

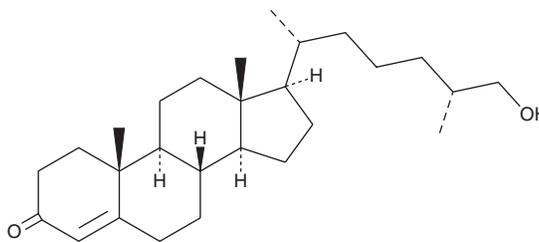
PRODUCT INFORMATION



27-hydroxy Cholestenone

Item No. 25550

CAS Registry No.: 56792-59-7
Formal Name: (25R)-26-hydroxy-cholest-4-en-3-one
MF: C₂₇H₄₄O₂
FW: 400.6
Purity: ≥95%
UV/Vis.: λ_{max}: 242 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

27-hydroxy Cholestenone is supplied as a crystalline solid. A stock solution may be made by dissolving the 27-hydroxy cholestenone in the solvent of choice, which should be purged with an inert gas. 27-hydroxy Cholestenone is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 27-hydroxy cholestenone in these solvents is approximately 2 and 0.1 mg/ml, respectively.

Description

27-hydroxy Cholestenone is an oxysterol.¹ Levels of 27-hydroxy cholestenone are increased in *Cyp46a*^{-/-} mouse brain. 27-hydroxy Cholestenone has been used as an intermediate in the synthesis of dafachronic acid ligands for the DAF-12 nuclear hormone receptor.²

References

1. Meljon, A., Wang, Y., and Griffiths, W.J. Oxysterols in the brain of the cholesterol 24-hydroxylase knockout mouse. *Biochem. Biophys. Res. Commun.* **446**(3), 768-774 (2014).
2. Sharma, K.K., Wang, Z., Motola, D.L., et al. Synthesis and activity of dafachronic acid ligands for the *C. elegans* DAF-12 nuclear hormone receptor. *Mol. Endocrinol.* **23**(5), 640-648 (2009).

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, 12/09/2022

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