

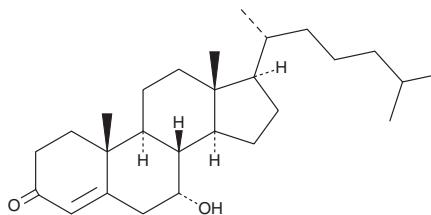
PRODUCT INFORMATION



7 α -hydroxy-4-Cholesten-3-one

Item No. 25442

CAS Registry No.: 3862-25-7
Formal Name: 7 α -hydroxy-cholest-4-en-3-one
MF: C₂₇H₄₄O₂
FW: 400.6
Purity: \geq 98%
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

7 α -hydroxy-4-Cholesten-3-one is supplied as a solid. A stock solution may be made by dissolving the 7 α -hydroxy-4-cholesten-3-one in the solvent of choice, which should be purged with an inert gas. 7 α -hydroxy-4-Cholesten-3-one is slightly soluble in methanol and chloroform.

Description

7 α -hydroxy-4-Cholesten-3-one is a metabolite of 7 α -hydroxy cholesterol and an intermediate in the biosynthesis of bile acids.¹ Plasma levels of 7 α -hydroxy-4-cholesten-3-one positively correlate with the rate of bile acid synthesis in humans and are increased in patients following ileal resection.²

References

1. Russell, D.W. The enzymes, regulation, and genetics of bile acid synthesis. *Annu. Rev. Biochem.* **72**, 137-174 (2003).
2. Axelson, M., Aly, A., and Sjövall, J. Levels of 7 α -hydroxy-4-cholesten-3-one in plasma reflect rates of bile acid synthesis in man. *FEBS Lett.* **239(2)**, 324-328 (1988).

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, 07/23/2018

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