

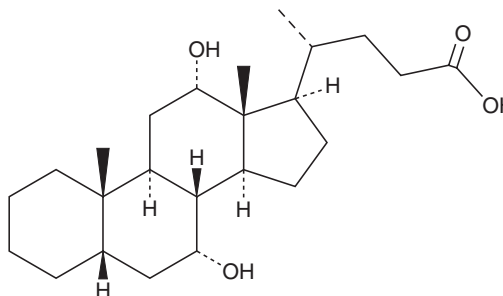
PRODUCT INFORMATION



Isodeoxycholic Acid

Item No. 29890

CAS Registry No.: 566-17-6
Formal Name: (5 β ,7 α ,12 α)-7,12-dihydroxy-
cholan-24-oic acid
Synonyms: 3-deoxyCA, 3-deoxy Cholic Acid,
3-DCA, iDCA, isoDCA,
7 α ,12 α -Dihydroxycholan-ic Acid,
7 α ,12 α -Dihydroxycholic Acid
MF: C₂₄H₄₀O₄
FW: 392.6
Purity: \geq 95%
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

Isodeoxycholic acid is supplied as a solid. A stock solution may be made by dissolving the isodeoxycholic acid in the solvent of choice, which should be purged with an inert gas. Isodeoxycholic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of isodeoxycholic acid in ethanol and DMSO is approximately 20 mg/ml and approximately 30 mg/ml in DMF.

Isodeoxycholic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, isodeoxycholic acid should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Isodeoxycholic acid has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Isodeoxycholic acid is a bile acid and an agonist of the farnesoid X receptor (FXR).¹ It induces FXR-mediated gene expression in a reporter assay using COS-7 cells when used at a concentration of 30 μ M. Isodeoxycholic acid also activates large-conductance calcium-activated potassium channels (K_{Ca}1.1/BK) in isolated rabbit superior mesenteric arteries.²

References

1. T., S., Tamehiro, N., Sato, Y., *et al.* The novel compounds that activate farnesoid X receptor: The diversity of their effects on gene expression. *J. Pharmacol. Sci.* **107(3)**, 285-294 (2008).
2. Dopico, A.M., Walsh, J.V., and Singer, J.J. Natural bile acids and synthetic analogues modulate large conductance Ca²⁺-activated K⁺ (BK_{Ca}) channel activity in smooth muscle cells. *J. Gen. Physiol.* **119(3)**, 251-273 (2002).

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