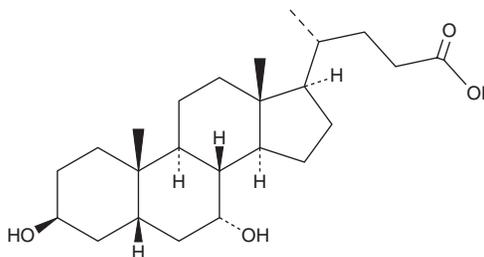
PRODUCT INFORMATION



Isochenodeoxycholic Acid

Item No. 44228

CAS Registry No.: 566-24-5
Formal Name: (5 β)-3 β ,7 α -dihydroxy-cholan-24-oic acid
Synonym: iso-CDCA
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

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

Description

Isochenodeoxycholic acid is a secondary bile acid.¹ It is produced from chenodeoxycholic acid (Item Nos. 10011286 | 35346) by gut bacteria.^{1,2}

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. Batta, A.K., Salen, G., and Shefer, S. Transformation of bile acids into iso-bile acids by *Clostridium perfringens*: Possible transport of 3 β -hydrogen via the coenzyme. *Hepatology* **5**(6), 1126-1131 (1985).

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, 10/08/2025

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