

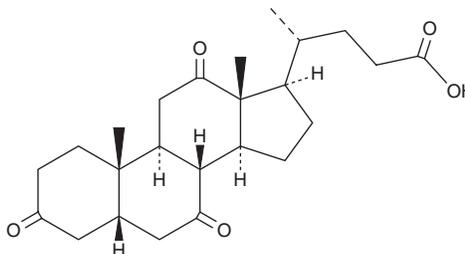
PRODUCT INFORMATION



Dehydrocholic Acid

Item No. 29176

CAS Registry No.: 81-23-2
Formal Name: (5 β)-3,7,12-trioxo-cholan-24-oic acid
Synonyms: NSC 8796, 3,7,12-Triketocholanoic Acid, Triketocholanoic Acid
MF: C₂₄H₃₄O₅
FW: 402.5
Purity: \geq 95%
Supplied as: A crystalline 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

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

Dehydrocholic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, dehydrocholic acid should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Dehydrocholic acid has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Dehydrocholic acid is a synthetic bile acid and a derivative of cholic acid (Item No. 20250).¹ It increases bile flow by 2.7-fold and decreases biliary levels of phospholipids, cholesterol, and bilirubin in conscious dogs when administered at a dose of 50 mg/kg.² Dehydrocholic acid (1 μ mol/min/0.1 kg) increases bile flow by 253% and decreases the secretion rate of phospholipids and cholesterol by 64 and 94%, respectively, in rats when infused intravenously at a dose of 2 μ mol/min/0.1 kg.³

References

1. Soloway, R.D., Hofmann, A.F., Thomas, P.J., *et al.* Triketocholanoic (dehydrocholic) acid. Hepatic metabolism and effect on bile flow and biliary lipid secretion in man. *J. Clin. Invest.* **52(3)**, 715-724 (1973).
2. Yanaura, S. and Ishikawa, S. Choleric properties of ursodeoxycholic acid and chenodeoxycholic acid in dogs. *Jpn. J. Pharmacol.* **28(3)**, 383-389 (1978).
3. Yousef, I.M., Mignault, D., Weber, A.M., *et al.* Influence of dehydrocholic acid on the secretion of bile acids and biliary lipids in rats. *Digestion* **45(1)**, 40-51 (1990).

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, 11/02/2022

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