

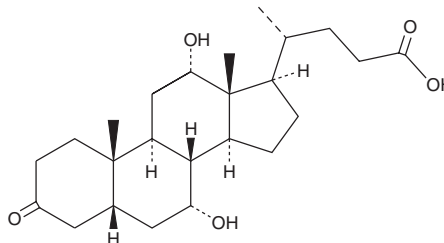
PRODUCT INFORMATION



3-Oxochohic Acid

Item No. 28403

CAS Registry No.: 2304-89-4
Formal Name: (5 β ,7 α ,12 α)-7,12-dihydroxy-3-oxo-
cholan-24-oic acid
Synonym: 3-Dehydrocholic Acid
MF: C₂₄H₃₈O₅
FW: 406.6
Purity: \geq 95%
UV/Vis.: λ_{max} : 284 nm
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

3-Oxochohic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 3-oxochohic acid in the solvent of choice, which should be purged with an inert gas. 3-Oxochohic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 3-oxochohic acid in these solvents is approximately 30 mg/ml.

3-Oxochohic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 3-oxochohic acid should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 3-Oxochohic 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

3-Oxochohic acid is a secondary bile acid and a metabolite of cholic acid (Item No. 20250) in *C. perfringens*, a bacteria found in the intestines of mammals and in the environment.^{1,2} 3-Oxochohic acid serum levels increase following ileal transposition surgery in rats.³

References

1. Wahlström, A., Sayin, S.I., Marschall, H.-I., *et al.* Intestinal crosstalk between bile acids and microbiota and its impact on host metabolism. *Cell Metab.* **24**(1), 41-50 (2016).
2. Macdonald, I.A., Forrest, T.P., Costain, G.A., *et al.* Identification of 7 α -, 12 α -dihydroxy-3-oxo cholanoic acid as the major degradation product from cholic by *C. perfringens*. *J. Steroid Biochem.* **9**(4), 353-358 (1978).
3. Yan, K., Chen, W., Zhu, H., *et al.* The changes of serum metabolites in diabetic GK rats after ileal transposition surgery. *Obes. Surg.* **29**(3), 882-890 (2019).

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, 11/11/2022

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