

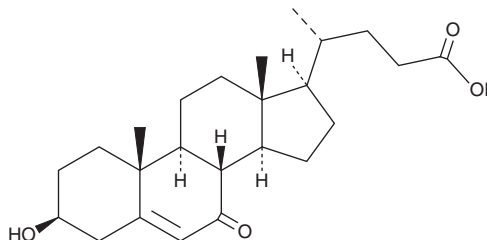
PRODUCT INFORMATION



3 β -OH-7-Oxocholeonic Acid

Item No. 29541

CAS Registry No.: 25218-38-6
Formal Name: (3 β)-3-hydroxy-7-oxo-chole-5-en-24-oic acid
Synonyms: 3 β -hydroxy-7-oxo-5-Cholenoic Acid, 3 β -hydroxy-7-Oxochole-5-enoic Acid
MF: C₂₄H₃₆O₄
FW: 388.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 238 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 β -OH-7-Oxocholeonic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 3 β -OH-7-oxocholeonic acid in the solvent of choice, which should be purged with an inert gas. 3 β -OH-7-Oxocholeonic acid is soluble in organic solvents such as ethanol, methanol, and acetonitrile. The solubility of 3 β -OH-7-oxocholeonic acid in these solvents is approximately 1 mg/ml.

Description

3 β -OH-7-Oxocholeonic acid is a bile acid.¹ It is also a metabolite of 7 β -hydroxy cholesterol (Item No. 20099) in rats. Conjugated forms of 3 β -OH-7-oxocholeonic acid have been found in the urine of patients with Neimann-Pick disease type C.^{2,3}

References

1. Norii, T., Yamaga, N., and Yamasaki, K. Metabolism of 7 β -hydroxycholesterol-4-¹⁴C in rat. *Steroids* **15**(3), 303-326 (1970).
2. Alvelius, G., Hjalmarson, O., Griffiths, W.J., *et al.* Identification of unusual 7-oxygenated bile acid sulfates in a patient with Niemann-Pick disease, type C. *J. Lipid Res.* **42**(10), 1571-1577 (2001).
3. Maekawa, M., Omura, K., Sekiguchi, S., *et al.* Identification of two sulfated cholesterol metabolites found in the urine of a patient with Niemann-Pick disease type C as novel candidate diagnostic markers. *Mass Spectrom. (Tokyo)* **5**(2), S0053 (2016).

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, 03/27/2025

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